

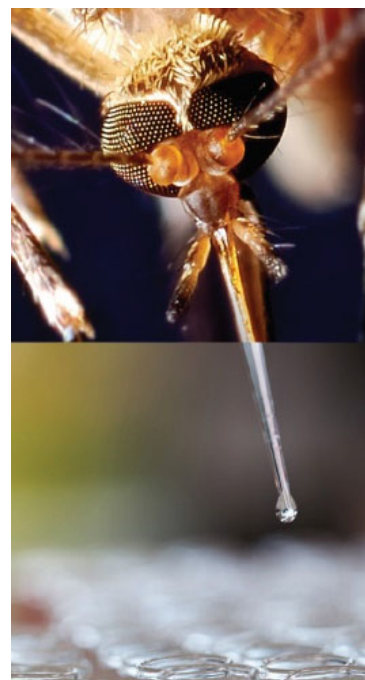
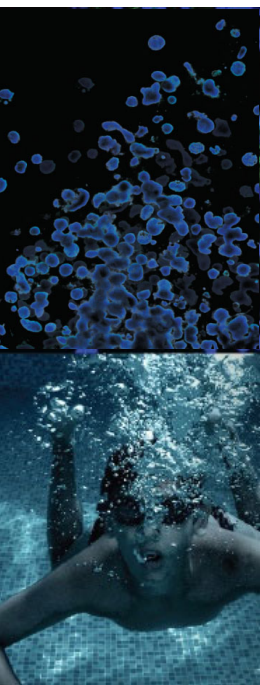


Extracellular Vesicles, HIV-1, and Neurospheres: Touching a Nerve

Heather Branscome, Ph.D.
Supervisor, Laboratory Operations, ATCC

Fatah Kashanchi, Ph.D.
Professor, George Mason University

Credible Leads to Incredible™

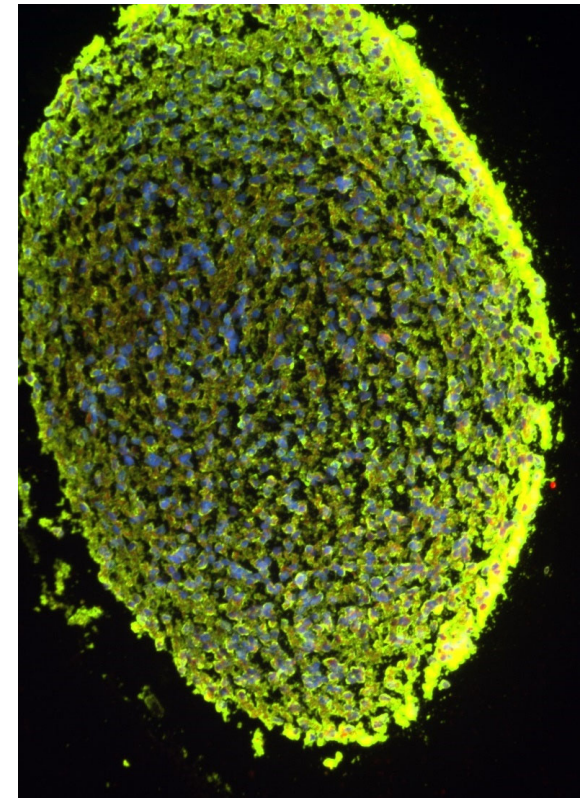


About ATCC

- Founded in 1925, ATCC is a non-profit organization with HQ in Manassas, VA, and an R&D and Services center in Gaithersburg, MD
- World's largest, most diverse biological materials and information resource for cell culture – the “*gold standard*”
- Innovative R&D company featuring gene editing, differentiated stem cells, advanced models
- cGMP biorepository
- Partner with government, industry, and academia
- Leading global supplier of authenticated cell lines, viral and microbial standards
- Sales and distribution in 150 countries, 19 international distributors
- Talented team of 450+ employees, over one-third with advanced degrees

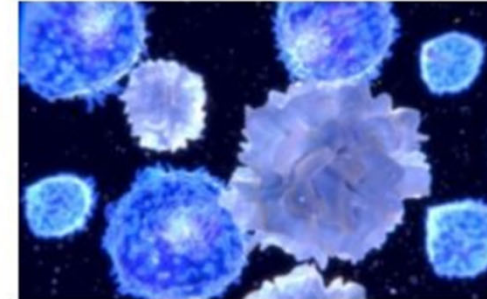
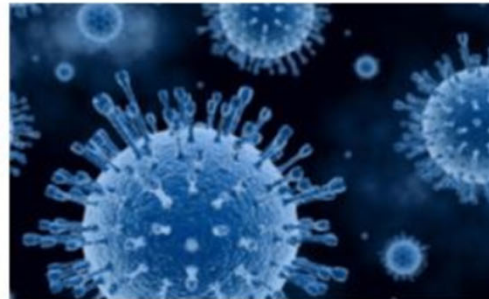
Overview

- Update on the most recent literature surrounding EVs from virally infected cells (“Damaging EVs”)
- Generation and infection of iPSC-derived neurospheres
- Effect of stem cell EVs (“Reparative EVs”) on HIV-1 infected neurospheres
- Summary



Neural progenitor cell-derived neurosphere

Meeting of the American Society for Intercellular Communication



It gives us great pleasure to announce the 1st meeting of the **American Society for Intercellular Communication (ASIC)**

The meeting will be held on **October 21st – October 23rd 2021**.

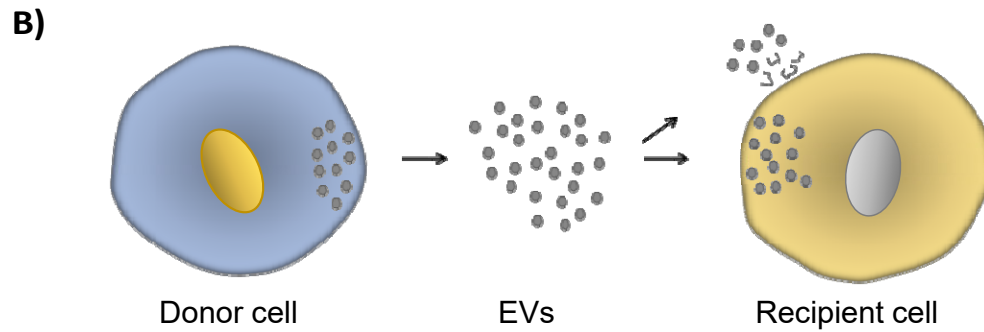
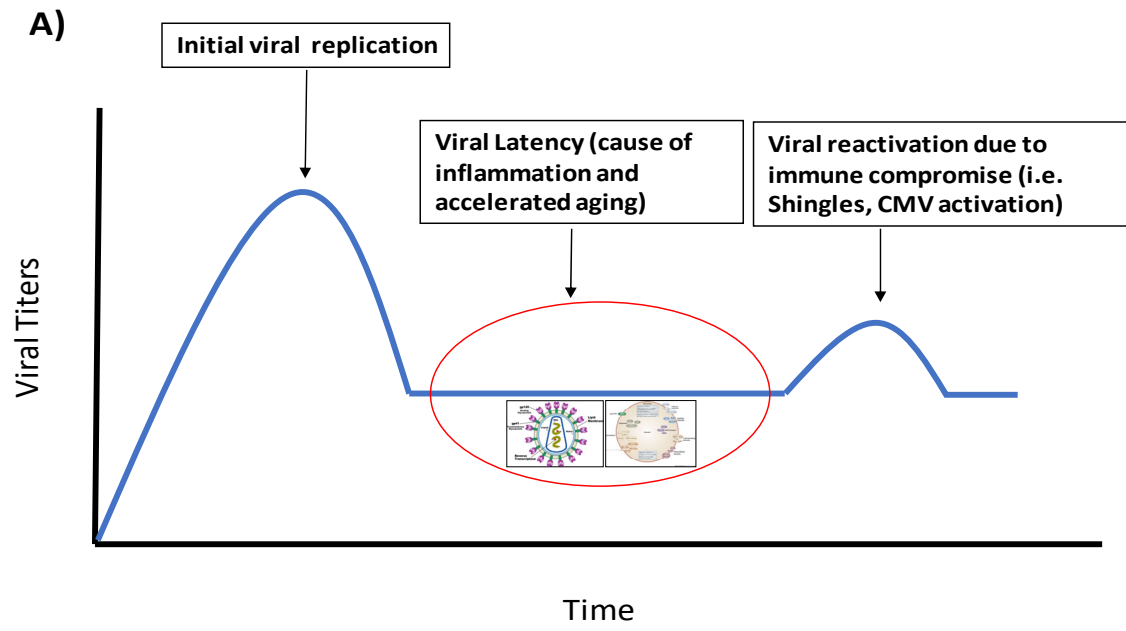
The meeting is organized by a group of investigators seeking to create a platform for informal exchange of ideas on emerging questions and cutting-edge developments in the field of Extracellular Vesicles (EVs), Extracellular Particles (EPs), and particulate carriers of extracellular RNA (exRNA) as biological mediators, regulators and diagnostic analytes. Our program of relatively short talks (15 and 30 minutes) encourage discussions that can take place after each scientific talk, where the audience and attendees can debate the concepts presented to foster a dynamic scientific exchange.

We invite you to join us for a 3-day meeting (Thursday afternoon, all day Friday and Saturday until noon; October 21st – October 23rd) at the Bolger Center in Potomac, Maryland. The program will include the functions of EVs, EPs, exRNA in cancer, CNS diseases, infections (both bacterial and viral), the mechanistic basis for their biogenesis and activities, and technological and diagnostic advances in the field. *In the interest of everyone's health and safety, vaccines are required to attend this meeting in person. We will also be following CDC guidelines for this meeting.*

Late Breaking Abstracts are due by September 30, 2021

Late Breaking Registration: September 30, 2021

EVs from latent cells



A)

Table 1. Baseline Characteristics Among 190 Study Subjects

Characteristic	Value
Age, years	51 (44–57)
Male sex, % of patients	92
CD4 ⁺ T-cell count, cells/mm ³	523 (249–728)
T cells expressing CD4, %	23 (16–32)
CD8 ⁺ T-cell count, cells/mm ³	844 (606–1185)
T cells expressing CD8, %	48 (37–57)
Nadir CD4 ⁺ T-cell count, cells/mm ³	113 (29–227)
Duration of HAART suppression, months	31 (14–66)
Agent included in HAART, no. of patients	
Protease inhibitor	124
NNRTI	73
Raltegravir	14
Maraviroc	1
Enfuvirtide	3

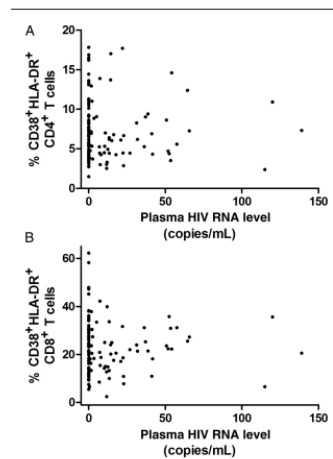


Figure 1. No associations between ultrasensitive plasma human immunodeficiency virus (HIV) RNA levels and T-cell activation. Ultrasensitive plasma HIV RNA levels were measured using the COBAS AmpliPrep/COBAS TaqMan HIV-1 test, version 2.0. All *P* values are >.40.

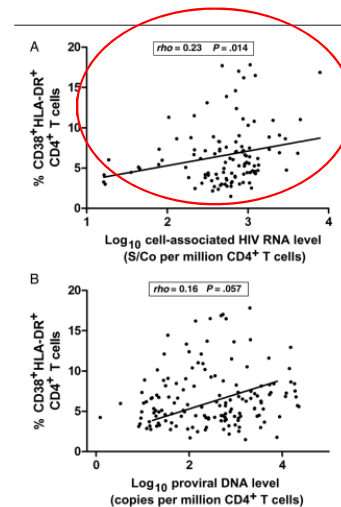


Figure 2. Cell-based measures of viral persistence are modestly associated with immune activation.

B)

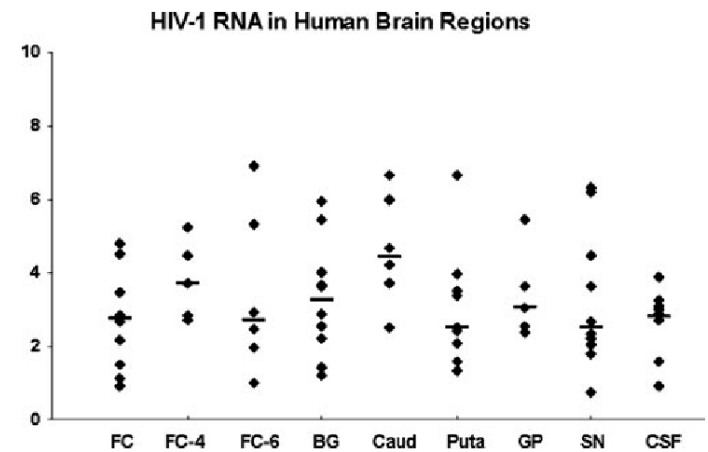
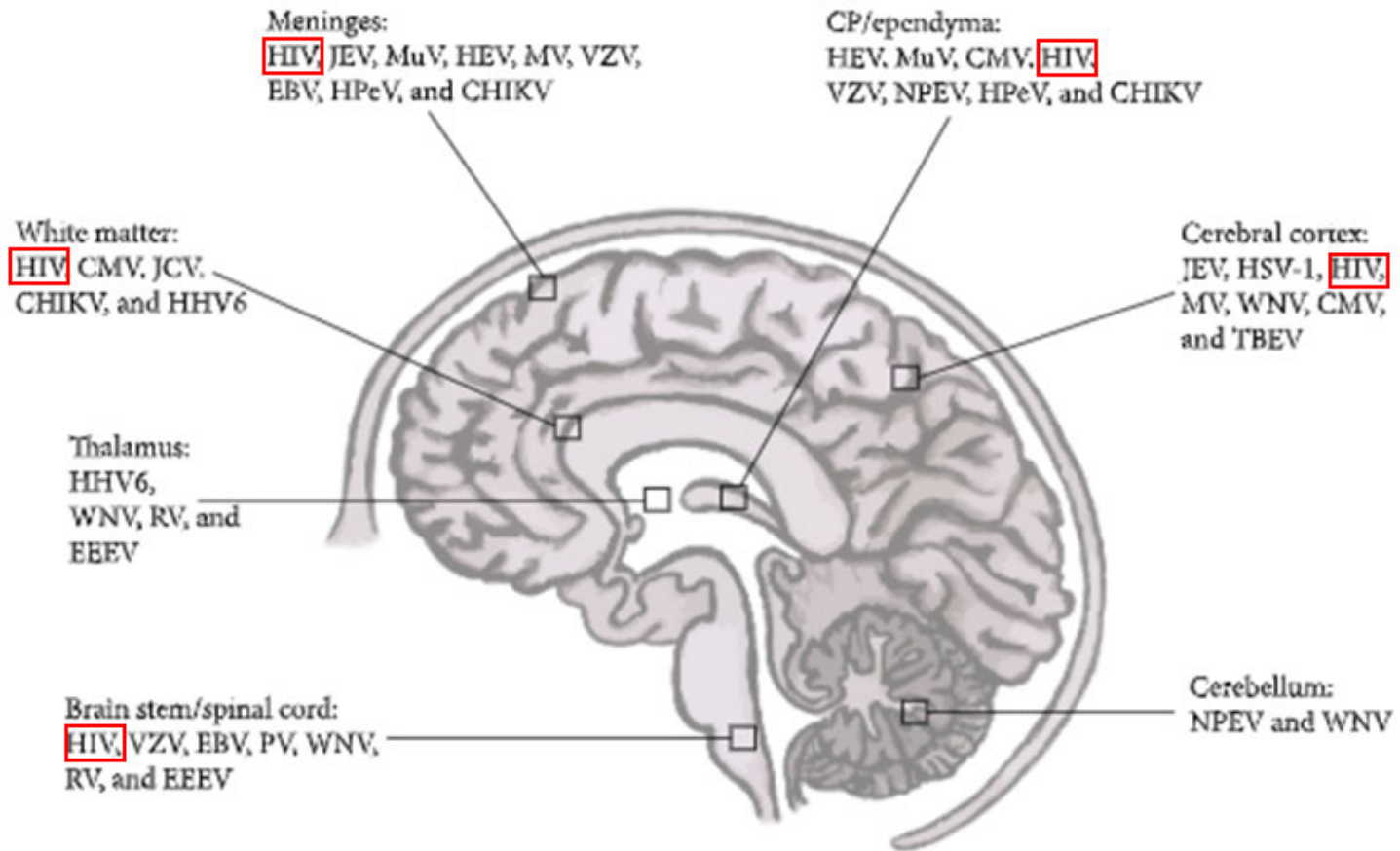


Figure 3 HIV-1 RNA load (\log_{10} copies/g tissue) in each region of HIV-1+ individuals is presented in a scatter plot. The horizontal bars represent the median value in each region. Viral load values were higher in caudate nucleus compared to that in the other regions (FC-4, FC-6, BG, putamen, globus pallidus, SN), and the lowest values were found in FC and CSF.

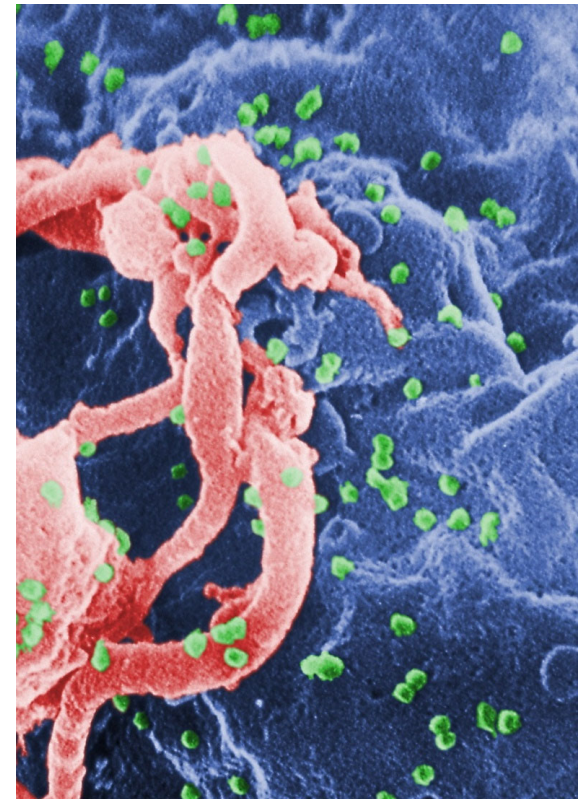
A) Hatano/Deeks et. al., JID. 2013
 B) Kumar et. al., J Neurovirol. 2007

CNS Viral Infections



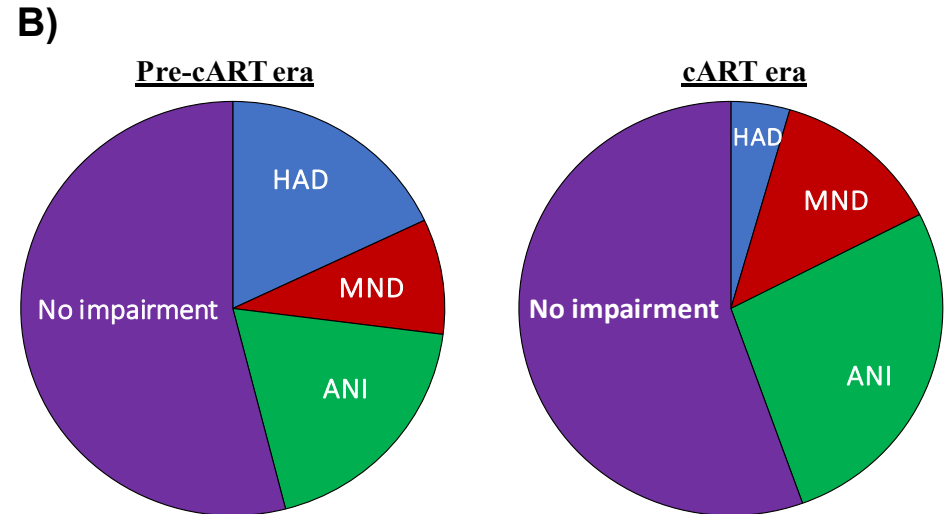
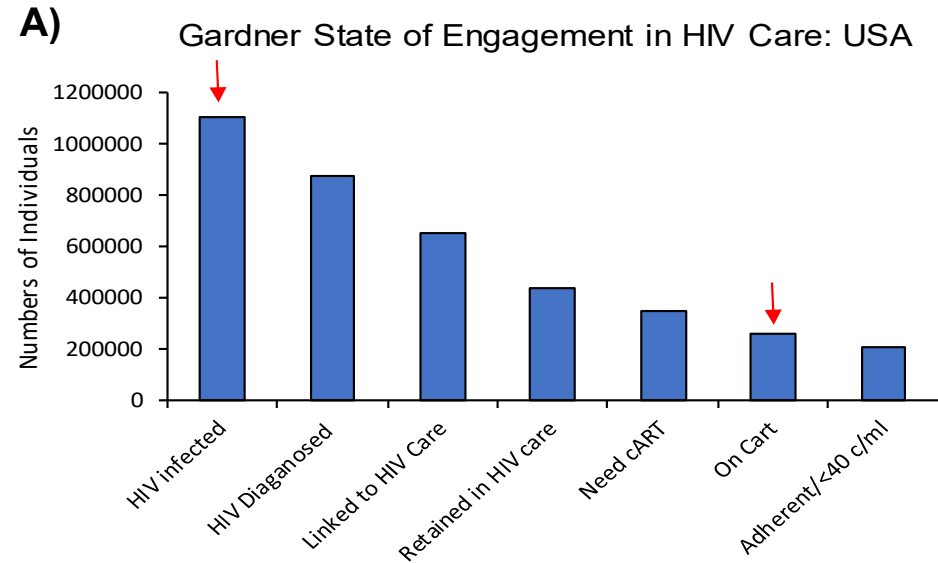
HIV-1 and HAND

- HIV-associated neurocognitive disorders (HAND) are comprised of a range of neurological dysfunctions which are commonly associated with HIV-1 infection.
- Despite cART, HAND still persists in HIV-1 patients (>50%).
- Exosomes secreted from HIV-1 infected cells under cART can damage CNS cells (damaging exosomes).
 - Damaging exosomes contain viral non-coding RNA (TAR) and other viral proteins (Tat, Nef, Env) that cause neuroinflammation.
- Damaging exosomes have also been isolated from 6 other CNS tropic viral infections including HTLV-1, Zika, RVFV, Ebola, HSV-1, and HHV-6A.

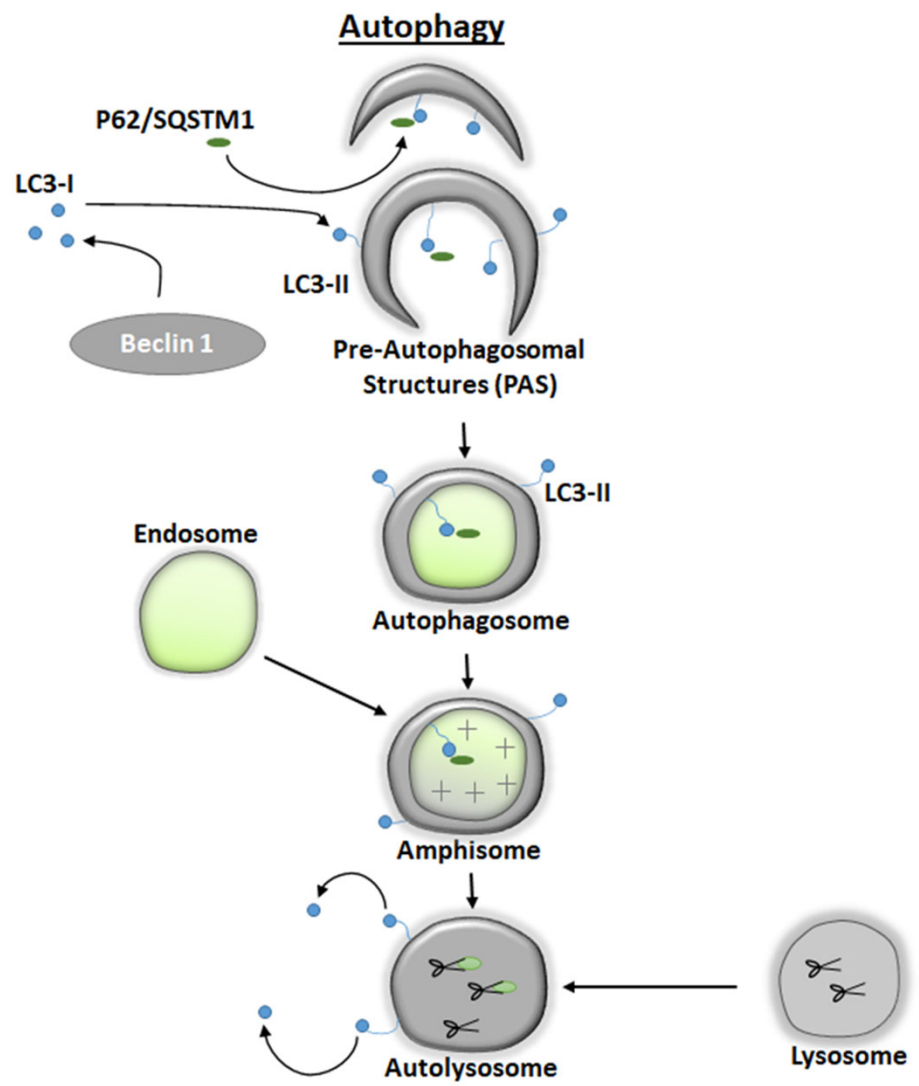


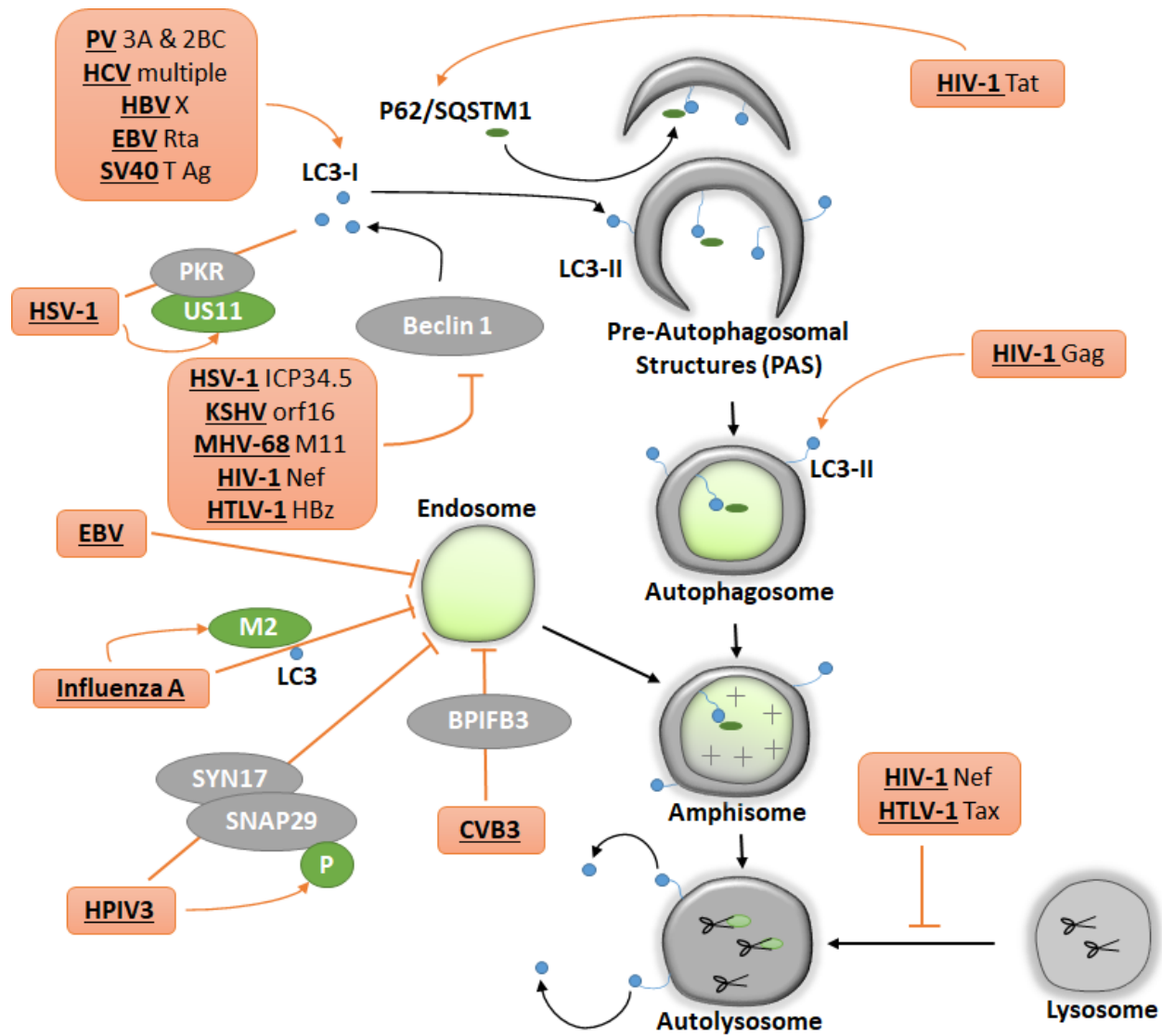
HIV-1 budding (in green) from cultured lymphocyte. Credit: C Goldsmith, P Feorino, EL Palmer, WR McManus, Centers for Disease Control and Prevention (CDC)

HIV-1 and HAND

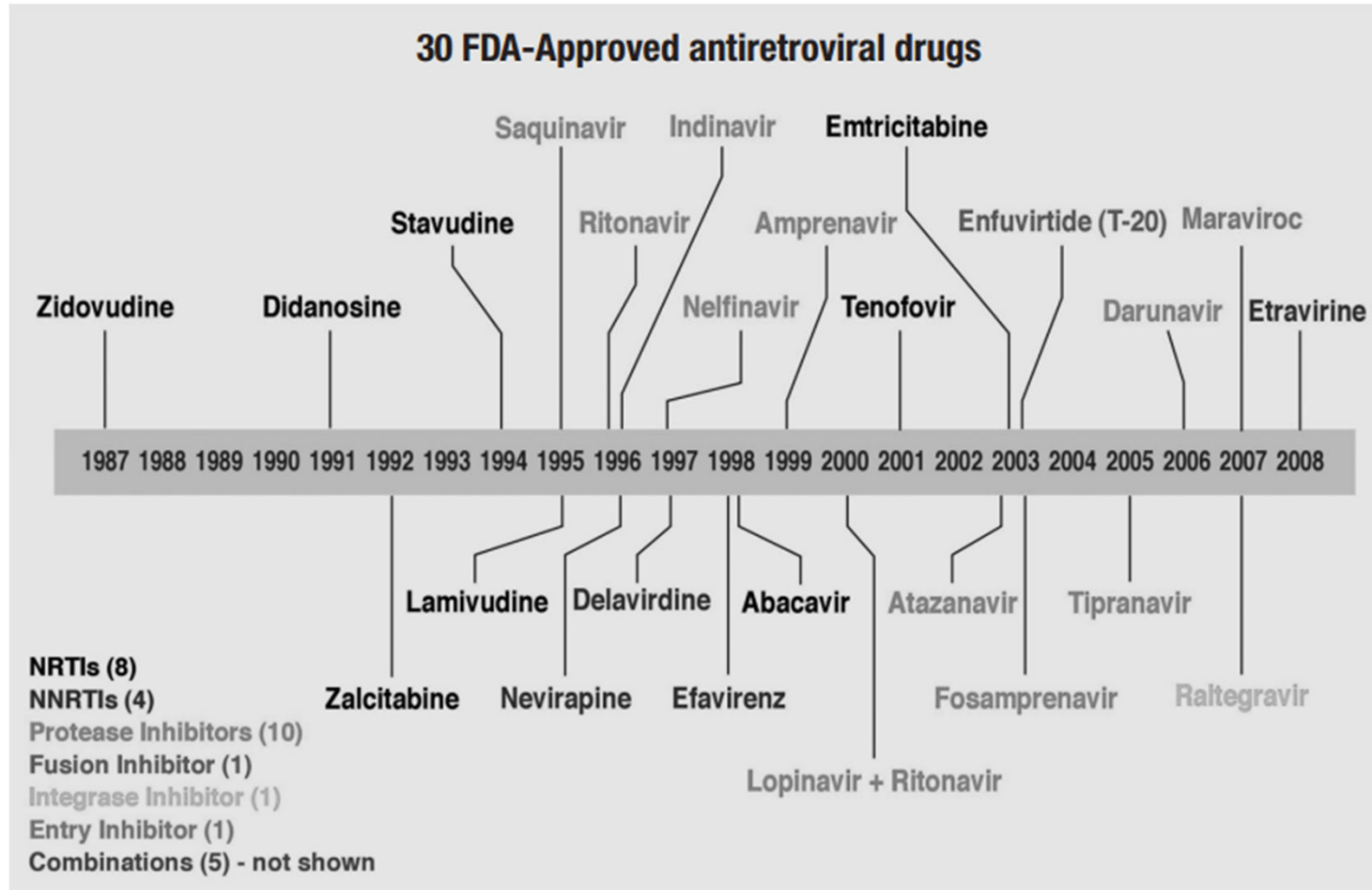


HIV-associated dementia (HAD), asymptomatic neurocognitive impairment (ANI), and mild neurocognitive disorder (ANI); from Justin McArthur at Hopkins

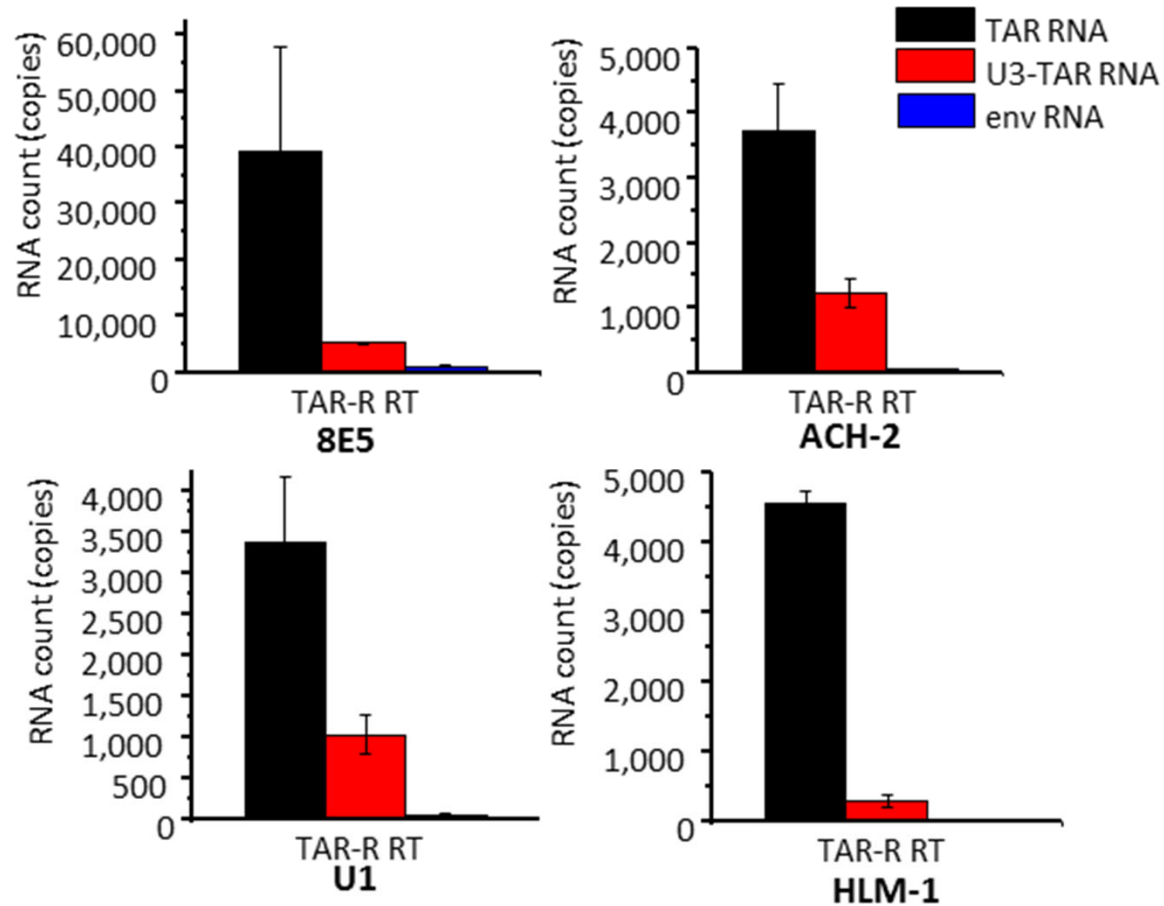




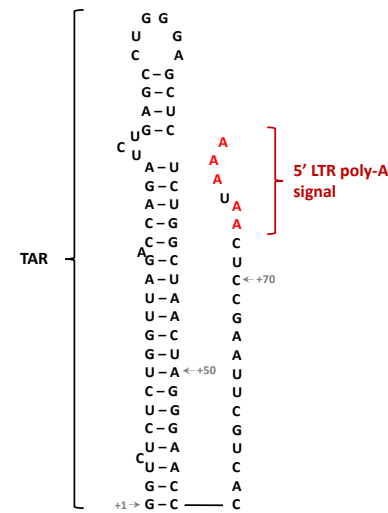
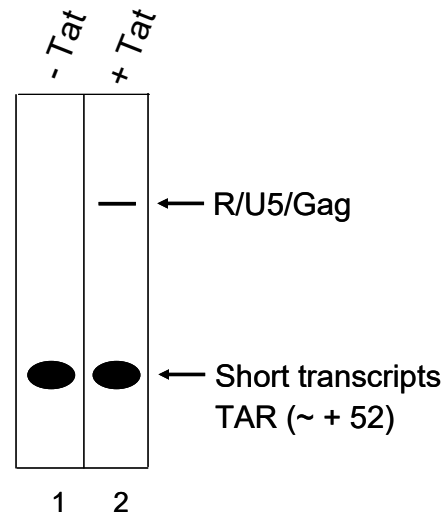
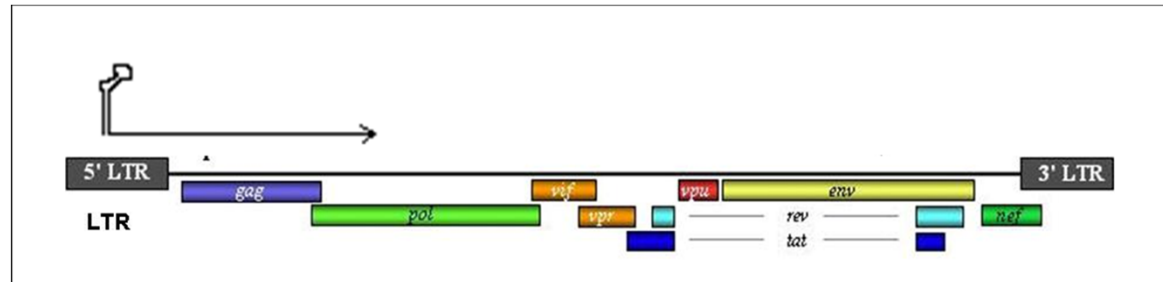
No FDA-approved transcription inhibitors



TAR RNA is present in culture supernatants of HIV-1 infected cells

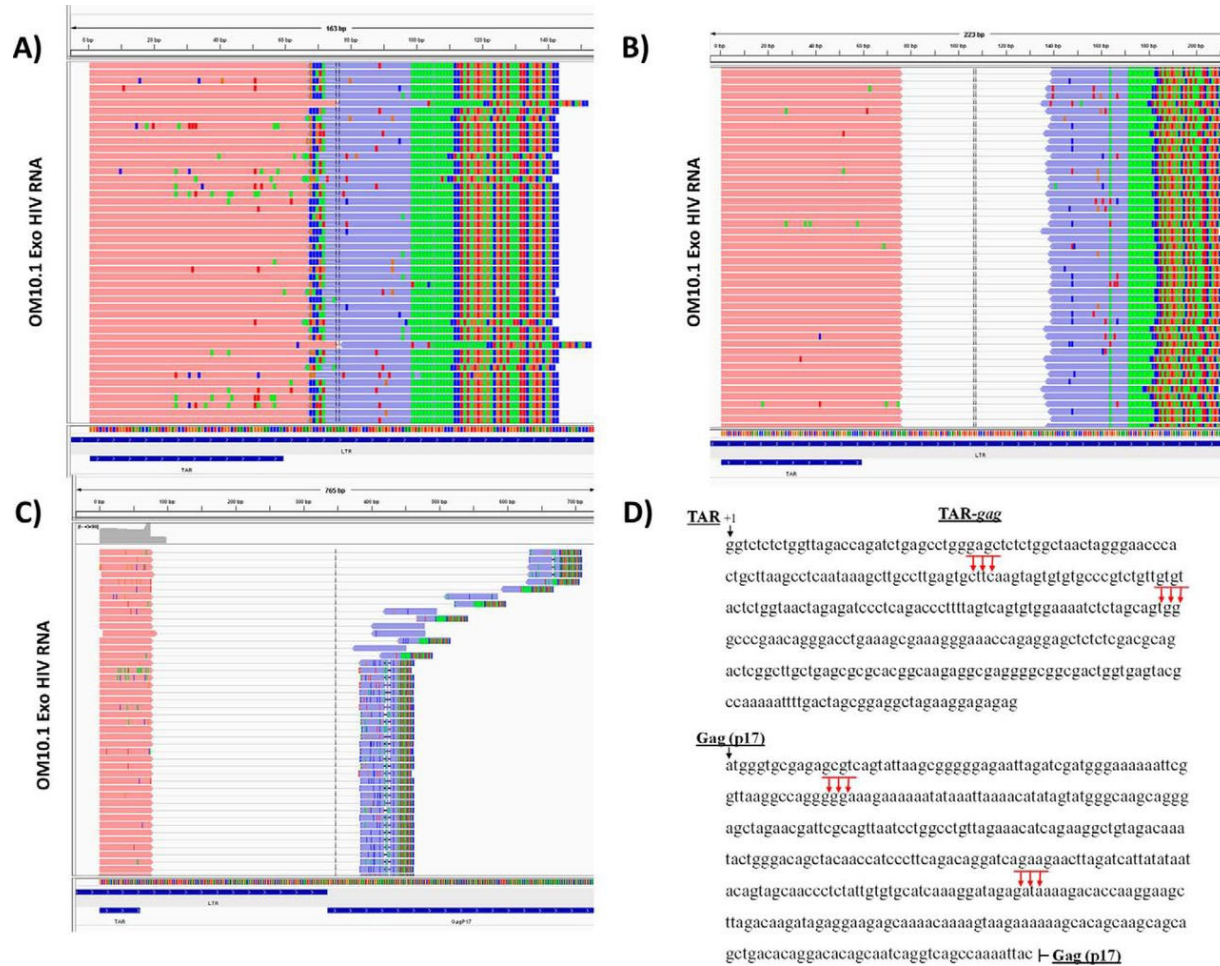


TAR RNA (short transcripts)



- More than 70 papers from late 80s to present
- Detected in vitro, in vivo and latent patient cells

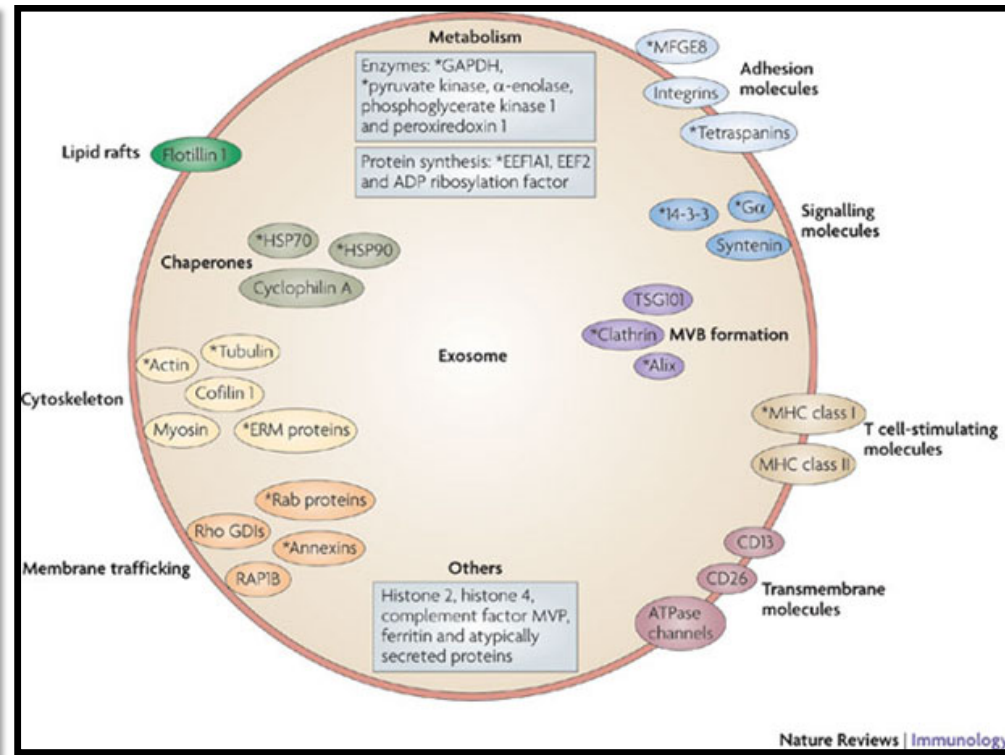
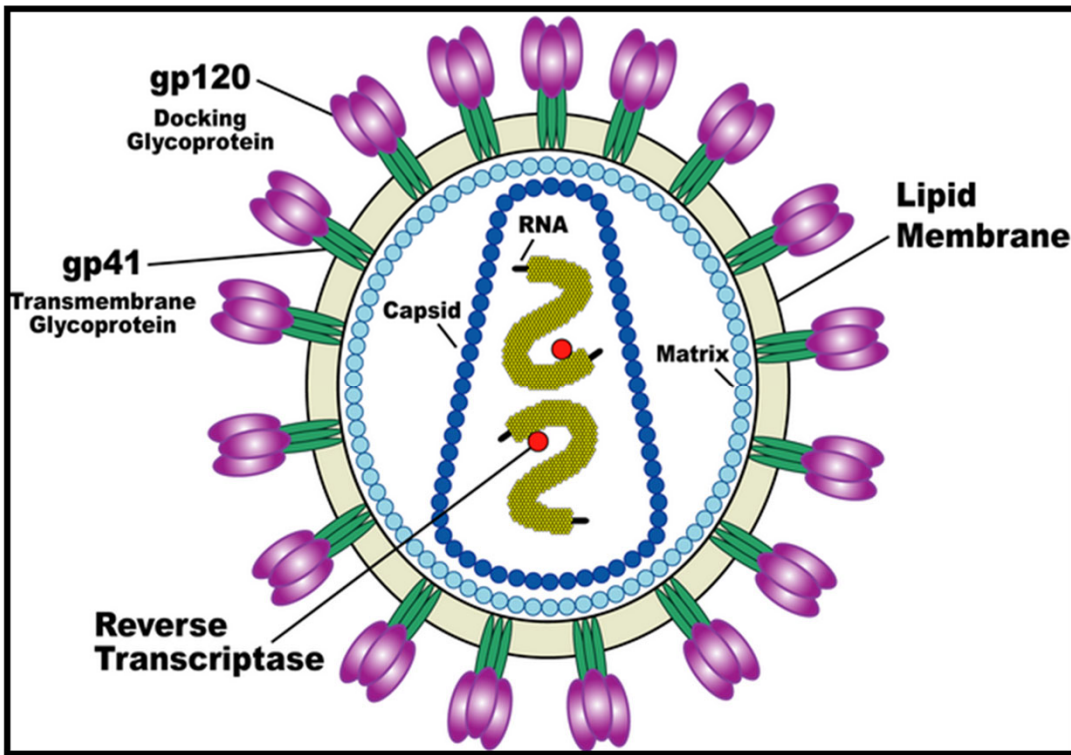
Sequence alignment of exosomal RNA using Integrated Genomics Viewer



Barclay et al., J Biol Chem. 2017.



Membrane vesicles as conveyors of immune responses



They et al., Nat Rev Immunol. 2009.



OPEN

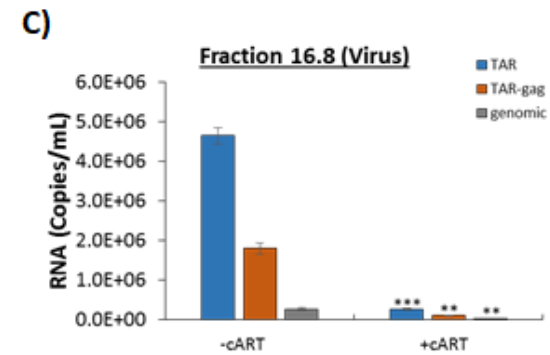
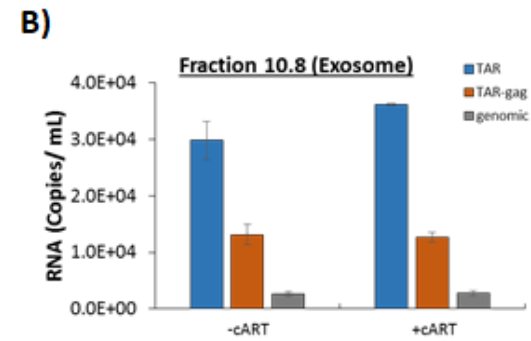
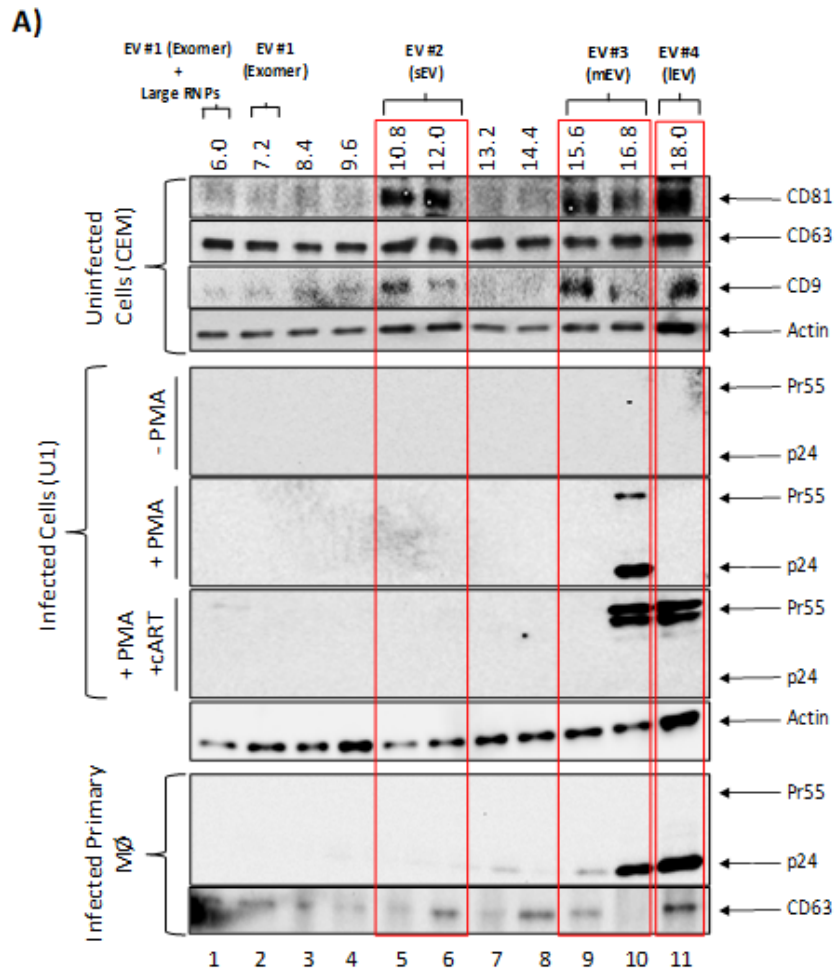
Antiretroviral Drugs Alter the Content of Extracellular Vesicles from HIV-1-Infected Cells

Received: 3 January 2018
Accepted: 1 May 2018
Published online: 16 May 2018

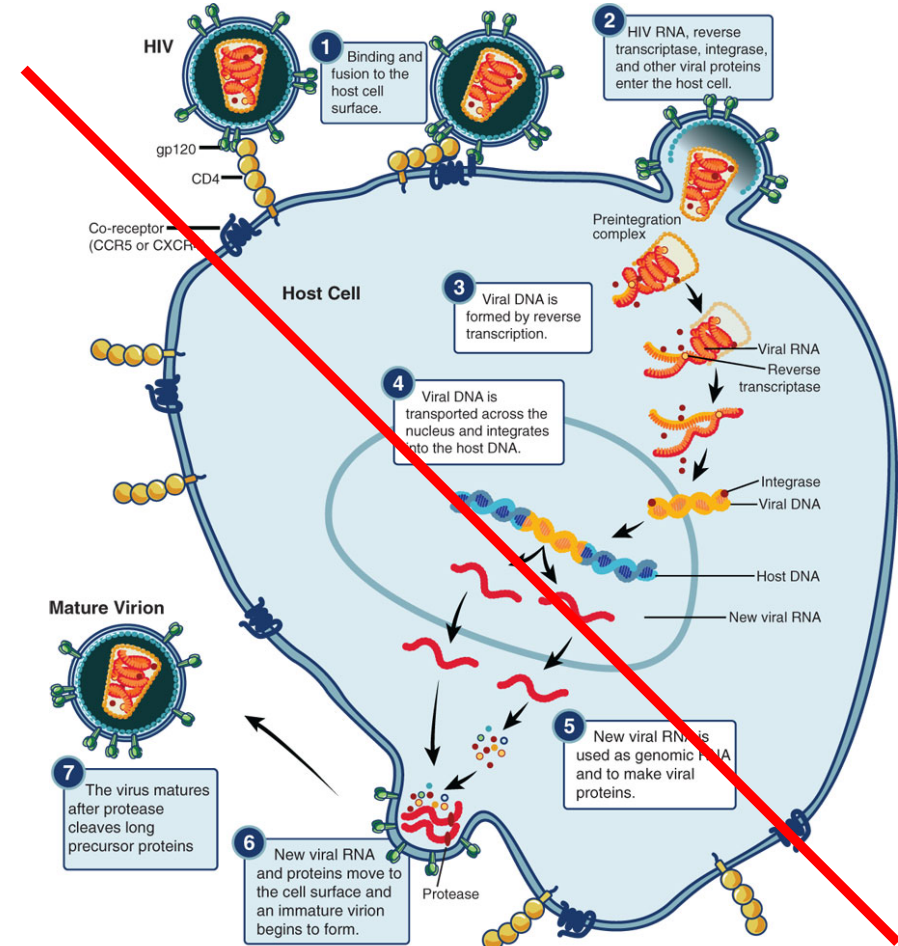
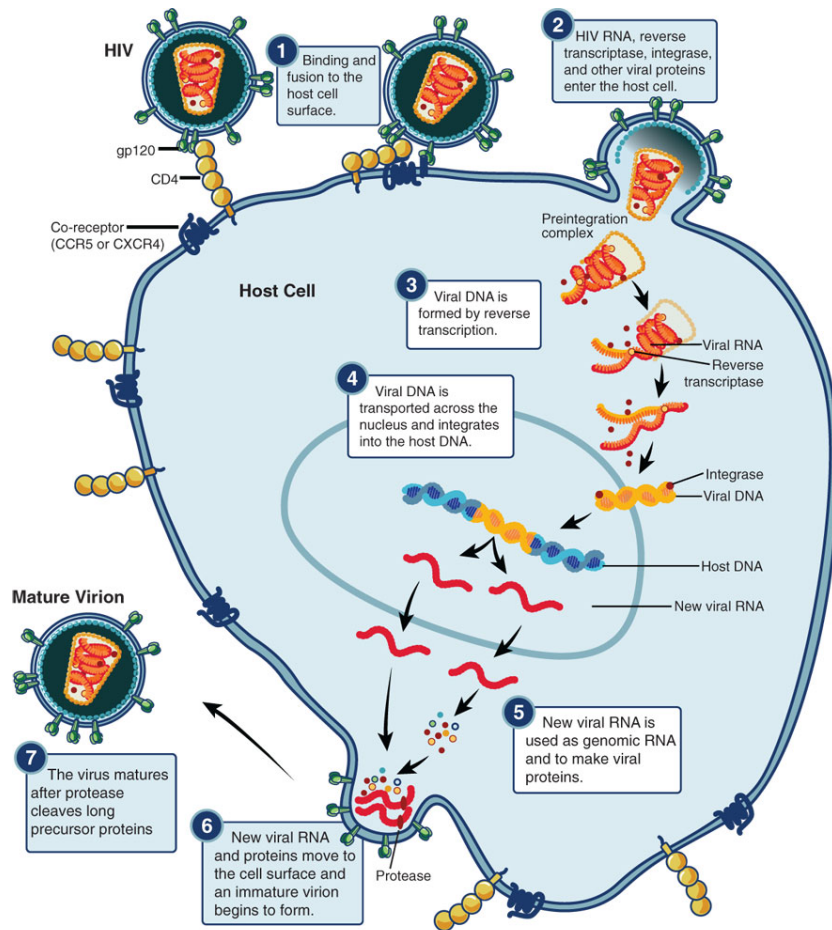
Catherine DeMarino¹, Michelle L. Pleet¹, Maria Cowen¹, Robert A. Barclay¹, Yao Akpamagbo¹, James Erickson¹, Nicaise Ndembu², Manhattan Charurat², Jibreel Jumare², Sunday Bwala³, Peter Alabi⁴, Max Hogan⁵, Archana Gupta⁵, Nicole Noren Hooten ⁶, Michele K. Evans⁶, Benjamin Lepene⁷, Weidong Zhou⁸, Massimo Caputi⁹, Fabio Romerio², Walter Royal 3rd¹⁰, Nazira El-Hage¹¹, Lance A. Liotta⁸ & Fatah Kashanchi¹

To date, the most effective treatment of HIV-1 is a combination antiretroviral therapy (cART), which reduces viral replication and reverses pathology. We investigated the effect of cART (RT and protease inhibitors) on the content of extracellular vesicles (EVs) released from HIV-1-infected cells. We have previously shown that EVs contain non-coding HIV-1 RNA, which can elicit responses in recipient cells. In this manuscript, we show that TAR RNA levels demonstrate little change with the addition of cART treatment in cell lines, primary macrophages, and patient biofluids. We determined possible mechanisms involved in the selective packaging of HIV-1 RNA into EVs, specifically an increase in EV-associated hnRNP A2/B1. More recent experiments have shown that several other FDA-approved drugs have the ability to alter the content of exosomes released from HIV-1-infected cells. These findings on cART-altered EV content can also be applied to general viral inhibitors (interferons) which are used to treat other chronic infections. Additionally, we describe unique mechanisms of ESCRT pathway manipulation by antivirals, specifically the targeting of VPS4. Collectively, these data imply that, despite antiretroviral therapy, EVs containing viral products are continually released and may cause neurocognitive and immunological dysfunction.

Isolation of EVs away from virus

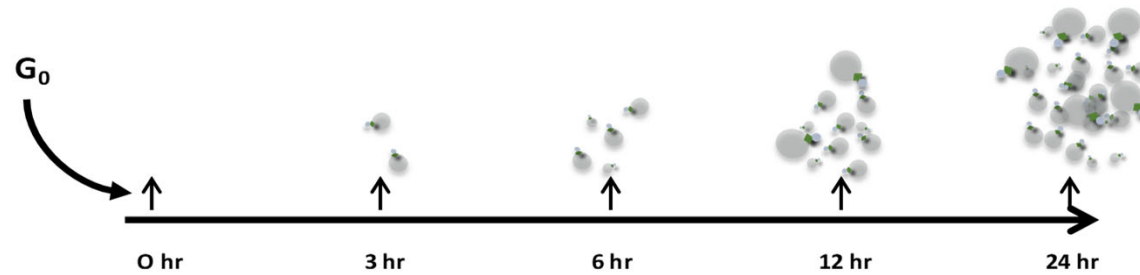


HIV life cycle

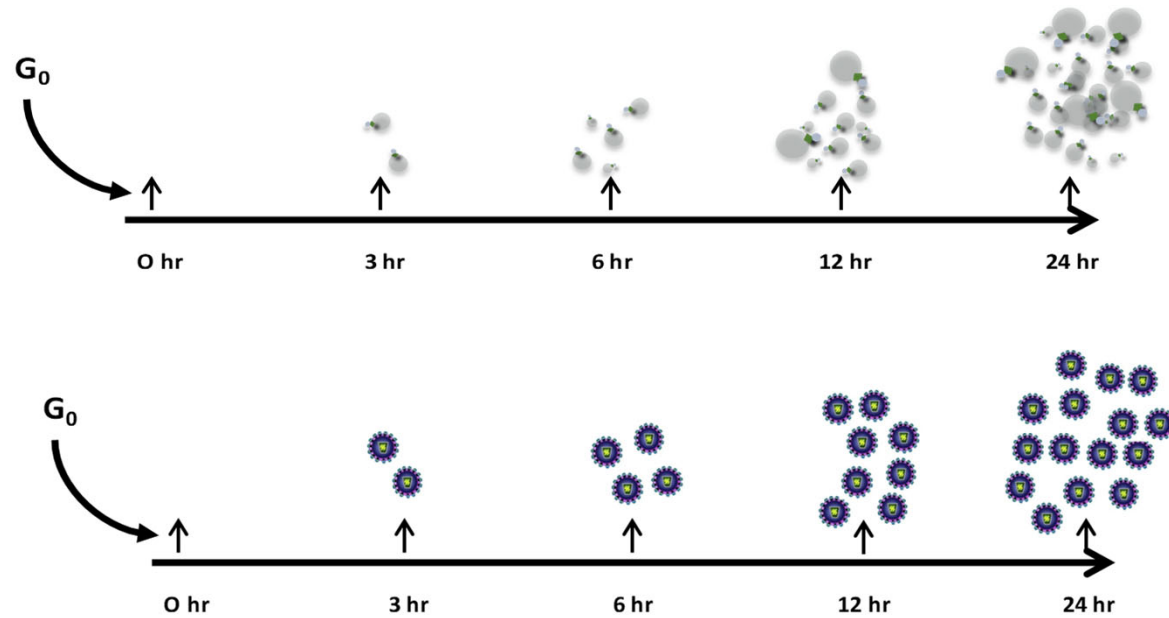


Source: J.L. Jameson, A.S. Fauci, D.L. Kasper, S.L. Hauser, D.L. Longo, J. Loscalzo: Harrison's Principles of Internal Medicine, 20th Edition Copyright © McGraw-Hill Education. All rights reserved.

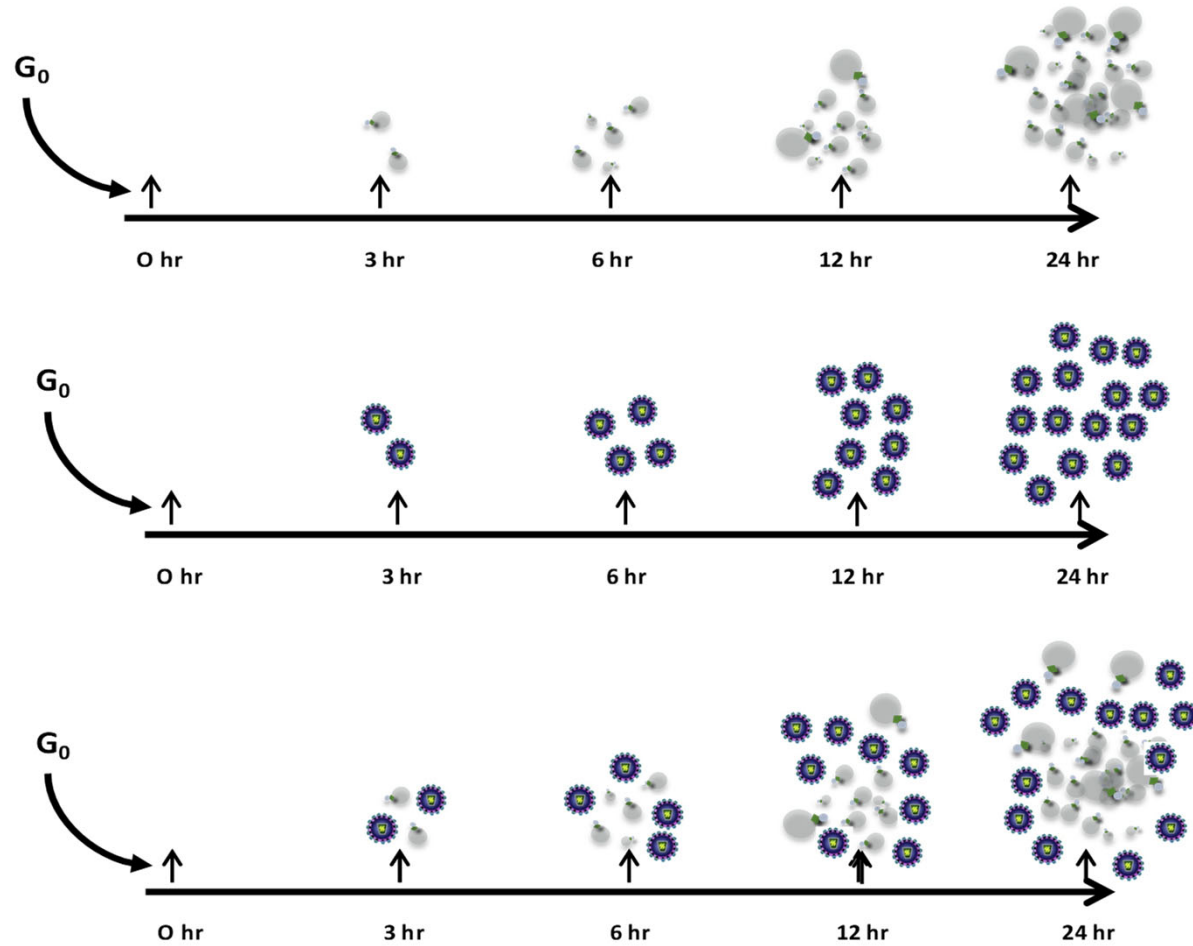
Question: Which particles are released first from infected cells- EVs or Viruses ?



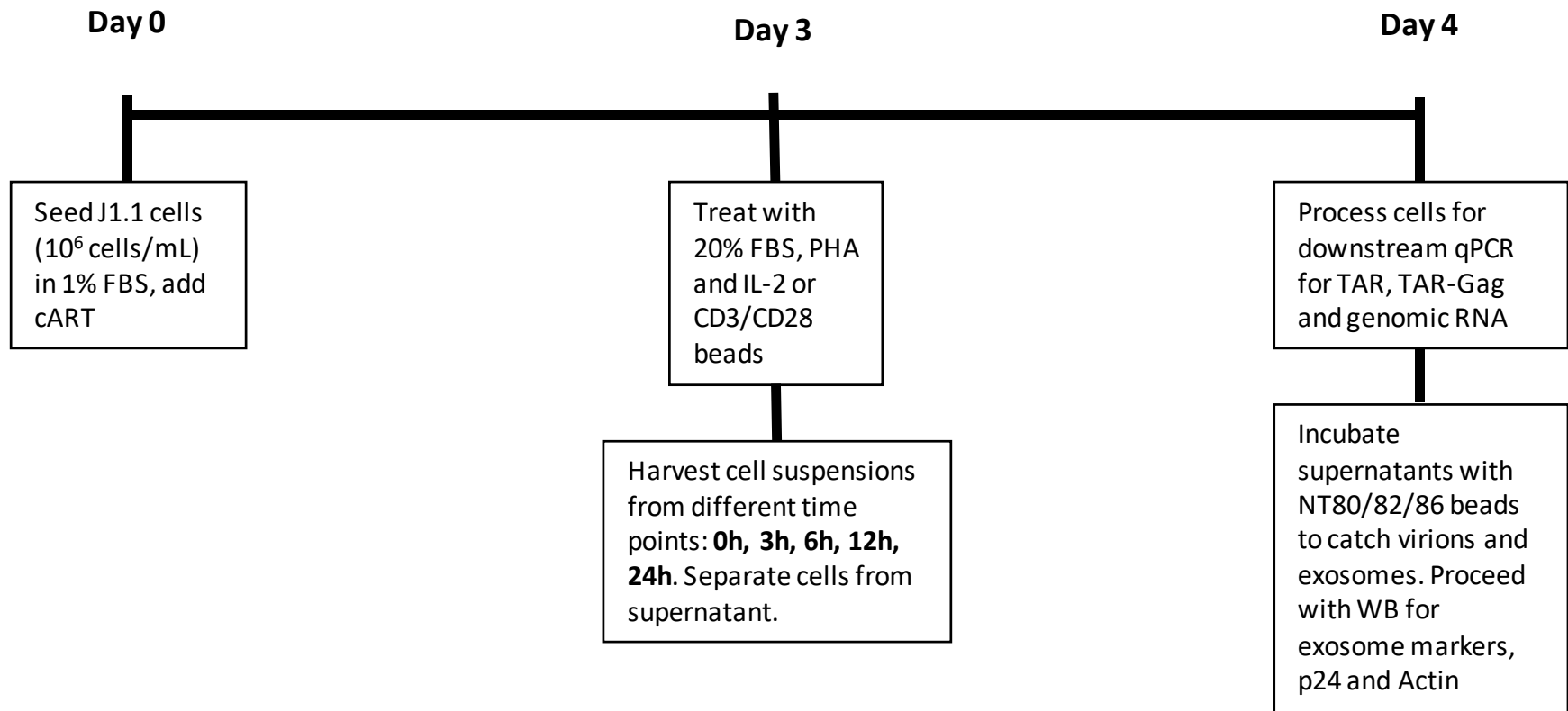
Question: Which particles are released first from infected cells- EVs or Viruses ?

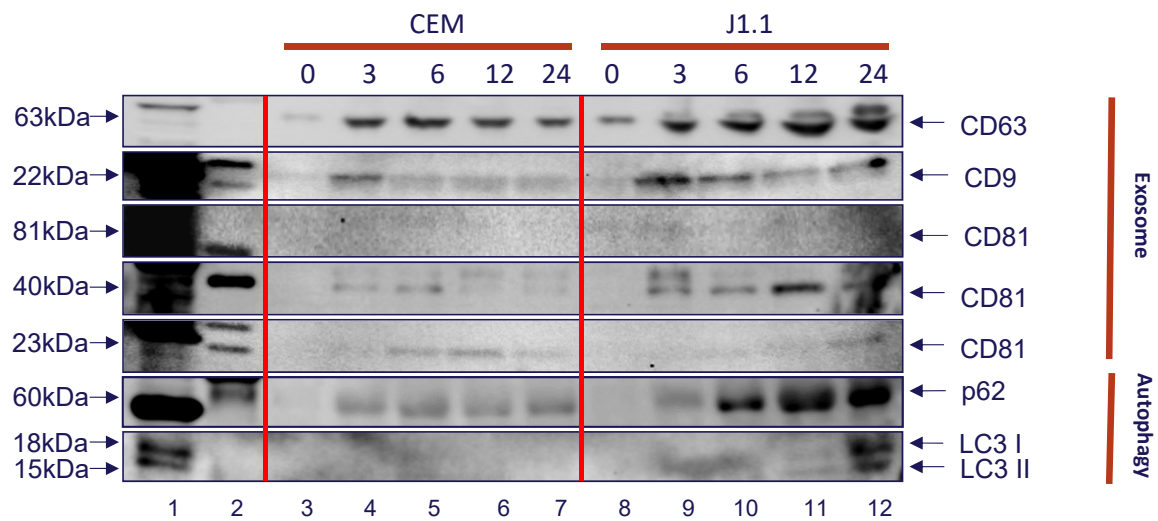


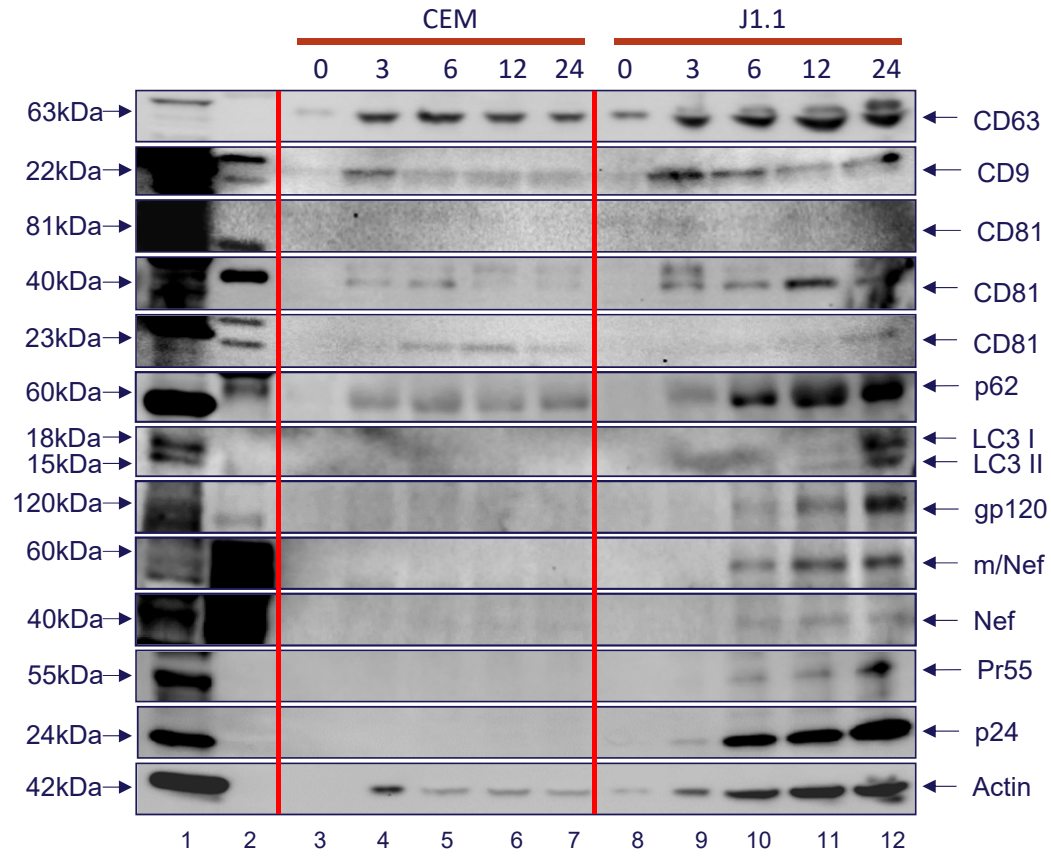
Question: Which particles are released first from infected cells- EVs or Viruses ?



Experimental procedure

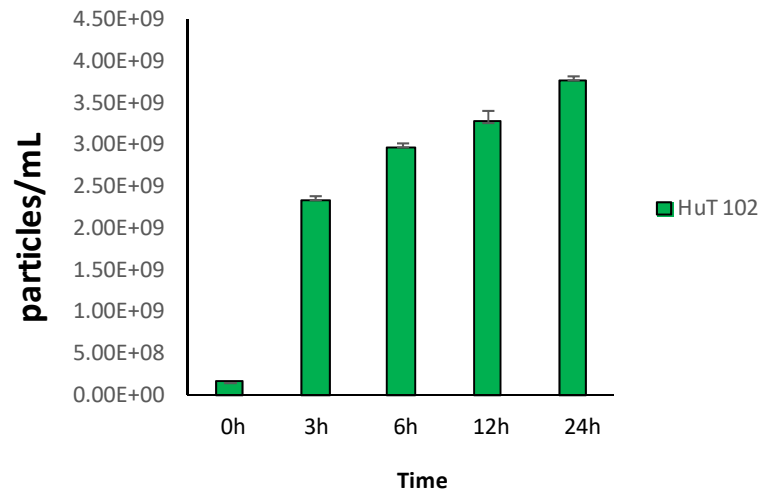
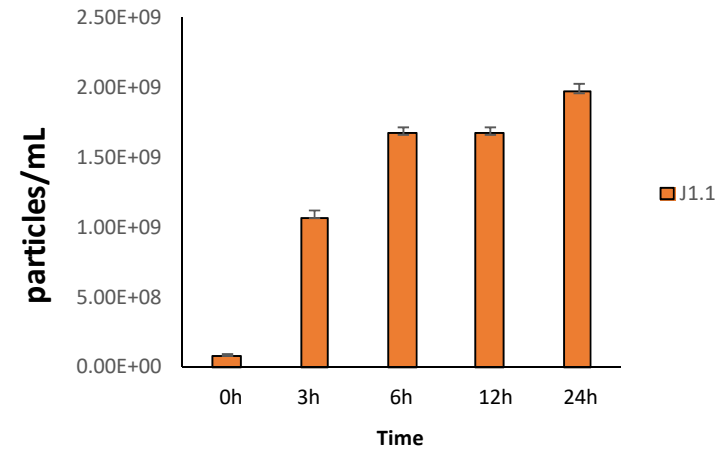
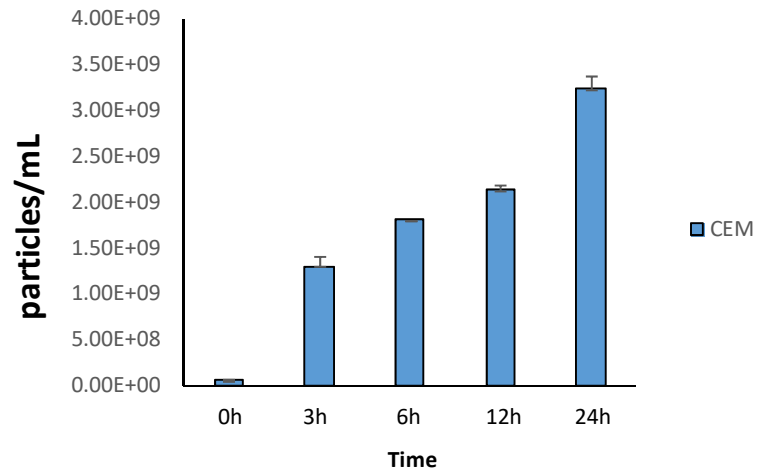




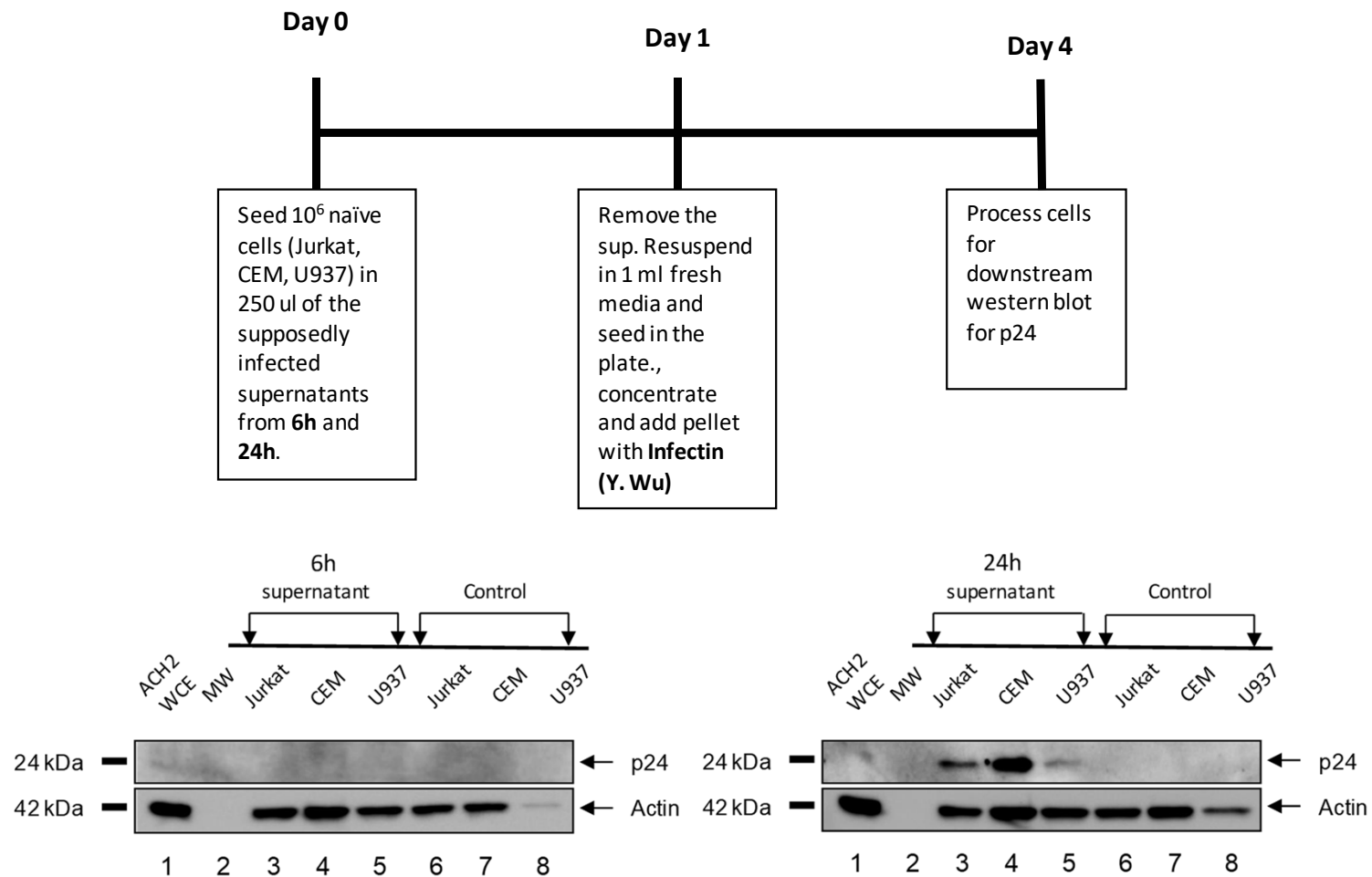


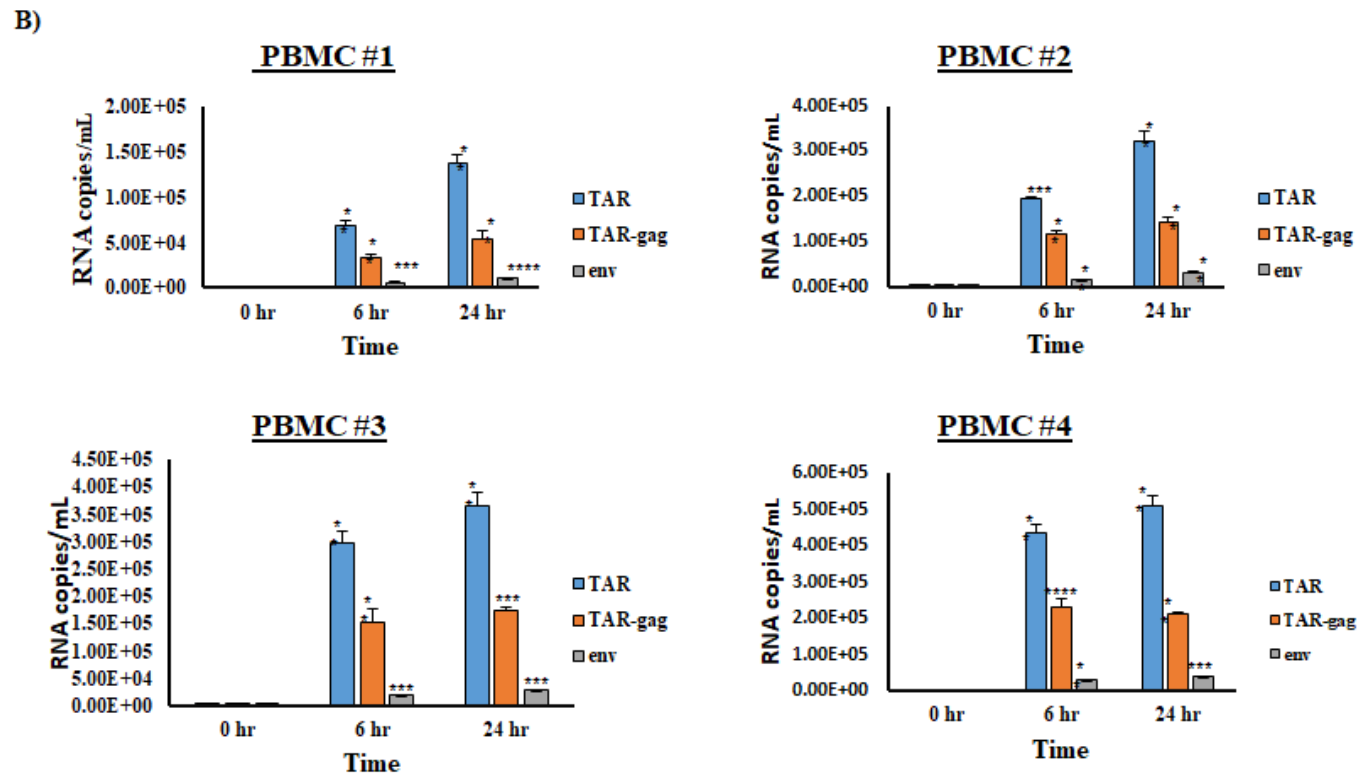
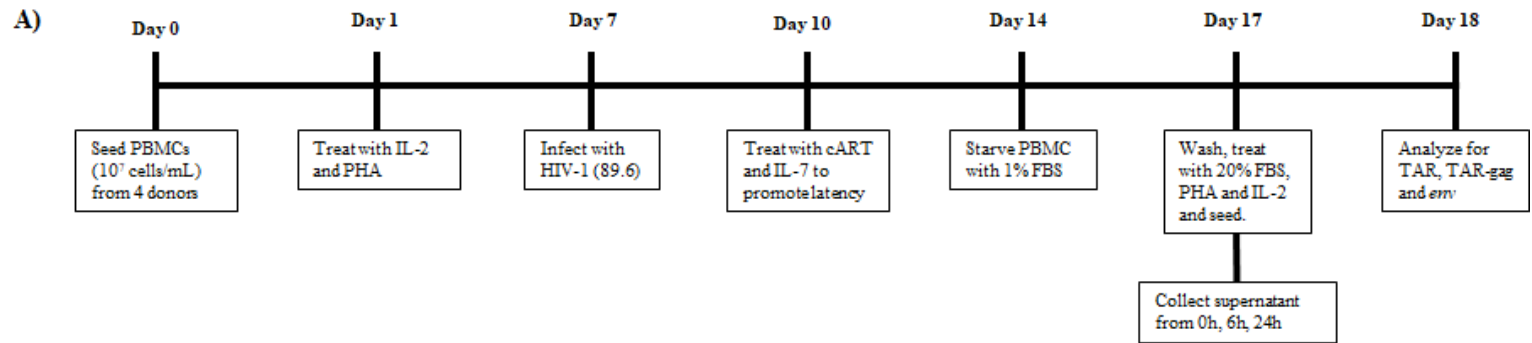
Exosome
Autophagy
Virus

Particle release over time

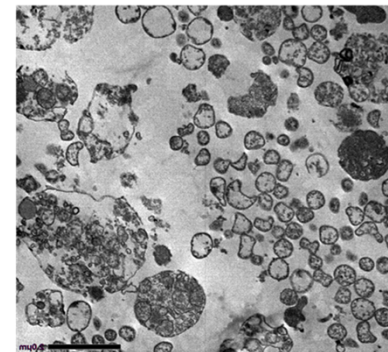
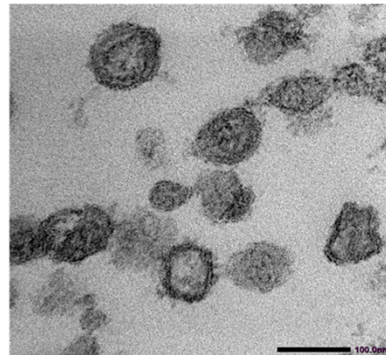
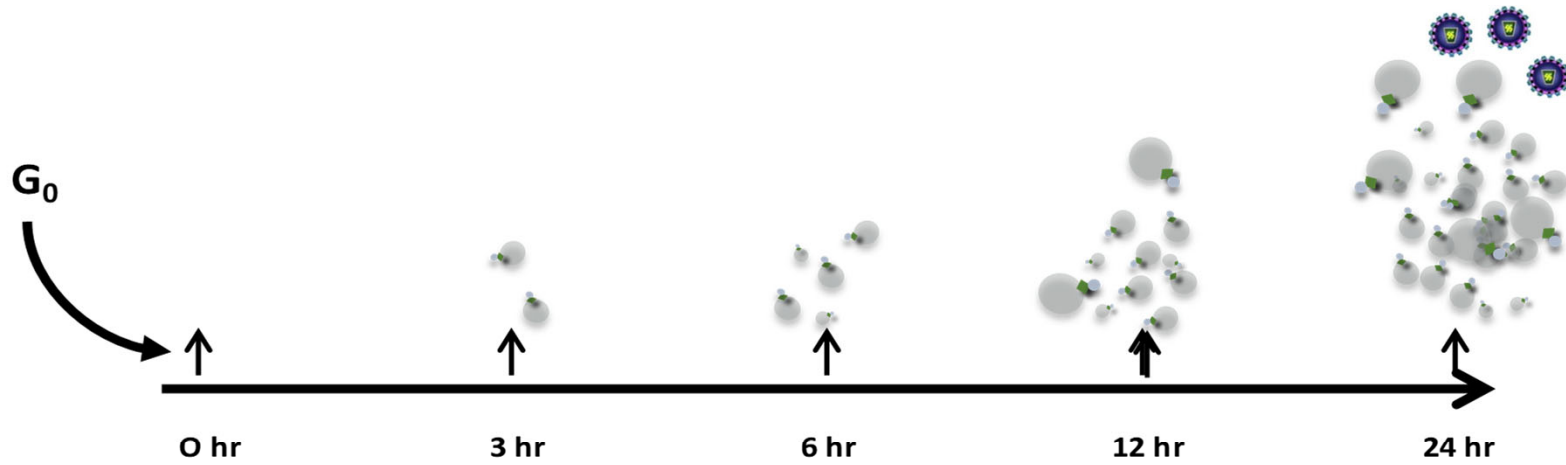


Virus rescue assay





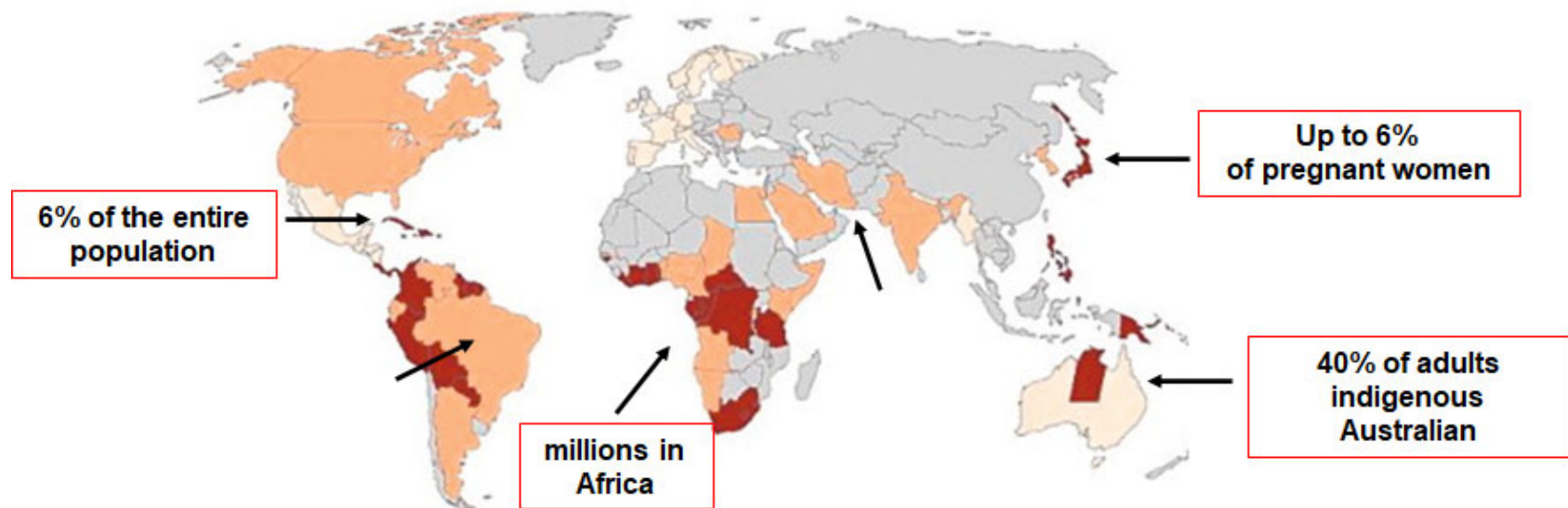
Answer: EVs come out of infected cells first and then a mixture of both EVs and viruses at a later time



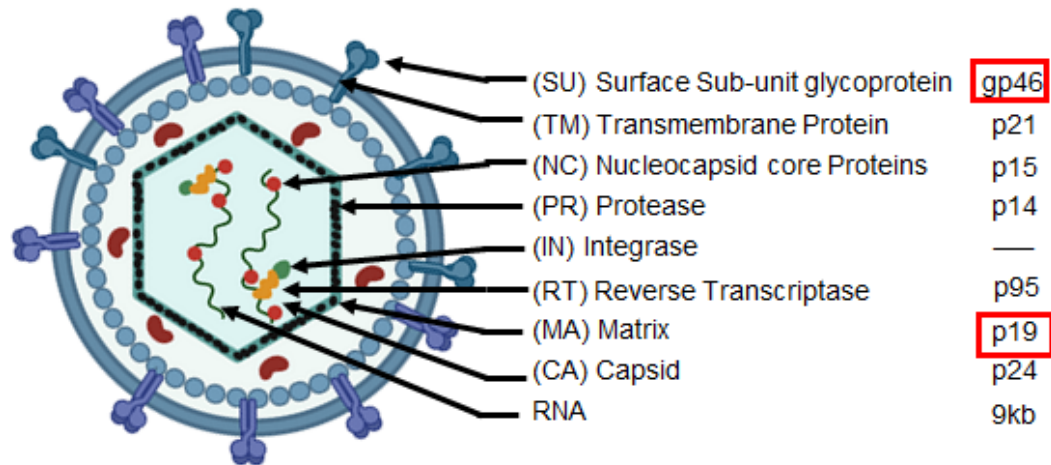
Human T-lymphotropic virus type 1 (HTLV-1)

Epidemiology

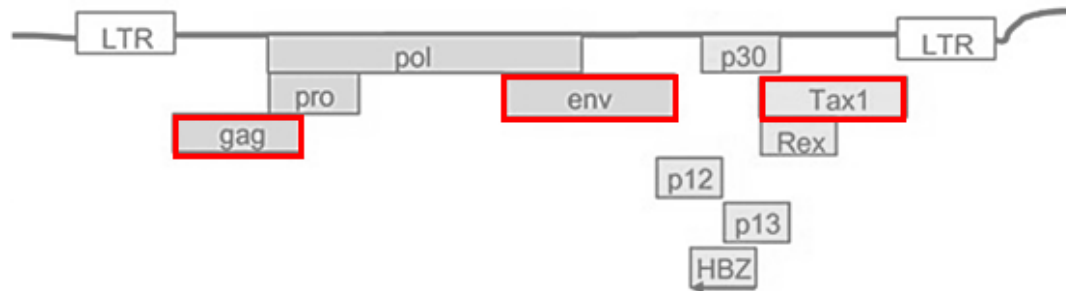
- Approximately **5-10 million** patients worldwide infected with HTLV-1.
- The true figure is greater as many cases are unreported from highly populated regions, such as **China, India, Northwest Africa (i.e., the Arab Maghreb), and East Africa.**
- It is estimated that the actual global prevalence may be as **high as 20 million** worldwide.



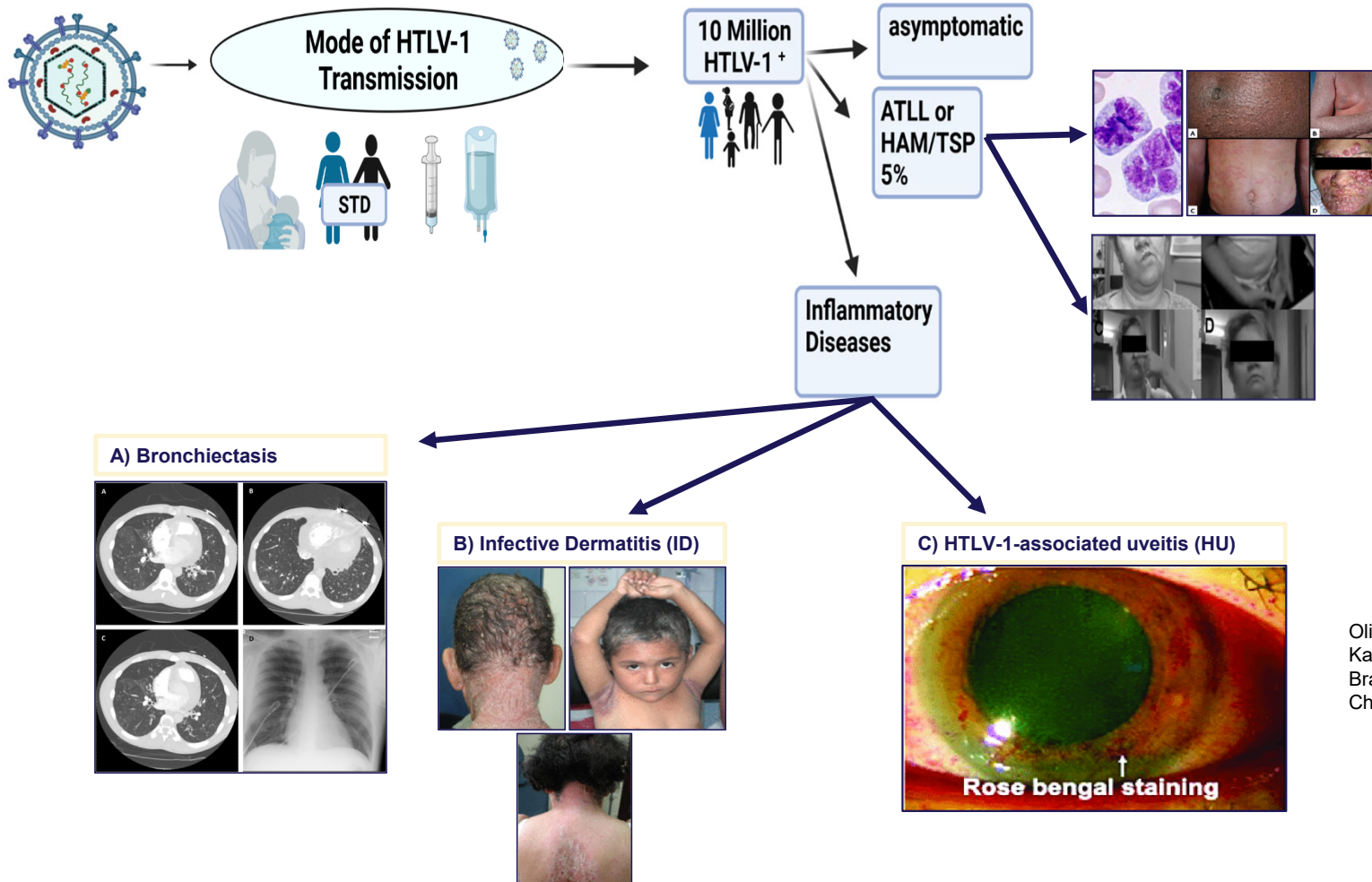
A) VIRION



B) HTLV-1 Genome Structure

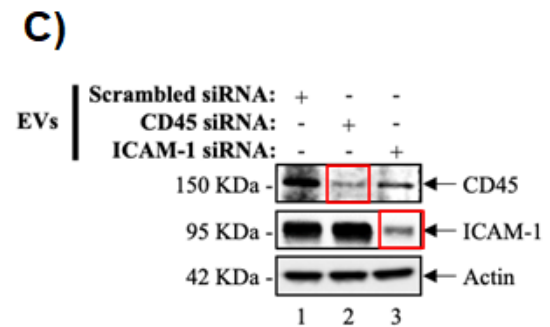
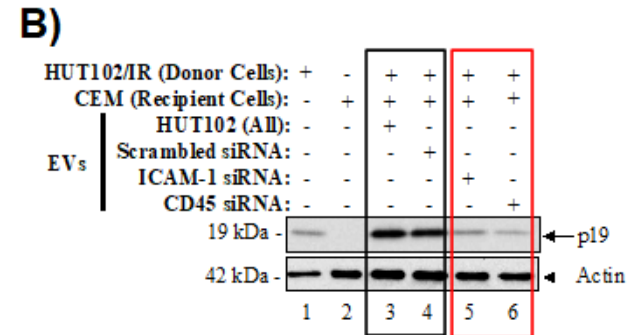
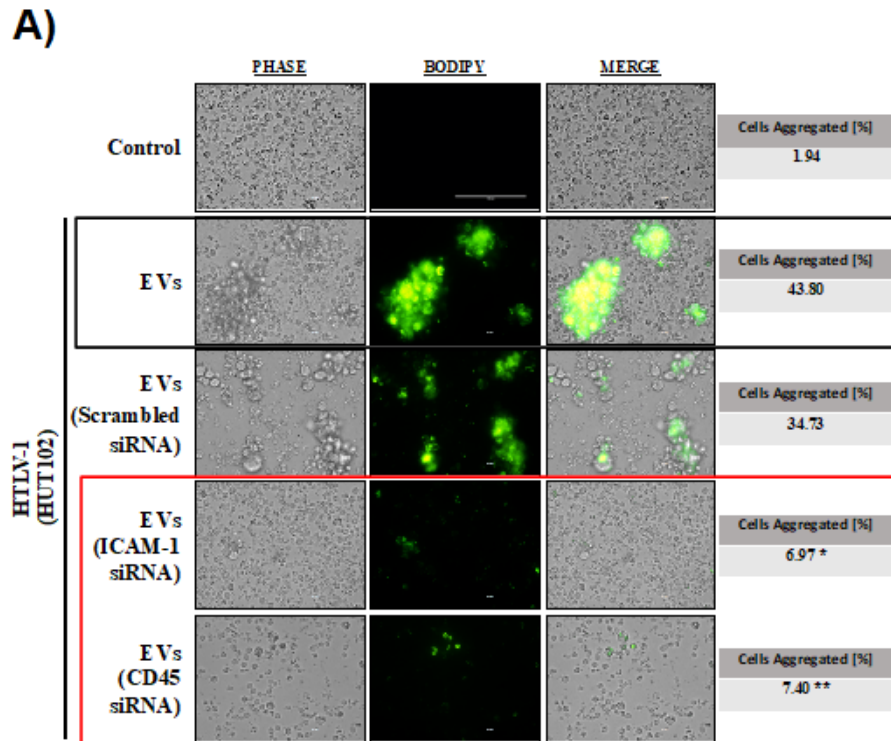


HTLV-1 overview

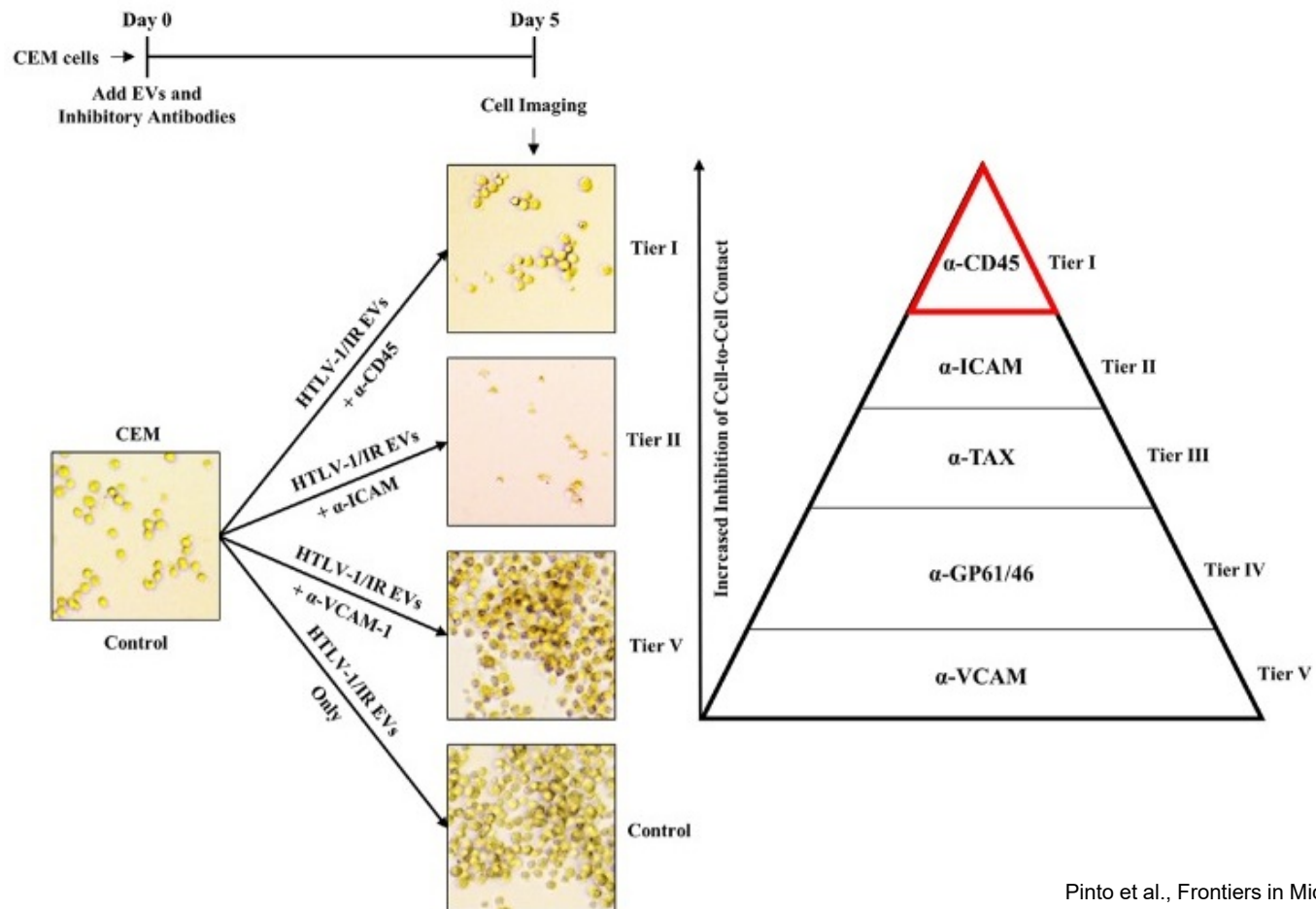


Oliveira PD et al., Rev Assoc Médica Bras. 2016.
 Kamoi K., Frontiers in Microbiology. 2020.
 Bravo FG., Sem Diagnostic Pathol. 2020.
 Chiong et al., IDCases, 19, e00714.

Inhibition of ICAM-1 and CD45 via small interfering RNA prevents cell-to-cell contact



Antibodies against specific cellular surface receptors inhibit cell-to-cell contact

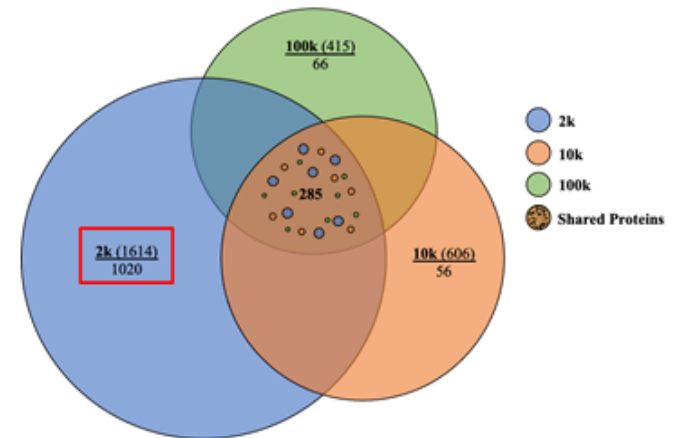


Proteomic analysis of the 2k, 10k, and 100k showed that HTLV-1 EVs carry viral/host proteins

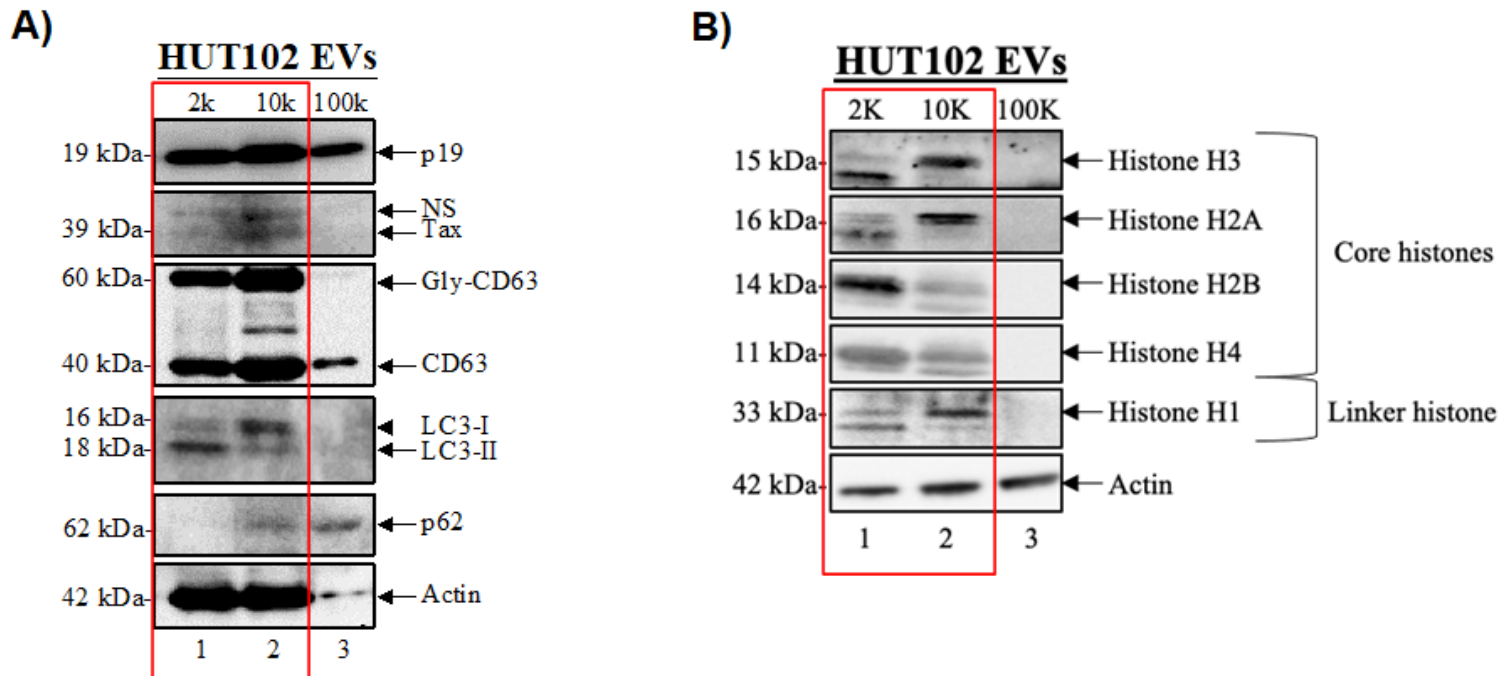
A)

Viral Proteins			
EVs	Reference	Hits	Sequence (Most Abundant Peptide)
2k	Polyprotein (Gag-Pro-Pol)	57	KPPPNQPF ^R
	Gag (p19)	15	IALETPVWIPINYSLLASLLPK
	Envelope	11	PYLGQSWTPYTGAVSSPYWK
	Tax	4	VIGSALQFLIP ^R
10k	Polyprotein (Gag-Pro-Pol)	28	KPPPNQPF ^R
	Gag (p19)	16	IALETPVWIPINYSLLASLLPK
	Gag-Pol	9	DQIYLNPSQVQSLVQL ^R
	Envelope	5	NGGGYY SASYSDPSL ^K
100k	Protease	5	NTSVLGAGGQTQDHF ^K
	Polyprotein (Gag-Pro-Pol)	30	DPSWASILQGLEEPPYHAFV ^E R
	Gag (p19)	12	IALETPVWIPINYSLLASLLPK; VQANNPQQQL ^R
Envelope	2	NGGGYY SASYSDPSL ^K	

B)

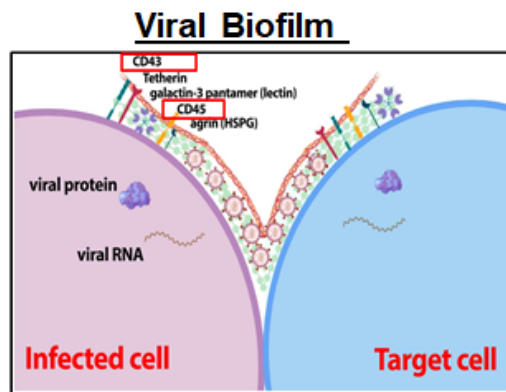
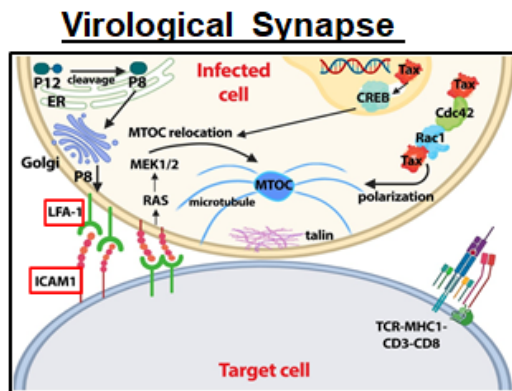


2k & 10k HTLV-1 EVs contain highest amount of viral/host proteins

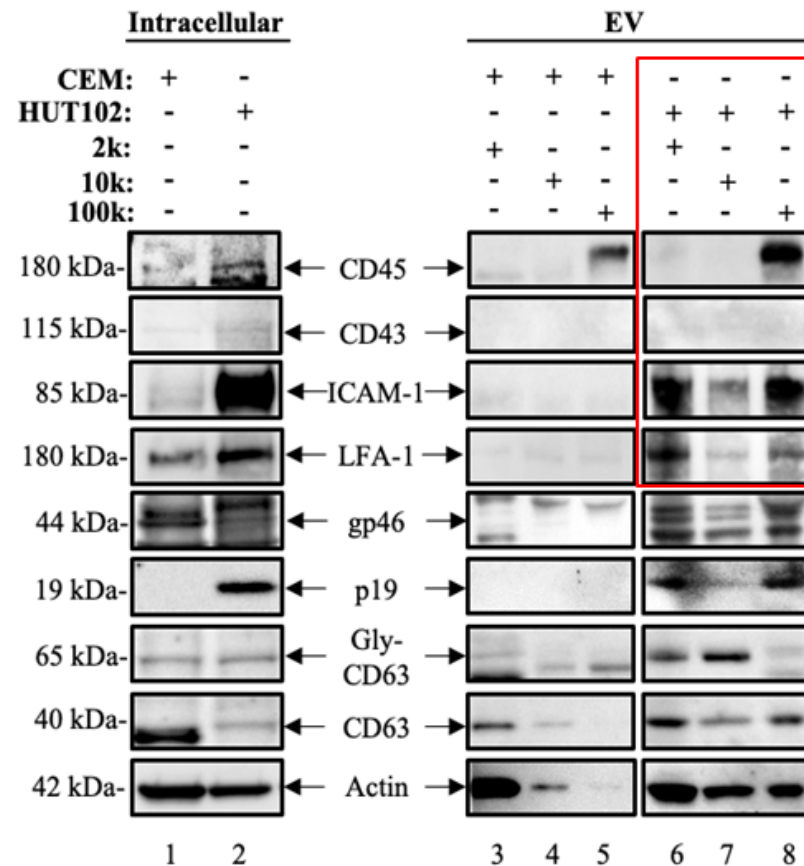


HTLV-1 EVs contain CD45, ICAM-1 and LFA-1 may contribute to enhance cell-to-cell contact

A)

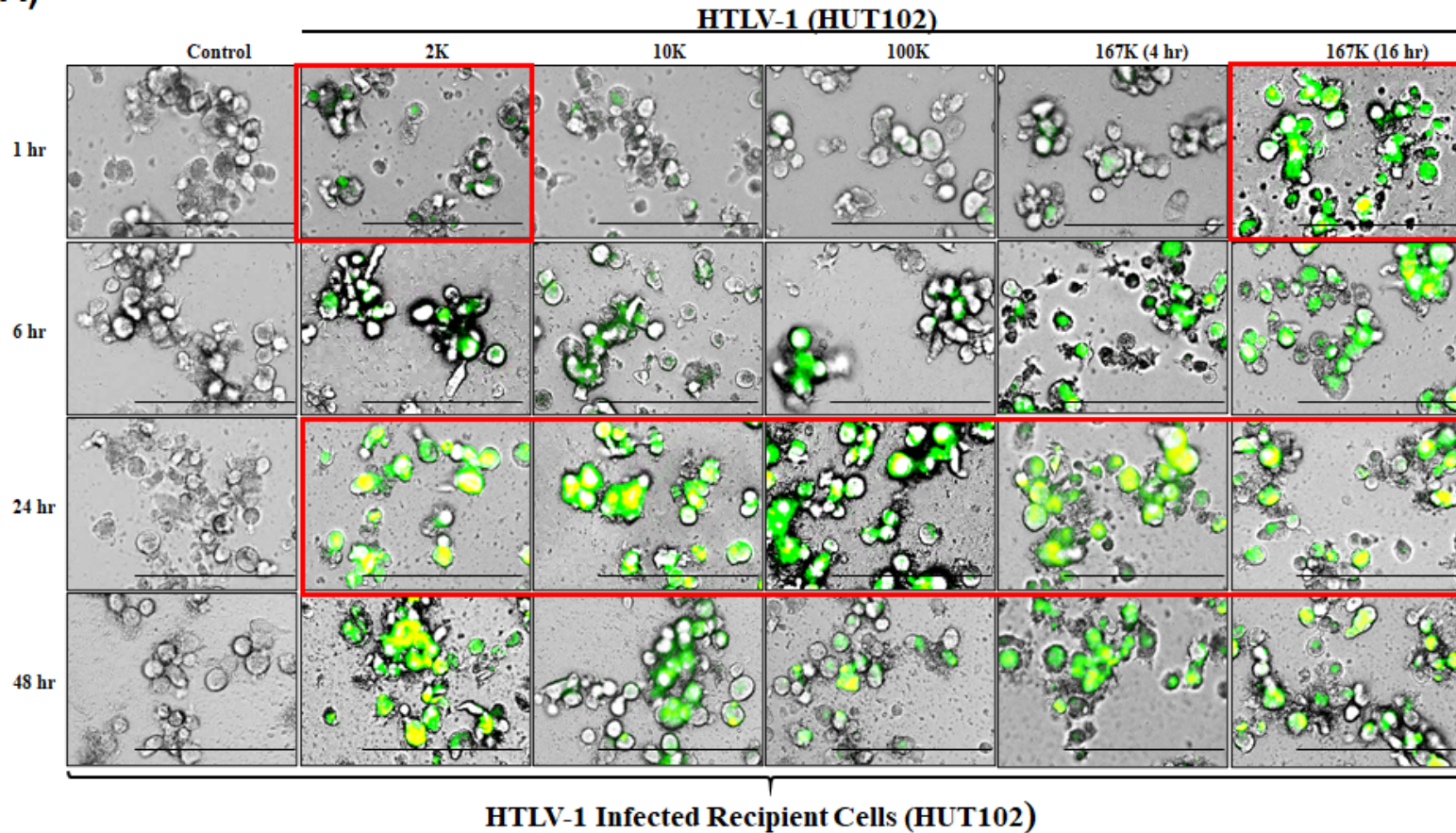


B)

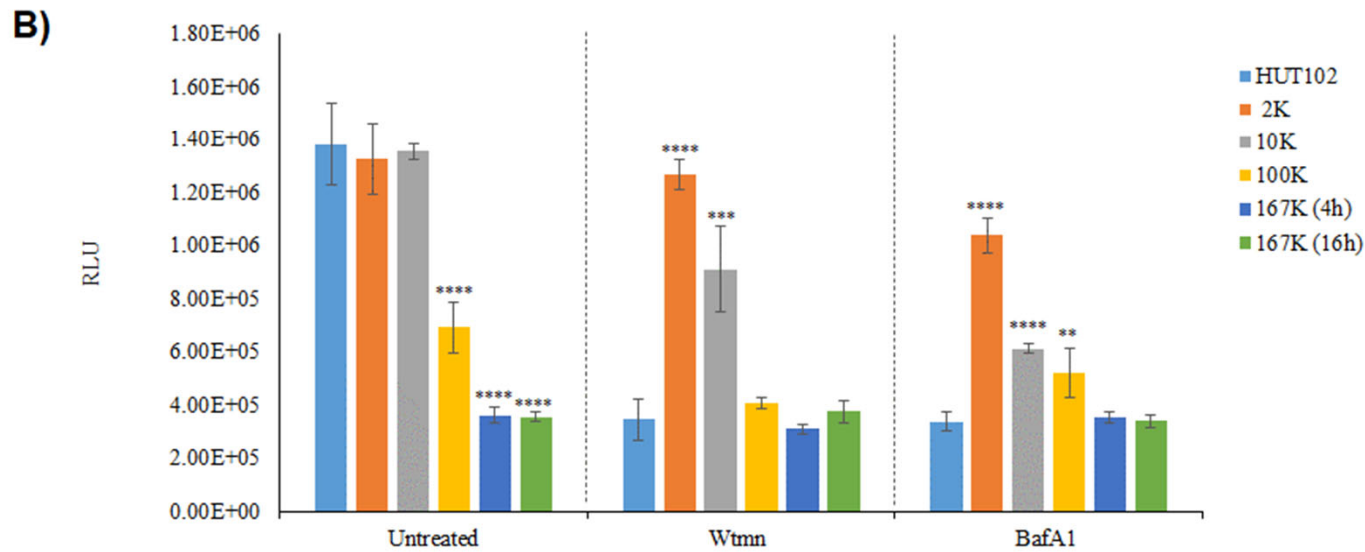


2K & 167K (16 hr) HTLV-1 EV uptake occurs before other EV subpopulations by the same cells

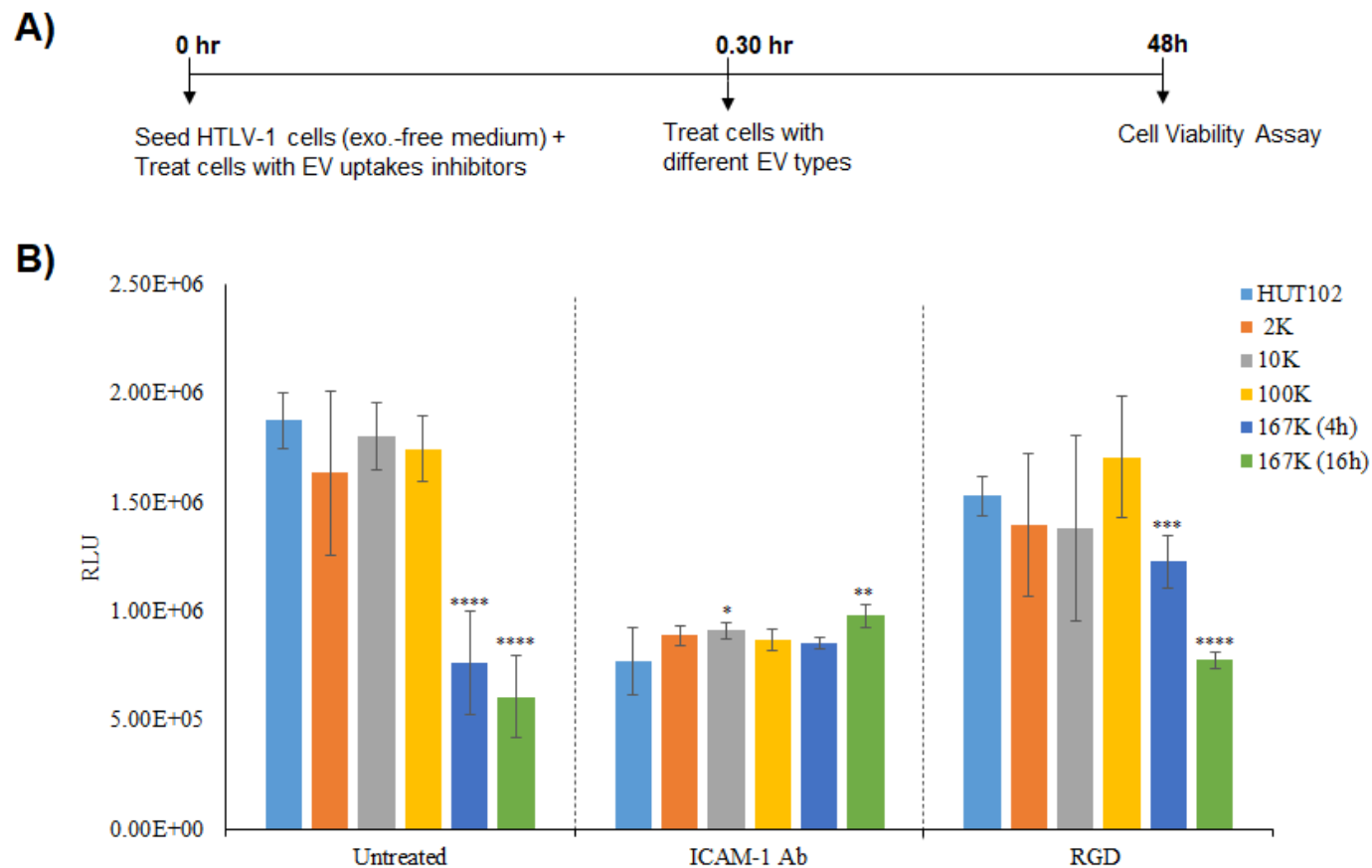
A)



Inhibition of micropinocytosis attenuates 10K and 100K EV internalization and decrease cell viability in HUT102 cells

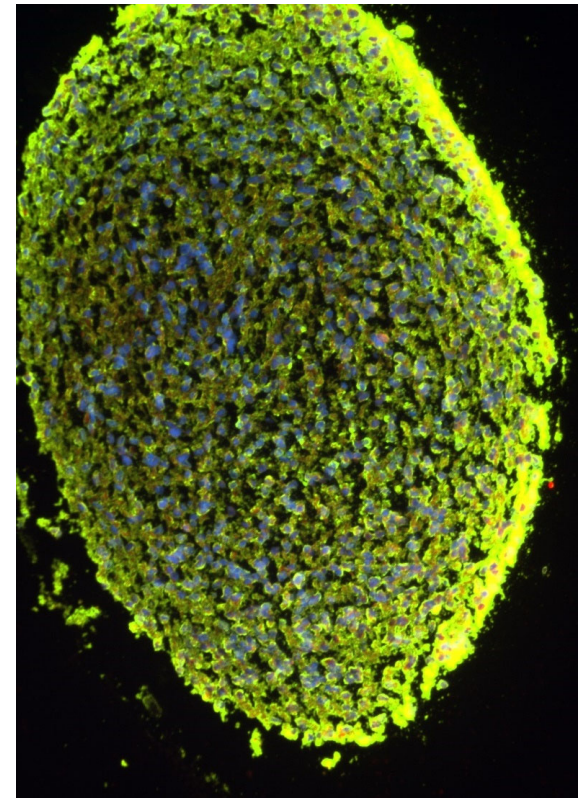


Inhibition of receptor-mediated endocytosis attenuates EV internalization and decrease cell viability in HUT102 cells



Overview

- Update on the most recent literature surrounding EVs from virally infected cells (“Damaging EVs”)
- Generation and infection of iPSC-derived neurospheres
- Effect of stem cell EVs (“Reparative EVs”) on HIV-1 infected neurospheres
- Summary



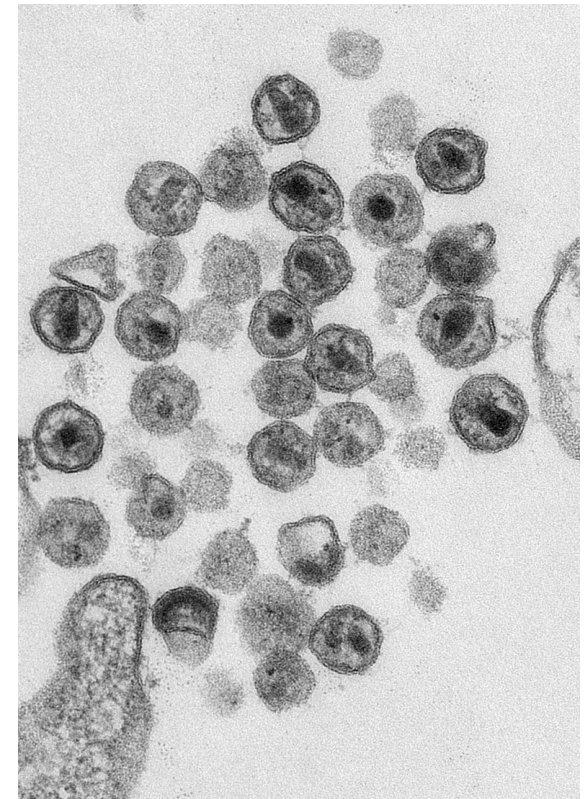
Background

Epidemiology

- HIV-associated neurocognitive disorders (HAND) are commonly associated with HIV-1 infection
- Despite combination antiretroviral therapy (cART) HAND persists in HIV-1 patients (>50%)

Significance

- There is an unmet need to develop novel platforms for CNS disease modeling and therapeutic intervention
- EVs from stem cells have demonstrated reparative properties in a wide range of pathologies, including those related to the CNS



HIV particles; Photo credit: M Metcalfe, T Hodge, CDC

REPORT

Zika virus impairs growth in human neurospheres and brain organoids

Patricia P. Garcez^{2,1,4}, Erick Correia Loliola^{1,1}, Rodrigo Madeiro da Costa^{1,1}, Luiza M. Higa^{3,1}, Pablo Trindade^{1,1}, Rodrigo Delvecchio³, Juliana Minardi Nascimento^{1,4}, Rodrigo Brindeiro³, Amilcar Tanuri³, Stevens K. Rehen^{1,2,*}

SCIENTIFIC REPORTS

OPEN

Zika virus disrupts molecular fingerprinting of human neurospheres

Received: 01 August 2016
Accepted: 09 December 2016
Published: 23 January 2017

Patricia P. Garcez^{2,3}, Juliana Minardi Nascimento^{1,1}, Janaina Mota de Vasconcelos³, Rodrigo Madeiro da Costa³, Rodrigo Delvecchio³, Pablo Trindade¹, Erick Correia Loliola¹, Luiza M. Higa¹, Juliana S. Cassoli¹, Gabriela Vitória¹, Patricia C. Sequeira¹, Jaroslaw Sochacki¹, Renato S. Aguiar¹, Hellen Thaís Fuzzi¹, Ana M. Bispo de Filippis¹, João Lúcio da Silva Gonçalves Vianez Júnior¹, Amilcar Tanuri¹, Daniel Martins-de-Souza¹ & Stevens K. Rehen^{1,2}

D'Aiuto et al. *Stem Cell Research & Therapy* (2018) 9:134
<https://doi.org/10.1186/s13287-018-0881-6>

Stem Cell Research & Therapy

RESEARCH

Open Access



Generation of three-dimensional human neuronal cultures: application to modeling CNS viral infections

Leonardo D'Aiuto¹, Jennifer Naciri¹, Nicholas Radio², Sessa Tekur², Dennis Clayton³, Gerard Apodaca³, Roberto Di Maio⁴, Yun Zhi⁵, Peter Dimitriou², Paolo Piazza⁶, Matthew Demers¹, Joel Wood¹, Charleen Chu^{4,9,12}, Jason Callio¹², Lora McClain¹³, Robert Yolken¹, James McNulty², Paul Kinchington², David Bloom¹⁰ and Vishwajit Nimgaonkar^{1,11}

Neuroinvasive potential of SARS-CoV-2 revealed in a human brain organoid model

Eric Song, Ca Zhang, Benjamin Israelow, Peiwen Liu, Orr-El Weizman, Felmet Liu, Yile Dai, Clara Szegedi-Buck, Yuki Yasumoto, Guilin Wang, Christopher Castaldi, Jaime Helkio, Evelyn Ng, John Wheeler, Mia Misdal Albjørn, Benjamin Fontes, Neal G. Ravindra, David Van Dijk, Sivakani Mane, Murat Gund, Aaron Ring, Craig B Wilton, Tomas L. Horvath, Angeliki Louvi, Shelli F. Firdad, Kaya Bilguvar, Akiko Iwasaki
doi: <https://doi.org/10.1101/2020.06.25.169946>

ACS Chemical
Neuroscience

pubs.acs.org/chemneuro

Viewpoint

iPSCs-Derived Platform: A Feasible Tool for Probing the Neurotropism of SARS-CoV-2

Xiao-Yuan Mao* and Wei-Lin Jin*

THE
EMBO
JOURNAL

SARS-CoV-2 targets cortical neurons of 3D human brain organoids and shows neurodegeneration-like effects

Anand Raman, Lisa Müller, Philipp Niklas Ostermann, Elie Gabriel, Pransy Abida-Islam, Andreas Müller-Schiffmann, Anujodhi Marappan, Olivier Goureau, Henning Gruell, Andreas Walker, Marcel André, Sandra Huska, Torsten Houwaart, Alexander Dilthey, Kai Wohlgemuth, Heymut Omran, Florian Klein, Dagma-Wieczorek, Orewin Adams, Jörg Timm, Carsten Koroh, Heiner Schaal, Jay Gopalakrishnan

Modeling HIV-1 neuropathogenesis using three-dimensional human brain organoids (hBORGs) with HIV-1 infected microglia

Roberta S. dos Reis¹, Shilpa Sant^{2,3,10}, Hannah Keeney¹, Marc C. E. Wagner¹ & Velupandi Ayyavoo^{1,11}

SCIENTIFIC
REPORTS
nature research



bioRxiv

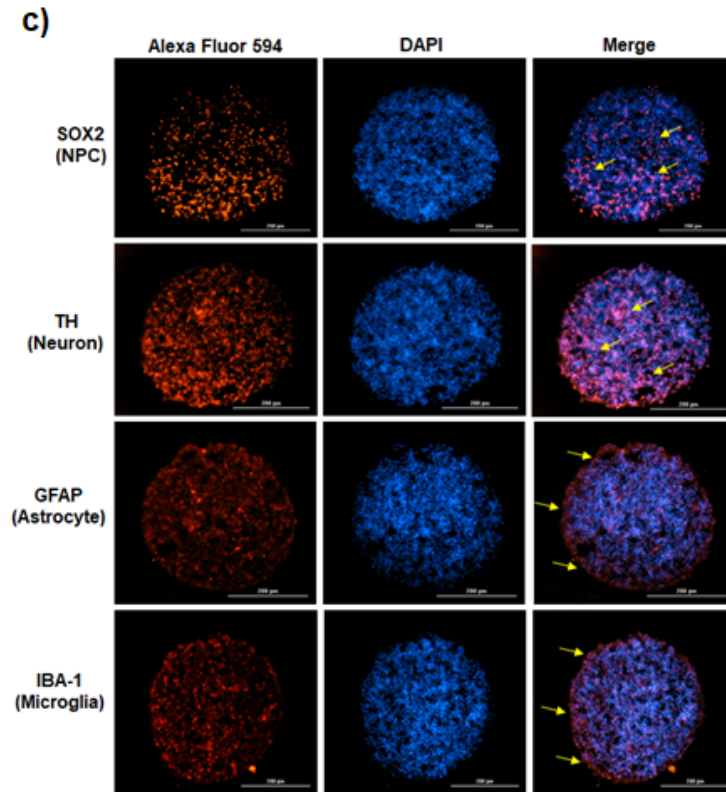
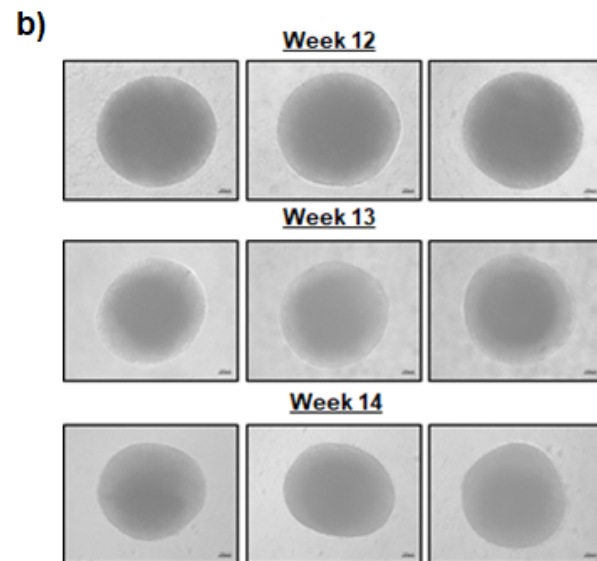
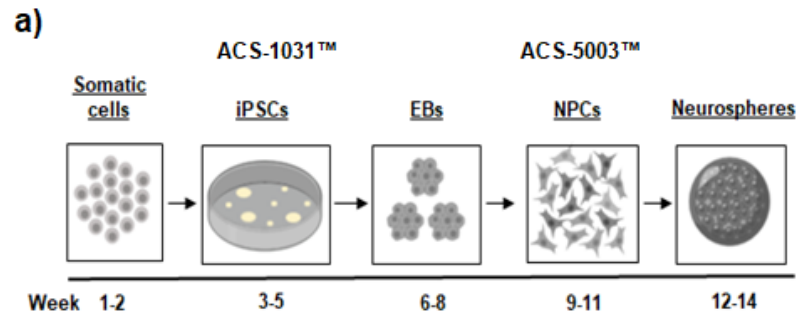
THE PREPRINT SERVER FOR BIOLOGY

Retroviral Infection of Human Neurospheres and Use of Stem Cell EVs to Repair Cellular Damage

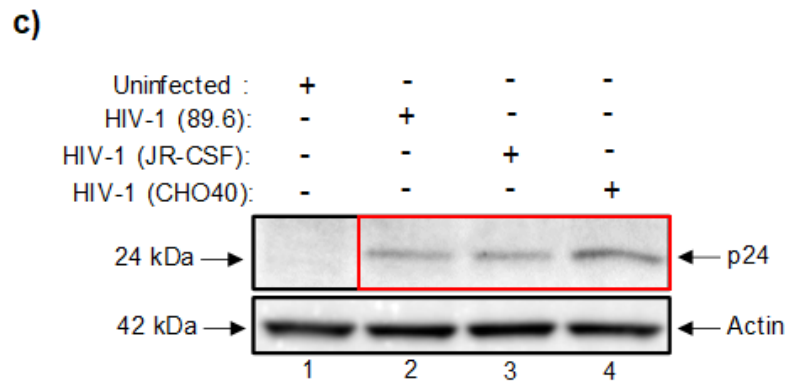
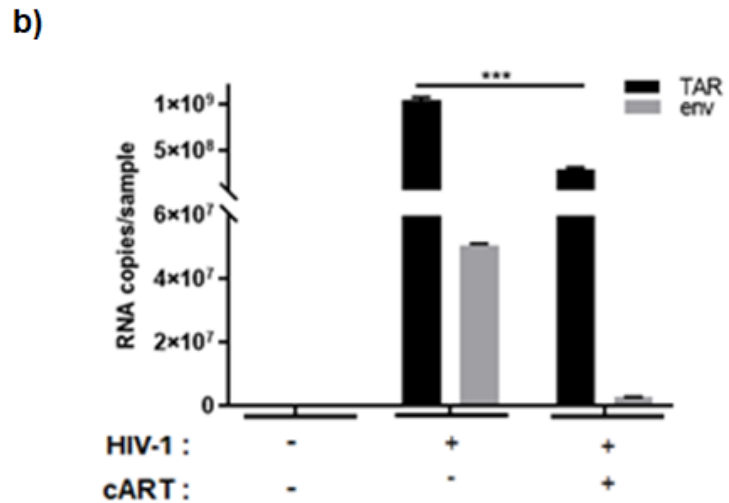
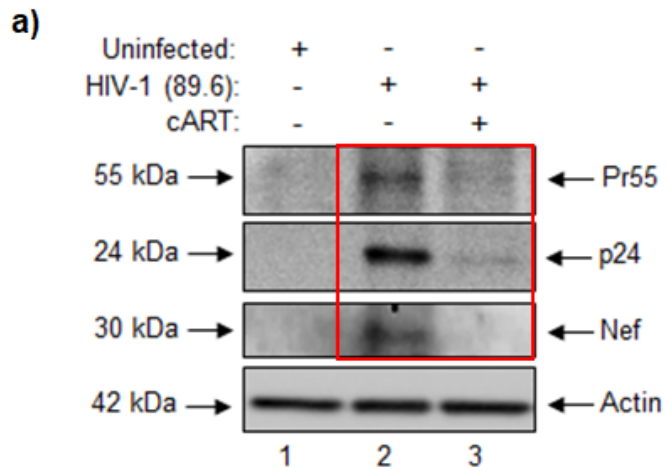
Heather Branscome, Dezhong Yin, Sheela Jacob, Maria Cowen, Yuriy Kim, Pooja Khatkar, James Erickson, Nazira El-Hage, Lance A. Liotta, Fatah Kashanchi

doi: <https://doi.org/10.1101/2020.12.31.424849>

Generation of neurospheres

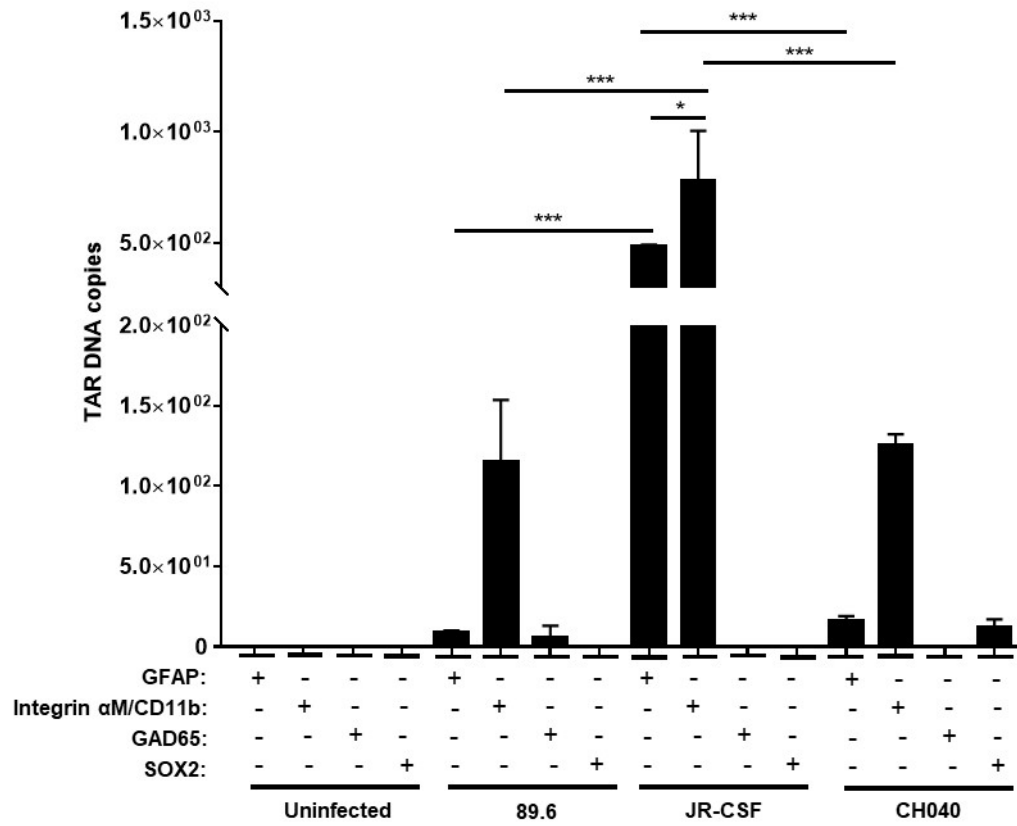


Infection of neurospheres

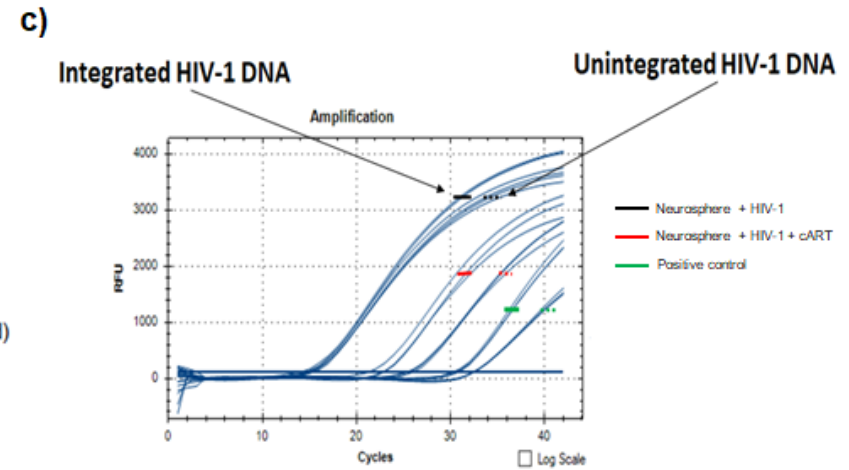
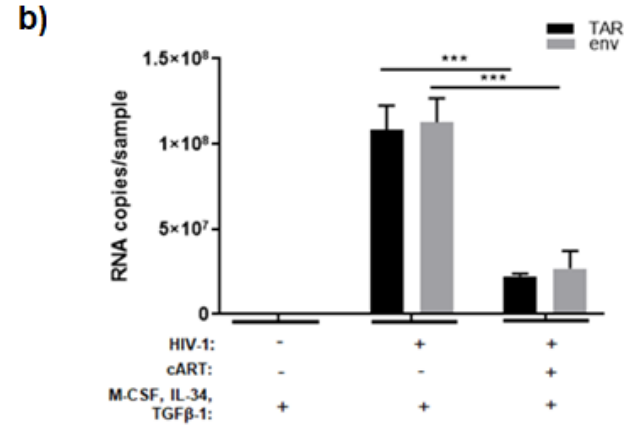
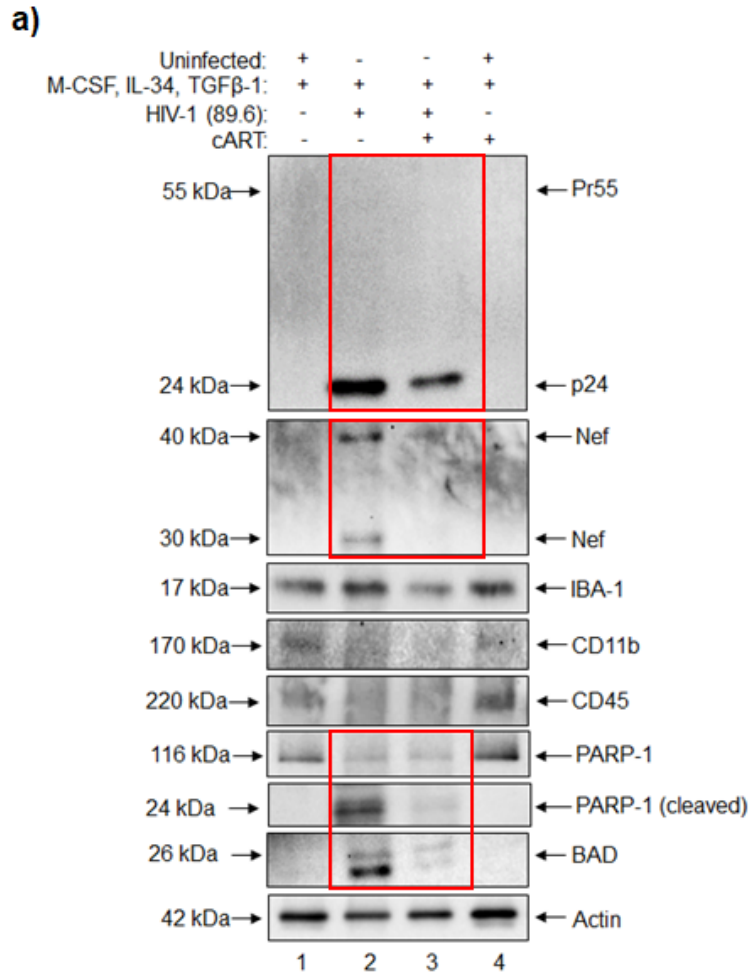


HIV-1 tropism

d)

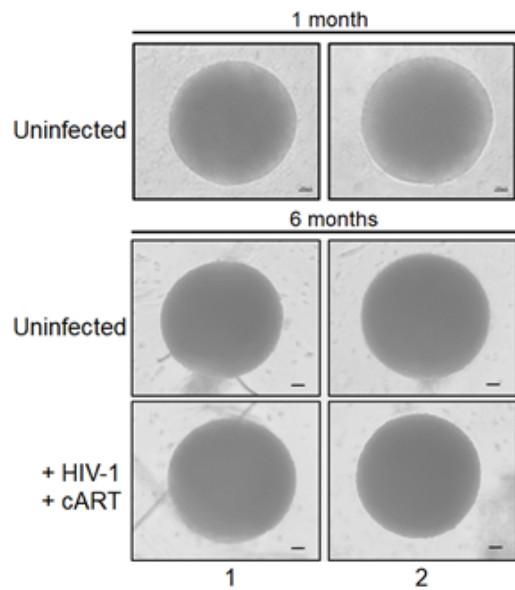


Infection of neurospheres (+ cytokines)

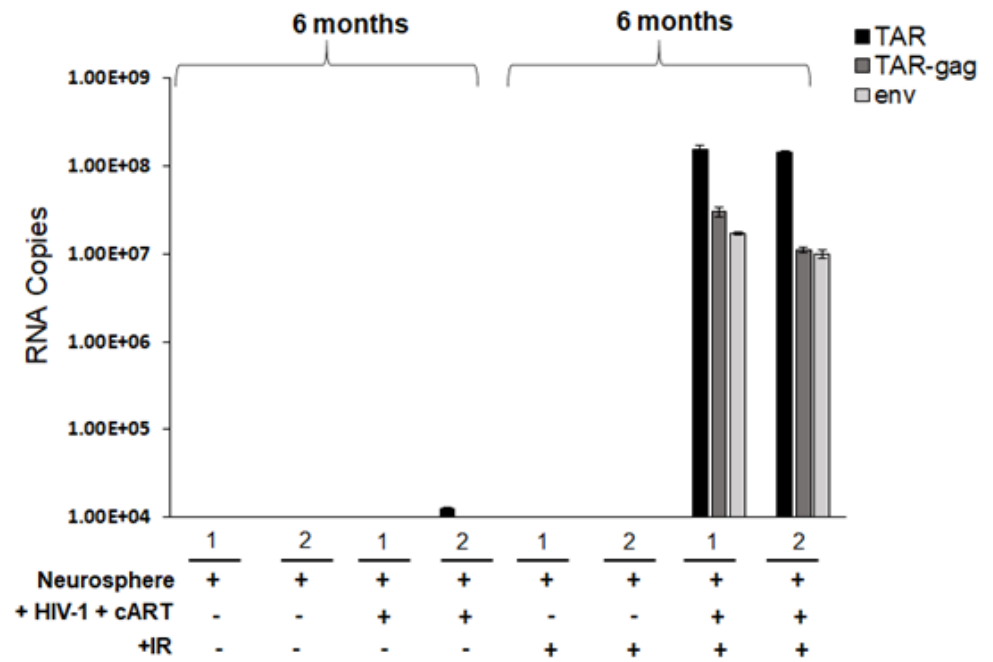


Infection of neurospheres (long-term)

a)

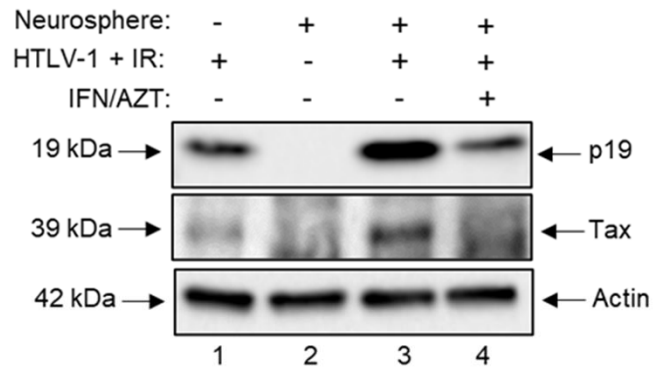


b)

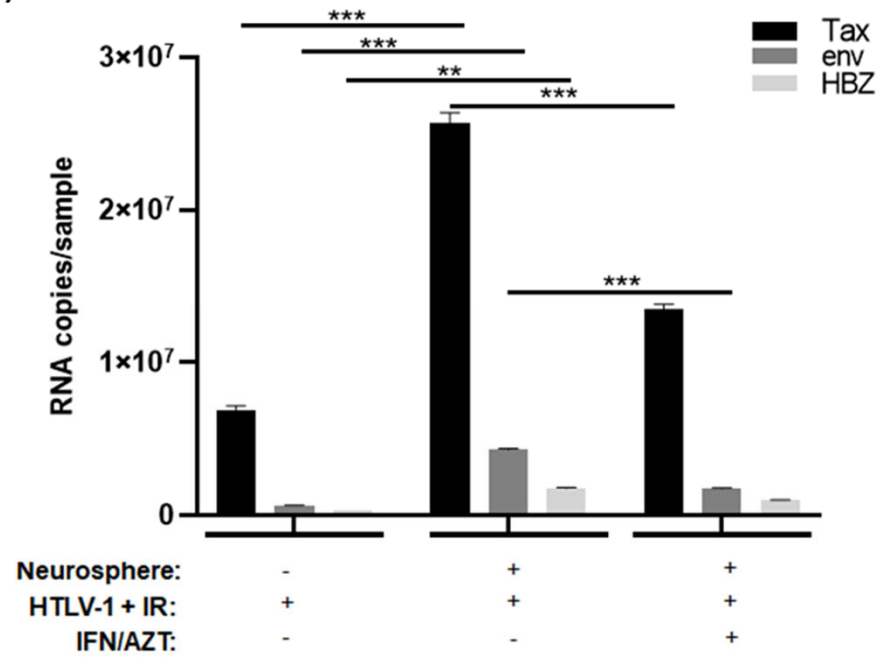


Infection of neurospheres (HTLV-1)

a)

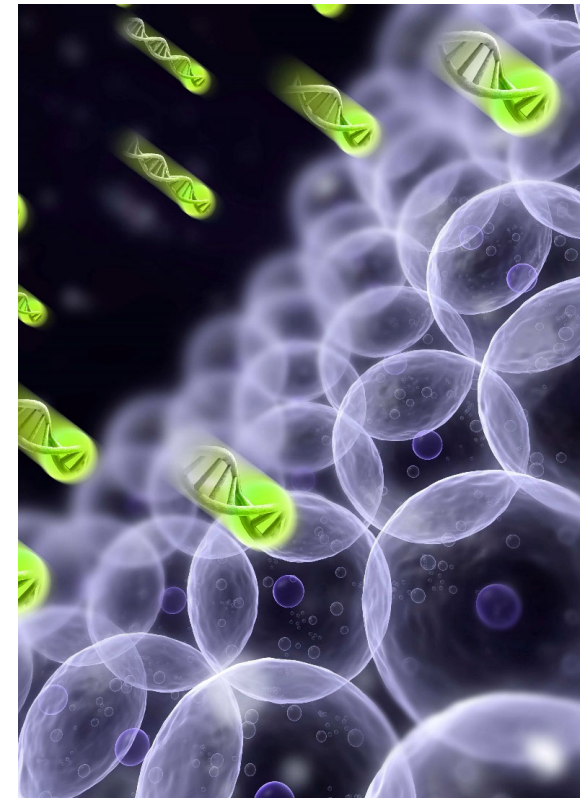


b)



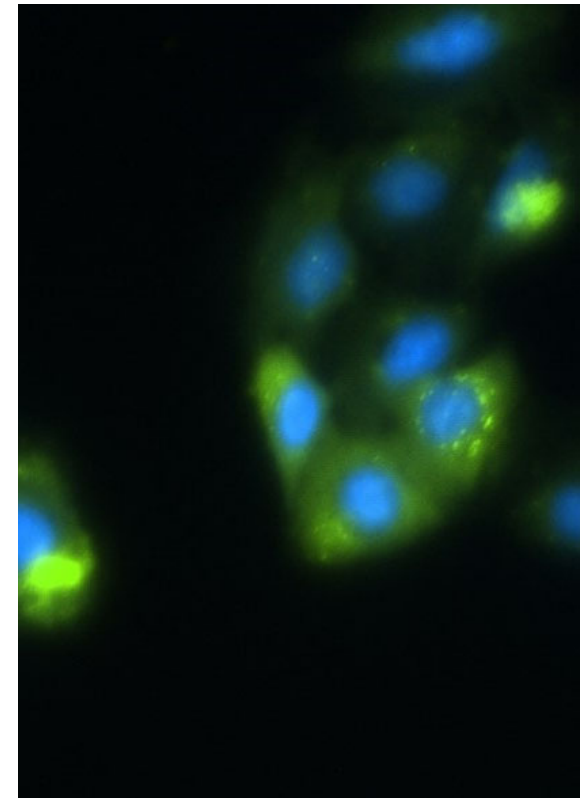
Stem cell EVs

- Contain various biological cargo (miRNAs, lncRNAs, proteins, cytokines) that can be transferred to recipient cells
- **Proposed to play a role in homeostasis through tissue repair, regeneration, and immunomodulation**
- Potential alternative to stem cell therapy due to higher potency, increased stability/shelf life, and lower immunogenicity
- Widely studied for reparative purposes (e.g., skin/wound healing, cardiac repair, **CNS-related pathologies**)



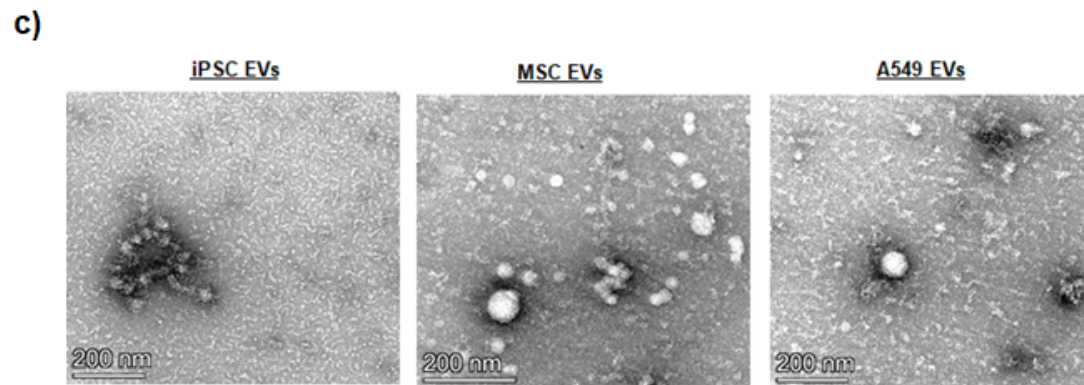
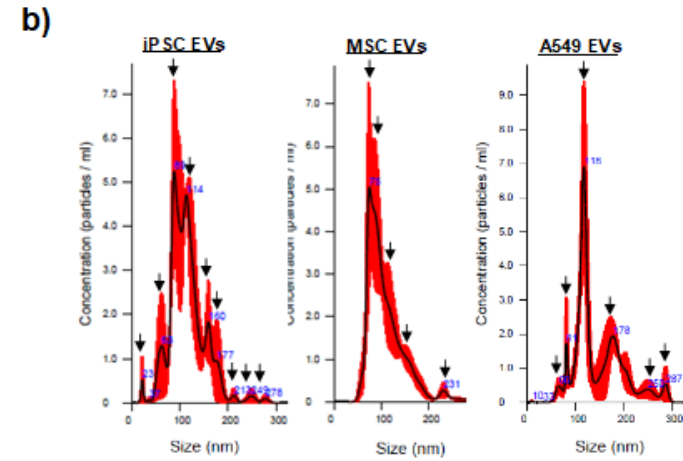
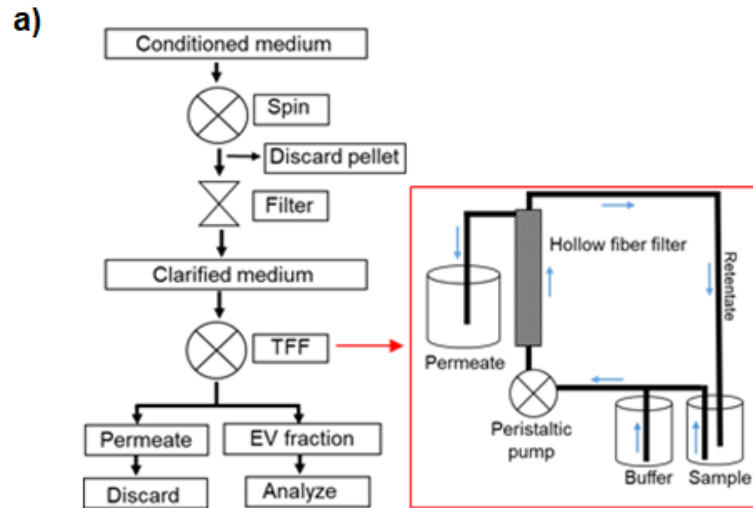
EVs and donor cells

- Mesenchymal Stem Cells (MSCs; ATCC® PCS-500-012™)
 - Human, normal
 - Bone-marrow derived
 - Authenticated for characteristic surface marker expression (CD90, CD73, CD105 positive; CD14, CD34, CD45 negative)
 - Multi-lineage differentiation potential (adipocyte, chondrocyte, osteocyte)
 - Induced Pluripotent Stem Cells (iPSCs; ATCC® ACS-1019™)
- Human, normal
 - Foreskin fibroblast-derived
 - Sendai virus reprogrammed
 - Authenticated for expression of stem cell markers (TRA-1-60, SSEA-4 positive; SSEA-1 negative)
 - Evaluated for pluripotency
- A549 Lung Carcinoma (ATCC® CCL-185™)
 - Control used for large-scale manufacturing and isolation
 - Equivalent to CCL-185-EXM™ exosomes (also: ATCC® SCRC-4000-EXM™, CRL-1435-EXM™, CCL-247-EXM™)



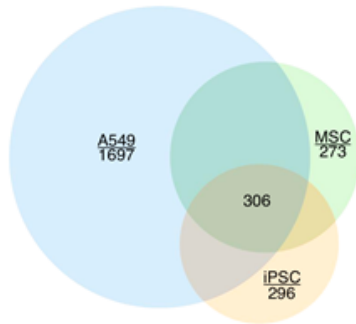
EVs synthesized in A549 cells

EV isolation and characterization

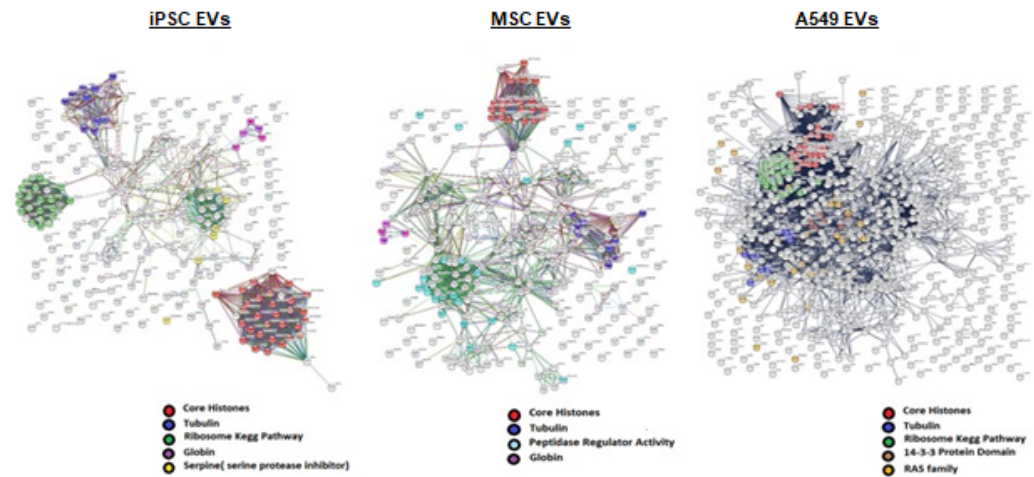


EV characterization

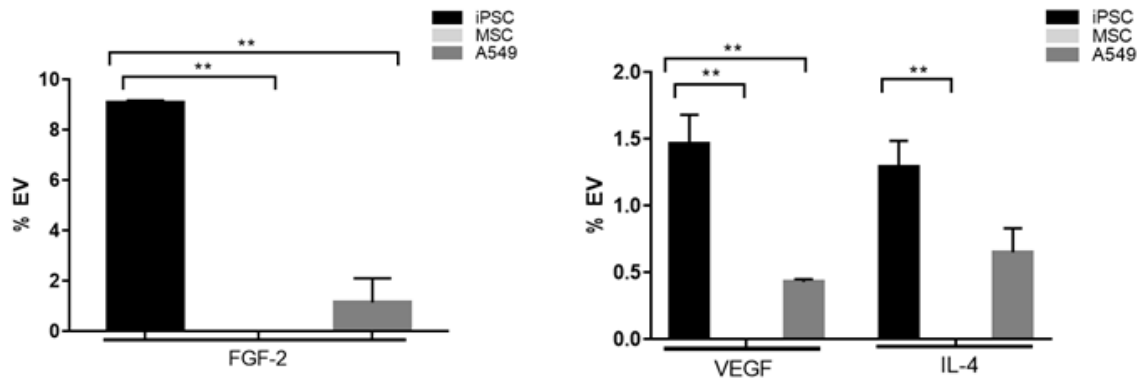
a)



b)

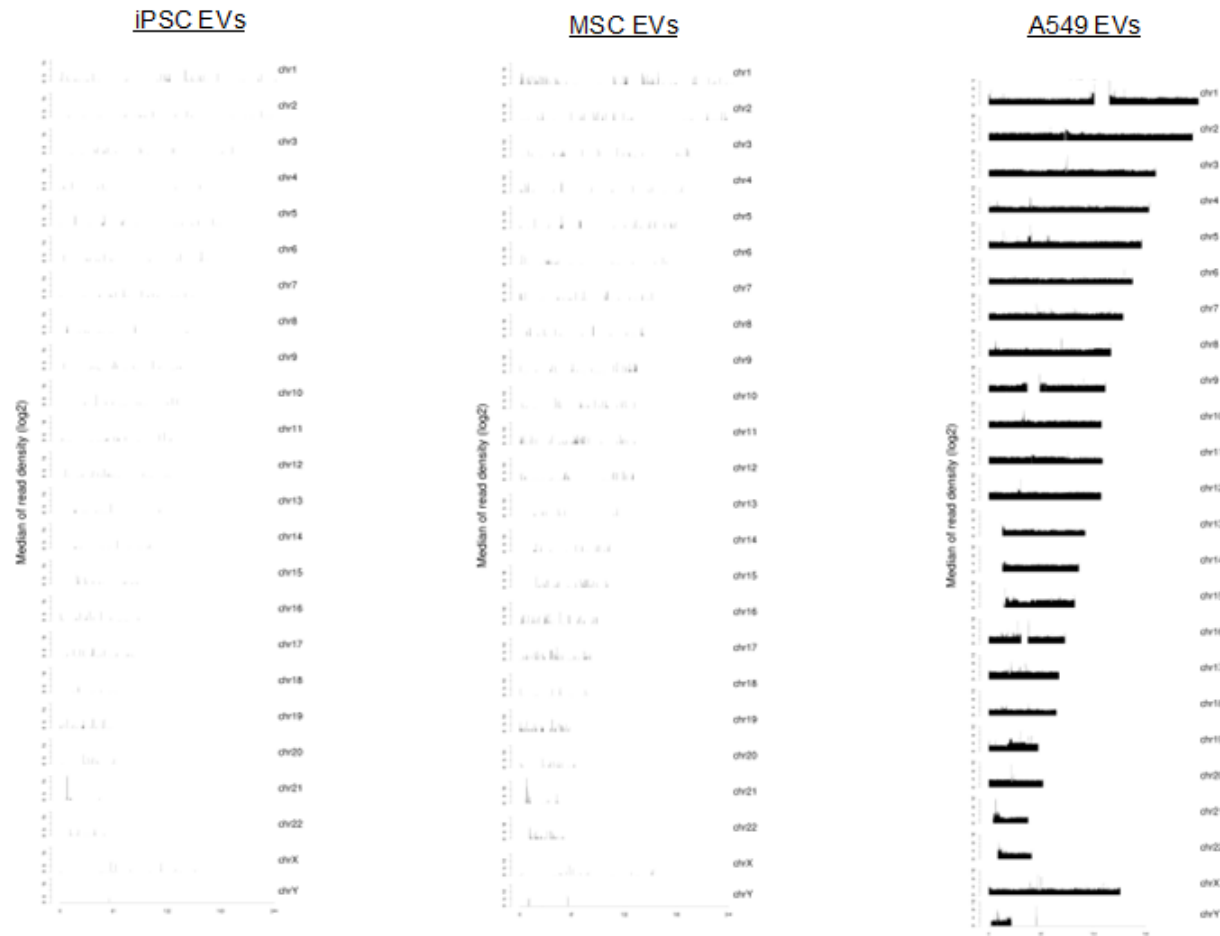


c)

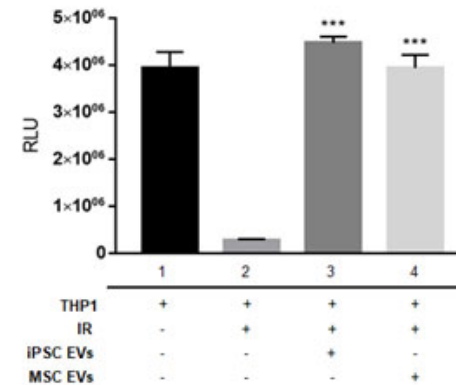
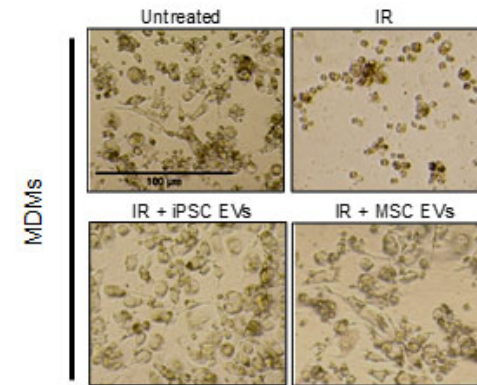
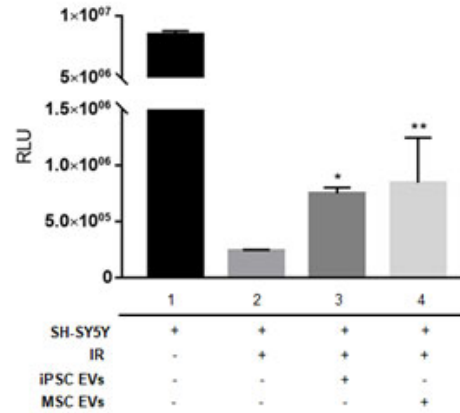
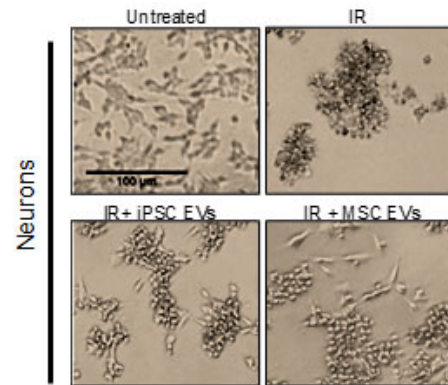
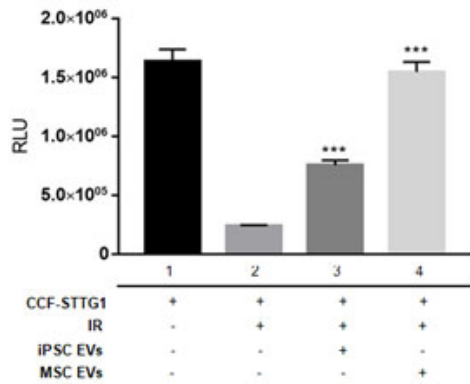
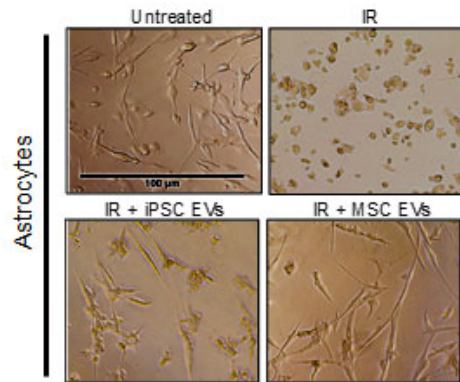


EV characterization

d)



EV function

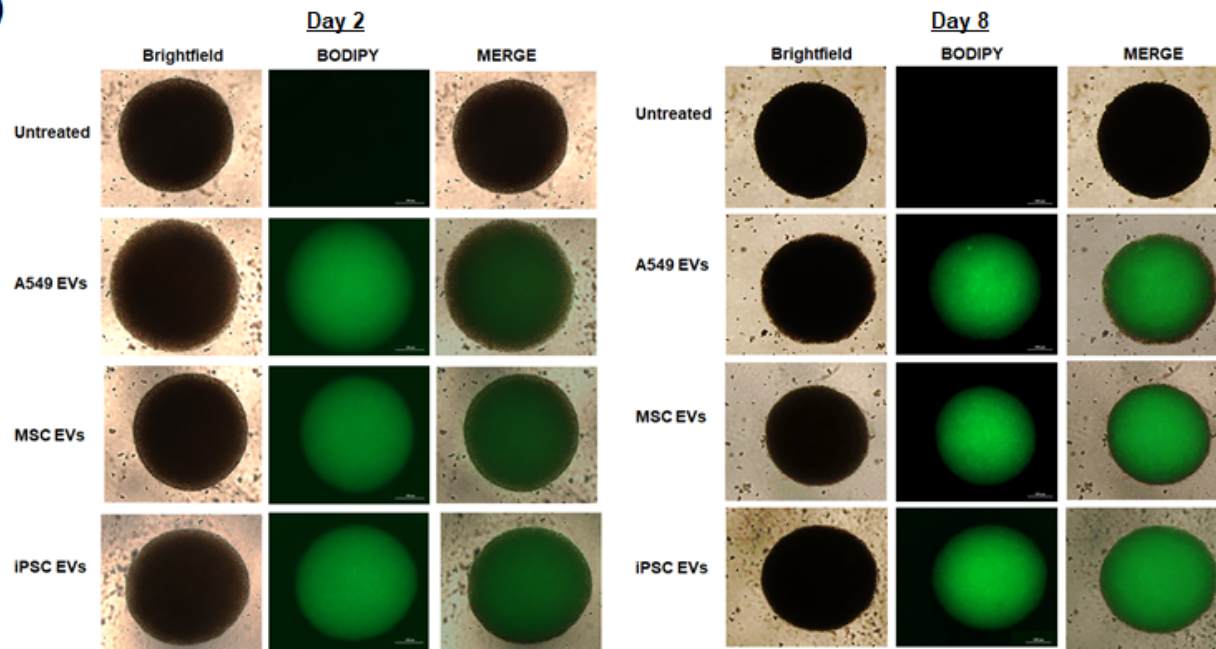


Stem Cell Extracellular Vesicles and Their Potential to Contribute to the Repair of Damaged CNS Cells

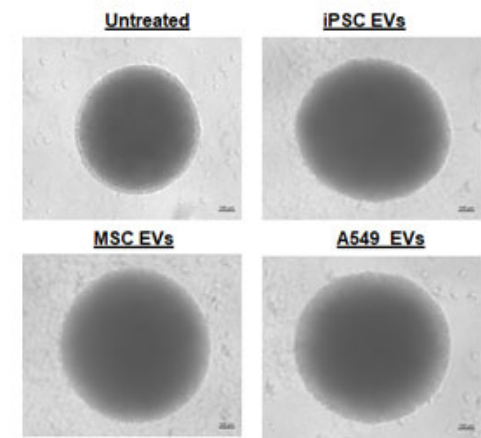
Heather Branscome^{1,2}, Siddhartha Paul¹, Pooja Khatkar¹, Yurly Kim¹, Robert A. Barclay¹, Daniel O. Pinto¹, Dezhong Yin¹, Weidong Zhou¹, Lance A. Liotta², Nazira El-Hage¹, Fatah Kashanchi^{1,7}

EV function

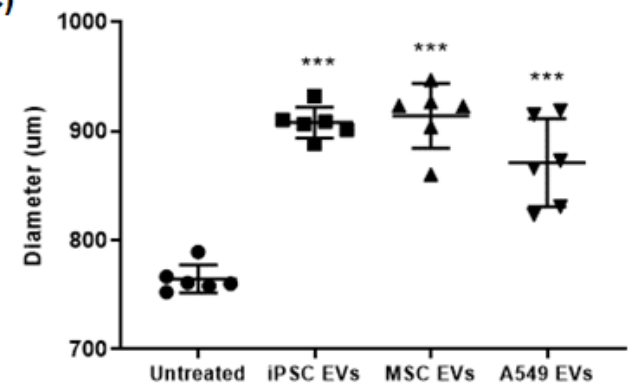
a)



b)

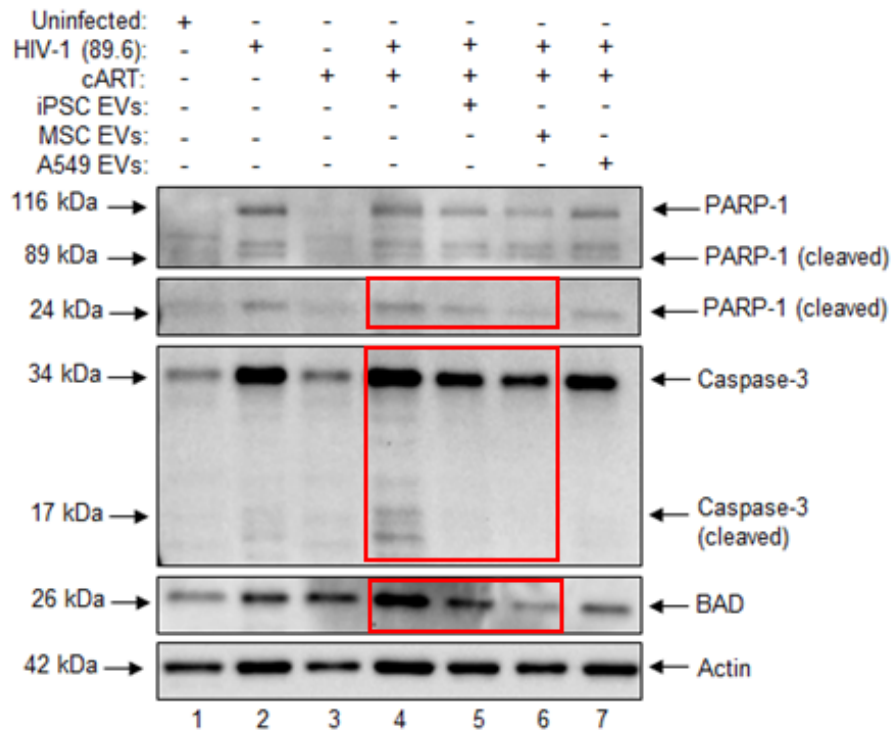


c)

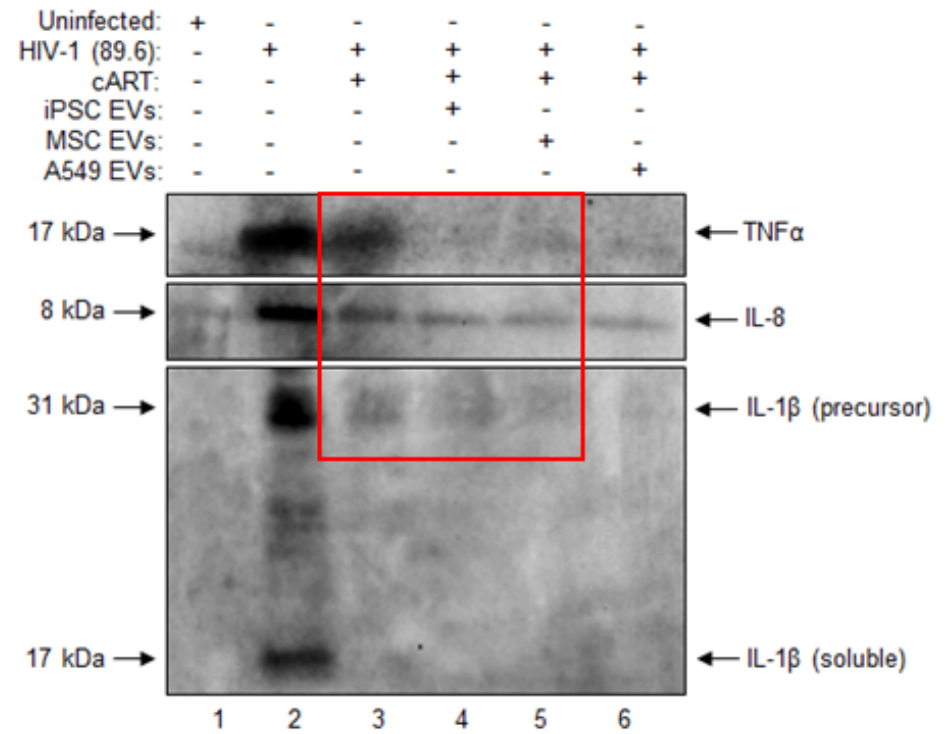


EV function

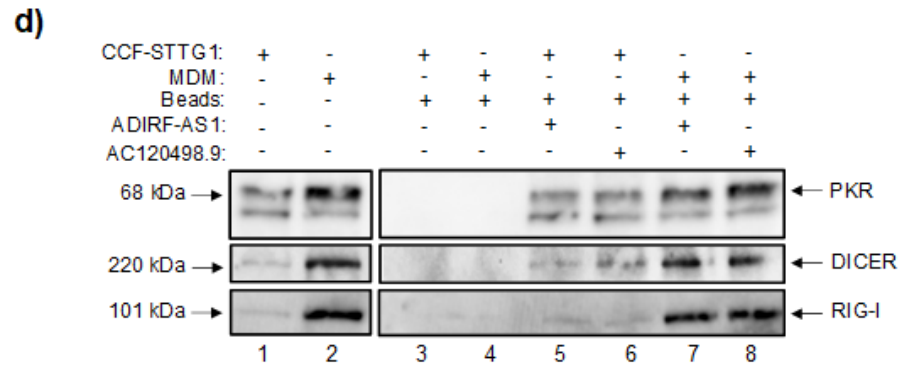
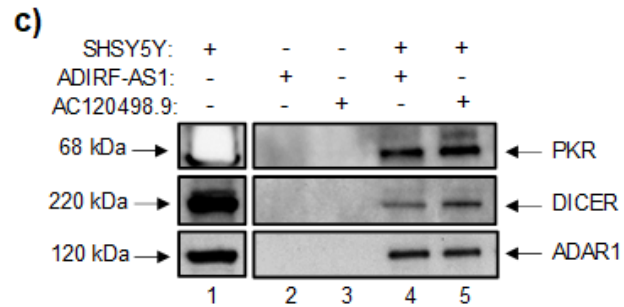
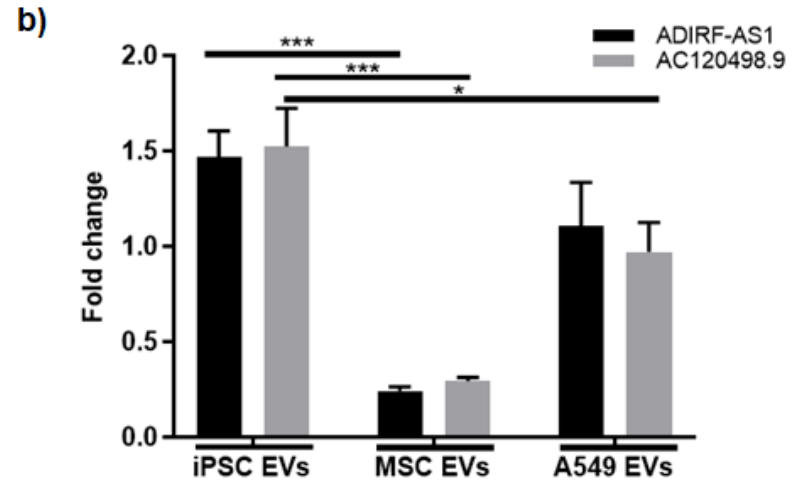
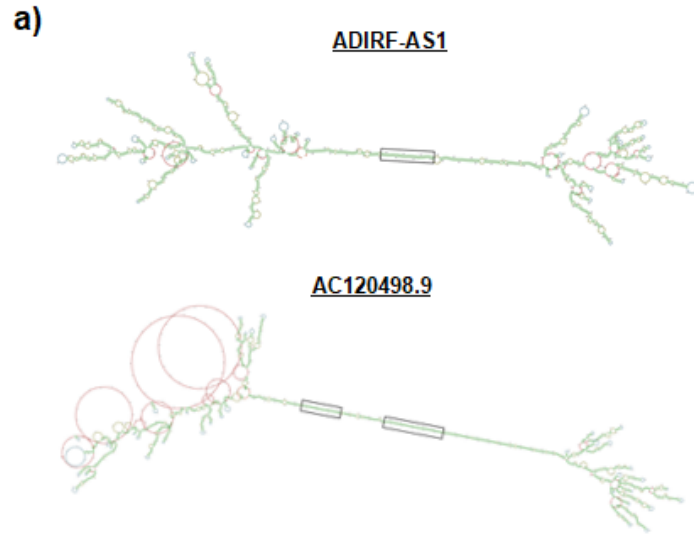
a)



b)

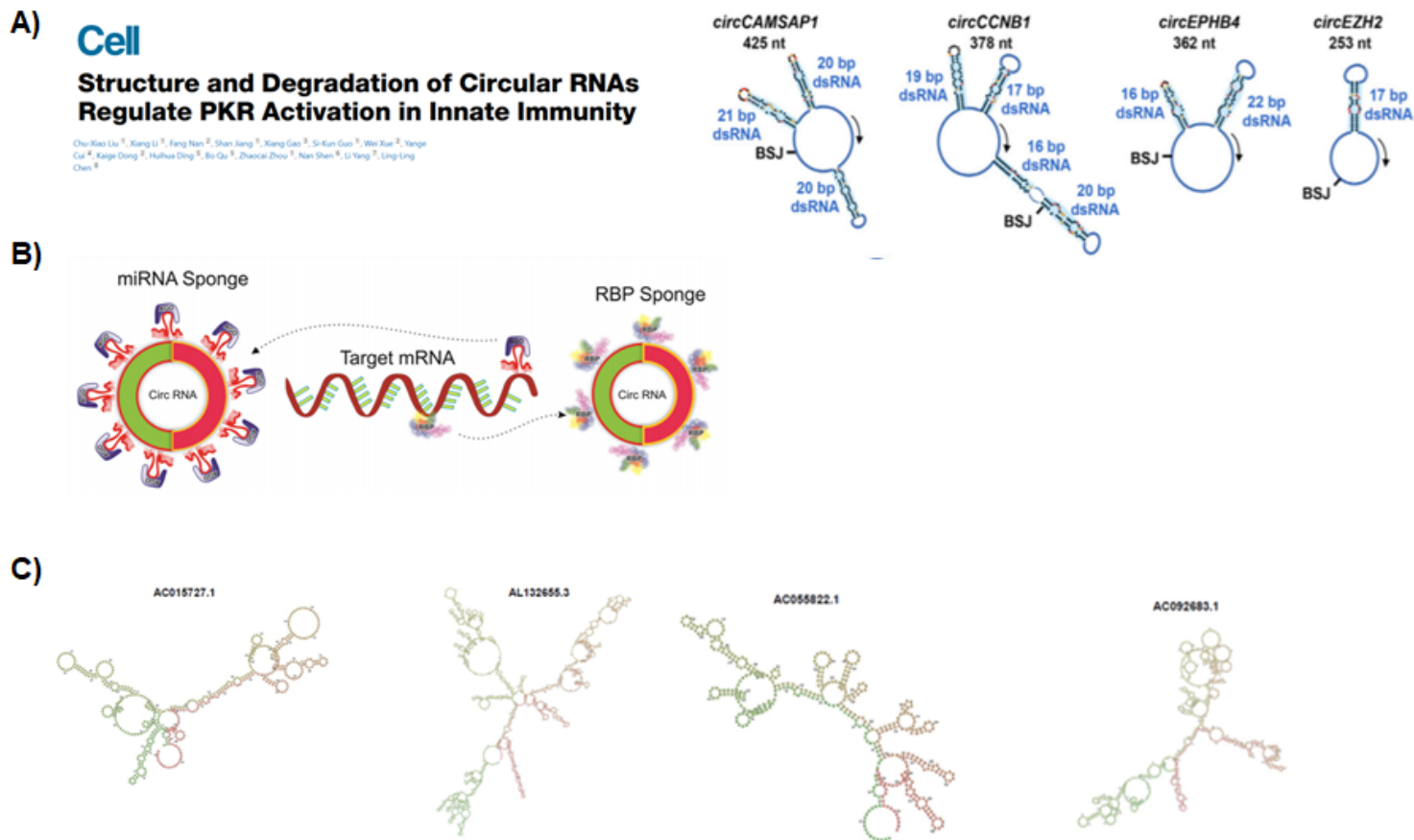


EV function



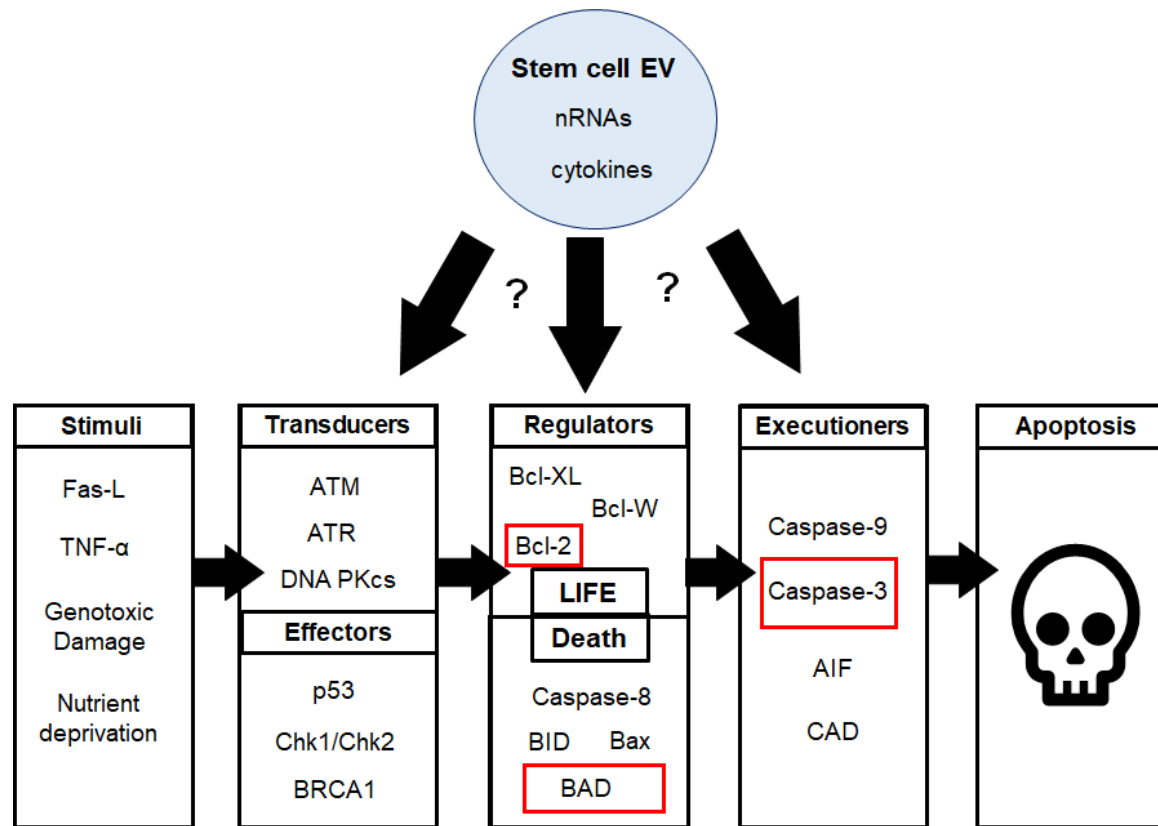
Future directions

- Further assessment of potential functional effects of EV-associated lncRNAs



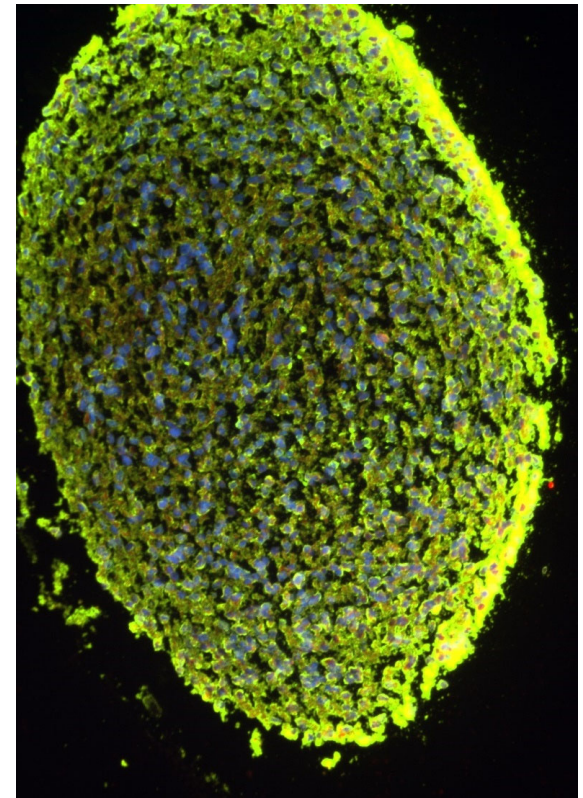
Future directions

- Better definitions of death and mechanisms of repair



Overview

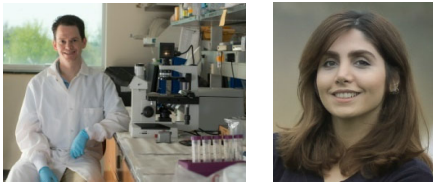
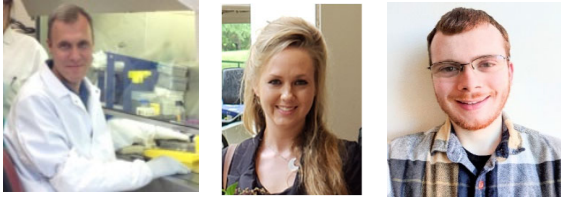
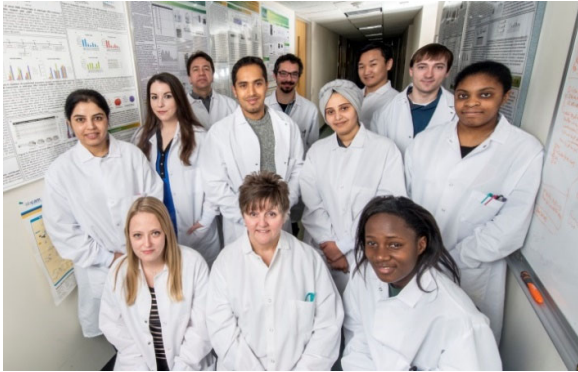
- Update on the most recent literature surrounding EVs from virally infected cells (“Damaging EVs”)
- Generation and infection of iPSC-derived neurospheres
- Effect of stem cell EVs (“Reparative EVs”) on HIV-1 infected neurospheres
- Summary



Summary

- HIV-1 TAR as non-coding RNA in EVs can be found in cell lines and patient samples.
- EVs are released first prior to viral release to potentially prime the neighboring uninfected cells.
- Early EV samples (i.e., 6 hours) contain p24, Nef, gp120 proteins, non-coding RNA, or coding RNAs; however, they are not infectious virions.
- The majority of the particles released (i.e., 24 hours) from infected activated cells are not infectious virions.
- NPC-derived 3-D neurospheres can be generated in a reproducible manner and can differentiate mature CNS cell types.
- 3-D neurospheres are permissive to retroviral infection.
- High yields of stem cell EVs can be recovered using advanced filtration methods.
- Stem cell EVs can exert functional effects in CNS-related cells in both 2-D and 3-D cultures.

Acknowledgements



ATCC

- Dr. Siddhartha Paul
- Dr. Dezhong Yin
- Dr. Sheela Jacob
- Dr. Mindy Goldsborough
- Dr. James Kramer
- James Fantuzzo
- Brian Shapiro
- Steve Budd
- Caitlyn Cabral
- Dong Kim

GMU

- Pooja Khatkar
- Yuriy Kim
- Fatemeh Dehbandi
- James Erickson
- Zachary Cuba
- Anastasia Williams
- Sebastian Molnar
- Michelle Pleet
- Maria Cowen
- Alex Barclay
- Catherine DeMarino
- Daniel Pinto
- Gwen Cox
- Gifty Mensah
- Sergey Iordanskiy
- Lance Liotta
- Y. Wu
- Kamel Khalili
- Gavin Sampey
- Rafal Kaminski
- Archana Gupta
- L. Shultz
- D. Price
- F. Romerio
- O. Kutsch
- V. Planelles
- Lena Al-Harhi
- Kevin Morris



THANK YOU

Credible Leads to Incredible™

