

#### Advanced models of Parkinson's disease

Brian Shapiro, Ph.D.

Technical Writer, Cell Biology, ATCC







## About ATCC

- Founded in 1925, ATCC is a non-profit organization with headquarters in Manassas, VA
- World's premiere biological materials resource and standards development organization
- ATCC collaborates with, and supports, the scientific community with industry-standard biological products and innovative solutions
- Strong team of 400+ employees; over onethird with advanced degrees







#### Agenda

#### Neural Progenitor Cells (NPCs) and Media

- Background information
- Differentiation potential of ATCC NPCs
- Toxicological studies
- Summary





## The current status of neurobiology research

- Primary cells from animals (mouse and rat neurons)
  - Not predictive
  - Donor variation
- Continuous cell lines (originally isolated from tumors)
  - Not normal
  - Not predictive
- Induced pluripotent stem cells (iPSCs; commercial or self-made)
  - Time and labor intensive
  - Often not validated for neural development



NPC-derived neurons





## What is needed?

- Biologically relevant models
- A true disease model
- Validated neural functioning
- Predictive for screening applications



NPC-derived neurons



# Neural progenitor cells (NPCs) - Neuronal differentiation



### Neural Progenitor Cells (NPCs) from ATCC



#### NPC-derived astrocytes

Value:

- Human models with no donor variation
- Live imaging is possible
- Cells exhibit full differentiation spectrum
- Complete system of cells and media will be available
   Key benefits:
  - More biologically relevant results/more predictive system
  - Markers allow for easy endpoint readout
  - Can differentiate to neuronal and glial cells
  - Easy to use and saves time







## ATCC<sup>®</sup> NPC offerings

ATCC <sup>®</sup> No.	Designation
ACS-3003	NPC Growth Kit – <i>add to DMEM/F12</i>
ACS-3004	NPC Dopaminergic Differentiation Kit – <i>add to DMEM/F12</i>
ACS-5001	NPCs derived from ATCC-DYS0530 Parkinson's Disease (ACS-1013) New!
ACS-5003	NPCs derived from ATCC-BXS0117 (ACS-1031)
ACS-5004	NPCs derived from ATCC-BYS0112 (ACS-1026)
ACS-5005	Neural Progenitor Cells derived from XCL-1 DCX-GFP (for late neuron differentiation)
ACS-5006	Neural Progenitor Cells derived from XCL-1 GFAP-Nanoluc <sup>®</sup> -Halotag <sup>®</sup> (for astrocyte differentiation)
ACS-5007	Neural Progenitor Cells derived from XCL-1 MAP2-Nanoluc <sup>®</sup> -Halotag <sup>®</sup> (for early neuron differentiation)
ACS-2103	Screening Fee – For Profit

ATCC<sup>®</sup> ACS-1026 – iPSC derived from bone marrow CD34+ cell from Caucasian male ATCC<sup>®</sup> ACS-1031 – iPSC derived from bone marrow CD34+ cell from Asian female



Reporter lines from iPSC derived from cord blood CD34+ from a Caucasian male (XL-1 iPSCs from NIH)



## QC testing of ATCC<sup>®</sup> NPCs

- Post-thaw cell viability: >80%
- Post-thaw viable cell number: >1x10<sup>6</sup> cells/vial
- Longevity: >15 PDLs or 5 passages
- NPC marker expression: Nestin<sup>+</sup>, Pax-6<sup>+</sup>, and Tra-I-60<sup>-</sup>
- Differentiation potential:
  - Tuj1<sup>+</sup> early neurons
  - TH<sup>+</sup> dopaminergic neurons
- Identity: STR profile matching parental iPSC line
- Sterility, mycoplasma, and viral panel testing: None detected



NPC-derived oligodentrocytes





#### Agenda

#### **NPCs and Media**

- Background information
- Differentiation potential of ATCC NPCs
- Toxicological studies
- Summary





# ATCC normal NPCs express NPC markers but **not** iPSC markers





## ATCC Parkinson's disease NPCs express NPC markers but **not** iPSC markers







## Dopaminergic neuron differentiation of NPCs





TH/DAPI



## Dopaminergic neuron differentiation of Parkinson's disease NPCs





## Astrocyte and oligodendrocyte differentiation

#### Astrocyte differentiation







#### **Oligodendrocyte differentiation**



ACS-5003





ACS-5001



## Dopaminergic neuron differentiation of NPC reporter lines



MAP2- NanoLuc<sup>®</sup>-HaloTag<sup>®</sup> (ACS-5007) DCX-GFP (ACS-5005) GFAP-NanoLuc<sup>®</sup>-HaloTag<sup>®</sup> (ACS-5006)



# <sup>■</sup>Expression of the luciferase reporter during dopaminergic neuron or astrocyte differentiation

Luciferase secretion during dopaminergic neuron differentiation of NanoLuc<sup>®</sup>-HaloTag<sup>®</sup> NPCs





Luciferase secretion during astrocyte differentiation of GFAP-NanoLuc<sup>®</sup>- HaloTag<sup>®</sup> NPCs





# Expression of the GFP or HaloTag<sup>®</sup> reporter during dopaminergic neuron or astrocyte differentiation



ATCC

18

## Development of ATCC's NPC expansion and dopaminergic differentiation media

NPCs cultured in company A NPC expansion media (top row) or ATCC NPC Growth Kit (bottom row) for 3 passages prior to differentiation using ATCC's NPC Dopaminergic Differentiation Kit

ATCC <sup>®</sup> No.	Designation
ACS-3003	NPC Growth Kit
ACS-3004	NPC Dopaminergic Differentiation Kit





## Expression of genes associated with the differentiation of NPCs

- TaqMan<sup>®</sup> primers were used to identify the presence of other types of neurons during dopaminergic neuron differentiation using ATCC<sup>®</sup> ACS-3004<sup>™</sup> media
- Dopaminergic neurons: TH, Nurr1, VMAT2, AADC
- Glutamatergic neurons: GLS2, vGLUT1,vGLUT2
- Gabaergic neurons: GABA (GABRB3)
- Motor neurons: EN1, LIM3, and Hb9
- Cholinergic neurons: ChAT







#### Early and dopaminergic neuron gene expression

Upregulation of early and dopaminergic neuron genes in ACS-5001, ACS-5003, and ACS-5007 NPCs during dopaminergic neuron differentiation



NPC-derived dopaminergic neurons



#### Expression of early neuron gene MAP2





### Expression of early neuron gene Tuj1





### Expression of dopaminergic neuron gene TH





### Expression of dopaminergic neuron gene Nurr1







#### Expression of VMAT2





#### Expression of DAT





27

#### Expression of AADC







#### Glutamatergic and GABAergic gene expression

Upregulation of glutamatergic and GABAergic neuron genes in ACS-5001, ACS-5003, and ACS-5007 NPCs during dopaminergic neuron differentiation



NPC-derived neurons





#### Expression of GLS2







#### Expression of vGLUT1







#### Expression of vGLUT2







#### Expression of GABA







## Motor and cholinergic gene expression

Upregulation of neuron genes in ACS-5001, ACS-5003, and ACS-5007 NPCs during dopaminergic neuron differentiation:

- Motor
  - LIM3
  - Hb9
  - EN1
- Cholinergic
  - ChAT



NPC-derived dopaminergic neurons



#### Expression of LIM3





#### **Expression of Hb9**





#### Expression of EN1





37

#### Expression of ChAT





### **Protein expression**

Confirmation of protein expression in ACS-5007 NPCs during dopaminergic differentiation by ICC



NPC-derived neurons



# Confirmation of dopaminergic neuronal-specific protein expression during differentiation by ICC







#### Agenda

#### **Neural Progenitor Cells and Media**

- Background information
- Differentiation potential of ATCC NPCs
- Toxicological studies
- Summary



NPC-derived astrocytes





### Neurotoxicity studies

#### Neurotoxicity studies with ACS-5001 and ACS-5003 NPCs

- Resazurin viability
  - Reliablue<sup>™</sup> (ATCC<sup>®</sup> 30-1014<sup>™</sup>)



NPC-derived oligodendrocytes



## ATCC

## Neurotoxicity studies – undifferentiated NPCs



## Neurotoxicity studies – undifferentiated NPCs





### Neurotoxicity studies – ACS-5001 and ACS-5003

Toxin	ACS-5001	ACS-5003
Amiodarone	Toxic	Toxic
Chlorhexidine	Toxic	Toxic
Cisplatin	Resistant	Weakly toxic
Piperine	Resistant	Resistant
Vincristine	Toxic	Toxic
Hydroxyurea	Resistant	Weakly Toxic
Paclitaxel	Toxic	Toxic





#### Agenda

#### **Neural Progenitor Cells and Media**

- Background information
- Differentiation potential of ATCC NPCs
- Toxicological studies
- Summary



NPC-derived astrocytes





### NPCs – Summary

- Cells and media with easy to use protocols
  - Expansion and Differentiation Medium
- Human model with no donor variation
  - Ability to expand and bank
- Differentiation across a wide spectrum of neural and glial lineages
  - Neurons
  - Astrocytes
  - Oligodendrocytes
- Live imaging of differentiation
  - GFP expression upon neural differentiation



NPC-derived neurons



## NPCs – Summary

- Our studies demonstrated that ATCC normal and PD NPCs have the potential to be differentiated into:
  - Dopaminergic neurons
  - GABAergic neurons
  - Glutamatergic neurons
  - Motor neurons
  - Cholinergic neurons

after treatment of NPCs with ATCC dopaminergic differentiation media

ATCC NPCs are suitable for drug screening applications



NPC-derived astrocytes



#### Disclaimers

© 2018 American Type Culture Collection. The ATCC trademark and trade name, and any other trademarks listed in this publication are trademarks owned by the American Type Culture Collection unless indicated otherwise. NanoLuc<sup>®</sup>, HaloTag<sup>®</sup> and Promega are registered trademarks or trademarks of Promega Corporation. TaqMan<sup>®</sup> is a registered trademark of Roche Molecular Systems, Inc.



NPC-derived astrocytes





## Thank you for joining today!

#### Visit us at SOT 18, booth 1422, March 11-15

- Exhibitor-hosted Session: March 14, 2018
   Mindy Goldsborough, Chief Science and
   Technology Officer and General Manager, ATCC
   Kevin Grady, Product Line Business Manager, ATCC
   Chaozhong Zou, Senior Scientist, ATCC
   Advanced In Vitro Solute Carrier Transporter
   Models for Renal Toxicity Studies and Screening
- Poster Presentations: March 12, 2018
   Chaozhong Zou, Senior Scientist, ATCC
   Advanced In Vitro Solute Carrier Transporter
   Models for Renal Toxicity Studies and Screening

Sheela Jacob, Scientist, ATCC

Comprehensive Gene Expression Analysis and Neurotoxicity Testing of Human iPSCs-Derived Neural Progenitor Cells and Neurons



Please email additional questions to: tech@atcc.org

