

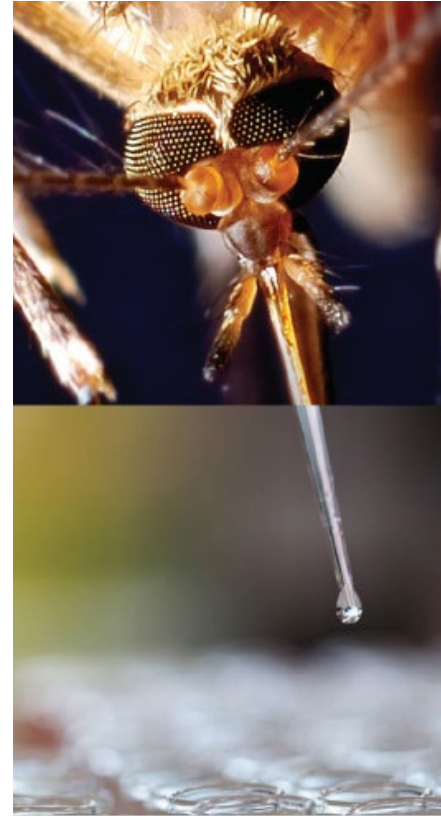
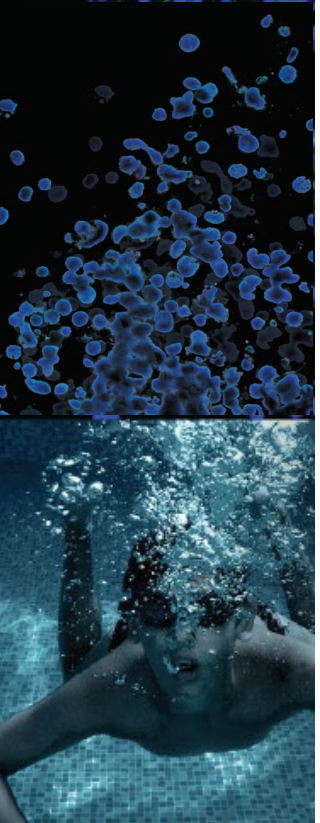


Does All Disease Begin in the Gut?

Monitoring the Barrier Function of an In Vitro Gut Mimic

Eline Geervliet, MSc
Application Scientist
Locsense B.V.

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 advanced models, differentiated stem cells, gene editing
- 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 550+ employees, over one-third with advanced degrees

Does all disease begin in the gut?

~Hippocrates

Monitoring the barrier function of an in
vitro gut mimic

Eline Geervliet, MSc
Application scientist at Locsense B.V.
www.Locsense.nl



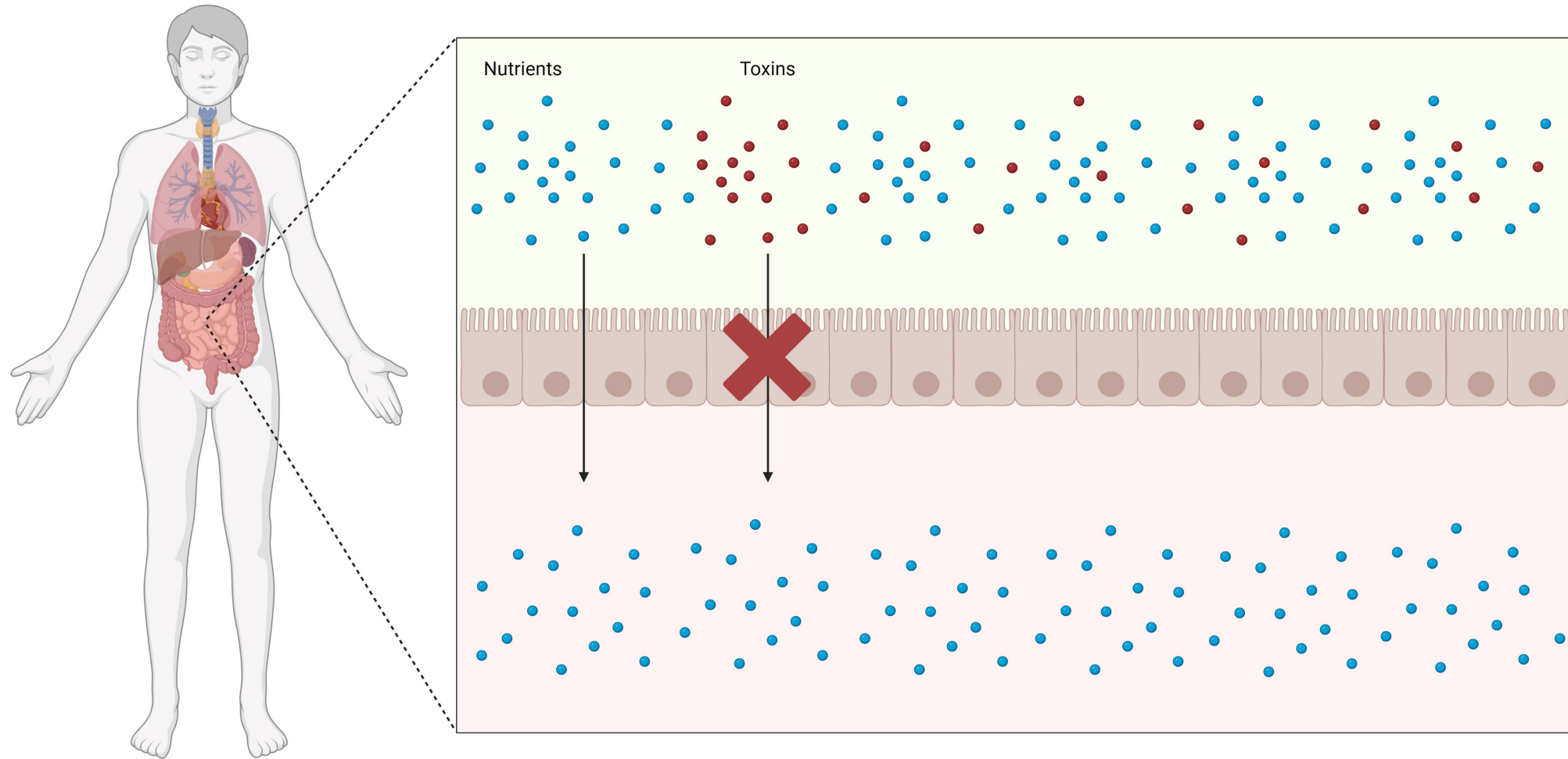
Index

- Function of the gut
- Gut mimic: *in vitro* translation
- Impedance spectroscopy & transepithelial electrical resistance (TEER)
- Interfering with the tight junction/barrier
- Measuring the tight junction/barrier
- Conclusions

Index

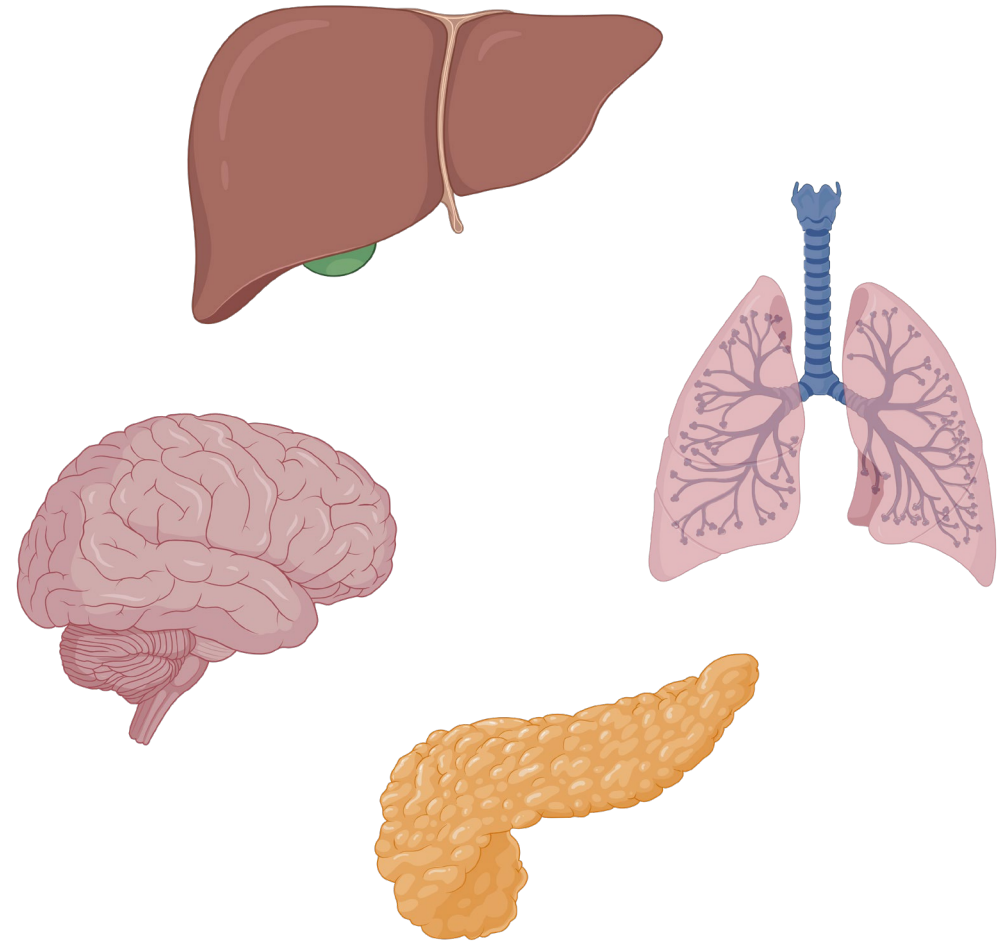
- Function of the gut
- Gut mimic: *in vitro* translation
- Impedance spectroscopy & transepithelial electrical resistance (TEER)
- Interfering with the tight junction/barrier
- Measuring the tight junction/barrier
- Conclusions

Function of the gut



Impaired gut barrier leads to:

- Liver disease [1]
- Brain damage [2]
- Lung disease [3]
- Pancreas [4]
- Etc.....



[1] A. Albillos et al. "The gut-liver axis in liver disease: Pathophysiological basis for therapy", *J. Hepatol* 2020

[2] J. Appleton et al. "The Gut-Brain Axis: Influence of Microbiota on Mood and Mental Health," *Integr Med* 2018

[3] L. Qu et al. "COPD and Gut-Lung Axis: How Microbiota and Host Inflammasome Influence COPD and Related Therapeutics," *Front Microbiol* 2022

[4] Z. Zhang, et al. "Intestinal homeostasis and inflammation: Gut microbiota at the crossroads of pancreas-intestinal barrier axis," *Eur J Immunol* 2022

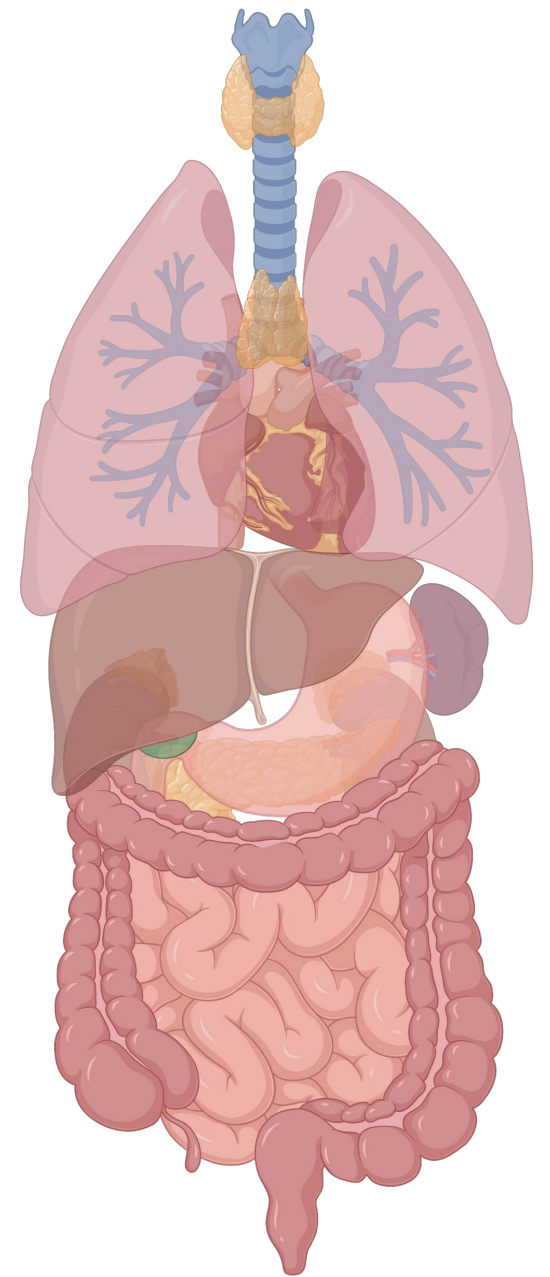
In vitro analysis

A representable *in vitro* gut mimic is necessary to investigate:

- which factors influence the barrier function
- the cross talk between the gut and other organs

Aim:

- Propose a representable gut mimic
- Evaluate the barrier function of the gut mimic

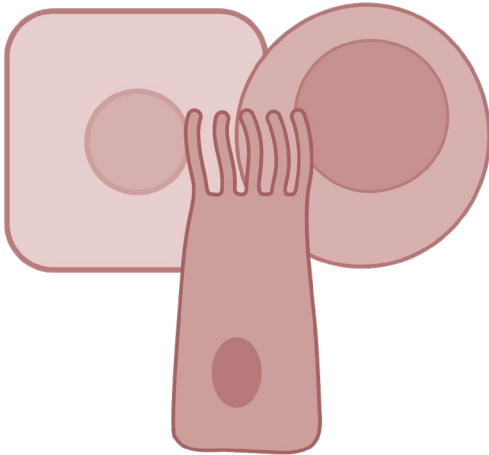


Index

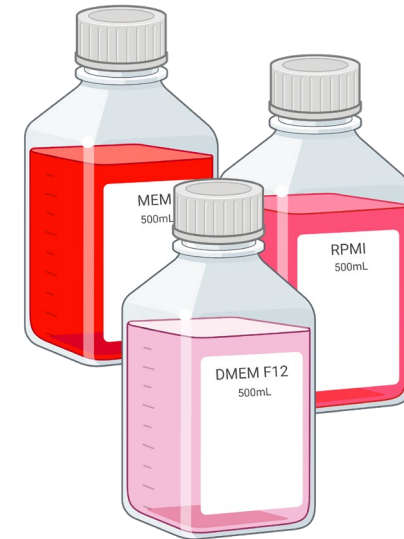
- Function of the gut
- Gut mimic: *in vitro* translation
- Impedance spectroscopy & transepithelial electrical resistance (TEER)
- Interfering with the tight junction/barrier
- Measuring the tight junction/barrier
- Conclusions

In vitro translation of the gut barrier

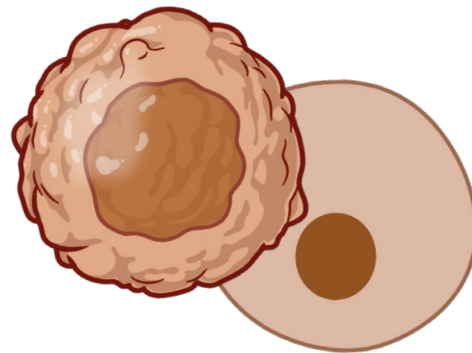
1. Cell type





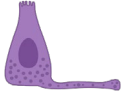




2. Cell medium

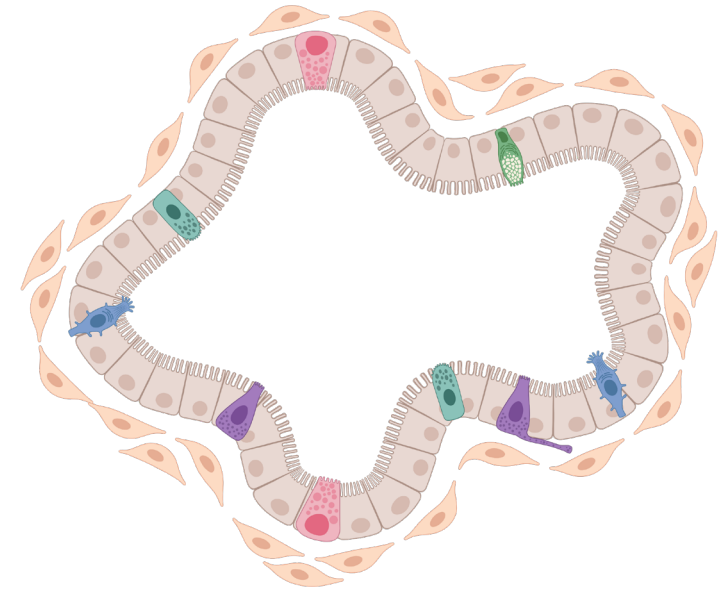


3. 2D/3D structure



Cell type [1, 2]

- Villi – differentiated compartment
 - Enterocytes – nutrient absorption 
 - Goblet cells – production of mucus 
 - Enteroendocrine cells – hormone production 
 - Cup cells – function not clear 
- Crypts – proliferative compartment
 - Paneth cells – regulate gut microbiota 
- Other
 - Microfold cells (M-cells) – contact with immune system 
 - Tuft cells – contact with immune system and taste perception 



[1] V. Bonis, et al. “The Intestinal Epithelium – Fluid Fate and Rigid Structure From Crypt Bottom to Villus Tip,” *Front. Cell Dev. Biol.* 2021

[2] F. Gerbe, et al. “The intestinal epithelium tuft cells: specification and function” *Cell Mol. Life Sci.* 2012

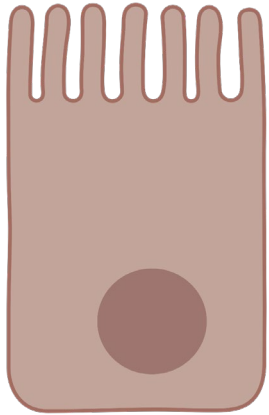
Cell type

Few - simple

Complex - many



Cell type



Caco-2 [1]
Enterocyte

Derived from the colon

ATCC[®] HTB-37[™]

20% FBS

90% cell number [1]



HT29-mtx [1]
Goblet-like cell

Derived from the colon

Mucus producing cells

10% FBS

10% cell number [1]

13[1] P. Hoffmann et al., "Caco-2/HT29-MTX co-cultured cells as a model for studying physiological properties and toxin-induced effects on intestinal cells," *PLoS One* 2021

Culturing medium

Cultured in DMEM-F12-GlutaMAX

Alternative DMEM-F12-Glutamine (ATCC[®] 30-2006[™])

1% Supplemented with non-essential amino acids

1% Pen/strep, removed during standard culture and monolayer formation

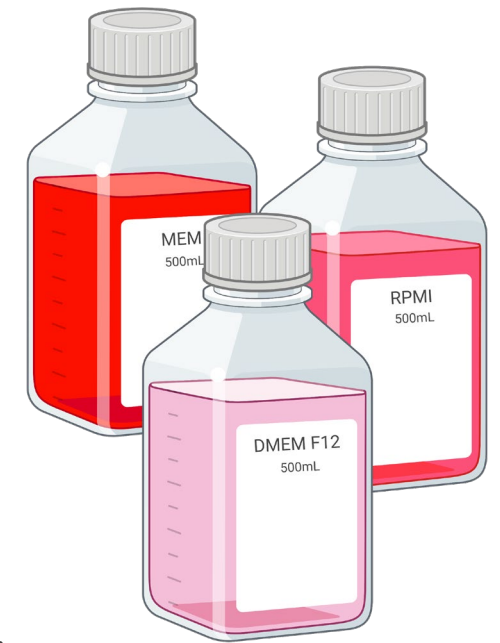
Penicillin-Streptomycin Solution (ATCC[®] 30-2300[™])

15% FBS, compromise

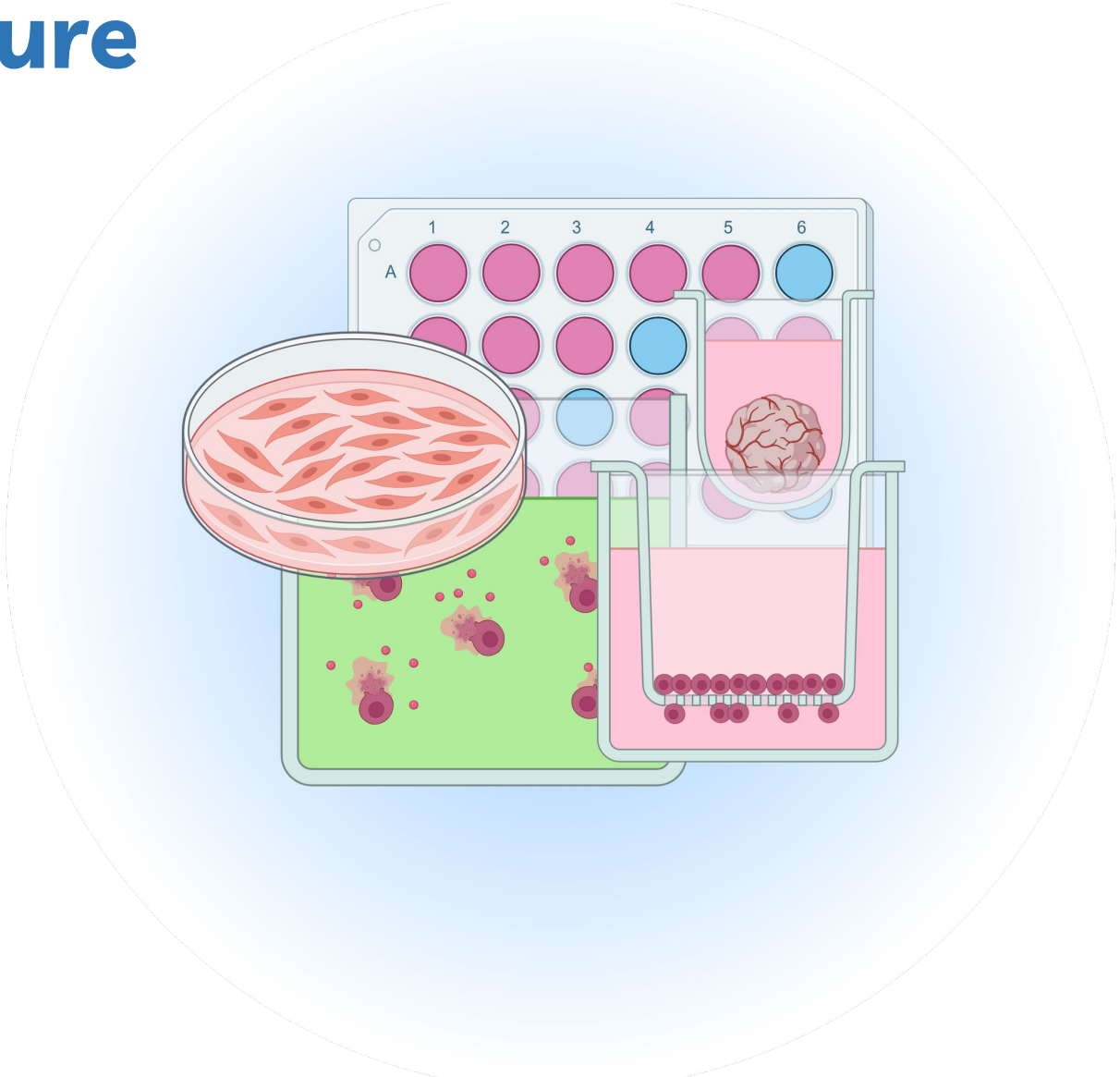
Fetal Bovine Serum (FBS) (ATCC[®] 30-2020[™])

Passaged via trypsin-EDTA

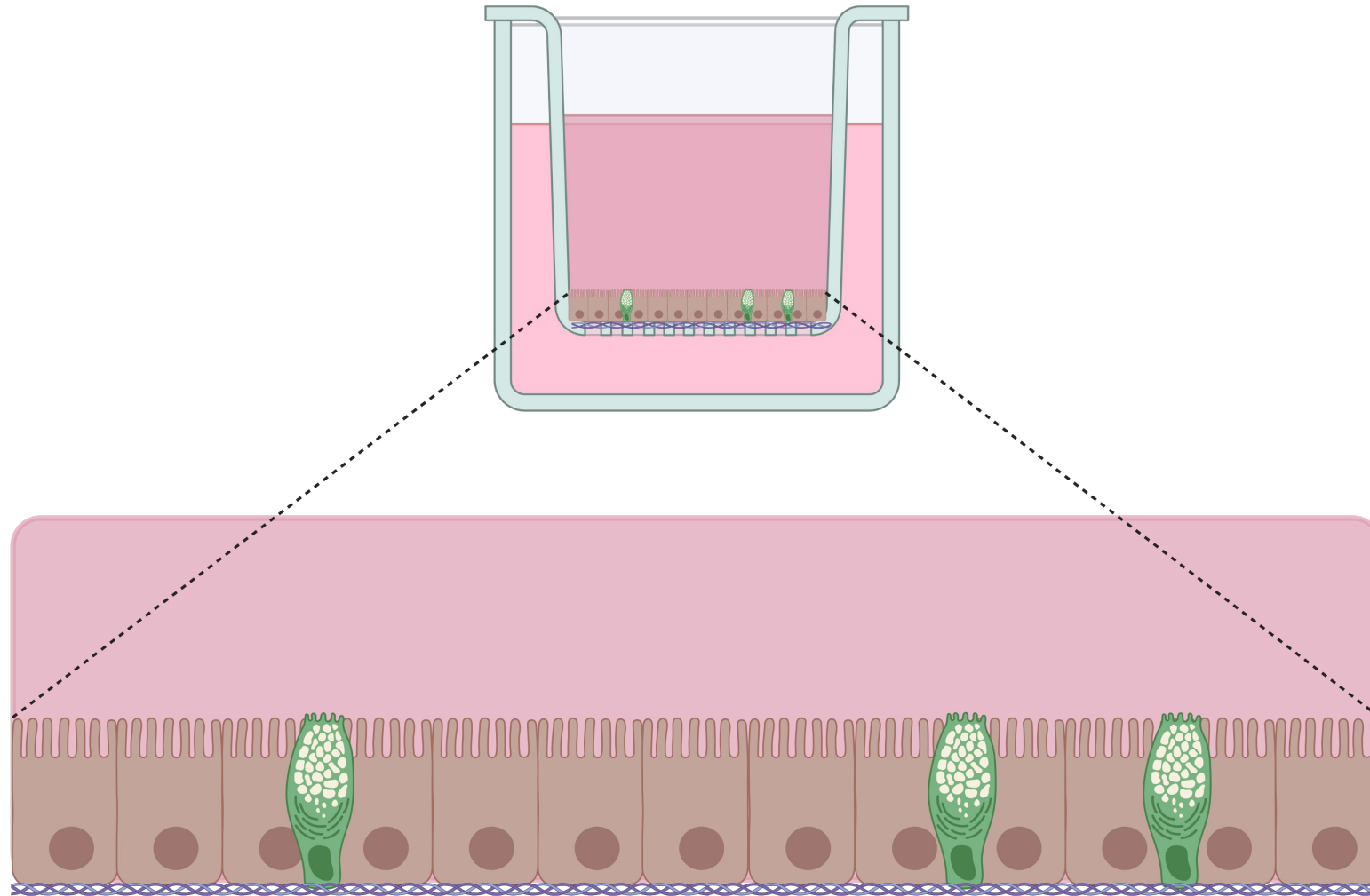
Trypsin-EDTA (ATCC[®] 30-2101[™])



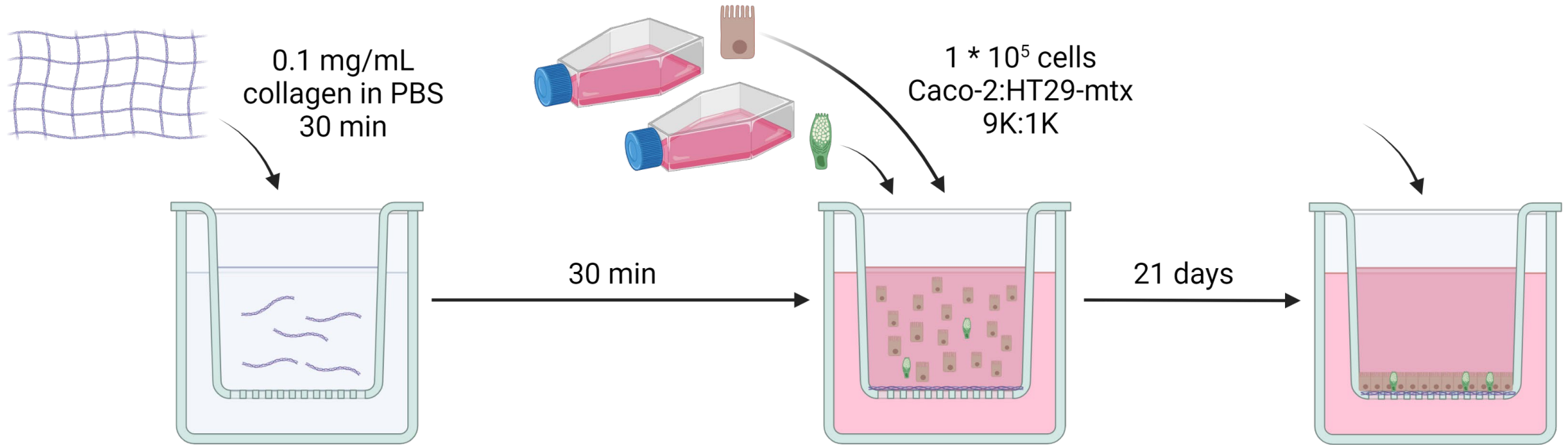
3D structure



Transwell



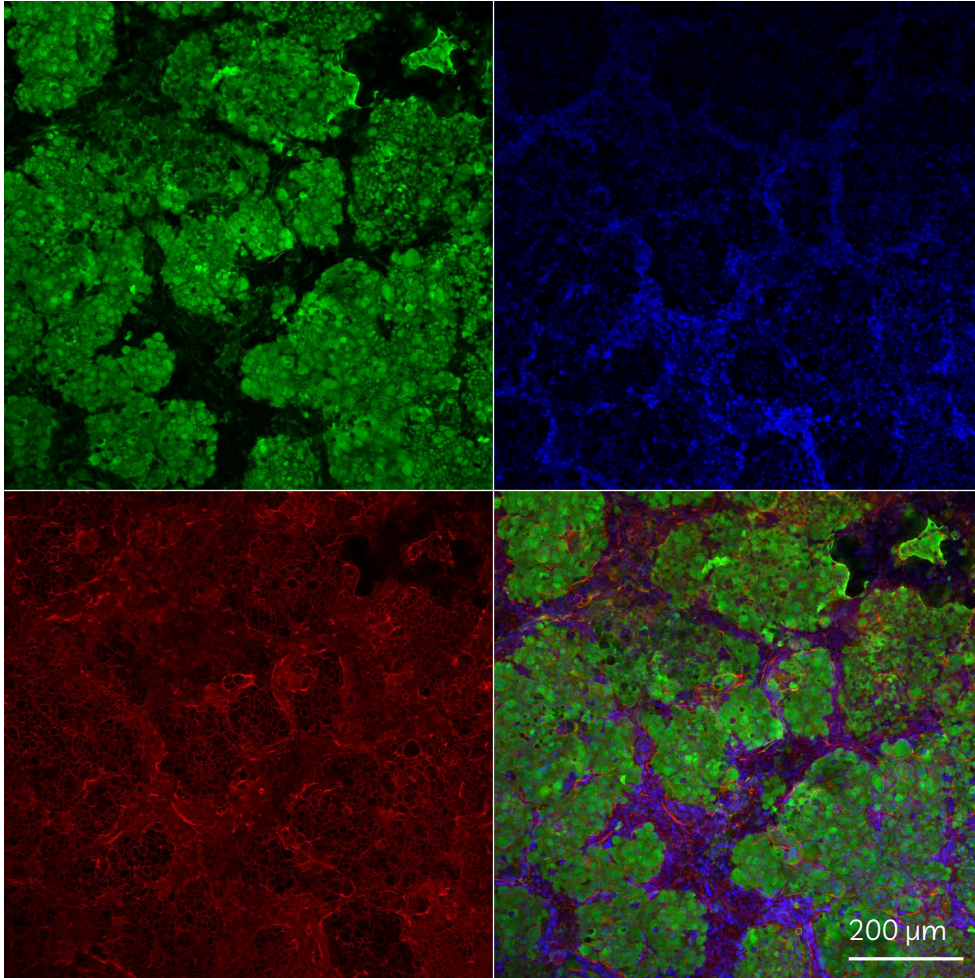
Complete culturing procedure



17 Prepared and performed by Lena Koch, Applied Stem Cell Technologies

Figures made with BioRender, Copyright (c) by Locsense BV All Rights Reserved.

Caco-2, HT29-mtx co-culture



Green = Wheat germ agglutinin (WGA) → mucus marker

Blue = DAPI → nucleus marker

Red = Actin → cytoskeleton marker

18 Prepared and performed by Lena Koch, Applied Stem Cell Technologies

Index

- Function of the gut
- Gut mimic: *in vitro* translation
- Impedance spectroscopy & transepithelial electrical resistance (TEER)
- Interfering with the tight junction/barrier
- Measuring the tight junction/barrier
- Conclusions

Impedance spectroscopy and Transepithelial electrical resistance (TEER)

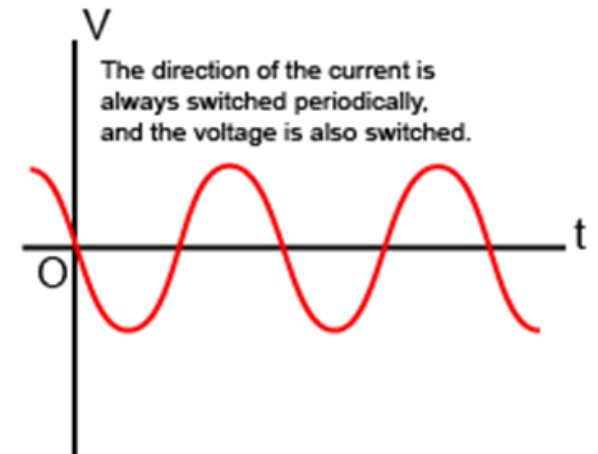
Impedance = The combined effect of **resistance** and **reactance** in a circuit [1]

All the obstacles that impede the current to flow to the other electrode

TEER = The measurement of **electrical resistance** across a cellular layer [1]

Tight junction resistance between the cells = barrier function

Alternating Current (AC)



Impedances = Resistance + Reactance

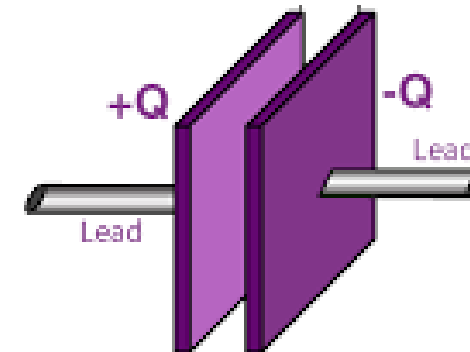
Resistance

Constant impedance of electrical current



Reactance - capacitance

Capability to store and release electrical charge



Resistance and capacitance

Resistance = swimming

It takes longer to get to the other side, but going is continuous

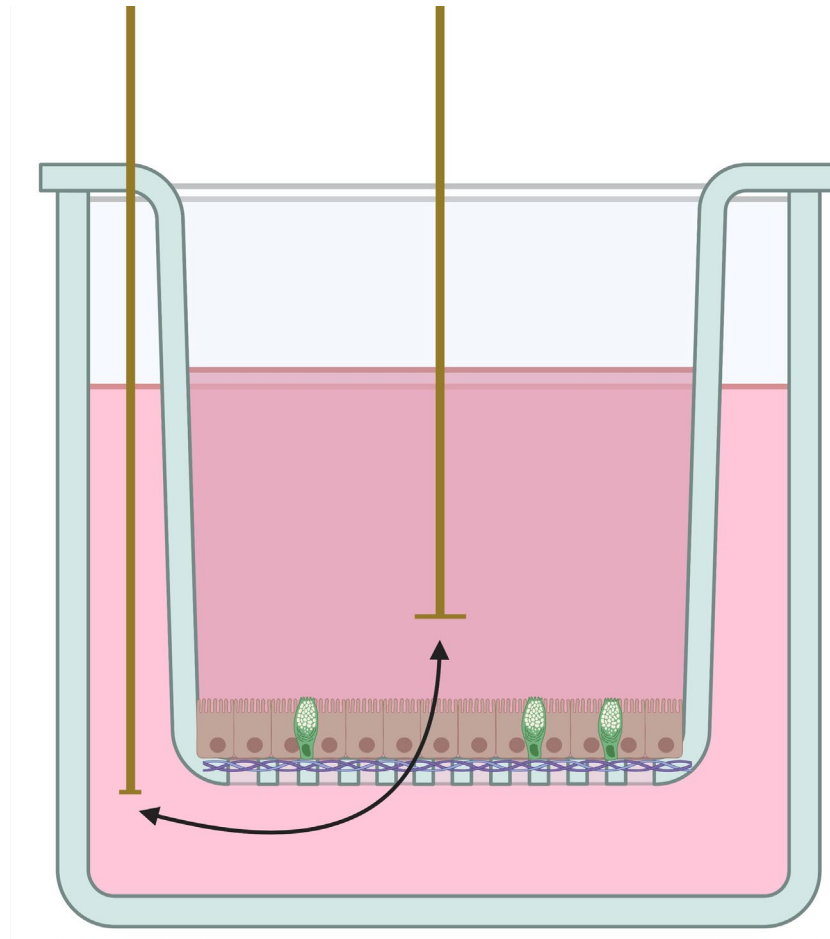


Capacitance = a ferry boat

It has to wait for other passengers, but it travels fast when it is going



Impedance measurement

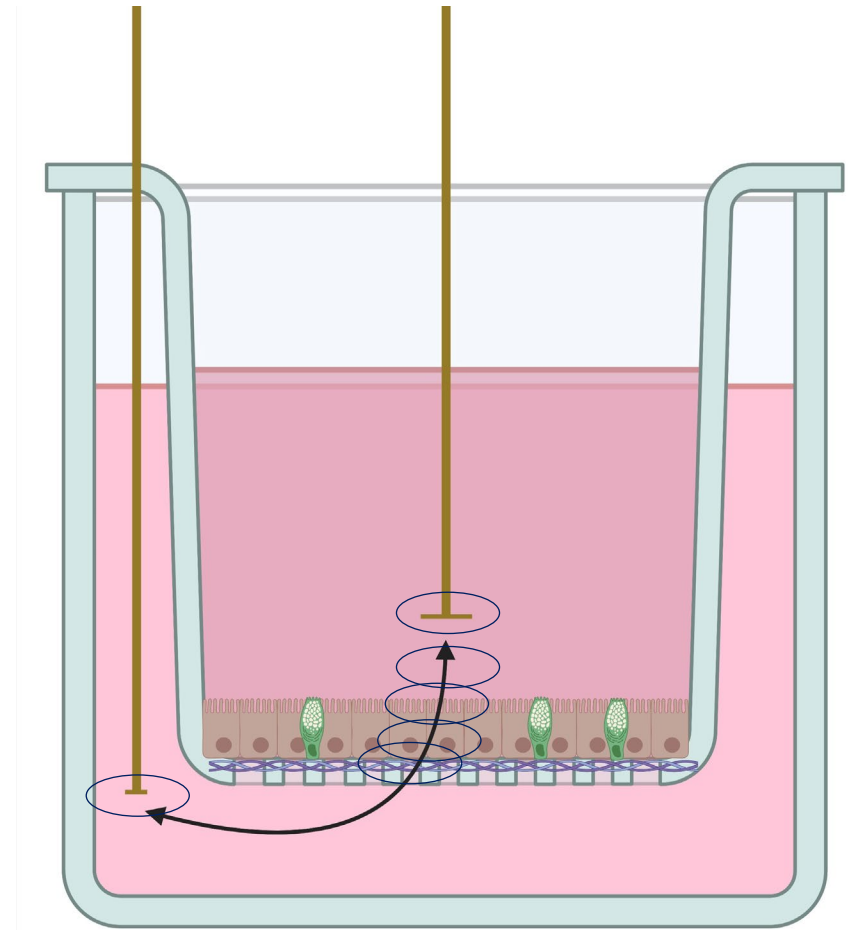




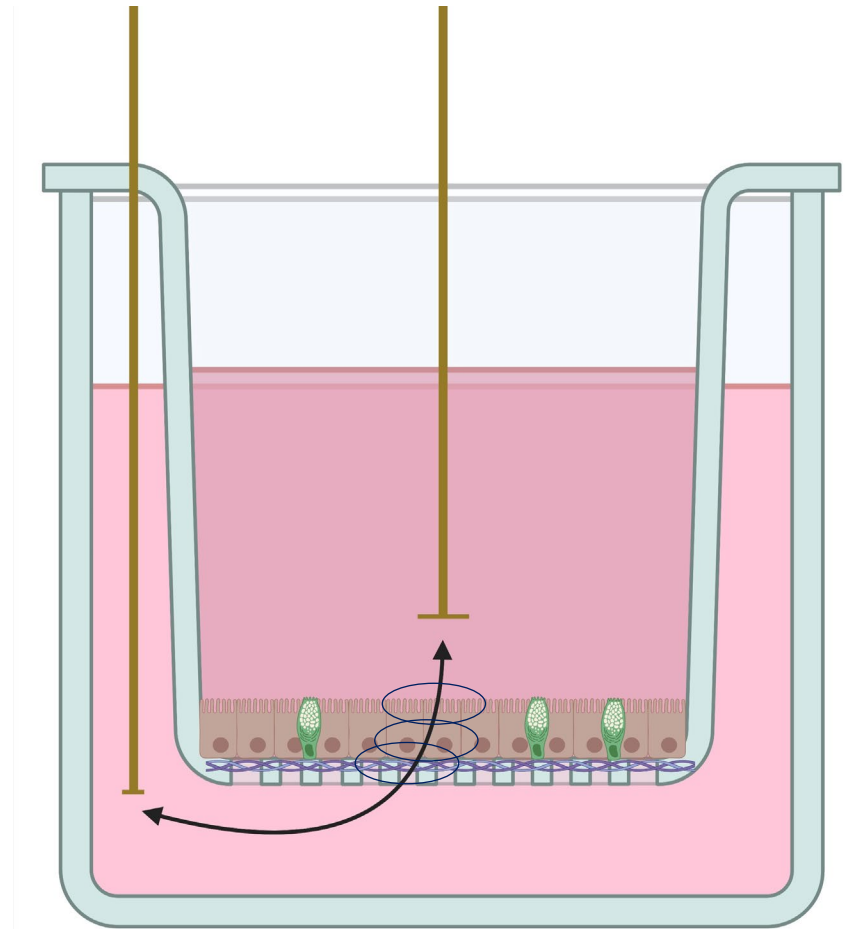
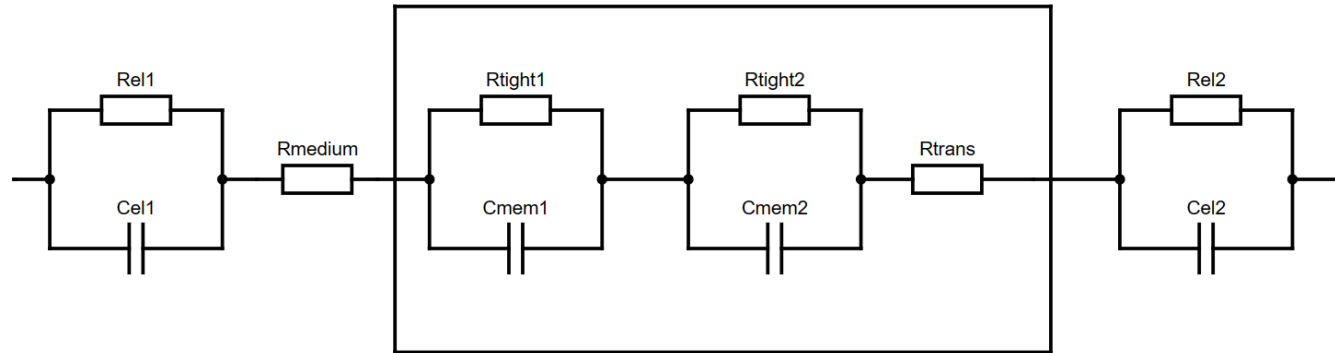
Circuit model fitting

Components of the circuit:

1. $R_{el1} + C_{el1}$ = Resistance + capacitance of electrode 1
2. R_{medium} = Resistance of the medium
3. $R_{tight1} + C_{mem1}$ = the tight junction resistance and membrane capacitance of the apical side
4. $R_{tight2} + C_{mem2}$ = the tight junction resistance and membrane capacitance of the basolateral side
5. R_{trans} = resistance of the coated transwell
6. $R_{el2} + C_{el2}$ = Resistance + capacitance of electrode 2



Transepithelial electrical resistance (TEER)



The cells influence

$R_{\text{tight1}} + C_{\text{mem1}}$ = the tight junction resistance and membrane capacitance of the apical side

$R_{\text{tight2}} + C_{\text{mem2}}$ = the tight junction resistance and membrane capacitance of the basolateral side

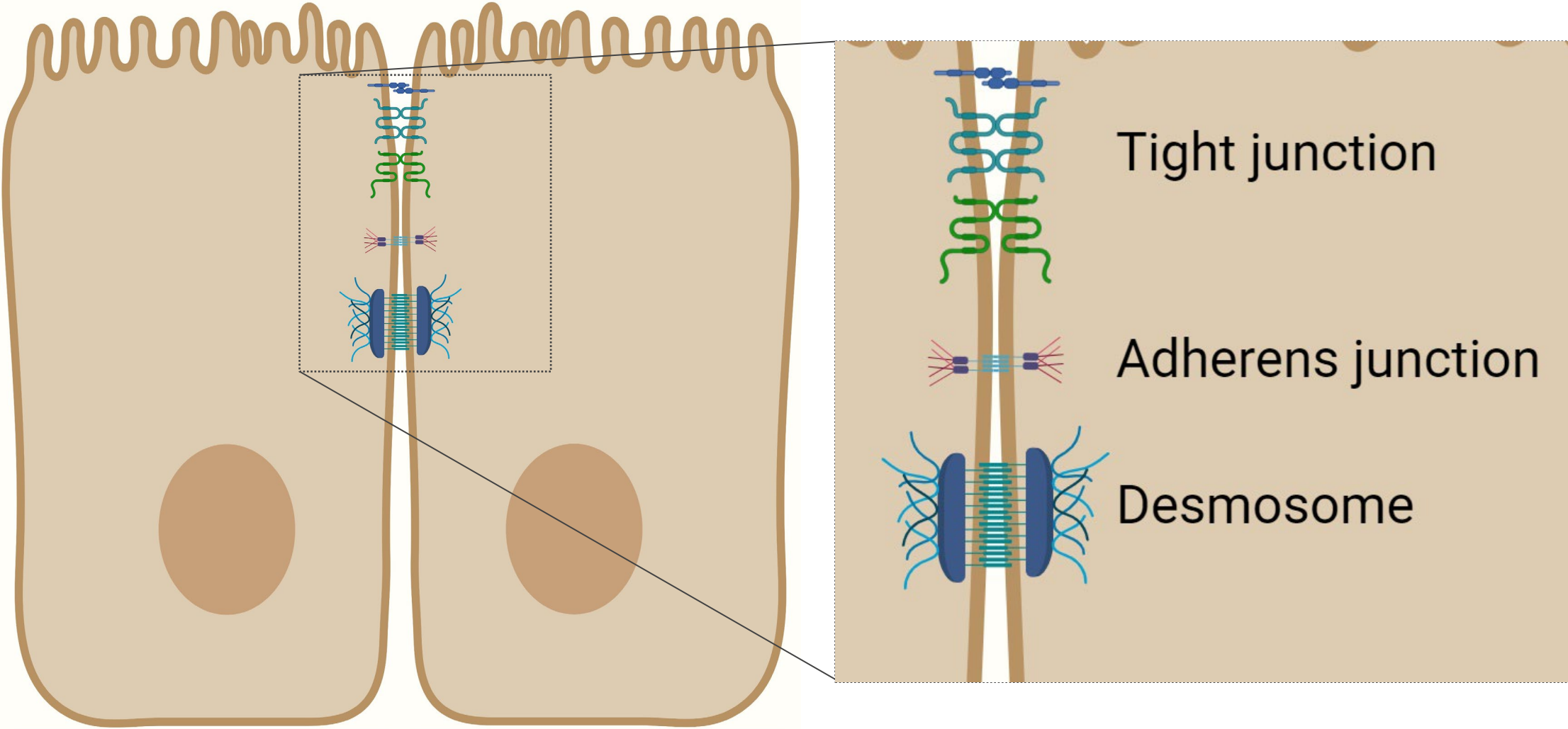
R_{trans} = resistance of the coated transwell

The total TEER = $R_{\text{tight1}} + R_{\text{tight2}}$

Index

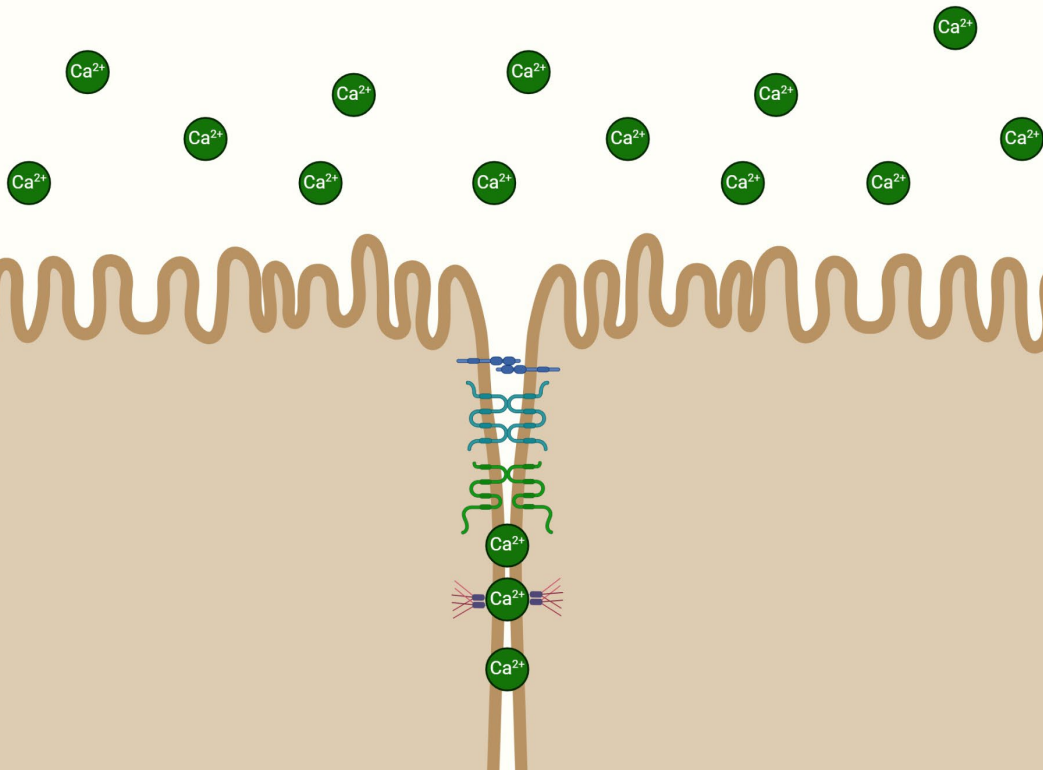
- Function of the gut
- Gut mimic: *in vitro* translation
- Impedance spectroscopy & transepithelial electrical resistance (TEER)
- Interfering with the tight junction/barrier
- Measuring the tight junction/barrier
- Conclusions

Tight junctions

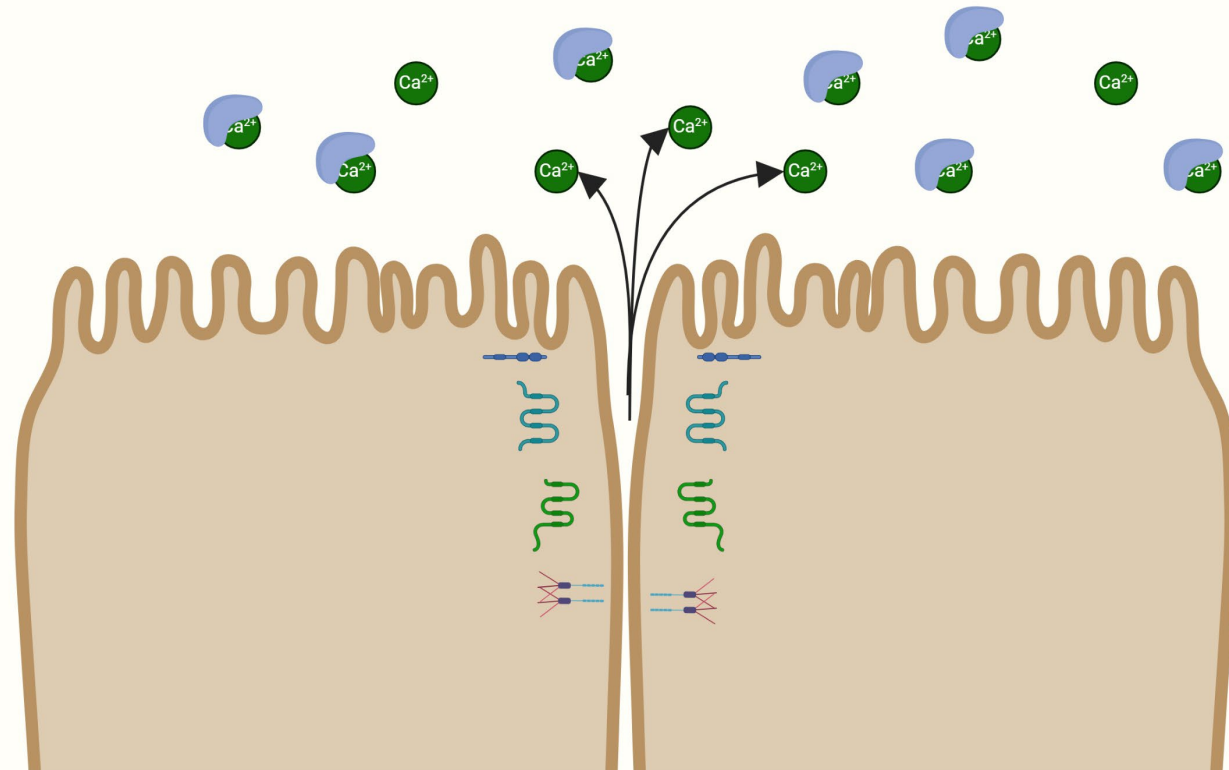


EGTA and tight junctions

High levels of free calcium



EGTA mediated depletion of free calcium



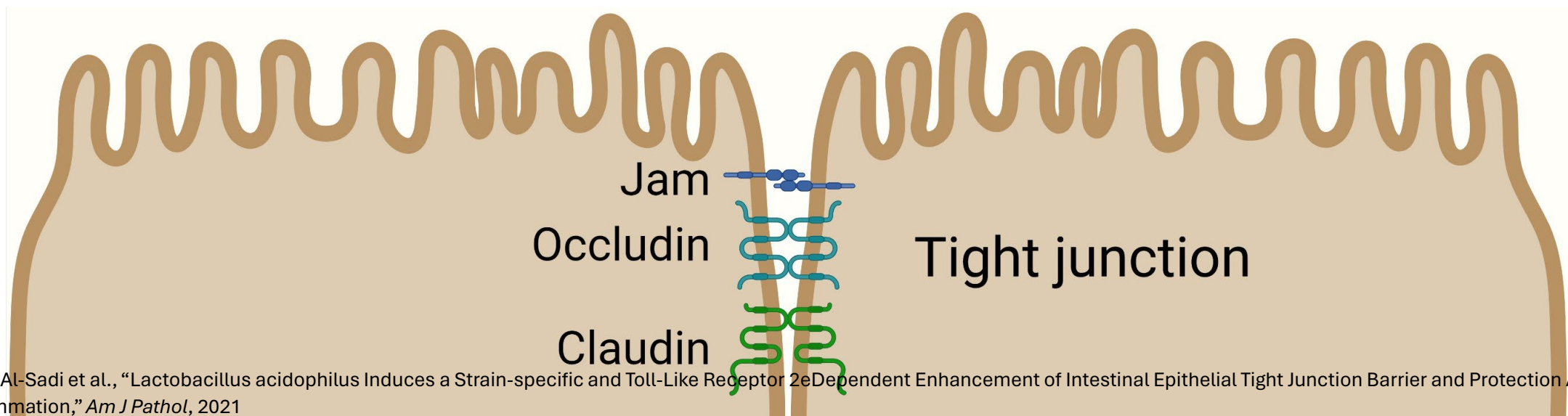
29

Lactobacillus acidophilus (LA) and tight junctions [1]

LA attaches to the apical side of the Enterocyte → ↑ TLR-2 membrane protein

TLR-2 membrane protein → ↑ interaction with TLR-1 and TLR-6

TLR2/TLR1 and TLR2/TLR6 → ↑ occludin

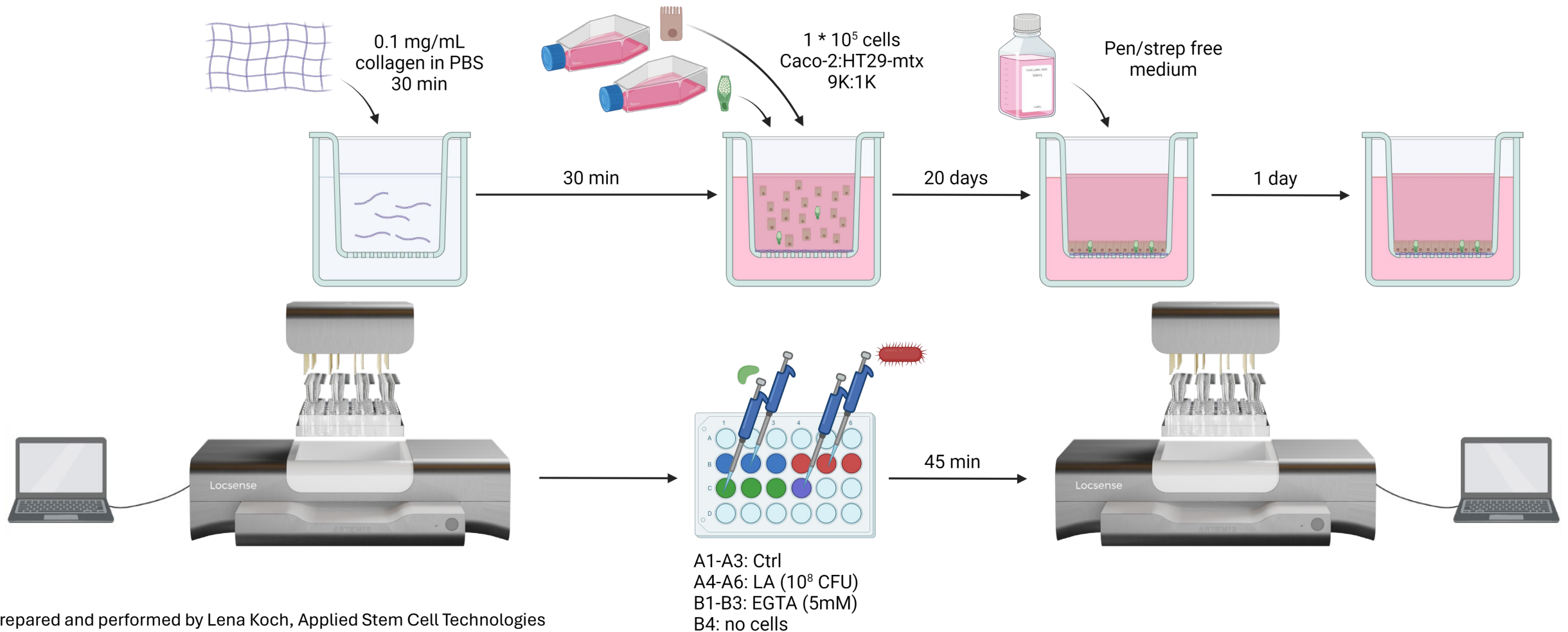


[1] R. Al-Sadi et al., "Lactobacillus acidophilus Induces a Strain-specific and Toll-Like Receptor 2-Dependent Enhancement of Intestinal Epithelial Tight Junction Barrier and Protection Against Intestinal Inflammation," *Am J Pathol*, 2021

Index

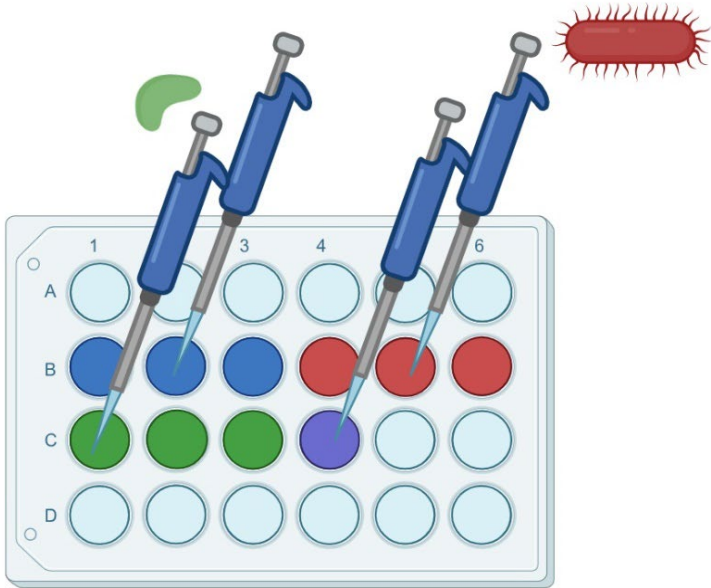
- Function of the gut
- Gut mimic: *in vitro* translation
- Impedance spectroscopy & transepithelial electrical resistance (TEER)
- Interfering with the tight junction/barrier
- Measuring the tight junction/barrier
- Conclusions

Complete procedure



32 Prepared and performed by Lena Koch, Applied Stem Cell Technologies

Conditions



A1-A3: Ctrl
A4-A6: LA (10^8 CFU)
B1-B3: EGTA (5mM)
B4: no cells

T=0 → Measurement 1

After measurement → LA is added to LA wells

T=24 → Measurement is performed

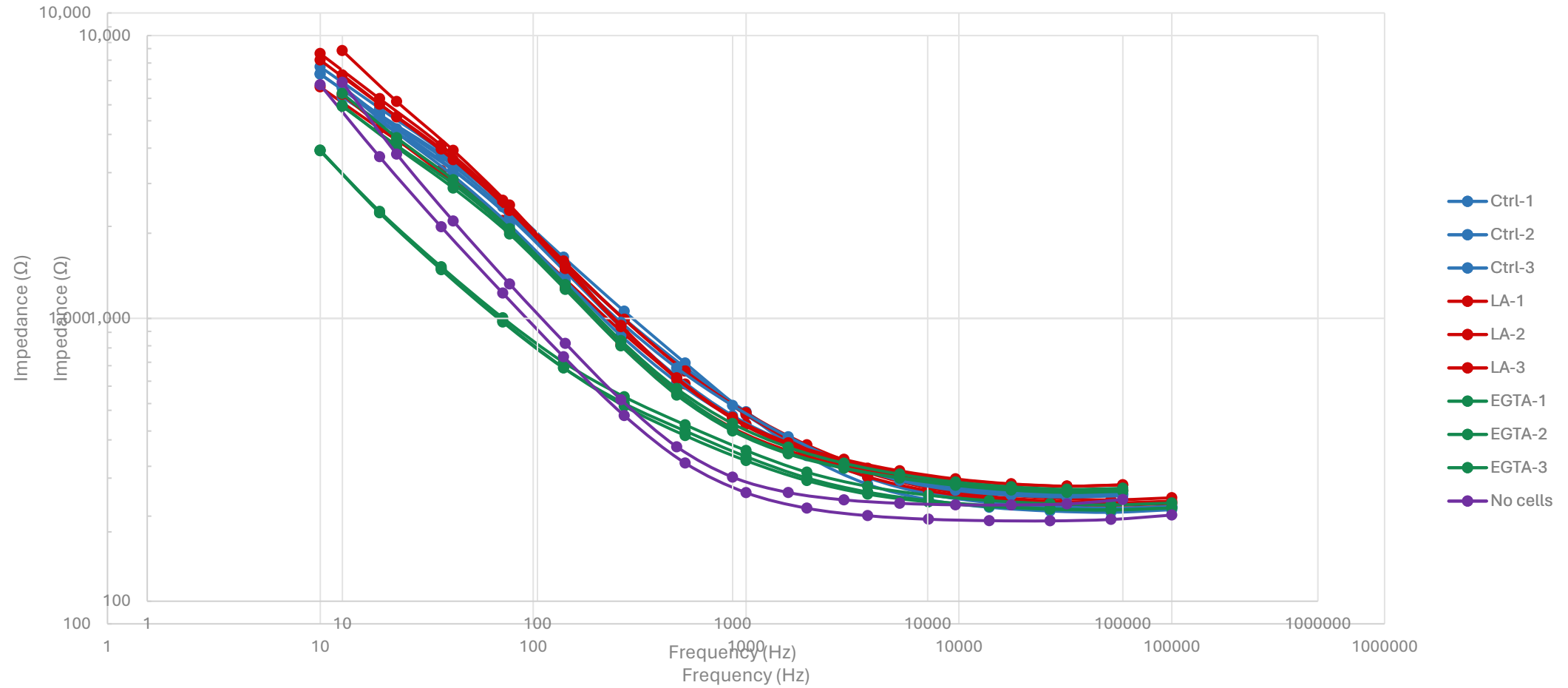
After T=24 → medium of EGTA wells is removed, EGTA in PBS is added for 45 min → old medium (+10% fresh) is added

T=25 → Measurement is performed

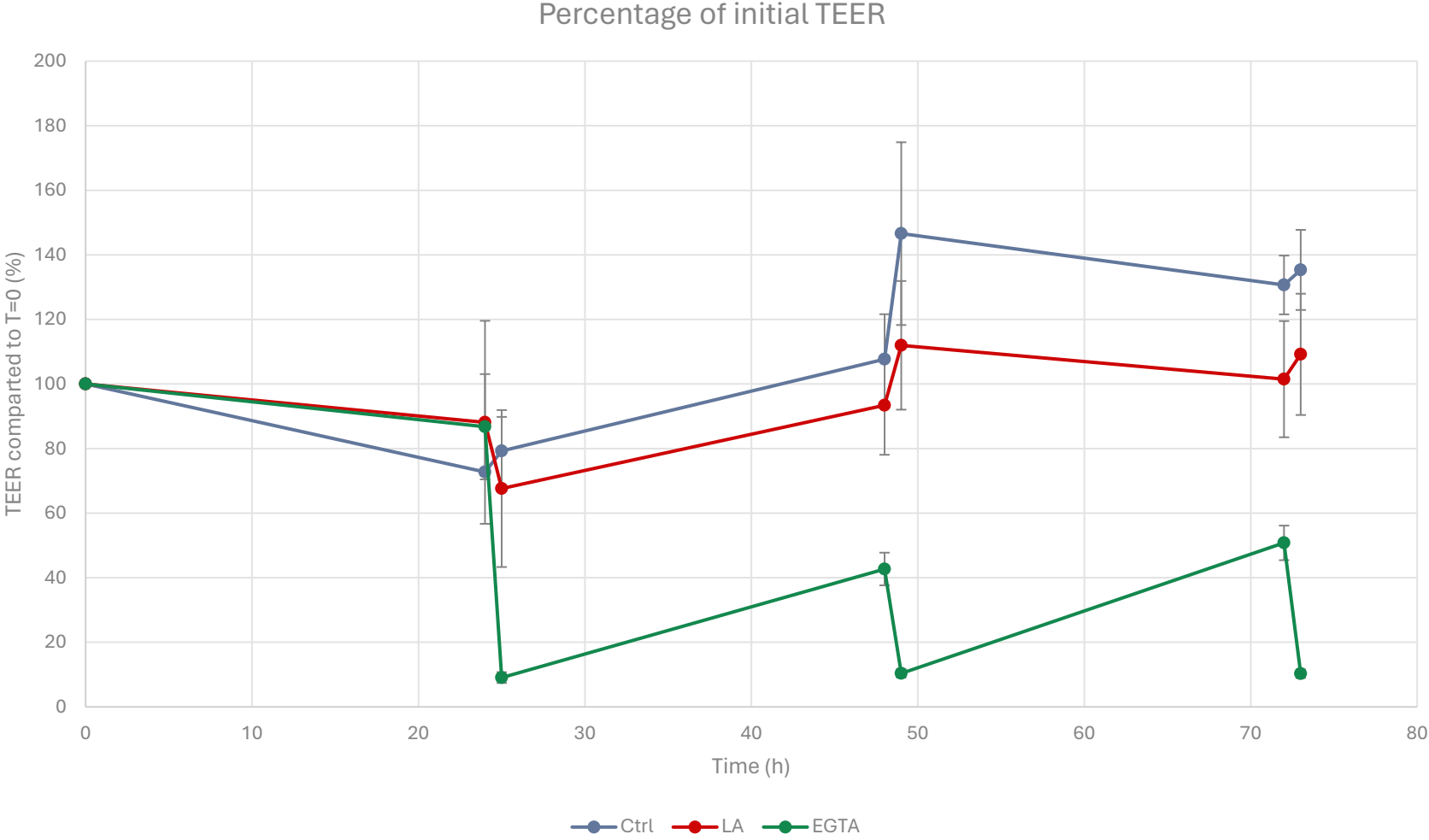
Repeated at T=48 + T=49 and T=72 + T=73

Raw impedance data

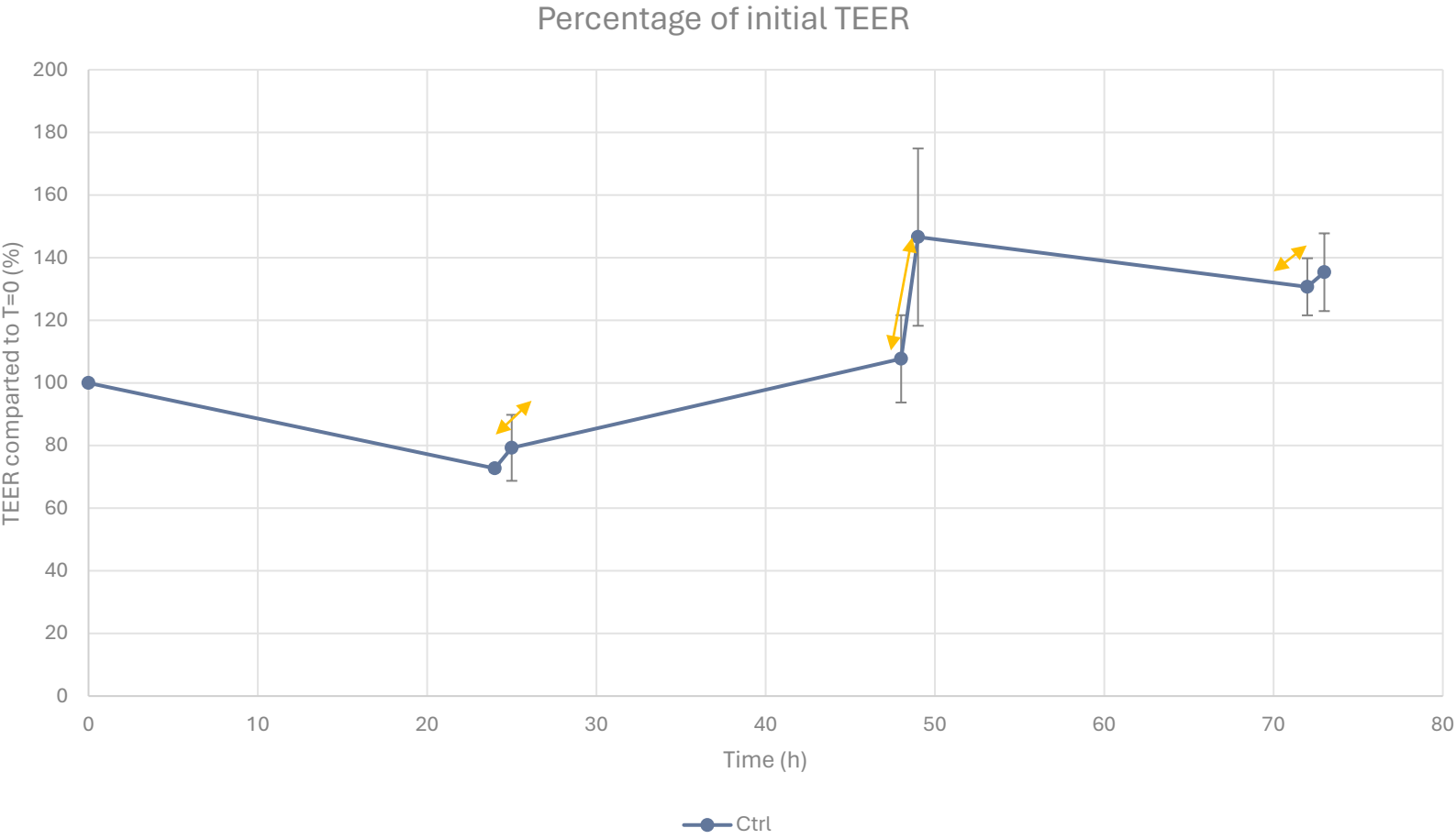
$\tau = 0.8$



TEER results



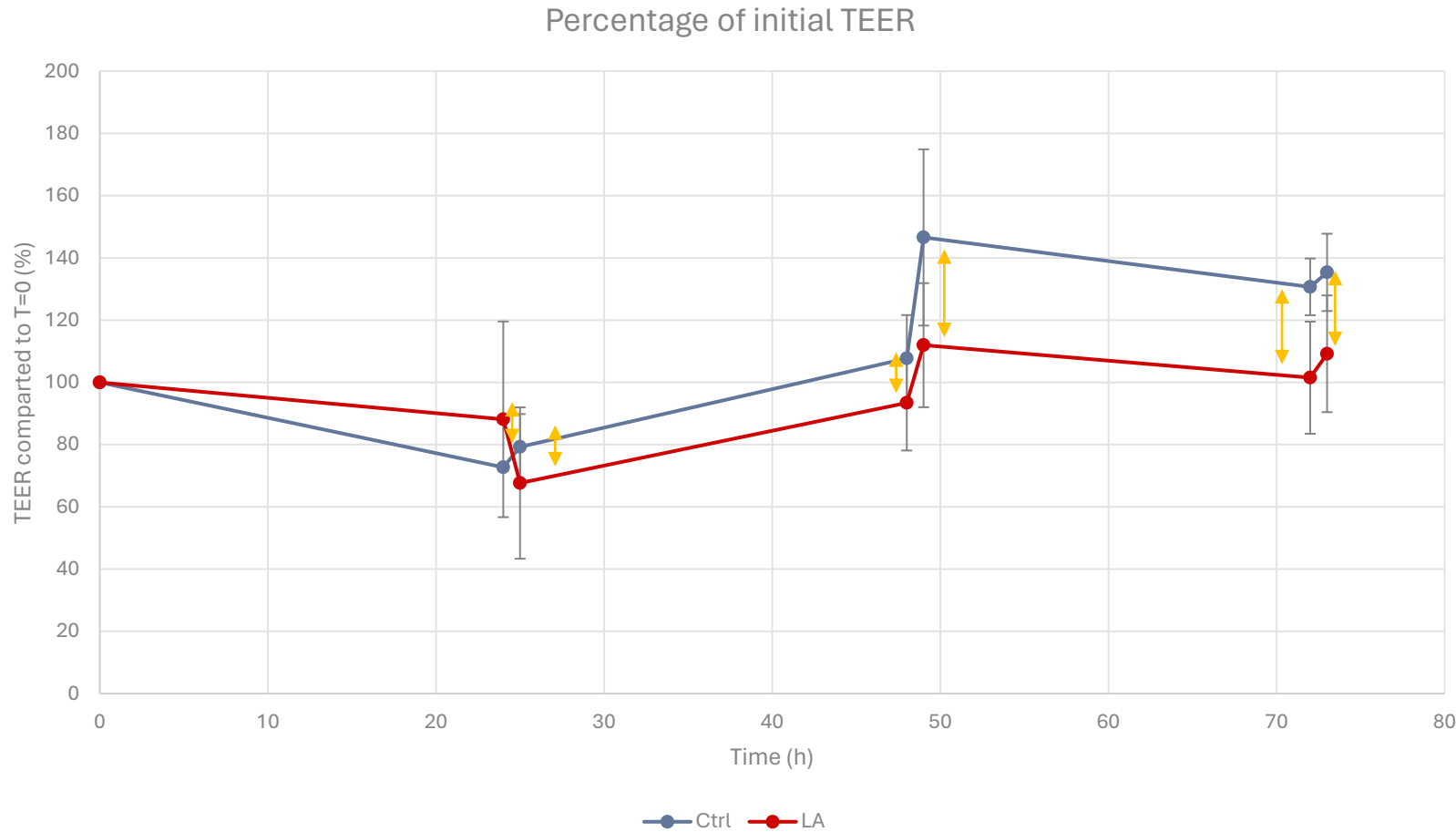
Temp. does not significantly change TEER



Difference between 2 measurements

| | Ctrl |
|-------|------|
| Day 1 | 6,5 |
| | ns |
| Day 2 | 38,9 |
| | ns |
| Day 3 | 4,7 |
| | ns |

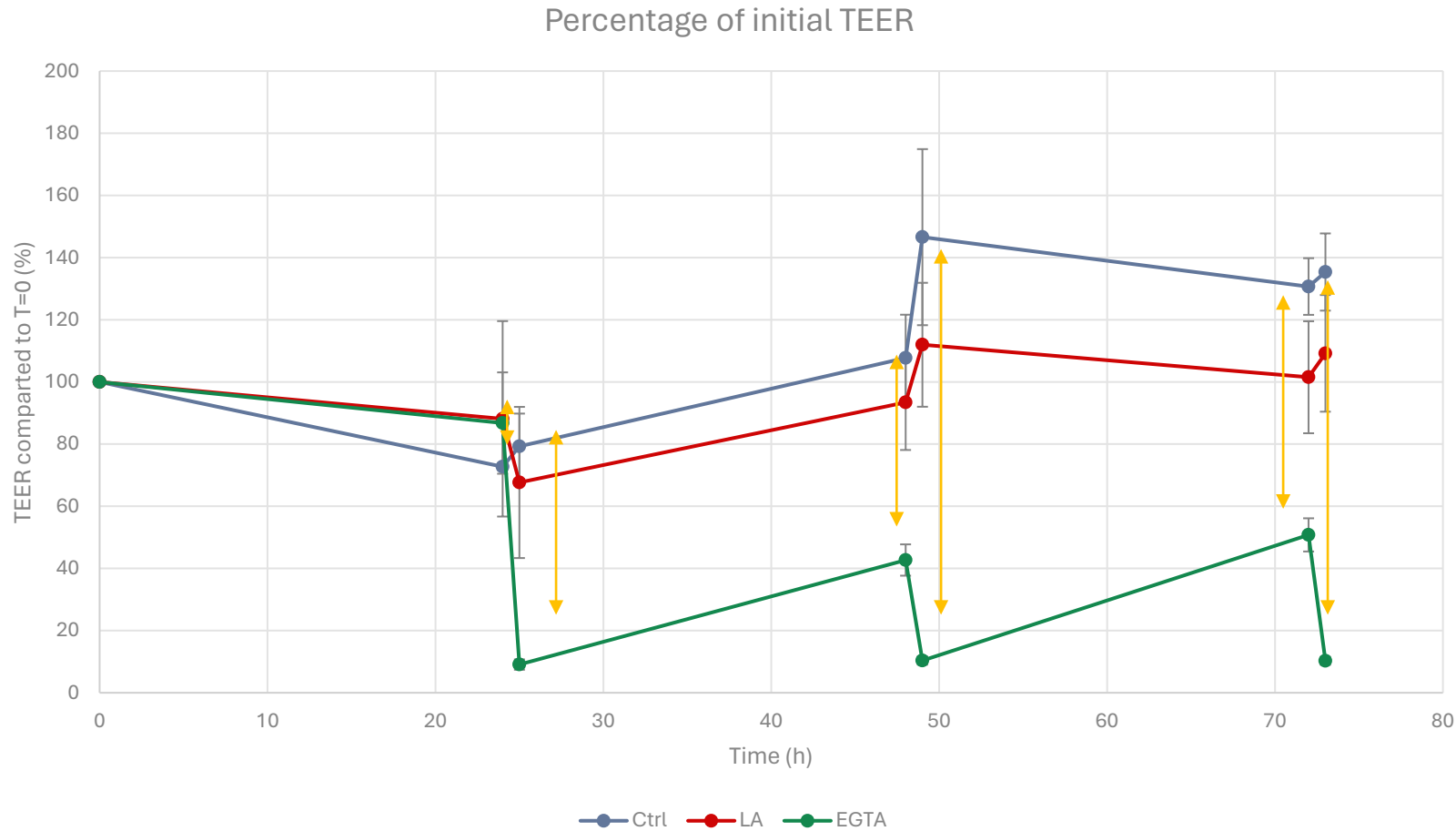
LA does not significantly change the TEER



Difference between Ctrl and LA

| Time (h) | LA | |
|----------|-------|----|
| 24 | 15,4 | ns |
| 25 | -11,6 | ns |
| 48 | -14,3 | ns |
| 49 | -34,6 | ns |
| 72 | -29,2 | ns |
| 73 | -26,2 | ns |

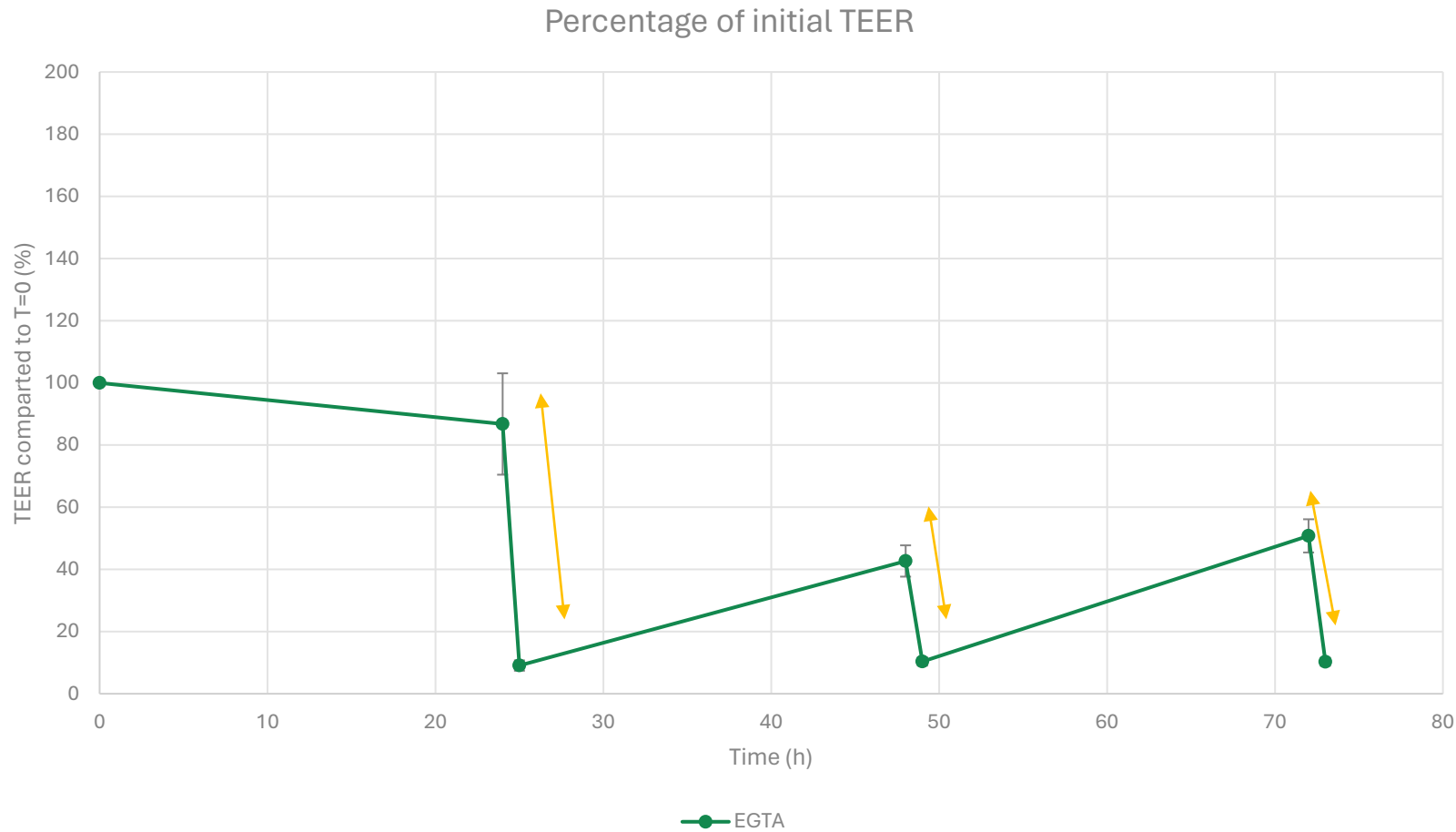
EGTA does significantly change the TEER



Difference between Ctrl and LA or EGTA

| Time (h) | LA | | EGTA | |
|----------|-------|--------------|--------|--------------|
| | Value | Significance | Value | Significance |
| 24 | 15,4 | ns | 14,0 | ns |
| 25 | -11,6 | ns | -70,2 | *** |
| 48 | -14,3 | ns | -65,0 | ** |
| 49 | -34,6 | ns | -136,2 | ** |
| 72 | -29,2 | ns | -79,9 | *** |
| 73 | -26,2 | ns | -125,1 | **** |

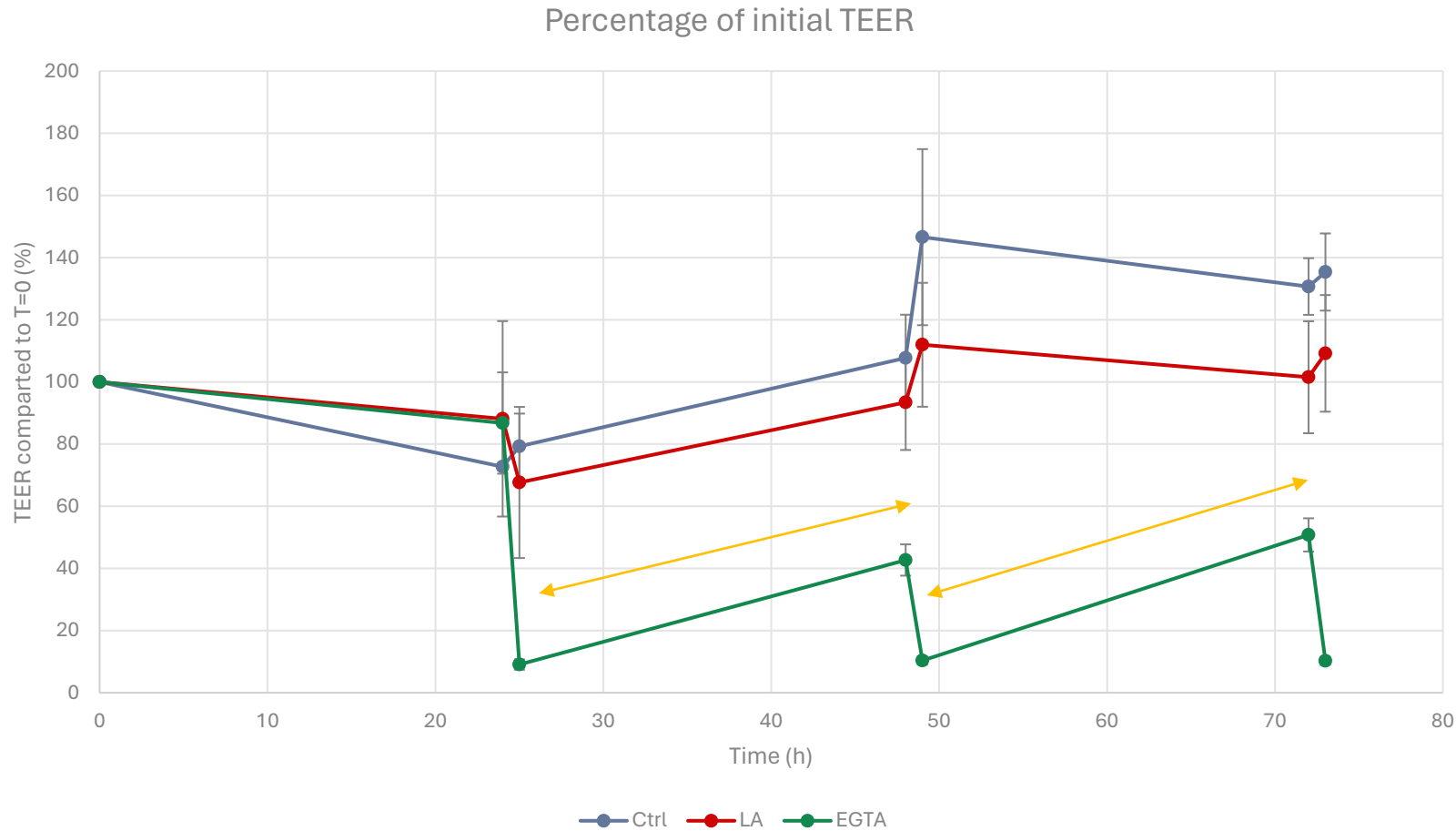
EGTA does significantly change TEER in 45 min



Difference between before and after EGTA addition

| | EGTA |
|-------|-------|
| Day 1 | -77,7 |
| | *** |
| Day 2 | -32,3 |
| | *** |
| Day 3 | -40,5 |
| | *** |

EGTA treatment is reversible



EGTA is capable of partial recovery after treatment

| | Difference | |
|-------|------------|----|
| 25-48 | 33,6 | ** |
| 49-72 | 40,4 | ** |

Index

- Function of the gut
- Gut mimic: *in vitro* translation
- Impedance spectroscopy & transepithelial electrical resistance (TEER)
- Interfering with the tight junction/barrier
- Measuring the tight junction/barrier
- Conclusions

Conclusions cell culturing

- A Caco-2 (ATCC[®] HTB-37[™]) HT29-mtx co-culture is a representable *in vitro* gut mimic
- 9:1 ratio
- Cultured in DMEM-F12-GlutaMAX culturing medium
 - 15% FBS, compromise
- 21-day pre-culture

Conclusions impedance spectroscopy

- Impedance spectroscopy is a non-destructive measurement technique
- Circuit model fitting enables to extract TEER data from complex impedance measurements
- The Artemis device
 - Automated measurement
 - User friendly
 - Time saving
 - Reduces cross contamination risks

Conclusions EGTA and LA

- Temperature influences, but changes the TEER non-significantly
- Lactobacillus acidophilus does not significantly influence the tight junction barrier
- EGTA successfully hampers the tight junctions
- EGTA barrier function impairment is reversible

Collaboration

This webinar was a collaboration between Locsense B.V. and Applied Stem Cell Technologies, University of Twente, the Netherlands

Work is conceptualized and performed by Lena Koch

Conceptualization and funding acquisition Kerensa Broersen

Artemis TEER/impedance spectroscopy instrument is provided by Locsense

THANKS

Follow Locsense on LinkedIn for updates



info@locsense.nl
www.locsense.nl
Institutenweg 25
7521PH Enschede



Questions?



For more information:

www.atcc.org/tox

<https://locsense.nl>

