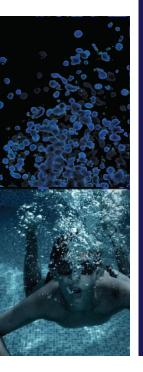


Best Practices for Cell Culture

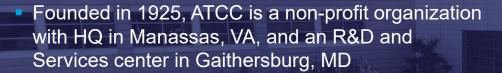
Credible Leads to Incredible™

Steven Budd, MS, MBA Product Line Business Specialist, ATCC









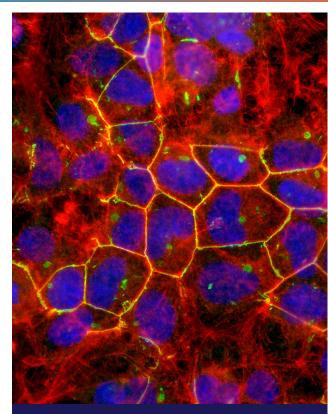
- World's largest, most diverse biological materials and information resource for cell culture – the "gold standard"
- Innovative R&D company featuring gene editing, microbiome, NGS, advanced models
- cGMP biorepository

- 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 onethird with advanced degrees



Agenda

- Cell and Media handling
 - Cryopreservation
 - Aseptic technique and contamination
 - Authentication
 - Mycoplasma effects and detection
- Workflow and how to choose the best model
- Applications
 - Transfection
 - Cell viability testing



HEK-293 (ATCC[®] CRL-1573[™])





Cell handling/media handling



Thawing cells

- Thaw in 37°C water bath for approximately 2 minutes with gentle agitation
- Spray vial with 70% ethanol
- Transfer to 10 mL centrifuge tube with 9 mL of appropriate growth media (10% FBS)
- *Centrifuge, resuspend in 2 mL of growth media
- Transfer to cell culture vessel

When bringing out of liquid nitrogen, thaw as quickly as possible

*For certain primary cells, centrifugation may be detrimental, refer to specific protocol







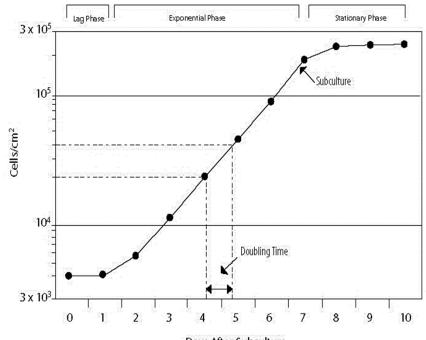
Cell expansion

- After thawing, cells should be plated in an appropriate cell culture vessel with complete media
- 24 hours after seeding, check for confluency
- Note, primary cells may take up to several days to reach 80% confluency for subculturing





Cell expansion



Days After Subculture **Figure 1**. Growth curve for cells grown in culture. Cells should be subcultured while still in the exponential phase.

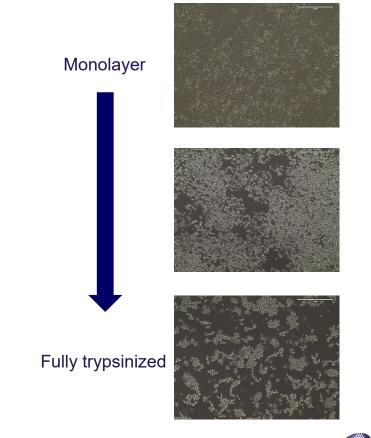


ATCC Animal Cell Culture Guide, 2014.

Trypsinization

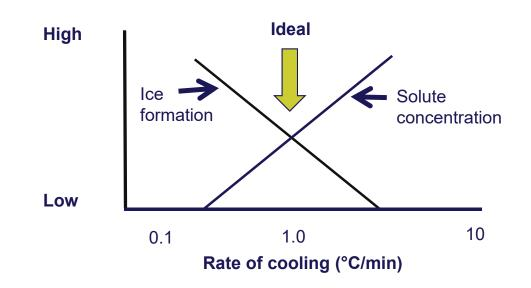
At 80% confluency (primary cells), cells can be passed using Trypsin-EDTA

- Using warm trypsin-EDTA for about 3-5 minutes, cells will detach with gentle agitation
- Trypsin-EDTA for Primary Cells (ATCC® PCS-999-003™) is a low concentration formula (.05% Trypsin and .002% EDTA) – necessary for primary cell survival
- A Trypsin Soybean Neutralizing Solution (ATCC® 30-2104[™]) is also needed to prevent cell damage





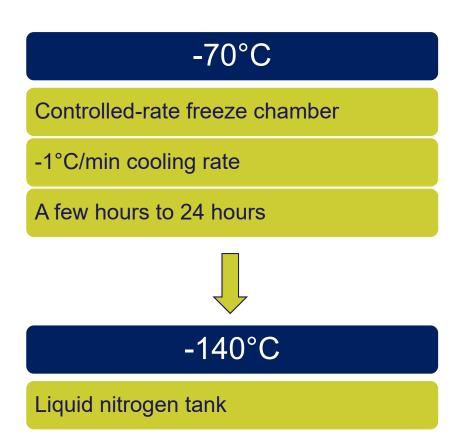
Cryopreservation



- High levels of ice formation and increased solute concentration have a negative impact on cell viability
- Optimal cooling rate for cell viability is 1 to 3°C/min



Freezing down cells







Low temperature storage



For the best security, always store your cells in liquid nitrogen freezers

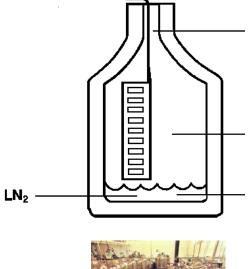




Low temperature storage

Mammalian cells

- Long-term storage should be below -140°C
- Vials should be stored in a liquid nitrogen unit above the volume of liquid at the bottom of the tank
- This temperature should be between -140°C and -180°C





ATCC[°]



Cell characterization

Characterizing cells

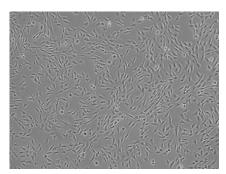
Cell count before plating
Calculating % viability

Morphology

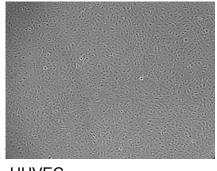
 Make sure the morphology is consistent with cell type

Doubling time

 Contamination from other cell types can affect growth rate







HUVEC



Contamination in cultures

Sources

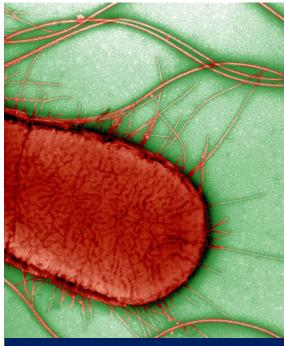
- Contaminated cell lines
- Improper aseptic technique

Types

- Microbial bacteria, mycoplasma, fungi, viruses
- Cellular cross contamination

Signs

- Turbid media
- Rapid decline in pH color change
- Morphological changes
- Filamentous structures
- Changes in physiological reponses



E. coli, image courtesy of David Gregory and Debbie Marshall

ATCC°

Bacterial and Fungal Contamination

Personnel and equipment

- Poor culturing practices
- Dust and aerosol

Contamination

- Aerosol dispersion of contaminated cell cultures
- Faulty laminar flow

Culture reagents

- Sera
- Media
- Reagents





Contamination prevention and aseptic technique

Good aseptic technique

- Make it difficult for microorganisms to invade culture vessels
 - -Sealed cultured vessels
 - -Vented cap flasks
- Disposable aspirators
 - -Cell culture hoods with good laminar flow
 - Do not use as a storage area!
- Spray media bottles/reagents with alcohol





Contamination prevention and aseptic technique

Use small volumes of reagents at a time

Aliquot stock solutions and reagents

Always wear clean lab coats and protective clothing

Use seed stocks

Create master stocks

Avoid using antibiotics in media!

- Can contribute to chronic contamination
- Rarely prevents contamination
- Toxic to cells





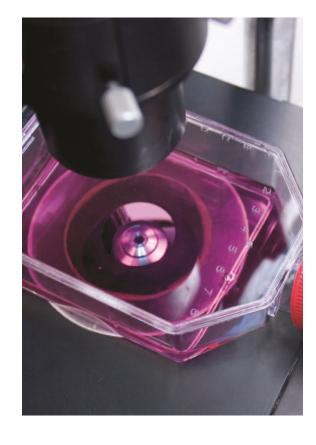
Mycoplasma contamination

Not easily detected

- Does not cause media turbidity
- Does not alter the pH of the media
- Few metabolic byproducts
- Cannot be detected by microscopy

Results in a number of deleterious effects

- Chromosomal aberrations
- Disruption of nucleic acid synthesis
- Changes in membrane antigenicity
- Inhibition of cell proliferation and metabolism
- Decreased transfection rates
- Changes in gene expression profiles
- Affects virus production
- Cell death



ΔΤCC



Mycoplasma Detection Kit

PCR-based kit

Detects any of the 60 most common mycoplasmas

Kit Components/Procedure

Buffer

- Mix of primers specific to mycoplasma
- Positive control
 - DNA from mycoplasma







Cell cross-contamination

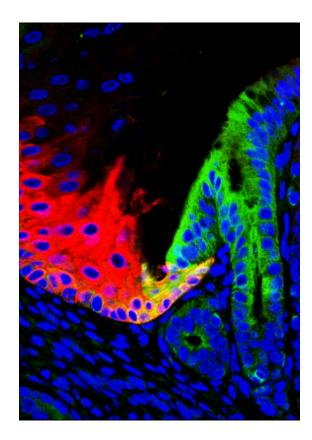
Leads to the replacement of the original cell line with the contaminant

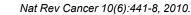
Causes

- Multiple cell lines under the hood at the same time
- Failure to change out pipettes
- Receiving cell lines from other labs

20% of scientific publications include misidentified cultures

50% of preclinical research is not reproducible





Short Tandem Repeats (STR)

- ATCC worked with NIST to pioneer STR profiling for human <u>and</u> mouse cell lines
- ATCC authentication services are simple and inexpensive, after placing your order:
 - Spot
 - Dry
 - Mail
 - Receive you results in three to five days

Report includes:

- Submitted & Matched Allele Calls
- Contamination check
- Comparative output for database comparison
- PDF of the submitted sample profile







Workflow and how to choose the best model



Media choices

Continuous (human and animal) cell lines – generally require a classical medium + 10% FBS

Specialized cell types (primary cells and stem cells require their own specially formulated media, specific to each cell type

Generally maintain cells in the recommended medium

If an application requires a different medium formulation, to transfer to new media:

- Use 1:1 mix (50% old, 50% new media)
- 1:2 mix
- 1:3 mix
- 1:7 mix

Heat inactivation of FBS? Not recommended



ATCC°



Pros and cons of different cell models

	Primary Cells	iPSC-derived	hTERT Immortalized	Continuous Cell Lines
Mimic <i>in vivo</i> tissue type	++++	++	+++	+
Genotypic stability	Diploid	Diploid	Diploid/ Near Diploid	Aneuploid
Proliferative capacity	+	+	+++	+++
Supply	+	+++	+++	+++
Inter-experimental reproducibility	+	+++	+++	+++
Cost	+	+	++	+++
Ease-of-use	+	++	++	+++
Predictability in studies	+++	++	+++	+



Primary cells as a control

Primary neonatal keratinocytes (ATCC® PCS-201-010™) differentiated into physiological epidermis

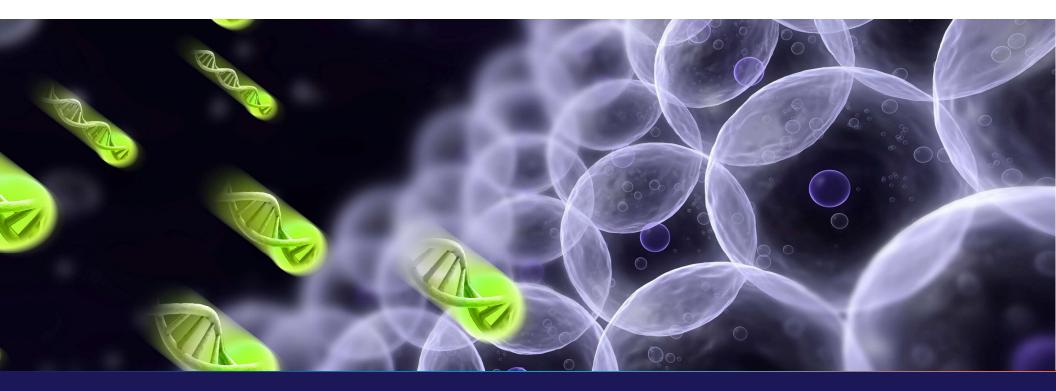
- Continuous cell lines are cells isolated from primary tissue (often a tumor) that have mutated to survive a "crisis"
- Continuous cell lines have deviated from original source

In every continuous cell line experiment, primary cells should be used as one of the controls



ATCC°

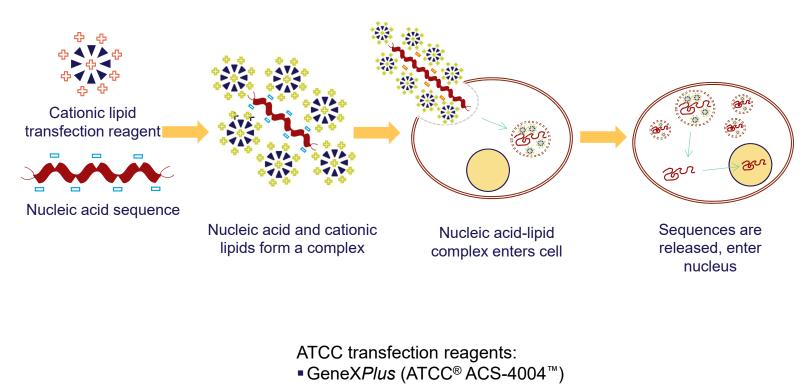




Transfection



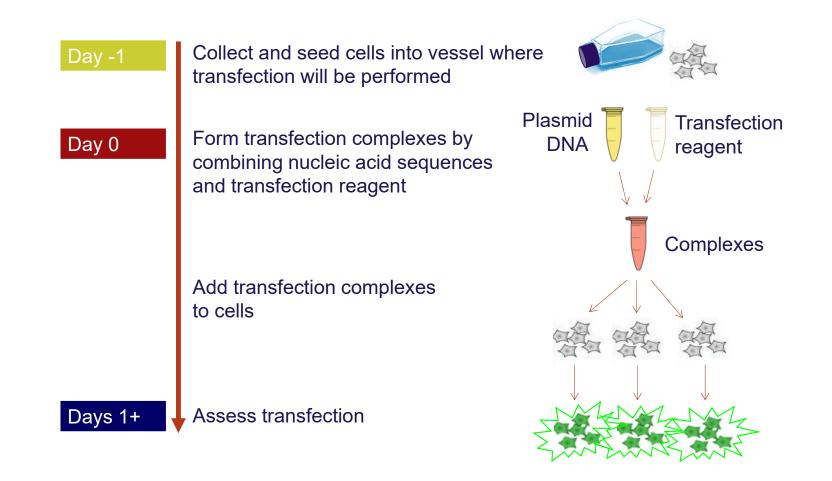
Mechanism of lipid-based transfection





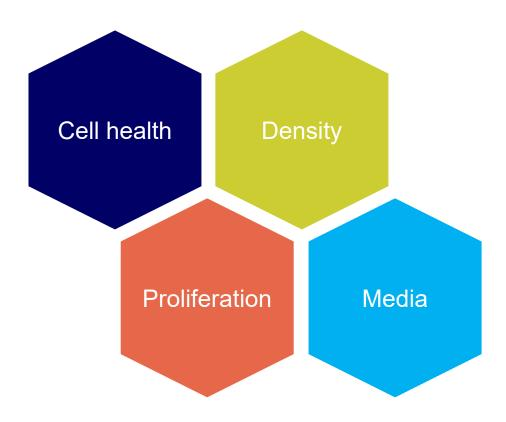
■ TransfeX[™] (ATCC[®] ACS-4005[™])

Typical transfection workflow



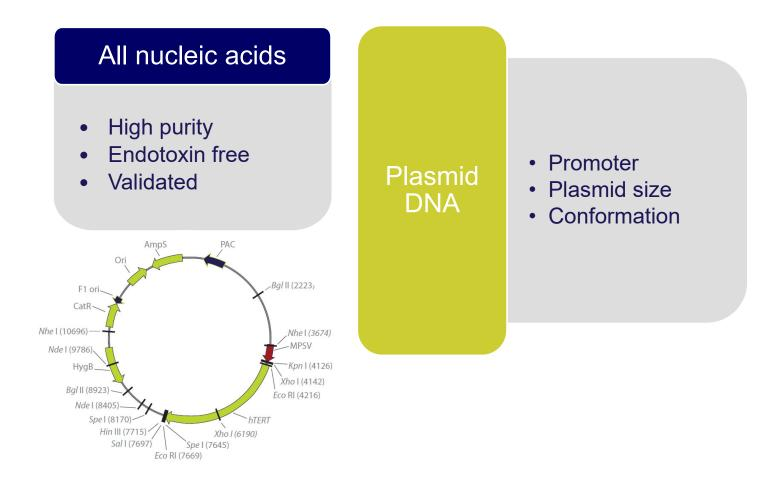


Cell culture conditions



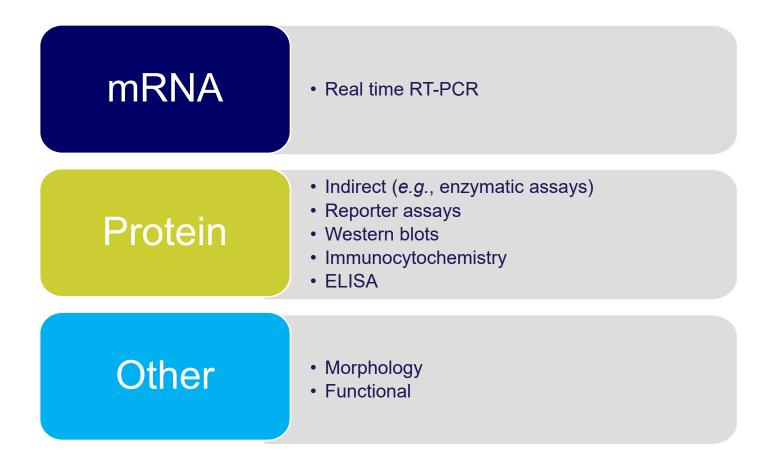


Nucleic acids



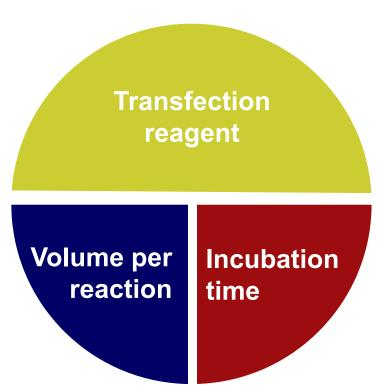


Transfection best practices – assay methods



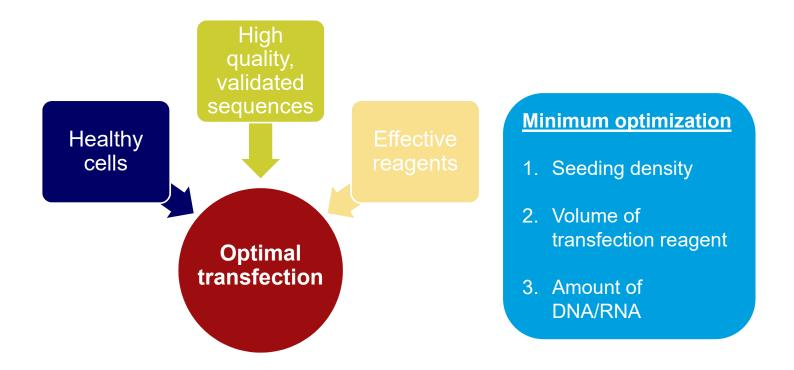


Transfection reagents

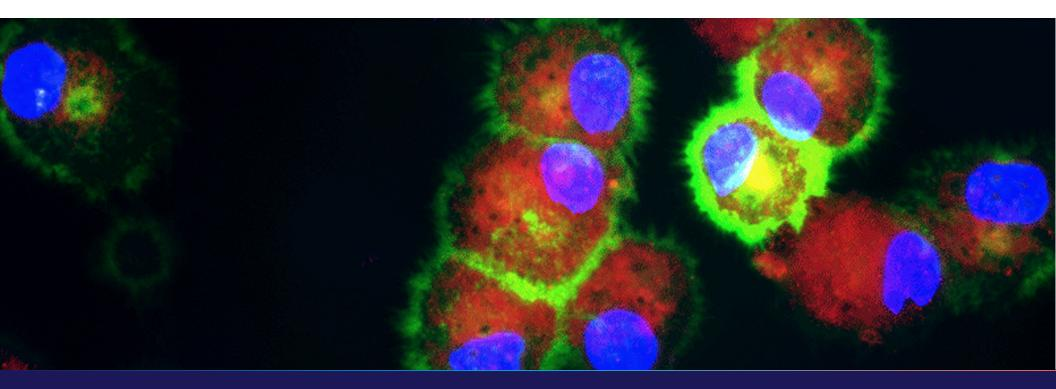




Best practices summary









Quantitative evaluation of cell proliferation rate and response to external factors that affect cell viability

- Commonly used for cytoxicity, high-throughput screening (e.g., drug development)
- Uses tetrazolium salts in a colorimetric method for evaluating cell populations

MTT Cell Proliferation Assay (ATCC[®] 30-1010K[™])

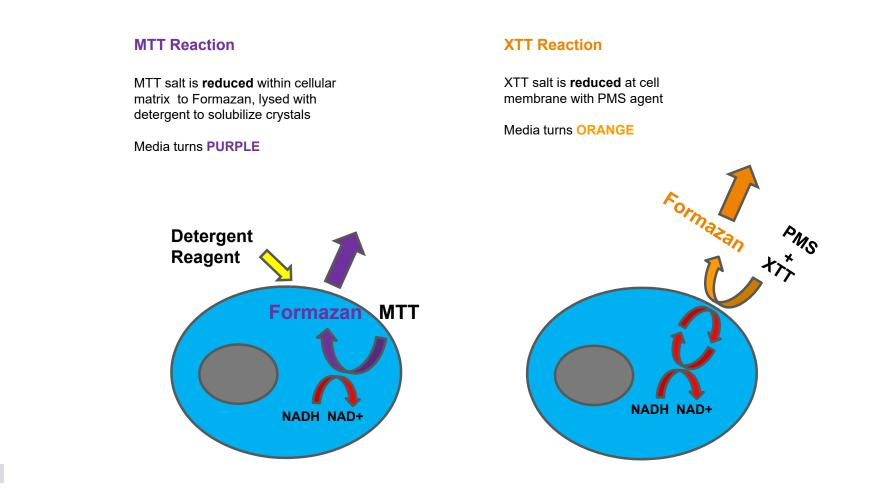
 Tetrazolium MTT (3-(4, 5-dimethylthiazolyl-2)-2, 5diphenyltetrazolium bromide)

XTT Cell Proliferation Assay (ATCC[®] 30- 1011K[™])

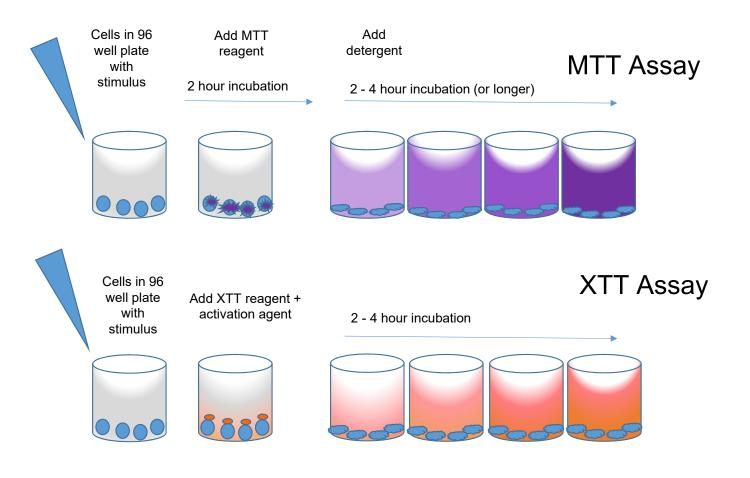
 Tetrazolium XTT (sodium 2,3,-bis(2-methoxy-4-nitro-5sulfophenyl)-5-[(phenylamino)-carbonyl]-2H-tetrazolium



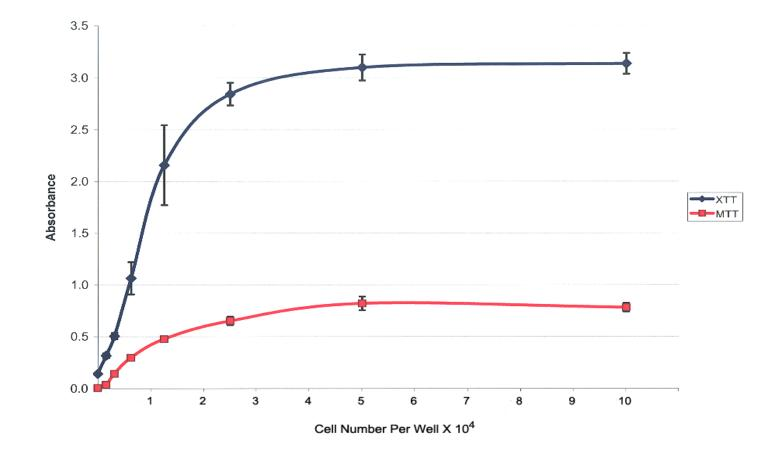
















Summary



Summary

Cell Handling/media handling

- Be sure to employ best practices to eliminate contamination and ensure optimal growth and storage
- Routinely authenticate your cells to ensure reliable results

Cell culture workflow

- Use cell lines for standardization and confirmation of each experiment; use primary cells after standardization to further validate the results with a biologically relevant model
- hTERT-immortalized cells provide continuous growth with near-primary cell performance

Transfection

 Effective lipid transfection starts with healthy cells, purified substrate, and optimal ratio of substrate and lipid

Viability assays

MTT, XTT assays can confirm cell growth characteristics







More webinars, coming soon!



iPSC-derived Primary Cells: Expand Your Cell-based Assays with an Unlimited, Biologically Relevant Source

Presented by Brian Shapiro, Ph.D.

12 ET, June 4, 2020

Understanding COVID-19: A Global Pandemic

Presented by Britany Tang

12 ET, June 11, 2020

Viral Metagenomics and the Use of Standards: From Biology to Clinical Applications

Presented by Tasha M. Santiago-Rodriguez, Ph.D.

12 ET, June 25, 2020

© 2020 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.

