

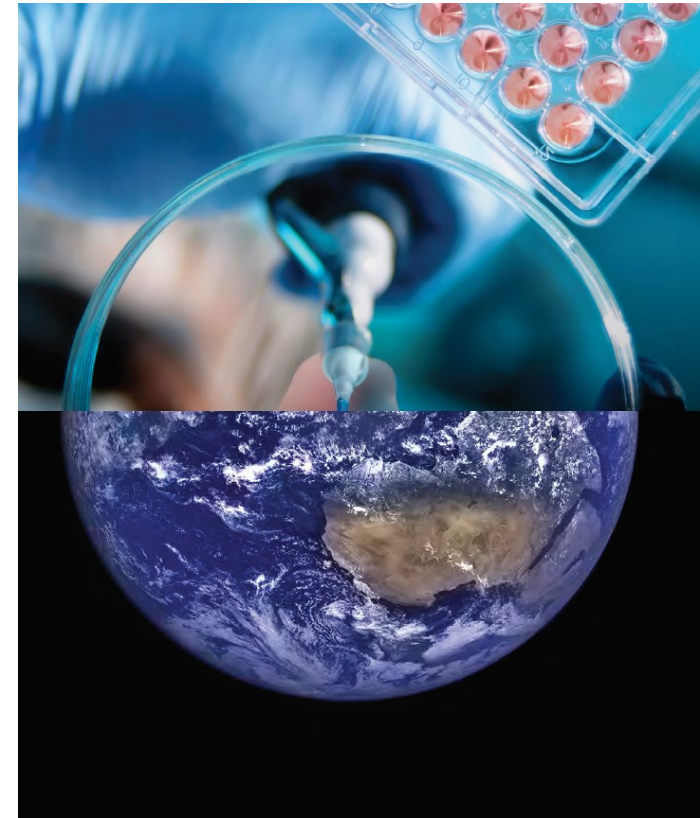
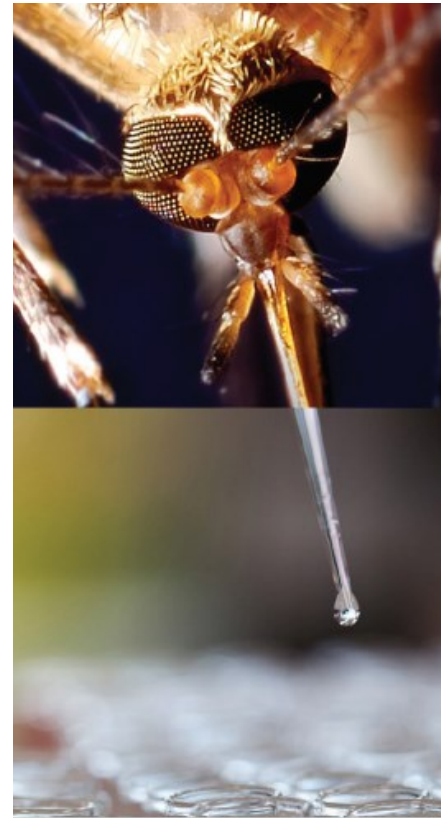
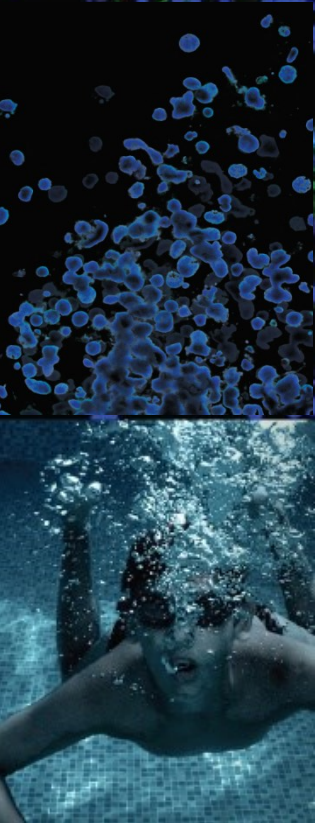


# Cell culture fundamentals: Your questions answered

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

Kevin Grady, BS  
*Senior Product Line Business Manager, ATCC*

Credible Leads to Incredible™



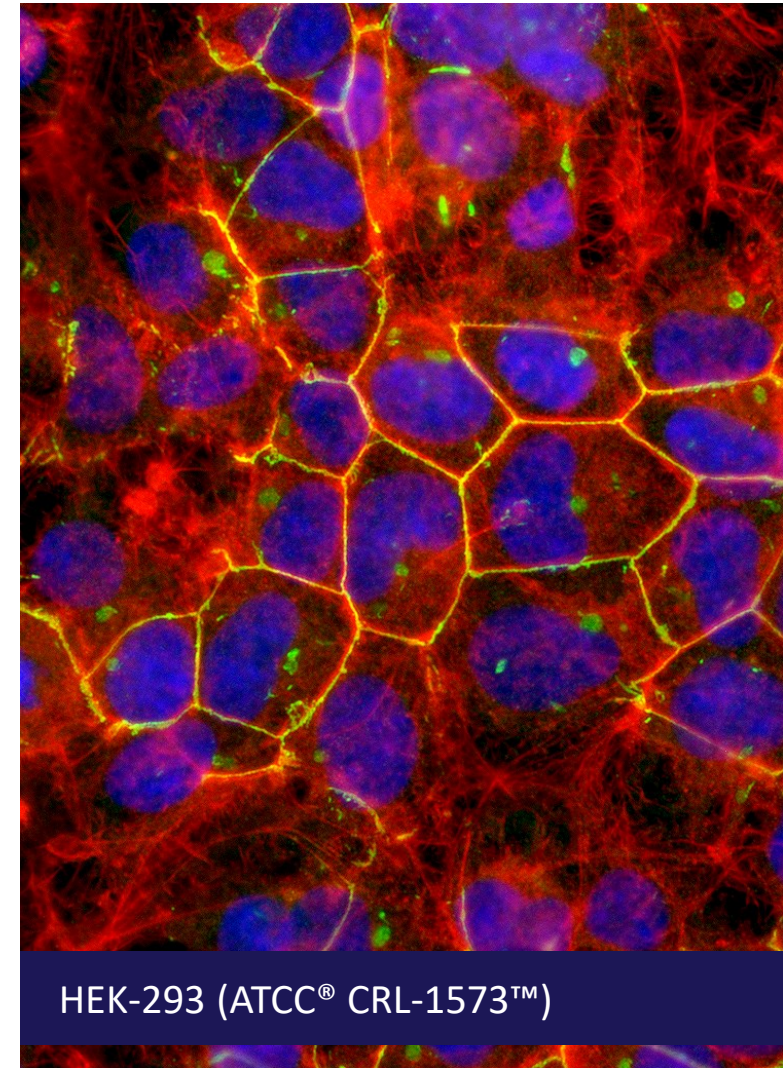
# ATCC – Credible leads to Incredible

- ATCC has provided credible biomaterials for over 90 years
- We continue to cultivate collaboration
  - Among scientists across disciplines
  - Essential for accelerating innovative research
  - Leading to incredible, high-impact results
- Our Cultivating Collaboration pledge: We bring scientists together to discuss
  - Breakthroughs in the state of science
  - Multidisciplinary approaches to key areas of research
  - Breaking the silos that impede research
- Our partnership with you, the scientific community, allows us all to reach the incredible

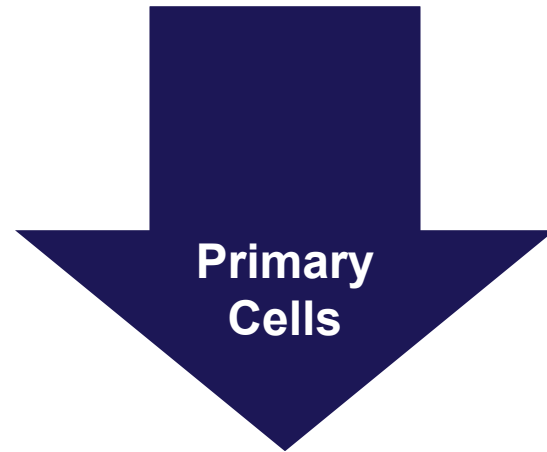


# Agenda

- Cell culture workflow - what cells to use
- Cell handling/media handling
- Viability assays
- Summary

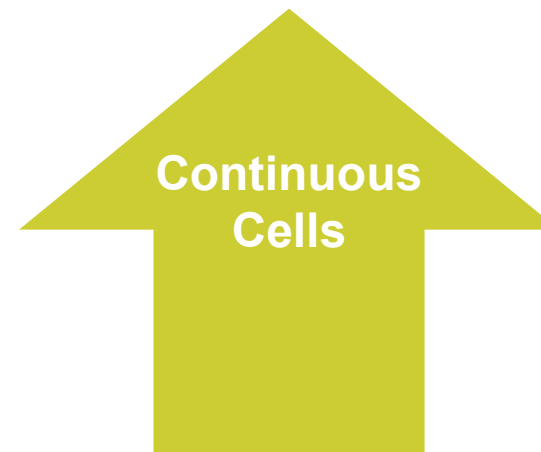


# Primary cells versus continuous cells



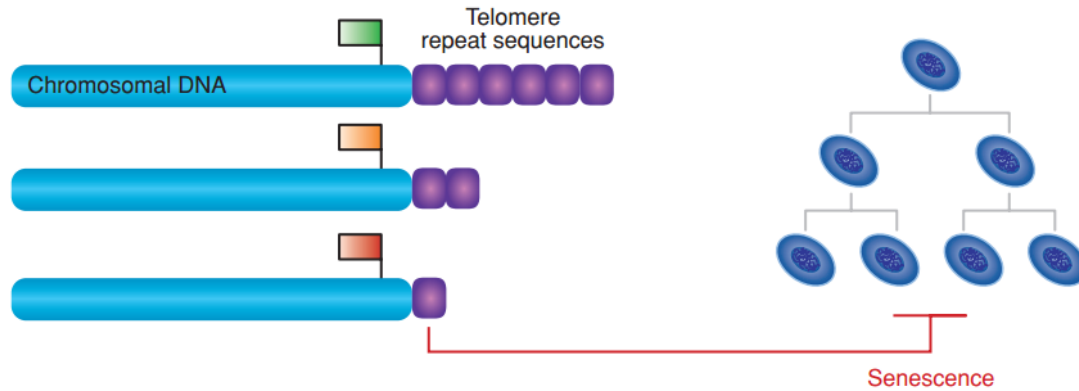
- Prepared directly from tissue
- Physiologically relevant
- Low risk for phenotypic or genotypic drift

- Easy to propagate in vitro
- Easy to generate large quantities of cells
- Inexpensive to maintain

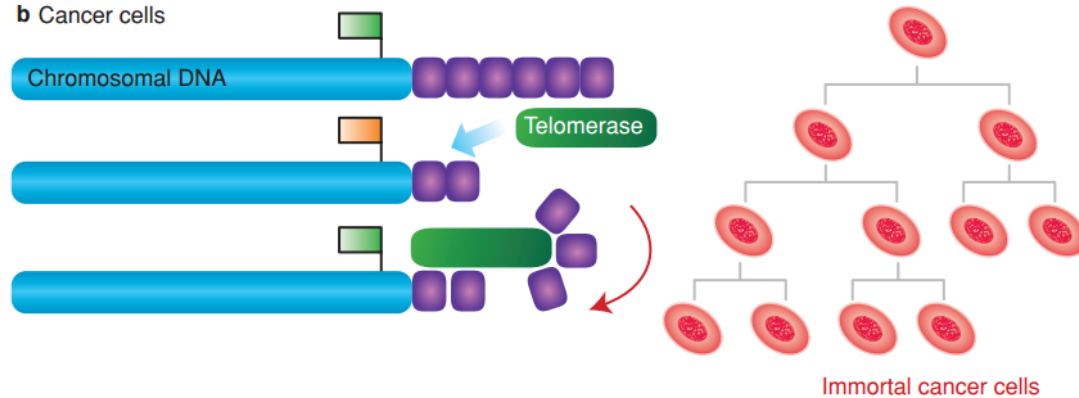


# hTERT-immortalized primary cells

a Normal somatic cells



b Cancer cells



- Bypass replicative senescence by telomerase
- Maintain primary cell function with the lifecycle of a continuous cell line

Regulation of telomere length in normal and cancer cells by telomerase

Expert Reviews in Molecular Medicine ©2002 Cambridge University Press

# Potential workflow situations: Standardization and validation

Use cell lines for standardization and confirmation of each experiment

- Large number of cells needed
- Samples with limited variability
- Generally easy to manipulate

Use primary cells after standardization to further validate the results

- Donor variability
- Biological relevance



# Potential workflow situations: High-throughput screening

## Screening work flow

- Initial screening in cell lines
  - Large number of cells needed
  - Samples with limited variability
- Next level in hTERT-immortalized primary cell lines
  - Large number of cells needed
  - Samples with limited variability
  - More physiologically relevant results
- Final screen in primary cells (**Results with the most biological relevance**)

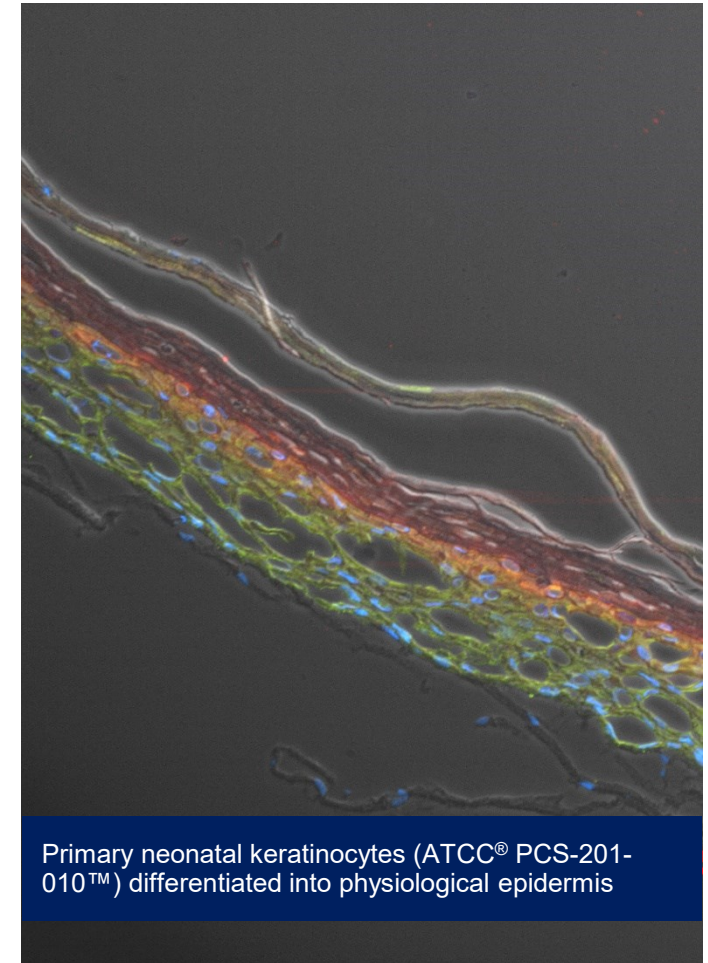


# Primary cells as a control

**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**







# 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**

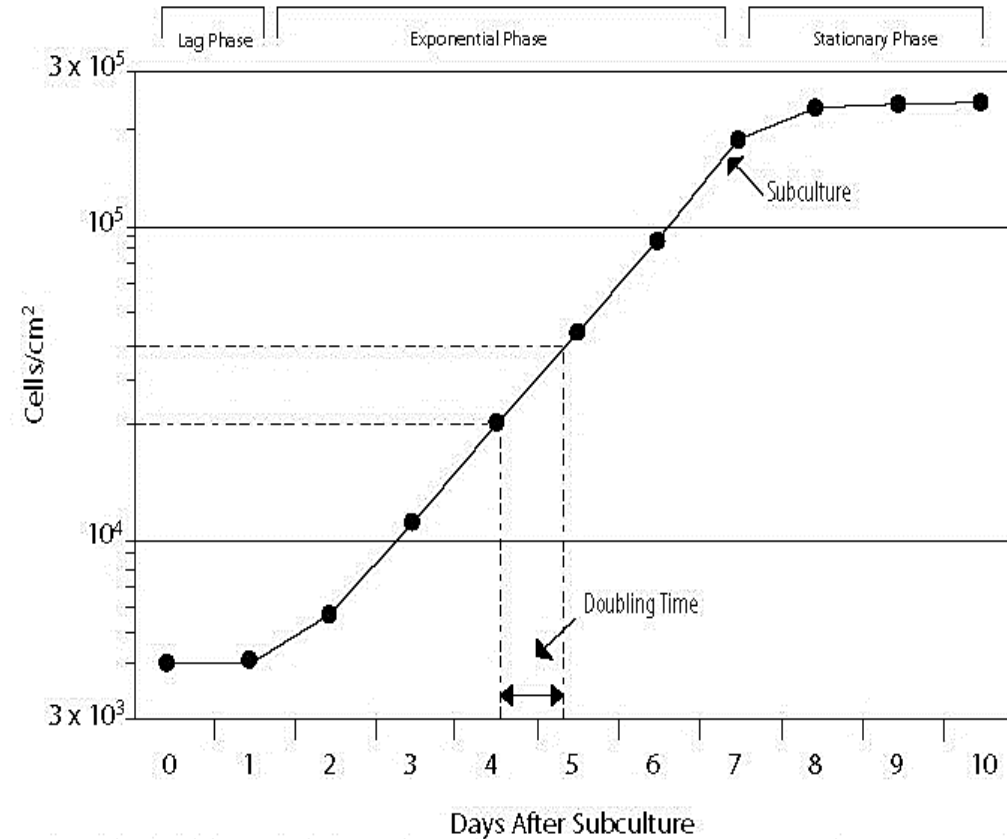


# Cell expansion

- 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



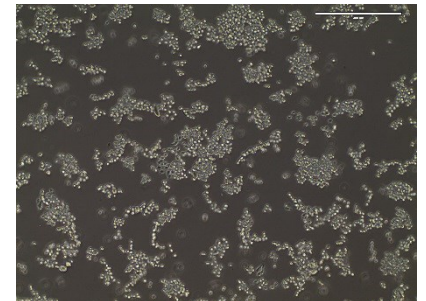
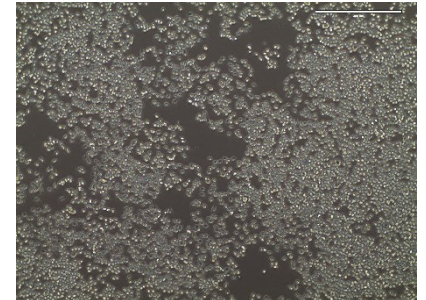
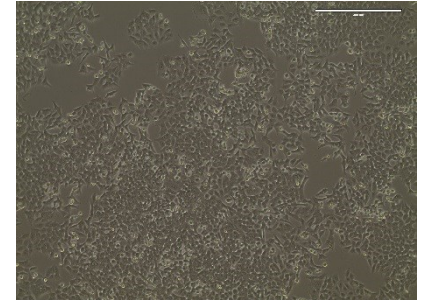
**Figure 1.** Growth curve for cells grown in culture. Cells should be subcultured while still in the exponential phase.

# Trypsinization

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

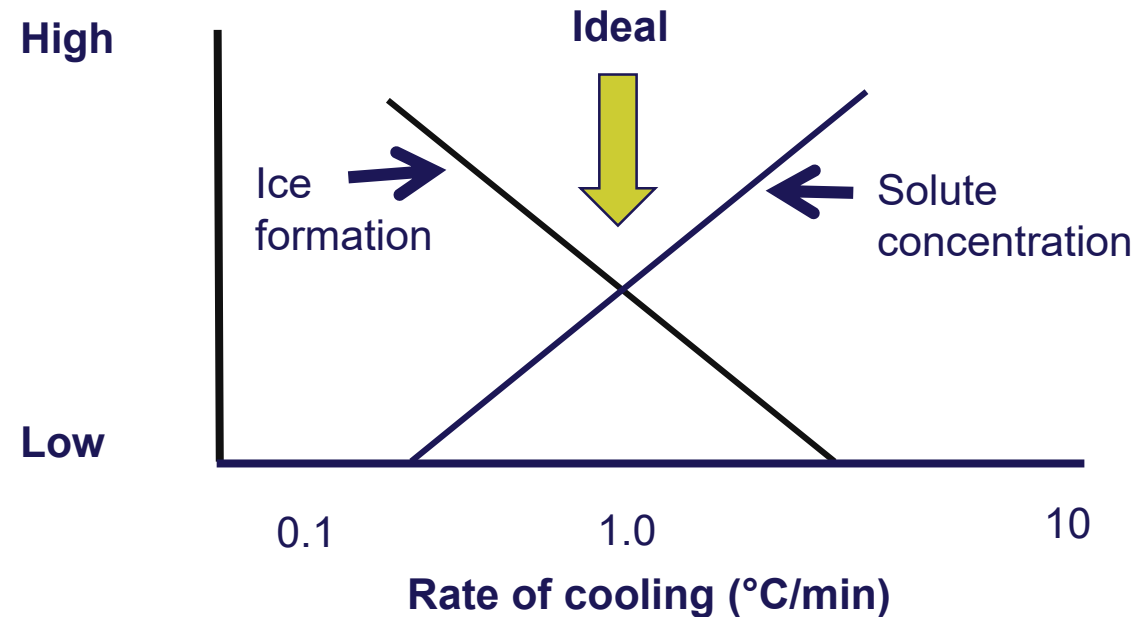
- Using warm trypsin-EDTA for about 3-5 minutes, cells will detach with gentle agitation
- ***Trypsin-EDTA for Primary Cells (ATCC® PCS-999-003™) is a low concentration formula (.05% Trypsin and .002% EDTA) – necessary for primary cell survival***
- A Trypsin Soybean Neutralizing Solution (ATCC® 30-2104™)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

Monolayer



Fully trypsinized

# 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

-70°C

Controlled-rate freeze chamber

-1°C/min cooling rate

A few hours to 24 hours



-140°C

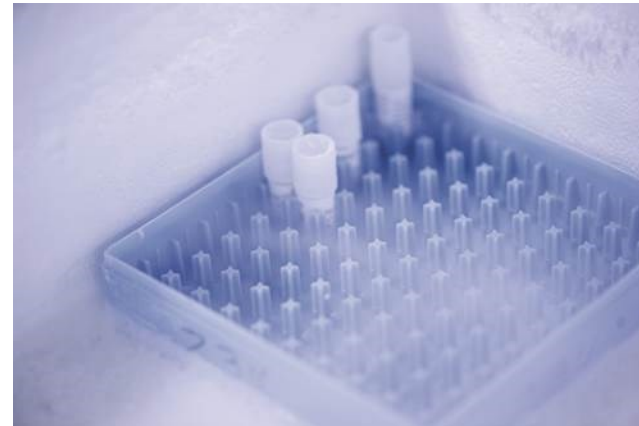
Liquid nitrogen tank



# Low temperature storage



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

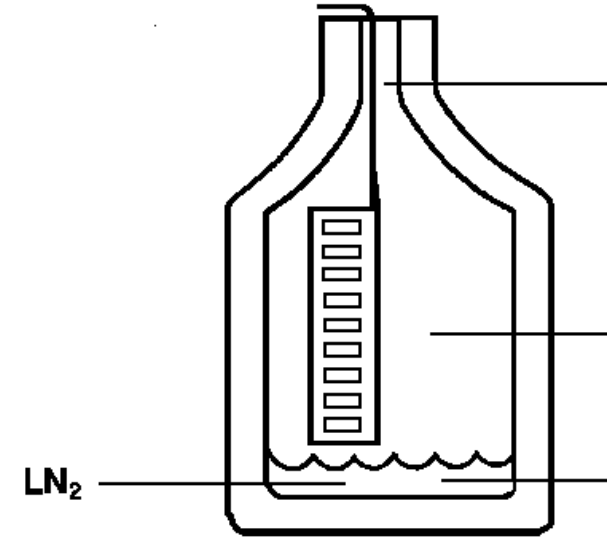




# Low temperature storage

## Mammalian cells

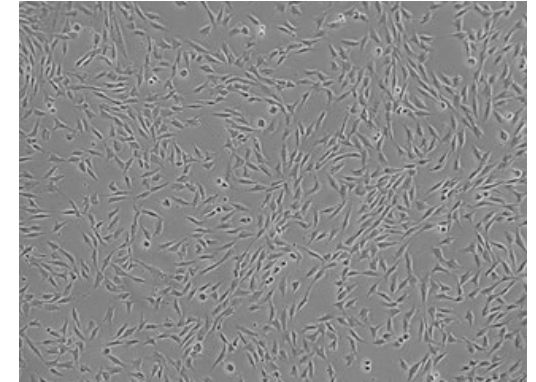
- Long-term storage should be below  $-140^{\circ}\text{C}$
- Vials should be stored in a liquid nitrogen unit **above** the volume of liquid at the bottom of the tank
- This temperature should be between  **$-140^{\circ}\text{C}$**  and  **$-180^{\circ}\text{C}$**



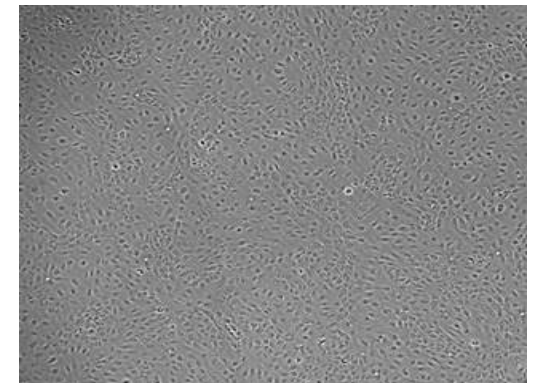
# Cell characterization

## 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



Fibroblasts



HUVEC

# Contamination

## Sources

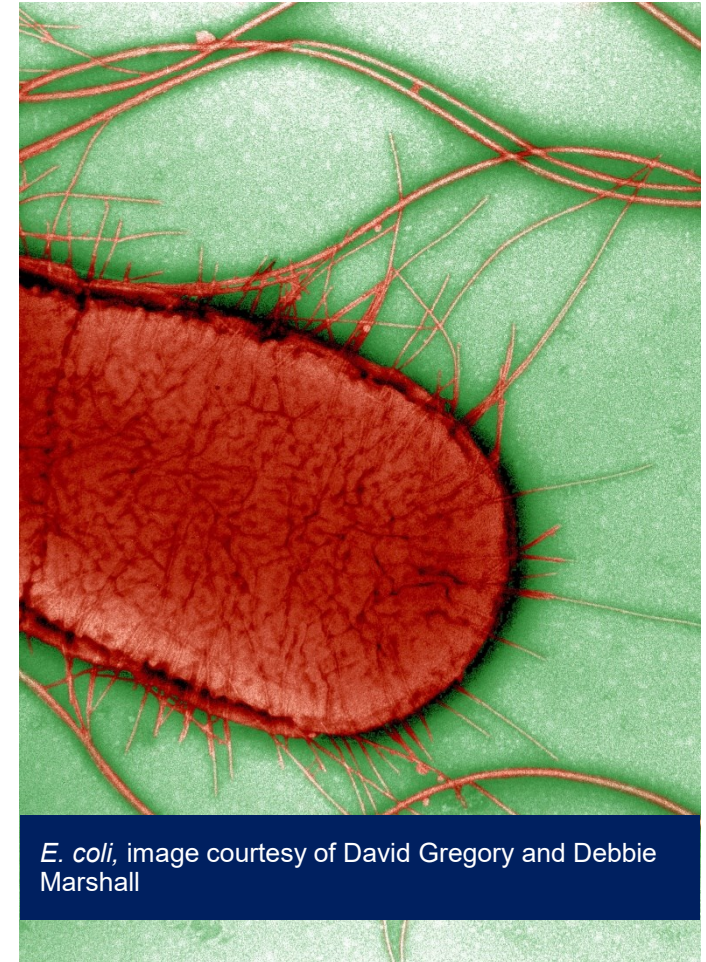
- Contaminated cell lines
- Improper aseptic technique

## Types

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

## Signs

- Turbid media
- Rapid decline in pH – color change
- Morphological changes
- Filamentous structures



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

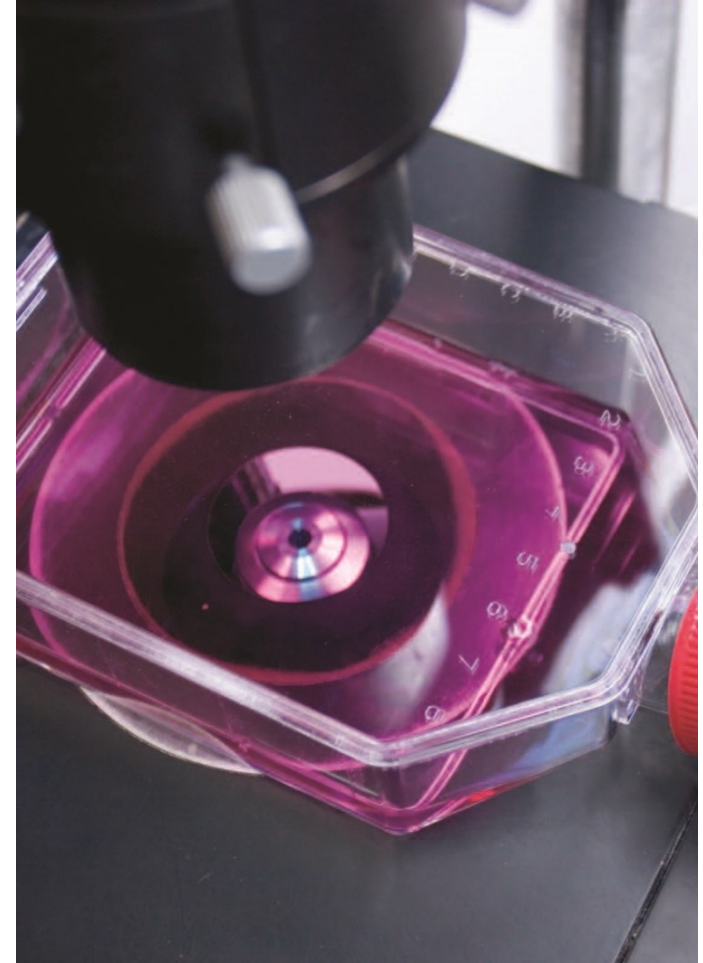
# Mycoplasma contamination

## 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



# Contamination

## Cross Contamination

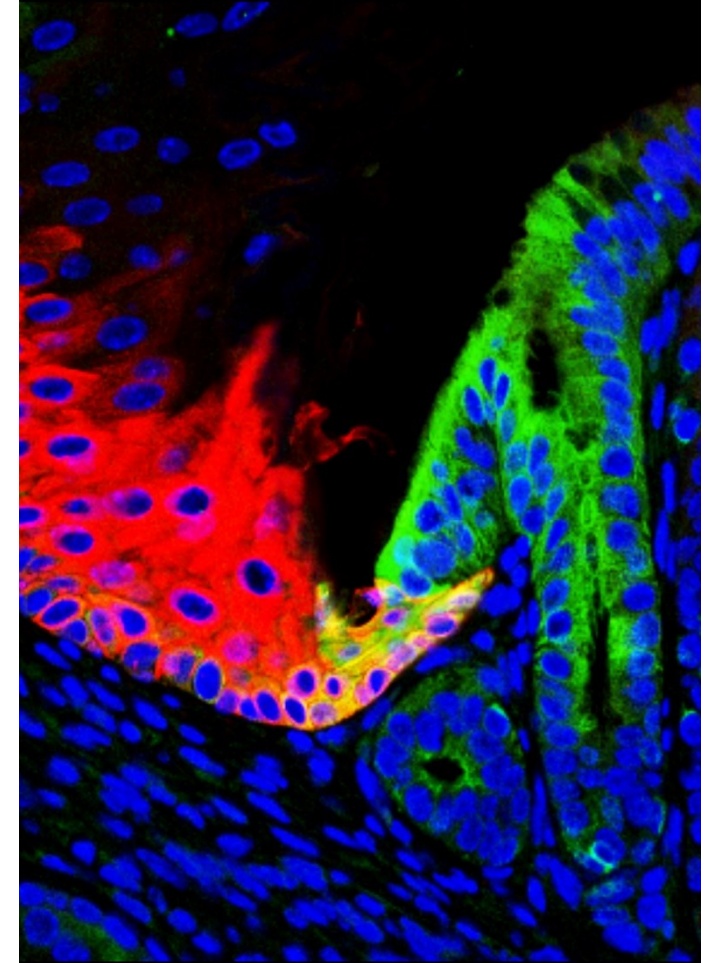
**Leads to the replacement of the original cell line with the contaminant**

### 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**



# Cell characterization

## Universal Mycoplasma Detection Kit

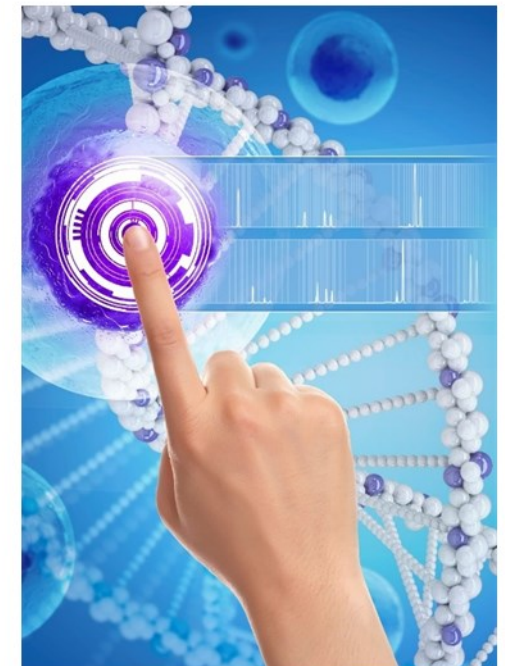
PCR-based kit (ATCC® 30-1012K™)

Detects any of the 60 most common mycoplasmas

## ATCC STR Profiling (Human and Mouse)

Ensures your cells are what you think they are

- STR profile of your cell line
- Comparison of your cells against ATCC STR Profile database at [www.atcc.org/str](http://www.atcc.org/str)
- Electropherograms supporting the allele calls at each locus
- Comprehensive interpretation of results



# Contamination

## Personnel and equipment

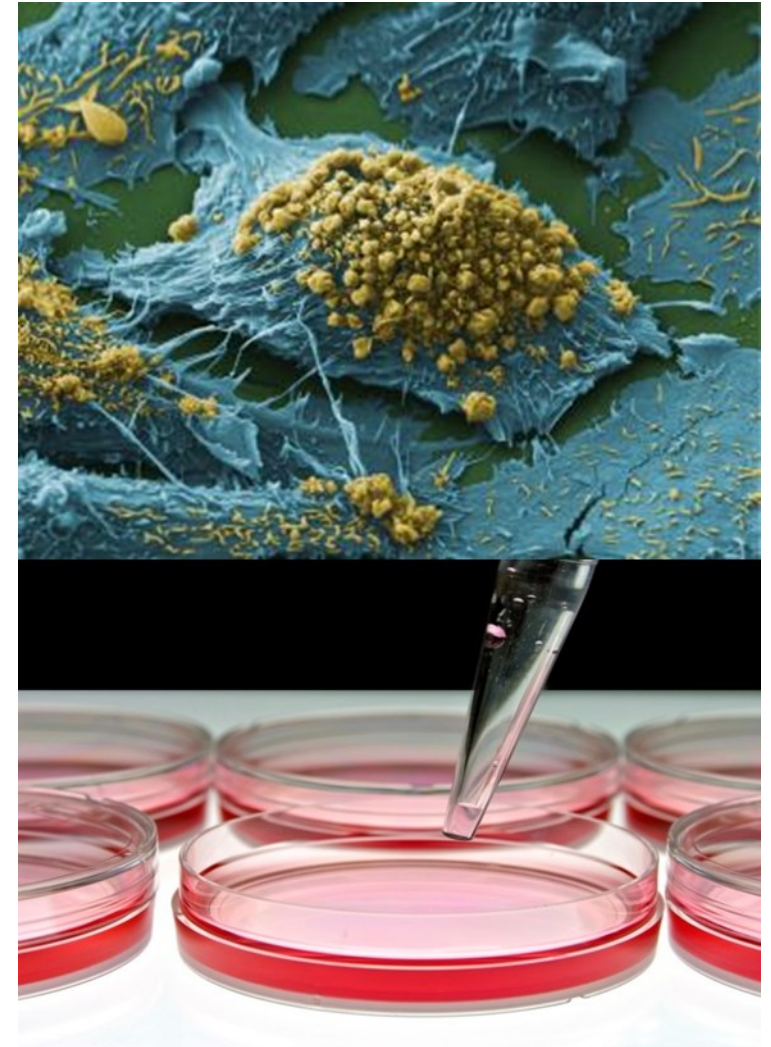
- Poor culturing practices
- Dust and aerosol

## 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



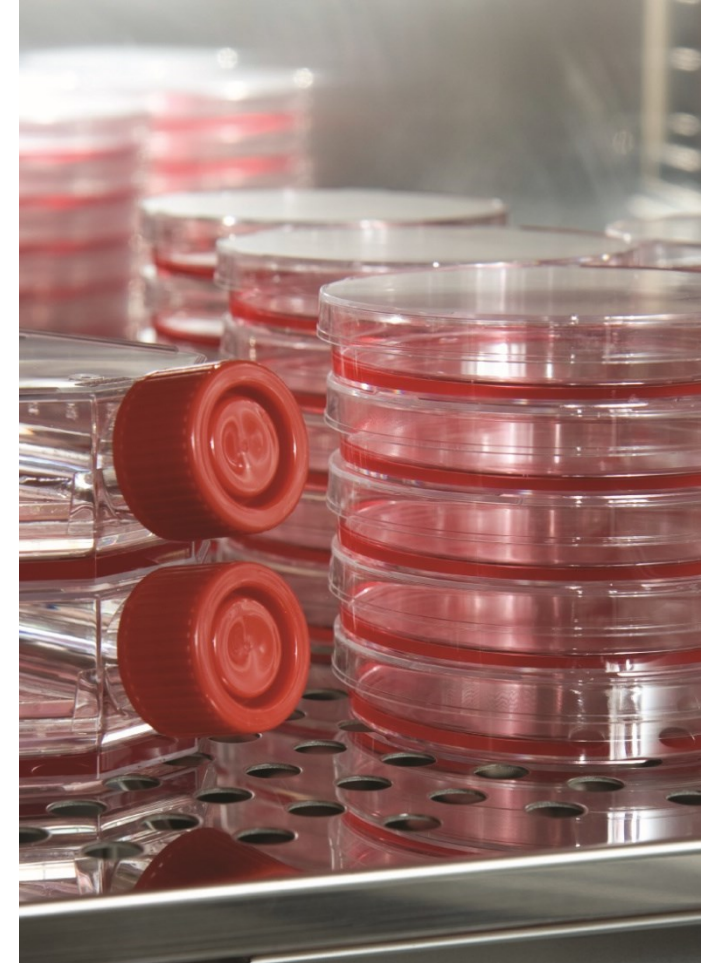
# Media choices

## Animal cell lines – media + 10% FBS

- Eagle's Minimum Essential Medium (EMEM; ATCC® 30-2003™)
- Dulbecco's Modified Eagle's Medium (DMEM; ATCC® 30-2002™)
- Iscove's Modified Dulbecco's Medium (IMDM; ATCC® 30-2005™)
- Kaighn's Modification of Ham's F-12 Medium (ATCC® 30-2004™)
- DMEM/ F12 Medium (ATCC® 30-2006™)
- McCoy's 5A (ATCC® 30-2007™)
- RPMI-1640 (ATCC® 30-2001™)
- Leibovitz's L-15 (ATCC® 30-2008™)

## Primary Cells – Primary Cell Basal Media and Growth Kits

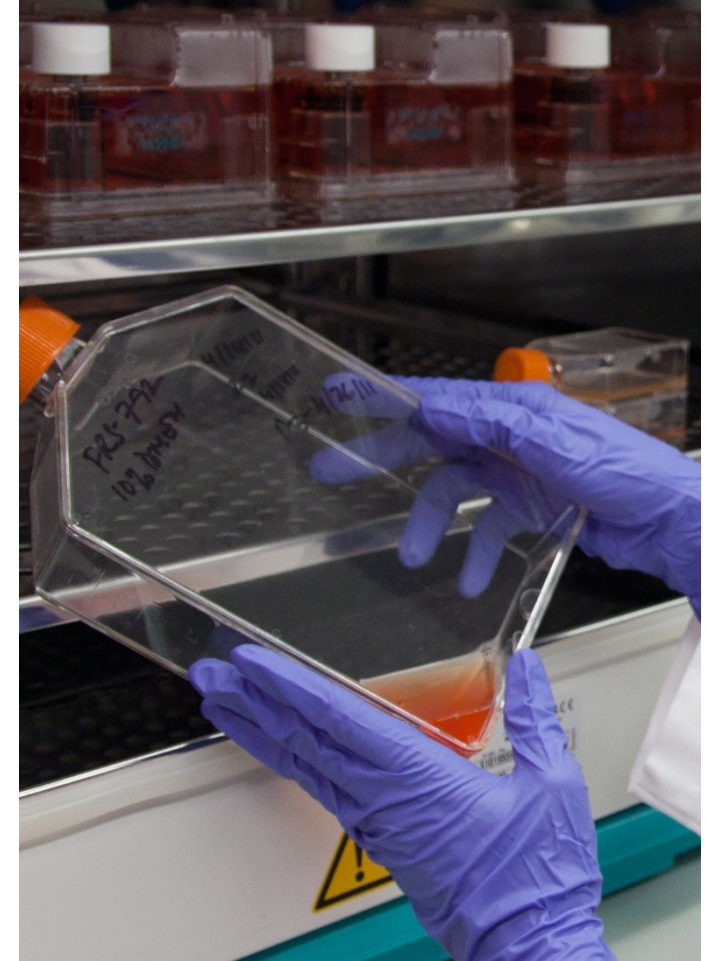
- Primary cells require their own specially formulated media, specific to each cell type

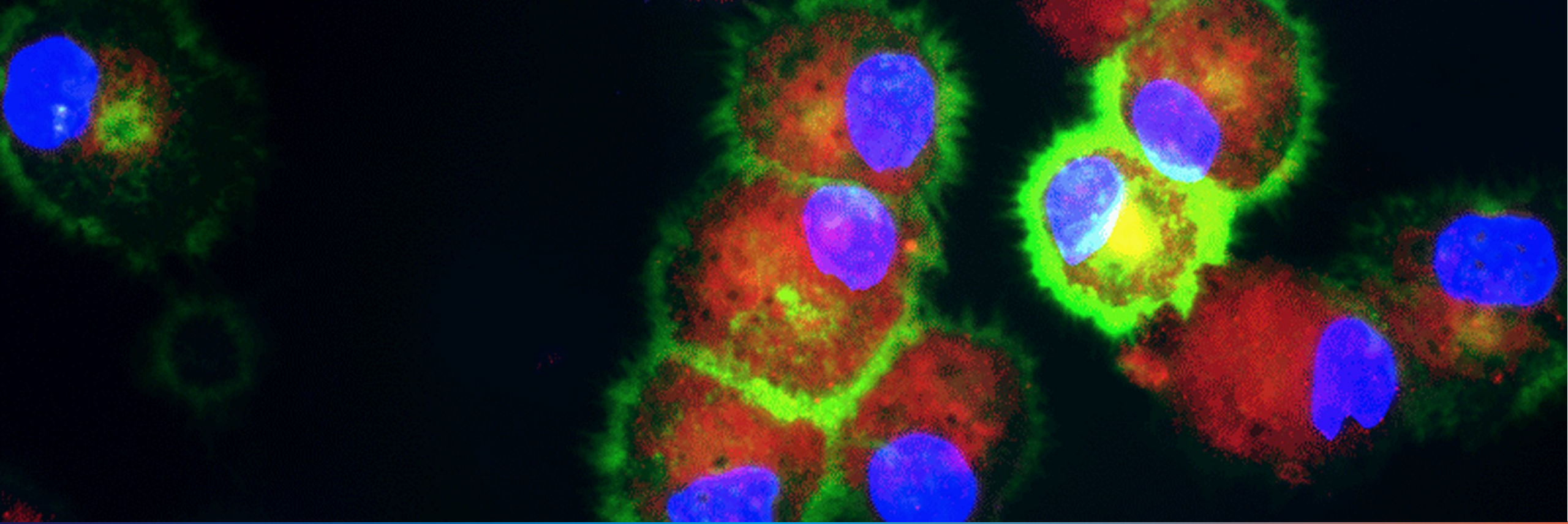


# Media choices

## Special notes:

- Maintain cells in the same media
- Vendor to vendor media variability
  - Possible osmotic shock
- When transferring 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***





# Viability assays

# Viability assays

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

- Commonly used for cytotoxicity, 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, 5-diphenyltetrazolium bromide)

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

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

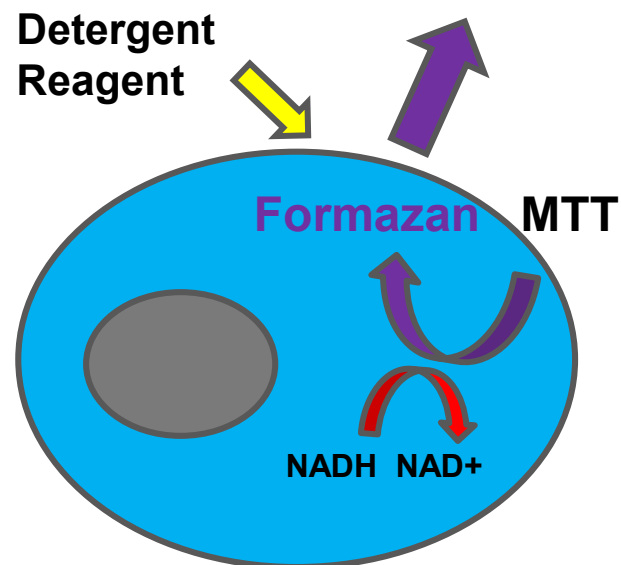


# Viability assays

## MTT Reaction

MTT salt is **reduced** within cellular matrix to Formazan, lysed with detergent to solubilize crystals

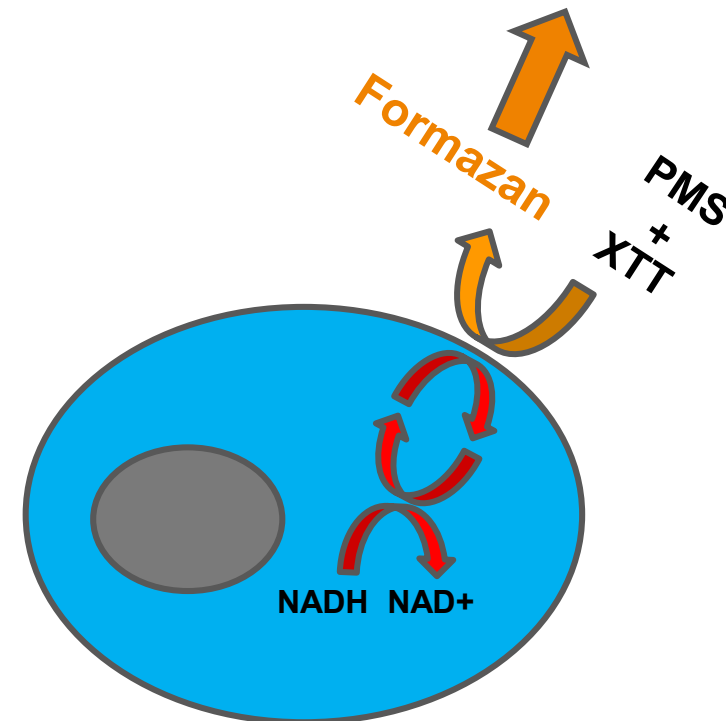
Media turns **PURPLE**



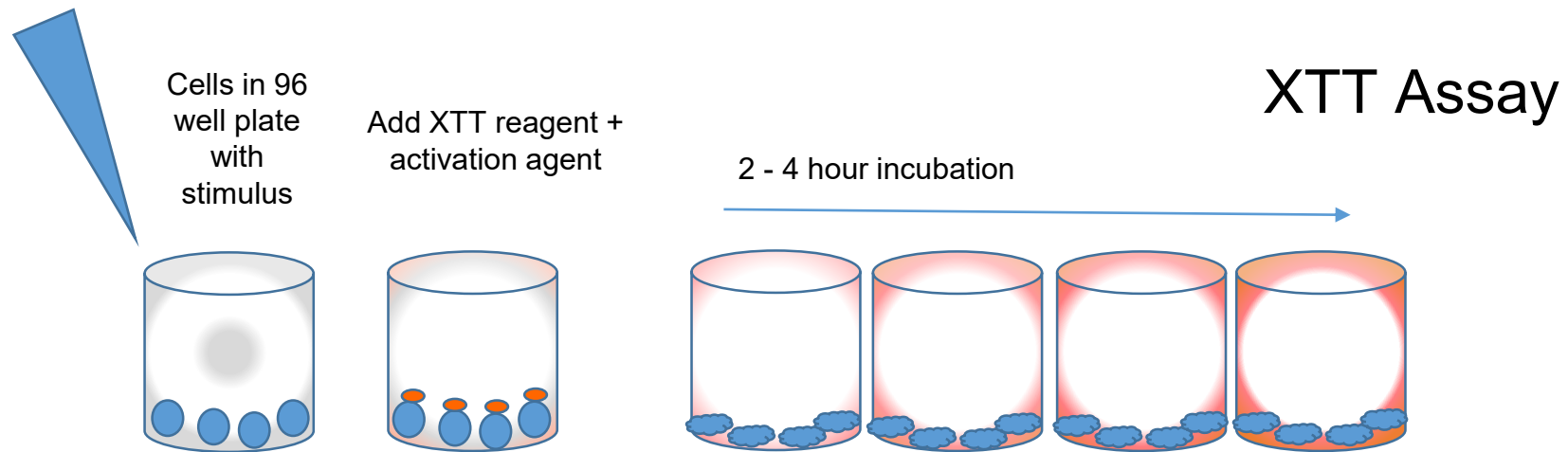
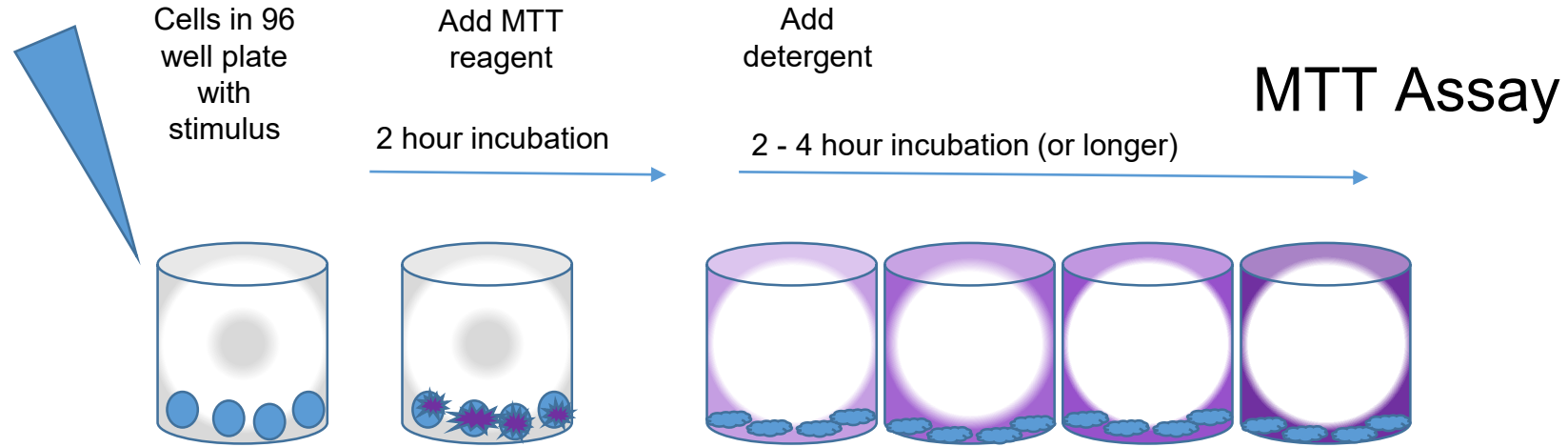
## XTT Reaction

XTT salt is **reduced** at cell membrane with PMS agent

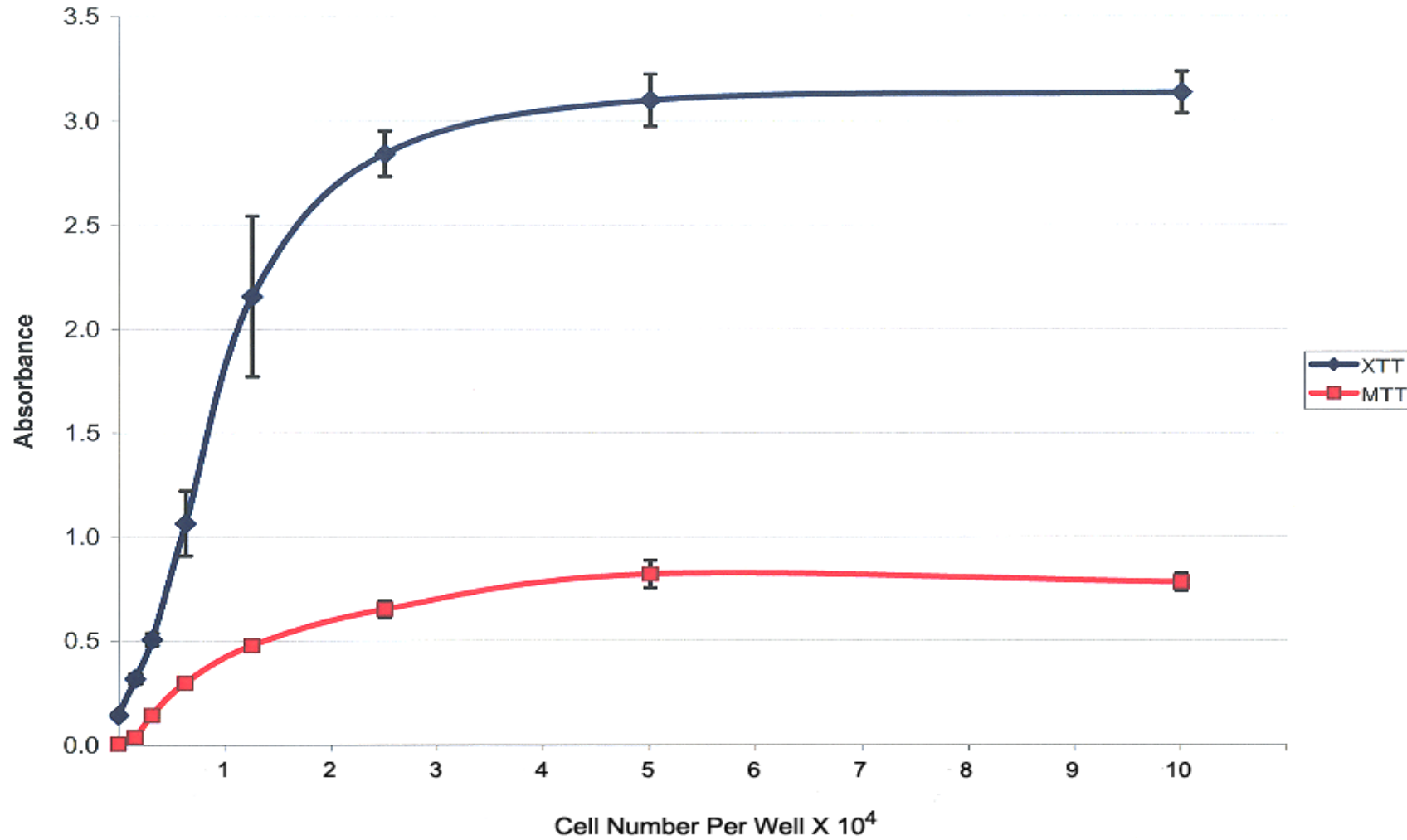
Media turns **ORANGE**



# Viability assays



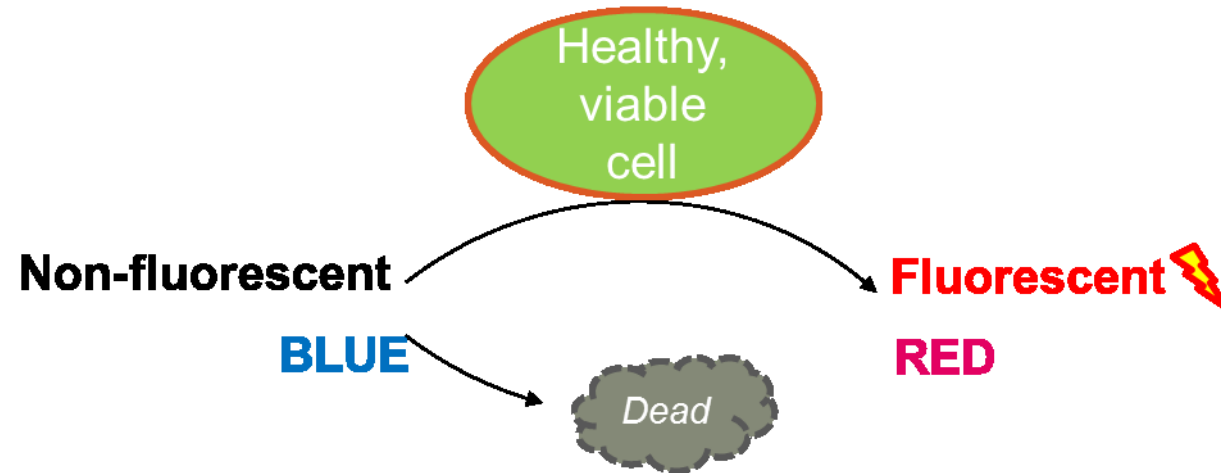
# Viability assays



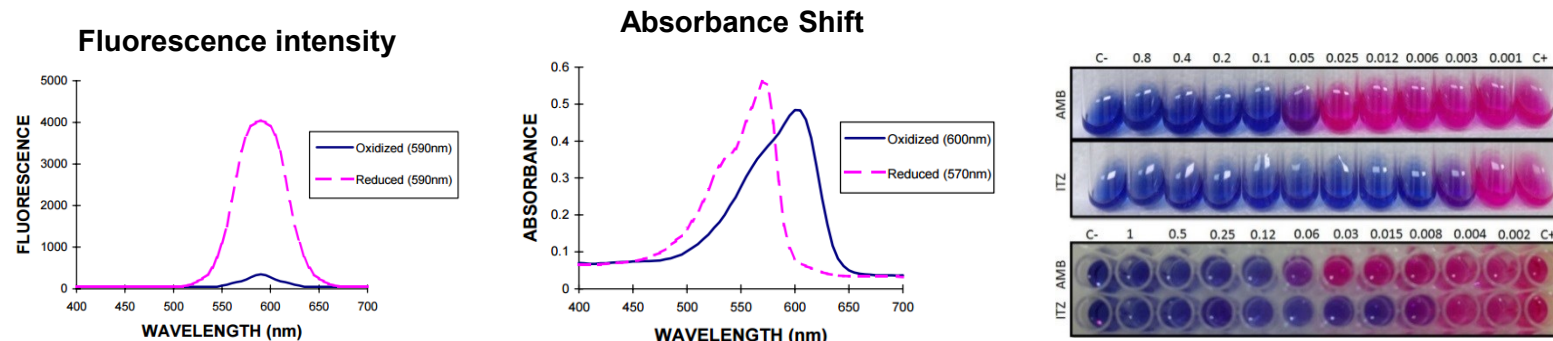


# Reliablue™ Cell Viability Reagent

Resazurin (7-Hydroxy-3H-phenoxazin-3-one 10-oxide) is a blue dye that is weakly fluorescent until reduced (redox) at which point it becomes pink and highly fluorescent.



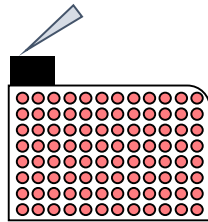
Resazurin is cell permeable but non-toxic and is metabolically reduced by living cells (but not dead cells or in the culture media) resulting in a change in absorbance and increase in fluorescence.



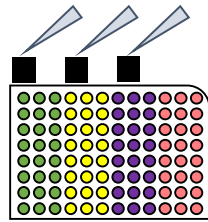
# Reliablue™ Cell Viability Reagent

Reliablue™ Reagent is supplied in a 10X ready-to-use format that can be added directly to cells, typically in multiwell plates. An overview of the workflow is shown below.

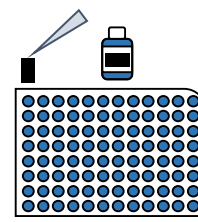
## Basic 4-Step Assay Workflow



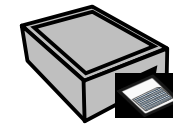
① Grow cells



② Add treatments



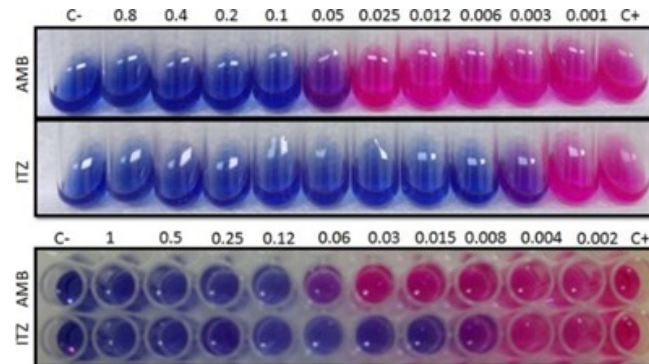
③ Add Reliablue



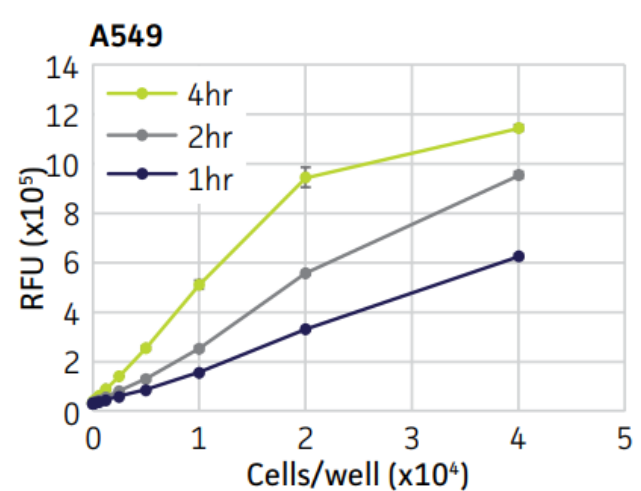
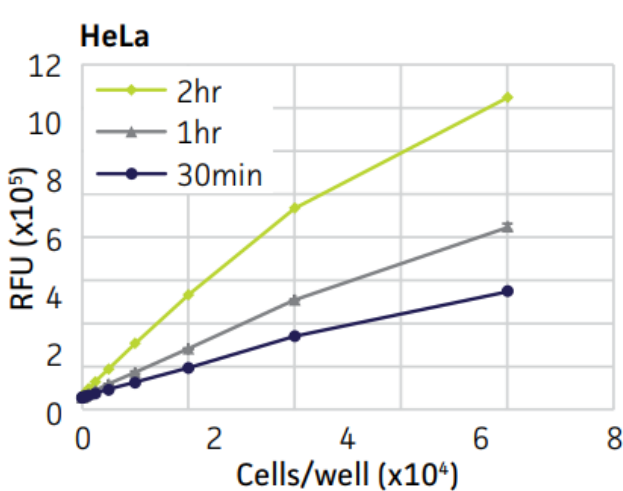
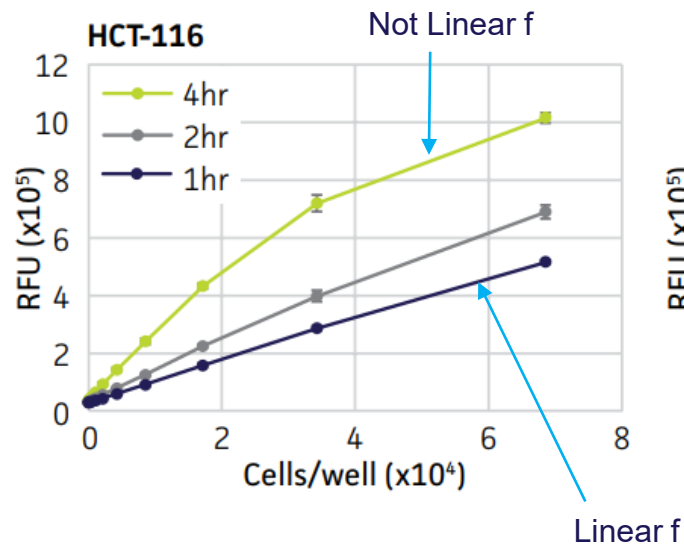
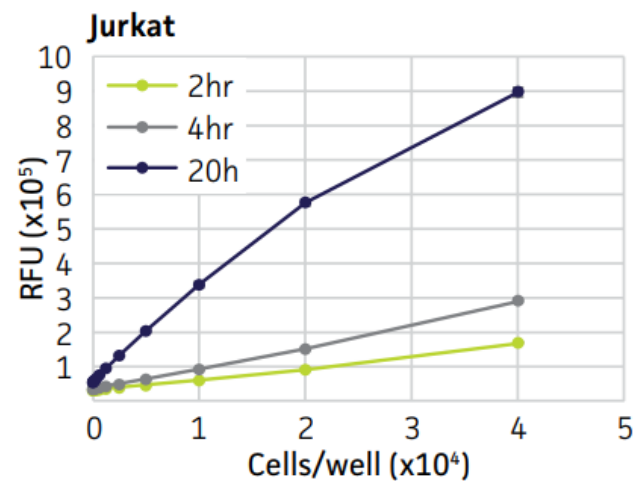
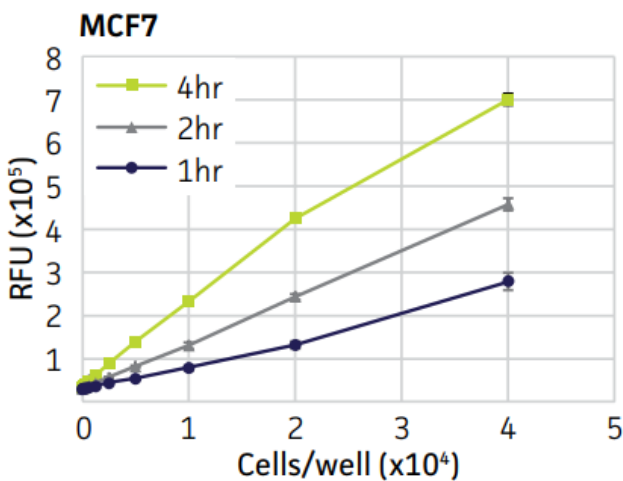
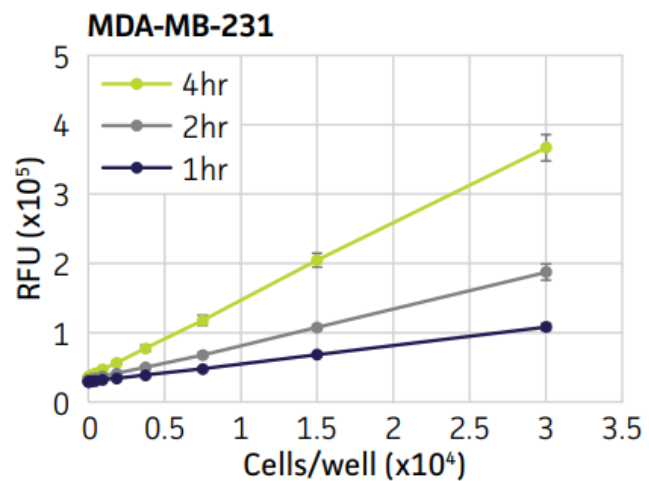
④ Read on plate-reader

Blue = Dead (or control)

Red = Alive and Actively Growing



# Reliablue™-derived viability curves





# Summary

# Summary

## Cell culture workflow

- Use cell lines for standardization and confirmation of each experiment
- Use primary cells after standardization to further validate the results
- hTERT-immortalized cells are the best of both worlds

## Cell Handling/media handling

- Be sure to employ best practices to eliminate contamination and ensure optimal growth and storage

## Viability assays

- MTT, XTT, and Reliablue can confirm cell growth characteristics



# Cultivating collaboration to support global health

## Upcoming webinars:

- Tick-borne Diseases: Developing Molecular Approaches to Detect Multiple Pathogens | August 15, 12:00 ET
- Making Sense Out of Microbiome Data – The Importance of Standards | August 22, 12:00 ET

[www.atcc.org/webinars](http://www.atcc.org/webinars)

