

# Cell Health and Viability

Steven Budd, M.S., M.B.A.  
*Product Line Business Specialist, 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 one-third with advanced degrees



Established partner to global researchers and scientists



# Outline

## Cell Health and Viability Topics

- Cell Culture
  - Media
  - Additives / Serum
- Cryopreservation / Post Thaw
- Cell Proliferation / Viability
  - MTT / XTT kits
  - Reliablue™ cell viability reagent
- Mycoplasma Effects / Detection



# Complete growth media

- Classical cell culture media
- Media ingredients
- Additives
- Animal sera



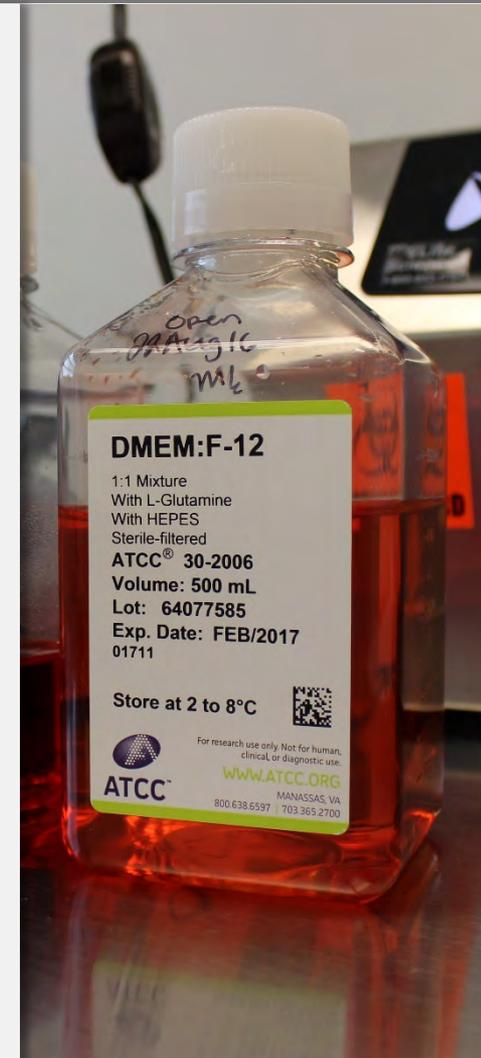
# Media choices

## 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 ingredients

## Sodium bicarbonate



## HEPES buffer

- Can buffer without CO<sub>2</sub> enrichment
- Good for working under the hood

## Phenol Red

- Monitors pH of media
- Yellow = acidic
- Purple = basic
- May mimic action of steroid hormones.

## Sodium Pyruvate

- Helps maintain metabolism



# Additives

## Nonessential Amino Acids

- Can be added to reduce the metabolic burden on cells

## L-Glutamine (ATCC® 30-2214™)

- Present in ATCC classical cell culture media
- Relatively stable in bottles kept at 4°C - 8°C
- **Glutamine degradation increases ammonia toxicity**
- **Generally not recommended to “spike” media with L-Glutamine**

## Antibiotics and Antimycotics

- Penicillin-Streptomycin, Gentamicin Sulfate
- Amphotericin B
- **Generally not recommended**



# Animal sera



## Fetal Bovine Serum (ATCC® 30-2020™)

### Fetal Bovine Serum, Embryonic Stem Cell Qualified (ATCC® SCRR-30-2020™)

- Very rich in growth factors, most common choice
- **Heat inactivation: Not Advised**

## Calf Bovine Serum (ATCC® 30-2030™)

- Lower concentrations of growth factors, good for contact inhibition studies

## Horse Serum (ATCC® 30-2040™)

- Collected from closed herds, lot-to-lot consistency, no bovine viruses

# Media usage considerations

**Maintain cells in the same media**

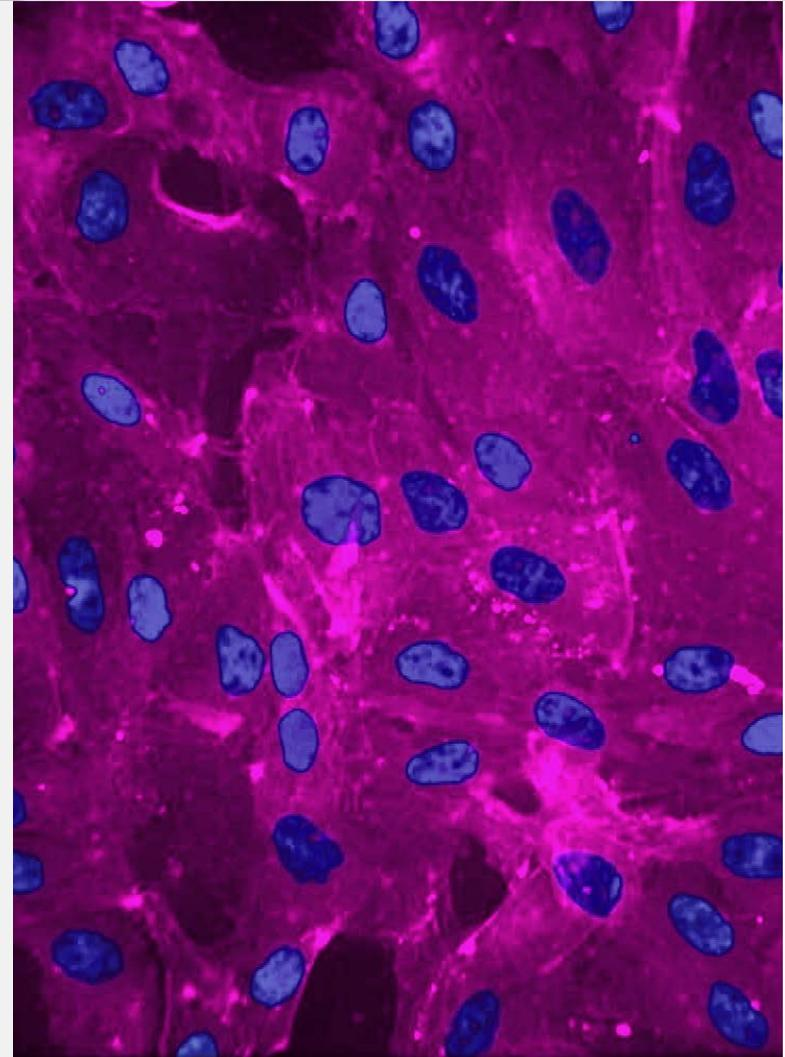
**Media variability**

- 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

**ATCC® Animal Cell Culture Guide**



# Outline

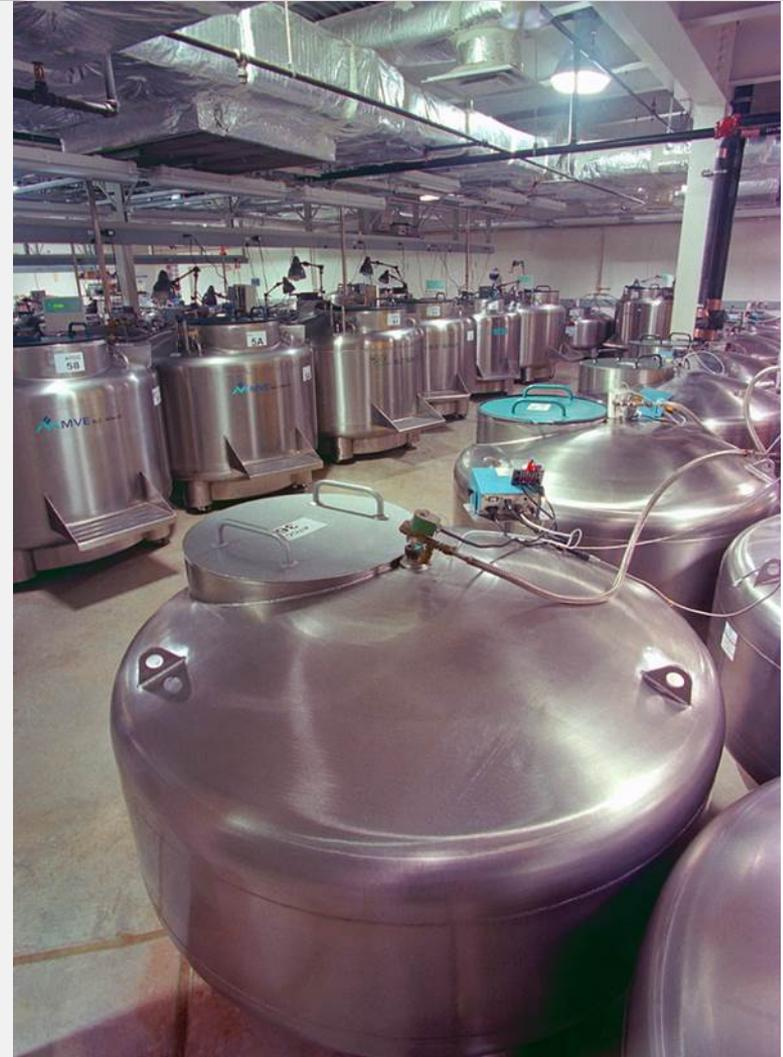
## Cell Health and Viability Topics

- Cell Culture
  - Media
  - Additives / Serum
- **Cryopreservation / Post Thaw**
- Cell Proliferation / Viability
  - MTT / XTT kits
  - Reliablue™ cell viability reagent
- Mycoplasma Effects / Detection

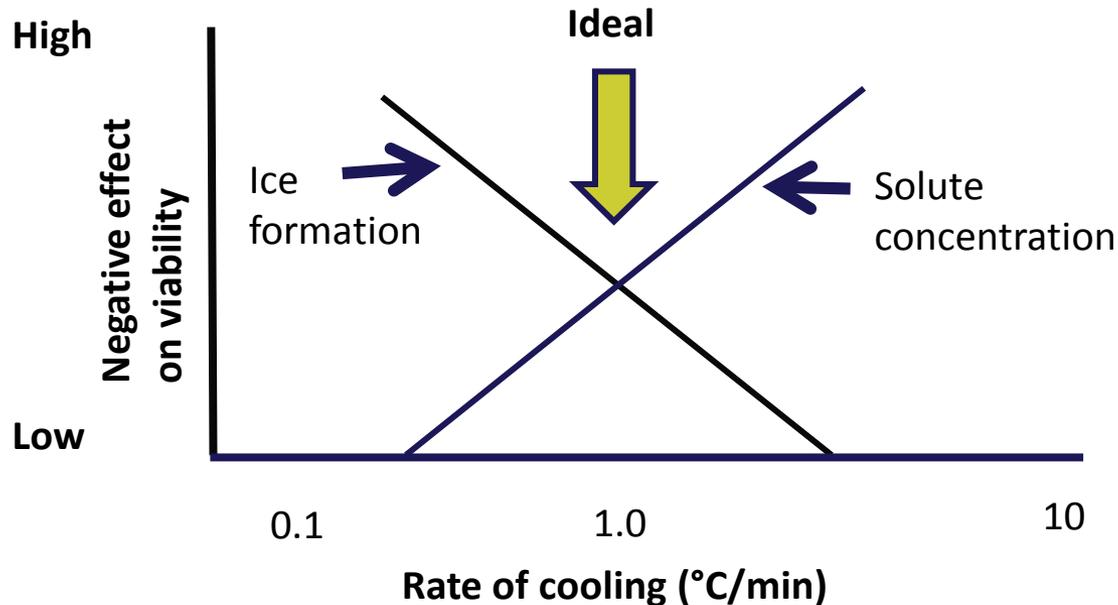


# Cryopreservation procedure

- Overview
- Cryoprotectants and media preparation
- Freezing cells in a controlled-rate chamber
- Long-term storage
- Thawing / Plating / Trypsinization



# Cryopreservation principles



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

# Cryoprotectants

Cell type	Cryoprotectant	Temperature	Number of cells
Animal cells	DMSO (5-10%) or Glycerol (5-10%)	-140°C	10 <sup>6</sup> to 10 <sup>7</sup> /mL
Bacteria	Glycerol (5-10%)	-80°C	10 <sup>7</sup> /mL
Yeast	Glycerol (10%)	-140°C	10 <sup>7</sup> /mL
Protozoa	DMSO (5-10%) or Glycerol (10-20%)	-140°C	10 <sup>5</sup> to 10 <sup>7</sup> /mL
Plant cells	DMSO (5-10%) and Glycerol (5-10%)	-140°C	3% to 20% cell volume
Animal viruses (free)	None	-80°C	NA
Animal viruses (infected cells)	DMSO (7%)	-10°C	10 <sup>6</sup> /mL

# Media preparation

## Classical cell culture media – DMEM, EMEM, RPMI-1640 (for suspension cells)

- 5-10% DMSO
- 20% fetal bovine serum (FBS) or bovine serum albumin (BSA)
  - Additional cryoprotectant properties
  - Necessary for post-thaw cell survival

## ATCC Serum-free Freezing Media (ATCC® 30-2600™)

- All in one media
- 10% DMSO with proteins and additives for cell survival



# Freezing cells

-70°C

Controlled rate freeze chamber

-1°C/min cooling rate

A few hours to 24 hours

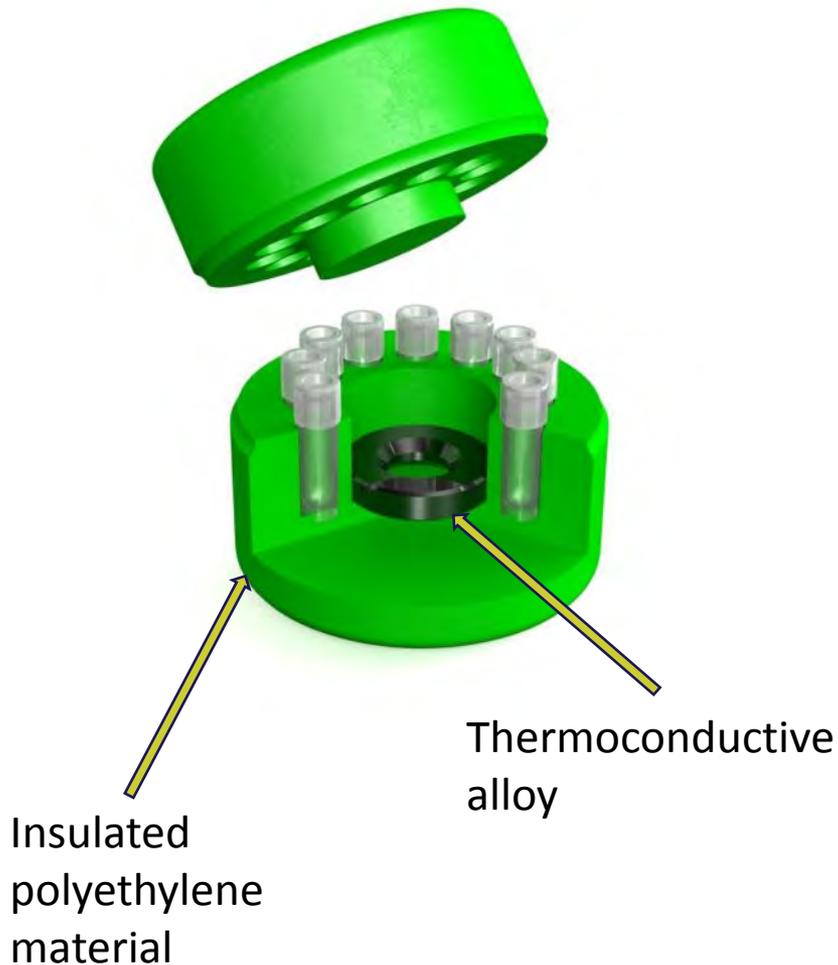


-140°C

Liquid nitrogen tank



# Freezing cells



## CoolCell® LX (ATCC® ACS-6000™)

- Reliable  $-1^{\circ}\text{C}/\text{min}$  cooling rate
- 4 hours in  $-70^{\circ}\text{C}$  freezer
- Comfortable to touch
- No alcohol use or maintenance
- Can be used with all cell types
  - Verified use with organoids

# Low temperature storage

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



# Low temperature storage

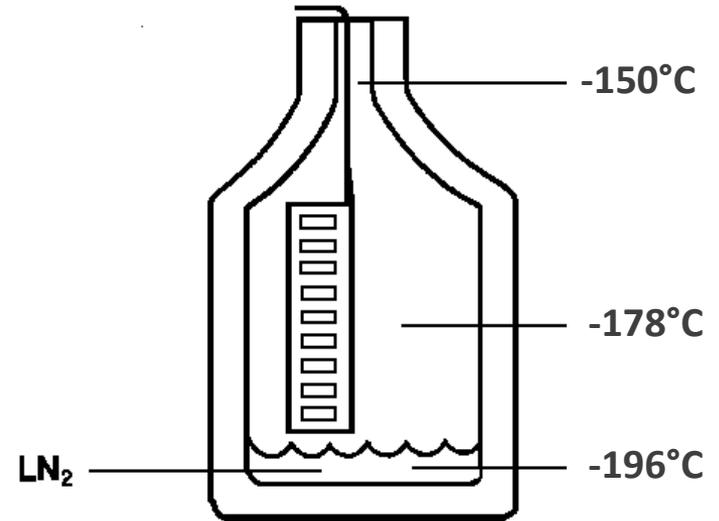
## Mammalian cells

Long-term storage should be below  $-140^{\circ}\text{C}$

- $-140^{\circ}\text{C}$  for an indefinite length of time
- $-80^{\circ}\text{C}$  for less than 1 year

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  $-180^{\circ}\text{C}$



# 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**

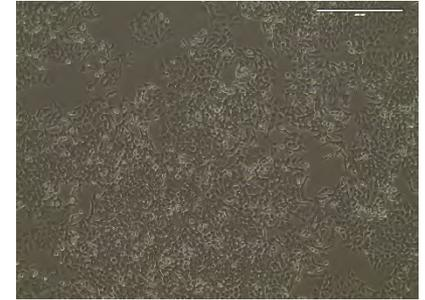


# Trypsinization

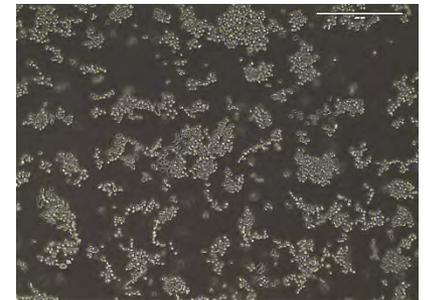
At 80% confluency (primary cells), cells can be passed using Trypsin-EDTA

- 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



# Outline

## Cell Health and Viability Topics

- Cell Culture
  - Media
  - Additives / Serum
- Cryopreservation / Post Thaw
- **Cell Proliferation / Viability**
  - MTT / XTT kits
  - **Reliablue™ cell viability reagent**
- Mycoplasma Effects / Detection

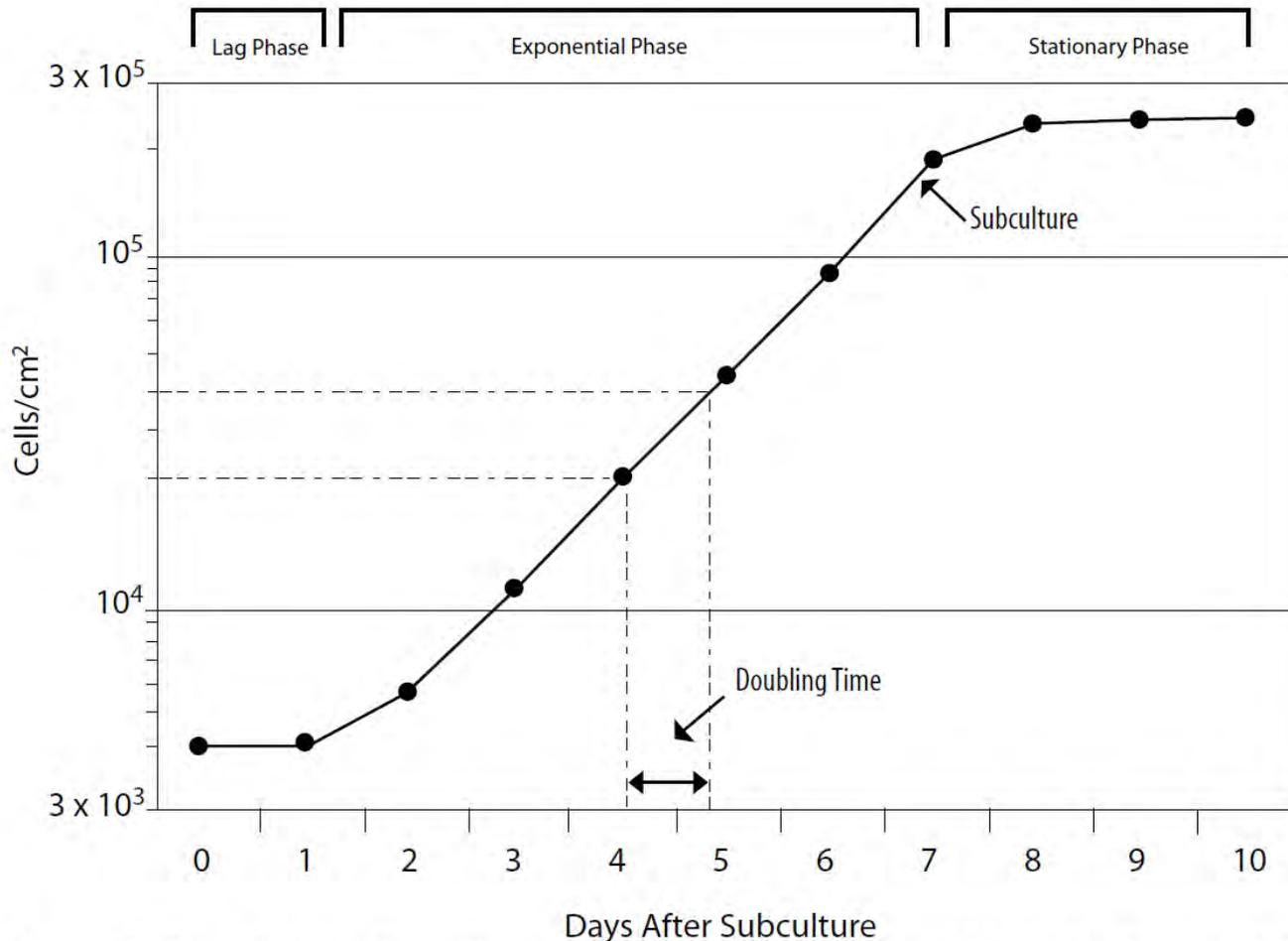


# Cell growth and propagation

- Population doubling level
- Measuring cell viability and growth
  - MTT Assay
  - XTT Assay
  - Reliablue™ Reagent



# Population doubling level



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

# Growth and viability

- Quantitative evaluation of **cell proliferation rate** and response to external factors that affect cell viability
- **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)
- **Reliablue™ Cell Proliferation Reagent (ATCC® 30-1014)**
  - Resazurin (7-Hydroxy-3H-phenoxazin-3-one 10-oxide)

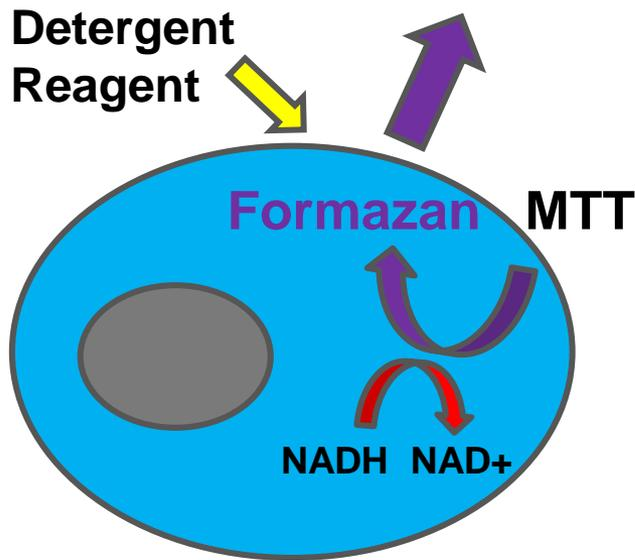
**New!**

# MTT / XTT

## MTT Reaction

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

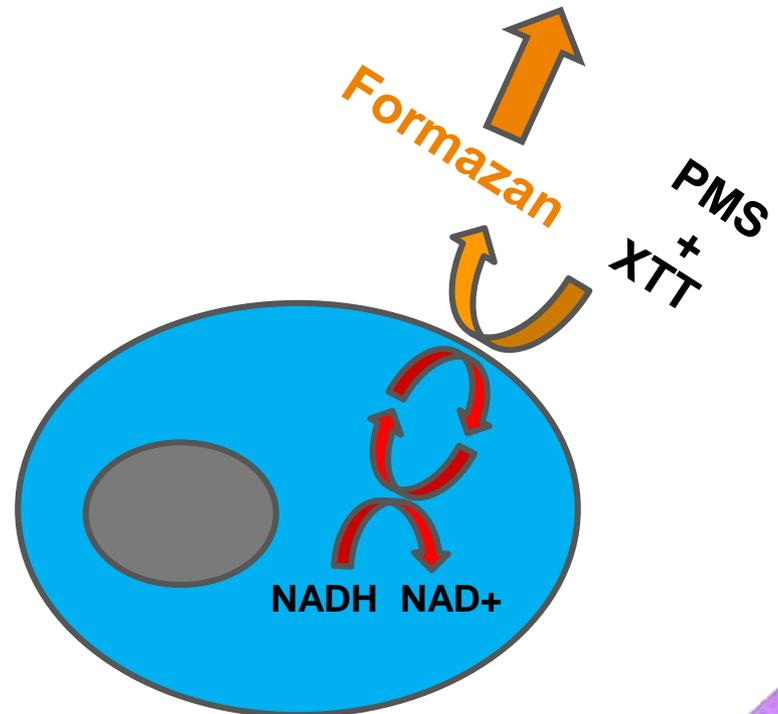
Media turns **PURPLE**



## XTT Reaction

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

Media turns **ORANGE**



# MTT / XTT

## Determining Optimal Cell Counts

- Plate, in triplicate, a serial dilution of  $1 \times 10^6$  to  $1 \times 10^3$  cells per mL (96-well plate)

### MTT Assay

Add MTT Reagent

Incubate 2-4 hours – add Detergent

Incubate 2-4 hours or overnight

### XTT Assay

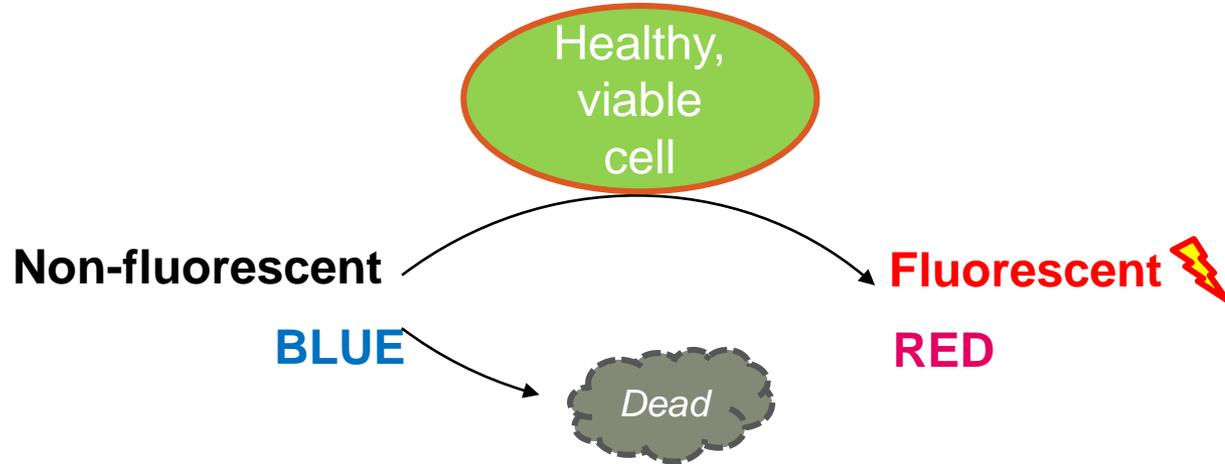
Add XTT Reagent + Activation Agent

Incubate 2-4 hours

- Determine optimal number of cells to use
- Repeat assay with experimental factors – compare absorbance at optimal cell volume
- Plots absorbance versus cell number

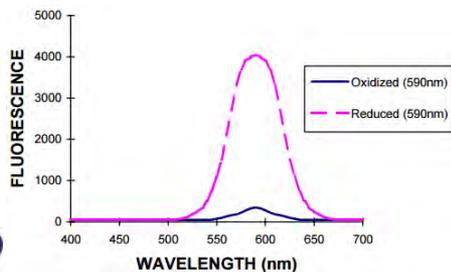
# Reliablue™ Cell Viability Reagent

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.

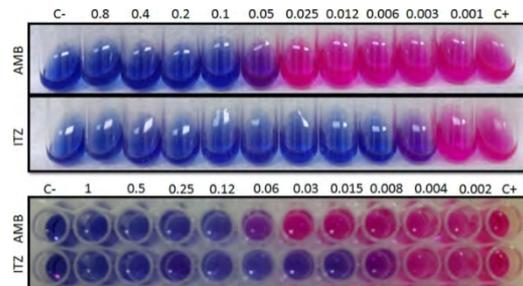
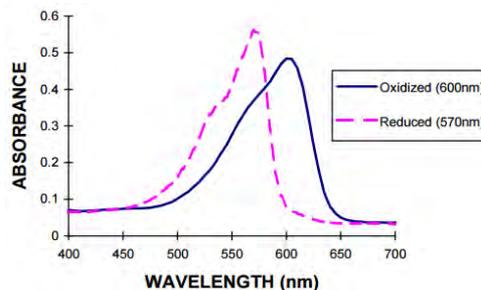


Resazurin is cell permeable but non-toxic and is metabolically reduced by living cells (but not dead cells or in the culture media) resulting in a change in absorbance and increase in fluorescence.

Fluorescence intensity



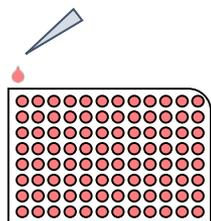
Absorbance Shift



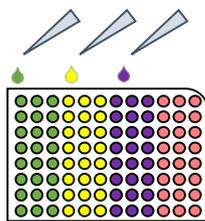
# Reliablue™ Cell Viability Reagent

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.

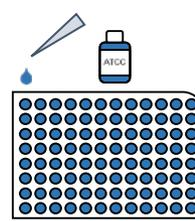
## Basic 4-Step Assay Workflow



① Grow cells



② Add treatments



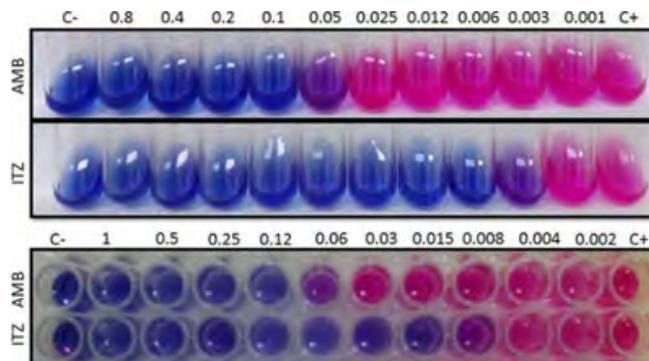
③ Add Reliablue



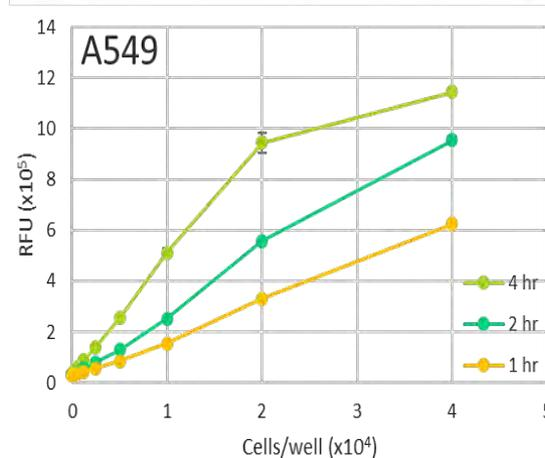
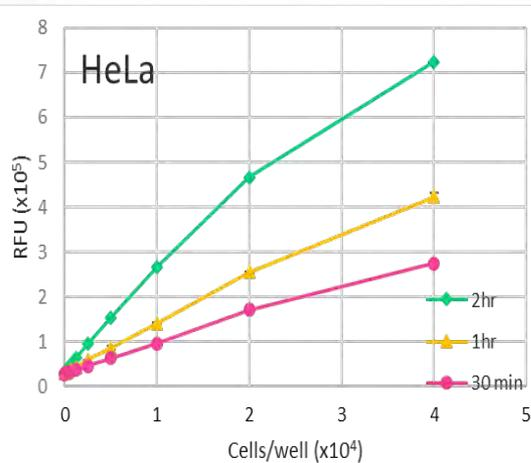
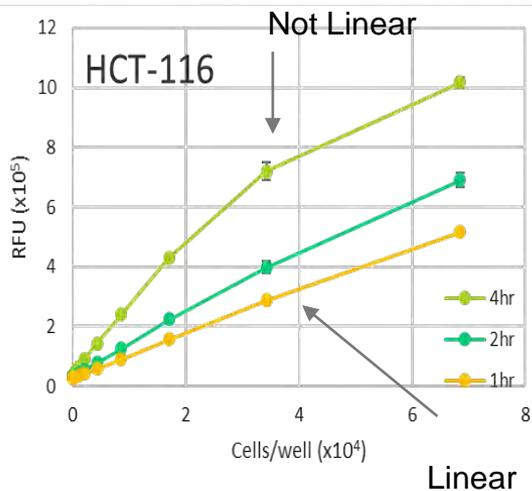
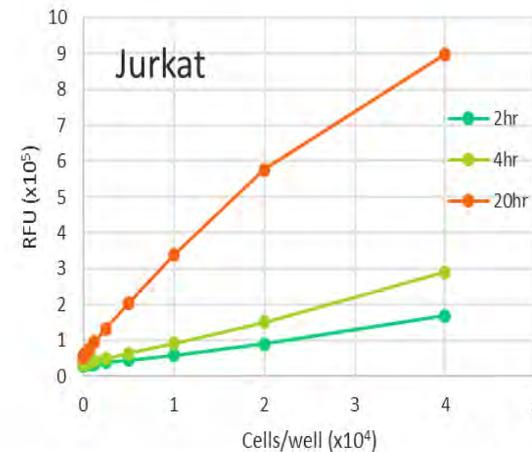
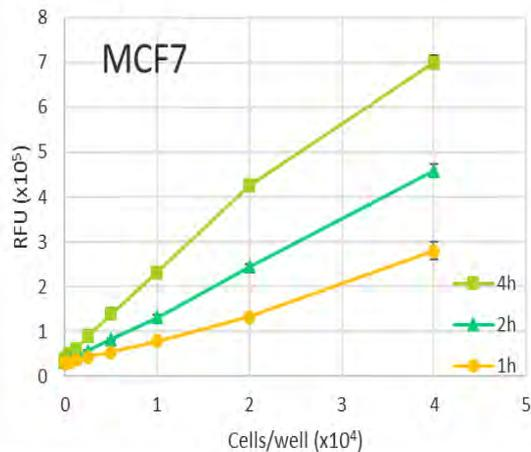
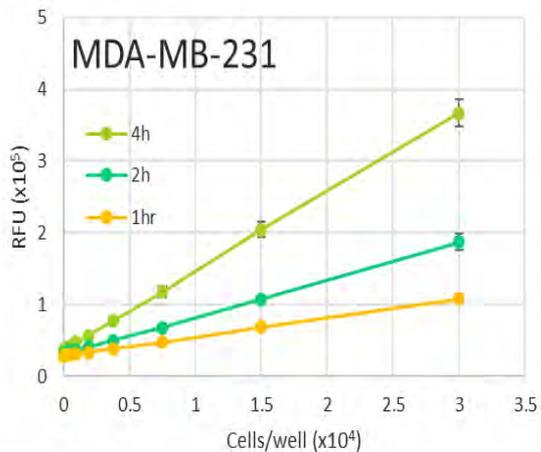
④ Read on plate-reader

Blue = Dead (or control)

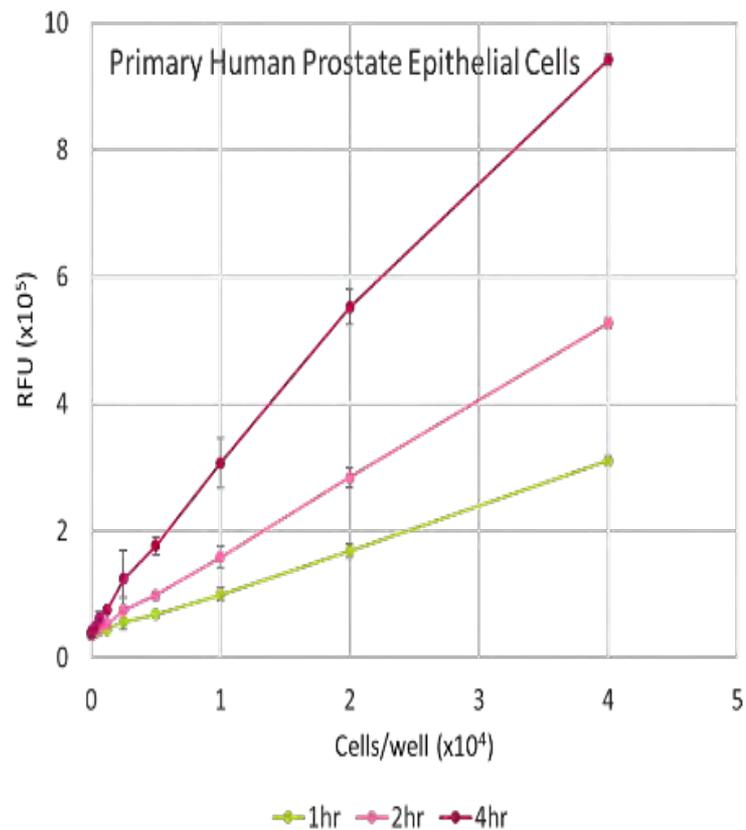
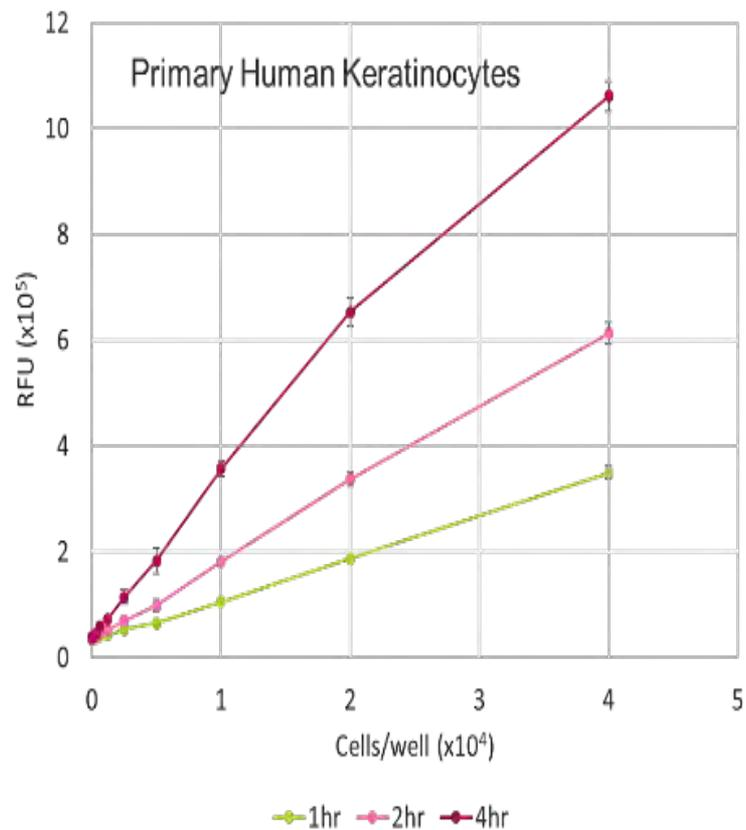
Red = Alive and Actively Growing



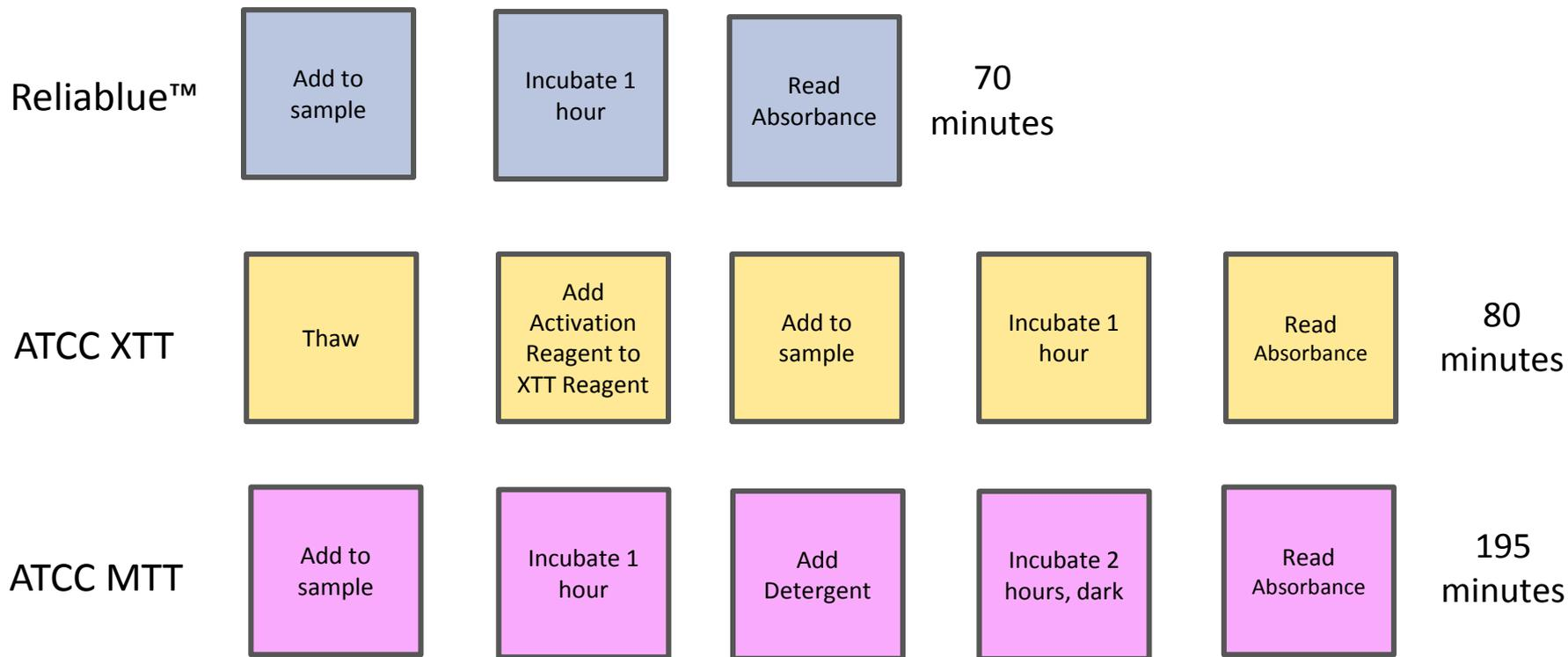
# Cell Tested Reliablue™



# Cell Tested Reliablue™



# Reliablue vs. MTT / XTT Kit



## Reliablue™

- Quick, one step, one reagent
- Nontoxic
- Inexpensive
- **HIGH-THROUGHPUT SCREENING**

# Outline

## Cell Health and Viability Topics

- Cell Culture
  - Media
  - Additives / Serum
- Cryopreservation / Post Thaw
- Cell Proliferation / Viability
  - MTT / XTT kits
  - Reliablue™ cell viability reagent
- Mycoplasma Effects / Detection



# Mycoplasma contamination

- Not easily detected, cannot be seen by microscopy
- 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
- Cell death

Mycoplasma (-)

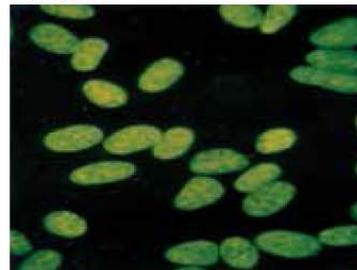


Direct culture method

Mycoplasma (+)

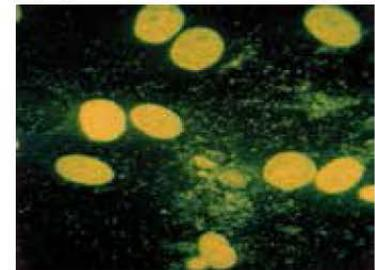


Mycoplasma (-)



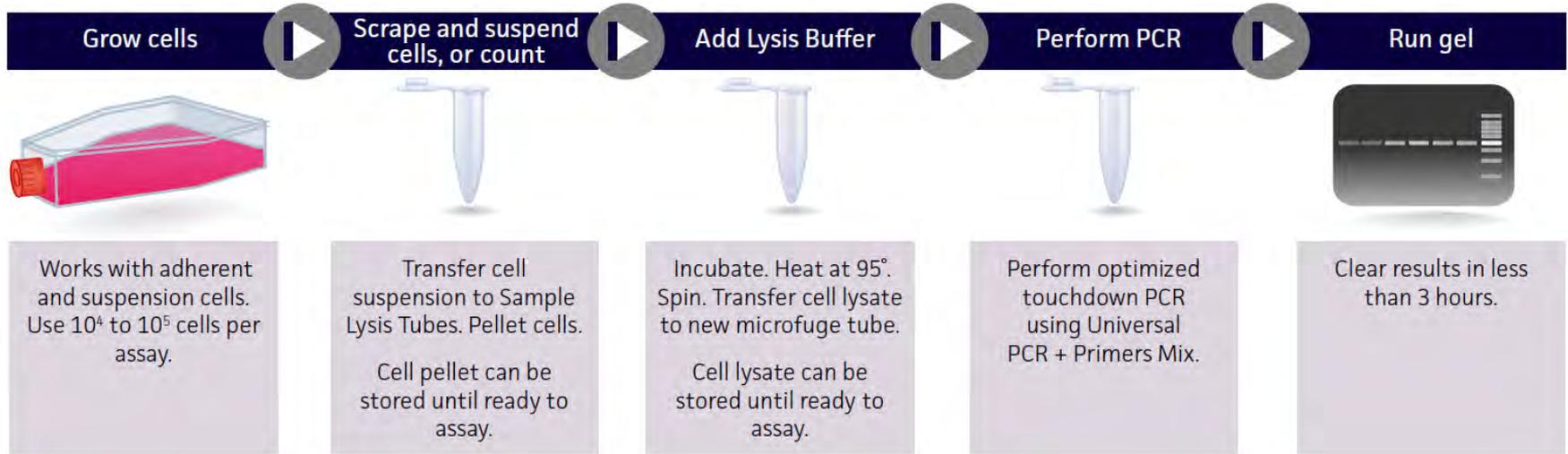
Hoechst DNA staining method

Mycoplasma (+)



# Universal Mycoplasma Detection Kit

## ATCC® 30-1012K™



- Detects over 60 species of *Mycoplasma*, *Acholeplasma*, *Spiroplasma*, and *Ureaplasma*
- All components for the PCR reaction are provided and optimized for amplification



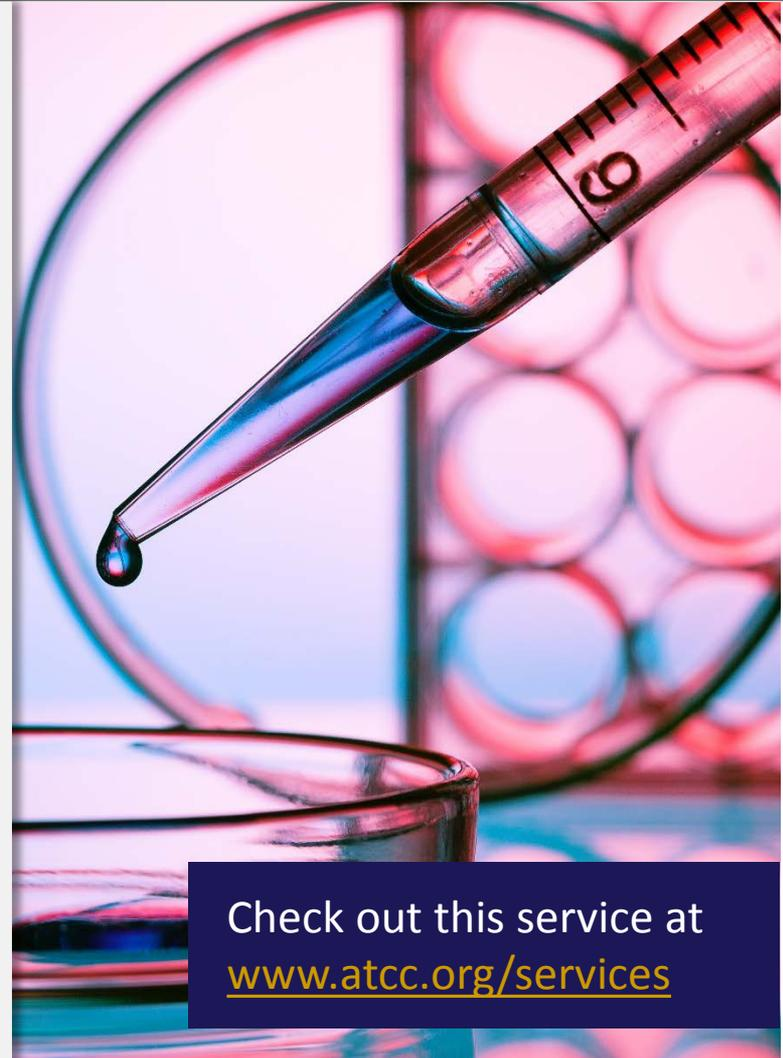
# Mycoplasma testing service

## Direct and indirect culture (bundled service)

- Direct culture – Uses both broth and agar
- Indirect culture – Hoechst DNA stain

## PCR-based testing

- Detection using the ATCC Universal Mycoplasma Detection Kit



Check out this service at  
[www.atcc.org/services](http://www.atcc.org/services)

# Summary points

## Cell culture

- Select appropriate media for your cells
- Understand issues/considerations for adding additional ingredients
- Consistently use same media whenever possible

## Cryopreservation

- Use a reliable rate-controlled cooler
- Keep mammalian cells at  $-140^{\circ}\text{C}$  for long term
- Keep cells in the vapor phase in liquid nitrogen tanks

## Measuring Cell Viability

- Importance of understanding growth rates in cell culture
- Use of viability assays and reagents in measuring proliferation

## Mycoplasma Detection

- Mycoplasma can cause cell death or inhibit proliferation and viability
- Routinely check for mycoplasma in cell cultures

# Disclaimers

© 2017 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.



# Thank you for joining today!

Register for more ATCC webinars at  
[www.atcc.org/webinars](http://www.atcc.org/webinars)

Learn more about cell health and viability at  
[www.atcc.org/cellhealth](http://www.atcc.org/cellhealth)

Learn more about ATCC services at  
[www.atcc.org/services](http://www.atcc.org/services)

