

From Curiosity to Breakthroughs: Accelerate Your Drug Development with Assay Ready Cells

Fatah Kashanchi, PhD, Professor, George Mason University
Heather Branscome, PhD, Senior Scientist, ATCC



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 new product formats that support drug development, safety testing, and high-throughput screening
- Partner with government, industry, and academia
- Leading global supplier of authenticated cell lines, viral and microbial standards
- Sales and distribution in 150 countries, 20 international distributors
- Talented team of 550+ employees, over one-third with advanced degrees



Outline



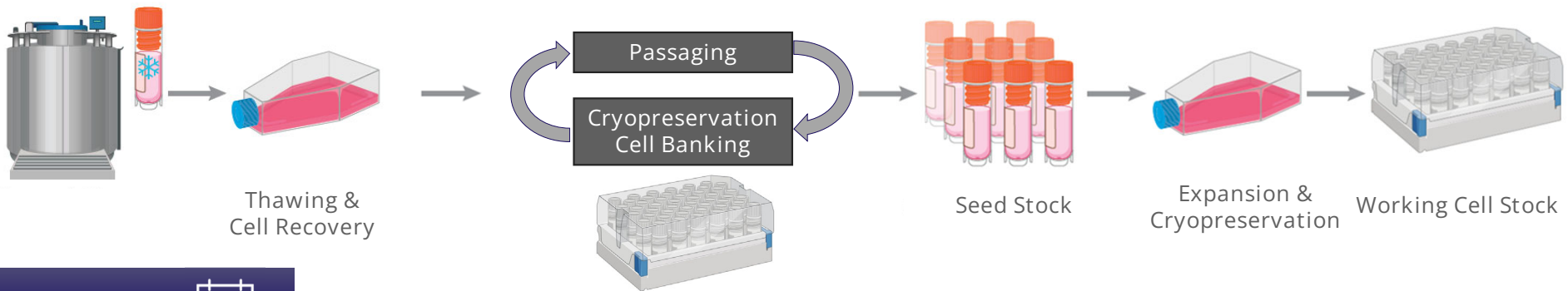
- ThawReady™ introduction
- ThawReady™ solution offerings
- Speaker introduction
- ThawReady™ applications



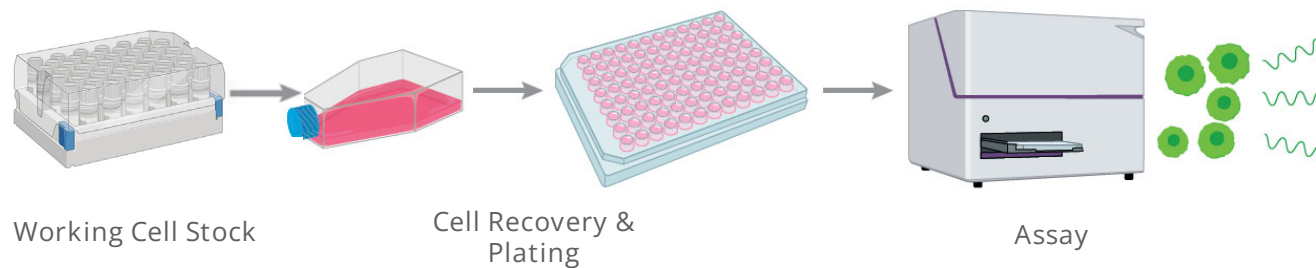
The Problem



Phase 1 Cell Banking



Phase 2 Assay Workflow



ThawReady[™]

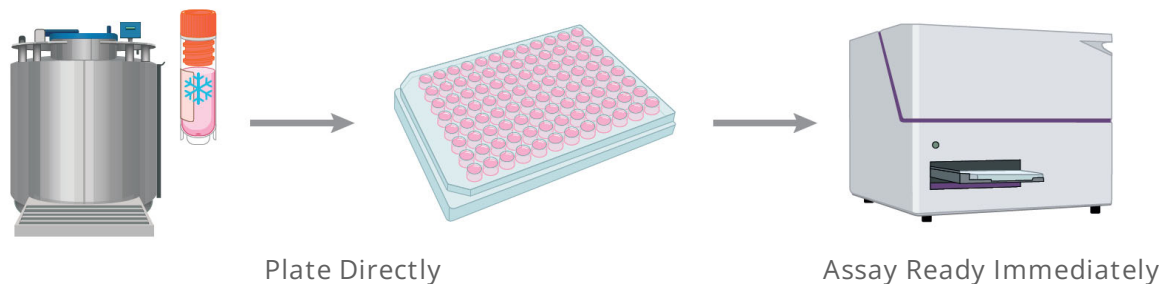
by ATCC



ThawReady™ Solutions



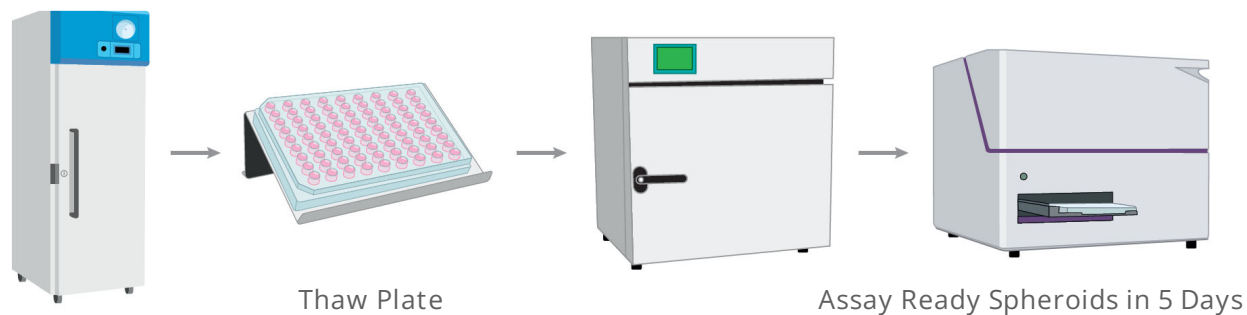
ATCC® ThawReady™ Assay Ready Cells – From frozen to data in one day



Product Format : Vials; Parental and reporter cell lines

ATCC® No.	Product Designation
TIB-202-NFkB-LUC2-AR™	ThawReady™ THP-1 NF-kB-Luc2
TIB-202-AR™	ThawReady™ THP-1

ATCC® ThawReady™ 3-D Spheroid Kits – A uniform spheroid in every well



Product Format : Pre-plated cells; 96-well plates

ATCC® No.	Product Designation
SCM-CCL-185-2PLT™	ThawReady™ A549 Spheroid Kit
SCM-CCL-247-2PLT™	ThawReady™ HCT 116 Spheroid Kit
SCM-HTB-133-2PLT™	ThawReady™ T47D Spheroid Kit

Our Speakers



Fatah Kashanchi, PhD,
Professor, Laboratory of Molecular
Virology, George Mason University



Heather Branscome, PhD,
Senior Scientist, ATCC

Application Data

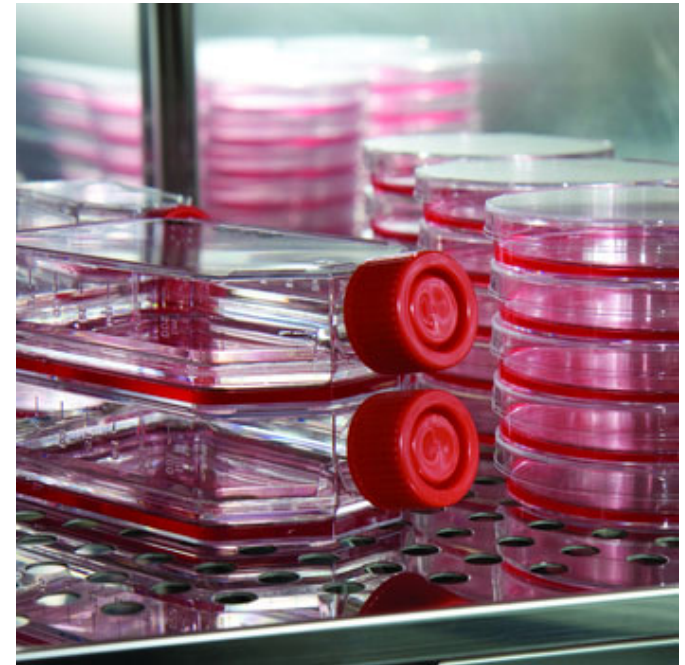
Heather Branscome, PhD



Overview



- Challenges with bioassays
- ThawReady™ overview
- Application data
 - Post-thaw recovery
 - Functionality testing using extracellular vesicles (EVs)
 - HIV-1 infection and drugs of abuse



Challenges with Bioassays



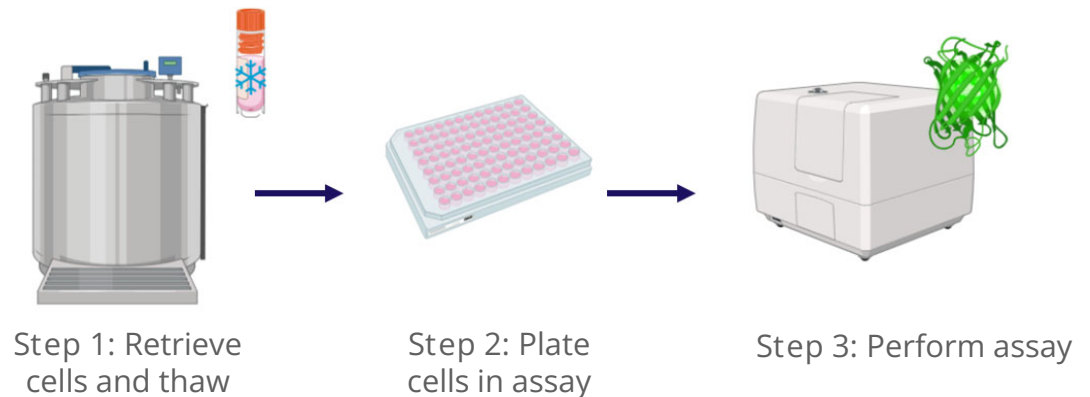
Upstream/downstream challenges

- Time consuming
- Cost and resource dependent
- Introduces variability and genetic drift
- Increased risk of contamination
- Post-thaw recovery period is required



ThawReady™ Solution

- Consumable format
- High reproducibility
- High functionality
- Unique cryobiology



ThawReady™ Overview

THP-1 (ATCC® TIB-202™)

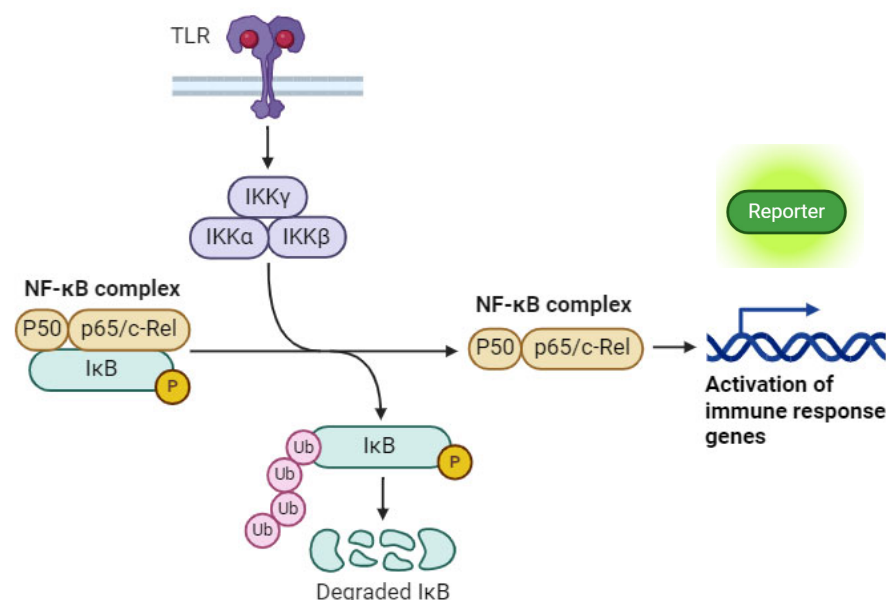
- Monocytic cells derived from the peripheral blood of an acute monocytic leukemia patient
- Versatile and essential for the biomedical research community
- Routinely used in research related to immune system, toxicology/drug development, and cell therapy

ThawReady™ THP-1 NF-κB-Luc2 (ATCC® TIB-202-NFκB-LUC2-AR™)

- Assay-ready, luciferase reporter cell line derived from the THP-1 parental cell line
- Animal by-product free
- Expresses the firefly luciferase gene (luc2) under the control of a NFκB promoter
- Useful for monitoring the activity of NFκB and cellular immune responses in vitro



ThawReady™
by ATCC

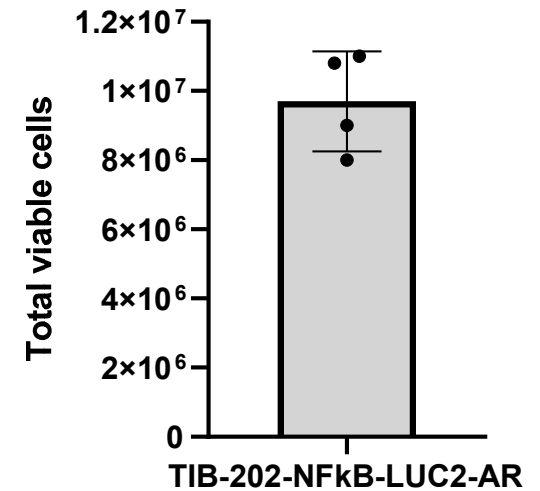
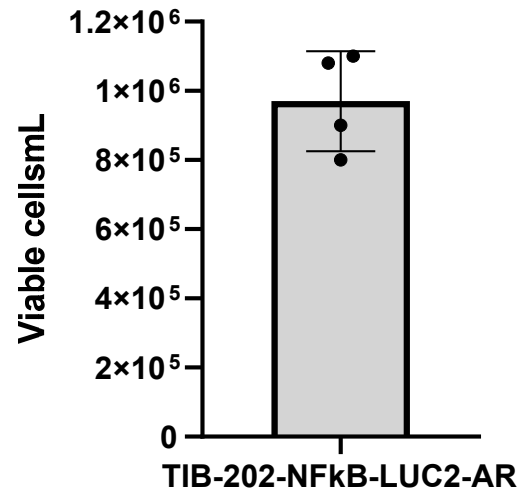
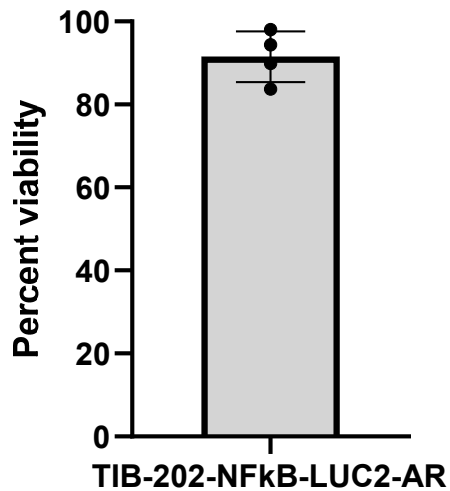


ThawReady™ THP-1 NF-kB-Luc2

ATCC® TIB-202-NFkB-LUC2-AR™



- Reproducible and consistent viability and cell yield
- Measure immediately post-thaw



ThawReady™ THP-1 NF-kB-Luc2

ATCC® TIB-202-NFkB-LUC2-AR™

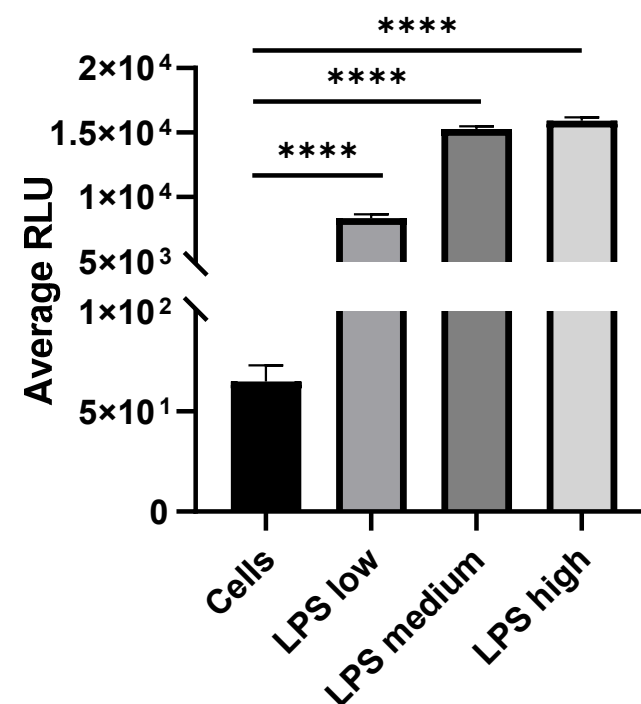


ThawReady™
by ATCC

LPS Challenge Assay

- Average post-thaw viability: 89.9%
- Seeding: Immediate post-thaw
- Seeding density: 50,000 viable cells
- Treatment:
 - LPS (*E. coli* O111:B4)
 - Stimulation period: 3 hours
 - Low dose: 1 µg/mL
 - Medium dose: 10 µg/mL
 - High dose: 25 µg/mL
- Assay/Instrument: Bright-Glo™ assay on BioTek Cytation®

Assayed
Directly
from
Frozen



Application Data: Extracellular Vesical (EV) Functionality

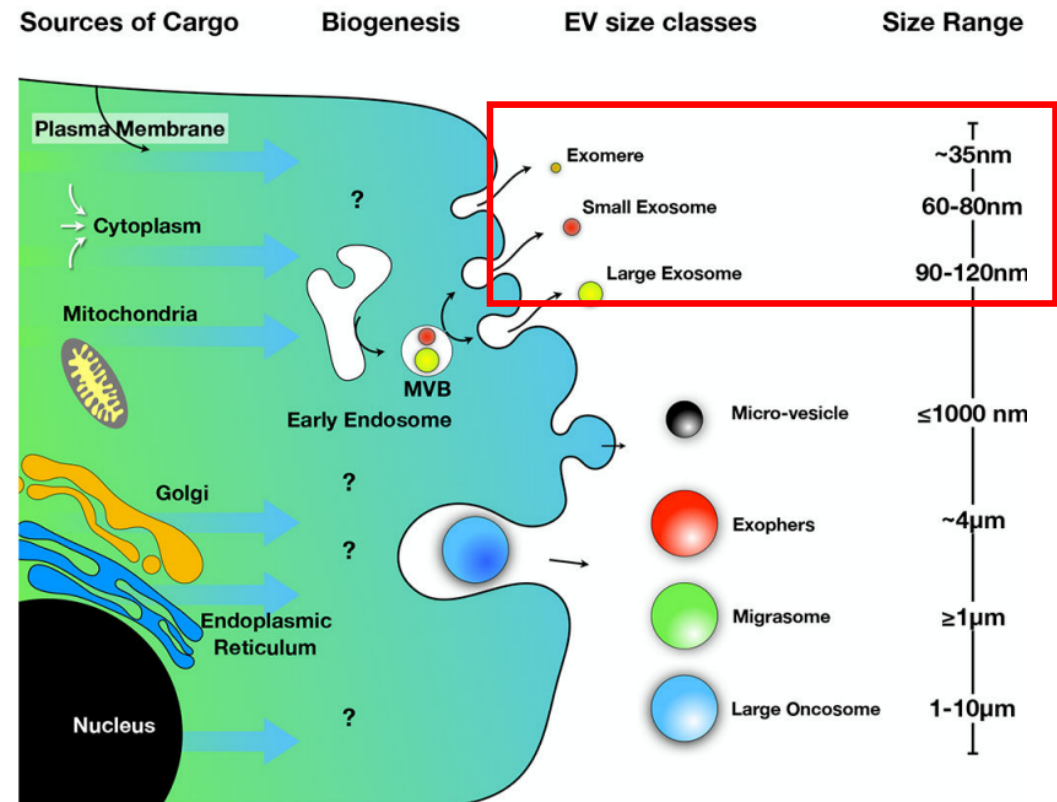


Application Data: EV Functionality



Extracellular Vesicles:

- Nano-sized particles (30 nm-1000 nm) released by all cell types
- Carry various components of the cytoplasm and cell membrane
- Mediate intercellular communication (physiological and pathological)
- Can be utilized as diagnostic markers (e.g., cancer EVs) or therapeutic tools (e.g., stem cell EVs)



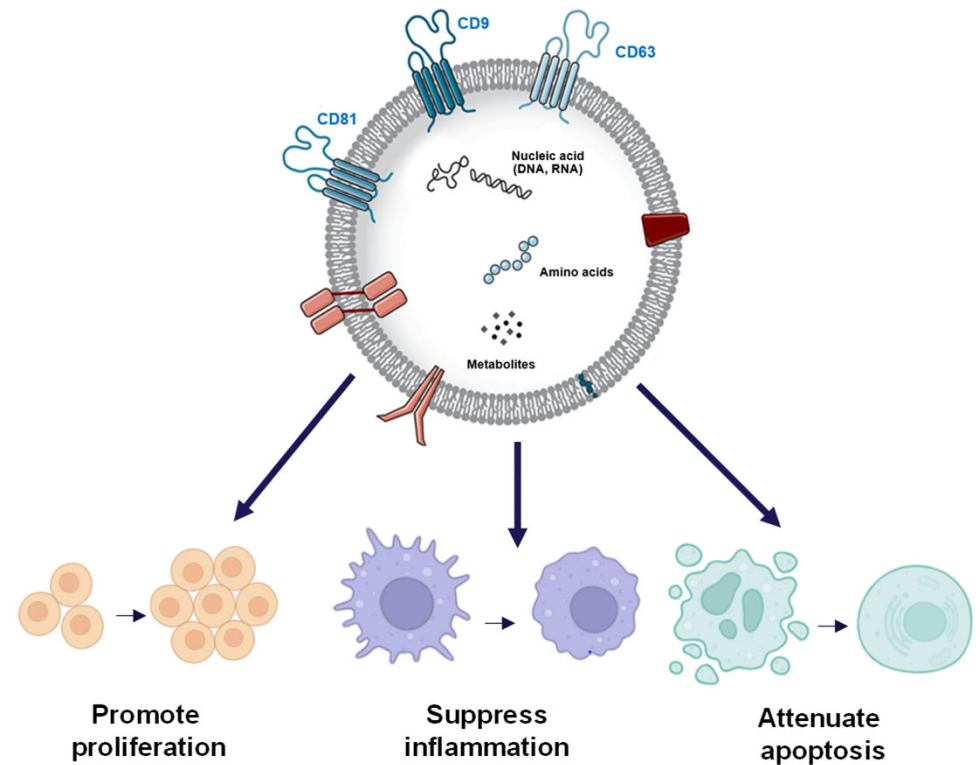
Zijlstra A, Di Vizio D. Nat Cell Biol 20(3): 228-230, 2018. PubMed: 29476154

Application Data: EV Functionality



Stem cell EVs:

- Contribute to tissue repair and regeneration
- Have reduced immunogenicity compared to parental stem cells
- Can be engineered to carry additional therapeutic agents to target damaged cells
- Can be engineered to carry additional therapeutic agents to target cells



Application Data: EV Functionality

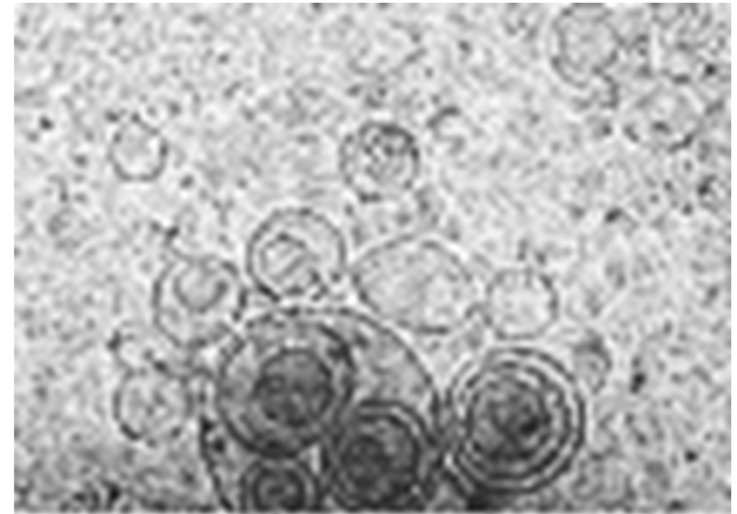


Donor cells:

- Mesenchymal Stem Cells (MSCs) (ATCC® PCS-500-012™)
- Induced Pluripotent Stem Cells (iPSCs) (ATCC® ACS-1019™)
- hTERT-immortalized MSCs (ATCC® SCRC-4000™)

In vitro functionality:

- Pro-migratory
 - Epithelial cells, endothelial cells, fibroblasts
- Pro-angiogenic
 - Mesenchymal stem cells and aortic endothelial cells (co-culture)
- Anti-apoptotic
 - Neurons (ATCC® CRL-2266™)
 - Astrocytes (ATCC® CRL-1718™)
 - Monocyte-derived macrophages (ATCC® TIB-202™)
 - Retinal epithelium (ATCC® CRL-2302™)
- Anti-inflammatory
 - Neurospheres (ATCC® ACS-5003™)



doi.org/10.1007/s11481-019-09865-y



doi.org/10.1038/s41598-022-05848-x



doi.org/10.3390/cells13100861

Application Data: EV Functionality



ATCC® No.	Parental Cell Designation	Model	Applications (pre-clinical)
SCRC-4000-EXM™	hTERT-immortalized adipose-derived mesenchymal stem cell (MSC)	Stem cell (non-cancer)	Therapeutics/ Drug delivery
CRL-1740-EXM™	LNCaP	Carcinoma, prostate	Diagnostics/biomarker discovery
CRL-1435-EXM™	PC-3	Adenocarcinoma, prostate	Diagnostics/biomarker discovery
CCL-185-EXM™	A549	Carcinoma, lung	Diagnostics/biomarker discovery
CCL-247-EXM™	HCT 116	Carcinoma, colorectal	Diagnostics/biomarker discovery

Application Data: ATCC® Stem Cell EVs

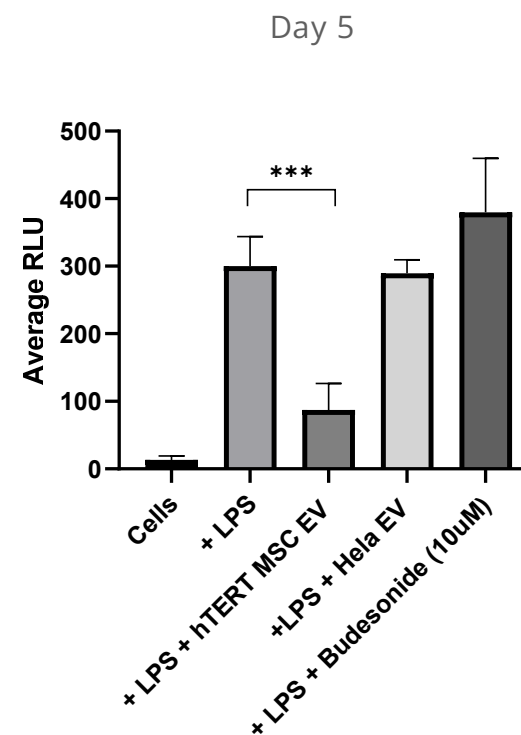
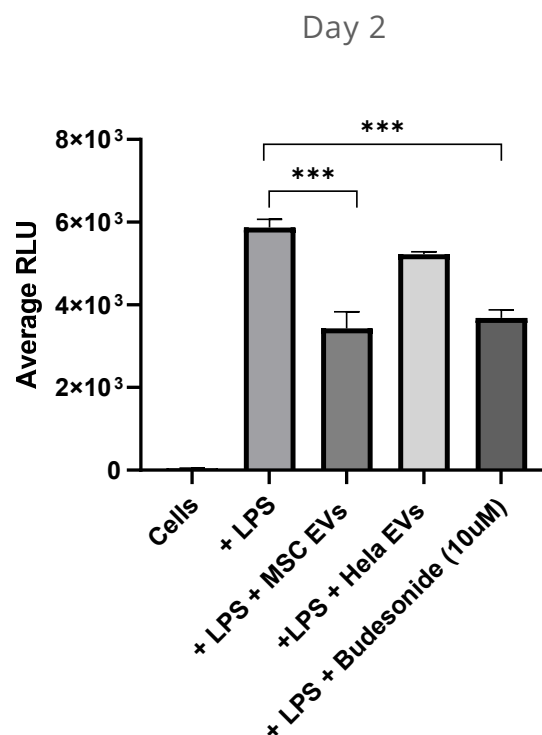
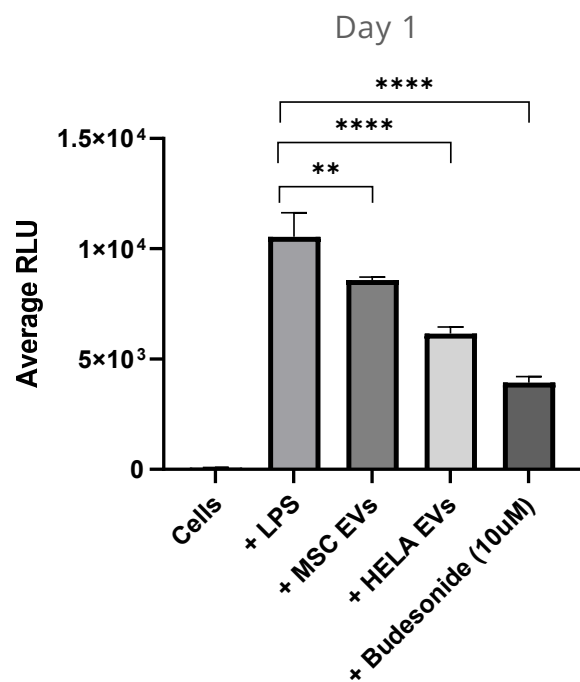
TIB-202-NFkB-LUC2-AR™ EV Function Assay



- Seeding: Immediate post-thaw
- Seeding density: 50,000 viable cells
- Treatment:
 - LPS: *E. coli* O111:B4
 - 3 hour stimulation; 10 µg/mL
 - LPS + stem cell EVs
 - LPS + cancer cell EVs
 - Positive control: Budesonide (10µm)
- Assay/Instrument: Bright-Glo™ assay/CellTiter-Glo® on BioTek Cytation®

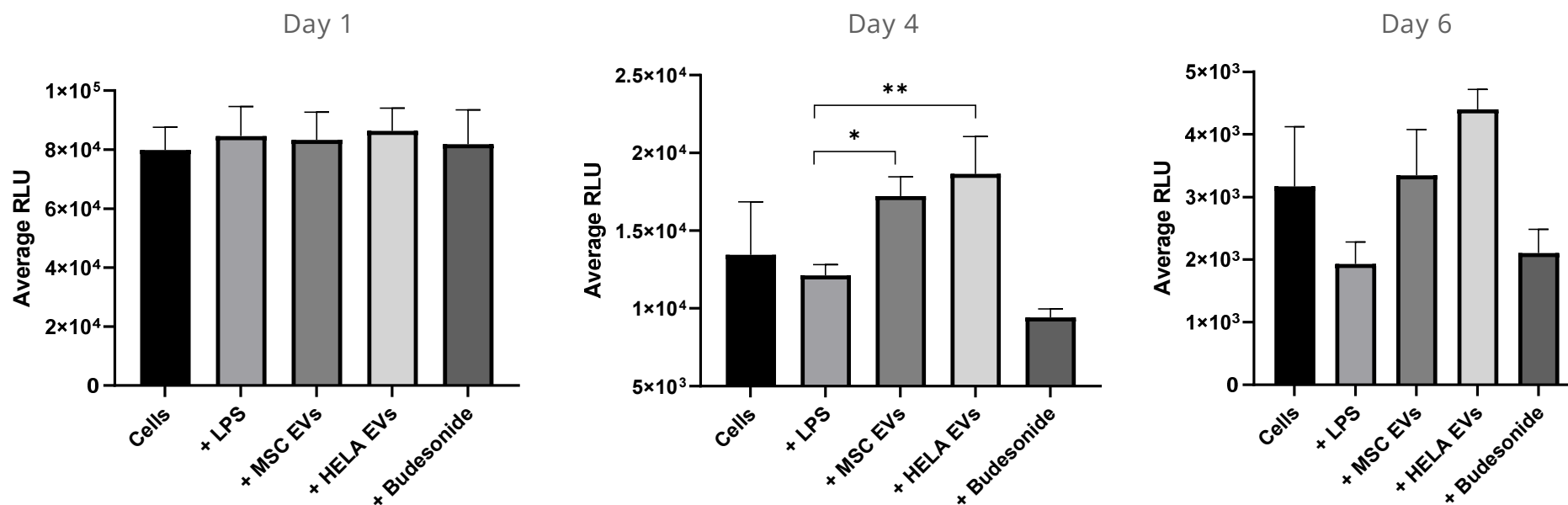
Application Data: ATCC® Stem Cell EVs

ThawReady™
by ATCC



Application Data: ATCC® Stem Cell EVs

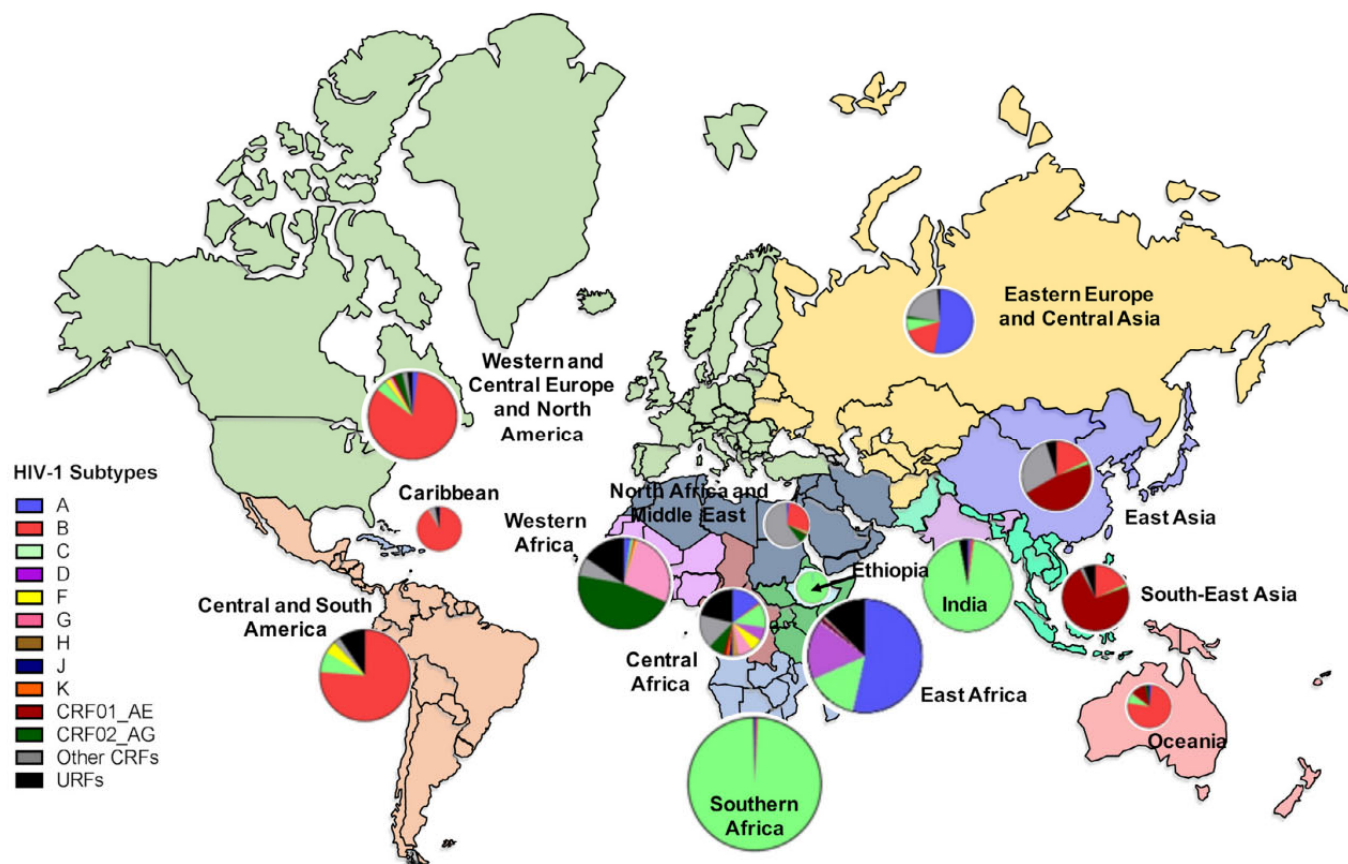
ThawReady™
by ATCC





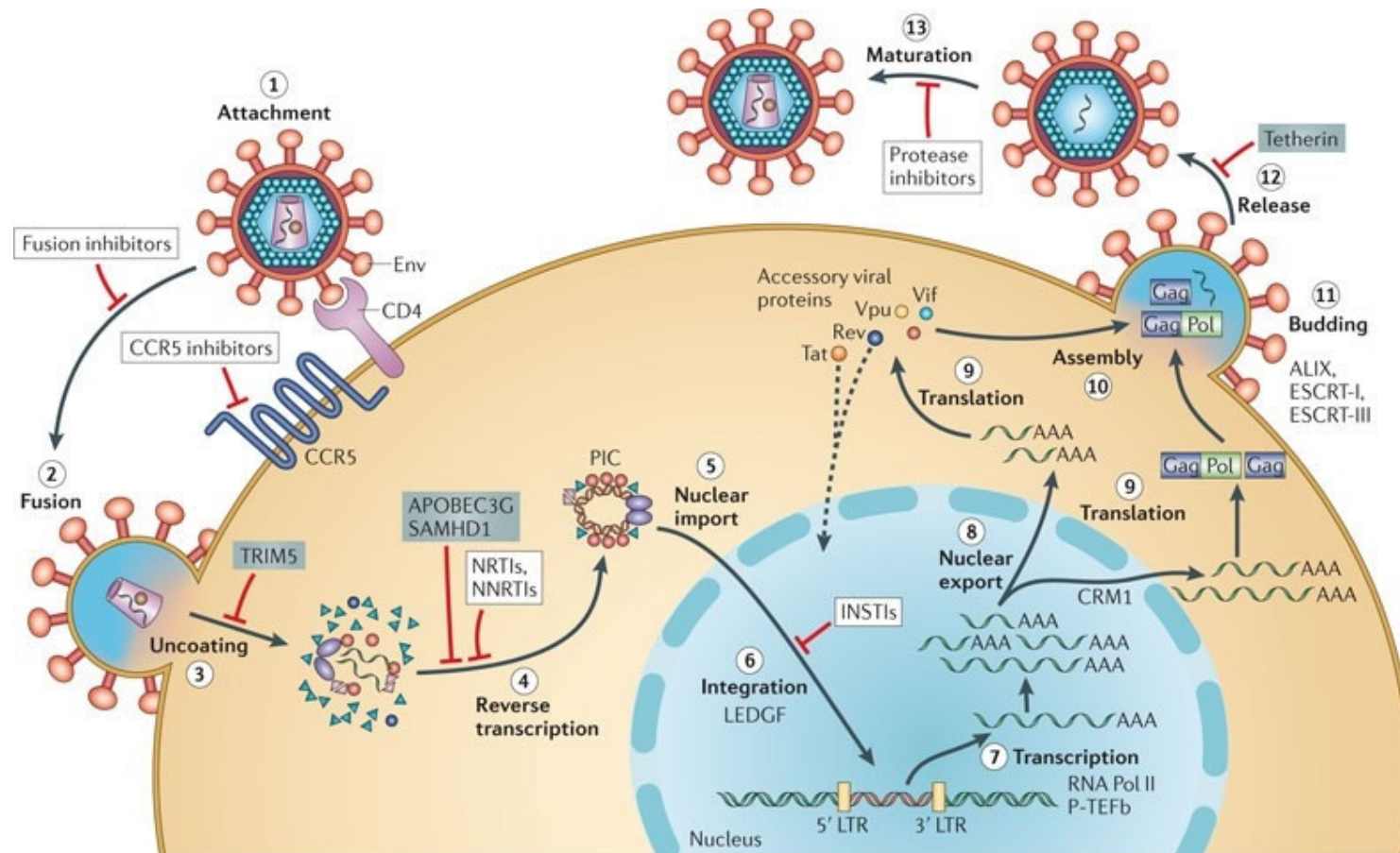
Application Data: HIV-1 Infection

Global prevalence of HIV-1



Gartner MJ, et al. EBioMedicine 53: 102682, 2020. PubMed: 32114391

Application Data: HIV-1 Infection



Engelman A, Cherepanov P. Nature Rev Microbiol 10: 279-290, 2012.

Application Data: HIV-1 Infection

ATCC® TIB-202-NFkB-LUC2-AR™ HIV-1 infection assay

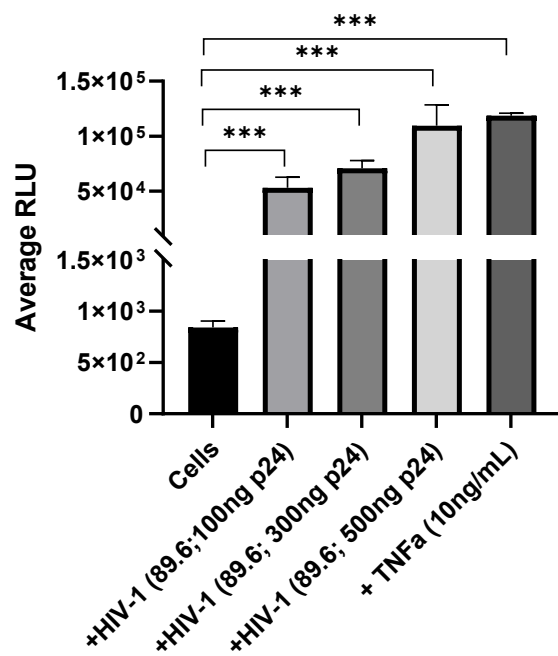
- Seeding: Immediate post-thaw
- Seeding density: 50,000 viable cells
- Treatment:
 - PMA (100 ng/mL)
 - HIV-1 dual tropic 89.6 (100, 300, 500 ng p24)
 - cART (5 mM)
 - TNFα (10 ng/mL)
 - CBD (5 uM)
 - HU308 (5 uM)
- Assay/Instrument: Bright-Glo™ assay/GloMax® (multiple reads over 10 minutes)



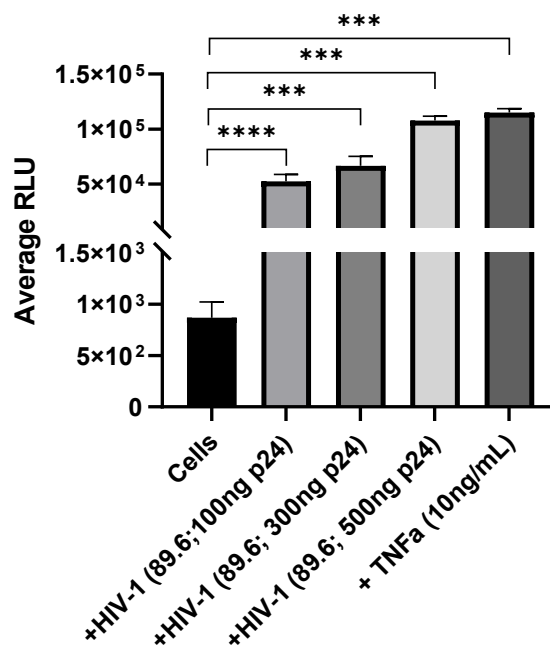
Application Data: HIV-1 Infection

ThawReady™
by ATCC

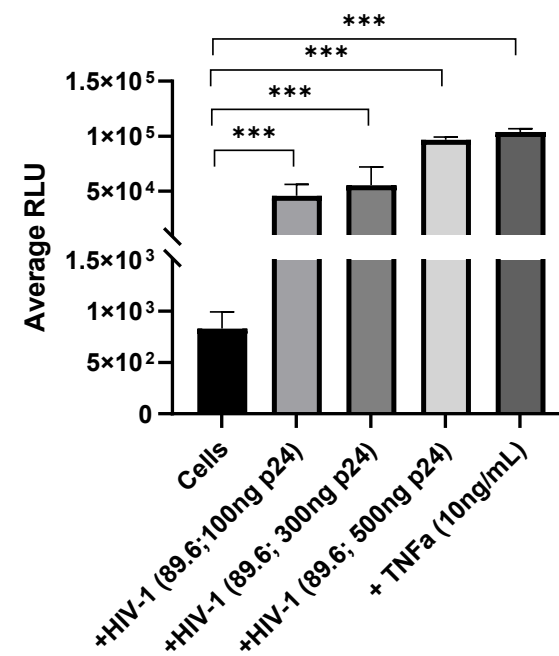
Assay 1



Assay 2



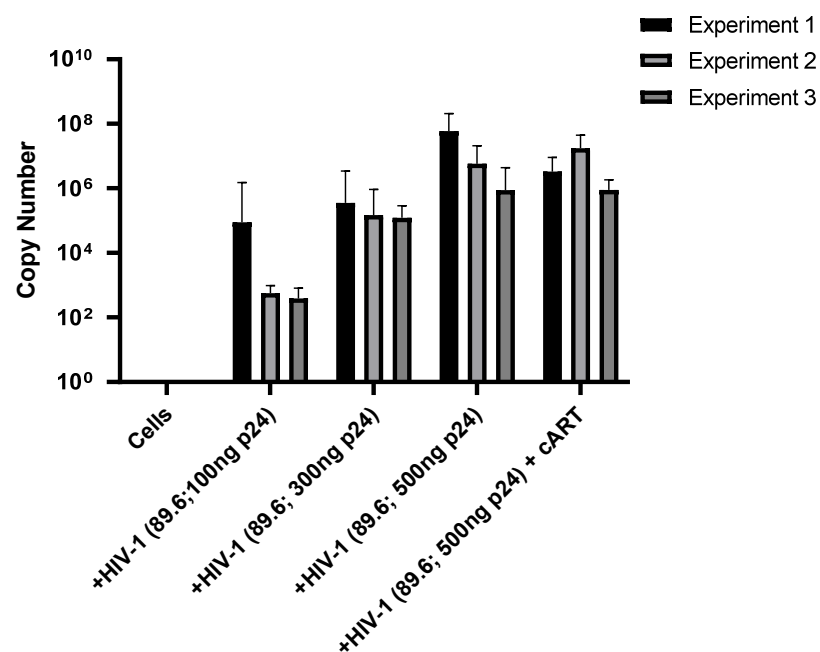
Assay 3



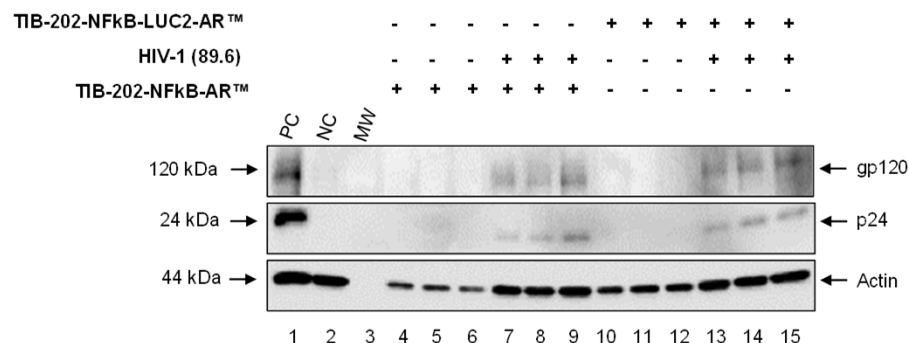
Application Data: HIV-1 Infection

ThawReady™
by ATCC

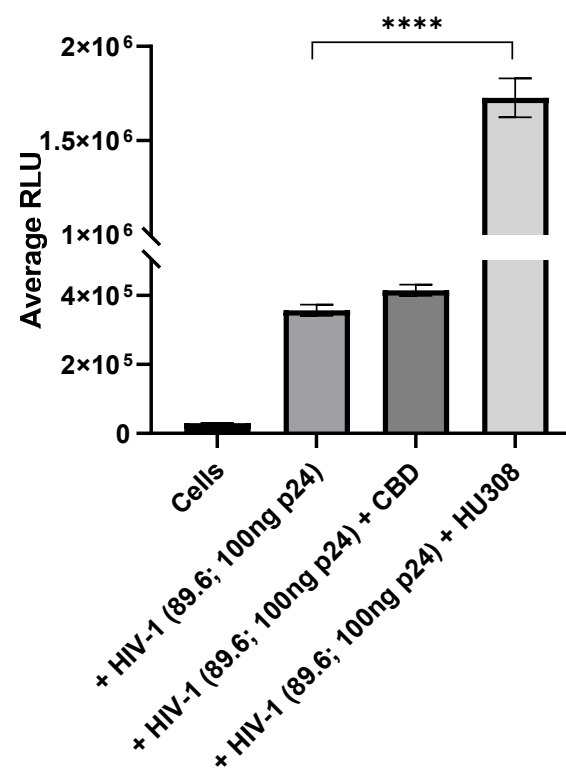
a)



b)

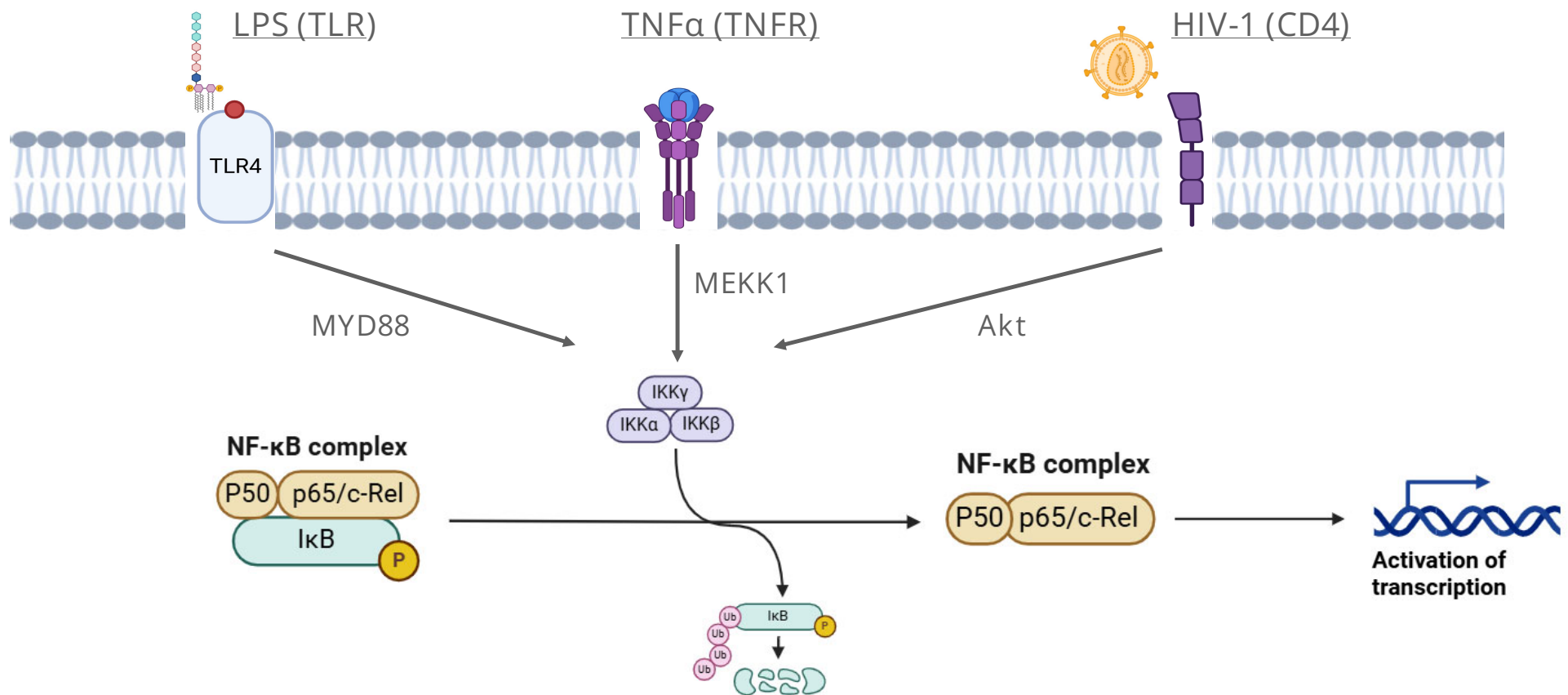


c)



Mechanism

Activation of NF κ B signaling via multiple stimulants/receptors



Summary



- ThawReady™ cells overcome many of the challenges associated with traditional bioassays
- ThawReady™ cells offer high viability, high cell yields, and high functionality in a reproducible manner
- ThawReady™ cells can be used in assays relating to cellular damage (e.g., cancer EVs, viral infection) or cellular repair (e.g., stem cell EVs, drug screening)
- ThawReady™ cells can be used to study various mechanisms of NFκB activation, including TLR4, TNFR, and viral attachment

Acknowledgements



ATCC

- Dr. Joseph Leonelli
- Becky Bradford
- Dr. Heather Couch
- Dr. Sujata Choudhury
- Dr. Fang Tian
- Zach Cuba
- Sarah Swineford
- Jessica Moore
- Dr. Utsav Sharma
- Dr. Brian Shapiro
- Dr. Nilay Chakraborty
- Tiffany Cato



GMU

- Dr. Yuriy Kim
- Kajal Patil
- Anastasia Williams
- Gwen Cox



Learn more about ThawReady™ cells



Visit us online at
www.atcc.org/thawready-cells



Listen to our podcast episode 26:
ThawReady™ – Transforming Cryopreservation for
Instant Cell Functionality
www.atcc.org/behindthebiology