

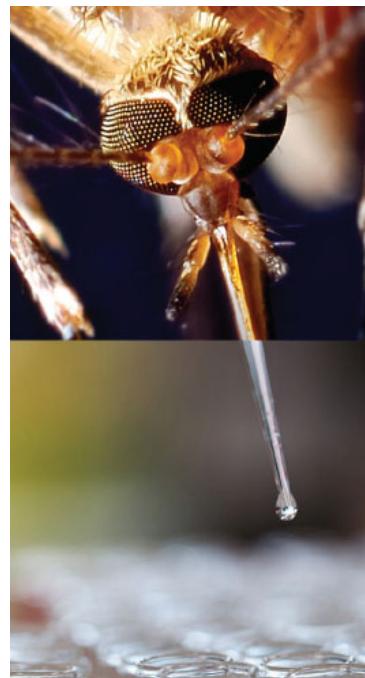
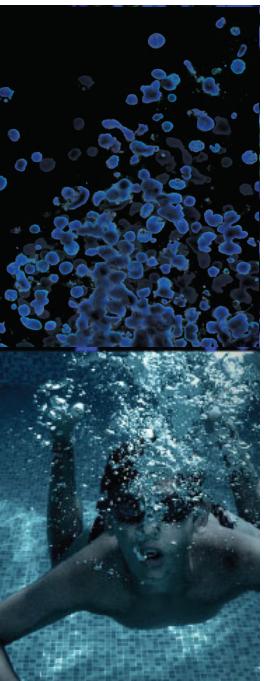


# Extracellular Vesicles: For Better and for Worse

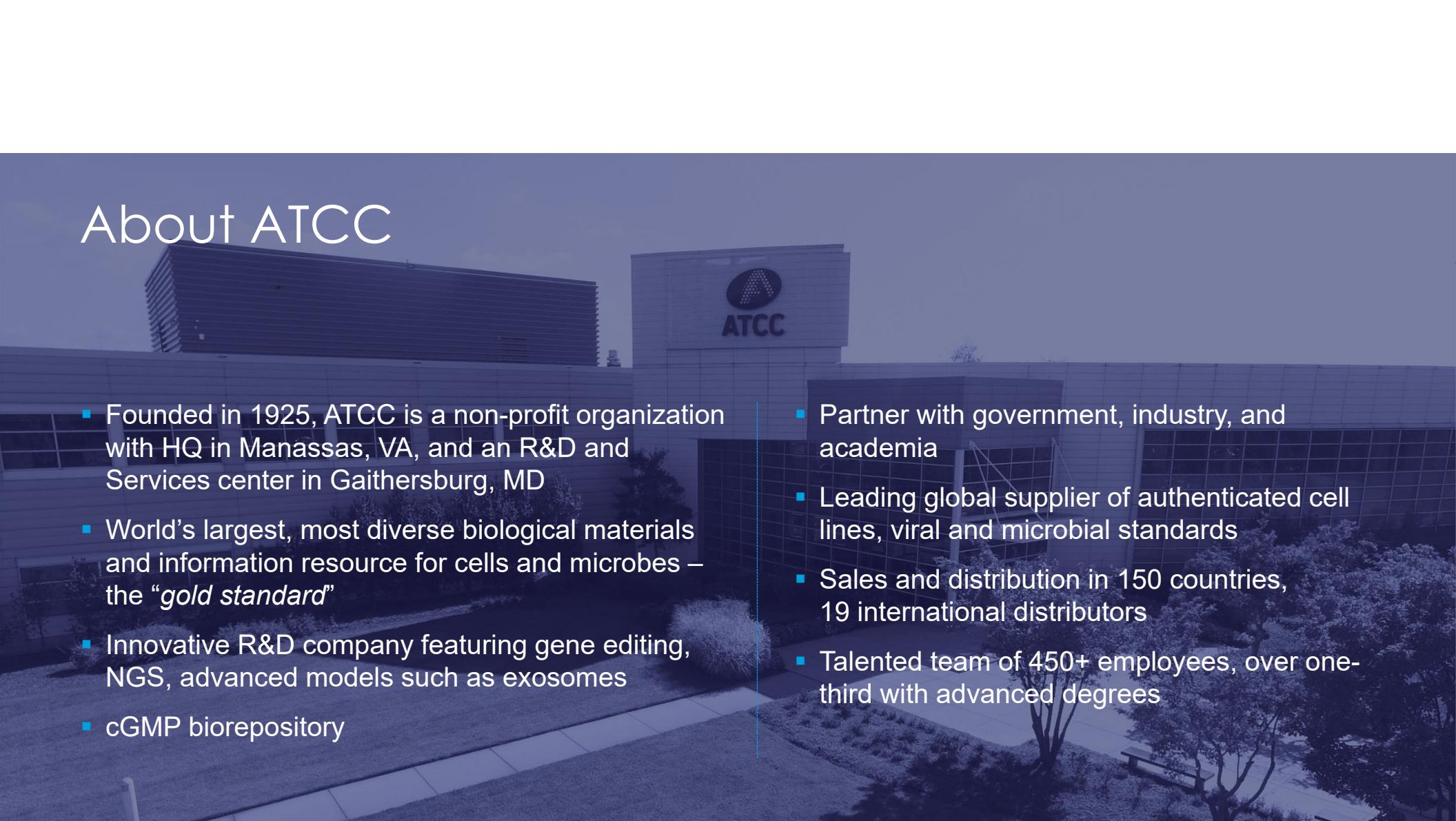
Heather Branscome, M.S.  
*Lead Biologist, ATCC*

Fatah Kashanchi, Ph.D.  
*Professor, George Mason University*

Credible Leads to Incredible™



# About ATCC

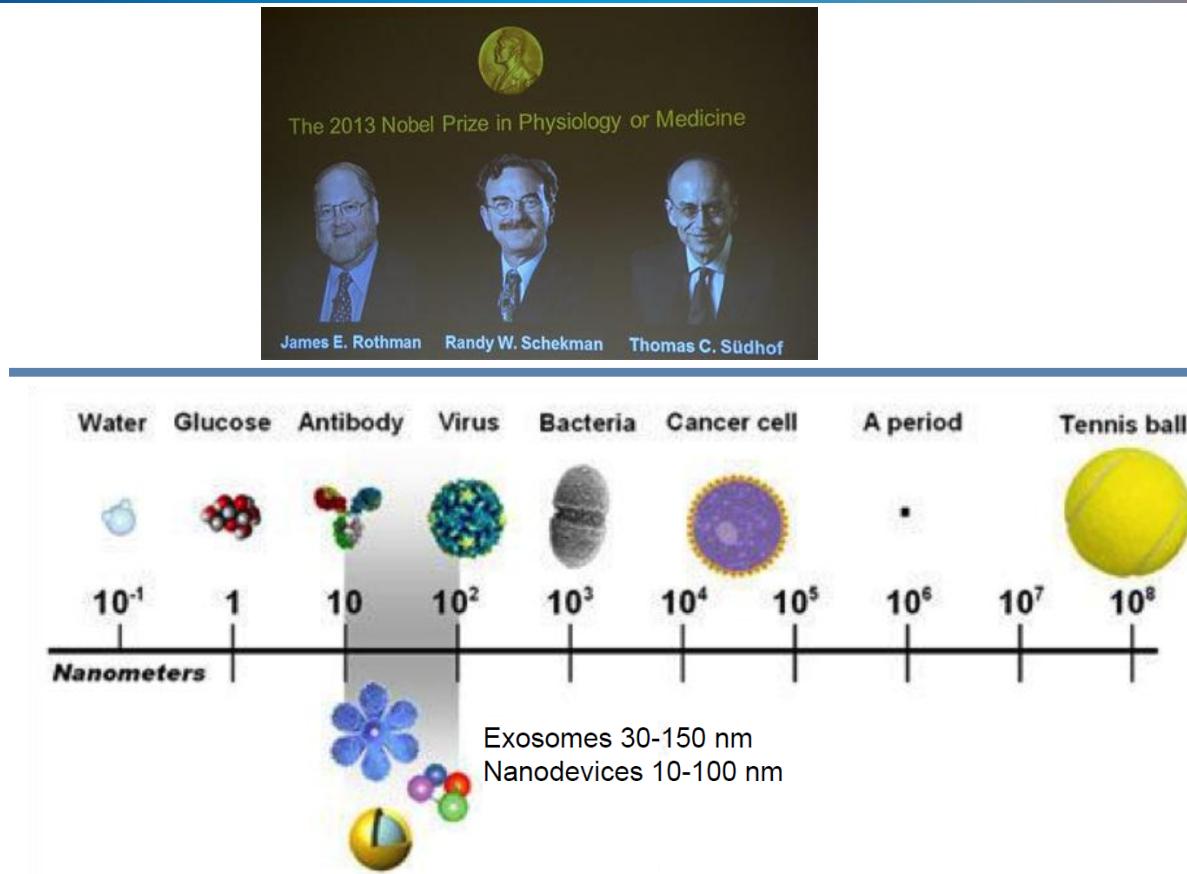
- 
- Founded in 1925, ATCC is a non-profit organization with HQ in Manassas, VA, and an R&D and Services center in Gaithersburg, MD
  - World's largest, most diverse biological materials and information resource for cells and microbes – the “*gold standard*”
  - Innovative R&D company featuring gene editing, NGS, advanced models such as exosomes
  - 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 one-third with advanced degrees

# Overview

---

- Overview of Extracellular Vesicles (EVs)
- Characterization and Analysis of “Damaging EVs”
- Characterization and Analysis of “Reparative EVs”
- Summary and Future Directions

# Overview

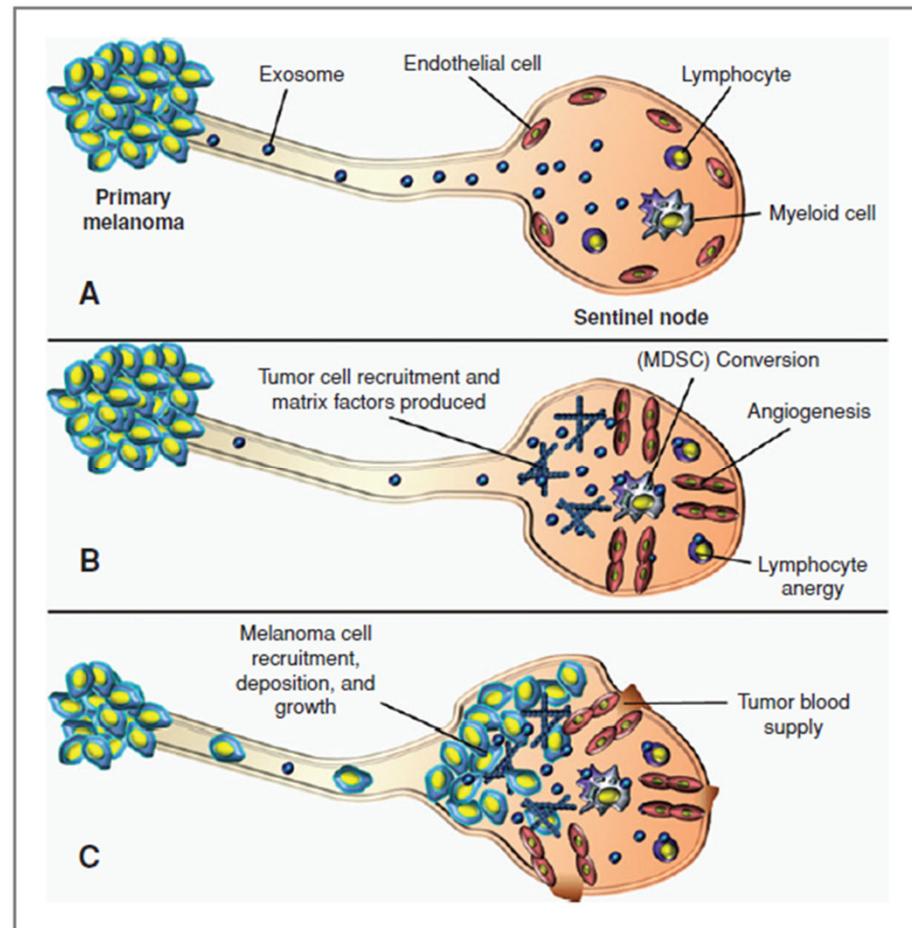
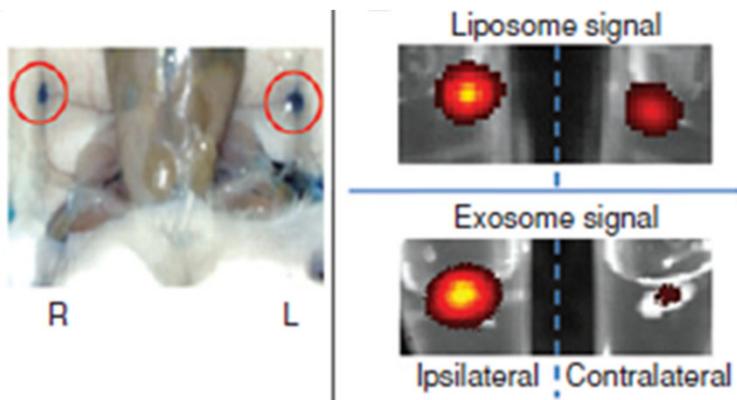


Internal volume of exosomes (diameter 30-150 nm) is  $\sim 4-1500 \times 10^{-24} \text{ m}^3$

The volume of one average 50 kDa protein or 100 nt RNA molecule is  $\sim 6 \times 10^{-26} \text{ m}^3$

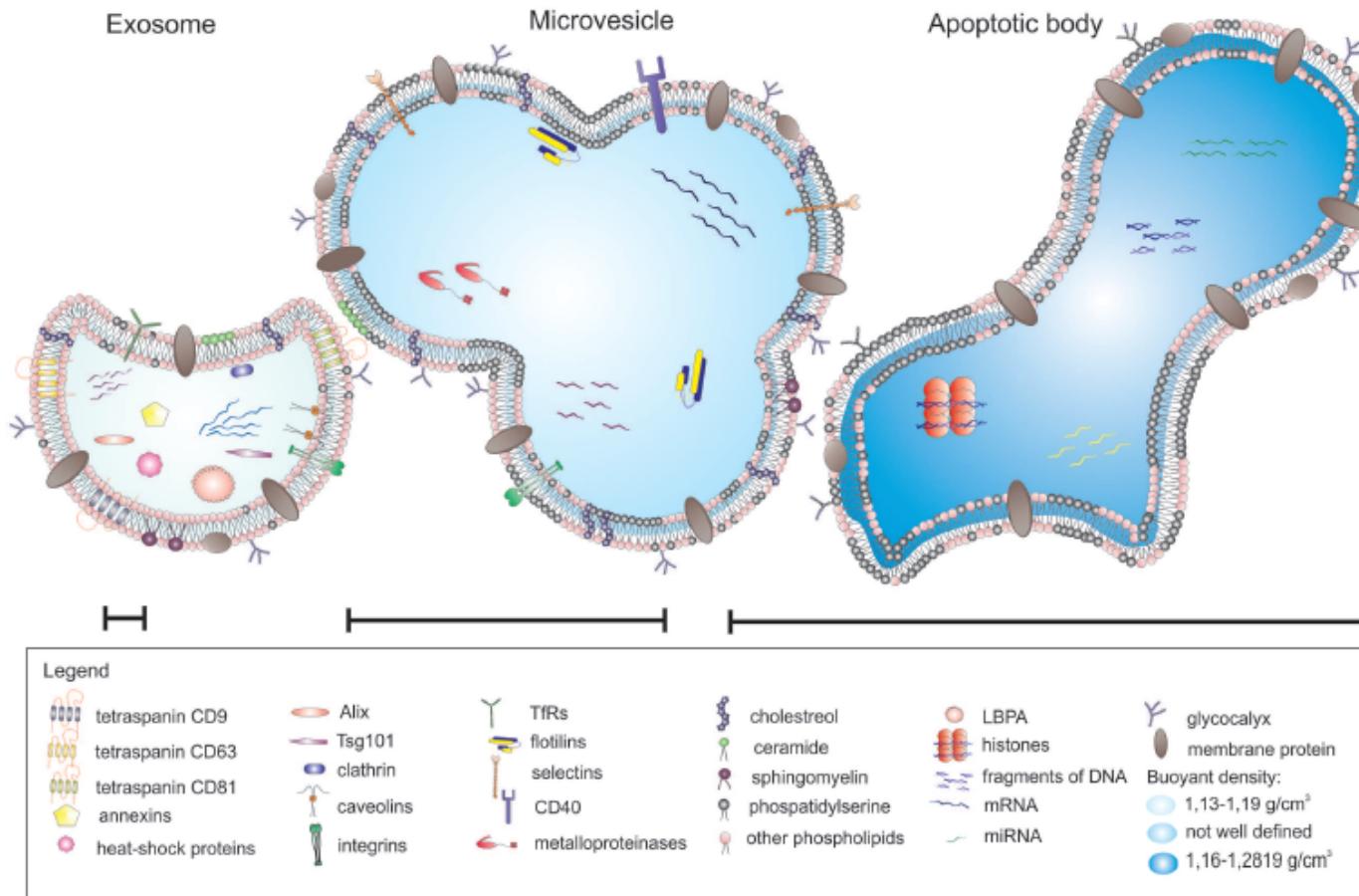
->Exosome theoretically can hold up  $\sim 70-25,000$  small proteins or short RNA

# Clinical Significance of EVs



Hood et al. *Cancer Res.* 2011

# Types of EVs



Koniusz et al. *Frontiers in Cellular Neuroscience*. 2016



# EVs (from uninfected cells)

## Exosomes

- Well defined shape under EM
- 50 to 100 nm
- Formed intracellularly
- Contain tetraspanin proteins: CD63, CD81, CD9
- Contain ESCRT-pathway proteins ALIX and TSG101

## Exosome-like vesicles

- 20 to 50 nm
- Originate from multivesicular bodies formed by other organelles (Golgi)
- Lack ESCRT proteins ALIX and TSG101

## Microvesicles (ectosomes)

- 100 to 1000 nm
- Originate from outward budding/fission of plasma membrane (exocytosis)
- Do not contain tetraspanin proteins
- Similar lipid composition to plasma membrane

## Apoptotic bodies

- 500 to 4000 nm
- Formed only during programmed cell death
- Contain DNA and histones

# EV Research and Infections: Tip of an Iceberg

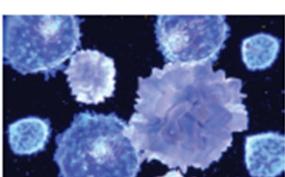
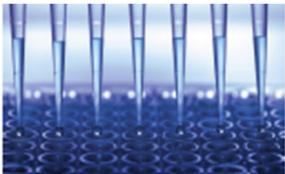
Baltimore groups are:

- I: dsDNA type (examples: Adenovirus, Herpesvirus, Poxvirus)
- II: ssDNA type (+)sense DNA (example: Parvovirus)
- III: dsRNA type (example: Reovirus)
- IV: (+)ssRNA type (+)sense RNA (examples: Picornavirus, Togavirus)
- V: (-)ssRNA type (-)sense RNA (examples: Orthomyxovirus, Rhabdovirus)
- VI: ssRNA-RT type (+)sense RNA with DNA intermediate to life-cycle (example: Retrovirus)
- VII: dsDNA-RT type (example: Hepadnavirus)

PubMed Search:	# of Manuscripts:
Viruses	~10 <sup>6</sup>
Intracellular Pathogens & Viruses (all DNA and RNA)	1,749
Intracellular Pathogens & Fungi (Candida, Cryptococcus, Aspergillus)	6,049
Intracellular Pathogens & Parasites (Leishmania, Malaria, Chagas, Sleeping sickness)	1,300
Intracellular Pathogens & Bacteria (TB, Listeria, Salmonella)	7,703

PubMed Search:	# of Manuscripts:
Parasites	239,892
Bacteria	~2.1 x 10 <sup>6</sup>
Fungi	~1.6 x 10 <sup>6</sup>

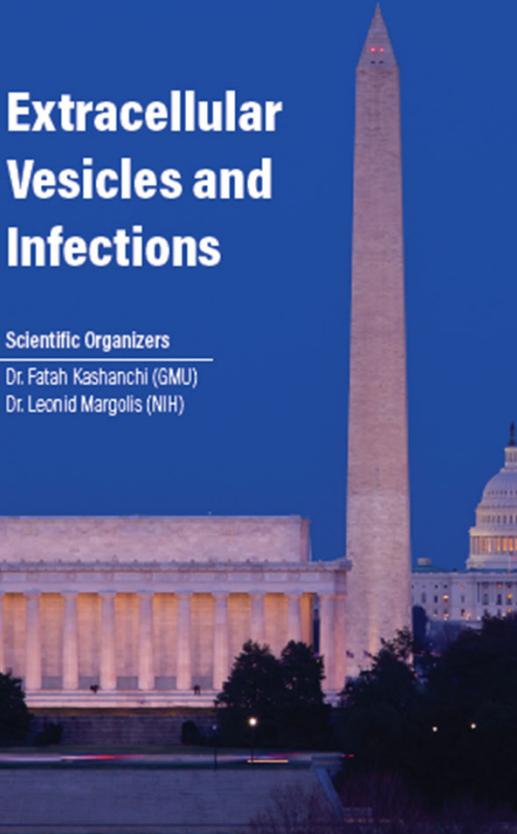




# Extracellular Vesicles and Infections

## Scientific Organizers

Dr. Fatah Kashanchi (GMU)  
Dr. Leonid Margolis (NIH)



**June 1 – 2, 2018**

Bolger Center in Potomac, Maryland



## Exosomes in Human Infectious Diseases Conference

05 Mar 2020 - 08 Mar 2020 | Nassau, Bahamas

### Important Dates

Early Bird	19 Sep 2019
Talk Submission	26 Sep 2019
Poster Submission	09 Jan 2020
Registration Deadline	10 Jan 2020

### Synopsis

Extracellular vesicles (EVs), which are generated by almost all living cells, are now considered to be an important system of cell-cell communication. A large volume of data has been published on EVs in different fields of biology and medicine, from immunology to marine biology. These data are discussed at large international meetings that gather thousands of researchers.

However, there is still a need for more-focused meetings at which the importance and reliability of the data should be evaluated by a narrow circle of experts. Here, we propose to organize a small conference that will focus on one of the most important aspects of EVs, their role in infectious human diseases. The aim of the proposed conference is to improve our understanding of the mechanisms of this involvement and to evaluate EVs as potential therapeutic targets.

Join us alongside chairs Leonid Margolis (NIH/NICHD), Fatah Kashanchi (George Mason University) and Graça Raposo (Institut Curie, CNRS)

### Plenary Speakers

Randy Schekman  
University of California, Berkeley  
Philip Stahl  
Washington University in St. Louis  
Yoel Sadovsky Magee  
Womens Research Institute  
Andrew Hill  
La Trobe University

Stephen Gould  
Johns Hopkins University  
Guillaume van Niel  
INSERM  
Genoveffa Franchini  
NIH

- \* 50% registration support for selected young investigators
- \* Poster Prizes Available

Registration rate includes 3 nights' fixed accommodation (05, 06 and 07 Mar 2020) and a 24-hour all-inclusive food and beverage package for the conference period.

[www.fusion-conferences.com](http://www.fusion-conferences.com)

# Overview

---

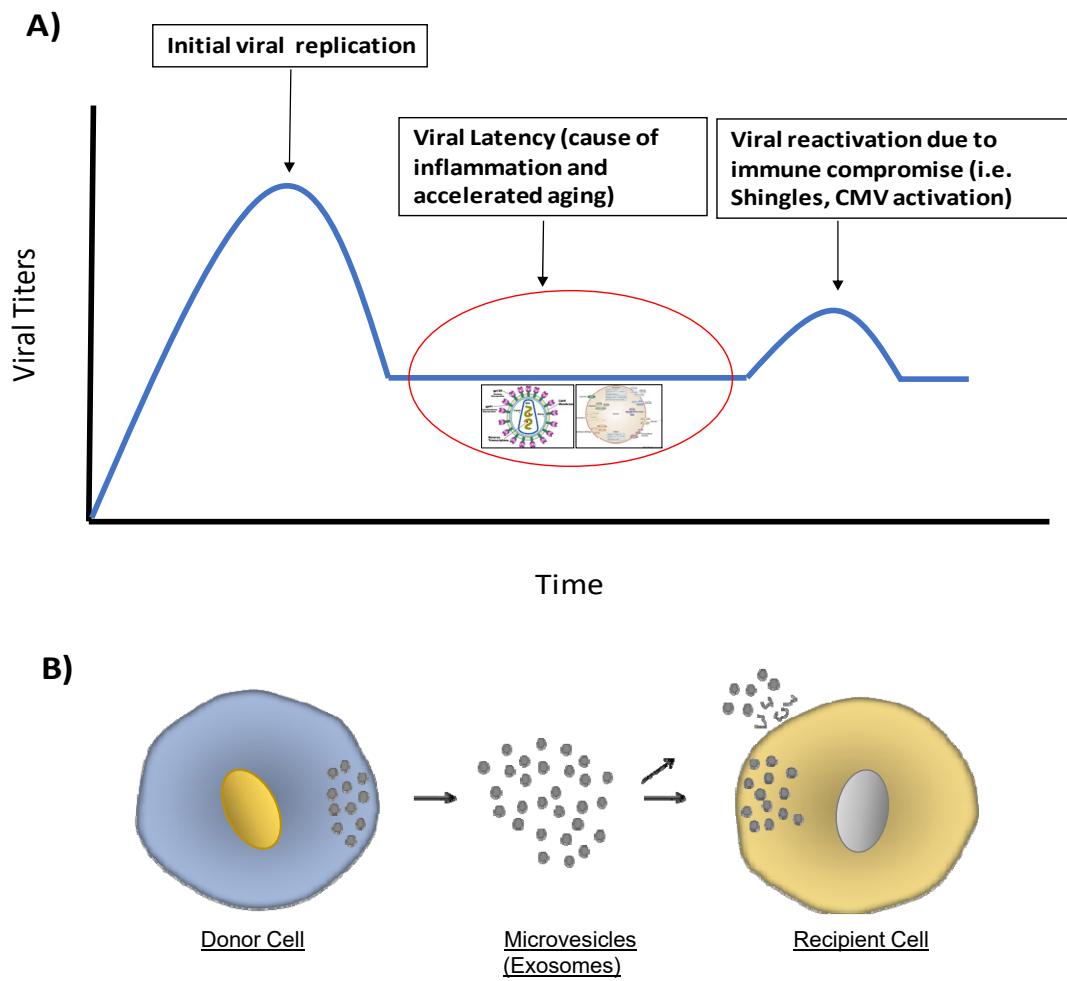
- Overview of Extracellular Vesicles (EVs)
- Characterization and Analysis of “Damaging EVs”
- Characterization and Analysis of “Reparative EVs”
- Summary and Future Directions

## Topics to Be Discussed

---

1. Background on viruses, autophagy and EVs  
(damaging EVs; diagnostics)
2. Isolation of EVs from latent HIV-1 and other virally infected cells (HTLV-1, RVFV, Ebola and Zika)

# Characterization and Analysis of “Damaging EVs”

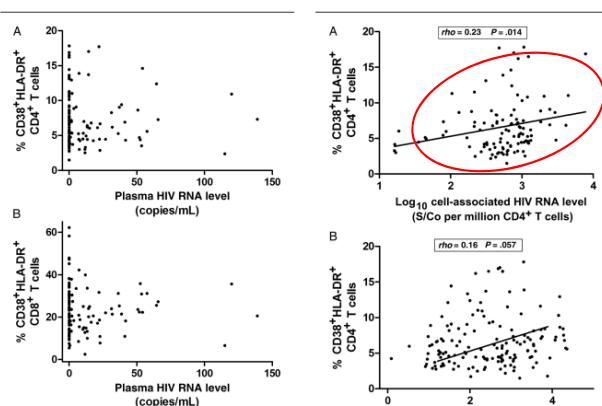


# Epidemiology

A)

**Table 1. Baseline Characteristics Among 190 Study Subjects**

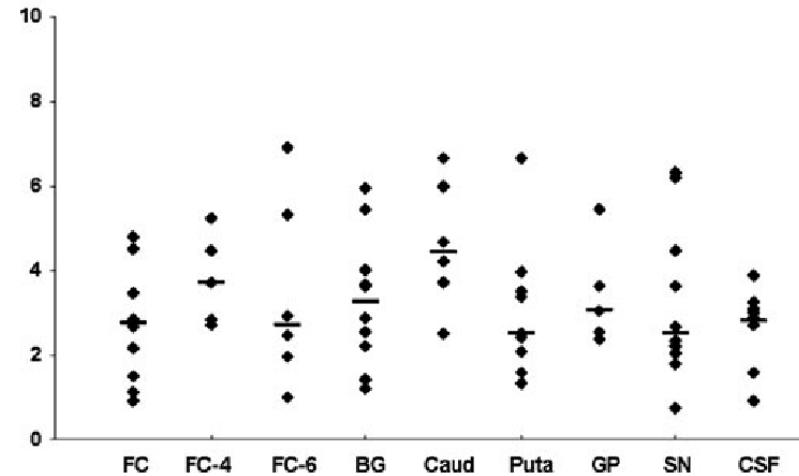
Characteristic	Value
Age, years	51 (44–57)
Male sex, % of patients	92
CD4 <sup>+</sup> T-cell count, cells/mm <sup>3</sup>	523 (249–728)
T cells expressing CD4, %	23 (16–32)
CD8 <sup>+</sup> T-cell count, cells/mm <sup>3</sup>	844 (606–1185)
T cells expressing CD8, %	48 (37–57)
Nadir CD4 <sup>+</sup> T-cell count, cells/mm <sup>3</sup>	113 (29–227)
Duration of HAART suppression, months	31 (14–66)
Agent included in HAART, no. of patients	
Protease inhibitor	124
NNRTI	73
Raltegravir	14
Maraviroc	1
Enfuvirtide	3



**Figure 1.** No associations between ultrasensitive plasma human immunodeficiency virus (HIV) RNA levels and T-cell activation. Ultrasensitive plasma HIV RNA levels were measured using the COBAS AmpliPrep/COBAS TaqMan HIV-1 test, version 2.0. All *P* values are >.40.

B)

**HIV-1 RNA in Human Brain Regions**

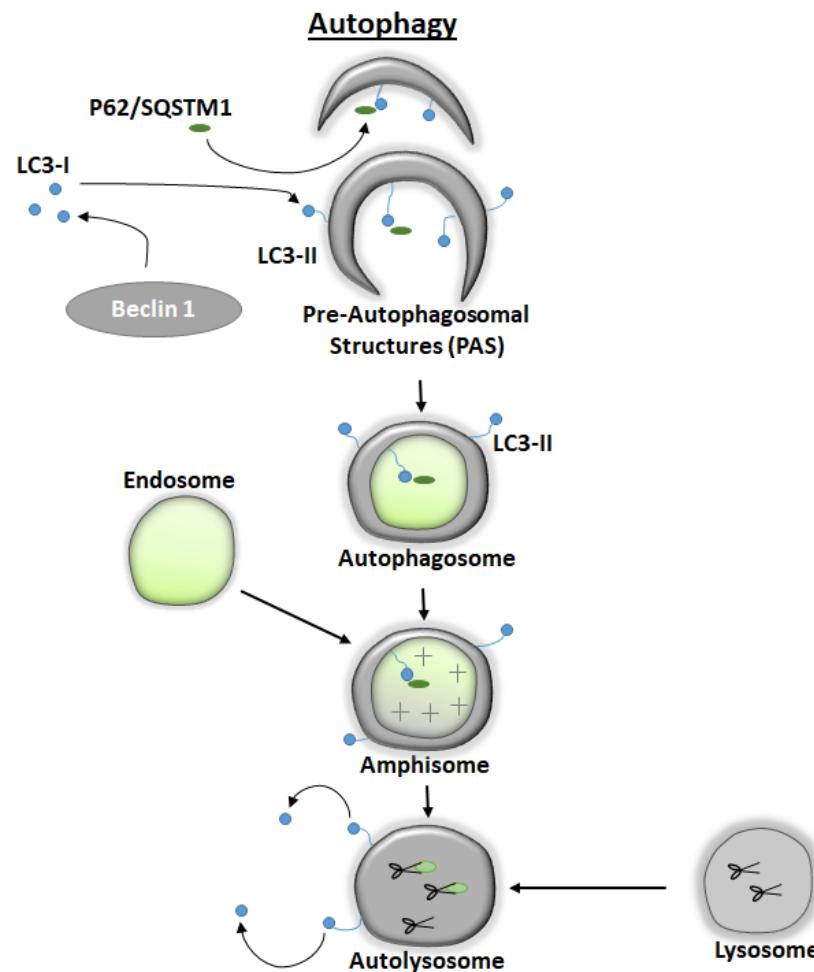


**Figure 1** HIV-1 RNA load ( $\log_{10}$  copies/g tissue) in each region of HIV-1+ individuals is presented in a scatter plot. The horizontal bars represent the median value in each region. Viral load values were higher in caudate nucleus compared to that in the other regions (FC-4, FC-6, BG, putamen, globus pallidus, SN), and the lowest values were found in FC and CSF.

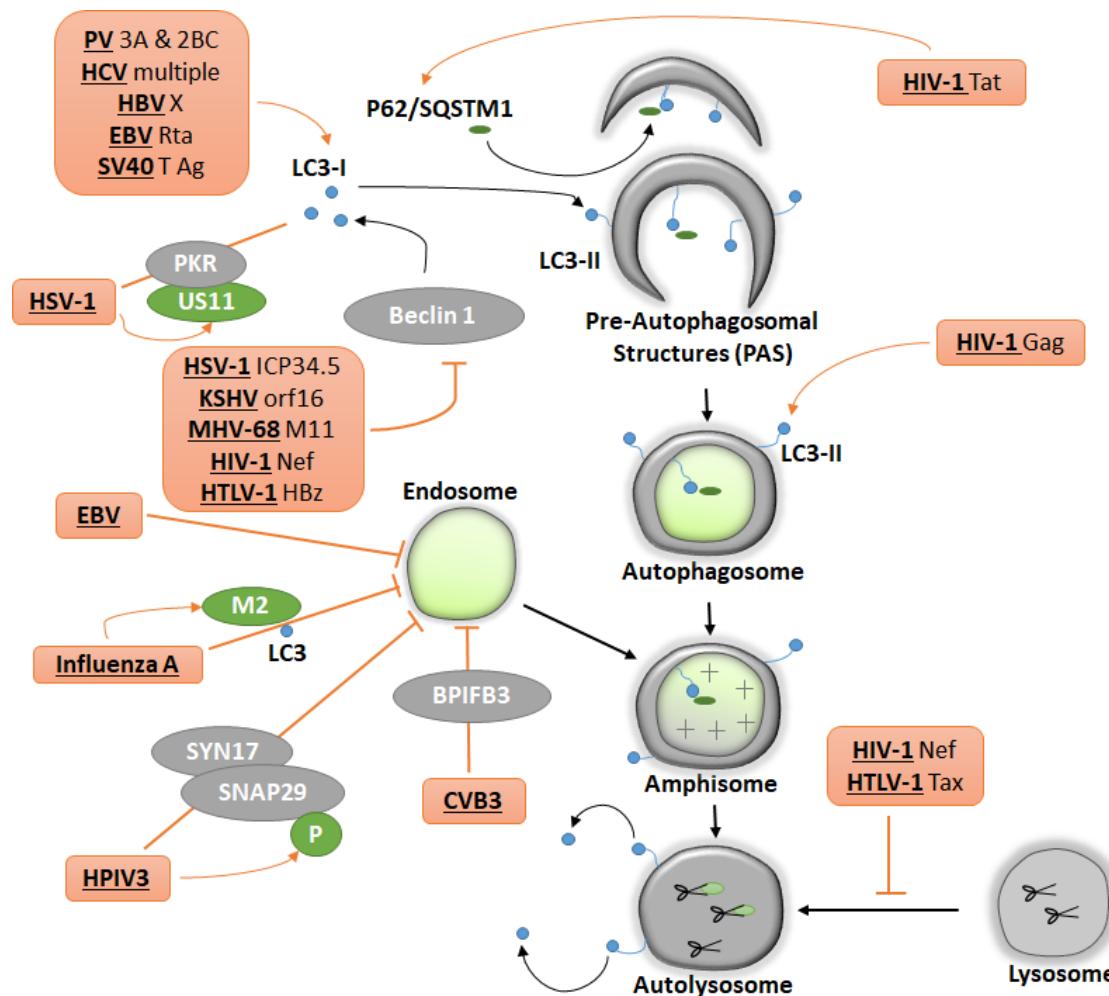
- A) Hatano et al. *JID*. 2013  
 B) Kumar et al. *J Neurovirol*. 2007



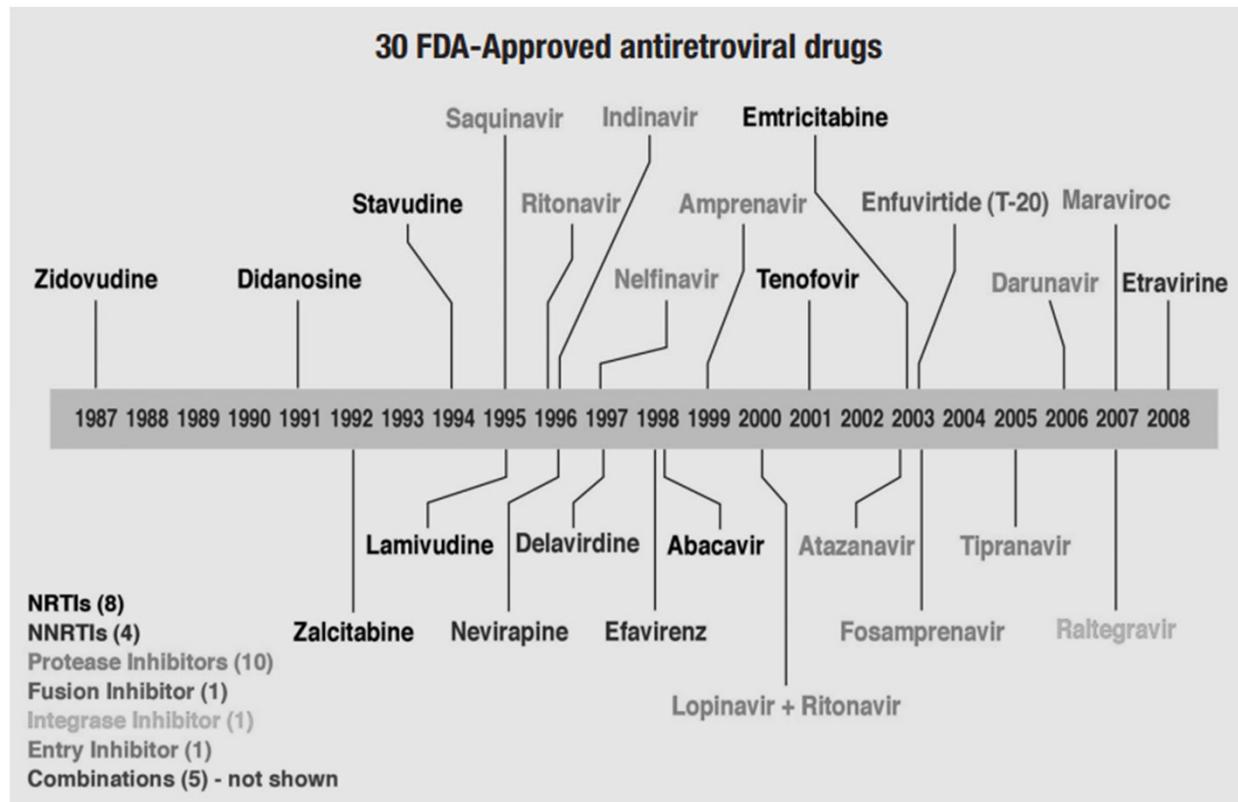
# Mechanism



# Mechanism



# Treatment



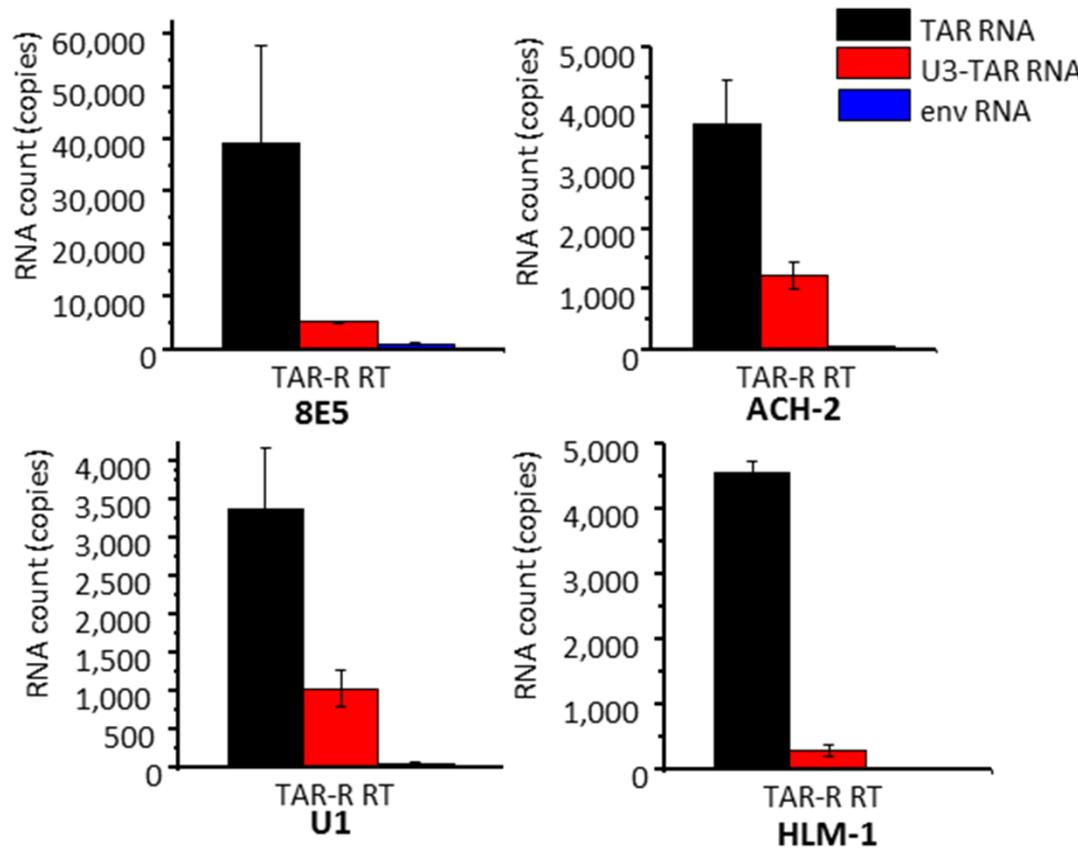
**No FDA-Approved Transcription Inhibitors**

Palmisano and Vella. Ann Ist Super Sanita. 2011



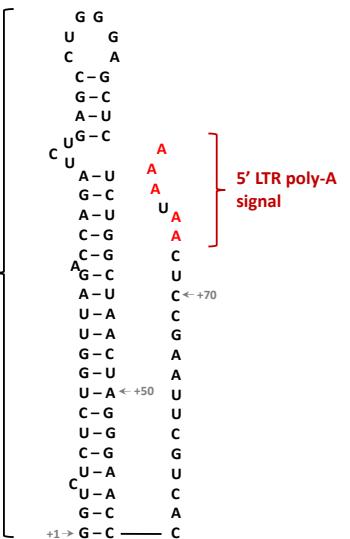
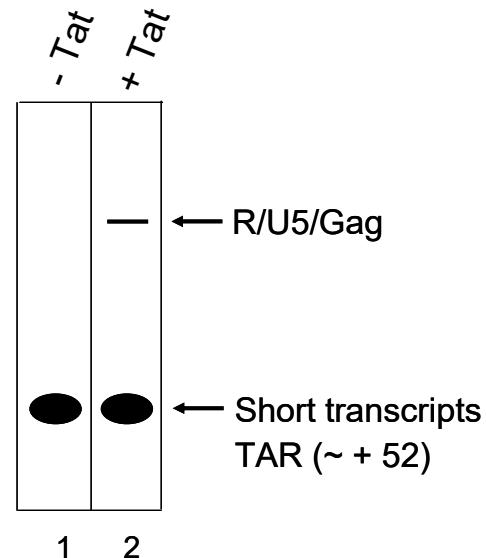
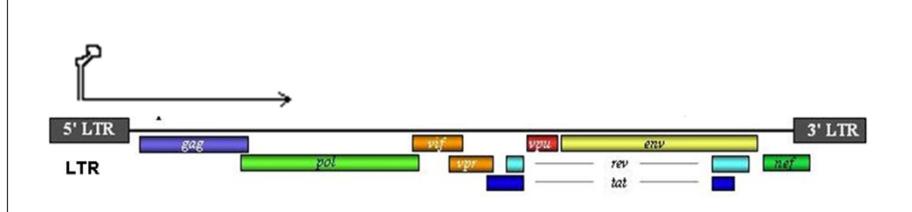
# Characterization and Analysis of “Damaging EVs”

## TAR RNA is present in culture supernatants of HIV-1 infected cells



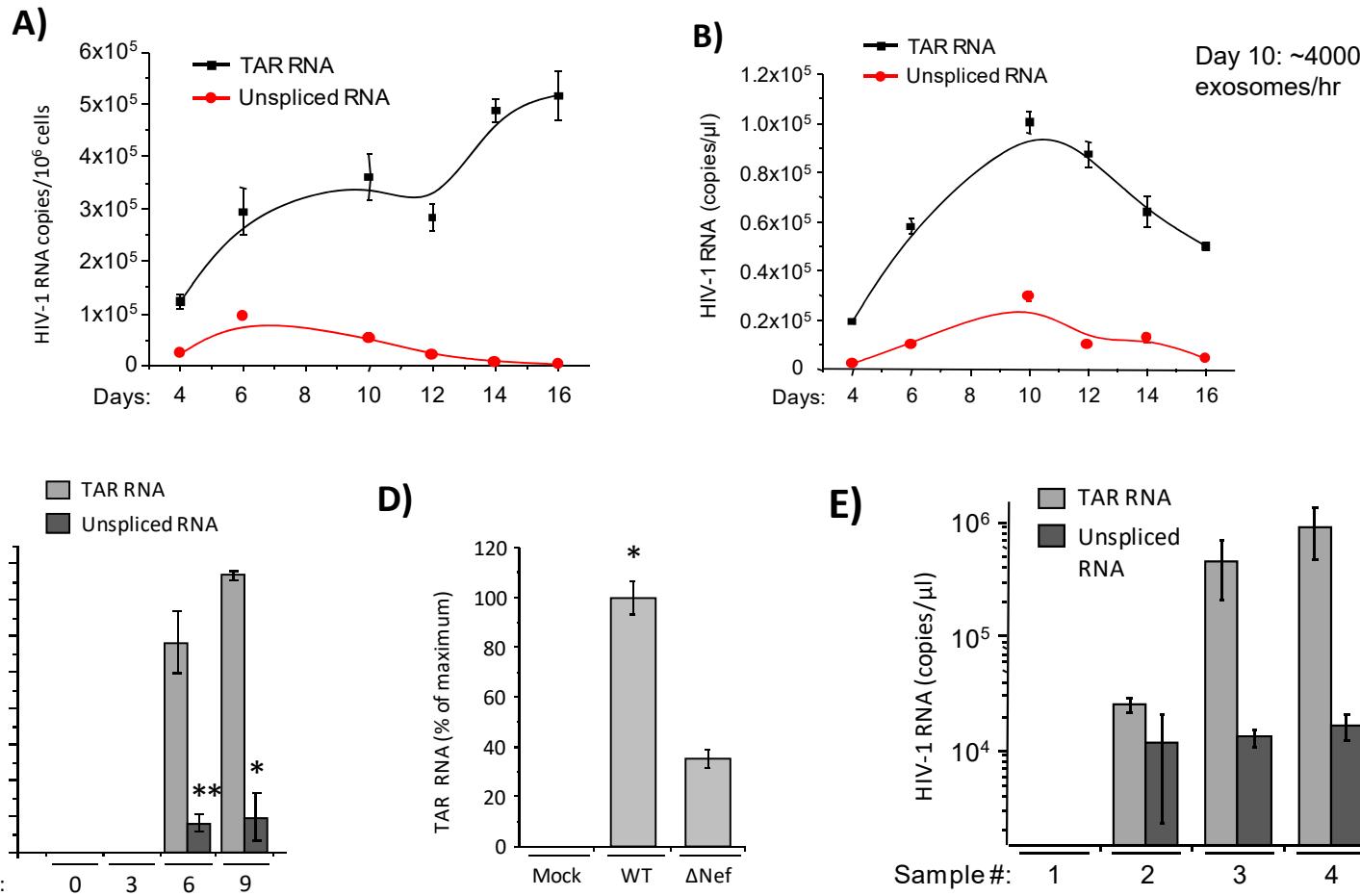
# Characterization and Analysis of “Damaging EVs”

## TAR RNA (short transcripts)

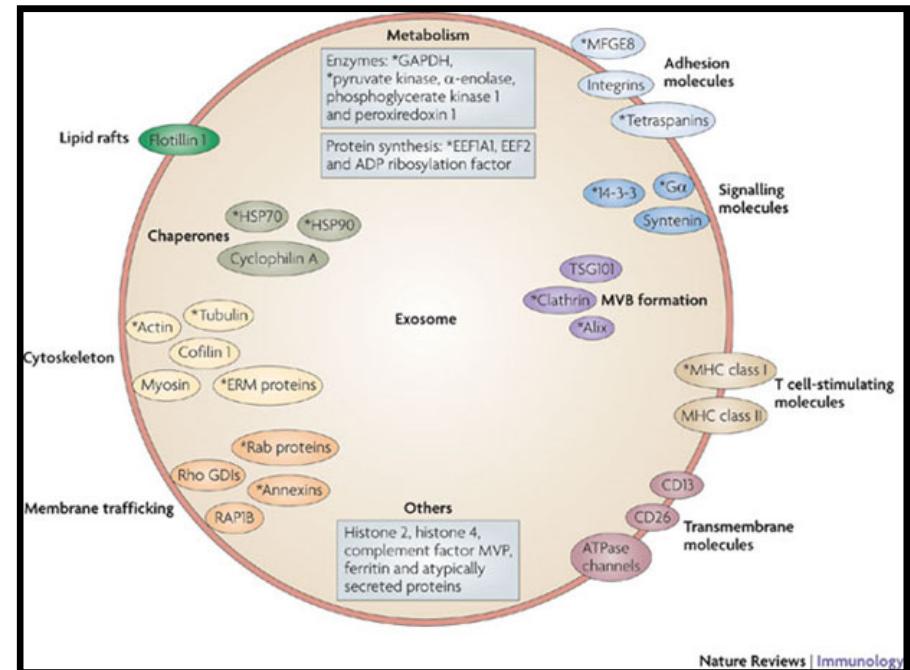
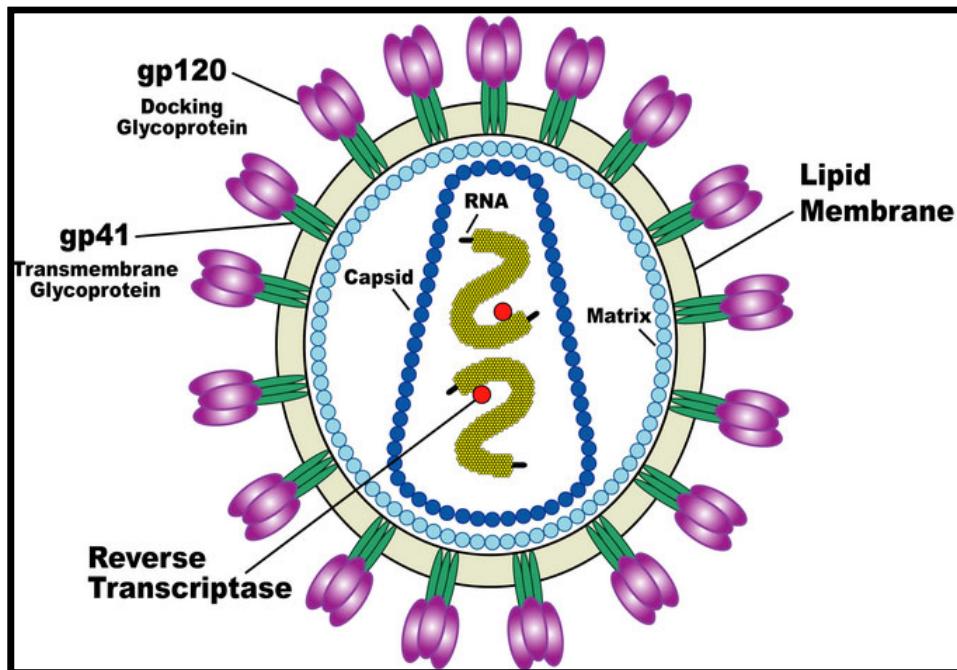


- More than 70 papers from late 80s to present
- Detected *in vitro*, *in vivo* and in latent patient cells

# Characterization and Analysis of “Damaging EVs”



# Characterization and Analysis of “Damaging EVs”



# Characterization and Analysis of “Damaging EVs”

## SCIENTIFIC REPORTS



Correction: Publisher Correction

OPEN

### Antiretroviral Drugs Alter the Content of Extracellular Vesicles from HIV-1-Infected Cells

Received: 3 January 2018

Accepted: 1 May 2018

Published online: 16 May 2018

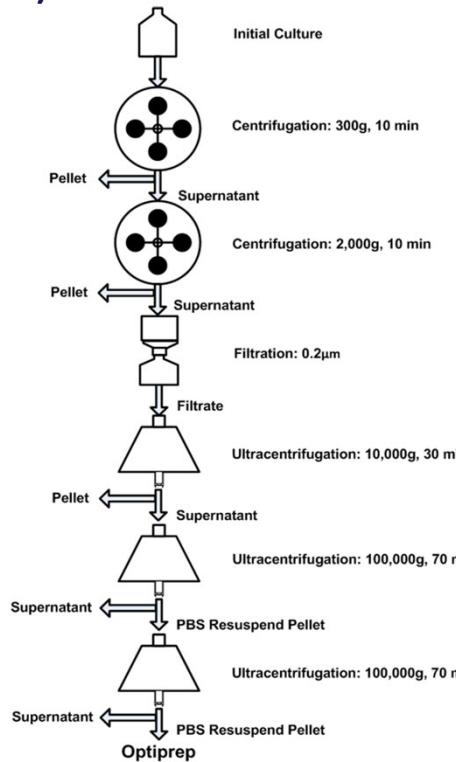
Catherine DeMarino<sup>1</sup>, Michelle L. Pleet<sup>1</sup>, Maria Cowen<sup>1</sup>, Robert A. Barclay<sup>1</sup>, Yao Akpamagbo<sup>1</sup>, James Erickson<sup>1</sup>, Nicaise Ndemb<sup>2</sup>, Manhattan Charurat<sup>2</sup>, Jibreel Jumare<sup>2</sup>, Sunday Bwala<sup>3</sup>, Peter Alabi<sup>4</sup>, Max Hogan<sup>5</sup>, Archana Gupta<sup>5</sup>, Nicole Noren Hooten<sup>1</sup>, Michele K. Evans<sup>6</sup>, Benjamin Lepene<sup>7</sup>, Weidong Zhou<sup>8</sup>, Massimo Caputi<sup>9</sup>, Fabio Romerio<sup>2</sup>, Walter Royal 3rd<sup>10</sup>, Nazira El-Hage<sup>11</sup>, Lance A. Liotta<sup>8</sup> & Fatah Kashanchi<sup>1</sup>

To date, the most effective treatment of HIV-1 is a combination antiretroviral therapy (cART), which reduces viral replication and reverses pathology. We investigated the effect of cART (RT and protease inhibitors) on the content of extracellular vesicles (EVs) released from HIV-1-infected cells. We have previously shown that EVs contain non-coding HIV-1 RNA, which can elicit responses in recipient cells. In this manuscript, we show that TAR RNA levels demonstrate little change with the addition of cART treatment in cell lines, primary macrophages, and patient biofluids. We determined possible mechanisms involved in the selective packaging of HIV-1 RNA into EVs, specifically an increase in EV-associated hnRNP A2/B1. More recent experiments have shown that several other FDA-approved drugs have the ability to alter the content of exosomes released from HIV-1-infected cells. These findings on cART-altered EV content can also be applied to general viral inhibitors (interferons) which are used to treat other chronic infections. Additionally, we describe unique mechanisms of ESCRT pathway manipulation by antivirals, specifically the targeting of VPS4. Collectively, these data imply that, despite antiretroviral therapy, EVs containing viral products are continually released and may cause neurocognitive and immunological dysfunction.

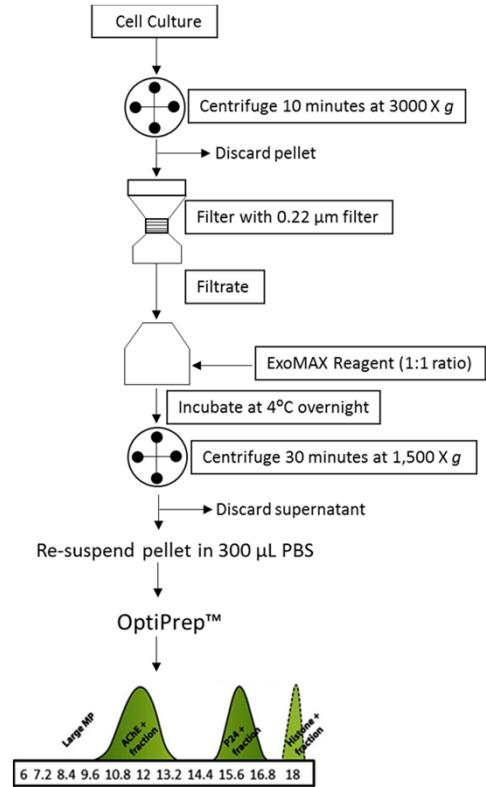


# Isolation

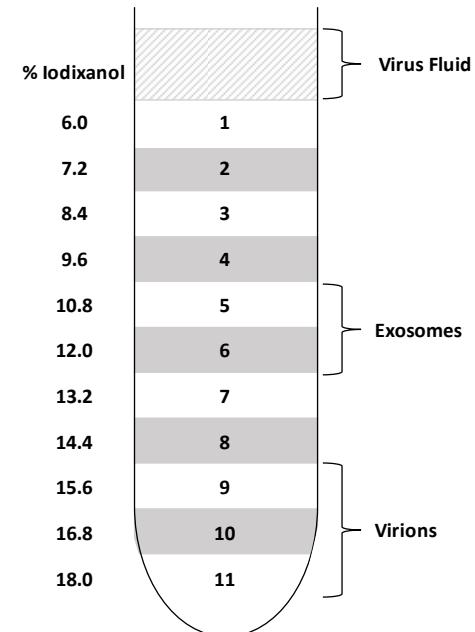
A)



B)

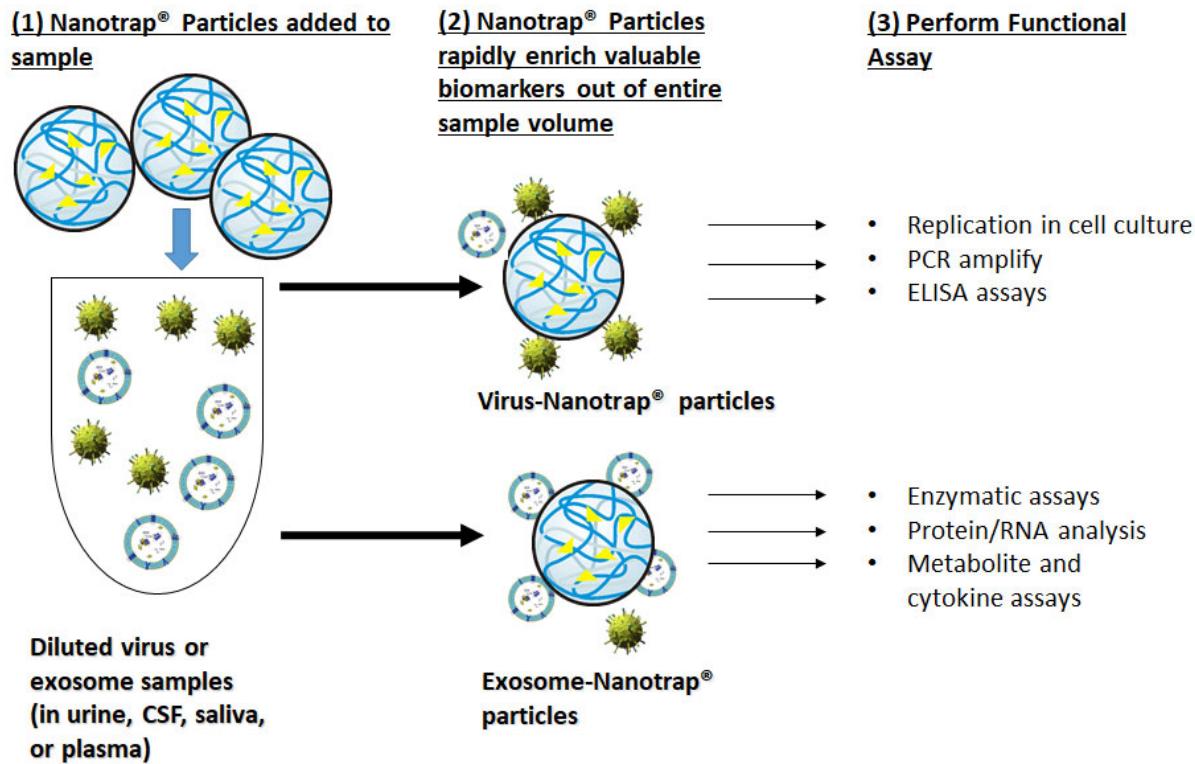


C)



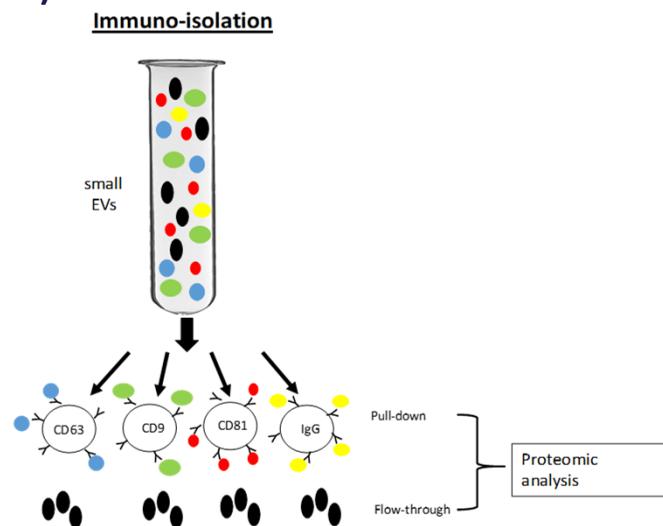
# Characterization and Analysis of “Damaging EVs”

## Using Nanotrap® particles to capture exosomes or viruses



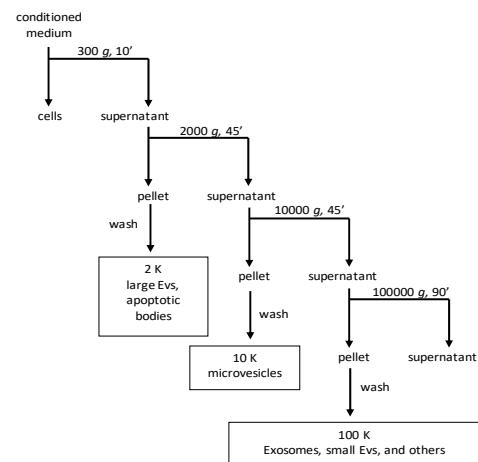
# Characterization and Analysis of “Damaging EVs”

A)

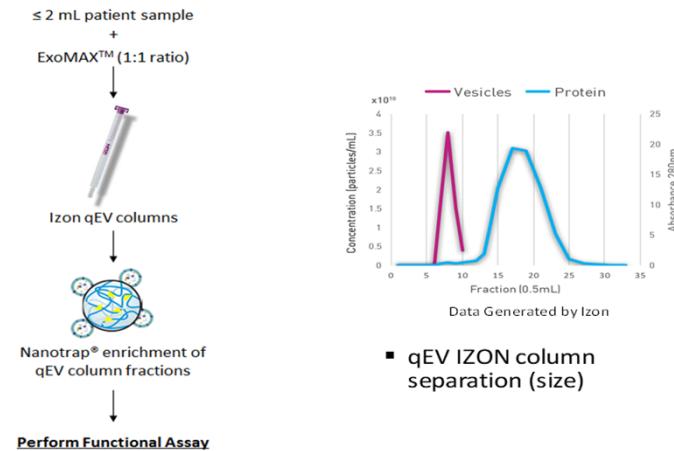


B)

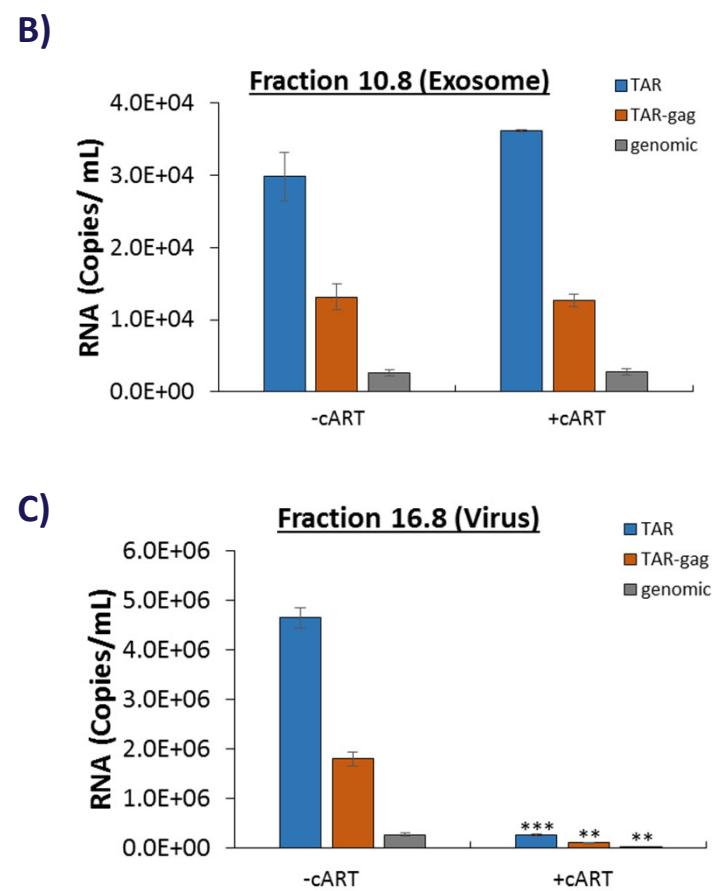
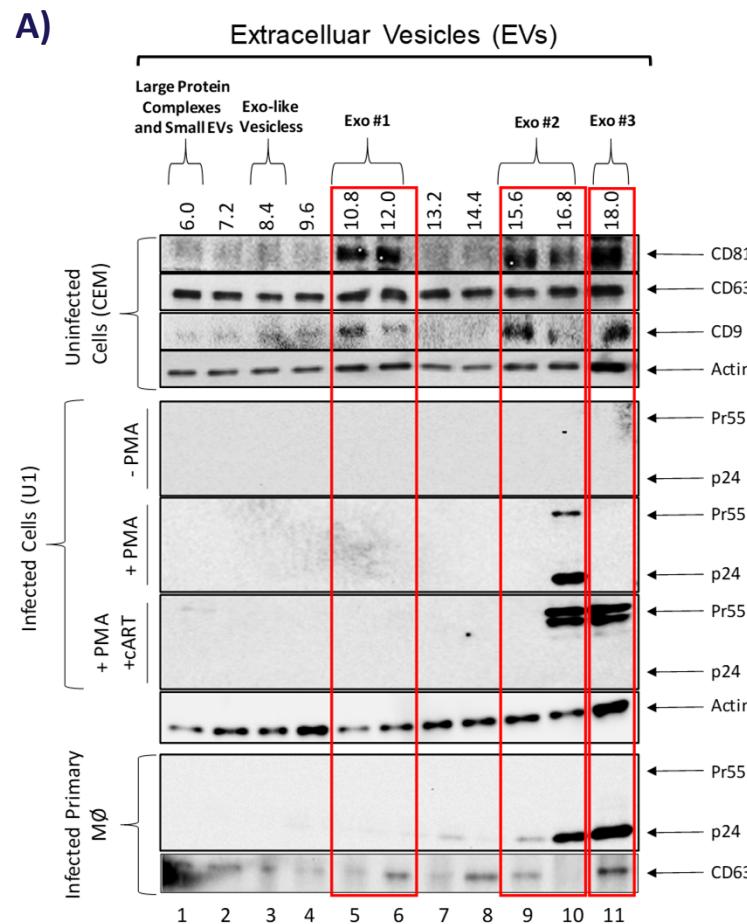
## Ultracentrifugation only for small samples



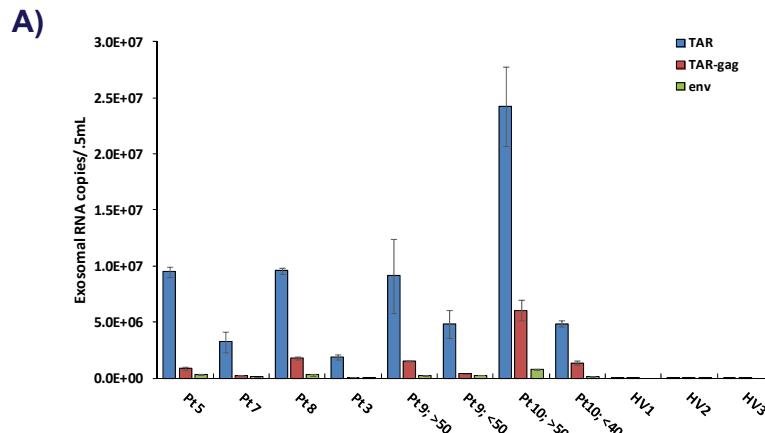
C)



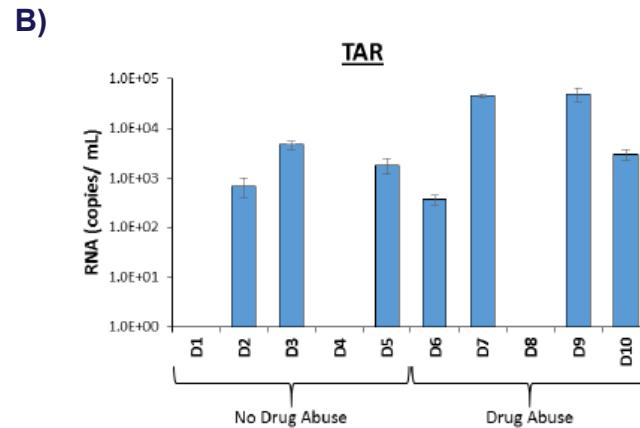
# Characterization and Analysis of “Damaging EVs”



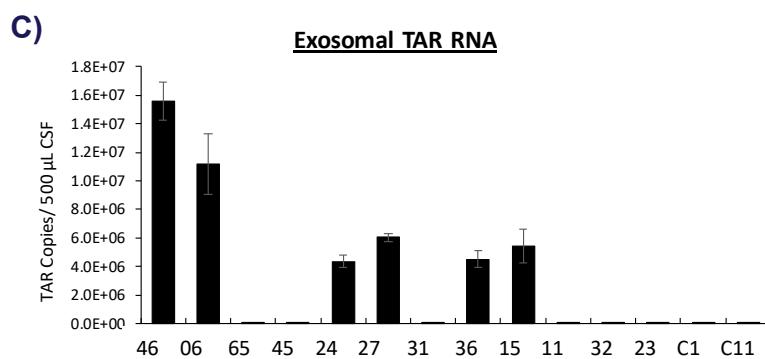
# Characterization and Analysis of “Damaging EVs”



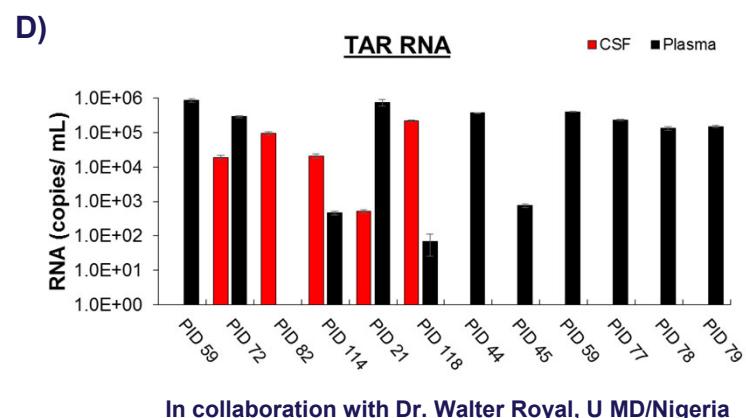
In collaboration with Dr. Cliff Lane's Lab, NIH



In collaboration with Dr. Nikki Noren Hooten, NIA



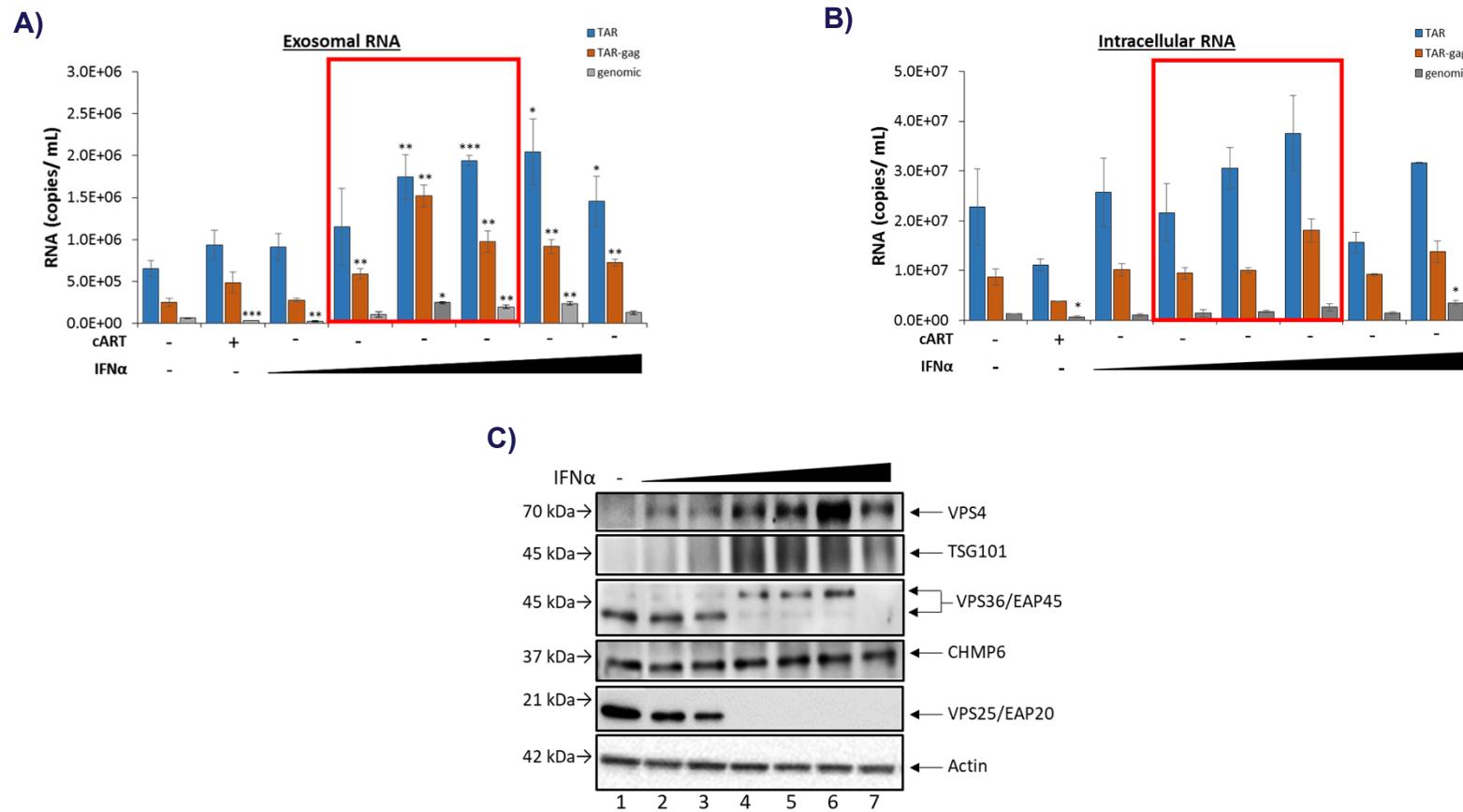
In collaboration with Dr. Avi Nath's Lab, NIH



In collaboration with Dr. Walter Royal, U MD/Nigeria

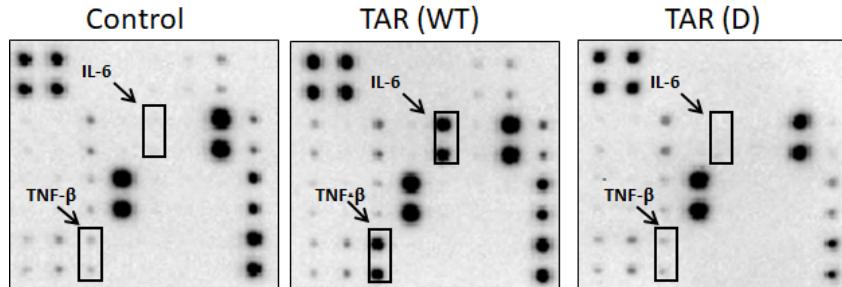
# Characterization and Analysis of “Damaging EVs”

## Interferon Alters EV Content Through the ESCRT Pathway

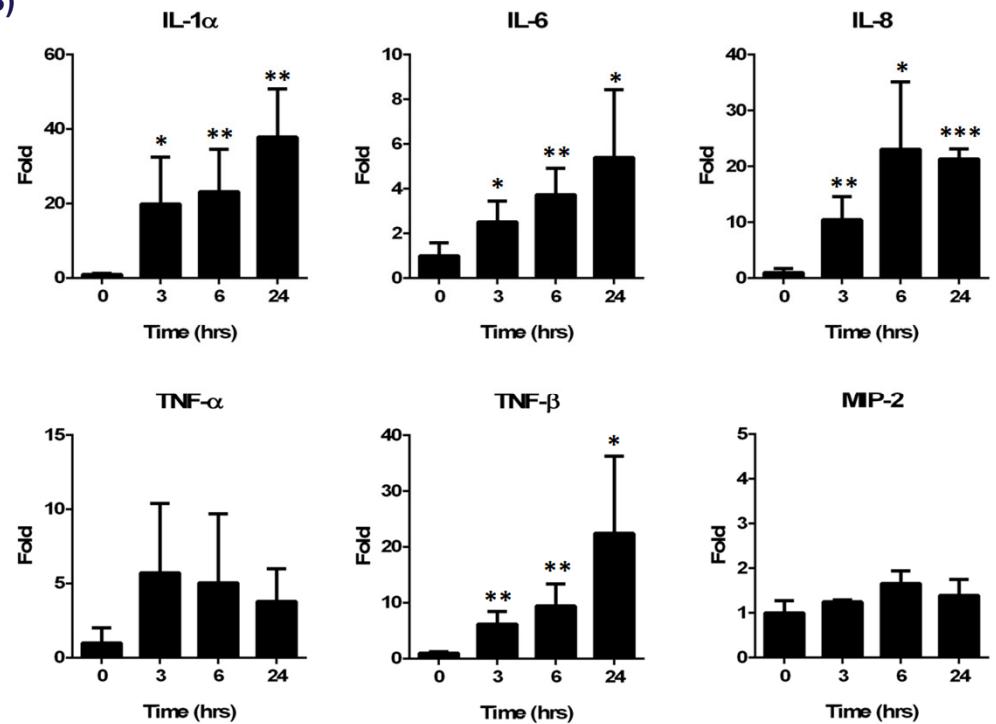


# Characterization and Analysis of “Damaging EVs”

A)



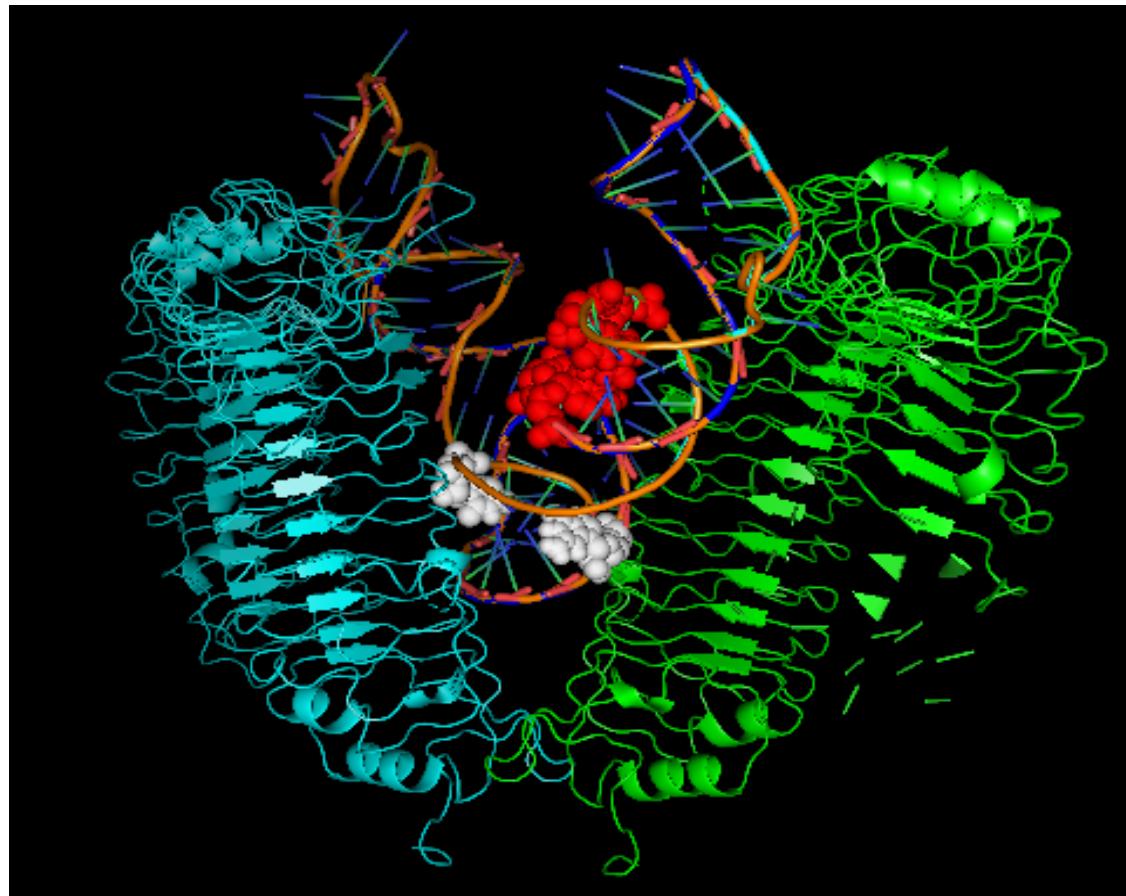
B)



\*P<0.05, \*\*P<0.01, \*\*\*P<0.001 vs. 0 h control

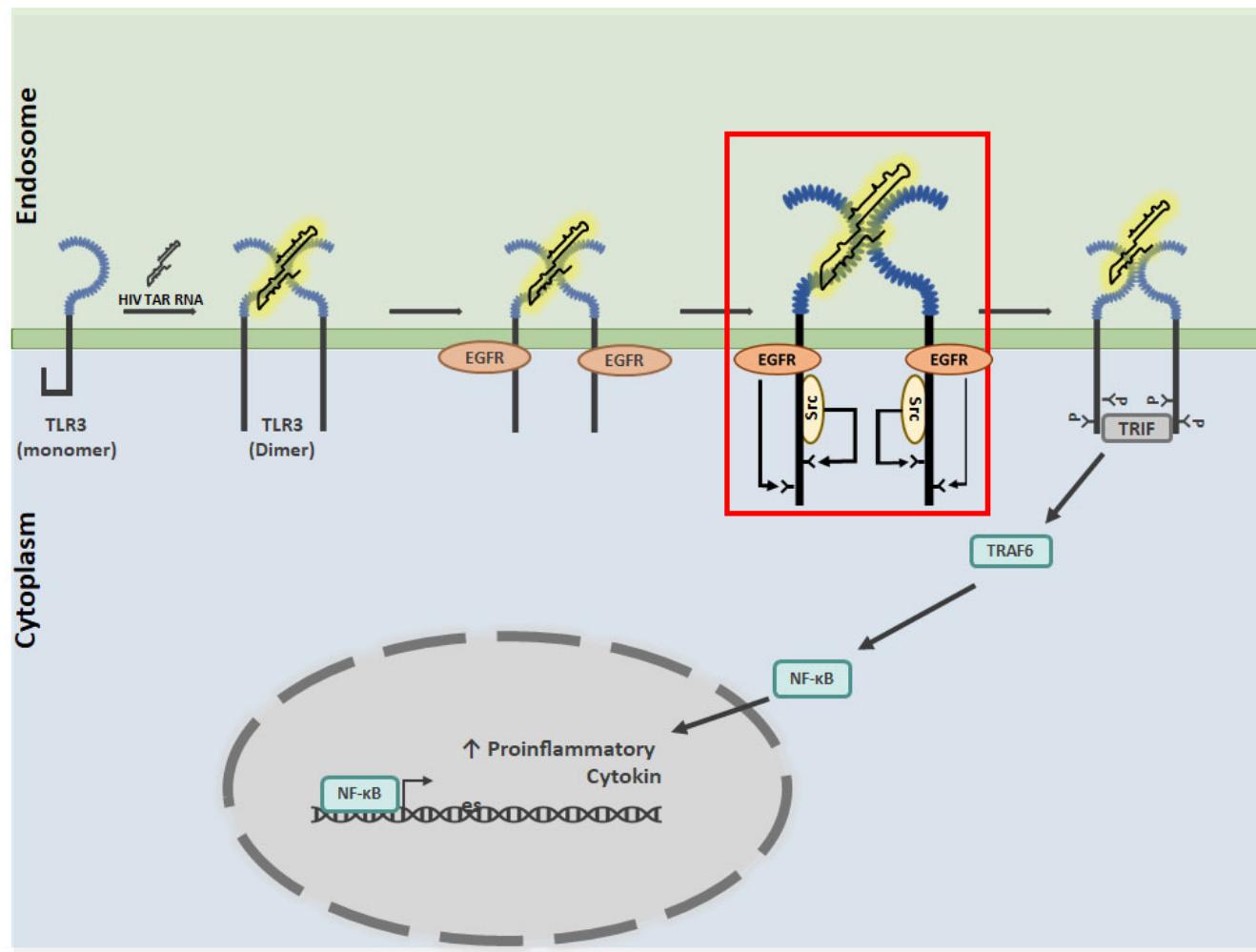


# Characterization and Analysis of “Damaging EVs”

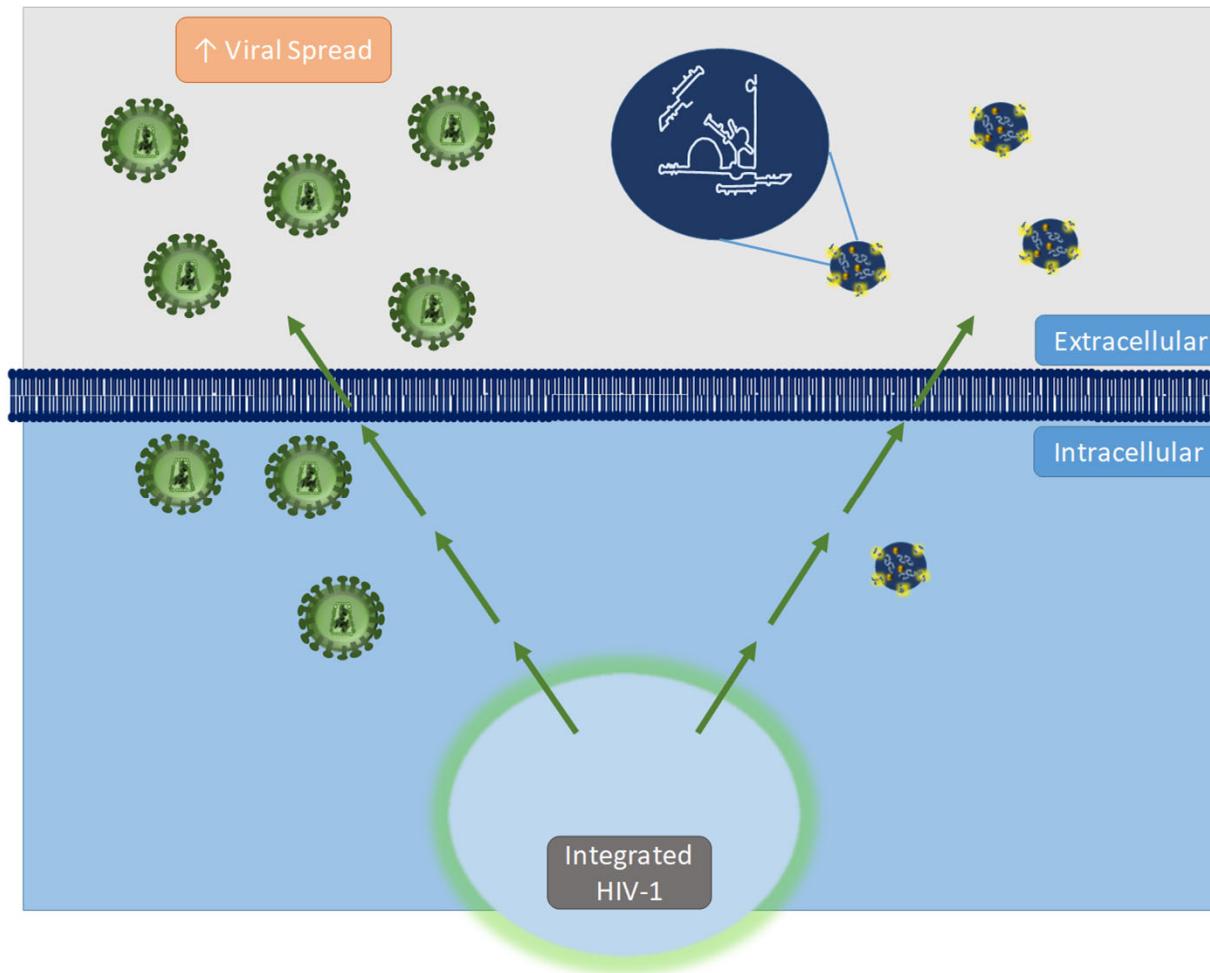


TAR RNA: TLR3 signaling complex.

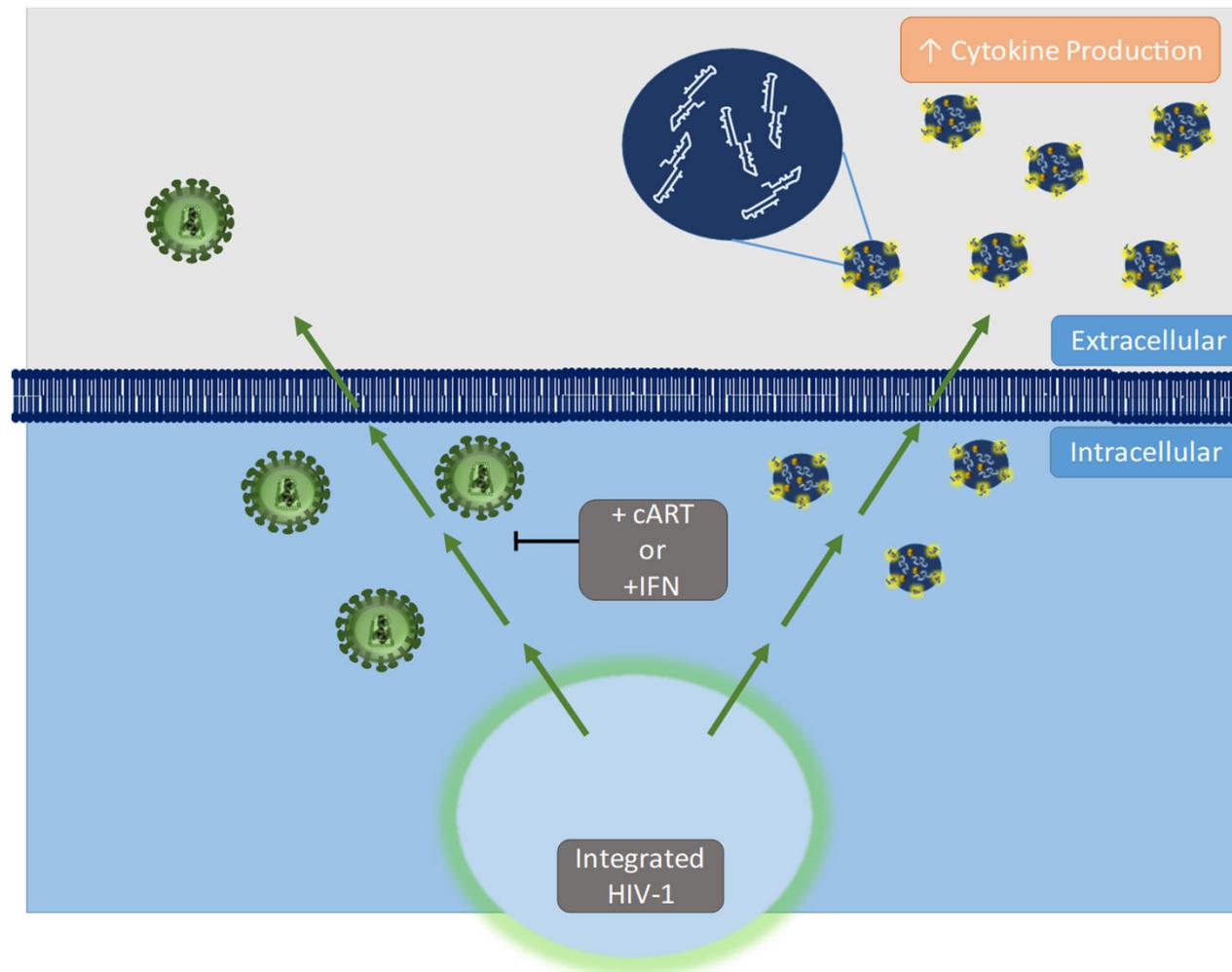
# Mechanism



# Mechanism



# Mechanism



# Overview

---

- Overview of Extracellular Vesicles (EVs)
- Characterization and Analysis of “Damaging EVs”
- **Characterization and Analysis of “Reparative EVs”**
- Summary and Future Directions

# Topics to Be Discussed

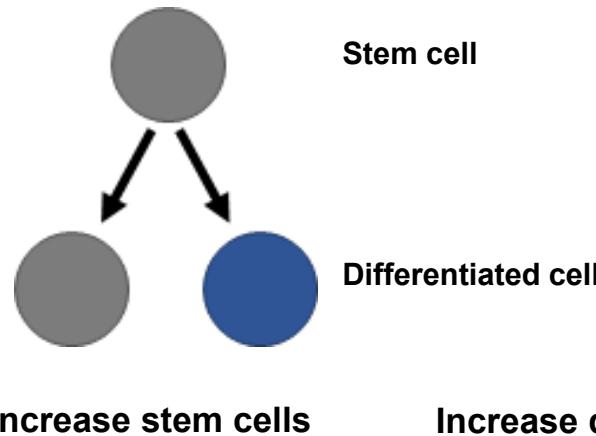
---

1. Background on stem cells and stem cell EVs  
**(reparative EVs)**
  
2. Functional effects of stem cell EVs
  - Wound healing and skin repair
  - Cardiac, eye repair
  - CNS repair (stroke, TBI, SCI, neurodegenerative disorders)

# Stem Cells

- **Defining characteristics**
  - Self-renewal
  - Differentiation potential
- **Types of stem cells**
  - Embryonic stem cells
  - Adult (somatic) stem cells
  - Induced pluripotent stem cells

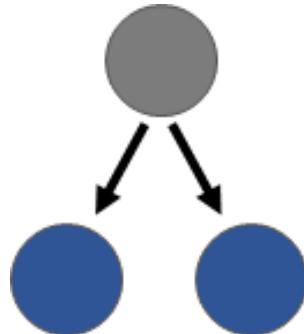
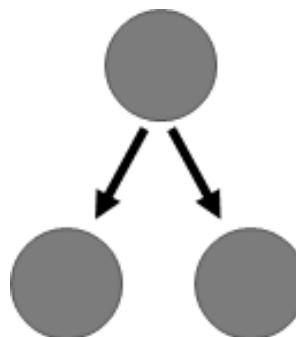
Maintain stem cell population



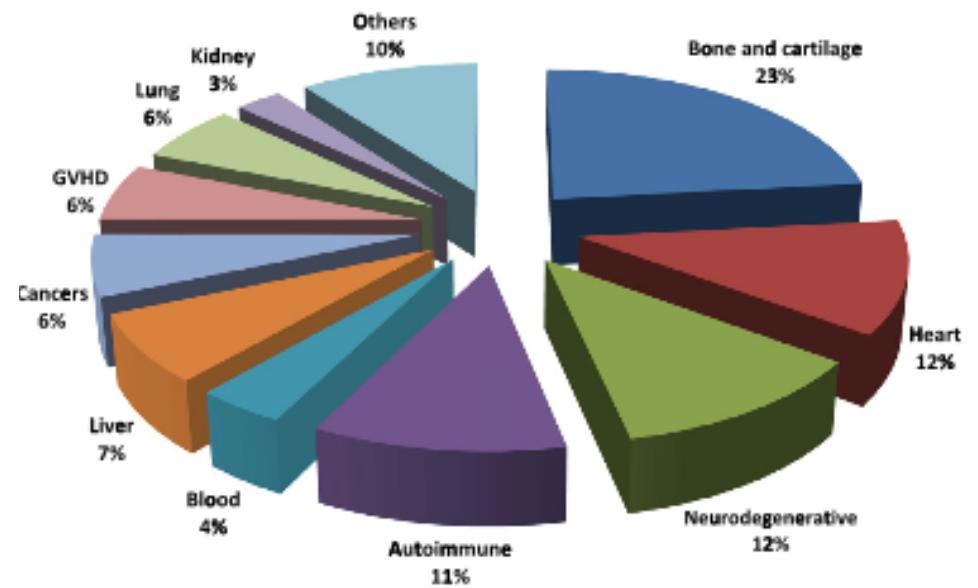
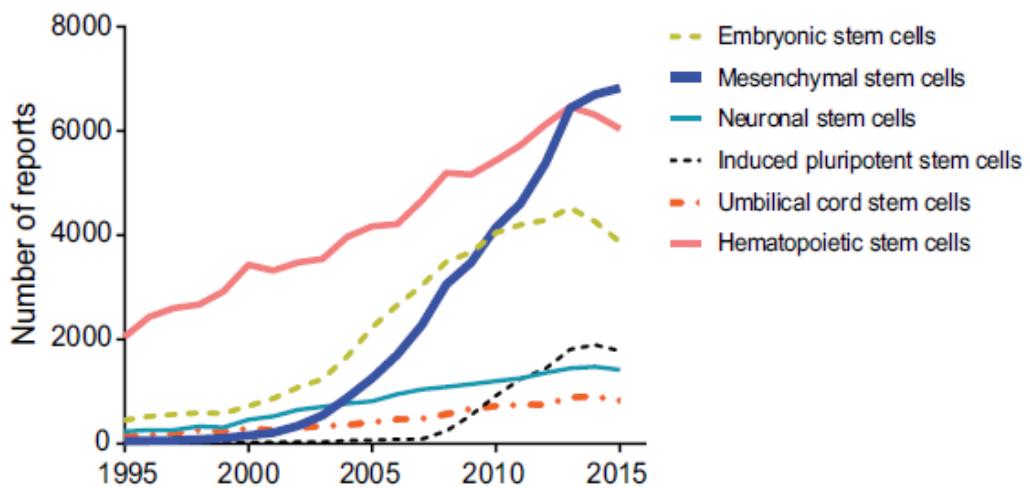
Stem cell

Increase stem cells

Increase differentiating cells



# Stem Cell Therapy



Samsonraj et al. *Stem Cells Translational Medicine*. 2017  
Marquez-Curtis et al. *Cryobiology*. 2015

# Stem Cell EVs

---

- Contain various biological cargo (miRNAs, lncRNAs, proteins, cytokines) that can be transferred to recipient cells
- Proposed to play a role in homeostasis through tissue repair, regeneration, and immunomodulation
- **Potential alternative to stem cell therapy due to higher potency, increased stability/shelf life, and lower immunogenicity**
- Widely studied in the context of skin/wound healing, angiogenesis, cardiac repair, and CNS-related pathologies

# Stem Cell EVs

Stem Cell Research & Therapy

RESEARCH

Open Access



Exosomes from acellular Wharton's jelly of the human umbilical cord promotes skin wound healing

Nazihah Bakhtyar<sup>1</sup>, Marc G. Jeschke<sup>1,2\*</sup>, Elaine Herer<sup>1,5</sup>, Mohammadali Sheikholeslam<sup>1,3</sup> and Saeid Amini-Nik<sup>1,3,4\*</sup>

**WJSC** World Journal of Stem Cells

Submit a Manuscript: <http://www.ifpublishing.com>  
DOI: 10.4237/wjsc.v10i8106  
ISSN 1946-0210 (online)  
REVIEW

Stem cell-derived exosomes - an emerging tool for myocardial regeneration

Erzsebet Lazar, Theodora Benedek, Szilámer Korodi, Nora Rat, Jocelyn Lo, Imre Benedek

J Trauma Acute Care Surg. 2019 Dec 5. doi: 10.1097/TA.0000000000002593. [Epub ahead of print]

**Early Single-Dose Treatment with Exosomes Provides Neuroprotection and Improves Blood-Brain Barrier Integrity in Swine Model of Traumatic Brain Injury and Hemorrhagic Shock.**

Williams AM<sup>1</sup>, Bhatti UF<sup>1</sup>, Brown JF<sup>1</sup>, Biesterveld BE<sup>1</sup>, Kathawate RG<sup>1</sup>, Graham NJ<sup>1</sup>, Chtraklin K<sup>1</sup>, Siddiqui AZ<sup>1</sup>, Dekker SE<sup>1</sup>, Andjelkovic A<sup>2</sup>, Higgins GA<sup>3</sup>, Buller B<sup>1</sup>, Alam HB<sup>1</sup>.

Neural Regen Res. 2019 Sep; 14(9): 1626–1634.  
doi: [10.4103/1673-5374.255978](https://doi.org/10.4103/1673-5374.255978)

PMCID: PMC6557105  
PMID: 31089063

Mesenchymal stem cell-derived exosomes promote neurogenesis and cognitive function recovery in a mouse model of Alzheimer's disease

Edwin E. Reza-Zaldivar,<sup>1</sup> Mercedes A. Hernández-Sapiens,<sup>1</sup> Yanet K. Gutiérrez-Mercado,<sup>1</sup> Sergio Sandoval-Ávila,<sup>1</sup> Ulises Gomez-Pinedo,<sup>2</sup> Ana L. Márquez-Aquirre,<sup>1</sup> Estefanía Vázquez-Méndez,<sup>1</sup> Eduardo Padilla-Camberos,<sup>1</sup> and Alejandro A. Canales-Aquirre<sup>1</sup>

Proc Natl Acad Sci U.S.A. 2017 Apr 25; 114(17): E3536–E3545.  
Published online 2017 Apr 10. doi: [10.1073/pnas.1703920114](https://doi.org/10.1073/pnas.1703920114)

PNAS Plus  
Neuroscience

PMCID: PMC5410779  
PMID: 28396435

Intranasal MSC-derived A1-exosomes ease inflammation, and prevent abnormal neurogenesis and memory dysfunction after status epilepticus

Qianfa Long,<sup>a,1,2</sup> Dinesh Upadhyaya,<sup>a,b,c,2,3</sup> Bharathi Hatiangady,<sup>a,b,c</sup> Dong-Ki Kim,<sup>a</sup> Su Yeon An,<sup>a</sup> Bing Shuai,<sup>a,b,c</sup> Darwin J. Prockop,<sup>a,c,d,4,5</sup> and Ashok K. Shetty,<sup>a,b,c,4,5</sup>

Hindawi  
Stem Cells International  
Volume 2019, Article ID 5738510, 9 pages  
<https://doi.org/10.1155/2019/5738510>

Research Article

**Mesenchymal Stem Cell-Derived Extracellular Vesicles for Corneal Wound Repair**

Hongyan Tao,<sup>1,2</sup> Xiaoniao Chen,<sup>2</sup> Hongmei Cao,<sup>1</sup> Lingyue Zheng,<sup>3</sup> Qian Li,<sup>2</sup> Kaiyue Zhang,<sup>1</sup> Zhibo Han,<sup>4,5</sup> Zhong-Chao Han,<sup>4,5</sup> Zhikun Guo,<sup>6</sup> Zongjin Li,<sup>7</sup> and Liqiang Wang<sup>2</sup>

Contents lists available at ScienceDirect  
 Materials Science & Engineering C  
journal homepage: [www.elsevier.com/locate/msec](http://www.elsevier.com/locate/msec)

hucMSC derived exosomes promote functional recovery in spinal cord injury mice via attenuating inflammation

Guodong Sun<sup>a,1</sup>, Guangqiang Li<sup>a,1</sup>, Dehai Li<sup>a</sup>, Wanjun Huang<sup>b</sup>, Renwen Zhang<sup>b</sup>, Hua Zhang<sup>b</sup>, Yuanyuan Duan<sup>a,b</sup>, Baocheng Wang<sup>a,b</sup>

Journal of Neuroimmune Pharmacology  
<https://doi.org/10.1007/s11481-019-09865-y>

ORIGINAL ARTICLE



**Stem Cell Extracellular Vesicles and their Potential to Contribute to the Repair of Damaged CNS Cells**

Heather Branscome,<sup>1,2</sup> Siddhartha Paul<sup>3</sup>, Pooja Khatkar<sup>1</sup>, Yuriy Kim<sup>1</sup>, Robert A. Barday<sup>1</sup>, Daniel O. Pinto<sup>1</sup>, Dezhong Yin<sup>2</sup>, Weidong Zhou<sup>4</sup>, Lance A. Liotta<sup>4</sup>, Nadia El-Hage<sup>5</sup>, Fatah Kashanchi<sup>1</sup>



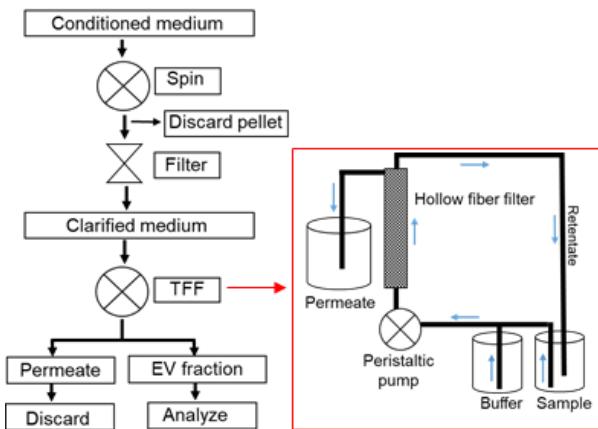
# EVs and EV Donor Cells

- **Mesenchymal Stem Cells (MSCs; ATCC® PCS-500-012™)**
  - Human, normal
  - Bone-marrow derived
  - Authenticated for characteristic surface marker expression (CD90, CD73, CD105 positive; CD14, CD34, CD45 negative)
  - Multi-lineage differentiation potential (adipocyte, chondrocyte, osteocyte)
- **Induced Pluripotent Stem Cells (iPSCs; ATCC® ACS-1019™)**
  - Human, normal
  - Foreskin fibroblast-derived
  - Sendai virus reprogrammed
  - Authenticated for expression of stem cell markers (TRA-1-60, SSEA-4 positive; SSEA-1 negative)
  - Evaluated for pluripotency
- **A549 Lung Carcinoma (ATCC® CCL-185™)**
  - Control used for large-scale manufacturing and isolation
  - Equivalent to CCL-185-EXM™ (also: ATCC® SCRC-4000-EXM™, CRL-1435-EXM™, CCL-247-EXM™)

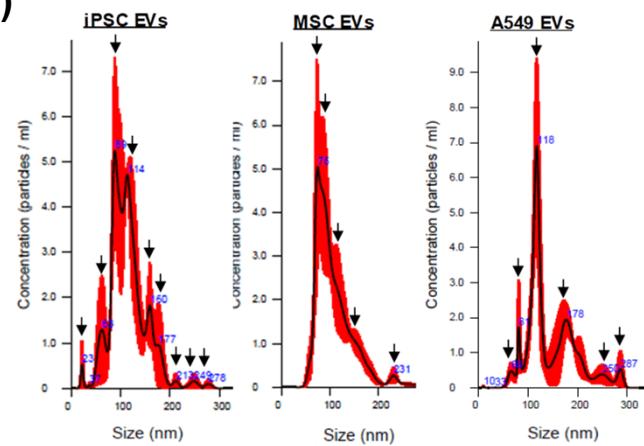


# EV Isolation and Characterization

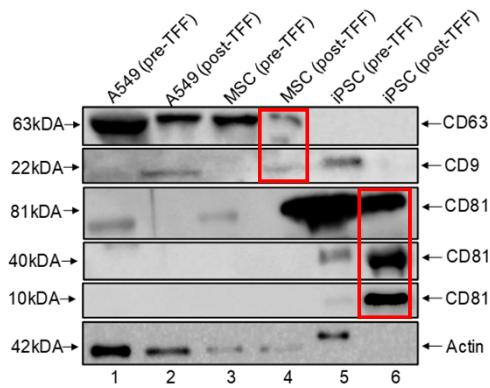
**A)**



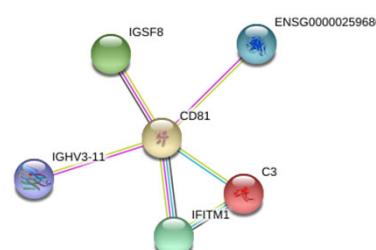
**B)**



**C)**

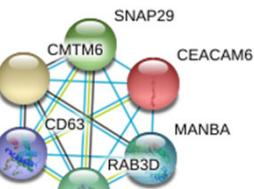


**D)**



**CD81**

**CD63**

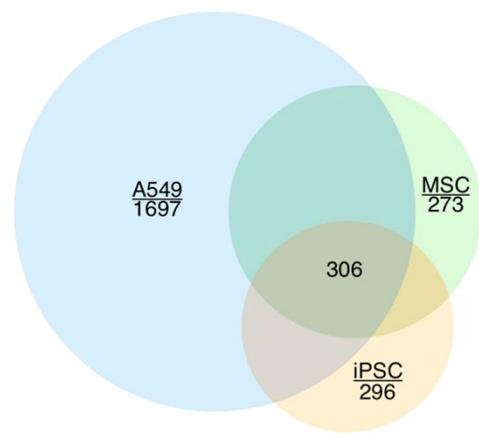


GeneCards.org

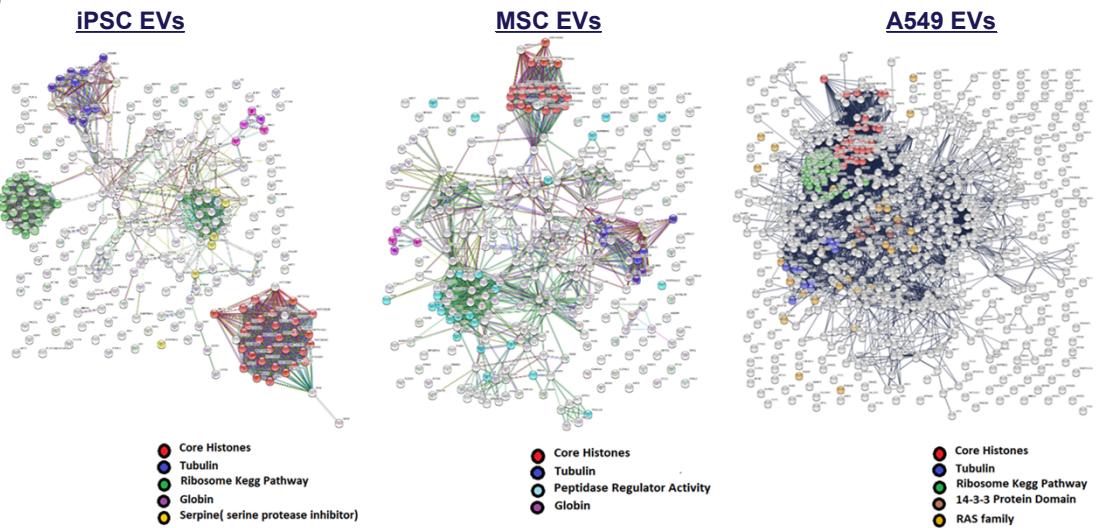


# EV Characterization

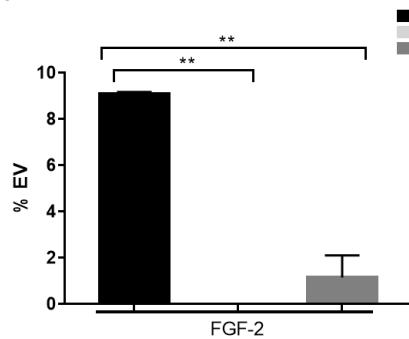
A)



B)



C)



D)

SCIENTIFIC REPORTS

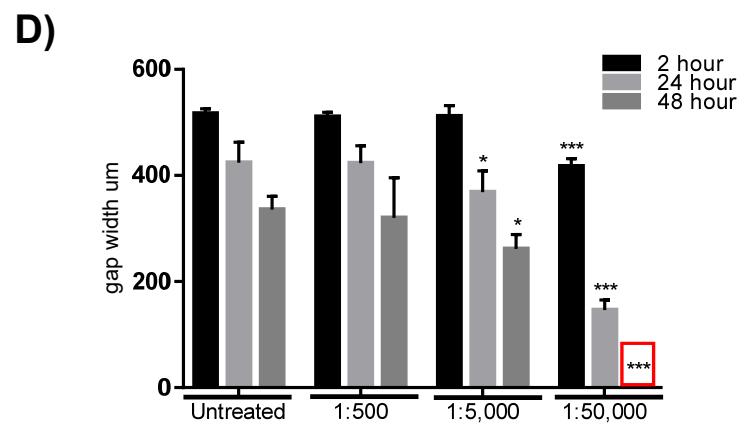
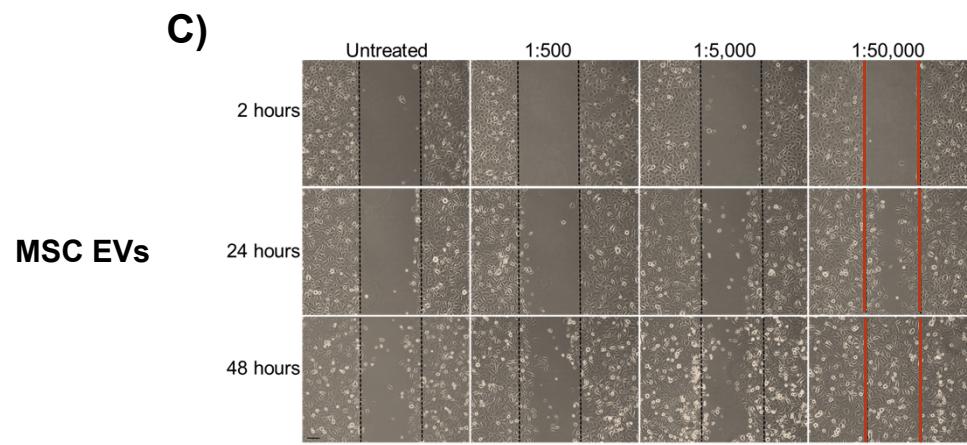
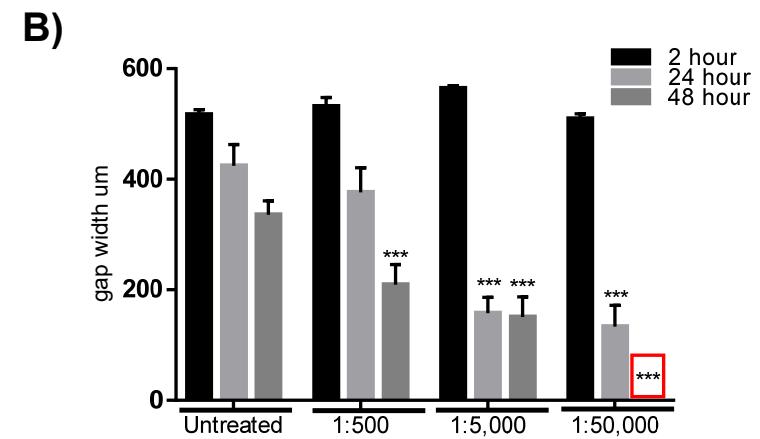
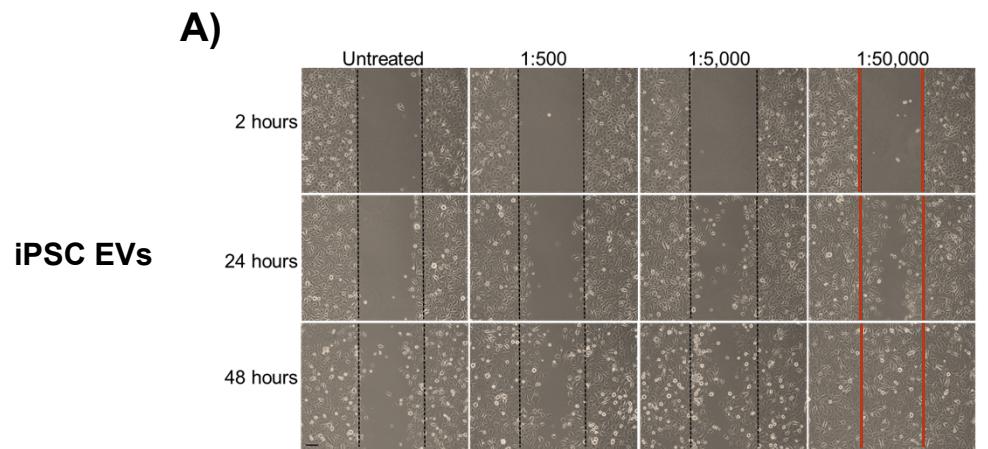
OPEN | A System of Cytokines Encapsulated in ExtraCellular Vesicles

Received: 25 January 2018  
Accepted: 14 May 2018  
Published online: 12 June 2018

Wendy Fitzgerald<sup>1</sup>, Michael L. Freeman<sup>2</sup>, Michael M. Lederman<sup>3</sup>, Elena Vasilieva<sup>4</sup>, Roberto Romero<sup>5</sup> & Leonid Margolis<sup>2</sup>



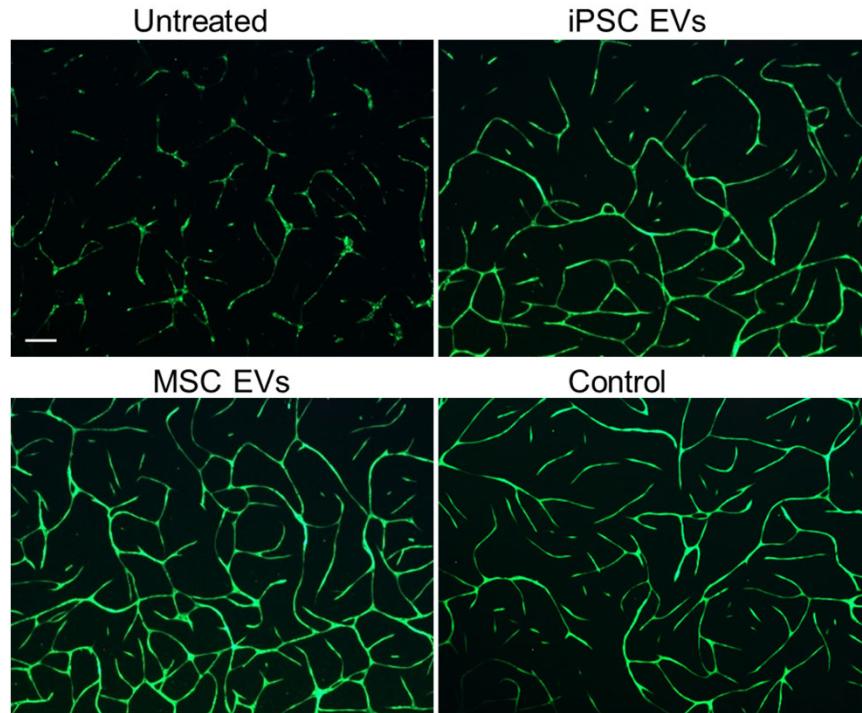
# EV Functionality



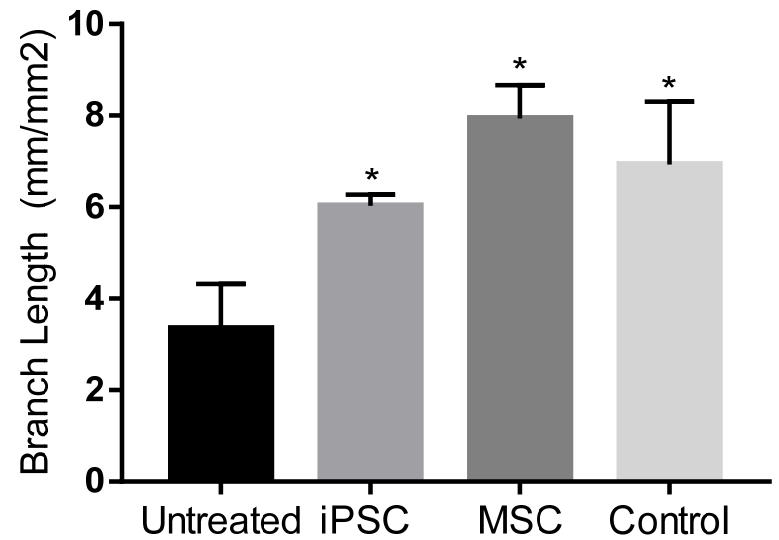
# EV Functionality

Angio-**Ready™** Angiogenesis Assay System (ATCC® ACS-2001-2™)

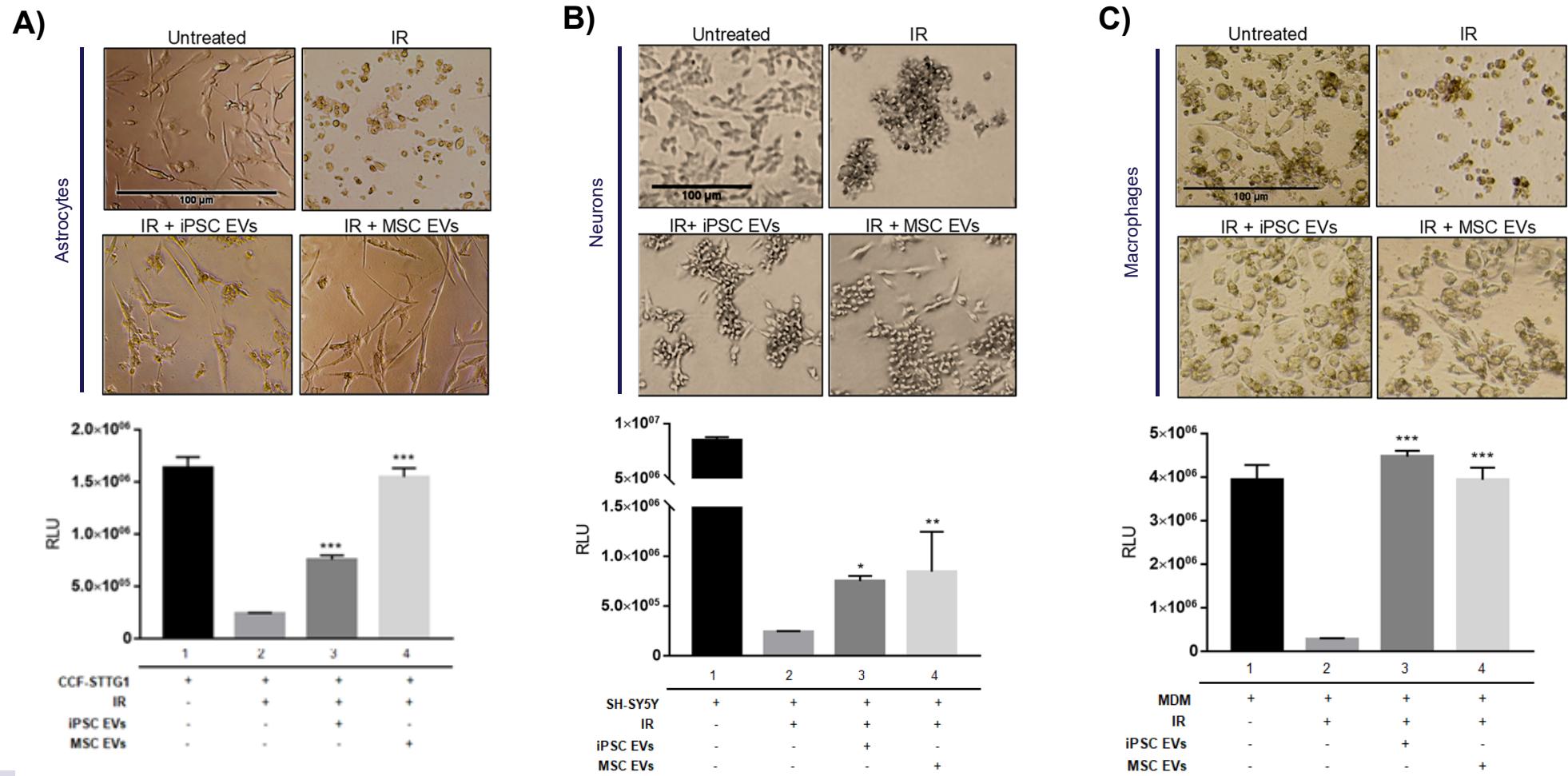
A)



B)

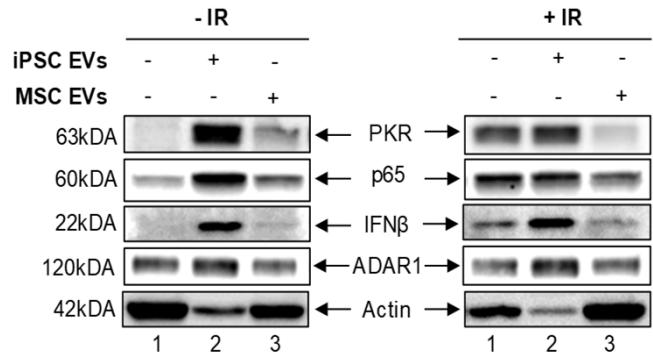


# EV Functionality

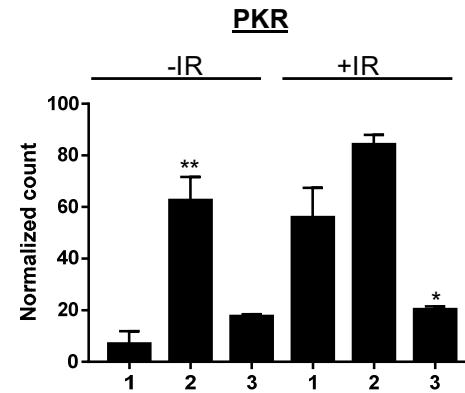


# EV Functionality

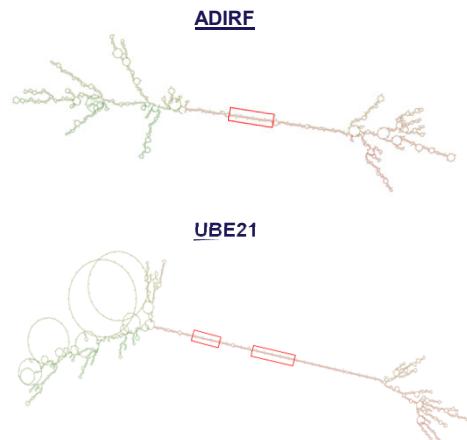
**A)**



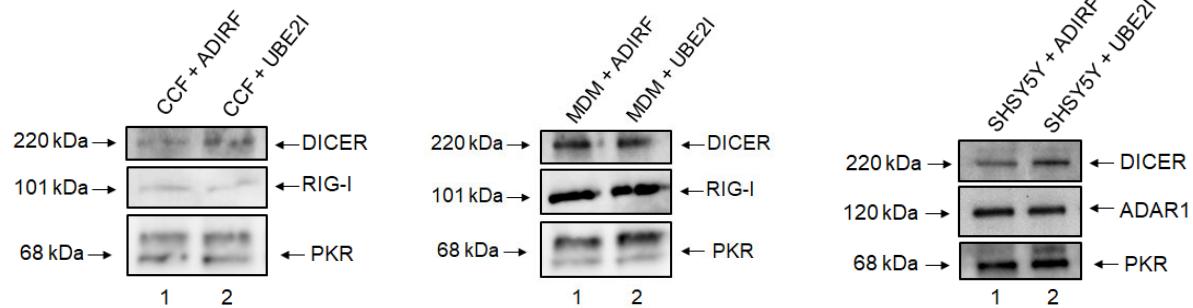
**B)**



**C)**



**D)**



# Overview

---

- Overview of extracellular Vesicles (EVs)
- Characterization and Analysis of “Damaging EVs”
- Characterization and Analysis of “Reparative EVs”
- Summary and Future Directions

# Summary

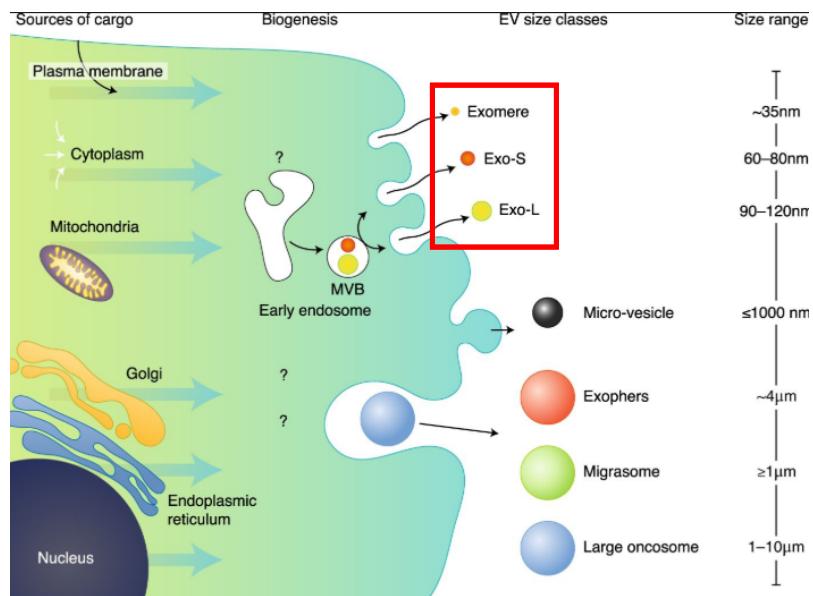
---

- Extracellular vesicles (EVs) such as exosomes are critical mediators of intercellular communication. The diverse biological cargo that is associated with these vesicles is believed to mediate the pleiotropic effects of EVs.
- Infected cells secrete “Damaging EVs” which contain viral non-coding RNAs and other viral proteins and these EVs can be separated from infectious virions.
- Stem cells secrete “Reparative EVs” which contain both non-coding RNAs and cytokines that can contribute to processes relevant to cellular repair.

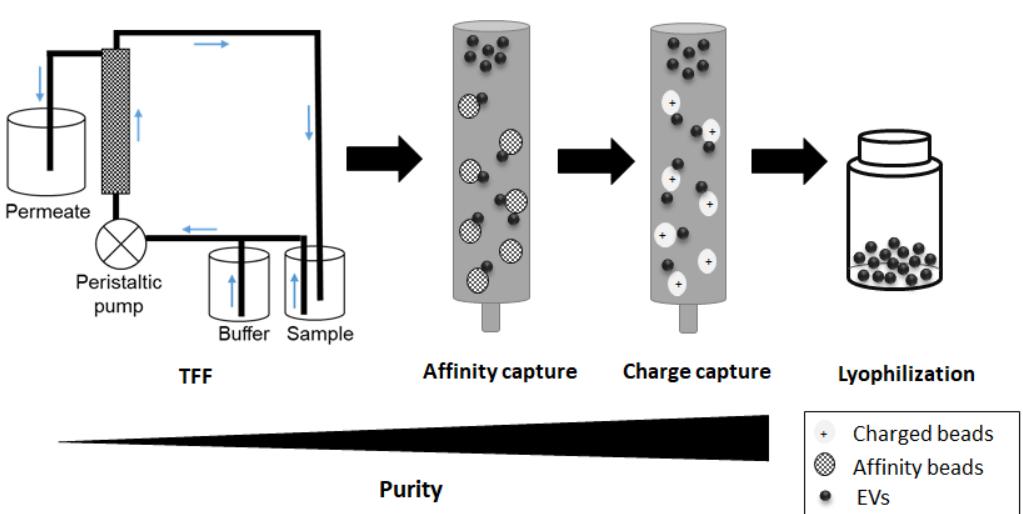
# Future Directions

## Further Purification of EV subpopulations

A)

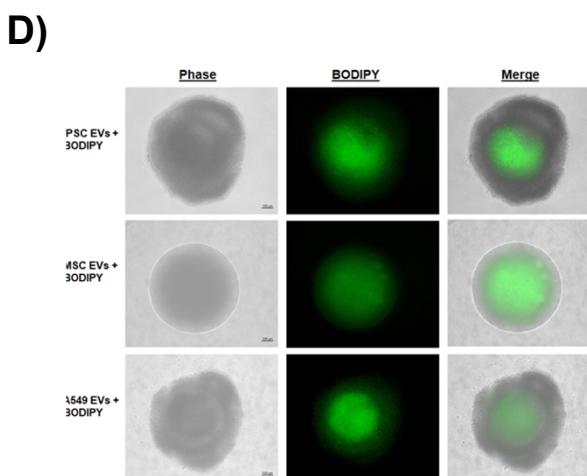
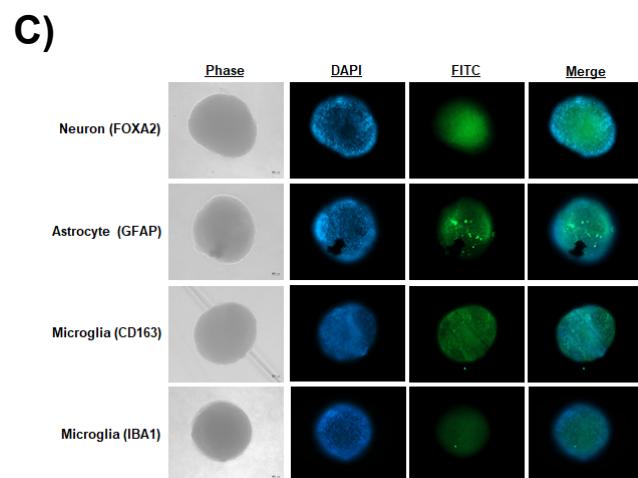
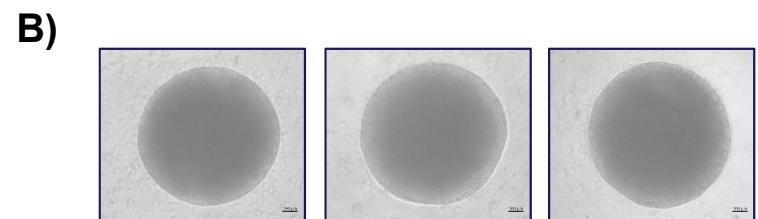
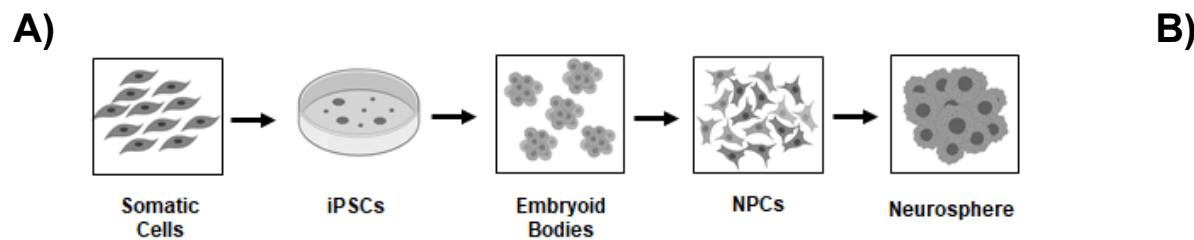


B)

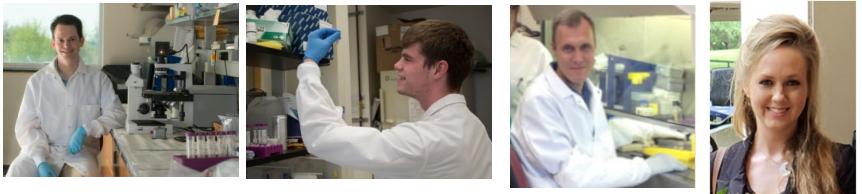
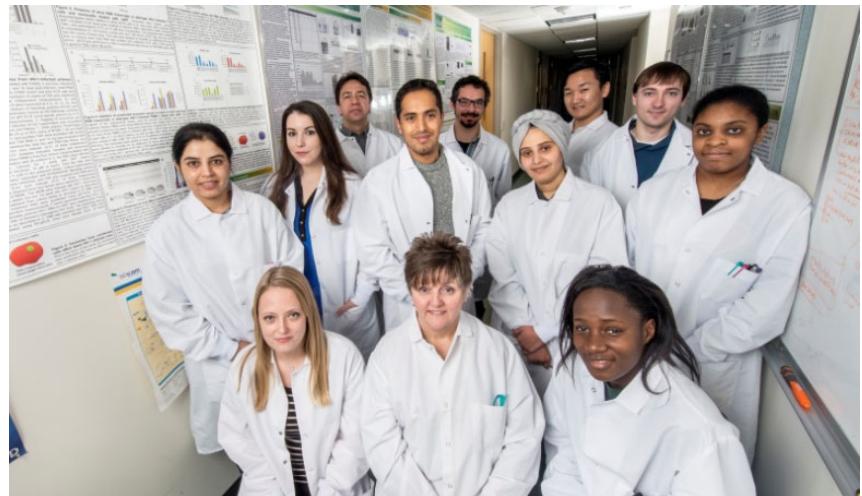


# Future Directions

Advanced functionality assays using neurospheres derived from ATCC ACS-5003™  
(unpublished data)



# Acknowledgements



## ATCC

- Dr. Siddhartha Paul
- Dr. Dezhong Yin
- Dr. Mindy Goldsborough
- Dr. James Kramer
- Dr. Katalin Kiss

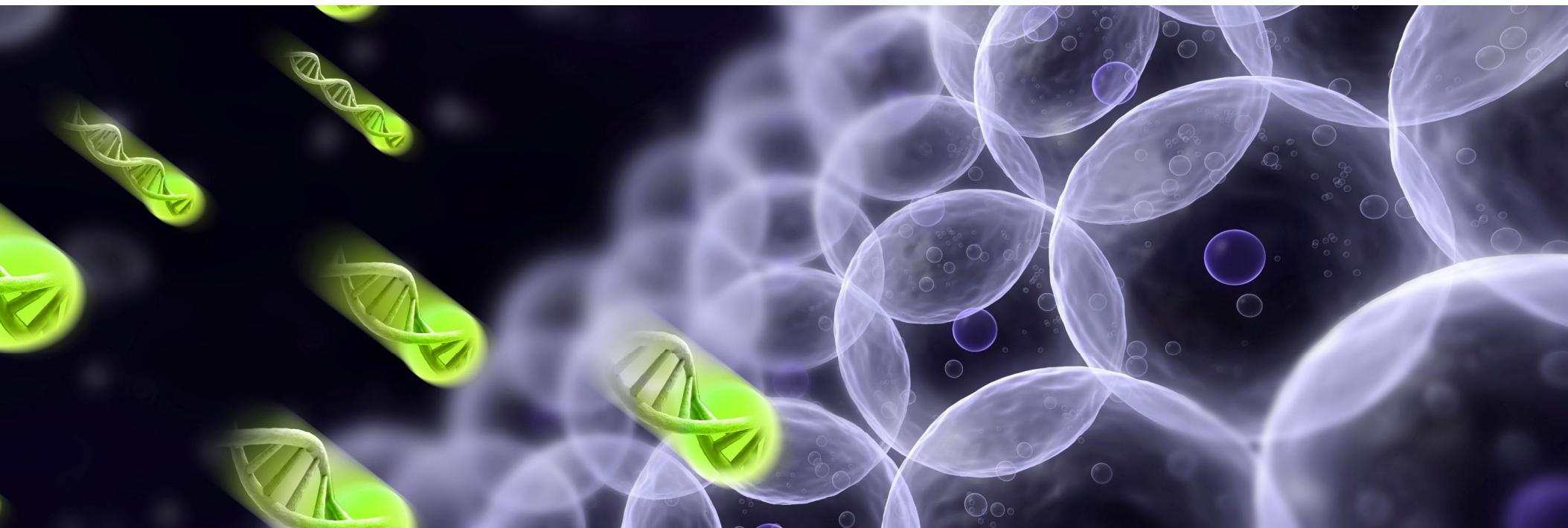
## GMU

- Dr. Fatah Kashanchi
- Dr. Lance Liotta
- Dr. Weidong Zhou
- Dr. Robert Barclay
- Dr. Catherine DeMarino
- Dr. Michelle Pleet
- Daniel Pinto
- Yuriy Kim
- Pooja Khatkar
- Gwen Cox

## Other

- Kamel Khalili (Temple)
- Rafal Kaminski (Temple)
- Archana Gupta (SBI)
- Ben Lepene (Ceres Nanosciences)
- Robbie Barbero (Ceres Nanosciences)
- L.Shultz (Jackson Lab)
- D. Price (Univ. of Iowa)
- F. Romerio (University of Maryland)
- O. Kutsch (UAB)
- V. Planelles (Utah)
- Lena Al-Harthi (Rush)
- Kevin Morris (City of Hope)





THANK YOU

# Cultivating collaboration to support global health

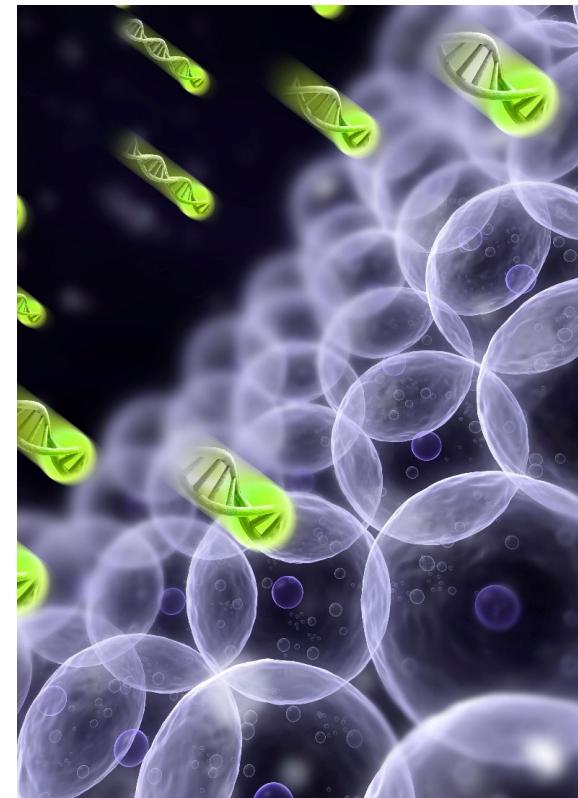
Learn more:

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

## Upcoming Webinar:

Antimicrobial Resistance: Arm Your Lab in the Fight Against Superbugs

- Presented by Christine Fedorchuk, Ph.D.
- February 27, 12:00 ET



© 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. Nanotrap is a trademark of Ceres Biosciences, Inc. OptiPrep is a trademark of Sigma Aldrich. qEV is a trademark of Izon Science, LTD. ExoMax is a trademark of System Biosciences.

