### TRANSFECTION REAGENTS; POWERFUL TOOLS TO ENABLE GENETIC MANIPULATION

James Clinton, Ph.D. Scientist, ATCC April 02, 2015







THE ESSENTIALS OF LIFE SCIENCE RESEARCH GLOBALLY DELIVERED\*

### 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 products and innovative solutions
- Strong team of 400+ employees; over one third with advanced degrees
- Broad range of biomaterials
  - Cell lines, iPSCs, primary cells, and hTERT immortalized cells
  - Bacteria, yeasts, protists, and viruses
  - Tumor cell panels
  - Media, sera, and reagents



Established partner to global researchers and scientists



2

# Introduction to Transfection

## **Best Practices**

# ATCC Transfection Reagents



### Introduction to transfection

- Method for introducing exogenous nucleic acid sequences into mammalian cells
- Widely used technique that has made expressing DNA or RNA in most types of cells relatively easy
- A variety of approaches have been developed for use across a range of applications
- No single approach will work for all conditions/cell types/applications





### **Applications**

- Gene function
- RNAi gene silencing
- Pathway analysis
- Functional screening
- Virus production
- Protein production
- Generation of stable cell lines
- Stem cell reprogramming
- Cell differentiation
- Genome editing with CRISPR/Cas9





### **Transfection methods**

**ATCC**°

Lipid	<ul> <li>Easy, most common method</li> <li>Variable efficiencies</li> <li>Will not work with all cell types</li> </ul>		
Viral	<ul> <li>Will transfect non-dividing cells</li> <li>Technically challenging, expensive</li> <li>Safety issues, immune response, mutagenesis</li> </ul>		
Electroporation	<ul> <li>Requires specialized equipment</li> <li>Cells must be in suspension</li> <li>Toxicity can be an issue</li> </ul>		
Physical	<ul> <li>Technically challenging, expensive</li> <li>Requires specialized equipment</li> <li>Works with non-nucleic acids; single cell transfection</li> </ul>		
Other	<ul> <li>Not common, may be technically challenging</li> <li>Non-lipid based chemicals</li> <li>Nanoparticles/ Laser/ Ultrasound/ Magnetic</li> </ul>		

### **Mechanism of lipid based transfection**





### **Typical transfection workflow**





Transfection

reagent

Complexes

### **Overexpression vs. knockdown**



Introduce foreign plasmid DNA/mRNA to induce expression of a desired transcript/protein



Utilize RNAi pathway to degrade or inhibit translation of mRNA transcripts and subsequently reduce the amount of protein



### **Transient vs. stable transfection**

#### Transient

- Foreign gene not integrated into genome
- Expression persists for limited time
- Foreign gene lost due to cell division, degradation, or other factors

#### Stable

- Initially a transient transfection
- Use co-expressed selection markers
- Long-term, only cells that have integrated the foreign gene persist



# Introduction to Transfection

## **Best Practices**

# ATCC Transfection Reagents



### **Transfection: best practices**





### **Cell culture conditions**





### Example: culture conditions can be critical

#### Primary Uterine Smooth Muscle Cells (SMCs; ATCC<sup>®</sup> No. PCS-460-011)

Phase Contrast

Transfected in complete growth media

10X

Transfected in differentiation media



ATCC



Transfection is completely inhibited

### **Nucleic acids**





### **Experimental design & execution**

Transfection Protocol	<ul> <li>Use master mixes</li> <li>Distribute complexes evenly</li> <li>Store DNA/RNA properly</li> </ul>			
Proper Controls	<ul> <li>Positive and negative control</li> <li>Transfected and un-transfected controls</li> </ul>			
Monitor Toxicity/ Off-target Effects	<ul> <li>Morphological changes</li> <li>Presence of vacuoles</li> <li>Changes in proliferation</li> </ul>			
Validate Results	<ul> <li>Multiple assays</li> <li>For siRNA: test multiple sequences</li> <li>For miRNA: increase &amp; suppress</li> </ul>			



### **Assay methods**





### Assay timing



**Time post-transfection** 



### **Transfection reagents considerations**

### **Ideal reagent**

- Effective in all cell types
- No optimization necessary
- No cytotoxicity

### Reality

- Effective in your cell type of interest
- Broad activity across culture conditions and protocols
- Minimal cytotoxicity
- An optimized protocol delivers desired expression



### **Transfection reagents**





### **Best practices summary**





# Introduction to Transfection

## **Best Practices**

ATCC Transfection Reagents



### **ATCC** transfection reagents overview

- Used for transfection of mammalian adherent and suspension cells
- Formulated for low cytotoxicity and high efficiency
- Produces high levels of gene expression (or inhibition)
- Suitable for both transient and stable transfection
- Sterility, purity, and performance tested
- Animal component-free

Reagent	ATCC <sup>®</sup> No.	Volume	Storage
GeneXPlus Transfection Reagent	ACS-4004	1 mL	-20°C
TransfeX <sup>™</sup> Transfection Reagent	ACS-4005	500 µL	4°C
siFEX™ RNAi Transfection Reagent	ACS-4006	500 µL	4°C

For more information on our transfection reagents: www.atcc.org/transfection



### **ATCC** transfection reagents selection guide

Reagent	Plasmid DNA	mRNA	siRNA & miRNA	Suspension	Hard to Transfect Cells			
GeneX <i>Plus</i>	$\checkmark \checkmark \checkmark$			$\checkmark$				
TransfeX™	<b>√ √ √</b>	√ √		~	<b>√ √ √</b>			
Newly released in 2015:								
siFEX™			<b>√ √ √</b>	<b>~</b>				



### Available GeneXPlus optimized protocols



Optimized protocol list as of April 2015

- Find the current list of available protocols at <u>www.atcc.org/transfection</u>
- Contact Technical Service at tech@atcc.org





#### SH-SY5Y (ATCC<sup>®</sup> No. CRL-2266<sup>™</sup>)





#### HEK 293T SF (ATCC<sup>®</sup> No. ACS-4500<sup>™</sup>)

THP-1 (ATCC<sup>®</sup> No. TIB-202<sup>™</sup>) **Phase Contrast** 

Transfection of suspension cells with GeneXPlus

GFP

### Available TransfeX<sup>™</sup> optimized protocols

#### Continuous

- A549
- HeLa
- LNCap
- MDA-MB-231
- HepG2
- Caco-2
- C2C12
- 3T3-L1
- CHO-K1
- HEK293
- HUV-EC-C
- MCF7
- NuLi-1
- TIME
- BT-142



- Bone marrowderived MSCs
- Adipose tissuederived MSCs
- Cord blood-derived MSCs
- Induced pluripotent stem cells (iPSCs)

#### **Primary**

- Dermal fibroblasts (DFs)
- DMEC
- HEMCs
- HUVECs
- RPTECs
- Large airway epithelial cells
- Large airway SMCs
- Uterine fibroblasts
- Uterine SMCs
- Prostate epithelial cells

Optimized protocol list as of April 2015

- Find the current list of available protocols at <u>www.atcc.org/transfection</u>
- Contact Technical Service at tech@atcc.org

27



# **ATCC**<sup>®</sup>

#### MCF-7 (ATCC<sup>®</sup> No. HTB-22<sup>™</sup>)



#### NuLi-1 (ATCC<sup>®</sup> No. CRL-4011<sup>™</sup>)

### A549 (ATCC<sup>®</sup> No. CCL-185<sup>™</sup>)

**Transfect continuous cells with TransfeX™** 



### Transfect primary cells with TransfeX<sup>™</sup>

**Phase Contrast** 

Primary DFs (ATCC<sup>®</sup> No. PCS-100-012)

#### Primary HUVECs (ATCC<sup>®</sup> No. PCS-100-010)

**Primary Uterine SMCs** 

(ATCC<sup>®</sup> No. PCS-460-011)

10X

GFP

ATCC<sup>°</sup>

### **Transfect stem cells with TransfeX™**

iPSCs (ATCC<sup>®</sup> No. ACS-1012<sup>™</sup>)



# Bone Marrow-derived MSCs (ATCC<sup>®</sup> No. PCS-500-012)







### **Transfection of mRNA with TransfeX™**

#### HeLa (ATCC<sup>®</sup> No. CCL-2<sup>™</sup>)





#### Primary DF



A549



### Available siFEX optimized protocols

#### Continuous

- TeloHAEC-GFP
- HepG2
- MCF-7
- A549
- MDA-MB-231
- HeLa
- MRC-5
- HEK293T/17
- HUV-EC-C
- 3T3-L1
- C2C12
- Caco2
- LNCap

#### Stem

- Adipose tissue derived MSCs
- Cord blood-derived MSCs

#### **Primary**

- Primary DFs
- Primary prostate epithelial cells

Optimized protocol list as of April 2015

- Find the current list of available protocols at <u>www.atcc.org/transfection</u>
- Contact Technical Service at tech@atcc.org



### **Transfection of fluorescently labeled siRNA**



### **Knockdown of constitutive GFP expression**

#### TeloHAEC-GFP (ATCC<sup>®</sup> No. CRL-4054<sup>™</sup>)

#### Untransfected



+Anti-GFP siRNA

For more information on TeloHAEC and TeloHAEC-GFP cells, visit the ATCC<sup>®</sup> Research page in our Learning Center at <u>www.atcc.org/LearningCenter</u>



### siFEX: performance tested



### **Transfection of pre-miRNA with siFEX**



Competing Product #2 Cell viability expressed in blue

- Control SIFEXTM
- HeLa cells in 24-well plate.
- Transfected with 20 nM hsa-miR-1 pre-miRNA
- Expression of PTK9 mRNA assessed 48 h post-transfection via RT-qPCR
- Results were calculated via ΔΔCT method, n=6 transfections, mean ± STD



### Summary

- Lipid-mediated transfection is a powerful tool and useful in a variety of applications
- ATCC offers a variety of transfection reagents suitable for the transfection of plasmid DNA, mRNA, and dsRNA into a variety of cell types
- ATCC transfection reagents have been performance tested to deliver high efficiency and low cytotoxicity
- ATCC offers optimized transfection protocols for dozens of cell types to help you achieve results faster
  - Continuous cell lines
  - Primary cells
  - Stem cells
  - Adding new protocols all the time



### Thank you!

Register for more webinars in the ATCC "*Excellence in Research*" webinar series at <u>www.atcc.org/webinars</u>.



May 21, 2015 10:00 AM ET or 3:00 PM ET Jodie Lee, M.S., *Lead Biologist*, ATCC Seeing is Believing – Reporter Labeled Microbial Control Strains



The ATCC<sup>®</sup> *"Excellence in Research"* webinar series returns in Fall 2015. Look for webinars starting in August at www.atcc.org/webinars.



Thank you for joining today! Please send additional questions to <u>tech@atcc.org</u>