

# Drug Transporter Studies: Lysosomal Trapping

Andrew G. Taylor, Ph.D.

Manager, Technical Support for Services

XenoTech

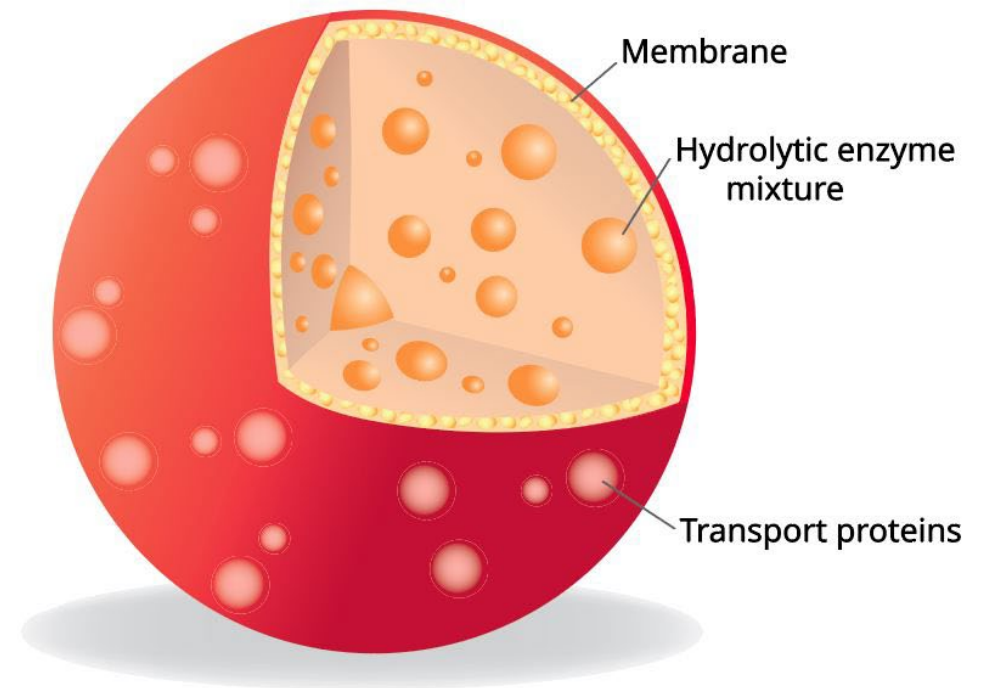




## What are Lysosomes?

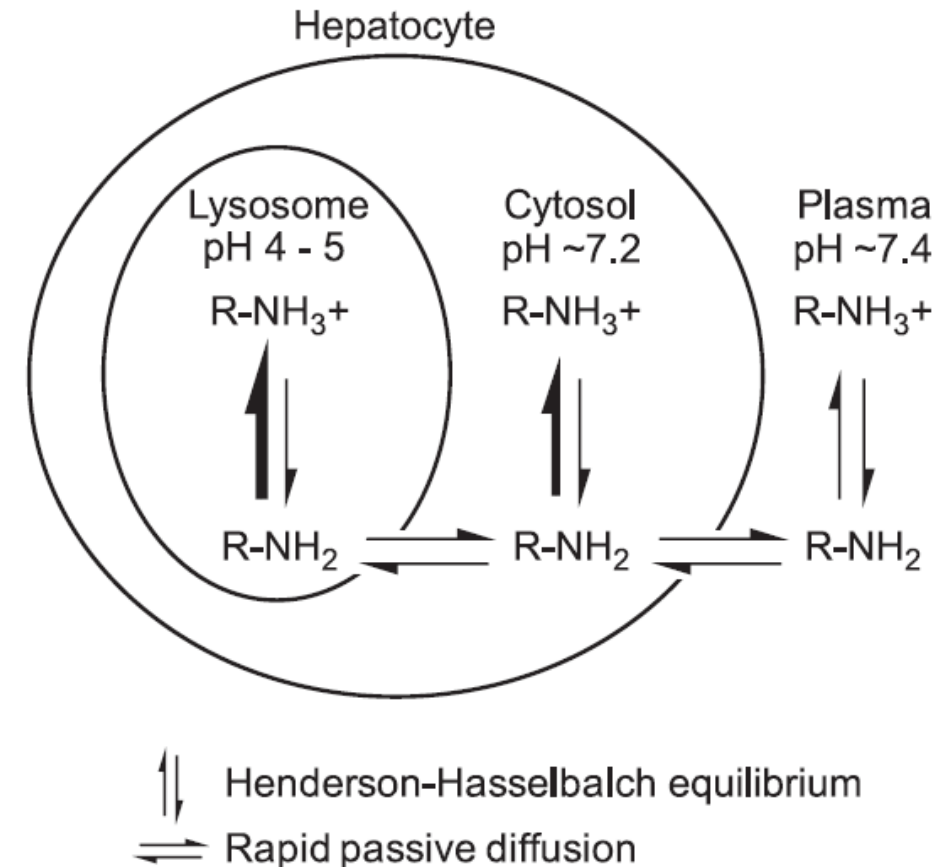
- Resident macrophage organelle in eukaryotic cells
- Acidic environment (pH ~4-5)
- Hydrolytic enzymes
- Break down larger molecules (e.g., proteins, polysaccharides, lipophilic compounds)
- ~1% of hepatic cell volume

## LYSOSOME



# Mechanism of Lysosomal Trapping

- Lysosomotropism (lysosomal trapping) is a physicochemical process
- Cationic amphiphilic compounds (many CNS and cardiovascular drugs)
- $\log P > 1$ ,  $pK_a > 6.5$
- Diffuse easily across membranes
- Become protonated (+ charge) in lysosome and become trapped



From Kazmi et al., 2013, DMD



## Lysosomal Trapping and DDI

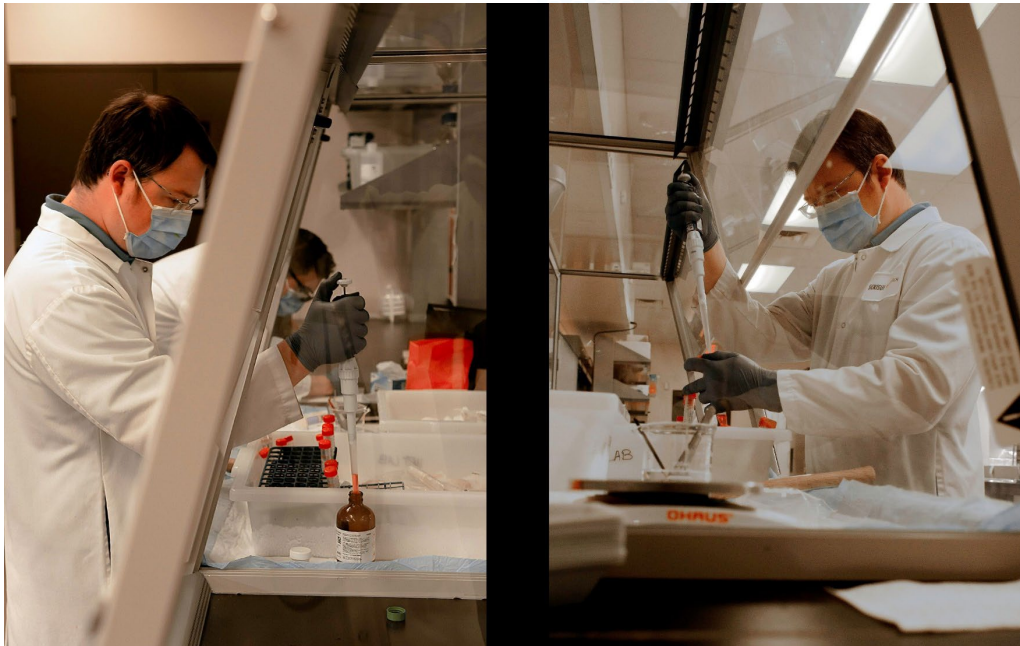


- Can lead to high organ-to-blood ratios, often mistaken for active drug transport
- Competition for lysosomal trapping, concomitant administration of lysosomotropic agents could lead to elevated drug exposure levels
- Accumulation of lipophilic amines can lead to drug-induced phospholipidosis due to decreased phospholipid catabolism



# Lysosomal Trapping Study Designs

Studies are conducted in Fa2N-4 cells (immortalized hepatocytes) that don't have a lot of drug metabolizing enzymes

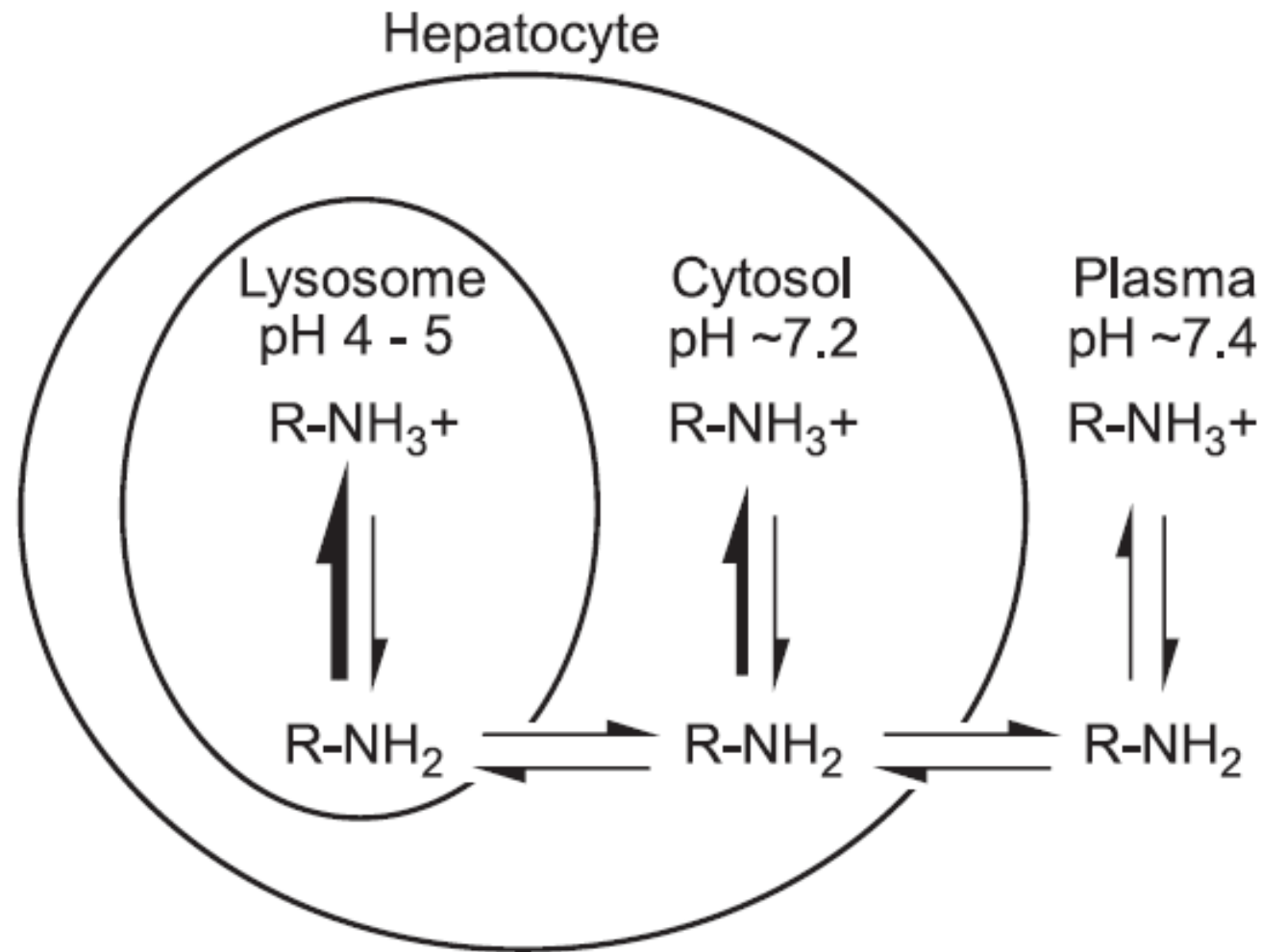


- 1) LysoTracker Inhibition Assay (screen)
  - LysoTracker Red used as a fluorescent probe (measured by fluorescence)
  - Six test article concentrations
- 2) Mechanistic Determination (definitive)
  - TA measured by LC/MS
  - 2 [TA], 2 time points, w/ and w/o NH<sub>4</sub>Cl
  - Positive control propranolol w/ and w/o NH<sub>4</sub>Cl
  - If worries of ammonium chloride interacting with TA, chloroquine can be used instead

A BioIVT Company

# Mechanistic Assay for Lysosomal Trapping

Without  $\text{NH}_4\text{Cl}$

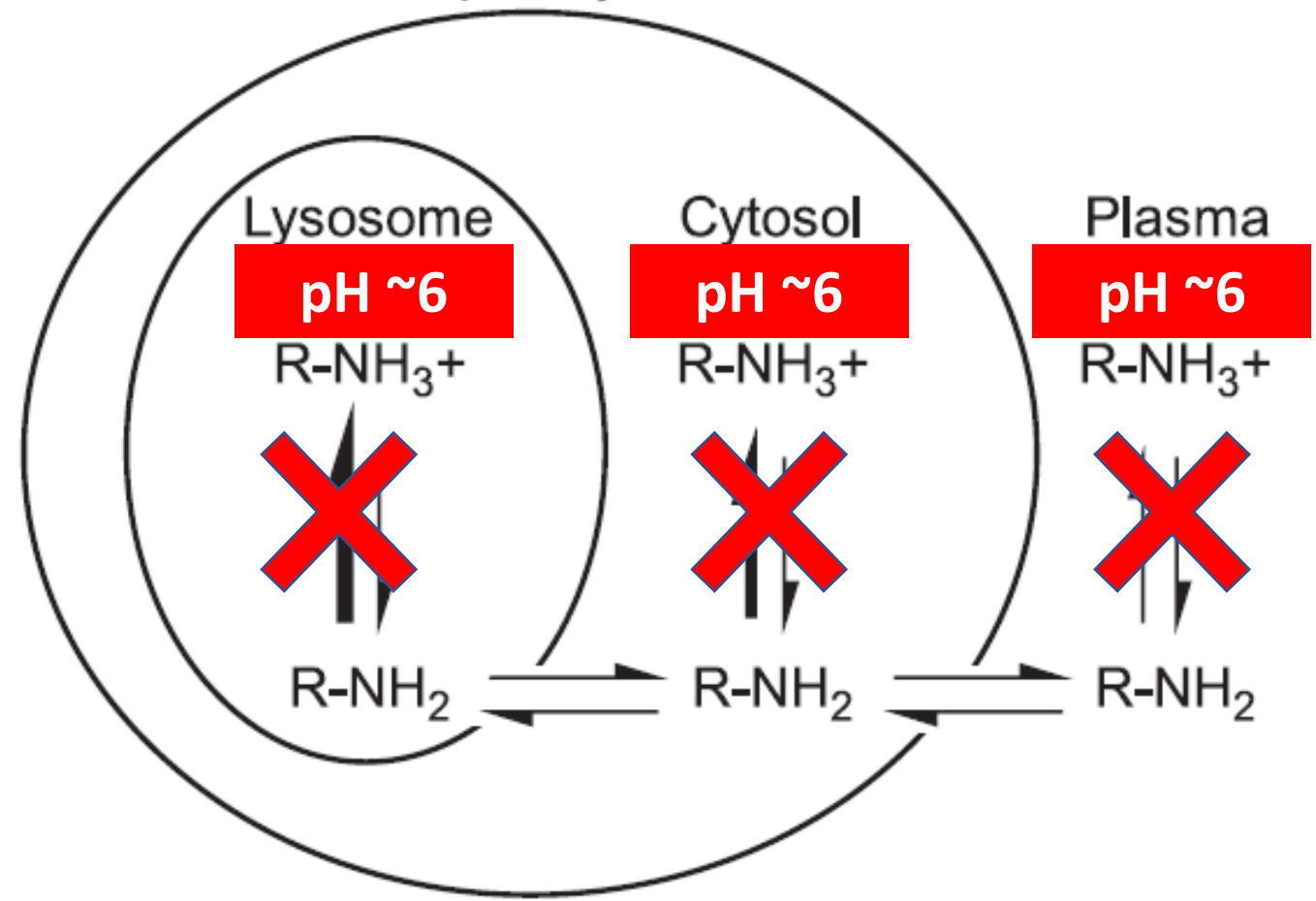


A BioIVT Company

# Mechanistic Assay for Lysosomal Trapping

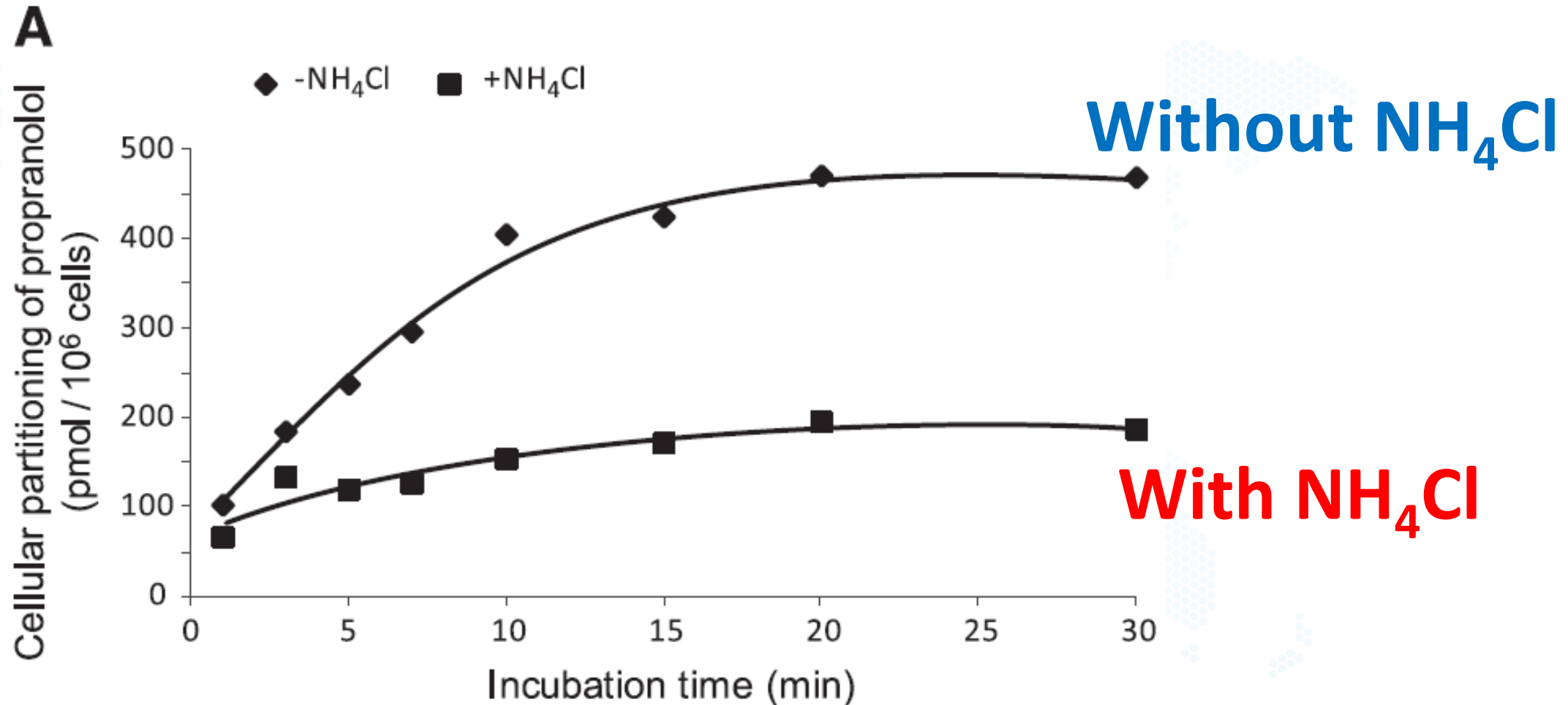
Hepatocyte

With  $\text{NH}_4\text{Cl}$



A BioIVT Company

# Mechanistic Assay for Lysosomal Trapping



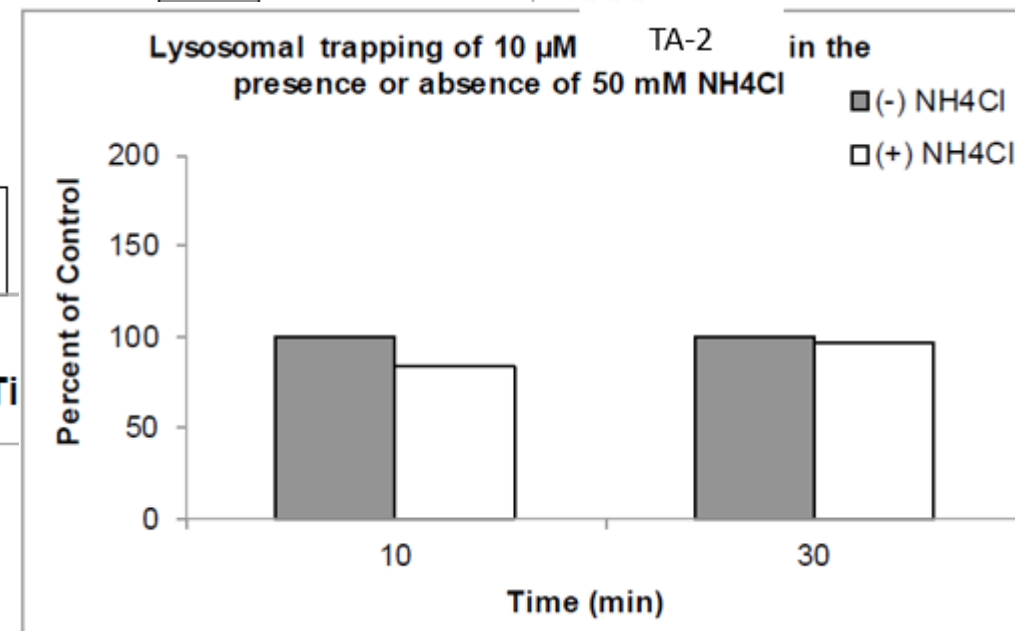
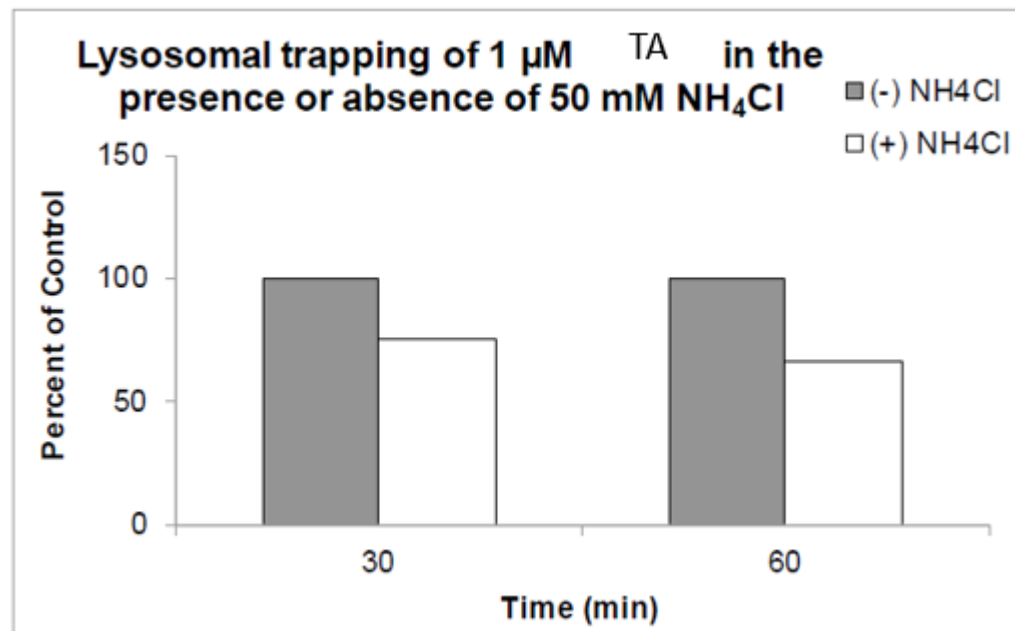
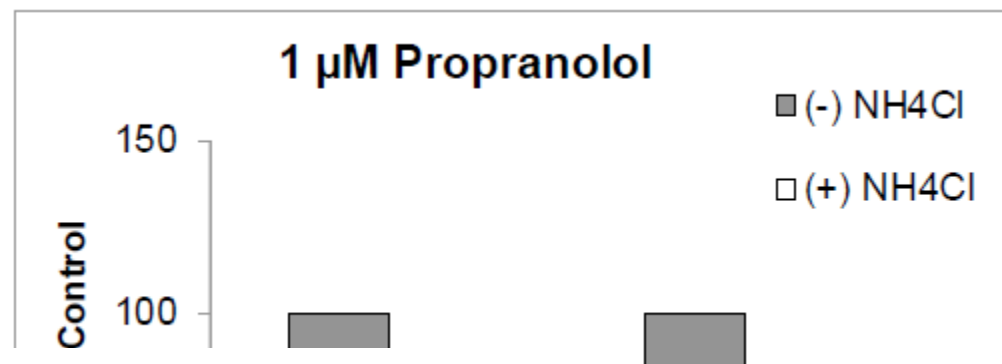
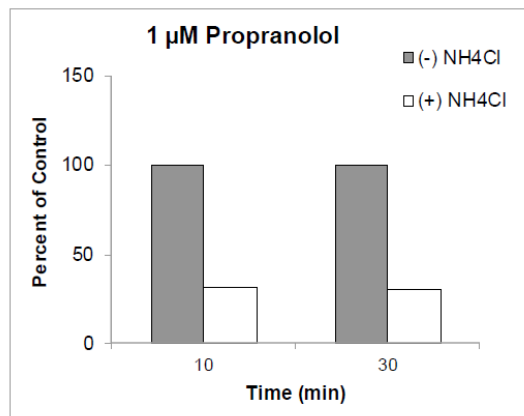
From Kazmi et al., 2013, DMD





A BioIVT Company

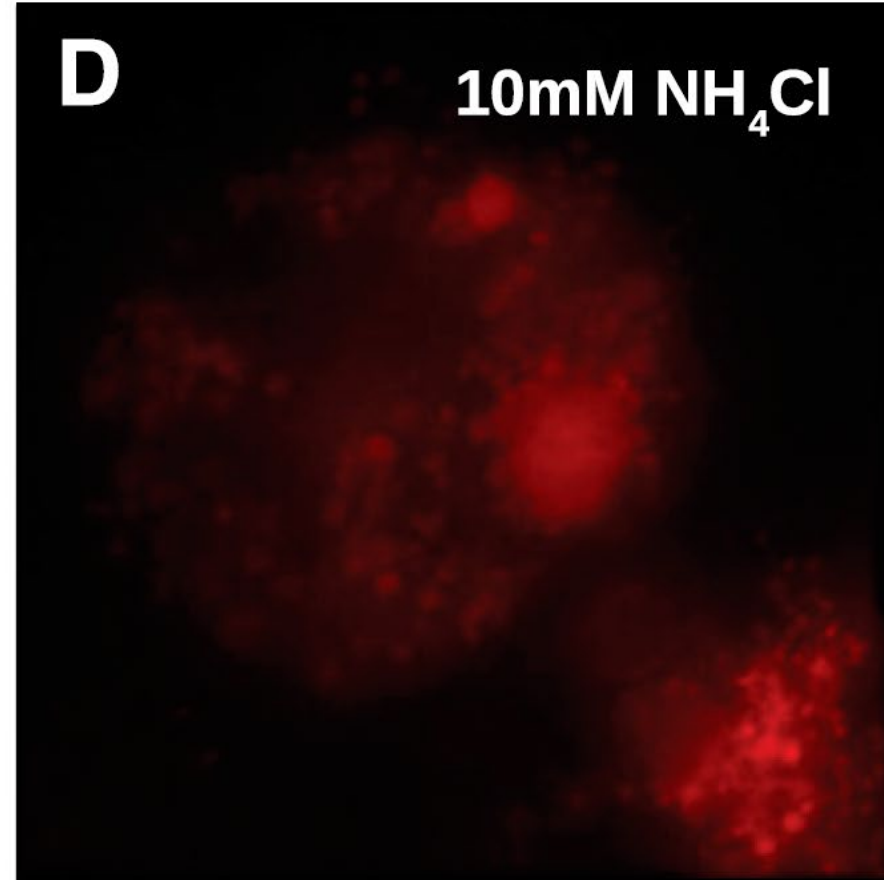
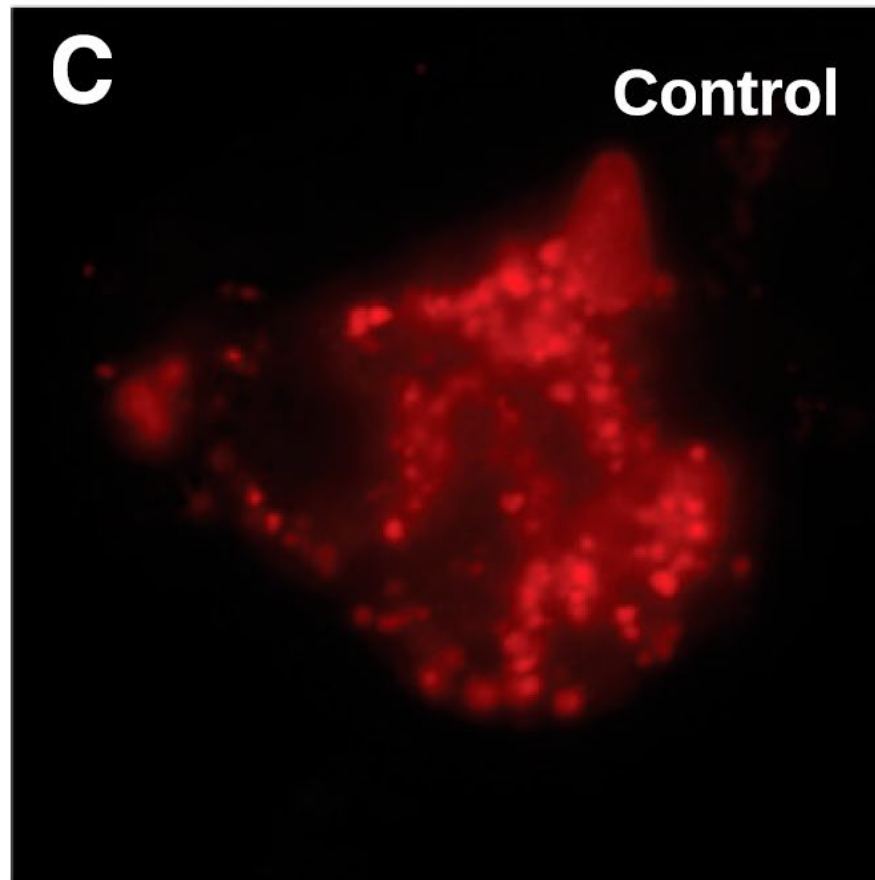
## Mechanistic Assay for Lysosomal Trapping



A BioIVT Company

# LysoTracker Assay for Lysosomal Trapping

Fa2N-4 Cells

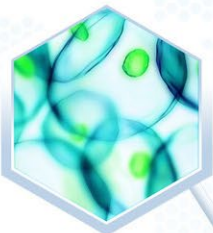


From Kazmi et al., 2013, DMD

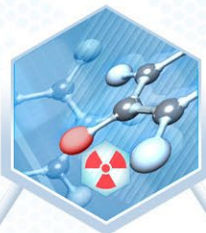
# XenoTech Products Available for Lysosomal Trapping

Test System	Details	Supporting Reagents
Immortalized human hepatocytes (Fa2N-4)	<ul style="list-style-type: none"> <li>Plated on 6, 12, 24, 48 or 96 well collagen-coated plate</li> <li>Consistent results, no lot-to-lot variability</li> </ul>	<ul style="list-style-type: none"> <li>Proprietary nutrient-rich media containing phenol red that is supplemented with serum (component B) prior to use</li> <li>Used for thawing, isolating, and seeding Fa2N-4 cells</li> <li>100 mL, 500 mL, or 1 L volumes</li> </ul>
Primary human hepatocytes	<ul style="list-style-type: none"> <li>Cryoplateable</li> <li>4 or 6 mil AMY</li> <li>Maintain enzyme and transporter activity</li> </ul>	<ul style="list-style-type: none"> <li>OptiTHAW</li> <li>OptiPLATE</li> <li>OptiCULTURE</li> </ul>

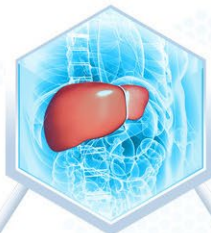
Cell & Tissue-Based Products



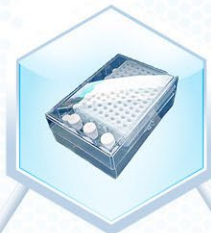
Radiolabeling



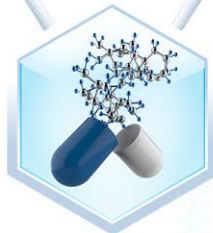
*in vitro* ADMET & Pharmacology



Metabolite ID & Production



Screening



API Manufacturing



*in vivo* ADMET & QWBA



Bioanalytical

**Thank you!**