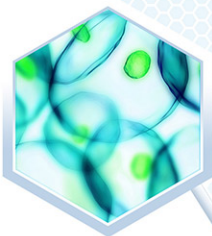
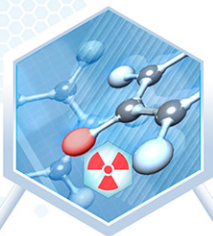


PROVEN GLOBAL CONTRACT RESEARCH EXPERTISE FROM DISCOVERY THROUGH CLINICAL SUPPORT

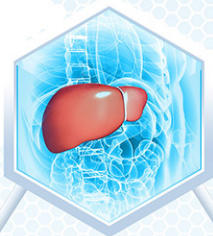
Cell & Tissue-Based Products



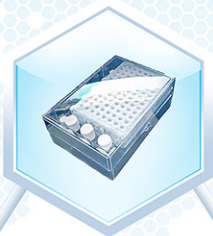
Radiolabeling



in vitro ADMET & Pharmacology



Metabolite ID & Production



Screening



API Manufacturing



in vivo ADMET & QWBA



Bioanalytical



EXPERTISE • EFFICIENCY • SUPPORT • PRECISION

Welcome to the webinar...

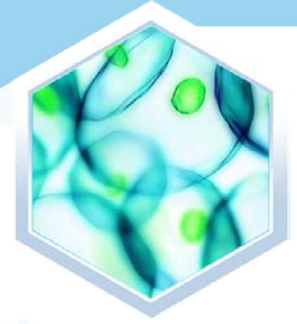
We will begin shortly

A BioIVT Company

PROVEN GLOBAL CONTRACT RESEARCH EXPERTISE
FROM DISCOVERY THROUGH CLINICAL SUPPORT

Investigation of freshly purified rat tritosomes and human hepatic lysosomes as an *in vitro* tool for characterization of biologic drugs

Chris Bohl**Research Scientist, Products R&D**
cbohl@xenotechllc.com



Lysosome Background

Discovered and named by Christian de Duve (Nobel Prize in 1974)

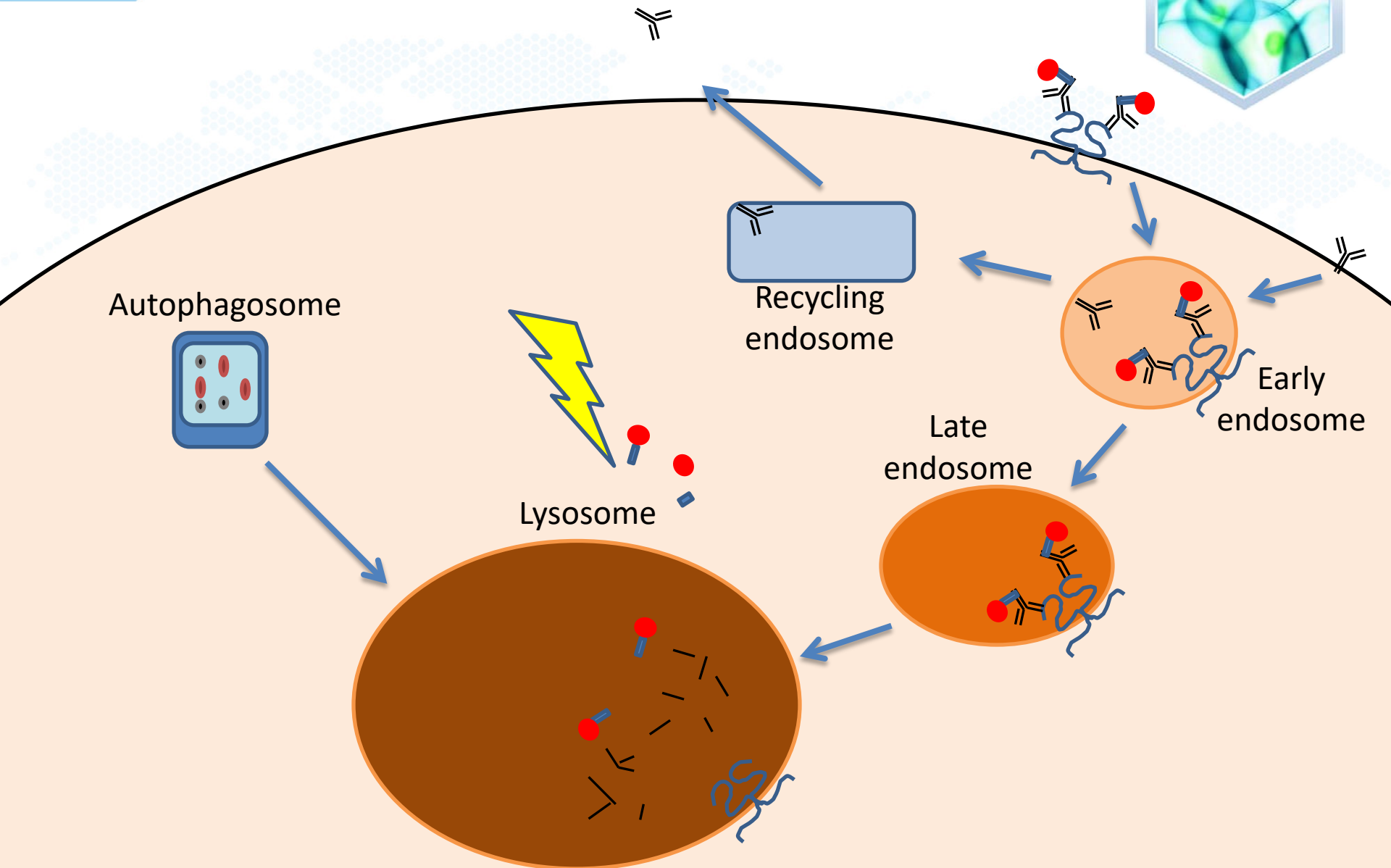
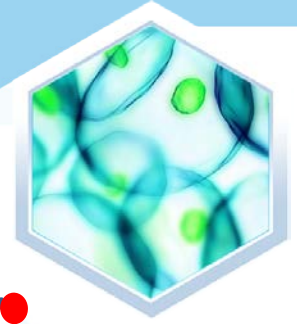
Lysosomes are a membrane bound, cellular organelle that is the sight of degradation/catabolism.

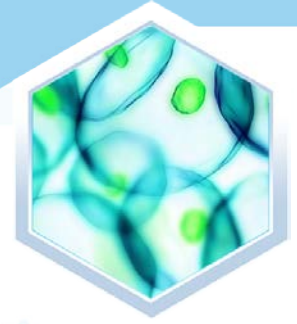
- extracellular substrates (endocytic pathway)
- intracellular substrates (autophagy pathway)

Contain a multitude of acidic hydrolytic enzymes and vary greatly in size.

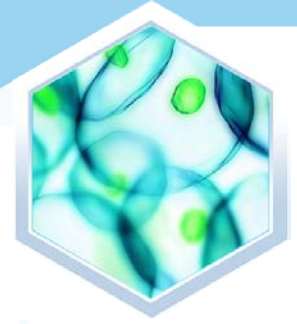
Lysosomes contain a variety of catabolic enzymes and are being designed as the first site of catabolism/activation of targeted biopharmaceuticals that enters cells through the endosomal-lysosomal pathway.

Isopycnic densities similar to mitochondria.





Rat Tritosomes



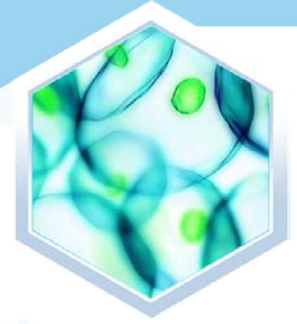
What are Tritosomes?

Tritosomes comes from the use of **Triton** WR 1339 (now called Tyloxapol) to modify **Lysosomes** density.

Method was first described in the 60's and was developed due to overlapping densities between lysosomes and mitochondria in sucrose (most common density gradient material at the time of tritosome development).

Tyloxapol is trafficked to the lysosomes and results in altered lysosomal lipid composition. Combined with the uptake and sequestration of the Tyloxapol, the Tritosomes have a lighter density than untreated lysosomes and allows improved separation from mitochondria using sucrose density gradients.

Common technique used in research to purify and enrich lysosomes.



What are Tritosomes?

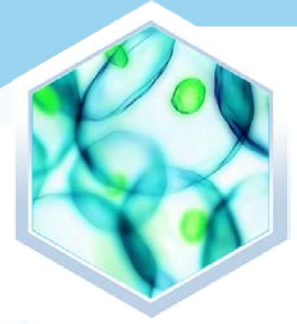
Highly purified lysosomes

High specific activity of lysosomal enzymes and activities

- Acid phosphatase
- Cathepsin B
- RNase

Simplified *in vitro* system

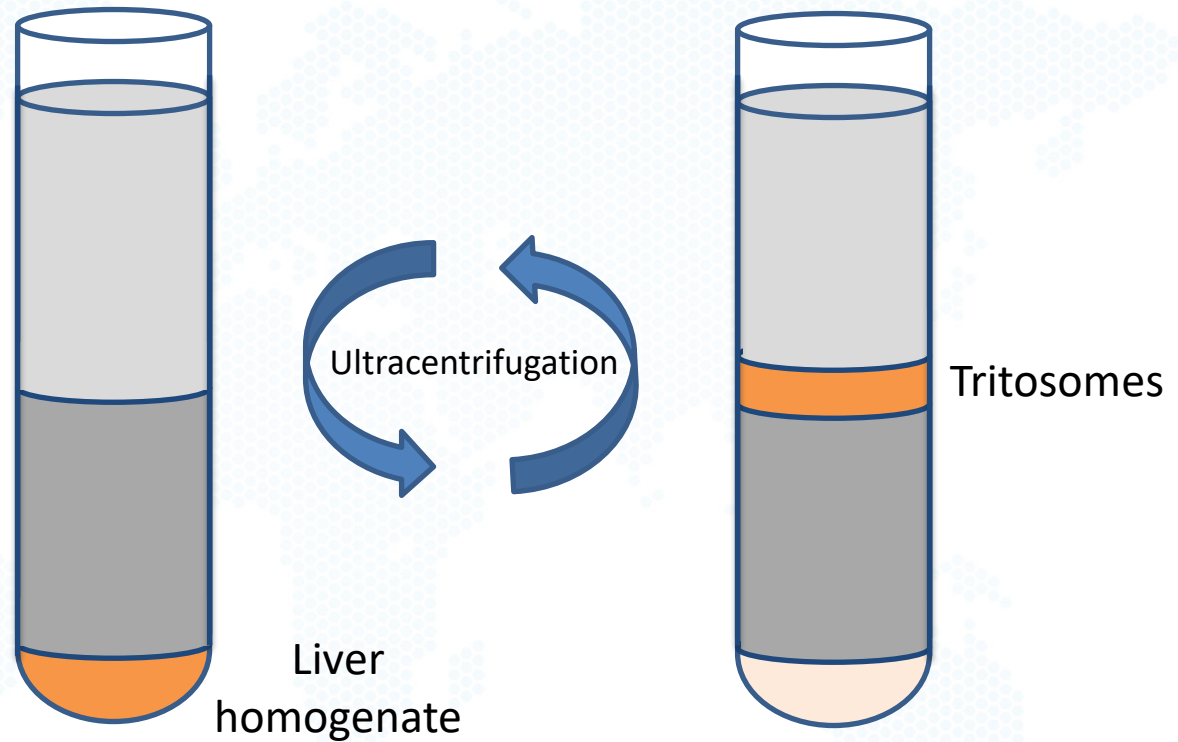
Tritosomes are disrupted and are ready- to – use
(EDTA and protease inhibitor free)

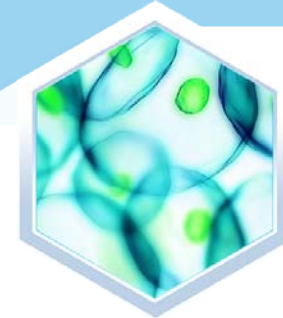


Tritosomes Production and Isolation



<http://www.criver.com>

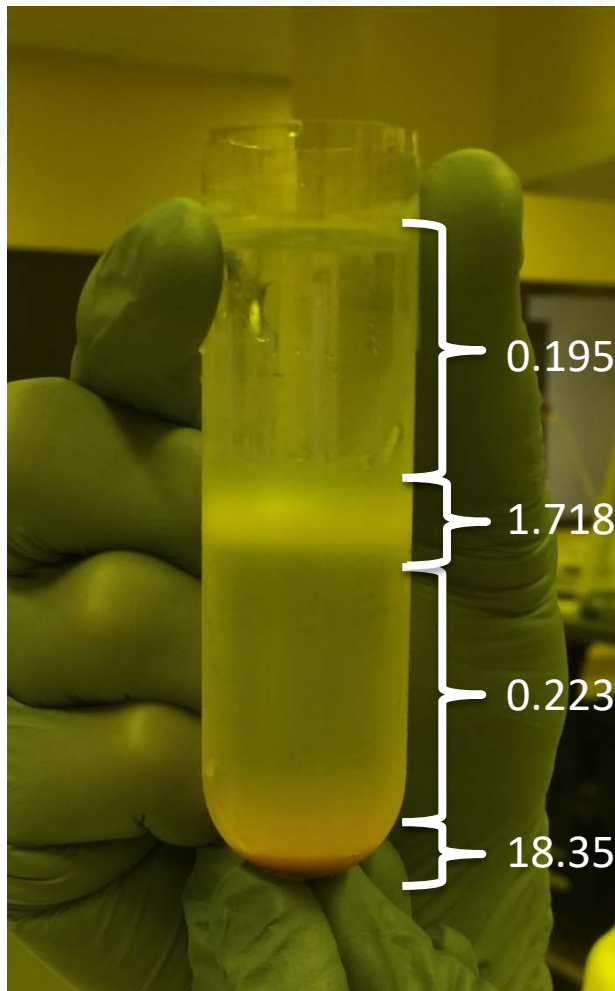




Tritosome Purity

Initial Liver Homogenate

34.05mg/ml Cat B = 1.43 U/mg COX = 0.0367 U/mg



0.195mg/ml Cat B = below detection level COX = 5.104U/mg

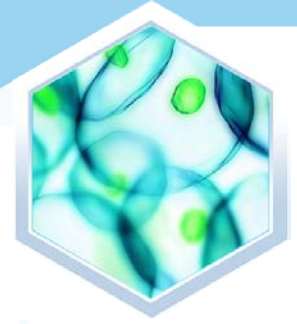
1.718mg/ml Cat B = 212.4 U/mg COX = 0.124 U/mg

0.223mg/ml Cat B = 98.9 U/mg COX = 38.906U/mg

18.35mg/ml Cat B = 5.82 U/mg COX = .0047U/mg

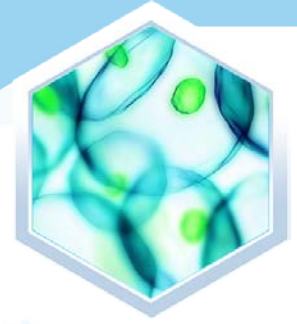
Cat B units = 1 μ M of AMC released / min

COX units = 1 μ M of Cytochrome C oxidized / min

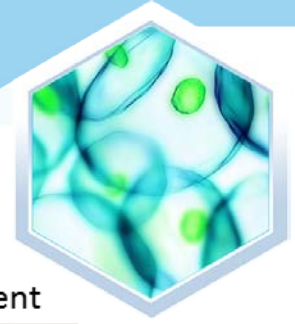


Conclusions

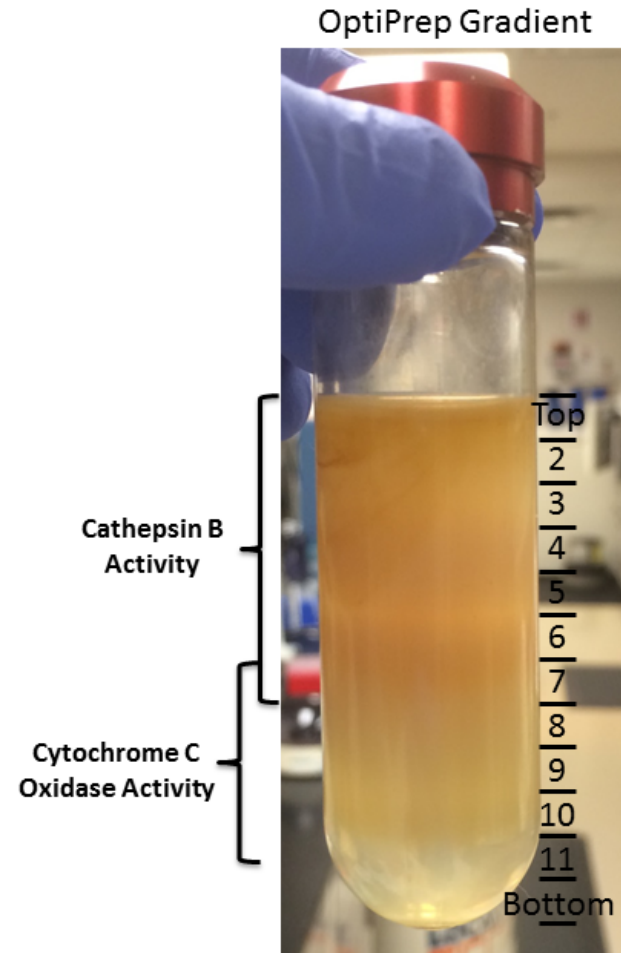
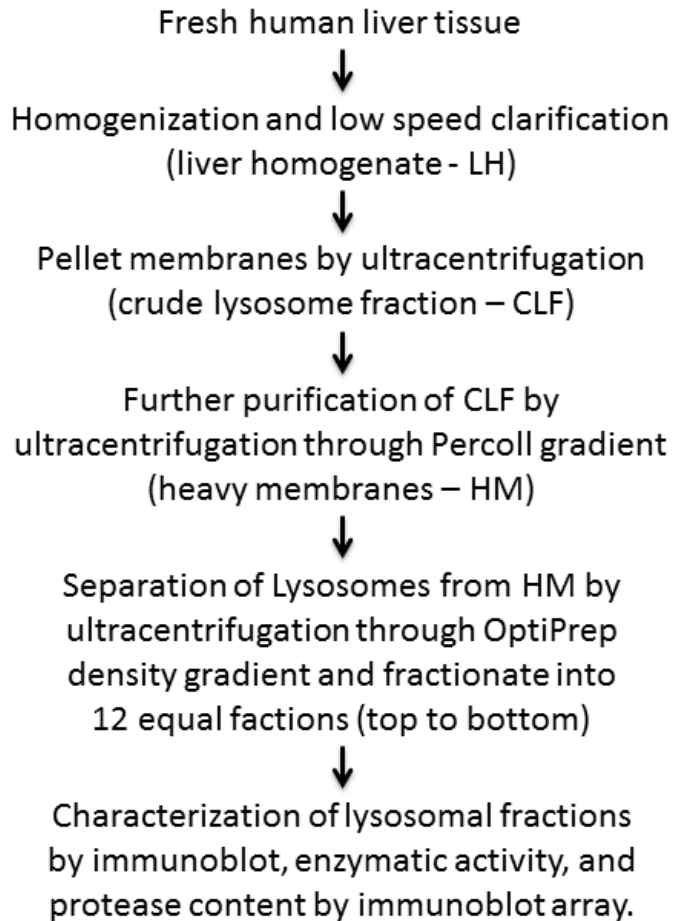
High specific activity for lysosomal enzymes (Acid phosphatase, cathepsin B, and nuclease) with minimal activity from enzymes associated with mitochondria.



Human Hepatic Lysosomes



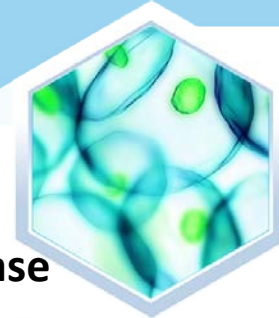
Lysosome Isolation



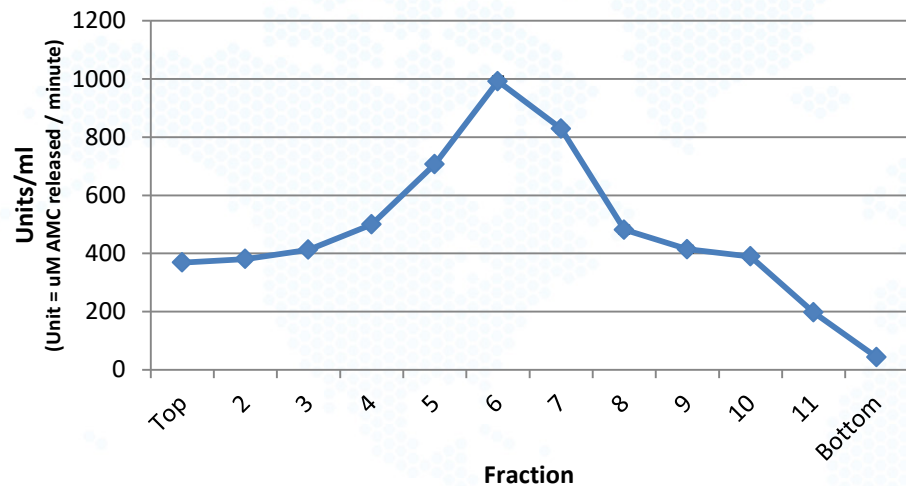


A BioIVT Company

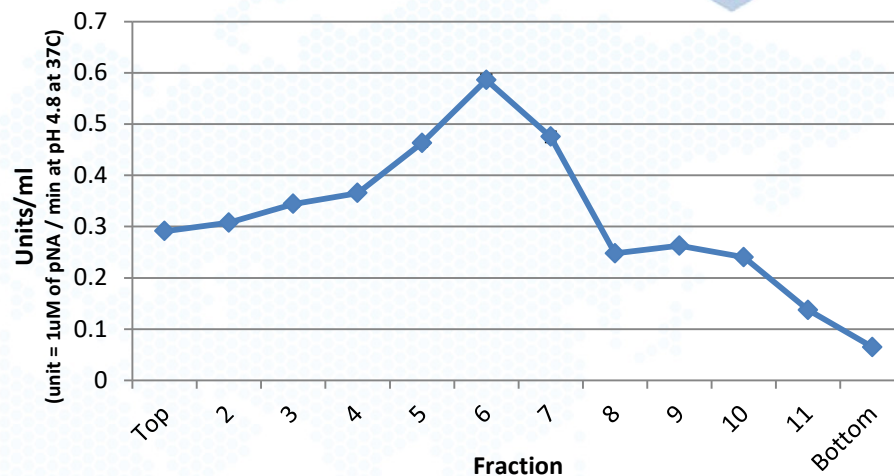
Identification of where the lysosome are in the gradient



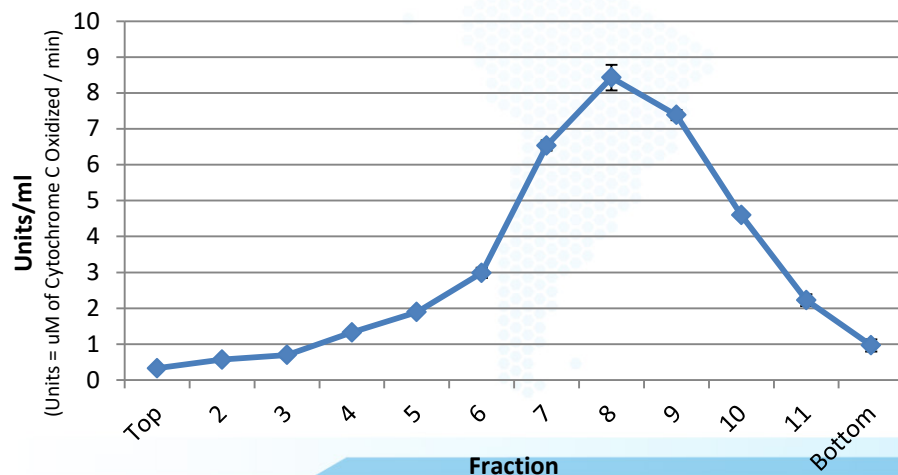
Cathepsin B



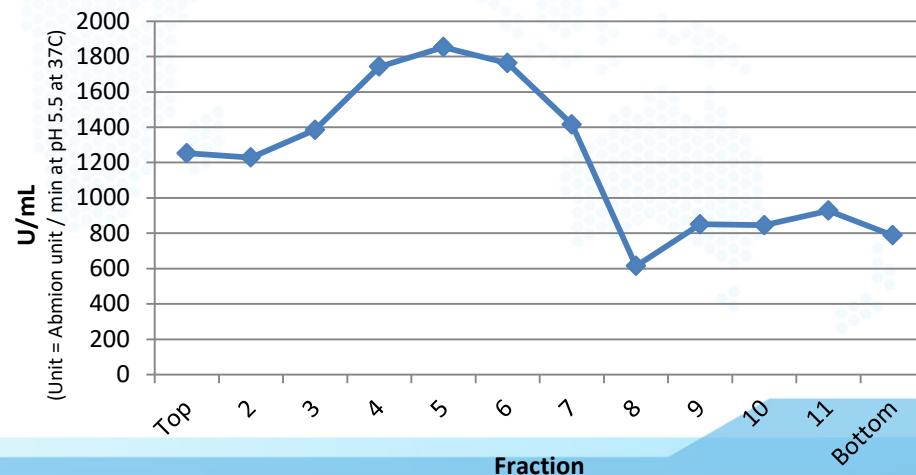
Acid Phosphatase



Cytochrome C Oxidase



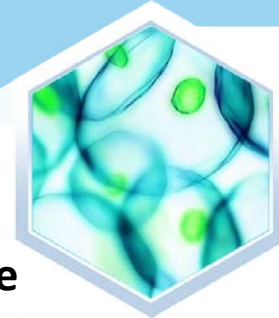
RNase



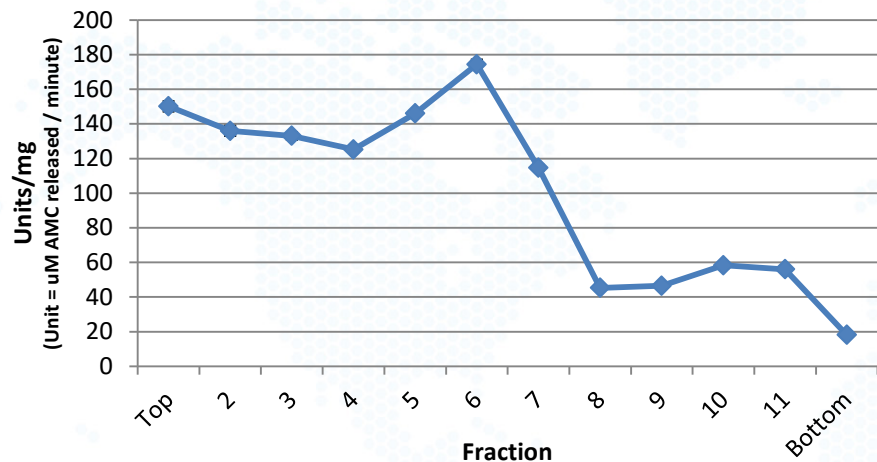


A BioIVT Company

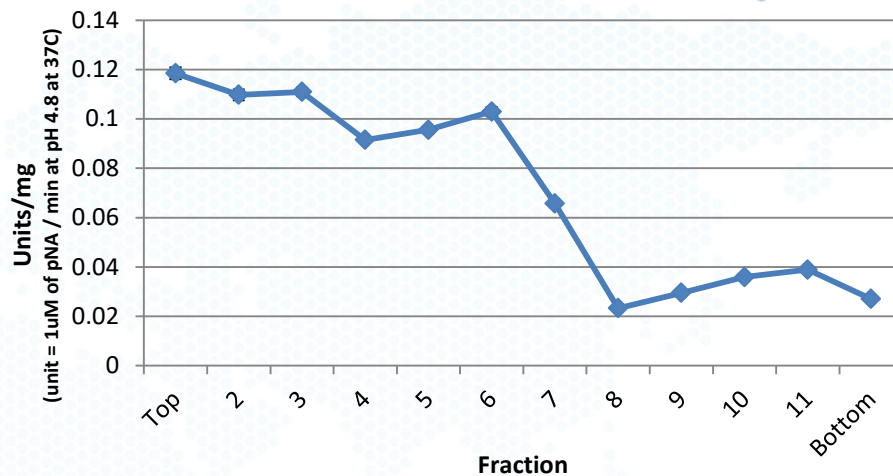
Identification of where the lysosome are in the gradient



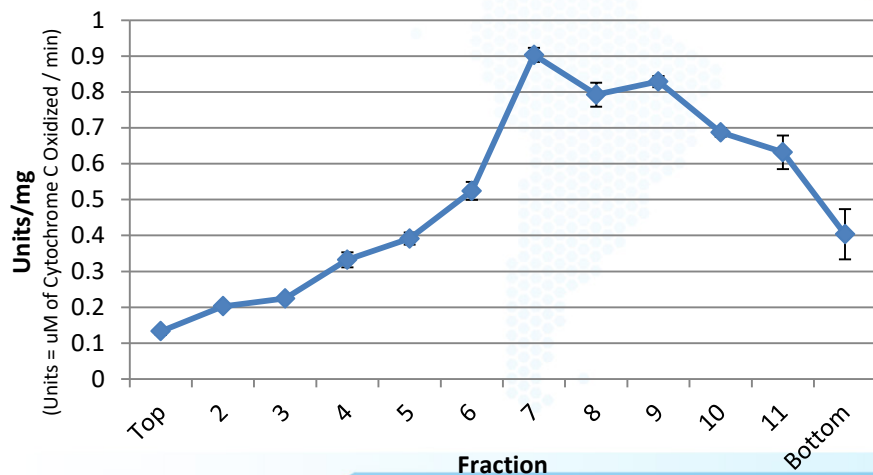
Cathepsin B



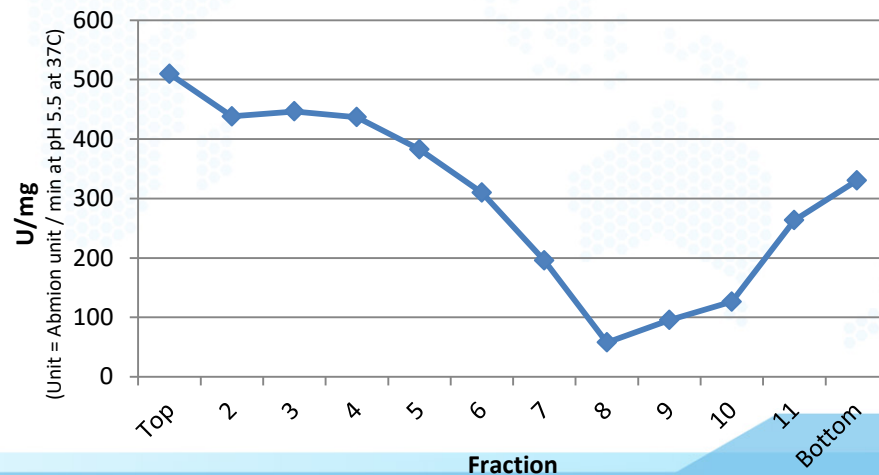
Acid Phosphatase

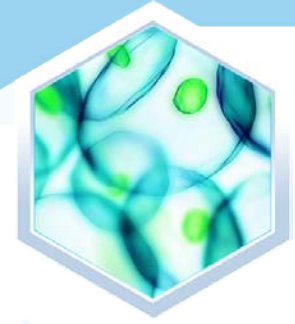


Cytochrome C Oxidase

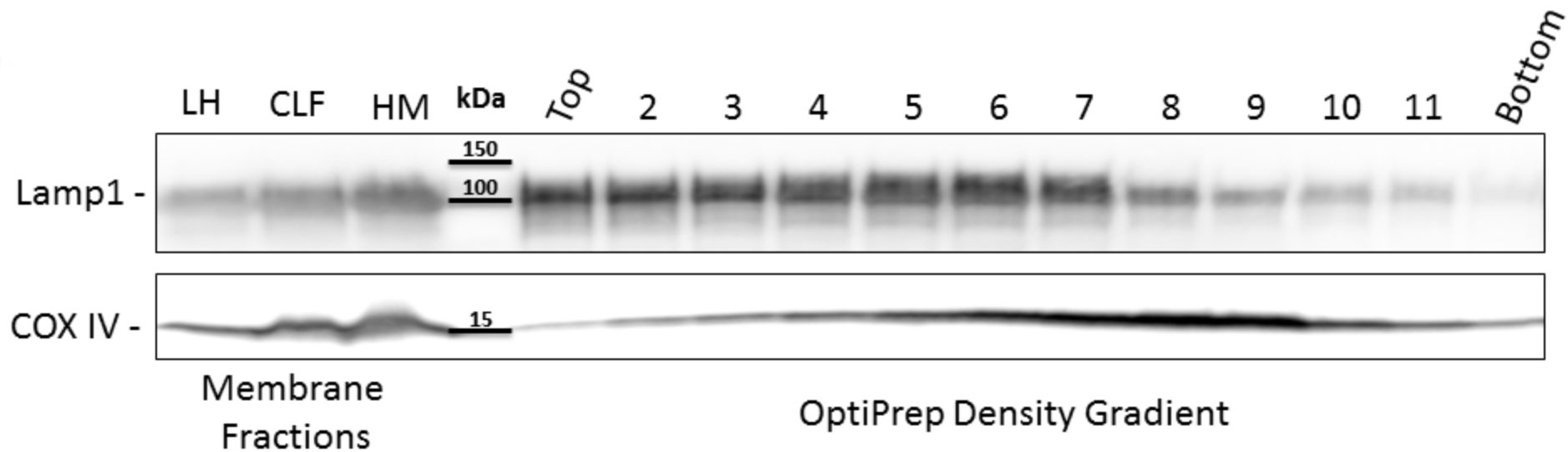


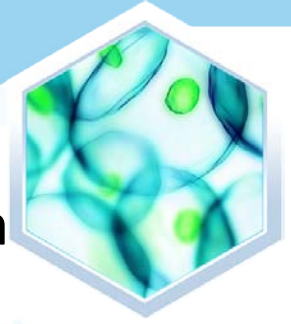
RNase



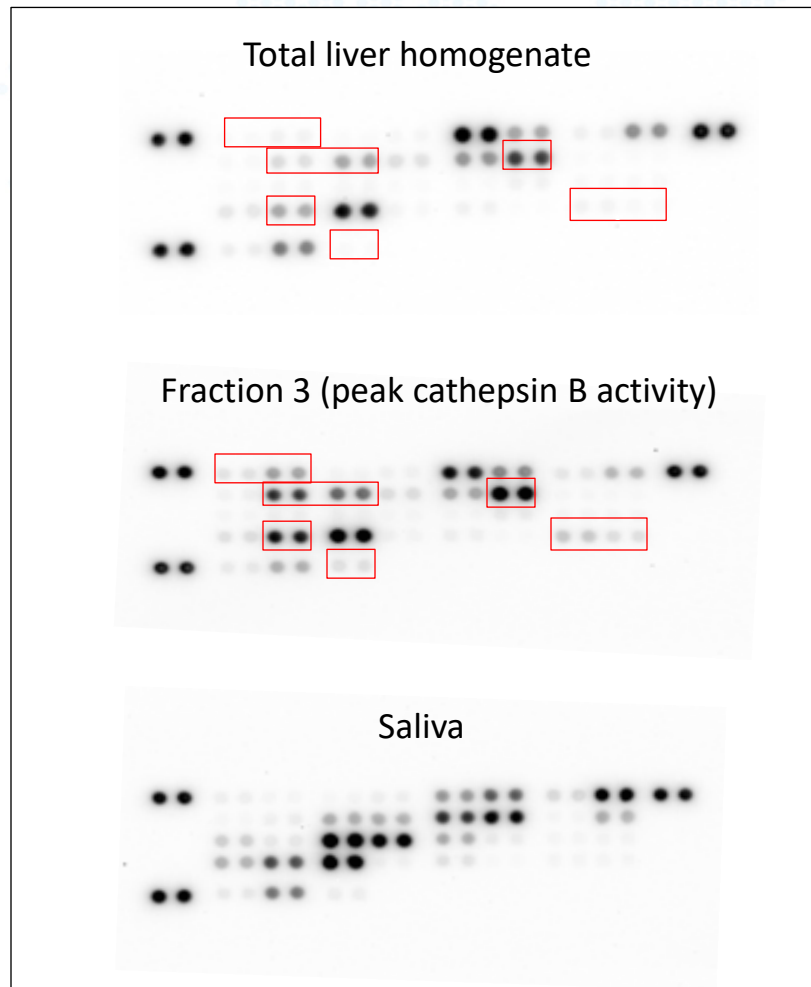


Identification of where the lysosome are in the gradient



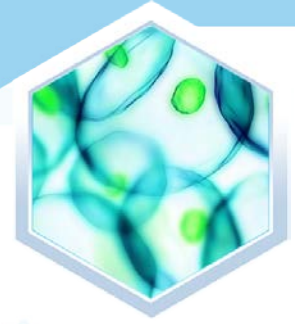


Proteases contained in fraction containing peak cathepsin



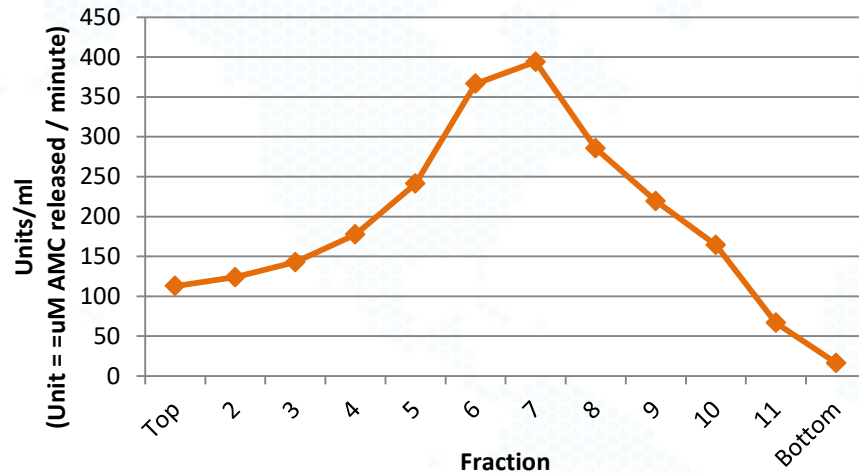
Cathepsin B-high activity fraction composed of disintegrin and metalloproteinase domain-containing proteins 8 and 9, cathepsins A, B, D, L, S and X/Z/P, dipeptidyl-peptidase 4, matrix metalloproteinases 8 and 9, neprilysin, presenilin and proteinase 3 (middle panel).

The gradient fractionation enriched cathepsins L and S, disintegrin and metalloproteinase domain-containing protein 8 and 9, dipeptidyl-peptidase 4, matrix metalloproteinase 8, Urokinase, neprilysin, and presenilin (boxed in red) but not the other proteases present in fraction 3.

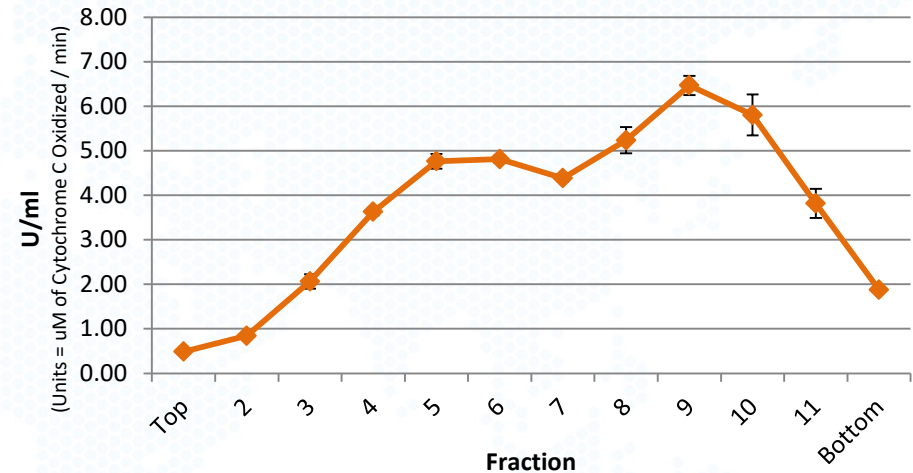


Donor/tissue variation

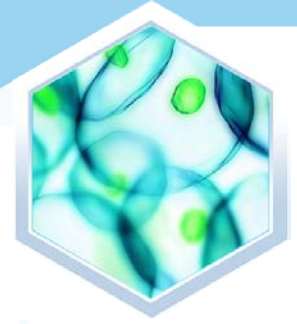
Cathepsin B



Cytochrome C Oxidase



We observed inter-donor variability in the separation of highest activities of the two enzymes

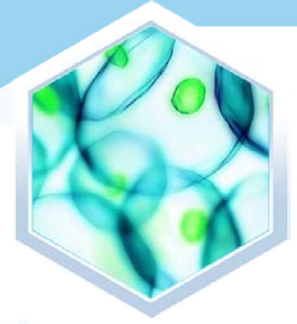


Conclusions

Purified human hepatic lysosomes show a 2-5 fold increase in lysosomal enzymatic activity per mg of total protein compared with initial liver homogenate.

Low contaminating activity from enzymes associated with mitochondria.

Serves as a *in vitro* reagent *with* a human matrix that compliments rat Tritosomes.



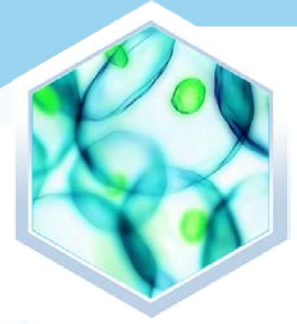
Further development of human hepatic lysosomes

Fractions 1 – 6 represent different densities, however they all contain significant amounts of Lamp1.

These fractions vary in density and likely vary in:

- enzymatic enzyme activity
- membrane composition
- Biological function

Are there different experimental conditions needed to assess catabolism with different classes of biopharmaceuticals?



What are the current uses?

In vitro diagnostic tool to conveniently and quickly evaluate potential changes in lysosomal stability due to targeted modifications of the biopharmaceutical/macromolecule during development

Help narrow and direct development tracks of biopharmaceutical

ADCs

siRNA/RNAi

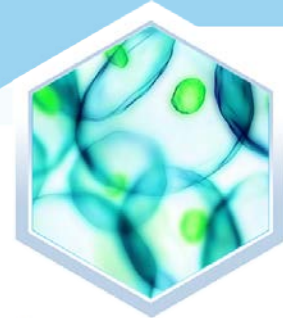
Biodegradable copolymers

Nonoparticles

**Pre-Register for the Upcoming Webinar on
Updated Lysosome Data, Lysosomal Trapping & More**
www.xenotech.com/scientific-resources/upcoming-webinars

View Our Other Available Webinars
www.xenotech.com/scientific-resources/webinar-series

Visit Our Lysosome & Tritosome Product Pages
www.xenotech.com/products/lysosomes-and-tritosomes
for more information or to purchase



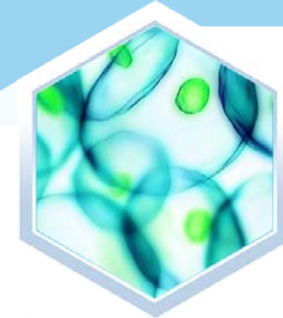
Available Product

H0610.L – Mixed Gender Human Liver Lysosomes, 0.25 mL

R0610.LT – Mixed Gender Rat Liver Tritosomes, 0.25 mL

Features & Benefits:

- Highly purified
- Characterized for lysosome specific enzymatic activity
- Less complex than *in vivo* models
- More representative than individually expressed/purified enzymes



Distributors

Europe



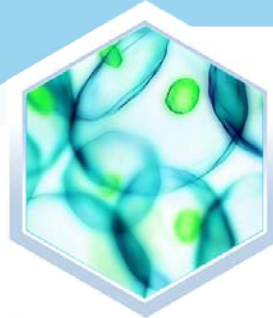
Jean-Francois Tetu, Ph.D.
Sales Manager – Cells and Cell-Based Assays
(T) +33 1 30 46 39 53
Jean-francois.tetu@tebu-bio.com
www.tebu-bio.com

Japan



Miki Fujishima
Sales & Marketing
Sekisui – Drug Development Solutions Center
(T) +81-3-3271-5634
smd-adme@sekisui.com

Complete list available online @ www.xenotech.com



2016-2017 Product Technical Guide



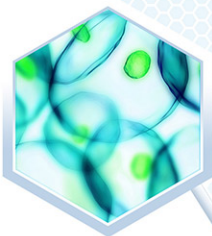
**Download the new
In Vitro Products & Reagents
Technology Guide**

Featuring...
a complete list of
XenoTech's products,
their uses & applications,
and protocols & handling
instructions.
([click here](#))

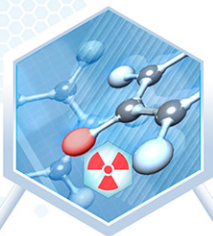
- Complete list of all XenoTech products
- Protocols, Applications, Handling Instruction

PROVEN GLOBAL CONTRACT RESEARCH EXPERTISE FROM DISCOVERY THROUGH CLINICAL SUPPORT

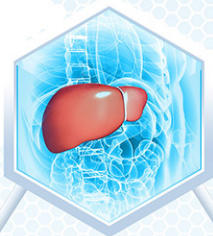
Cell & Tissue-Based Products



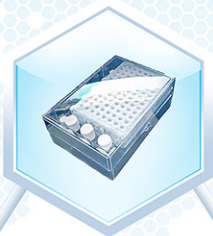
Radiolabeling



in vitro ADMET & Pharmacology



Metabolite ID & Production



Screening



API Manufacturing



in vivo ADMET & QWBA



Bioanalytical



EXPERTISE • EFFICIENCY • SUPPORT • PRECISION

Thank You!
XenoTech, LLC
913-438-7450
info@xenotechllc.com