



**XenoTech Offers Rat Tritosomes, lysosomes isolated from animals dosed with Tyloxapol (Triton WR 1339).**

## Tritosomes

Rat hepatic tritosomes are lysosomes that have been isolated from the livers of animals dosed with Tyloxapol (Triton WR 1339), a nonionic surfactant. Tyloxapol is taken up by hepatocytes through endocytosis and is trafficked to the lysosomal compartment. Tyloxapol-containing lysosomes exhibit decreased density that allows their efficient isolation from organelles that have overlapping densities with naive lysosomes (Leighton et al., 1968; Trouet, 1974). Density of lysosomes isolated from the livers of rats dosed with Tyloxapol also decreases due to an accumulation of very-low-density lipoproteins elevated in the serum of the animal in response to the administration of the detergent (Hayashi et al., 1982). This hyperlipidemic effect of Tyloxapol found its application in screening for hypolipidemic drugs in animals (Schurr et al., 1972). Tritosomes are well-established test system for study lysosomal composition, function, and diseases (Maguire et al., 1983; Lubke et al., 2009; Della Valle et al., 2011). More recently they have gained popularity as a reagent to assess catabolism of small molecular moieties conjugated to nano-particles, polymers, as well as, antibody-drug conjugates that are processed in the cell along the endosomal-lysosomal pathway (Hreczuk-Hirst et al., 2001; Malugin et al., 2007; Lim et al., 2013; Barrett et al., 2014).

Sekisui XenoTech rat tritosomes are prepared according to a discontinuous sucrose-gradient centrifugation method, based on Bagshaw et al., (2003). Human hepatic microsomes prepared from fresh human livers are a companion custom product to the rat tritosomes. Human lysosomal fractions are characterized for acid phosphatase and cathepsin B activities.

## References

- Bagshaw RD, Pasternak SH, Mahuran DJ, and Callahan JW (2003) Nicastrin is a resident lysosomal membrane protein. *Biochem Biophys Res Commun* **300**:615-618.
- Barrett SE, Abrams MT, Burke R, Carr BA, Crocker LS, Garbaccio RM, Howell BJ, Kemp EA, Kowtoniuk RA, Latham AH, Leander KR, Leone AM, Patel M, Pechenov S, Pudvah NT, Riley S, Sepp-Lorenzino L, Walsh ES, Williams JM, and Colletti SL (2014) An in vivo evaluation of amphiphilic, biodegradable peptide copolymers as siRNA delivery agents. *International journal of pharmaceutics* **466**:58-67.
- Della Valle MC, Sleat DE, Zheng H, Moore DF, Jadot M, and Lobel P (2011) Classification of subcellular location by comparative proteomic analysis of native and density-shifted lysosomes. *Molecular & cellular proteomics : MCP* **10**:M110 006403.
- Hayashi H, Shitara M, and Yamasaki F (1982) The origin of lipid accumulated in liver lysosomes after administration of triton WR-1339. *Journal of biochemistry* **92**:1585-1590.
- Hreczuk-Hirst D, Chicco D, German L, and Duncan R (2001) Dextrins as potential carriers for drug targeting: tailored rates of dextrin degradation by introduction of pendant groups. *International journal of pharmaceutics* **230**:57-66.
- Leighton F, Poole B, Beaufay H, Baudhuin P, Coffey JW, Fowler S, and De Duve C (1968) The large-scale separation of peroxisomes, mitochondria, and lysosomes from the livers of rats injected with triton WR-1339. Improved isolation procedures, automated analysis, biochemical and morphological properties of fractions. *The Journal of cell biology* **37**:482-513.
- Lim EK, Jang E, Lee K, Haam S, and Huh YM (2013) Delivery of cancer therapeutics using nanotechnology. *Pharmaceutics* **5**:294-317.
- Lubke T, Lobel P, and Sleat DE (2009) Proteomics of the lysosome. *Biochim Biophys Acta* **1793**:625-635.
- Maguire GA, Docherty K, and Hales CN (1983) Sugar transport in rat liver lysosomes. Direct demonstration by using labelled sugars. *Biochem J* **212**:211-218.
- Malugin A, Kopeckova P, and Kopecek J (2007) Liberation of doxorubicin from HPMA copolymer conjugate is essential for the induction of cell cycle arrest and nuclear fragmentation in ovarian carcinoma cells. *Journal of controlled release : official journal of the Controlled Release Society* **124**:6-10.
- Schurr PE, Schultz JR, and Parkinson TM (1972) Triton-induced hyperlipidemia in rats as an animal model for screening hypolipidemic drugs. *Lipids* **7**:68-74.
- Trouet A (1974) Isolation of modified liver lysosomes. *Methods Enzymol* **31**:323-329.

## Packaging & Characterization

**Product ID - R0610.LT**

**Protein Concentration - 2.7 mg/mL**

**Volume - 0.25 mL**

**Pool Size - 60 animals**

Tritosomes are characterized for acid phosphatase and RNase activity.

Sekisui XenoTech also offers **K5200** 10x catabolism buffer that has been formulated and optimized to extract the most *in vitro* catabolic performance from isolated human lysosomes and rat tritosomes (figure 1 of lysosome flyer). All of the catabolic data presented by Sekisui XenoTech utilized this buffer and is crucial for full catabolic activity.



**Band of discontinuous sucrose-gradient isolated rat tritosomes**

