

Lysosome Isolation Kit ab234047

12 References 2 Images

Overview

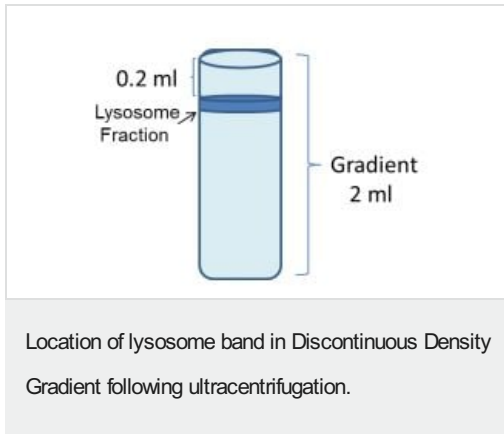
Product name	Lysosome Isolation Kit
Sample type	Tissue, Adherent cells, Suspension cells
Product overview	Lysosome Isolation Kit (ab234047) provides a procedure for isolating an enriched or purified lysosomal fraction from animal tissues and cultured cells by differential centrifugation followed by density gradient centrifugation.
Notes	<p>This product is manufactured by BioVision, an Abcam company and was previously called K235 Lysosome Purification Kit. K235-50 is the same size as the 50 test size of ab234047.</p> <p>Lysosomes are membrane-bound cell organelles found in most animal cells (except red blood cells). Lysosomes are spherical vesicles containing hydrolytic enzymes capable of breaking down different kinds of biomolecules, including proteins, nucleic acids, carbohydrates, lipids, and cellular debris. They are known to contain over 50 different enzymes, which are active under acidic conditions. Thus, lysosomes act as the waste disposal system of the cells by digesting unwanted materials in the cytoplasm.</p>

Properties

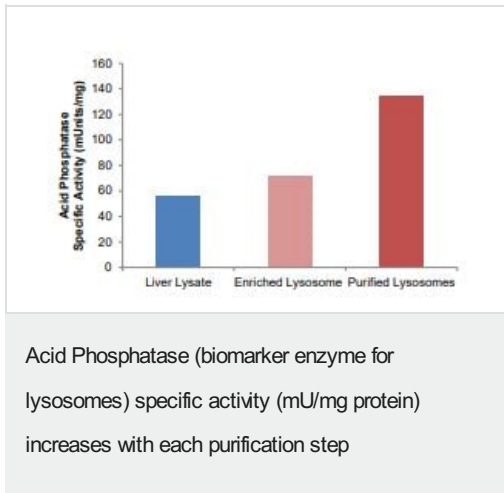
Storage instructions Store at -20°C. Please refer to protocols.

Components	50 tests
Lysosome Enrichment Buffer	1 x 100ml
Density Gradient Media II	1 x 85ml
Lysosome Isolation Buffer	1 x 25ml
Protease Inhibitor Cocktail II	1 x 1ml

Images



Tentative Lysosome band location in gradient density.



Data is normalized by the amount of protein used per well.

Please note: All products are "FOR RESEARCH USE ONLY. NOT FOR USE IN DIAGNOSTIC PROCEDURES"

Our Abpromise to you: Quality guaranteed and expert technical support

- Replacement or refund for products not performing as stated on the datasheet
- Valid for 12 months from date of delivery
- Response to your inquiry within 24 hours
- We provide support in Chinese, English, French, German, Japanese and Spanish
- Extensive multi-media technical resources to help you
- We investigate all quality concerns to ensure our products perform to the highest standards

If the product does not perform as described on this datasheet, we will offer a refund or replacement. For full details of the Abpromise, please visit <https://www.abcam.com/abpromise> or contact our technical team.

Terms and conditions

- Guarantee only valid for products bought direct from Abcam or one of our authorized distributors