

Human LAMP2 ELISA Kit ab277449

1 Image

Overview

Product name	Human LAMP2 ELISA Kit				
Detection method	Colorimetric				
Precision	Intra-assay				
	Sample	n	Mean	SD	CV%
	Overall				< 10%
	Inter-assay				
	Sample	n	Mean	SD	CV%
	Overall				< 12%
Sample type	Cell culture supernatant, Serum, Plasma				
Assay type	Sandwich (quantitative)				
Sensitivity	0.3 ng/ml				
Range	0.307 ng/ml - 75 ng/ml				
Recovery	Sample specific recovery				
	Sample type		Average %		Range
	Serum		107.2		86% - 129%
	Plasma		105.1		93% - 117%
	Cell culture media		113		103% - 130%
Assay duration	Multiple steps standard assay				
Species reactivity	Reacts with: Human				
Product overview	Human LAMP2 ELISA Kit (ab277449) is an <i>in-vitro</i> enzyme-linked immunosorbent assay for the quantitative measurement of Human LAMP2 in serum, plasma and cell culture supernatants.				

This assay employs an antibody specific for Human LAMP2 coated on a 96-well plate. Standards and samples are pipetted into the wells and Human LAMP2 present in a sample is bound to the

wells by the immobilized antibody. The wells are washed, and biotinylated anti-Human LAMP2 antibody is added. After washing away unbound biotinylated antibody, HRP-conjugated streptavidin is pipetted to the wells. The wells are again washed, a TMB substrate solution is added to the wells and color develops in proportion to the amount of Human LAMP2 bound. The Stop Solution changes the color from blue to yellow, and the intensity of the color is measured at 450 nm.

Platform Pre-coated microplate (12 x 8 well strips)

Properties

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

Components	1 x 96 tests
150X HRP-Streptavidin Concentrate	1 x 200µl
20X Wash Buffer	1 x 25ml
5X Assay Diluent B	1 x 15ml
5X Assay Diluent D	1 x 15ml
Biotinylated Anti-Human LAMP2 Antibody	2 vials
Anti-Human LAMP2 coated Microplate	1 unit
Human LAMP2 standard protein (Lyophilized)	2 vials
Stop Solution	1 x 8ml
TMB One-Step Substrate Reagent	1 x 12ml

Function Implicated in tumor cell metastasis. May function in protection of the lysosomal membrane from autodigestion, maintenance of the acidic environment of the lysosome, adhesion when expressed on the cell surface (plasma membrane), and inter-and intracellular signal transduction. Protects cells from the toxic effects of methylating mutagens.

Tissue specificity Isoform LAMP-2A is highly expressed in placenta, lung and liver, less in kidney and pancreas, low in brain and skeletal muscle. Isoform LAMP-2B is highly expressed in skeletal muscle, less in brain, placenta, lung, kidney and pancreas, very low in liver.

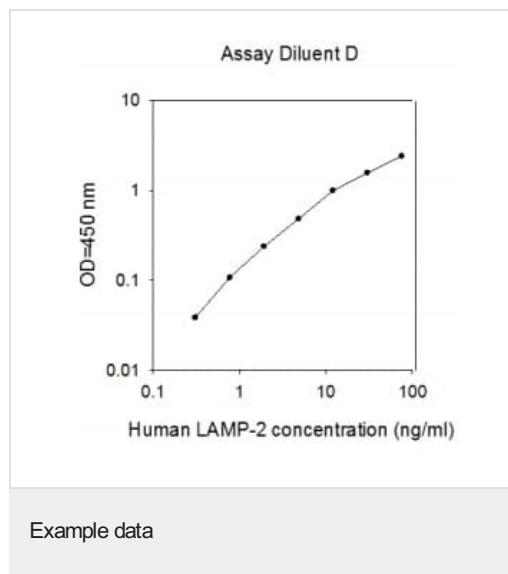
Involvement in disease Defects in LAMP2 are the cause of Danon disease (DAND) [MIM:300257]; also known as glycogen storage disease type 2B (GSD2B). DAND is a lysosomal glycogen storage disease characterized by the clinical triad of cardiomyopathy, vacuolar myopathy and mental retardation. It is often associated with an accumulation of glycogen in muscle and lysosomes.

Sequence similarities Belongs to the LAMP family.

Post-translational modifications O- and N-glycosylated; some of the 16 N-linked glycans are polylactosaminoglycans.

Cellular localization Cell membrane. Endosome membrane. Lysosome membrane. This protein shuttles between lysosomes, endosomes, and the plasma membrane.

Images



These standard curves are for demonstration only. A standard curve must be run with each assay.

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