

Product datasheet

Human LYVE ELISA Kit ab100585

★★★★★ 1 Abreviews 3 Images

Overview

Product name	Human LYVE ELISA Kit
Detection method	Colorimetric
Sample type	Cell culture supernatant, Serum, Plasma
Assay type	Sandwich (quantitative)
Sensitivity	< 8 pg/ml
Range	8.19 pg/ml - 2000 pg/ml
Recovery	78 %

Sample specific recovery

Sample type	Average %	Range
Cell culture supernatant	76	69% - 85%
Serum	76.38	67% - 87%
Plasma	79.07	68% - 86%

Assay duration Multiple steps standard assay

Species reactivity **Reacts with:** Human

Product overview Abcam's Human LYVE ELISA (Enzyme-Linked Immunosorbent Assay) kit is an *in vitro* enzyme-linked immunosorbent assay for the quantitative measurement of Human LYVE in serum, plasma, and cell culture supernatants.

This assay employs an antibody specific for Human LYVE coated on a 96-well plate. Standards and samples are pipetted into the wells and LYVE present in a sample is bound to the wells by the immobilized antibody. The wells are washed and biotinylated anti-Human LYVE 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 LYVE bound. The Stop Solution changes the color from blue to yellow, and the intensity of the color is measured at 450 nm.

Notes Optimization may be required with urine samples.

Platform Microplate

Properties

Storage instructions

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

Components	1 x 96 tests
200X HRP-Streptavidin Concentrate	1 x 200µl
20X Wash Buffer	1 x 25ml
5X Assay Diluent B	1 x 15ml
Assay Diluent A	1 x 30ml
Biotinylated anti-Human LYVE	2 vials
LYVE Microplate (12 x 8 wells)	1 unit
Recombinant Human LYVE Standard (lyophilized)	2 vials
Stop Solution	1 x 8ml
TMB One-Step Substrate Reagent	1 x 12ml

Function

Ligand-specific transporter trafficking between intracellular organelles (TGN) and the plasma membrane. Plays a role in autocrine regulation of cell growth mediated by growth regulators containing cell surface retention sequence binding (CRS). May act as a hyaluronan (HA) transporter, either mediating its uptake for catabolism within lymphatic endothelial cells themselves, or its transport into the lumen of afferent lymphatic vessels for subsequent re-uptake and degradation in lymph nodes.

Tissue specificity

Mainly expressed in endothelial cells lining lymphatic vessels.

Sequence similarities

Contains 1 Link domain.

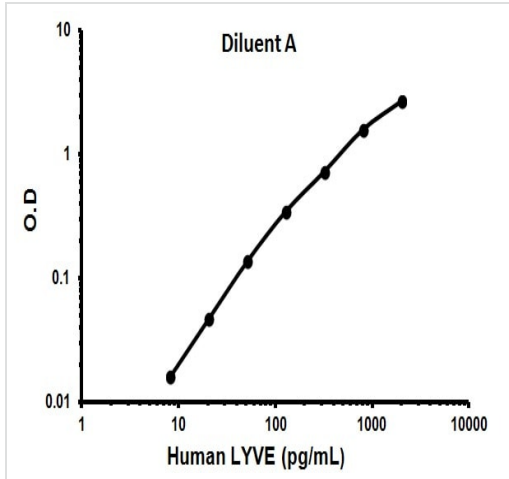
Post-translational modifications

O-glycosylated.

Cellular localization

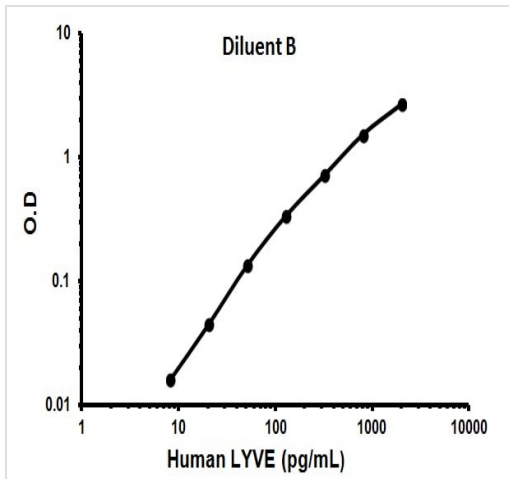
Membrane. Localized to the plasma membrane and in vesicles near extranuclear membranes which may represent trans-Golgi network (TGN) and endosomes/prelysosomal compartments. Undergoes ligand-dependent internalization and recycling at the cell surface.

Images



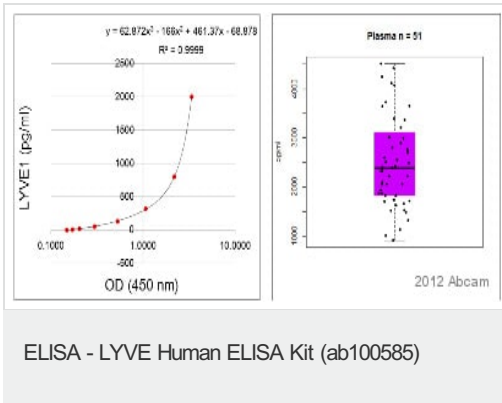
Representative Standard Curve using ab100585.

Typical Standard Curve



Representative Standard Curve using ab100585.

Typical Standard Curve



ab100585 was used to detect LYVE in Human heparin plasma samples according to the kit protocol. 10 µL sample was used per well.

This image is courtesy of an anonymous Abreview

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