

Human Legumain ELISA Kit ab267603

1 Image

Overview

Product name	Human Legumain 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	28.7 pg/ml				
Range	28.67 pg/ml - 7000 pg/ml				
Recovery	Sample specific recovery				
	Sample type		Average %		Range
	Serum		103.4		69% - 129%
	Plasma		109.1		82% - 144%
	Cell culture media		124.4		118% - 133%
Assay duration	Multiple steps standard assay				
Species reactivity	Reacts with: Human				
Product overview	Human Legumain ELISA Kit is designed for the quantitative determination of Legumain in cell culture supernatants, plasma and serum samples.				

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

the wells by the immobilized antibody. The wells are washed and biotinylated anti-human Legumain 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 Legumain bound. The Stop Solution changes the color from blue to yellow, and the intensity of the color is measured at 450 nm.

This ELISA kit shows no cross-reactivity with the following cytokines tested: human 2B4 (CD244), ADAM-9, Angiopoietin-2, APRIL, BMP-2, BMP-9, C5a, Cathepsin L, CD200, CD97, Chemerin, DcR3, FABP2, FAP, FGF-19, Galectin-3, HGFR, IFN-alpha / betaR2, IGF-2, IGF-2 R, IL-1 R6 (IL-1 Rrp2), IL-24, IL-33 (IL-1 F11), Kallikrein 14.

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
200X HRP-Streptavidin Concentrate	1 x 200µl
20X Wash Buffer	1 x 25ml
5X Assay Diluent	1 x 15ml
Anti-Human Legumain coated Microplate (12 x 8 wells)	1 unit
Biotinylated Anti-Human Legumain Detection Antibody	2 vials
Human Legumain Standard (Lyophilized)	2 vials
Stop Solution	1 x 8ml
TMB Substrate Solution	1 x 12ml

Function Has a strict specificity for hydrolysis of asparaginyl bonds. Can also cleave aspartyl bonds slowly, especially under acidic conditions. May be involved in the processing of proteins for MHC class II antigen presentation in the lysosomal/endosomal system.

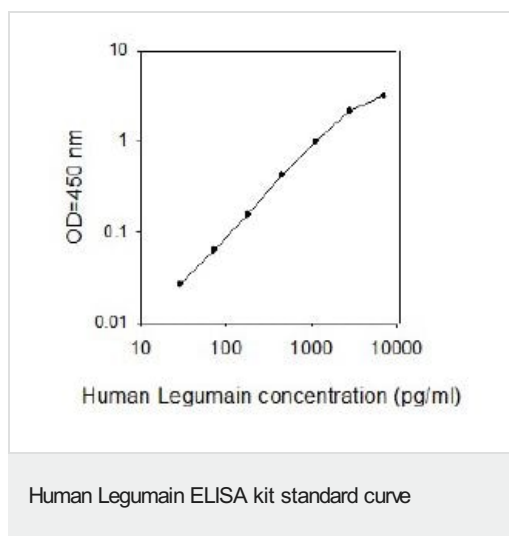
Tissue specificity Ubiquitous. Particularly abundant in kidney, heart and placenta.

Sequence similarities Belongs to the peptidase C13 family.

Post-translational modifications Glycosylated.

Cellular localization Lysosome.

Images



This standard curve is 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