

Human ULBP3 ELISA Kit ab267617

1 Image

Overview

Product name	Human ULBP3 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	1.3 pg/ml				
Range	1.229 pg/ml - 300 pg/ml				
Recovery	Sample specific recovery				
	Sample type		Average %		Range
	Serum		81.54		74% - 92%
	Plasma		77.77		72% - 84%
	Cell culture media		77.14		69% - 80%
Assay duration	Multiple steps standard assay				
Species reactivity	Reacts with: Human				
Product overview	Human ULBP3 ELISA Kit is designed for the quantitative determination of ULBP3 in cell culture supernatants, plasma and serum samples.				

This assay employs an antibody specific for human ULBP3 coated on a 96-well plate. Standards and samples are pipetted into the wells and ULBP3 present in a sample is bound to the wells by

the immobilized antibody. The wells are washed and biotinylated anti-human ULBP3 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 ULBP3 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 AMIGO, Aminopeptidase LRAP (ERAP2), Amnionless, Arylsulfatase A, Bcl-w, CD109, CD157, CD34, CD83, CLEC1, CLEC10A (CD301), CMG-2, CREG, Cystatin SN, Cytokeratin-8, Dectin-1 (CLEC7A), Desmocollin-3, Endoglycan, Galectin-4, HAPLN1, Jagged 1, Langerin (CD207), Lumican, Matriptase (ST14), MEP1B, Nectin-3, OX40 (TNFRSF4), OX40 Ligand (TNFSF4), p27, Pappalysin2, Plexin B3, Plexin D1, proGRP, PSA-total, Reg1B, RGM-A, ROBO2, Spinesin, TWEAK R (TNFRSF12).

Platform Microplate

Properties

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

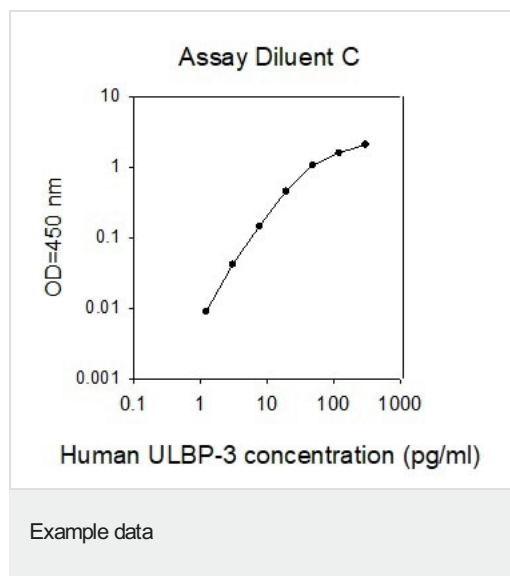
Components	1 x 96 tests
20X Wash Buffer	1 x 25ml
400X HRP-Streptavidin Concentrate	1 x 200µl
5X Assay Diluent B	1 x 15ml
Anti-Human ULBP3 coated Microplate (12 x 8 wells)	1 unit
Assay Diluent C	1 x 30ml
Biotinylated Anti-Human ULBP3 Detection Antibody	2 vials
Human ULBP3 Standard (Lyophilized)	2 vials
Stop Solution	1 x 8ml
TMB Substrate Solution	1 x 12ml

Function Ligand for the KLRK1/NKG2D receptor, together with at least ULBP1 and ULBP2. ULBPs activate multiple signaling pathways in primary NK cells, resulting in the production of cytokines and chemokines. Binding of ULBPs ligands to KLRK1/NKG2D induces calcium mobilization and activation of the JAK2, STAT5, ERK and PI3K kinase/Akt signal transduction pathway. Has lower affinity for KLRK1/NKG2D compared to ULBP1 and ULBP2 and induces weaker signaling responses than does ULBP2 or ULBP1.

Sequence similarities Belongs to the MHC class I family.

Cellular localization Cell membrane.

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