

Mouse LBP ELISA Kit ab213876

5 References 1 Image

Overview

Product name	Mouse LBP ELISA Kit			
Detection method	Colorimetric			
Precision	Intra-assay			
	Sample	n	Mean	SD
	1	16	3.313ng/ml	0.237
	2	16	12.997ng/ml	0.832
	3	16	39.089ng/ml	2.111
	Inter-assay			
	Sample	n	Mean	SD
	1	24	3.385ng/ml	0.244
	2	24	14.212ng/ml	1.151
	3	24	39.495ng/ml	2.804
Sample type	Cell culture supernatant, Serum, Cell Lysate, Hep Plasma, EDTA Plasma			
Assay type	Sandwich (quantitative)			
Sensitivity	< 50 pg/ml			
Range	1.56 ng/ml - 100 ng/ml			
Assay time	3h 30m			
Assay duration	Multiple steps standard assay			
Species reactivity	Reacts with: Mouse			
Product overview	The Mouse LBP Enzyme-Linked Immunosorbent Assay (ELISA) kit (ab213876) is designed for the quantitative detection of Mouse LBP in cell culture supernatants, cell lysates, serum and plasma (heparin, EDTA).			

The ELISA kit is based on standard sandwich enzyme-linked immunosorbent assay technology. A

polyclonal antibody from goat specific for LBP has been pre-coated onto 96-well plates. Standards (Expression system for standard: NSO; Immunogen sequence: G25-V481) and test samples are added to the wells, a biotinylated detection polyclonal antibody from goat specific for LBP is added subsequently and then followed by washing with PBS or TBS buffer. Avidin-Biotin-Peroxidase Complex was added and unbound conjugates were washed away with PBS or TBS buffer. HRP substrate TMB was used to visualize HRP enzymatic reaction. TMB was catalyzed by HRP to produce a blue color product that changed into yellow after adding acidic stop solution. The density of yellow is proportional to the Mouse LBP amount of sample captured in plate.

Notes Lipopolysaccharide binding protein is a protein that in humans is encoded by the LBP gene. This gene is mapped to 20q11.23. LBP is a soluble acute-phase protein that binds to bacterial lipopolysaccharide (or LPS) to elicit immune responses by presenting the LPS to important cell surface pattern recognition receptors called CD14 and TLR4. It is present in the cerebrospinal fluid of patients with pneumococcal meningitis. The protein encoded by this gene is involved in the acute phase immunologic response to gram-negative bacterial infections. LBP is made in the liver during the acute phase of infections and is thought to function as a carrier for LPS and to help control LPS dependent monocyte responses.

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

Properties

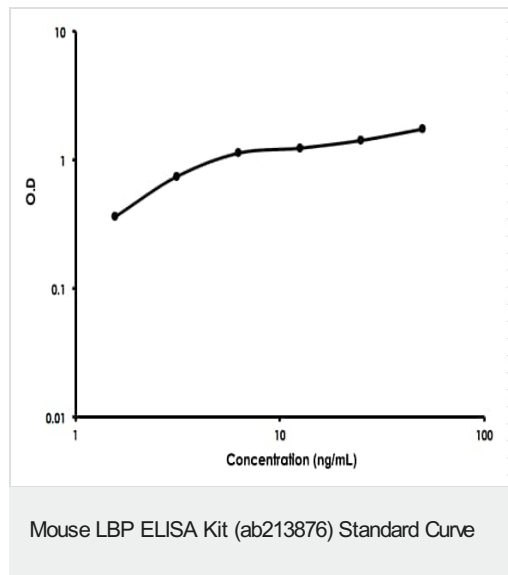
Storage instructions Store at -20°C. Please refer to protocols.

Components	Identifier	1 x 96 tests	1 x 96 tests
ABC Diluent Buffer	Blue Cap	1 x 12ml	1 x 12ml
Adhesive Plate Seal		4 units	4 units
Antibody Diluent Buffer	Green Cap	1 x 12ml	1 x 12ml
Anti-Mouse LBP coated Microplate (12 x 8 wells)		1 unit	1 unit
Avidin-Biotin-Peroxidase Complex (ABC)		1 x 100µl	1 x 100µl
Biotinylated anti- Mouse LBP antibody		1 x 100µl	1 x 100µl
Lyophilized recombinant Mouse LBP standard		2 vials	2 vials
Sample Diluent Buffer	Green Cap	1 x 30ml	1 x 30ml
TMB Color Developing Agent	Black Cap	1 x 10ml	1 x 10ml
TMB Stop Solution	Yellow Cap	1 x 10ml	1 x 10ml
Wash Buffer (25X)		1 x 20ml	1 x 20ml

Function Binds to the lipid A moiety of bacterial lipopolysaccharides (LPS), a glycolipid present in the outer membrane of all Gram-negative bacteria. The LBP/LPS complex seems to interact with the CD14 receptor.

Sequence similarities Belongs to the BPI/LBP/Plunc superfamily. BPI/LBP family.

Images



Mouse LBP ELISA Kit (ab213876) Standard Curve

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