

Product datasheet

Mouse Cathepsin B ELISA Kit ab119585

★★★★★ 1 Abreviews 1 References 1 Image

Overview

Product name	Mouse Cathepsin B ELISA Kit
Detection method	Colorimetric
Sample type	Cell culture supernatant, Tissue Extracts
Assay type	Sandwich (quantitative)
Sensitivity	< 5 pg/well
Range	156 pg/well - 10000 pg/well
Assay duration	Multiple steps standard assay
Species reactivity	Reacts with: Mouse Does not react with: Human

Product overview

Abcam's Mouse Cathepsin B *in vitro* ELISA (Enzyme-Linked Immunosorbent Assay) kit is designed for the accurate quantitative measurement of Mouse Cathepsin B in cell culture supernatants, serum and plasma (heparin, EDTA).

A Cathepsin B specific rat monoclonal antibody has been precoated onto 96-well plates. Standards and test samples are added to the wells and incubated. A biotinylated detection polyclonal antibody from goat, specific for Cathepsin B is then added followed by washing with PBS or TBS buffer. Avidin-Biotin-Peroxidase Complex is added and unbound conjugates are washed away with PBS or TBS buffer. TMB is then used to visualize the HRP enzymatic reaction. TMB is catalyzed by HRP to produce a blue color product that changes into yellow after adding acidic stop solution. The density of yellow coloration is directly proportional to the mouse Cathepsin B amount of sample captured in plate.

Platform

Microplate

Properties

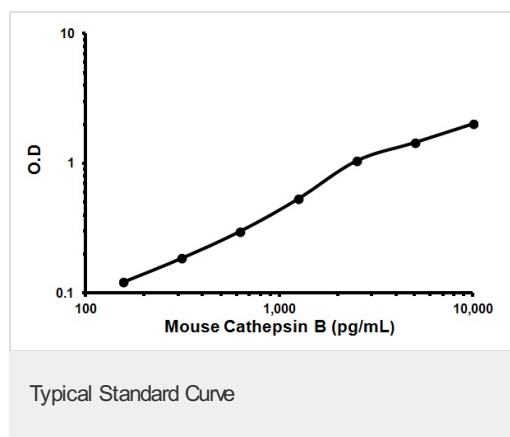
Storage instructions Store at -20°C. Please refer to protocols.

Components	Identifier	1 x 96 tests
ABC Diluent Buffer	Blue Cap	1 x 12ml
Antibody Diluent Buffer	Green Cap	1 x 12ml
Anti-mouse Cathepsin B antibody Microplate (12 x 8 wells)		1 unit

Components	Identifier	1 x 96 tests
Avidin-Biotin-Peroxidase Complex (ABC)		1 x 130µl
Biotinylated anti-mouse Cathepsin B antibody		1 x 130µl
Lyophilized recombinant mouse Cathepsin B standard		2 x 10ng
Plate Seal		1 x 4 units
Sample Diluent Buffer	Green Cap	1 x 30ml
TMB Color Developing Agent	Amber Bottle	1 x 10ml
TMB Stop Solution	Yellow Cap	1 x 10ml

Function	Thiol protease which is believed to participate in intracellular degradation and turnover of proteins. Has also been implicated in tumor invasion and metastasis.
Sequence similarities	Belongs to the peptidase C1 family.
Cellular localization	Lysosome. Melanosome. Identified by mass spectrometry in melanosome fractions from stage I to stage IV.

Images



Representative Standard Curve Using ab119585.

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