

Mouse Cathepsin B ELISA Kit ab119585

★★★★★ [1 Abreviews](#) [1 References](#) [1 Image](#)

Overview

Product name Mouse Cathepsin B ELISA Kit

Detection method Colorimetric

Precision Intra-assay

Sample	n	Mean	SD	CV%
1	16	242pg/ml	14.76	6.1%
2	16	1828pg/ml	100.54	5.5%
3	16	4749pg/ml	370.42	7.8%

Inter-assay

Sample	n	Mean	SD	CV%
1	24	229pg/ml	13.96	6.1%
2	24	1916pg/ml	128.37	6.7%
3	24	4534pg/ml	439.79	9.7%

Sample type Cell culture supernatant, Serum, Cell Lysate, Hep Plasma, EDTA Plasma

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.

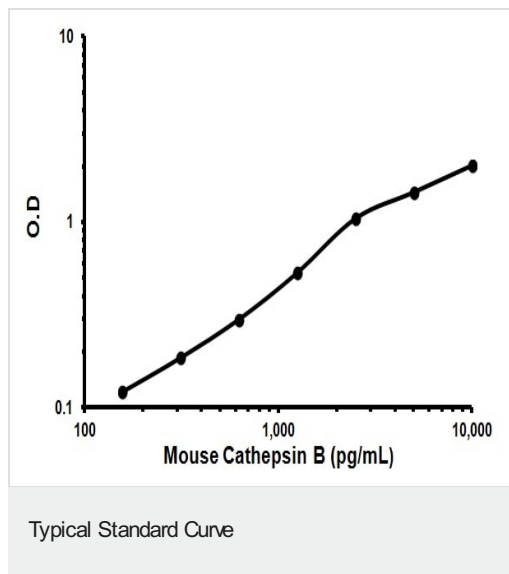
Components	Identifier	1 x 96 tests	1 x 96 tests
ABC Diluent Buffer	Blue Cap	1 x 12ml	1 x 12ml
Antibody Diluent Buffer	Green Cap	1 x 12ml	1 x 12ml
Anti-mouse Cathepsin B antibody Microplate (12 x 8 wells)		1 unit	1 unit
Avidin-Biotin-Peroxidase Complex (ABC)		1 x 100µl	1 x 100µl
Biotinylated anti-mouse Cathepsin B antibody		1 x 100µl	1 x 100µl
Lyophilized recombinant mouse Cathepsin B standard		2 x 10ng	2 x 10ng
Plate Seal		1 x 4 units	1 x 4 units
Sample Diluent Buffer	Green Cap	1 x 30ml	1 x 30ml
TMB Color Developing Agent	Amber Bottle	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 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