

Mouse Alkaline Phosphatase ELISA Kit ab285274

[1 Image](#)

Overview

Product name Mouse Alkaline Phosphatase ELISA Kit

Detection method Colorimetric

Precision

Intra-assay

Sample	n	Mean	SD	CV%
General				< 8%

Inter-assay

Sample	n	Mean	SD	CV%
General				< 10%

Sample type Cell culture supernatant, Serum, Plasma

Assay type Sandwich (quantitative)

Sensitivity < 18.75 pg/ml

Range 31.25 pg/ml - 2000 pg/ml

Assay duration Multiple steps standard assay

Species reactivity **Reacts with:** Mouse

Product overview The Mouse ALP ELISA kit (ab285274, E4573) is a sandwich ELISA assay for the quantitative measurement of ALP in mouse serum, plasma and cell culture supernatants. The density of color is proportional to the amount of ALP captured from the samples.

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

Properties

Storage instructions Store at +4°C. Please refer to protocols.

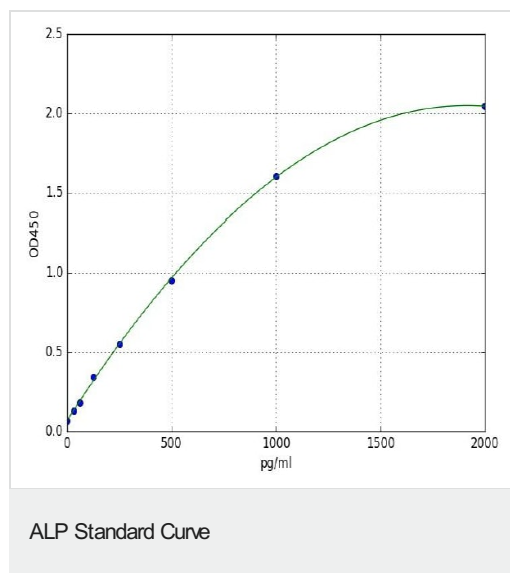
Components	96 tests
25X Wash buffer	1 x 30ml

Components	96 tests
Antibody dilution buffer	1 x 10ml
Biotin/Detection antibody (Concentrated)	1 x 120µl
HRP-Streptavidin Conjugate (SABC)	1 x 120µl
Lyophilized Standard	1 x 2 vials
Micro ELISA Plate	1 unit
Plate Sealer	1 x 5 units
SABC dilution buffer	1 x 10ml
Sample / Standard Dilution Buffer	1 x 20ml
Stop Solution	1 x 10ml
TMB Substrate	1 x 10ml

Sequence similarities Belongs to the alkaline phosphatase family.

Cellular localization Cell membrane.

Images



Standard curve analysis. For reference use only.

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