

Mouse CD30 ELISA Kit ab113341

2 Images

Overview

Product name	Mouse CD30 ELISA Kit
Detection method	Colorimetric
Sample type	Cell culture supernatant, Serum, Plasma
Assay type	Sandwich (quantitative)
Sensitivity	< 10 pg/ml
Range	12.29 pg/ml - 3000 pg/ml
Recovery	76 %

Sample specific recovery

Sample type	Average %	Range
Cell culture supernatant	77.29	69% - 87%
Serum	75.99	68% - 84%
Plasma	74.88	67% - 83%

Assay duration Multiple steps standard assay

Species reactivity **Reacts with:** Mouse

Product overview Abcam's CD30 Mouse ELISA (Enzyme-Linked Immunosorbent Assay) kit is an *in vitro* enzyme-linked immunosorbent assay for the quantitative measurement of mouse CD30 in serum, plasma and cell culture supernatants.

This assay employs an antibody specific for mouse CD30 coated on a 96- well plate. Standards and samples are pipetted into the wells and CD30 present in a sample is bound to the wells by the immobilized antibody. The wells are washed and biotinylated anti-mouse CD30 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 CD30 bound. The Stop Solution changes the color from blue to yellow, and the intensity of the color is measured at 450 nm.

Platform Microplate

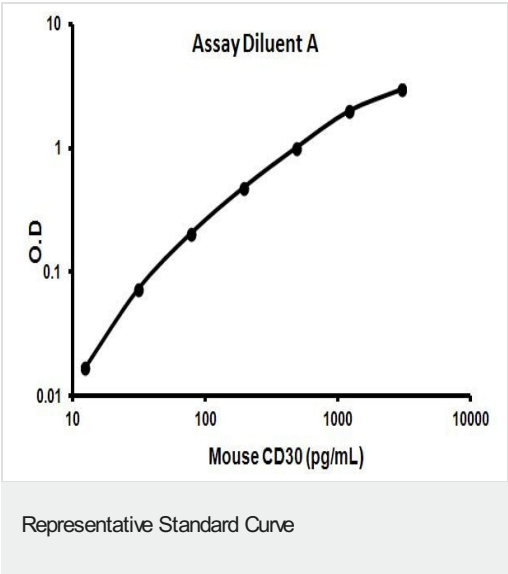
Properties

Storage instructions

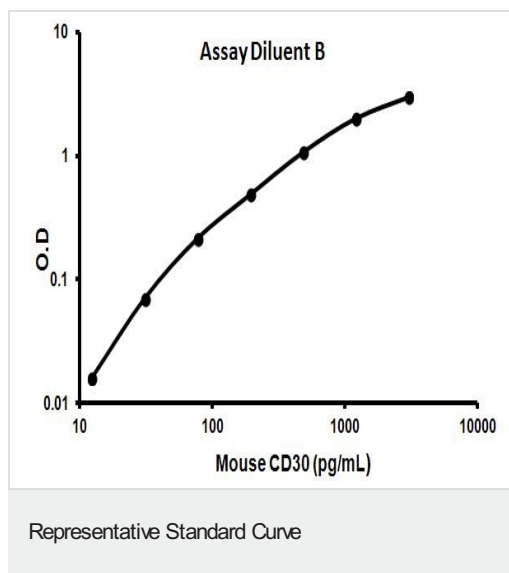
Store at -20°C. Please refer to protocols.

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
Assay Diluent A	1 x 30ml
Biotinylated anti-Mouse CD30	2 vials
CD30 Microplate (12 strips x 8 wells)	1 unit
Recombinant Mouse CD30 Standard (lyophilized)	2 vials
Stop Solution	1 x 8ml
TMB One-Step Substrate Reagent	1 x 12ml

Images



Representative Standard Curve using ab113341



Representative Standard Curve using ab113341

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