

Product datasheet

Mouse Flt 3 Ligand ELISA Kit ab113342

[2 Images](#)

Overview

Product name	Mouse Flt 3 Ligand ELISA Kit
Detection method	Colorimetric
Sample type	Cell culture supernatant, Serum, Plasma
Assay type	Sandwich (quantitative)
Sensitivity	< 15 pg/ml
Range	20 pg/ml - 5000 pg/ml
Recovery	87 %

Sample specific recovery

Sample type	Average %	Range
Cell culture supernatant	90.49	67% - 103%
Serum	91.81	83% - 99%
Plasma	77.99	69% - 85%

Assay duration Multiple steps standard assay

Species reactivity **Reacts with:** Mouse

Product overview

Abcam's Flt3 Ligand Mouse ELISA (Enzyme-Linked Immunosorbent Assay) kit is an *in vitro* enzyme-linked immunosorbent assay for the quantitative measurement of mouse Flt3 Ligand in serum, plasma and cell culture supernatants.

This assay employs an antibody specific for mouse Flt3 Ligand coated on a 96-well plate. Standards and samples are pipetted into the wells and Flt3 Ligand present in a sample is bound to the wells by the immobilized antibody. The wells are washed and biotinylated anti-mouse Flt3 Ligand 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 Flt3 Ligand bound. The Stop Solution changes the color from blue to yellow, and the intensity of the color is measured at 450 nm.

Tested applications **Suitable for:** Sandwich ELISA

Platform Microplate

Properties

Storage instructions

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

Components	1 x 96 tests
200X HRP-Streptavidin Concentrate	1 x 200µl
20X Wash Buffer Concentrate	1 x 25ml
5X Assay Diluent B	1 x 15ml
Assay Diluent A	1 x 30ml
Biotinylated anti-mouse Flt 3 Ligand	2 vials
Flt 3 Ligand Microplate (12 x 8 wells)	1 unit
Recombinant Mouse Flt 3 Ligand Standard (lyophilized)	2 vials
Stop Solution	1 x 8ml
TMB One-Step Substrate Reagent	1 x 12ml

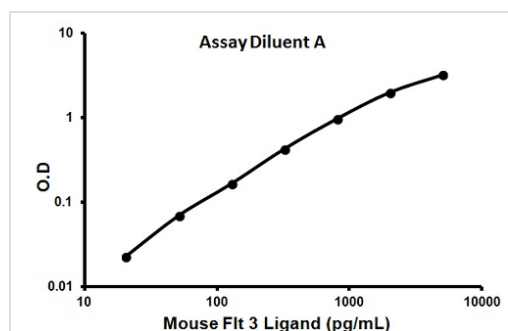
Applications

Our [Abpromise guarantee](#) covers the use of **ab113342** in the following tested applications.

The application notes include recommended starting dilutions; optimal dilutions/concentrations should be determined by the end user.

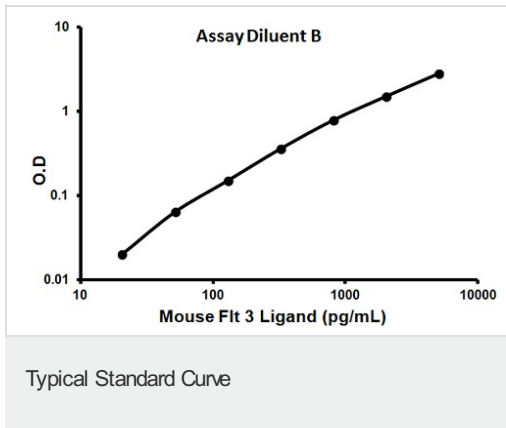
Application	Abreviews	Notes
Sandwich ELISA		Use at an assay dependent concentration.

Images



Typical Standard Curve

Representative standard curve using
ab113342



Representative standard curve using
ab113342

Please note: All products are "FOR RESEARCH USE ONLY AND ARE NOT INTENDED FOR DIAGNOSTIC OR THERAPEUTIC USE"

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