

## Product datasheet

### Mouse Flt-3 ELISA Kit ab256395

Recombinant SimpleStep ELISA®

[5 Images](#)

#### Overview

**Product name** Mouse Flt-3 ELISA Kit

**Detection method** Colorimetric

**Precision**

Intra-assay

Sample	n	Mean	SD	CV%
Extract	8			3%

Inter-assay

Sample	n	Mean	SD	CV%
Extract	7			2.9%

**Sample type** Cell Lysate

**Assay type** Sandwich (quantitative)

**Sensitivity** 0.915 ng/ml

**Range** 1.56 ng/ml - 100 ng/ml

**Recovery**

Sample specific recovery

Sample type	Average %	Range
Cell Lysate	109	97% - 119%

**Assay time** 1h 30m

**Assay duration** One step assay

**Species reactivity** **Reacts with:** Mouse

**Product overview**

Mouse Flt-3 ELISA Kit (ab256395) is a single-wash 90 min sandwich ELISA designed for the quantitative measurement of Flt-3 protein in cell lysate. It uses our proprietary SimpleStep ELISA® technology. Quantitate Mouse Flt-3 with 0.915 ng/ml sensitivity.

SimpleStep ELISA® technology employs capture antibodies conjugated to an affinity tag that is recognized by the monoclonal antibody used to coat our SimpleStep ELISA® plates. This

approach to sandwich ELISA allows the formation of the antibody-analyte sandwich complex in a single step, significantly reducing assay time. See the SimpleStep ELISA® protocol summary in the image section for further details. Our SimpleStep ELISA® technology provides several benefits:

- Single-wash protocol reduces assay time to 90 minutes or less
- High sensitivity, specificity and reproducibility from superior antibodies
- Fully validated in biological samples
- 96-wells plate breakable into 12 x 8 wells strips

A 384-well SimpleStep ELISA® microplate (**ab203359**) is available to use as an alternative to the 96-well microplate provided with SimpleStep ELISA® kits.

**Notes** Flt-3, also known as CD135, is a receptor type tyrosine kinase involved in the differentiation of hematopoietic cells. Binding of its ligand, FLT3L induces phosphorylation cascades through multiple kinases, including STAT5A and MAPK1. Flt-3 is used to differentiate multipotent progenitor cells and hematopoietic stem cells. Mutations in this gene are linked to acute myelogenous leukemia and its levels are negatively associated with survival. Given Flt-3's importance in this disease, it is currently being investigated as a therapeutic target.

Abcam has not and does not intend to apply for the REACH Authorisation of customers' uses of products that contain European Authorisation list (Annex XIV) substances.

It is the responsibility of our customers to check the necessity of application of REACH Authorisation, and any other relevant authorisations, for their intended uses.

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

## Properties

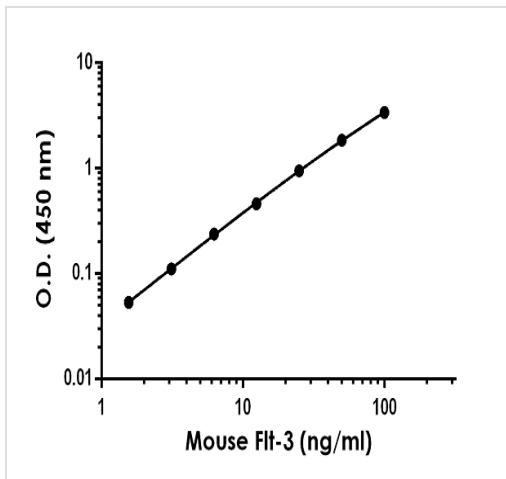
**Storage instructions** Store at +4°C. Please refer to protocols.

Components	1 x 96 tests
10X Mouse Flt-3 Capture Antibody	1 x 600µl
10X Mouse Flt-3 Detector Antibody	2 vials
10X Wash Buffer PT (ab206977)	1 x 20ml
50X Cell Extraction Enhancer Solution (ab193971)	1 x 1ml
5X Cell Extraction Buffer PTR (ab193970)	1 x 10ml
Antibody Diluent CPR2	1 x 6ml
Mouse Flt-3 Lyophilized Recombinant Protein	2 vials
Plate Seals	1 unit
Sample Diluent NS (ab193972)	1 x 12ml
SimpleStep Pre-Coated 96-Well Microplate (ab206978)	1 unit

Components	1 x 96 tests
Stop Solution	1 x 12ml
TMB Development Solution	1 x 12ml

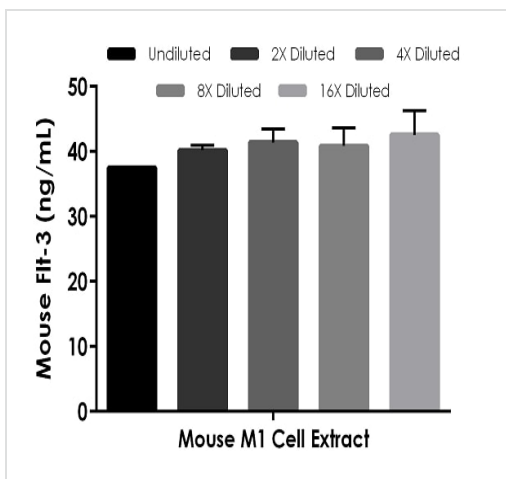
**Sequence similarities**      Belongs to the IFIT family.  
 Contains 8 TPR repeats.

**Images**



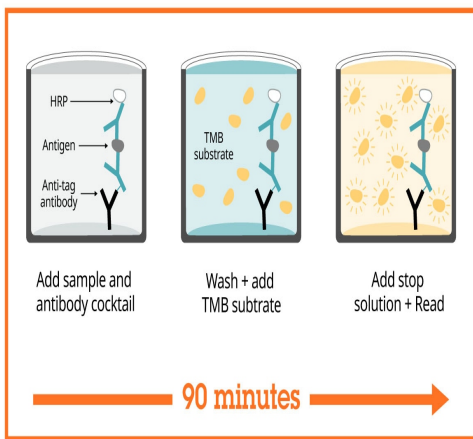
Example of mouse Flt-3 standard curve in 1X Cell Extraction Buffer PTR.

The Flt-3 standard curve was prepared as described in Section 10. Raw data values are shown in the table. Background-subtracted data values (mean +/- SD) are graphed.



Interpolated concentrations of native Flt-3 in mouse M1 cell extract based on a 200 µg/mL extract load.

The concentrations of Flt-3 were measured in duplicate and interpolated from the Flt-3 standard curve and corrected for sample dilution. The interpolated, dilution factor corrected, values are plotted (mean +/- SD, n=2). The mean Flt-3 concentration was determined to be 40.5 ng/mL.



Sandwich ELISA - Mouse Fit-3 ELISA Kit  
(ab256395)

SimpleStep ELISA technology allows the formation of the antibody-antigen complex in one single step, reducing assay time to 90 minutes. Add samples or standards and antibody mix to wells all at once, incubate, wash, and add your final substrate. See protocol for a detailed step-by-step guide.

Powered by  
recombinant antibodies

- Research with confidence**  
Consistent and reproducible results
- Long-term and scalable supply**  
Recombinant technology
- Success from the first experiment**  
Confirmed specificity
- Ethical standards compliant**  
Animal-free production

Sandwich ELISA - Mouse Fit-3 ELISA Kit  
(ab256395)

To learn more about the advantages of recombinant antibodies see [here](#).

### Get more done with SimpleStep ELISA



**Easy to use**  
Single-wash 90-minute protocol



**Flexible**  
Matched antibody pairs available



**Precision antibodies**  
High sensitivity, specificity and reproducibility



**Scalable**  
Now in 10-pack and 384-well formats

Sandwich ELISA - Mouse Fit-3 ELISA Kit  
(ab256395)

To learn more about the advantages of SimpleStep ELISA<sup>®</sup> kits see [here](#).

**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