

Human c-Kit ELISA Kit (CD117) ab45924

1 References 1 Image

Overview

Product name Human c-Kit ELISA Kit (CD117)

Detection method Colorimetric

Precision	Intra-assay				
	Sample	n	Mean	SD	CV%
	Sample A	8	5.629ng/ml	0.34	6%
	Sample B	8	1.231ng/ml	0.06	4.8%

	Inter-assay				
	Sample	n	Mean	SD	CV%
	Sample A	8	5.44ng/ml	0.43	7.9%
	Sample B	8	1.314ng/ml	0.1	7.6%

Sample type Cell culture supernatant, Serum, Plasma

Assay type Sandwich (quantitative)

Sensitivity < 0.095 ng/ml

Range 0.312 ng/ml - 10 ng/ml

Assay time 1h 45m

Assay duration Multiple steps standard assay

Species reactivity **Reacts with:** Human

Product overview Abcam's c-Kit Human *in vitro* ELISA (Enzyme-Linked Immunosorbent Assay) kit is designed for the quantitative measurement of c-Kit in Human serum, plasma, buffered solutions or supernatants.

A monoclonal antibody specific for c-Kit has been coated onto the wells of the microtiter strips provided. Samples, including standards of known c-Kit concentrations, control specimens or unknowns are pipetted into these wells. During the first incubation, the standards or samples and a biotinylated monoclonal antibody specific for c-Kit are simultaneously incubated. After washing, the enzyme Streptavidin-HRP, that binds the biotinylated antibody is added, incubated and

washed. A TMB substrate solution is added which acts on the bound enzyme to induce a colored reaction product. The intensity of this colored product is directly proportional to the concentration of c-Kit present in the samples.

This kit will recognize both endogenous and recombinant Human c-Kit.

Notes The standard is affinity purified.

Platform Microplate

Properties

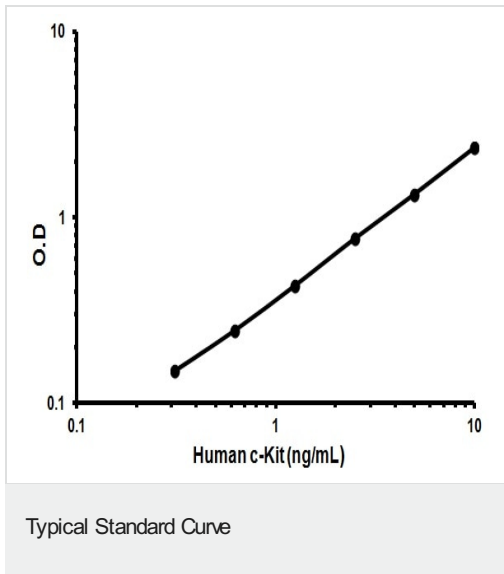
Storage instructions Store at +4°C. Please refer to protocols.

Components	Identifier	1 x 96 tests	2 x 96 tests
10X Standard Diluent Buffer	Black	1 x 15ml	1 x 25ml
200X Wash Buffer	White	1 x 10ml	2 x 10ml
Biotinylated anti c-Kit		1 x 0.4ml	2 x 0.4ml
Biotinylated Antibody Diluent	Red	1 x 7.5ml	1 x 13ml
Chromogen TMB Substrate Solution		1 x 11ml	1 x 24ml
c-Kit Control	Silver	2 vials	4 vials
c-Kit Microplate (12 x 8 well strips)		1 unit	2 units
c-Kit Standard (Lyophilized)	Yellow	2 vials	4 vials
HRP Diluent	Red	1 x 12ml	1 x 23ml
Plastic Plate Covers		1 x 2 units	1 x 2 units
Stop Reagent	Black	1 x 11ml	2 x 11ml
Streptavidin-HRP		2 x 5µl	4 x 5µl

Relevance KIT encodes the human homolog of the proto-oncogene c-kit. c-kit was first identified as the cellular homolog of the feline sarcoma viral oncogene v-kit. KIT is a type 3 transmembrane receptor for MGF (mast cell growth factor, also known as stem cell factor). Mutations in KIT are associated with gastrointestinal stromal tumors, mast cell disease, acute myelogenous leukemia, and piebaldism.

Cellular localization Cell Membrane; single-pass type I membrane protein

Images



Representative Standard Curve using ab45924.

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