

Human c-Kit ELISA Kit (CD117) ab45924

1 References 画像数 1

製品の概要

製品名	Human c-Kit ELISA Kit (CD117)				
検出方法	Colorimetric				
再現性	Intra-Assay (同時再現性)				
	サンプル	N	平均値	SD	CV%
	Sample A	8	5.629ng/ml	0.34	6%
	Sample B	8	1.231ng/ml	0.06	4.8%
	Inter-Assay (日差再現性)				
	サンプル	N	平均値	SD	CV%
	Sample A	8	5.44ng/ml	0.43	7.9%
	Sample B	8	1.314ng/ml	0.1	7.6%
サンプルの種類	Cell culture supernatant, Serum, Plasma				
アッセイタイプ	Sandwich (quantitative)				
検出感度	< 0.095 ng/ml				
検出範囲	0.312 ng/ml - 10 ng/ml				
全工程の試験時間	1h 45m				
ステップ	Multiple steps standard assay				
種交差性	交差種: Human				
製品の概要	Abcam's c-Kit Human <i>in vitro</i> ELISA (Enzyme-Linked Immunosorbent Assay) kit is designed for				

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.

#### 特記事項

The standard is affinity purified.

#### 試験プラットフォーム

Microplate

#### 製品の特性

#### 保存方法

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

内容	ラベル	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

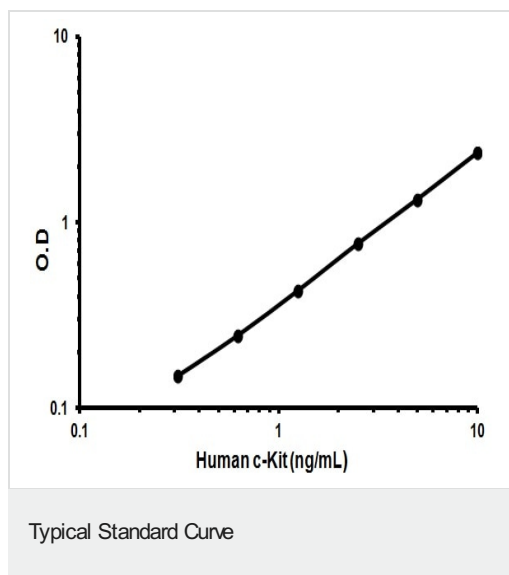
#### 関連性

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.

#### 細胞内局在

Cell Membrane; single-pass type I membrane protein

#### 画像



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.co.jp/abpromise> or contact our technical team.

#### Terms and conditions

---

- Guarantee only valid for products bought direct from Abcam or one of our authorized distributors