

# Human TREM1 ELISA Kit ab100659

[1 References](#) [2 图像](#)

### 概述

产品名称	人TREM1 ELISA试剂盒
检测方法	Colorimetric
样品类型	Cell culture supernatant, Serum, Plasma
检测类型	Sandwich (quantitative)
灵敏度	< 70 pg/ml
范围	0.069 ng/ml - 50 ng/ml
回收率	> 100 %

特定样本回收率

样品类型	平均%	范围
Cell culture supernatant	124.4	106% - 134%
Serum	125.5	91% - 145%
Plasma	116	94% - 115%

**实验步骤** Multiple steps standard assay

**种属反应性** 与反应: Human

**产品概述** Abcam's TREM1 Human ELISA (Enzyme-Linked Immunosorbent Assay) kit is an *in vitro* enzyme-linked immunosorbent assay for the quantitative measurement of Human TREM1 in serum, plasma and cell culture supernatants. (Human TREM1 concentration is pretty low in normal serum, it may not be detected in this assay).

This assay employs an antibody specific for Human TREM1 coated on a 96-well plate. Standards and samples are pipetted into the wells and TREM1 present in a sample is bound to the wells by the immobilized antibody. The wells are washed and biotinylated anti-Human TREM1 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 TREM1 bound. The Stop Solution changes the color from blue to yellow, and the intensity of the color is measured at 450 nm.

Get higher sensitivity in only 90 minutes with Human soluble TREM1 ELISA Kit ([ab185987](#)) from

our SimpleStep ELISA® range.

**说明**

Optimization may be required with urine samples.

**平台**

Microplate

**性能**

---

**存放说明**

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

组件	1 x 96 tests
20X Wash Buffer	1 x 25ml
5X Assay Diluent B	1 x 15ml
500X HRP-Streptavidin Concentrate	1 x 200µl
Assay Diluent A	1 x 30ml
Biotinylated anti-Human TREM1	2 vials
Recombinant Human TREM1 Standard (lyophilized)	2 vials
Stop Solution	1 x 8ml
TMB One-Step Substrate Reagent	1 x 12ml
TREM1 Microplate (12 x 8 wells)	1 unit

**相关性**

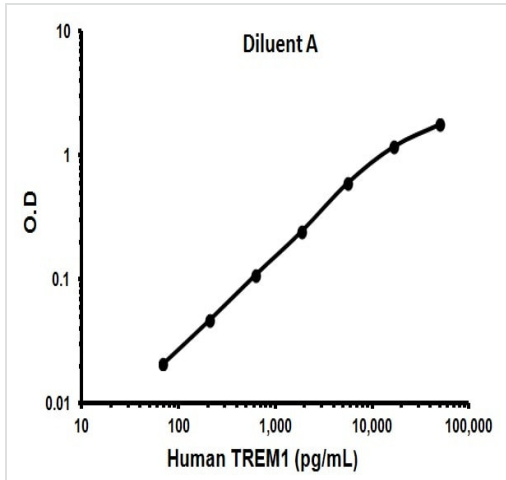
TREM1 is also known as triggering receptor expressed on myeloid cells 1, and triggering receptor expressed on monocytes 1. TREM1 is a Type I membrane protein that contains an immunoglobulin-like V-type domain. An alternatively spliced protein variant may be secreted. TREM1 is highly expressed on peripheral blood myeloid cells (particularly mature monocytes and granulocytes); TREM1 expression can be further upregulated by bacteria, fungi and lipopolysaccharide. TREM1 has been shown to interact with the adaptor protein DAP12 to stimulate neutrophil and monocyte-mediated inflammatory responses through the triggering and release of pro-inflammatory cytokines and chemokines. TREM1 is thought to amplify inflammatory responses to fungal and bacterial infections and potentiate septic shock.

**细胞定位**

Secreted and Cell membrane.

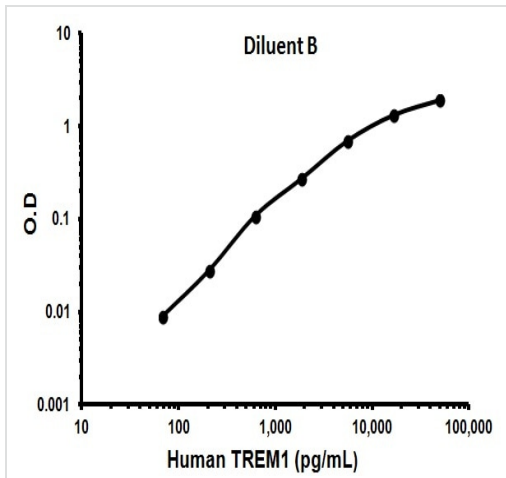
**图片**

---



Representative standard curve using ab100659

Typical standard curve



Representative standard curve using ab100659

Typical standard curve

**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.cn/abpromise> or contact our technical team.

### **Terms and conditions**

---

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