

Mouse Thrombopoietin ELISA Kit ab100748

[2 References](#) [2 图像](#)

概述

产品名称	小鼠Thrombopoietin 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	< 30 pg/ml
范围	27.4 pg/ml - 20000 pg/ml
回收率	94 %

特定样本回收率

样品类型	平均%	范围
Cell culture supernatant	96.23	83% - 103%
Serum	93.45	82% - 103%
Plasma	94.28	84% - 104%

实验步骤 Multiple steps standard assay

种属反应性 与反应: Mouse

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

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

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
300X HRP-Streptavidin Concentrate	1 x 200µl
5X Assay Diluent B	1 x 15ml
Assay Diluent A	1 x 30ml
Biotinylated anti-Mouse Thrombopoietin	2 vials
Recombinant Mouse Thrombopoietin Standard (lyophilized)	2 vials
Stop Solution	1 x 8ml
Thrombopoietin Microplate (12 x 8 wells)	1 unit
TMB One-Step Substrate Reagent	1 x 12ml

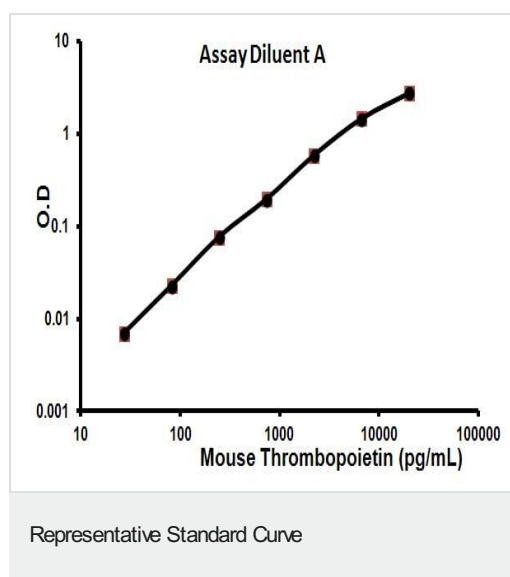
相关性

Thrombopoietin is a humoral growth factor that is necessary for megakaryocyte proliferation and maturation, as well as for thrombopoiesis. This protein is the ligand for MLP/cMPL, the product of myeloproliferative leukemia virus oncogene. Alternatively spliced transcript variants encoding distinct isoforms have been reported.

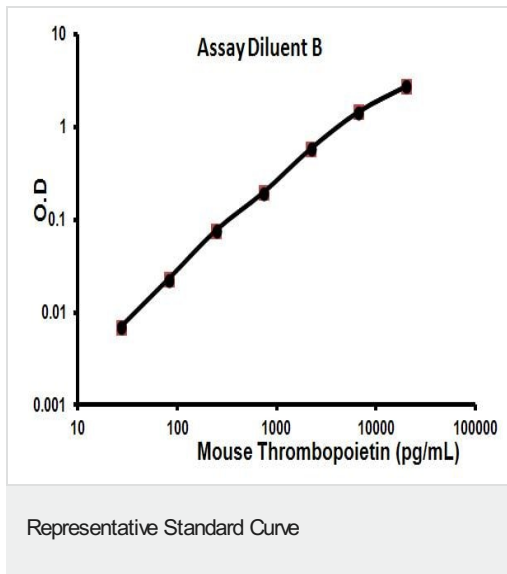
细胞定位

Secreted

图片



Representative Standard Curve using ab100748



Representative Standard Curve using ab100748

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