

Mouse IL-28 ELISA Kit ab100708

1 References 画像数 1

製品の概要

製品名	Mouse IL-28 ELISA Kit
検出方法	Colorimetric
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	< 2 pg/ml
検出範囲	2.46 pg/ml - 600 pg/ml
添加回収試験	88 %

特定サンプルでの回収試験

サンプルの種類	平均 %	測定範囲
Cell culture supernatant	74.3	67% - 87%
Serum	109.8	98% - 118%
Plasma	76.8	67% - 93%

ステップ Multiple steps standard assay

種交差性 交差種: Mouse

製品の概要

Abcam's IL-28 Mouse ELISA (Enzyme-Linked Immunosorbent Assay) kit is an *in vitro* enzyme-linked immunosorbent assay for the quantitative measurement of Mouse IL-28 in serum, plasma and cell culture supernatants.

This assay employs an antibody specific for IL-28 coated on a 96-well plate. Standards and samples are pipetted into the wells and IL-28 present in a sample is bound to the wells by the immobilized antibody. The wells are washed and biotinylated anti-Mouse IL-28 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 IL-28 bound. The Stop Solution changes the color from blue to yellow, and the intensity of the color is measured at 450 nm.

試験プラットフォーム Microplate

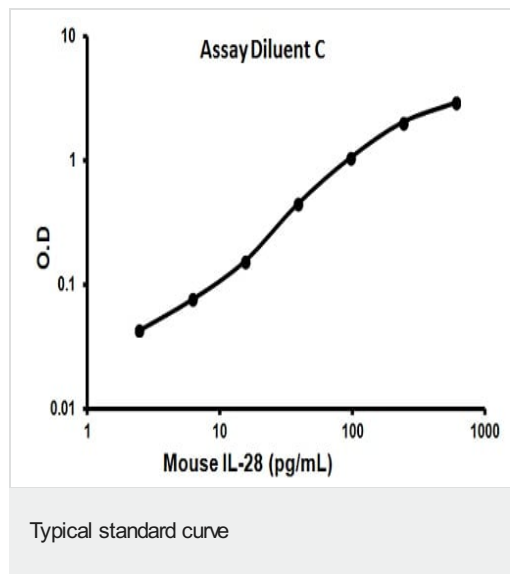
製品の特性

保存方法

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

内容	1 x 96 tests
200X HRP-Streptavidin Concentrate	1 x 200µl
20X Wash Buffer	1 x 25ml
5X Assay Diluent B	1 x 15ml
Assay Diluent C	1 x 30ml
Biotinylated anti-Mouse IL-28	2 vials
IL-28 Microplate (12 x 8 wells)	1 unit
Recombinant Mouse IL-28 Standard (lyophilized)	2 vials
Stop Solution	1 x 8ml
TMB One-Step Substrate Reagent	1 x 12ml

画像



Representative standard curve using ab100708

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