

Mouse IL-12 p40 + IL-12 p70 ELISA Kit ab100699

3 References 画像数 2

製品の概要

製品名	Mouse IL-12 p40 + IL-12 p70 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	< 5 pg/ml
検出範囲	2.74 pg/ml - 2000 pg/ml
添加回収試験	94 %

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	95.25	84% - 103%
Serum	93.29	82% - 102%
Plasma	94.52	83% - 103%

ステップ Multiple steps standard assay

種交差性 交差種: Mouse

製品の概要 Abcam's IL-12 p40 + IL-12 p70 Mouse ELISA (Enzyme-Linked Immunosorbent Assay) kit is an *in vitro* enzyme-linked immunosorbent assay for the quantitative measurement of Mouse IL-12 p40 + IL-12 p70 in serum, plasma and cell culture supernatants.

This assay employs an antibody specific for Mouse IL-12 p40 + IL-12 p70 coated on a 96-well plate. Standards and samples are pipetted into the wells and IL-12 p40 + IL-12 p70 present in a sample is bound to the wells by the immobilized antibody. The wells are washed and biotinylated anti-IL-12 p40 + IL-12 p70 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 bound. The Stop Solution changes the color from blue to yellow, and the intensity of the color is measured at 450 nm.

特記事項 The standard is affinity purified.

試験プラットフォーム Microplate

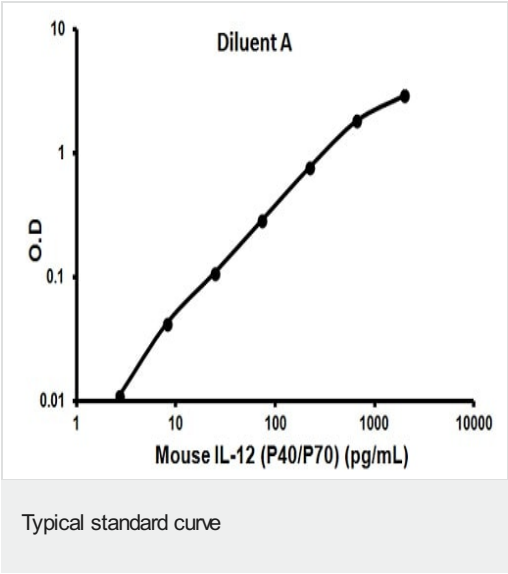
製品の特性

保存方法Store at -20°C. Please refer to protocols.

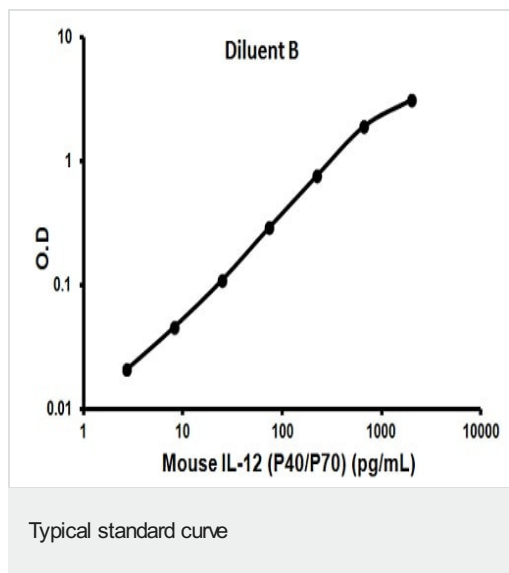
内容	1 x 96 tests
20X Wash Buffer	1 x 25ml
400X HRP-Streptavidin Concentrate	1 x 200µl
5X Assay Diluent B	1 x 15ml
Assay Diluent A	1 x 30ml
Biotinylated anti-Mouse IL-12 p40/p70	2 vials
IL-12 p40/p70 Microplate (12 x 8 wells)	1 unit
Recombinant Mouse IL-12 p40/p70 Standard (lyophilized)	2 vials
Stop Solution	1 x 8ml
TMB One-Step Substrate Reagent	1 x 12ml

関連性IL-12 is a heterodimeric cytokine comprised of p35 and p40 subunits. IL-12 plays a central role in mediating cell mediated immunity, promoting differentiation of CD4+ T cells to the Th1 subset and of CD8+ T cells into mature cytotoxic T lymphocytes (CTLs). IL-12 is a potent stimulator of NK cells. IL-12 therapy has been suggested as a method of enhancing cytotoxic anti tumour immune responses.

画像



Representative standard curve using ab100699



Representative standard curve using ab100699

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