

Human IL-17A Receptor ELISA Kit ab100558

画像数 2

製品の概要

製品名	Human IL-17A Receptor 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	< 10 pg/ml
検出範囲	12.35 pg/ml - 3000 pg/ml
添加回収試験	81 %

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	81.54	70% - 92%
Serum	87.64	83% - 104%
Plasma	73.27	67% - 87%

ステップ Multiple steps standard assay

種交差性 交差種: Human

製品の概要 Abcam's IL-17A Receptor Human ELISA (Enzyme Linked Immunosorbent Assay) kit is an *in vitro* enzyme-linked immunosorbent assay for the quantitative measurement of Human IL-17A in serum, plasma, and cell culture supernatants.

This assay employs an antibody specific for Human IL-17A coated on a 96-well plate. Standards and samples are pipetted into the wells and IL-17A present in a sample is bound to the wells by the immobilized antibody. The wells are washed and biotinylated anti-Human IL-17A 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-17A 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 Optimization may be required with urine samples.

試験プラットフォーム Microplate

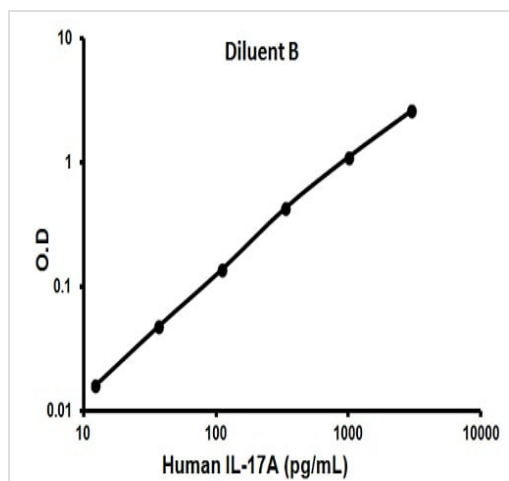
製品の特性

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-Human IL-17A	2 vials
IL-17A Microplate (12 x 8 wells)	1 x 200µl
Recombinant Human IL-17A Standard (lyophilized)	2 vials
Stop Solution	1 x 8ml
TMB One-Step Substrate Reagent	1 x 12ml

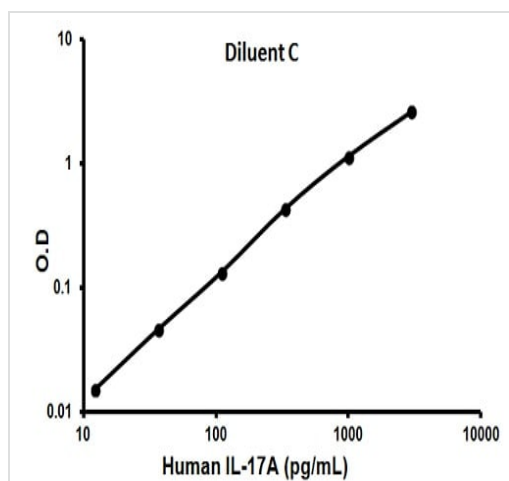
機能	Receptor for IL17A and IL17F. Binds its IL17A ligand with low affinity, suggesting that additional components are involved in IL17A-induced signaling.
組織特異性	Widely expressed.
関連疾患	Defects in IL17RA are the cause of familial candidiasis type 5 (CANDF5) [MIM:613953]. CANDF5 is a rare disorder with altered immune responses and impaired clearance of fungal infections, selective against Candida. It is characterized by persistent and/or recurrent infections of the skin, nails and mucous membranes caused by organisms of the genus Candida, mainly Candida albicans.
配列類似性	Contains 1 SEFIR domain.
翻訳後修飾	Glycosylated.
細胞内局在	Membrane.

画像



Representative standard curve using ab100558

Typical standard curve



Representative standard curve using ab100558

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.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