

Human IL-1 R4 ELISA Kit (ST2) ab100563

2 References 2 图像

概述

产品名称	人IL-1 R4 ELISA试剂盒(ST2)
检测方法	Colorimetric
样品类型	Cell culture supernatant, Serum, Plasma
检测类型	Sandwich (quantitative)
灵敏度	< 2 pg/ml
范围	1.65 pg/ml - 1200 pg/ml
回收率	97 %

特定样本回收率

样品类型	平均%	范围
Cell culture supernatant	98.43	91% - 105%
Serum	109.6	98% - 117%
Plasma	84.06	67% - 102%

检测时间	5h 00m
实验步骤	Multiple steps standard assay
种属反应性	与反应: Human
产品概述	Human IL-1 R4 ELISA Kit (ST2) (ab100563) is an <i>in vitro</i> enzyme-linked immunosorbent assay for the quantitative measurement of Human IL-1 R4 in serum, plasma, and cell culture supernatants.

This assay employs an antibody specific for Human IL-1 R4 coated on a 96-well plate. Standards and samples are pipetted into the wells and IL-1 R4 present in a sample is bound to the wells by the immobilized antibody. The wells are washed and biotinylated anti-Human IL-1 R4 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-1 R4 bound. The Stop Solution changes the color from blue to yellow, and the intensity of the color is measured at 450 nm.

说明	Optimization may be required with urine samples.
----	--

平台

Microplate

性能

存放说明

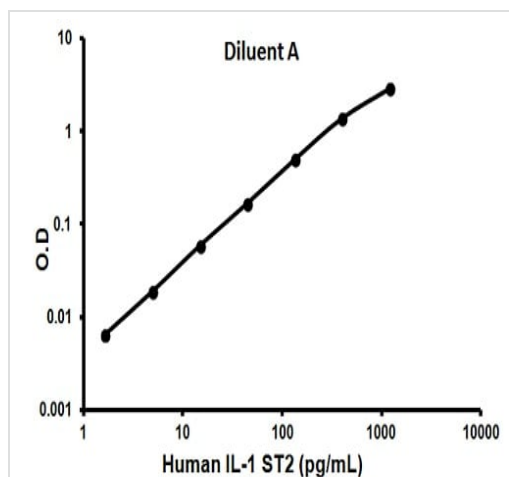
Store at -20°C. Please refer to protocols.

组 件	1 x 96 tests
20X Wash Buffer	1 x 25ml
300X HRP-Streptavidin Concentrate	1 x 200µl
5X Assay Diluent B	1 x 15ml
Assay Diluent A	1 x 30ml
Biotinylated anti-Human IL-1 R4	2 vials
IL-1 R4 Microplate (12 x 8 wells)	1 unit
Recombinant Human IL-1 R4 Standard (lyophilized)	2 vials
Stop Solution	1 x 8ml
TMB One-Step Substrate Reagent	1 x 12ml

相关性

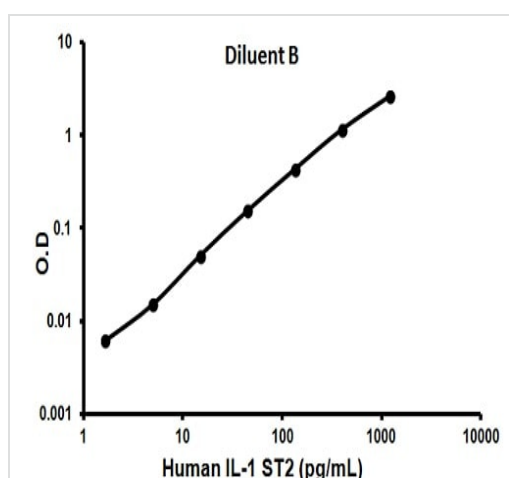
ST2 is a member of a superfamily containing the Interleukin 1 (IL1) receptor and the Toll-like receptors (TLRs). The TLRs are signaling molecules that recognize different microbial products during infection and serve as an important link between the innate and adaptive immune responses. ST2 was originally identified as a protein whose production was stimulated by various proliferation-inducing agents such as PDGF and FGF. More recently, it has been shown to negatively regulate IL1 receptor and Toll-like receptor 4 (TLR4) signaling and to maintain endotoxin tolerance. It has been suggested that the inhibition of TLR4 signaling occurs through the association and sequestering of TLR adaptor molecules such as MyD88 and TIRAP.

图片



Representative standard curve using ab100563

Typical standard curve



Representative standard curve using ab100563

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