

Human SAA ELISA Kit (SAA1) ab100635

13 References 画像数 1

製品の概要

製品名	Human SAA ELISA Kit (SAA1)
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	< 0.5 ng/ml
検出範囲	0.41 ng/ml - 300 ng/ml
添加回収試験	100 %

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	93.75	86% - 102%
Serum	114.8	105% - 124%
Plasma	112.1	83% - 137%

ステップ	Multiple steps standard assay
種交差性	交差種: Human
製品の概要	Human SAA ELISA (Enzyme-Linked Immunosorbent Assay) kit is an <i>in vitro</i> enzyme-linked immunosorbent assay for the quantitative measurement of human SAA in serum, plasma and cell culture supernatants.

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

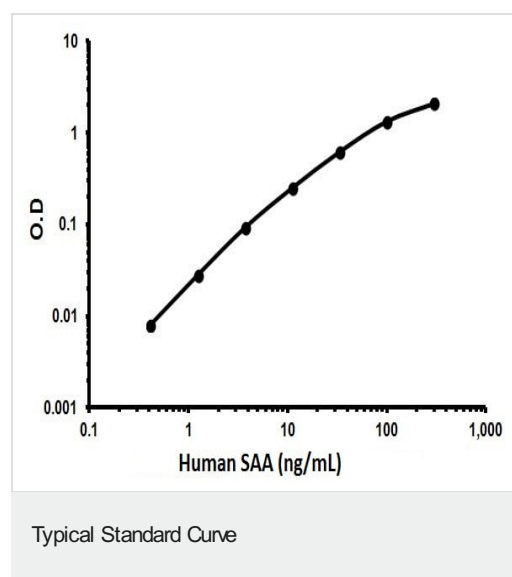
特記事項	Optimization may be required with urine samples. ab100635 cross reacts with SAA2.
試験プラットフォーム	Microplate

## 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
5X Assay Diluent B	1 x 15ml
800X HRP-Streptavidin Concentrate	1 x 200µl
Assay Diluent C	1 x 30ml
Biotinylated anti-Human SAA	2 vials
Recombinant Human SAA Standard (lyophilized)	2 vials
Serum Amyloid A Microplate (12 x 8 wells)	1 unit
Stop Solution	1 x 8ml
TMB One-Step Substrate Reagent	1 x 12ml

## 画像



Representative standard curve using ab100635.

**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