

Human Alpha 1 Acid Glycoprotein / AGP ELISA Kit (AGP) ab108852

4 References [画像数 1](#)

製品の概要

製品名 Human Alpha 1 Acid Glycoprotein / AGP ELISA Kit (AGP)

検出方法 Colorimetric

再現性 Intra-Assay (同時再現性)

サンプル	N	平均値	SD	CV%
Overall				5.1%

Inter-Assay (日差再現性)

サンプル	N	平均値	SD	CV%
Overall				9.9%

サンプルの種類 Cell culture supernatant, Saliva, Milk, Urine, Serum, Plasma, Cerebral Spinal Fluid

アッセイタイプ Sandwich (quantitative)

検出感度 = 1.6 ng/ml

検出範囲 3.125 ng/ml - 200 ng/ml

添加回収試験 = 96 %

全工程の試験時間 4h 00m

ステップ Multiple steps standard assay

種交差性 交差種: Human

製品の概要 Abcam's Alpha 1 Acid Glycoprotein / AGP Human *in vitro* ELISA (Enzyme-Linked Immunosorbent Assay) kit is designed for the quantitative measurement of AGP levels in cell culture supernatant, saliva, milk, and urine.

An Alpha 1 Acid Glycoprotein / AGP specific antibody has been precoated onto 96-well plates and blocked. Standards or test samples are added to the wells and subsequently an Alpha 1 Acid Glycoprotein / AGP specific biotinylated detection antibody is added and then followed by washing with wash buffer. Streptavidin-Peroxidase Conjugate is added and unbound conjugates are washed away with wash buffer. TMB is then used to visualize Streptavidin-Peroxidase enzymatic reaction. TMB is catalyzed by Streptavidin-Peroxidase to produce a blue color product

that changes into yellow after adding acidic stop solution. The density of yellow coloration is directly proportional to the amount of Alpha 1 Acid Glycoprotein / AGP captured in plate.

The entire kit may be stored at -20°C for long term storage before reconstitution - Avoid repeated freeze-thaw cycles.

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
100X Streptavidin-Peroxidase Conjugate	1 x 80µl
10X Diluent N Concentrate	1 x 30ml
20X Wash Buffer Concentrate	2 x 30ml
50X Biotinylated Human Alpha 1 Acid Glycoprotein / AGP Antibody	1 x 120µl
Alpha 1 Acid Glycoprotein / AGP Microplate (12 x 8 well strips)	1 unit
Alpha 1 Acid Glycoprotein / AGP Standard	1 vial
Chromogen Substrate	1 x 8ml
Sealing Tapes	3 units
Stop Solution	1 x 12ml

機能

Appears to function in modulating the activity of the immune system during the acute-phase reaction.

組織特異性

Expressed by the liver and secreted in plasma.

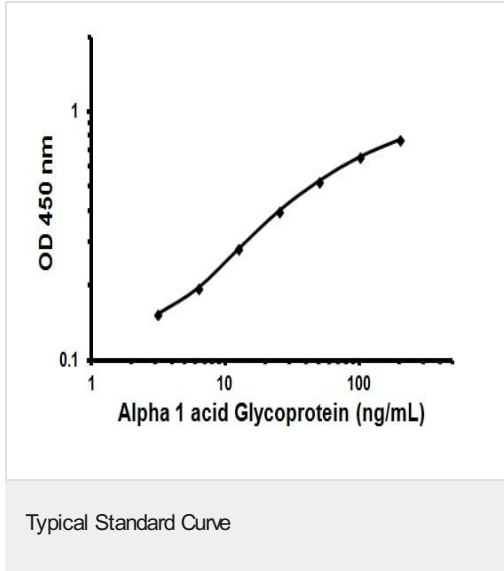
配列類似性

Belongs to the calycin superfamily. Lipocalin family.

細胞内局在

Secreted.

画像



Representative Standard Curve using ab108852.

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