

Mouse Alpha 1 Antitrypsin ELISA Kit ab205088

4 References 画像数 1

医薬用外毒物

医薬用外劇物

製品の概要

製品名 Mouse Alpha 1 Antitrypsin ELISA Kit

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%
	Serum				< 10%
	Inter-Assay (日差再現性)				
	サンプル	N	平均値	SD	CV%
	Serum				< 10%

サンプルの種類 Urine, Serum, Hep Plasma, EDTA Plasma, Cit plasma

アッセイタイプ Sandwich (quantitative)

検出感度 1.492 ng/ml

検出範囲 6.25 ng/ml - 200 ng/ml

添加回収試験 > 85 %

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
Serum	> 85	6.25ng/ml - 200ng/ml

全工程の試験時間 2h 10m

ステップ Multiple steps standard assay

種交差性 交差種: Mouse

製品の概要 Alpha 1 Antitrypsin *in vitro* ELISA (Enzyme-Linked Immunosorbent Assay) kit is designed for the quantitative measurement of Alpha 1 Antitrypsin protein in mouse biological samples.

特記事項 In this assay the A1AT present in samples reacts with the anti-A1AT antibodies which have been adsorbed to the surface of polystyrene microtitre wells. After the removal of unbound proteins by washing, anti-A1AT antibodies conjugated with horseradish peroxidase (HRP), are added. These

enzyme-labeled antibodies form complexes with the previously bound A1AT . Following another washing step, the enzyme bound to the immunosorbent is assayed by the addition of a chromogenic substrate, 3,3',5,5'-tetramethylbenzidine (TMB). The quantity of bound enzyme varies directly with the concentration of A1AT in the sample tested; thus, the absorbance, at 450 nm, is a measure of the concentration of A1AT in the test sample. The quantity of A1AT in the test sample can be interpolated from the standard curve constructed from the standards, and corrected for sample dilution.

試験プラットフォーム

Pre-coated microplate (12 x 8 well strips)

製品の特性

保存方法

Store at +4°C. Please refer to protocols.

内容	1 x 96 tests
100X HRP Conjugated Enzyme Antibody	1 x 150µl
20X Wash Buffer Concentrate	1 x 50ml
5X Diluent Concentrate	1 x 50ml
Anti-Mouse A1AT ELISA Microplate	1 unit
Chromogen Substrate Solution	1 x 12ml
Mouse A1AT Calibrator (lyophilized)	1 vial
Stop Solution	1 x 12ml

機能

Inhibitor of serine proteases. Its primary target is elastase, but it also has a moderate affinity for plasmin and thrombin. Irreversibly inhibits trypsin, chymotrypsin and plasminogen activator. The aberrant form inhibits insulin-induced NO synthesis in platelets, decreases coagulation time and has proteolytic activity against insulin and plasmin.
Short peptide from AAT: reversible chymotrypsin inhibitor. It also inhibits elastase, but not trypsin. Its major physiological function is the protection of the lower respiratory tract against proteolytic destruction by human leukocyte elastase (HLE).

組織特異性

Ubiquitous. Expressed in leukocytes and plasma.

関連疾患

Alpha-1-antitrypsin deficiency

配列類似性

Belongs to the serpin family.

ドメイン

The reactive center loop (RCL) extends out from the body of the protein and directs binding to the target protease. The protease cleaves the serpin at the reactive site within the RCL, establishing a covalent linkage between the carboxyl group of the serpin reactive site and the serine hydroxyl of the protease. The resulting inactive serpin-protease complex is highly stable.

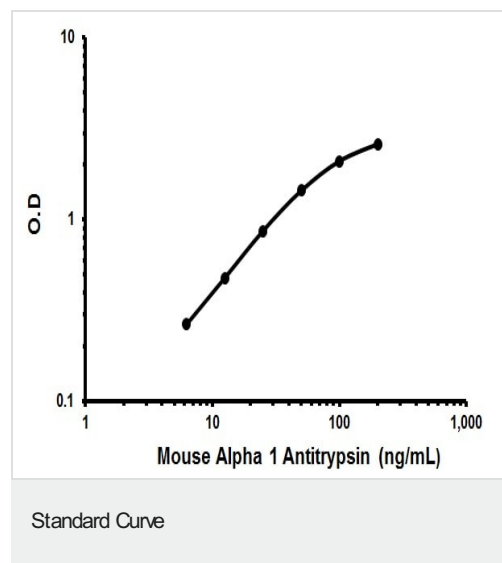
翻訳後修飾

N-glycosylated. Differential glycosylation produces a number of isoforms. N-linked glycan at Asn-107 is alternatively di-antennary, tri-antennary or tetra-antennary. The glycan at Asn-70 is di-antennary with trace amounts of tri-antennary. Glycan at Asn-271 is exclusively di-antennary. Structure of glycans at Asn-70 and Asn-271 is Hex5HexNAc4. The structure of the antennae is Neu5Ac(alpha1-6)Gal(beta1-4)GlcNAc attached to the core structure Man(alpha1-6)[Man(alpha1-3)]Man(beta1-4)GlcNAc(beta1-4)GlcNAc. Some antennae are fucosylated, which forms a Lewis-X determinant.
Proteolytic processing may yield the truncated form that ranges from Asp-30 to Lys-418.

細胞内局在

Secreted. Endoplasmic reticulum. The S and Z allele are not secreted effectively and accumulate intracellularly in the endoplasmic reticulum and Secreted, extracellular space, extracellular matrix.

画像



Representative standard curve using ab205088 Mouse Alpha 1 Antitrypsin ELISA Kit.

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