

Human Complement C3a des Arg ELISA Kit ab133037

5 References 画像数 1

製品の概要

製品名	Human Complement C3a des Arg 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%
	Low conc.	16	1.86ng/ml		8.7%
	Medium conc.	16	3.4ng/ml		9.8%
	High conc.	16	12.3ng/ml		11.1%
	Inter-Assay (日差再現性)				
	サンプル	N	平均値	SD	CV%
	Low		0.753ng/ml		5.7%
	Medium		1.529ng/ml		16.9%
	High		3.491ng/ml		28.6%
サンプルの種類	Plasma				
アッセイタイプ	Competitive				
検出感度	0.12 ng/ml				
検出範囲	0.313 ng/ml - 20 ng/ml				
添加回収試験	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	
	Plasma		94.8	% - %	
全工程の試験時間	3h 00m				
ステップ	Multiple steps standard assay				
種交差性	交差種: Human				

製品の概要

Abcam's Complement C3a des Arg Human *in vitro* competitive ELISA (Enzyme-Linked Immunosorbent Assay) kit is for the accurate quantitation of Human Complement C3a des Arg in plasma.

A goat anti-rabbit IgG antibody has been precoated onto 96-well plates. Standards or test samples are added to the wells, along with an alkaline phosphatase (AP) conjugated- Complement C3a des Arg antigen and a polyclonal rabbit antibody specific to Complement C3a des Arg. After incubation the excess reagents are washed away. pNpp substrate is added and after a short incubation the enzyme reaction is stopped and the yellow color generated is read at 405 nm. The intensity of the yellow coloration is inversely proportional to the amount of Complement C3a des Arg captured in the plate.

特記事項

The Human C3a des Arg molecule is one of three activation fragments formed from the activation of the complement cascade. C3a des Arg is formed from C3a via carboxypeptidase cleavage of the C-terminal arginine group. The structure of Human C3a des Arg was first reported in 1975 by Hugli. Human C3a des Arg contains 77 amino acids with 6 cysteines involved in disulfide bridges between residues 22-49, 23-56 and 36-57. The C terminal end of C3a des Arg in Human, porcine, rat, mouse and guinea pig is identical. C3a des Arg is a highly cationic molecule containing no carbohydrate. X-ray crystal data shows the N and C terminal residues to have highly flexible helical structures. C3a is one of the most potent constrictors of smooth muscle cells, and guinea pig airways are hyperresponsive to C3a when pretreated with histamine. The long term study of liver and other transplant recipients for both C3a des Arg and C4a des Arg may be useful in assessing a number of pathological conditions. The use of potent protease inhibitors, such as Futhan, in conjunction with EDTA, may allow complement activation factors to be quantitated specifically via inhibition of non-specific protease formation of C3a des Arg.

Cross Reactivity

Compound	Cross Reactivity
Human Complement C3a des Arg	100%
Human Complement C3	1.28%
Human Complement C4a des Arg	0.40%
Human Complement C4	0.31%
Human Complement C5 des Arg	0.10%
Human Complement C5	0.05%

Abcam has not and does not intend to apply for the REACH Authorisation of customers' uses of products that contain European Authorisation list (Annex XIV) substances.

It is the responsibility of our customers to check the necessity of application of REACH Authorisation, and any other relevant authorisations, for their intended uses.

試験プラットフォーム

Microplate

製品の特性

保存方法

Please refer to protocols.

内容	1 x 96 tests
20X Wash Buffer Concentrate	1 x 30ml
Assay Buffer 10 Concentrate	1 x 15ml
Complement Reagent A	1 x 15ml
Complement Reagent B	1 x 30ml
Goat anti-rabbit IgG Microplate (12 x 8 wells)	1 unit
Human Complement C3a des Arg EIA Antibody	1 x 6ml
Human Complement C3a des Arg EIA Conjugate	1 x 6ml
Human Complement C3a des Arg Standard	2 x 500ng
Plate Sealer	2 units
pNpp Substrate	1 x 23ml
Stop Solution	1 x 5ml

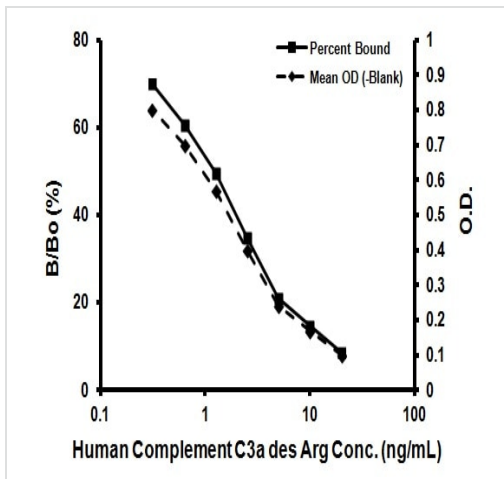
関連性

C3 plays a central role in the activation of the complement system. Its processing by C3 convertase is the central reaction in both classical and alternative complement pathways. After activation C3b can bind covalently, via its reactive thioester, to cell surface carbohydrates or immune aggregates. Derived from proteolytic degradation of complement C3, C3a anaphylatoxin is a mediator of local inflammatory process. It induces the contraction of smooth muscle, increases vascular permeability and causes histamine release from mast cells and basophilic leukocytes.

細胞内局在

Secreted

画像



Representative Standard Curve using ab133037

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