

Human Perforin ELISA Kit (PRF1) ab46068

7 References 画像数 3

製品の概要

製品名 Human Perforin ELISA Kit (PRF1)

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%
A	8	416pg/ml	20	5%
B	8	674pg/ml	25	4%
C	8	984pg/ml	24	3%

Inter-Assay (日差再現性)

サンプル	N	平均値	SD	CV%
A	9	580pg/ml	15	3%
B	9	665pg/ml	31	5%
C	9	1004pg/ml	45	5%

サンプルの種類 Cell culture supernatant

アッセイタイプ Sandwich (quantitative)

検出感度 < 40 pg/ml

検出範囲 62.5 pg/ml - 2000 pg/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
Cell culture supernatant	71	% - %

全工程の試験時間 2h 45m

ステップ Multiple steps standard assay

種交差性 交差種: Human

製品の概要

Abcam's Perforin (PRF1) Human *in vitro* ELISA (Enzyme-Linked Immunosorbent Assay) kit is designed for the quantitative measurement of Perforin in supernatant and buffered solutions.

A monoclonal antibody specific for Perforin has been coated onto the wells of the microtiter strips provided. Samples, including standards of known Perforin concentrations, control specimens or unknowns are pipetted into these wells. During the first incubation, the standards or samples and a biotinylated monoclonal antibody specific for Perforin are simultaneously incubated. After washing, the enzyme Streptavidin-HRP, that binds the biotinylated antibody is added, incubated and washed. A TMB substrate solution is added which acts on the bound enzyme to induce a colored reaction product. The intensity of this colored product is directly proportional to the concentration of Perforin present in the samples.

This kit will recognize both endogenous and recombinant Human Perforin.

Please note this kit is not suitable for use with serum and plasma.

試験プラットフォーム

Microplate

製品の特性

保存方法

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

内容	ラベル	2 x 96 tests	1 x 96 tests	1 x 96 tests	2 x 96 tests
1X Standard Diluent Buffer		2 x 25ml	1 x 25ml	1 x 25ml	2 x 25ml
200X Wash Buffer	White	2 x 10ml	1 x 10ml	1 x 10ml	2 x 10ml
Biotinylated anti PRF1	Red	2 x 0.4ml	1 x 0.4ml	1 x 0.4ml	2 x 0.4ml
Biotinylated Antibody Diluent	Red	1 x 13ml	1 x 7.5ml	1 x 7.5ml	1 x 13ml
Chromogen TMB Substrate Solution		1 x 24ml	1 x 11ml	1 x 11ml	1 x 24ml
HRP Diluent	Red	1 x 23ml	1 x 12ml	1 x 12ml	1 x 23ml
PRF1 Microplate (12 x 8 well strips)		2 units	1 unit	1 unit	2 units
PRF1 Standard (Lyophilized)	Yellow	4 vials	2 vials	2 vials	4 vials
Stop Reagent	Black	2 x 11ml	1 x 11ml	1 x 11ml	2 x 11ml
Streptavidin-HRP		4 x 5µl	2 x 5µl	2 x 5µl	4 x 5µl

機能

Plays a key role in secretory granule-dependent cell death, and in defense against virus-infected or neoplastic cells. Plays an important role in killing other cells that are recognized as non-self by the immune system, e.g. in transplant rejection or some forms of autoimmune disease. Can insert into the membrane of target cells in its calcium-bound form, oligomerize and form large pores. Promotes cytolysis and apoptosis of target cells by facilitating the uptake of cytotoxic granzymes.

関連疾患

Defects in PRF1 are the cause of hemophagocytic lymphohistiocytosis familial type 2 (FHL2) [MIM:603553]; also known as HPLH2. Familial hemophagocytic lymphohistiocytosis (FHL) is a

genetically heterogeneous, rare autosomal recessive disorder. It is characterized by immune dysregulation with hypercytokinemia and defective natural killer cell function. The clinical features of the disease include fever, hepatosplenomegaly, cytopenia, hypertriglyceridemia, hypofibrinogenemia, and neurological abnormalities ranging from irritability and hypotonia to seizures, cranial nerve deficits, and ataxia. Hemophagocytosis is a prominent feature of the disease, and a non-malignant infiltration of macrophages and activated T lymphocytes in lymph nodes, spleen, and other organs is also found.

配列類似性

Belongs to the complement C6/C7/C8/C9 family.

Contains 1 C2 domain.

Contains 1 EGF-like domain.

Contains 1 MACPF domain.

ドメイン

The C2 domain mediates calcium-dependent binding to lipid membranes. A subsequent conformation change leads to membrane insertion of beta-hairpin structures and pore formation. The pore is formed by transmembrane beta-strands.

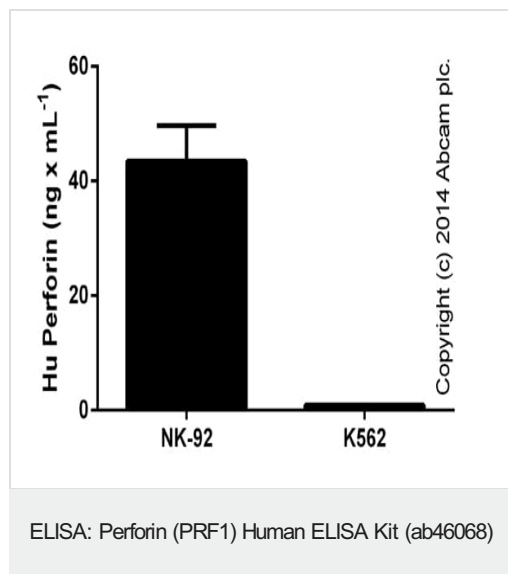
翻訳後修飾

N-glycosylated.

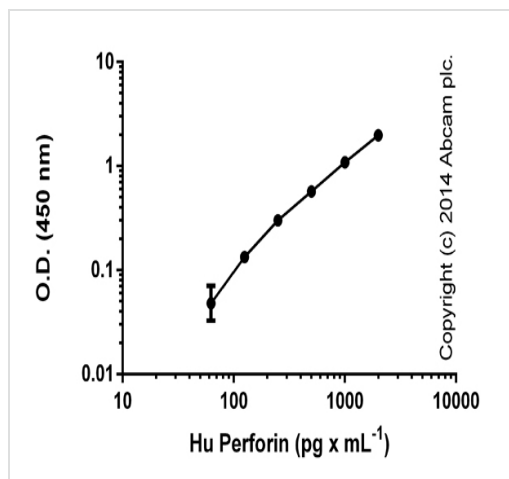
細胞内局在

Cytoplasmic granule lumen. Secreted. Cell membrane. Endosome lumen. Stored in cytoplasmic granules of cytolytic T-lymphocytes and secreted into the cleft between T-lymphocyte and target cell. Inserts into the cell membrane of target cells and forms pores. Membrane insertion and pore formation requires a major conformation change. May be taken up via endocytosis involving clathrin-coated vesicles and accumulate in a first time in large early endosomes.

画像

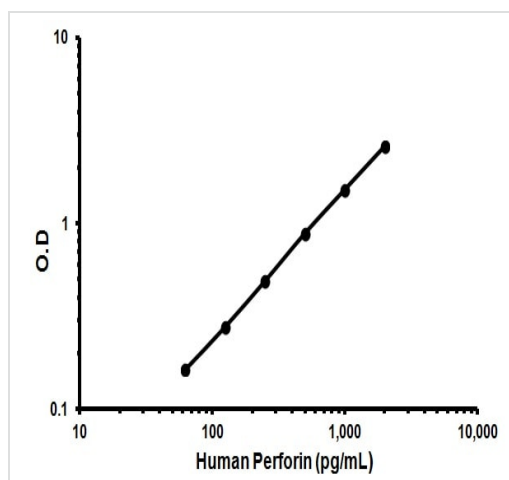


Perforin measured in cell culture supernatants (tested at 1/1-1/100 dilution range; duplicates +/- SD).



Standard curve with background signal subtracted (duplicates; +/- SD).

ELISA: Perforin (PRF1) Human ELISA Kit (ab46068)



Representative standard curve using ab46068

Typical data

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