

Human Perforin ELISA Kit (PRF1) ab46068

7 References 3 图像

概述

产品名称 人Perforin ELISA试剂盒(PRF1)

检测方法 Colorimetric

| 精确度 | 批次内 | | | | |
|-----|-----|---|----------|----|-----|
| | 样品 | n | Mean | SD | CV% |
| | A | 8 | 416pg/ml | 20 | 5% |
| | B | 8 | 674pg/ml | 25 | 4% |
| | C | 8 | 984pg/ml | 24 | 3% |

| | 批次间 | | | | |
|--|-----|---|-----------|----|-----|
| | 样品 | n | Mean | SD | CV% |
| | A | 9 | 580pg/ml | 15 | 3% |
| | B | 9 | 665pg/ml | 31 | 5% |
| | C | 9 | 1004pg/ml | 45 | 5% |

样品类型 Cell culture supernatant

检测类型 Sandwich (quantitative)

灵敏度 < 40 pg/ml

范围 62.5 pg/ml - 2000 pg/ml

| 回收率 | 特定样本回收率 | | |
|-----|--------------------------|-----|-------|
| | 样品类型 | 平均% | 范围 |
| | Cell culture supernatant | 71 | % - % |

检测时间 2h 45m

实验步骤 Multiple steps standard assay

种属反应性 与反应: Human

| | | | | | |
|--------------------------------------|--|--------------|--------------|--------------|--------------|
| 产品概述 | <p>Abcam's Perforin (PRF1) Human <i>in vitro</i> ELISA (Enzyme-Linked Immunosorbent Assay) kit is designed for the quantitative measurement of Perforin in supernatant and buffered solutions.</p> <p>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.</p> <p>This kit will recognize both endogenous and recombinant Human Perforin.</p> <p>Please note this kit is not suitable for use with serum and plasma.</p> | | | | |
| 平台 | Microplate | | | | |
| 性能 | | | | | |
| 存放说明 | Store at +4°C. Please refer to protocols. | | | | |
| 组件 | 标识符 | 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 |
| 功能 | <p>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.</p> | | | | |
| 疾病相关 | <p>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</p> | | | | |

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.

序列相似性

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

Contains 1 C2 domain.

Contains 1 EGF-like domain.

Contains 1 MACPF domain.

结构域

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.

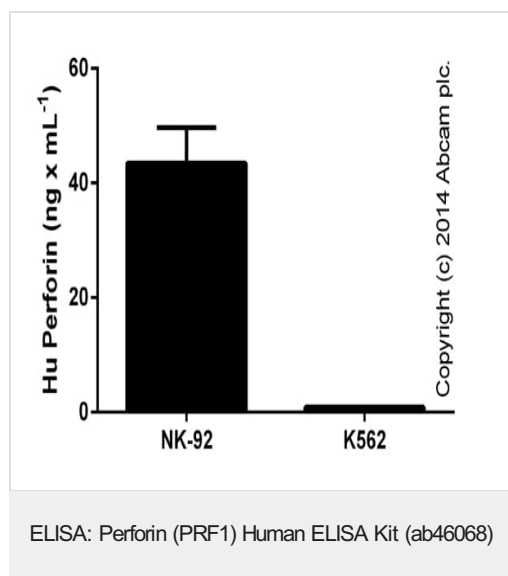
翻译后修饰

N-glycosylated.

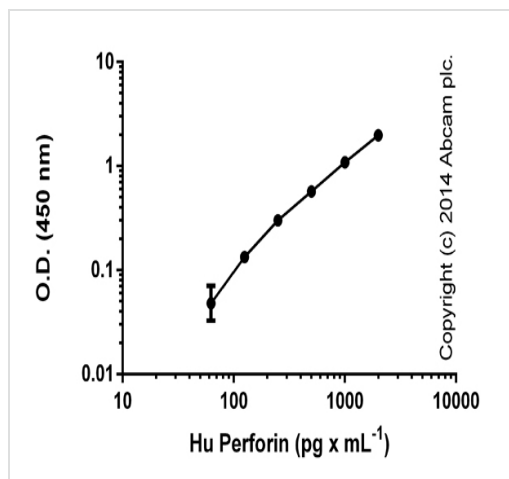
细胞定位

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.

图片

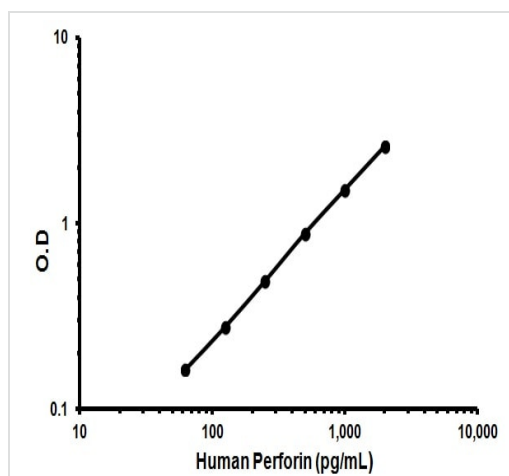


Perforin measured in cell culture supernatants (tested at 1/1-1/100 dilution range; duplicates +/- SD).



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.cn/abpromise> or contact our technical team.

Terms and conditions

- Guarantee only valid for products bought direct from Abcam or one of our authorized distributors