

Human Granulysin ELISA Kit ab213787

[1 References](#) [1 Image](#)

Overview

Product name Human Granulysin ELISA Kit

Detection method Colorimetric

Precision

Intra-assay

Sample	n	Mean	SD	CV%
1	16	57pg/ml	4.16	= 7.3%
2	16	258pg/ml	12.64	= 4.9%
3	16	911pg/ml	60.12	= 6.6%

Inter-assay

Sample	n	Mean	SD	CV%
1	24	56pg/ml	4.14	= 7.4%
2	24	266pg/ml	13.3	= 5%
3	24	859pg/ml	62.7	= 7.3%

Sample type Cell culture supernatant, Serum, Cell Lysate, Hep Plasma, EDTA Plasma

Assay type Sandwich (quantitative)

Sensitivity < 10 pg/ml

Range 31.2 pg/ml - 2000 pg/ml

Assay time 3h 30m

Assay duration Multiple steps standard assay

Species reactivity **Reacts with:** Human

Product overview Human Granulysin ELISA Kit (ab213787) is designed for the quantitative measurement of Human Granulysin in cell culture supernatants, cell lysates, serum and plasma (heparin, EDTA).

The ELISA kit is based on standard sandwich enzyme-linked immunosorbent assay technology. A monoclonal antibody from mouse specific for Granulysin has been precoated onto 96-well plates.

Standards (Expression system for standard: NSO; Immunogen sequence: M1-L145) and test samples are added to the wells, a biotinylated detection polyclonal antibody from goat specific for Granulysin is added subsequently and then followed by washing with PBS or TBS buffer. Avidin-Biotin-Peroxidase Complex was added and unbound conjugates were washed away with PBS or TBS buffer. HRP substrate TMB was used to visualize HRP enzymatic reaction. TMB was catalyzed by HRP to produce a blue color product that changed into yellow after adding acidic stop solution. The density of yellow is proportional to the Human Granulysin amount of sample captured in plate.

Notes Granulysin is a substance released by cytotoxic T cells (CD8) when they are attached to infected body cells. The product of this gene is a member of the saposin-like protein (SAPLIP) family. It is mapped to 2p11.2. Granulysin functions to create holes in the target cell membrane and destroy it. It is able to induce apoptosis in target cells and also has antimicrobial action. This gene is expressed in cytolytic granules with perforin, a pore forming protein, and granzymes that are also involved in cytolysis. In addition to it, Granulysin is broadly anti-microbial, killing microbes that cause, for example, tuberculosis and malaria, and can destroy some tumors. A series of peptides generated from the amino acid sequence of Granulysin are potential antibiotics. It has been found that secretory Granulysin is a key molecule responsible for the disseminated keratinocyte death in SJS/TEN.

Platform Pre-coated microplate (12 x 8 well strips)

Properties

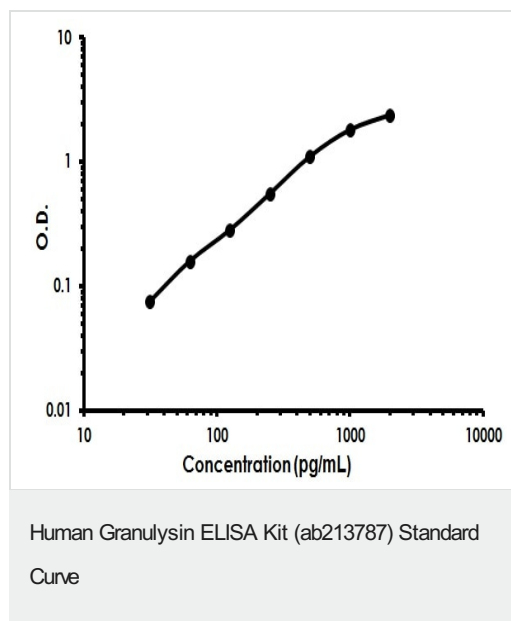
Storage instructions Store at -20°C. Please refer to protocols.

Components	Identifier	1 x 96 tests	1 x 96 tests
ABC Diluent Buffer	Blue Cap	1 x 12ml	1 x 12ml
Adhesive Plate Seal		4 units	4 units
Antibody Diluent Buffer	Green Cap	1 x 12ml	1 x 12ml
Anti-Human Granulysin Coated Microplate (12 x 8 wells)		1 unit	1 unit
Avidin-Biotin-Peroxidase Complex (ABC)		1 x 100µl	1 x 100µl
Biotinylated Anti-Human Granulysin Antibody		1 x 100µl	1 x 100µl
Lyophilized Recombinant Human Granulysin Standard		2 vials	2 vials
Sample Diluent Buffer	Green Cap	1 x 30ml	1 x 30ml
TMB Color Developing Agent	Black Cap	1 x 10ml	1 x 10ml
TMB Stop Solution	Yellow Cap	1 x 10ml	1 x 10ml
Wash Buffer (25X)		1 x 20ml	1 x 20ml

Function Antimicrobial protein that kills intracellular pathogens. Active against a broad range of microbes, including Gram-positive and Gram-negative bacteria, fungi, and parasites. Kills Mycobacterium tuberculosis.

Tissue specificity	Expressed in natural killer and T-cells.
Sequence similarities	Contains 1 saposin B-type domain.
Post-translational modifications	A 9 kDa form is produced by proteolytic processing of a 15 kDa protein.
Cellular localization	Secreted. Located in the cytotoxic granules of T-cells, which are released upon antigen stimulation.

Images



Human Granulysin ELISA Kit (ab213787) 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.com/abpromise> or contact our technical team.

Terms and conditions

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