

Granzyme B Inhibitor Screening Assay Kit ab157405

1 Image

Overview

Product name	Granzyme B Inhibitor Screening Assay Kit
Detection method	Fluorescent
Assay type	Quantitative
Product overview	<p>In Abcam's Granzyme B Inhibitor Screening Assay Kit, active human Granzyme B enzyme hydrolyzes the specific substrate to release the quench of fluorescent group, which can be detected fluorometrically at Ex/Em = 380/500 nm. In the presence of potent Granzyme B Inhibitor, the hydrolyzation of substrate will be impeded. The Kit provides a rapid, simple, sensitive and reliable test suitable for high throughput screening of Granzyme B inhibitors or characterize/study Granzyme B inhibitors. Inhibitor control, Ac-IEPD-CHO is included to compare the efficacy of test inhibitors.</p>
Notes	<p>This product is manufactured by BioVision, an Abcam company and was previously called K169 Granzyme B Inhibitor Screening Kit (Fluorometric). K169-100 is the same size as the 100 test size of ab157405.</p> <p>Granzyme B (GZMB, EC number 3.4.21.79) also known as Granzyme-2 is a serine protease stored in the granules of activated cytotoxic T cells and NK cells. Upon target cell contact, Granzyme B is directionally exocytosed and with the assistance of perforin enters the target cell. With its unique substrate specificity (Granzyme B prefers an aspartic acid residue at the P1 site of its substrates), Granzyme B processes and activates various pro-caspases thereby inducing apoptosis in the target cell.</p> <p>The standard is affinity purified.</p>
Platform	Microplate reader

Properties

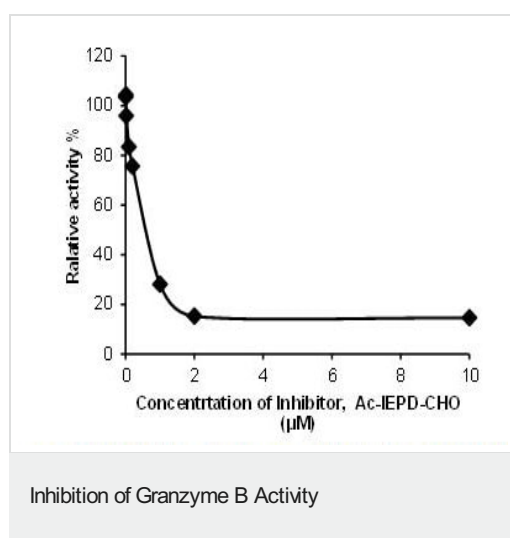
Storage instructionsStore at -20°C. Please refer to protocols.

Components	100 tests
Granzyme B Assay Buffer	1 x 25ml
Granzyme B Enzyme	1 vial
Granzyme B Substrate	1 x 0.5ml

Components	100 tests
Inhibitor Control (Ac-IEPD-CHO)	1 x 20µl

Function	This enzyme is necessary for target cell lysis in cell-mediated immune responses. It cleaves after Asp. Seems to be linked to an activation cascade of caspases (aspartate-specific cysteine proteases) responsible for apoptosis execution. Cleaves caspase-3, -7, -9 and 10 to give rise to active enzymes mediating apoptosis.
Sequence similarities	Belongs to the peptidase S1 family. Granzyme subfamily. Contains 1 peptidase S1 domain.
Cellular localization	Cytoplasmic granule. Cytoplasmic granules of cytolytic T-lymphocytes and natural killer cells.

Images



Inhibition of Granzyme B Activity by Granzyme B Inhibitor, Ac-IEPD-CHO. Assay was performed following Kit protocol.

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

