

Granzyme B Activity Assay Kit (Fluorometric) ab157403

1 References 1 图像

概述

产品名称	Granzyme B Activity Assay试剂盒(Fluorometric)
检测方法	Fluorescent
样品类型	Cell Lysate
检测类型	Enzyme activity (quantitative)
种属反应性	与反应: Mammals, Other species
产品概述	<p>In Abcam's Granzyme B Activity Assay Kit (Fluorometric), 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. This Assay Kit is rapid, reliable and sensitive and can detect Granzyme B activity in various biological samples.</p>
说明	<p>This product is manufactured by BioVision, an Abcam company and was previously called K168 Granzyme B Activity Fluorometric Assay Kit. K168-100 is the same size as the 100 test size of ab157403.</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>
平台	Microplate reader

性能

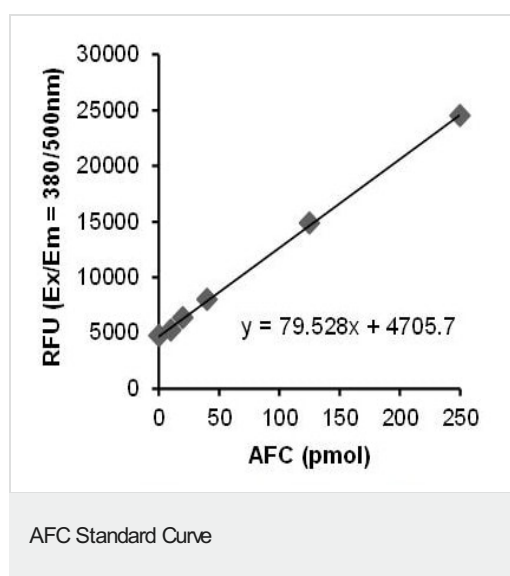
存放说明Store at -20°C. Please refer to protocols.

组件	100 tests
AFC Standard	1 x 0.1ml
Granzyme B Assay Buffer	1 x 25ml
Granzyme B Positive Control	1 vial

组件	100 tests
Granzyme B Substrate	1 x 0.5ml

功能	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.
序列相似性	Belongs to the peptidase S1 family. Granzyme subfamily. Contains 1 peptidase S1 domain.
细胞定位	Cytoplasmic granule. Cytoplasmic granules of cytolytic T-lymphocytes and natural killer cells.

图片



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