

Chitotriosidase Assay Kit ab241009

3 Images

Overview

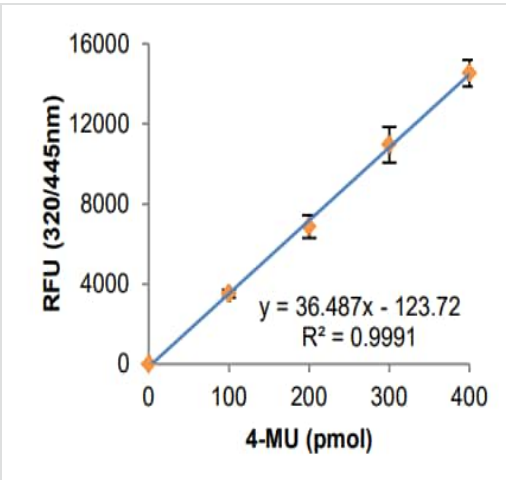
Product name	Chitotriosidase Assay Kit
Detection method	Fluorescent
Product overview	Chitotriosidase Assay Kit (ab241009) utilizes a fluorogenic substrate that can be hydrolyzed by chitinase and a set of proprietary assay buffers that can distinguish specific CHIT1 activity from other hydrolases including Acidic Mammalian Chitinase. The kit provides a simple, specific, sensitive assay that can detect as low as 0.2 mU/mL of CHIT1 in a variety of biological samples.
Notes	This product is manufactured by BioVision, an Abcam company and was previously called K512 Chitotriosidase Activity Assay Kit (Fluorometric). K512-100 is the same size as the 100 test size of ab241009.
Platform	Microplate reader

Properties

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

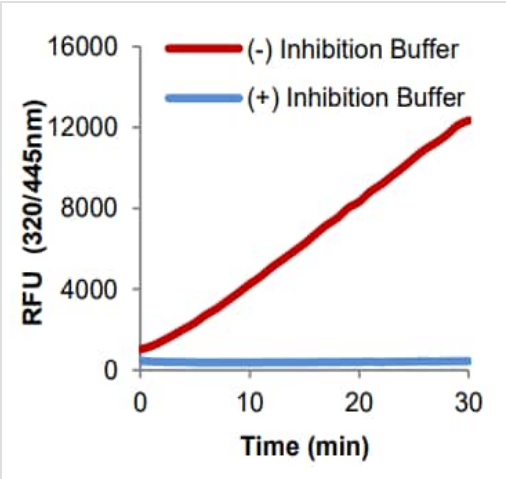
Components	100 tests
4-Methylumbelliferone Standard	1 x 35µl
Assay Buffer LI	1 x 25ml
CHIT1 Inhibition Buffer	1 x 18ml
Chitinase Substrate	1 x 25µl
Chitotriosidase	1 vial

Function	Degrades chitin, chitotriose and chitobiose. May participate in the defense against nematodes and other pathogens. Isoform 3 has no enzymatic activity.
Tissue specificity	Detected in spleen. Secreted by cultured macrophages.
Sequence similarities	Belongs to the glycosyl hydrolase 18 family. Chitinase class II subfamily. Contains 1 chitin-binding type-2 domain.
Cellular localization	Secreted. Lysosome. A small proportion is lysosomal.



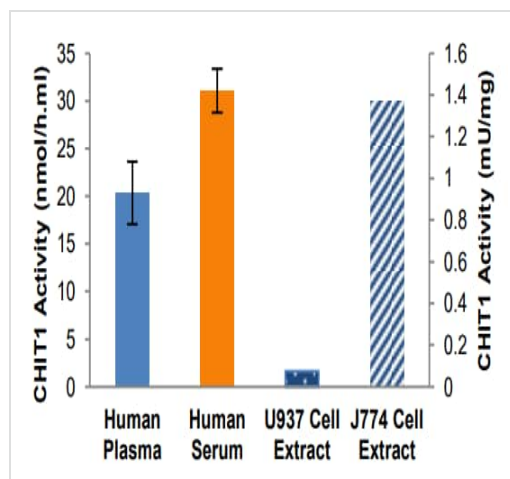
4-Methylumbelliferon Standard Curve.

4-Methylumbelliferon Standard Curve.



Measurement of purified Human Chitotriosidase activity with or without Inhibition Buffer

Measurement of purified Human Chitotriosidase activity with or without Inhibition Buffer



Measurement of CHIT1 activity in Human Plasma (4 μ L), Human Serum (4 μ L), U937 Cell Extract (10 μ g) and J774 Cell Extract (10 μ g). All assays were performed following kit protocol.

Measurement of CHIT1 activity

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