

Chymotrypsin Assay Kit (Fluorometric) ab234051

1 References 画像数 6

製品の概要

製品名	Chymotrypsin Assay Kit (Fluorometric)
検出方法	Fluorescent
サンプルの種類	Cell Lysate, Tissue Lysate
種交差性	交差種: Mammals, Other species
製品の概要	Chymotrypsin Assay Kit (Fluorometric) (ab234051) uses a synthetic fluorogenic substrate, enabling kinetic measurement of chymotrypsin activity in cell and tissue lysates. A chymotrypsin activator cleaves chymotrypsinogen to form active chymotrypsin, which then hydrolyzes the non-fluorescent substrate to release a stable fluorophore. The kit includes a selective chymotrypsin inhibitor that can be used to measure specific chymotrypsin activity in samples containing non-specific proteases and endopeptidases that may also metabolize the substrate. The assay can detect as low as 0.01 mU of Chymotrypsin.
特記事項	<p>This product is manufactured by BioVision, an Abcam company and was previously called K352 Chymotrypsin Activity Assay Kit (Fluorometric). K352-100 is the same size as the 100 test size of ab234051.</p> <p>Chymotrypsin (EC 3.4.21.1) is a key serine protease involved in dietary protein digestion in mammals. It is primarily produced by the pancreas, but may be expressed in other tissues, including spleen and liver. In the pancreas, chymotrypsin is initially expressed as the inactive proenzyme chymotrypsinogen, which is cleaved by other proteases to chymotrypsin, its active form. Chymotrypsin specifically cleaves peptide bonds at the C-terminal end of bulky hydrophobic or aromatic amino acids (such as tyrosine, tryptophan or phenylalanine).</p>
試験プラットフォーム	Microplate reader

製品の特性

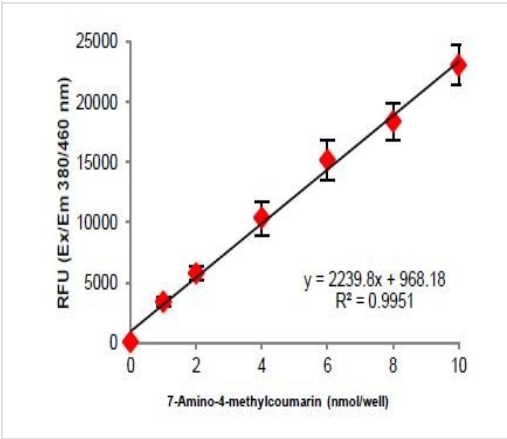
保存方法 Store at -20°C. Please refer to protocols.

内容	100 tests
AMC Standard	1 x 100µl
Assay Buffer I	1 x 25ml
Chymotrypsin Activator	1 vial

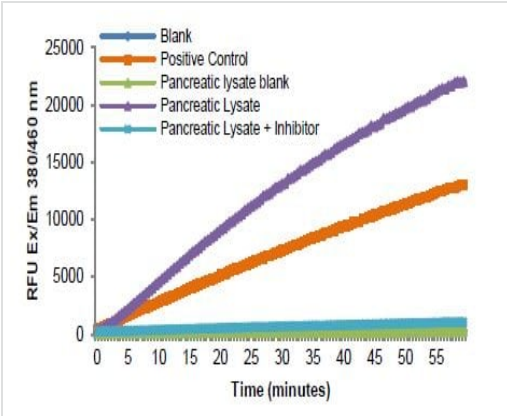
内容	100 tests
Chymotrypsin Inhibitor I	1 x 80µl
Chymotrypsin Positive Control	1 vial
Chymotrypsin Substrate	1 x 200µl

機能	Has chymotrypsin-type protease activity and hypocalcemic activity.
組織特異性	Pancreas.
関連疾患	Variations in CTRC influence susceptibility to chronic pancreatitis [MM:167800]. Chronic pancreatitis is a persistent inflammatory disorder characterized by permanent destruction of the pancreatic parenchyma.
配列類似性	Belongs to the peptidase S1 family. Elastase subfamily. Contains 1 peptidase S1 domain.

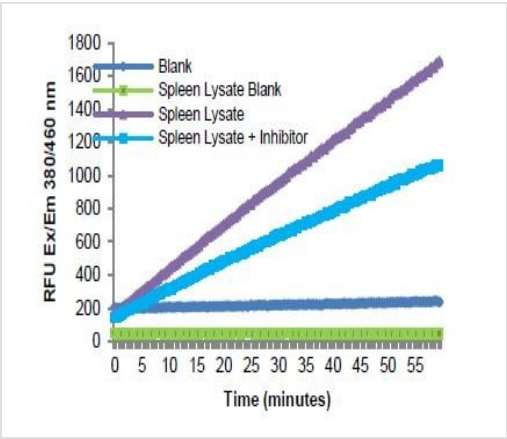
画像



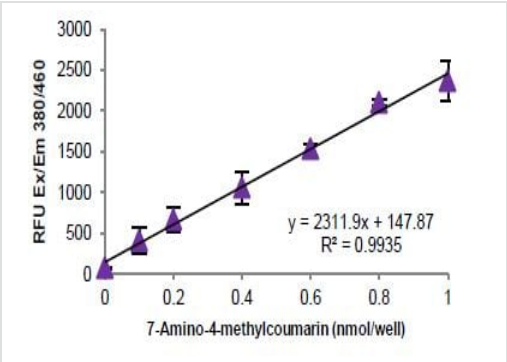
Coumarin standard curves showing a range of 0-10 nmol/well.



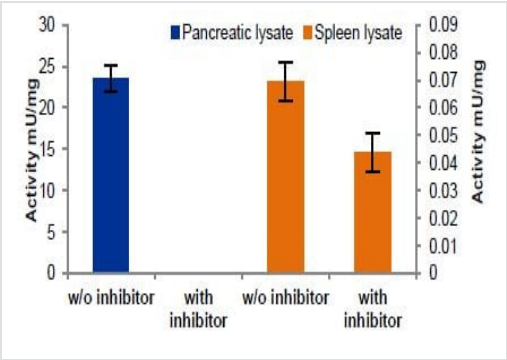
Reaction kinetics for positive control and rat pancreatic lysate (9 µg).



Reaction kinetics for rat spleen lysate (160 µg).

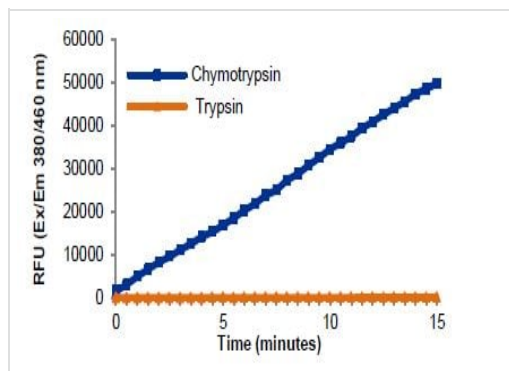


Coumarin standard curves showing a range of 0-1 nmol/well.



Chymotrypsin activity with and without inhibitor in rat pancreas and rat spleen.

The presence of non-specific chymotrypsin-like proteases in spleen leads to some activity in presence of the selective chymotrypsin inhibitor.



The substrate is cleaved by chymotrypsin but not trypsin, making the assay free from trypsin interference.

Reaction kinetics using substrate in the presence of Chymotrypsin or Trypsin.

**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.co.jp/abpromise> or contact our technical team.

### Terms and conditions

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