

## Product datasheet

# Enterokinase Activity Assay Kit (Fluorometric) ab170964

[2 Images](#)

### Overview

<b>Product name</b>	Enterokinase Activity Assay Kit (Fluorometric)
<b>Detection method</b>	Fluorescent
<b>Sample type</b>	Purified protein
<b>Assay type</b>	Enzyme activity (quantitative)
<b>Sensitivity</b>	>= 1 $\mu$ U
<b>Species reactivity</b>	<b>Reacts with:</b> Mammals, Other species
<b>Product overview</b>	Abcam's Enterokinase Activity Assay Kit (Fluorometric) utilizes a peptide substrate containing the Enterokinase recognition sequence along with a fluorescent label 'AFC'. Enterokinase catalyzes the cleavage of this substrate and releases the AFC molecule, which can be easily quantified by measuring its fluorescence at Ex/Em = 380/500 nm. This assay kit is simple and rapid and can detect Enterokinase activity as low as 1 mU.
<b>Notes</b>	<p>This product is manufactured by BioVision, an Abcam company and was previously called K758 Enteropeptidase/Enterokinase Activity Fluorometric Assay Kit. K758-100 is the same size as the 100 test size of ab170964.</p> <p>Enterokinase (Enteropeptidase, EC 3.4.21.9) is a serine protease involved in activation of trypsinogen to trypsin, which in turn results in the activation of various digestive enzymes. It recognizes a highly specific amino acid sequence 'DDDDK' and cleaves after the lysine residue. High specific activity of Enterokinase has been utilized in cleaving a variety of native or fusion protein tags containing the above recognition motif.</p>
<b>Platform</b>	Microplate reader

### Properties

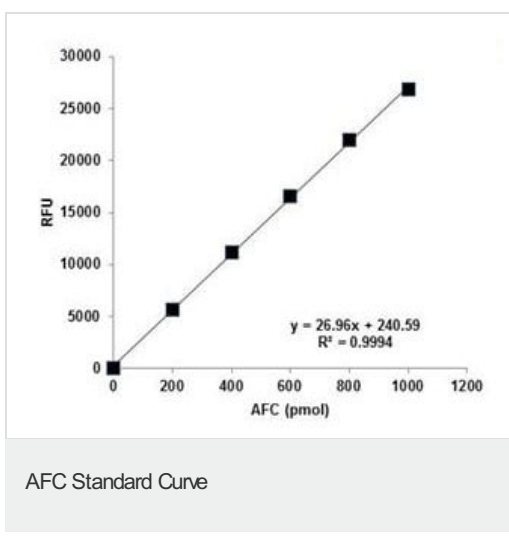
**Storage instructions** Store at -20°C. Please refer to protocols.

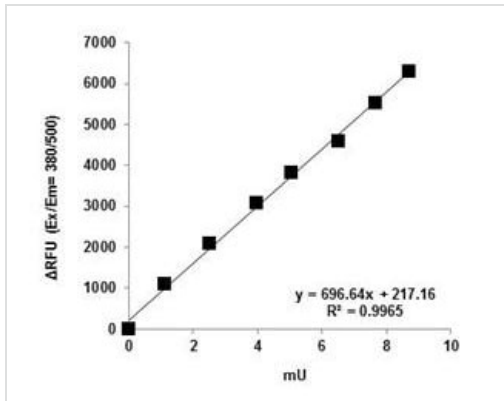
Components	Identifier	100 tests
AFC Standard	Yellow	1 x 100 $\mu$ l
Enteropeptidase Assay Buffer	WM	1 x 20ml
Enteropeptidase Substrate	Red	1 x 0.2ml

Components	Identifier	100 tests
Human Enteropeptidase	Green	1 x 50µl

<b>Function</b>	Responsible for initiating activation of pancreatic proteolytic proenzymes (trypsin, chymotrypsin and carboxypeptidase A). It catalyzes the conversion of trypsinogen to trypsin which in turn activates other proenzymes including chymotrypsinogen, procarboxypeptidases, and proelastases.
<b>Tissue specificity</b>	Intestinal brush border.
<b>Involvement in disease</b>	Defects in TMPRSS15 are a cause of enterokinase deficiency (ENTKD) [MIM:226200]; a life-threatening intestinal malabsorption disorder characterized by diarrhea and failure to thrive.
<b>Sequence similarities</b>	Belongs to the peptidase S1 family. Contains 2 CUB domains. Contains 2 LDL-receptor class A domains. Contains 1 MAM domain. Contains 1 peptidase S1 domain. Contains 1 SEA domain. Contains 1 SRCR domain.
<b>Post-translational modifications</b>	The chains are derived from a single precursor that is cleaved by a trypsin-like protease.
<b>Cellular localization</b>	Membrane.

**Images**





Measurement of Enterokinase Levels

**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