

# Enterokinase cleavage kit ab207001

## 2 图像

### 概述

产品名称	Enterokinase cleavage试剂盒
产品概述	<p>Abcam's Enterokinase cleavage kit (ab207001) efficiently remove tags from recombinant fusion proteins containing accessible enteropeptidase-specific recognition sequence.</p> <p>Enteropeptidase (enterokinase, EC 3.4.21.9) is a serine protease involved in activation of trypsinogen to trypsin. It recognizes a highly specific amino acid sequence 'DDDDK' and cleaves after the lysine (K) residue. The high specific activity of enteropeptidase is also used to cleave native and fusion proteins containing this recognition motif, and is often used to remove the tag from a recombinant protein. Abcam's Enterokinase cleavage kit contains a highly active light chain fragment of human enteropeptidase. Our pure enzyme is an excellent tool to obtain a wild type protein sequence from a fusion protein that contains the enteropeptidase recognition sequence. This kit is sufficient for cleaving at least 5 mg of target protein. The residual enteropeptidase left in the reaction mix will not interfere with most experiments that the target protein will be used in. Following cleavage of the target protein, the enteropeptidase may be removed using nickel or cobalt beads if required. However, this is only possible if the target protein contains a histidine residue. If no histidine residue is present, enteropeptidase cannot be removed.</p>
说明	<p>This product is manufactured by BioVision, an Abcam company and was previously called K760 Enteropeptidase/Enterokinase Cleavage Kit. K760-100 is the same size as the 100 test size of ab207001.</p>
经测试应用	适用于: Purification

### 性能

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

组件	100 tests
Cleavage Control Protein	1 vial
Enteropeptidase Assay Buffer	1 x 20ml
Human Enteropeptidase II	1 vial

## 应用

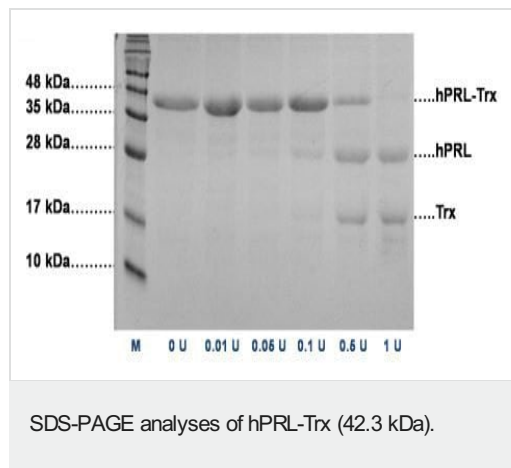
### The Abpromise guarantee

**Abpromise™** 承诺保证使用ab207001于以下的经测试应用

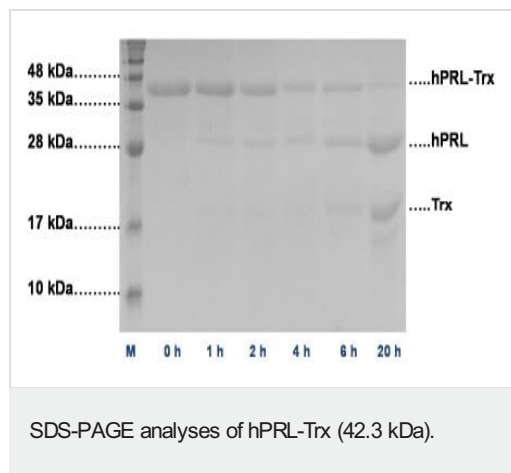
“应用说明”部分 下显示的仅为推荐的起始稀释度;实际最佳的稀释度/浓度应由使用者检定。

应用	Ab评论	说明
Purification		Use at an assay dependent concentration.

## 图片



SDS-PAGE analysis of the cleavage of hPRL-Trx using different amounts of enteropeptidase. hPRL-Trx (50 µg) was digested into its individual protein fragments hPRL (25.3 kDa) and Trx (17 kDa) using different amounts (0.01-1 U) of enteropeptidase at room temperature for 20 hours.



SDS-PAGE analysis of the cleavage of hPRL-Trx by enteropeptidase over time: Analysis of hPRL-Trx digestion at different time points at room temperature using 1 U of enteropeptidase.

**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