

| | |
|---------------------|---------------|
| E535-31H-5 | 5 µg |
| E535-31H-10 | 10 µg |
| E535-31H-100 | 100 µg |
| E535-31H-500 | 500 µg |

EK (enterokinase), Active

Recombinant swine protein expressed in yeast cells

Catalog # E535-31H

Lot # L2146-3

Product Description

Recombinant swine EK (800-end) was expressed in yeast cells using an N-terminal His tag. The enzyme commission number is EC 3.4.21.9.

Gene Aliases

Enteropeptidase; TMPRSS15; PRSS7; Serine protease 7; ENTK; MGC133046.

Formulation

Recombinant protein stored in 50mM Tris-HCl, pH 8, 200mM NaCl, 30% glycerol.

Storage and Stability

Store product at -20°C for up to 1 year. Aliquot enzymes to avoid freeze / thaw cycles.

Digestion Conditions

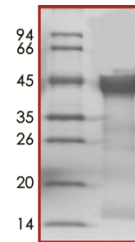
Catalytic pH Range: 4.5 ~ 9.5

Catalytic Temperature Range: 4 ~ 45°C

Scientific Background

Enterokinase (EK) is a member of the trypsin family of serine proteases, which cleaves proteins following the Lys at the Asp-Asp-Asp-Asp-Lys recognition sequence. In the stomach, EK converts trypsinogen into trypsin. EK has high specificity and high hydrolytic efficiency, making it a widely used biochemical tool for cleaving recombinant fusion proteins. EK can tolerate a variety of detergents and some denaturing agents. EK's activity is reduced by > 250 mM NaCl, > 2 M urea, > 20 mM, β-ME, > 0.1% SDS and > 250 mM imidazole.

Purity

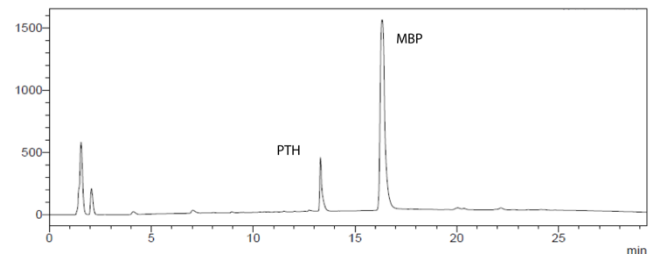


← Enterokinase
The purity of EK (enterokinase) was determined to be **>85%** by densitometry, approx. MW **44kDa**.

Specific Activity

The specific activity of EK was determined to be **>10,000 units/mg** as per the activity assay protocol.

Sample Data:



HPLC results for digestion of the parathyroid hormone - maltose binding protein (MBP-PTH) fusion protein by Enterokinase at an enzyme : substrate mass ratio of 1:250 for 8 hr at 25°C. The digestion products are: PTH and MBP.

EK (enterokinase), Active

Recombinant swine protein expressed in yeast cells

| | |
|--------------------|--|
| Catalog Number | E535-31H |
| Specific Activity | 10,000 units/mg |
| Lot # | L2146-3 |
| Purification | Affinity purification; his tagged |
| Purity | >85% |
| Concentration | 0.1 mg/ml |
| Stability | 1 yr at -20°C from date of shipment |
| Storage & Shipping | Store as supplied at -20°C for up to 1 year. |

To place your order, please contact us by phone 1-(604)-232-4600, fax 1-604-232-4601 or by email: orders@signalchem.com
www.signalchem.com

FOR IN VITRO RESEARCH PURPOSES ONLY. NOT INTENDED FOR USE IN HUMAN OR ANIMALS.

Assay Protocol

Reaction Components

Active EK Protease (Catalog #: E535-31H)

Active EK protease.

EK Digestion Buffer (User Prepared)

20 mM Tris-HCl pH 8, 200 mM NaCl, 2mM CaCl₂

Reaction Optimization Protocol

The following conditions may be different for different proteins. Optimize the protocol for each specific protein.

Step 1. Dissolve fusion protein in EK Digestion Buffer

Step 2. Add EK to fusion protein solution at a mass ratio of 1:50 and mix

Step 3. Incubate reaction mixture at room temperature for 2-16 hours

Step 4. Monitor the cleavage products by SDS-PAGE. Run an undigested sample of the target fusion protein as a control

Activity Definition (units/mg)

SignalChem's Enterokinase activity is defined by the following:

One unit equals the amount of EK required to produce 1 nmole of trypsin from trypsinogen per minute at pH 8.0 at 25 °C.

To place your order, please contact us by phone 1-(604)-232-4600, fax 1-604-232-4601 or by email: orders@signalchem.com
www.signalchem.com

FOR IN VITRO RESEARCH PURPOSES ONLY. NOT INTENDED FOR USE IN HUMAN OR ANIMALS.