

HtrA Protein

Recombinant bacterial protein expressed in E. coli cells

Catalog # H531-30CH

Lot # Z1336-3

Product Description

Recombinant streptococcus HtrA (33-393) protein was expressed in bacterial cells using an N-terminal His-tag. The gene is cloned from bacterial strain ATCC 700669 and the protein accession number is [B8ZQD0](#).

Gene Aliases

SPN23F22720

Formulation

Recombinant protein stored in 50mM Tris-HCl, pH 8.0, 300mM NaCl, 150mM imidazole and 10% glycerol.

Storage and Stability

Store product at -70°C. For optimal storage, aliquot target into smaller quantities after centrifugation and store at recommended temperature. For most favorable performance, avoid repeated handling and multiple freeze/thaw cycles.

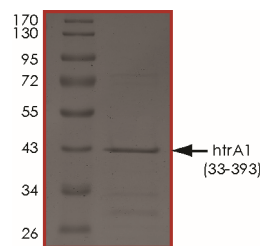
Scientific Background

HtrA (also designated as DegP) is a trypsin-like serine protease which contains PDZ domain and involves in heat-shock response, chaperone function, and apoptosis (1). HtrA is an error sensor that specifically represses genetic competence when the overall level of biosynthetic errors in cellular proteins is low (2).

References

1. Ge X. et.al: DegP primarily functions as a protease for the biogenesis of β -barrel outer membrane proteins in the Gram-negative bacterium Escherichia coli. FEBS J. 2013 Dec 26. doi: 10.1111/febs.12701.
2. Cassone M. et.al: The HtrA protease from Streptococcus pneumoniae digests both denatured proteins and the competence-stimulating peptide. J Biol Chem. 2012 Nov 9;287(46):38449-59.

Purity



The purity of HtrA was determined to be **>85%** by densitometry. Approx. MW **41 kDa**.

HtrA Protein

Recombinant bacterial protein expressed in E. coli cells

Catalog #	H531-30CH
Lot #	Z1336-3
Purity	>85%
Concentration	0.1 µg/µl
Stability	1yr at -70°C from date of shipment
Storage & Shipping	Store product at -70°C. For optimal storage, aliquot target into smaller quantities after centrifugation and store at recommended temperature. For most favorable performance, avoid repeated handling and multiple freeze/thaw cycles. Product shipped on dry ice.

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.