

HtrA (S234A) Protein

Recombinant bacterial protein expressed in *E. coli* cells

Catalog # H531-36CH

Lot # Z1324-1

Product Description

Recombinant streptococcus HtrA (S234A) (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](#). This protein has a serine protease dead mutation.

Gene Aliases

DegP, DegP/HtrA

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.

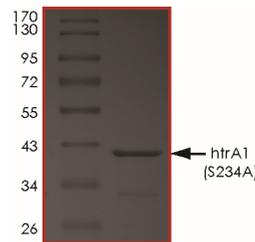
Scientific Background

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). Analysis of strains with the htrA (S234A) catalytic site mutation showed that the proteolytic activity of HtrA selectively repressed competence when translational fidelity was high but not when accuracy was 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. Stevens KE. et.al: Competence in *Streptococcus pneumoniae* is regulated by the rate of ribosomal decoding errors. MBio. 2011 Sep 20;2(5). pii: e00071-11.

Purity



The purity of HtrA (S234A) was determined to be **>95%** by densitometry. Approx. MW **40 kDa**.

HtrA (S234A) Protein

Recombinant bacterial protein expressed in *E. coli* cells

Catalog #	H531-36CH
Lot #	Z1324-1
Purity	>95%
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.