

HTRA4 Protein

Recombinant human protein expressed in E. coli cells

Catalog # H534-31H

Lot # Y972-4

Product Description

Recombinant human HTRA4 (159-376) protein was expressed in E. coli cells using an N-terminal His-tag. The protein accession number is [NM_153692](#). This protein contains a catalytic domain of serine protease HTRA4.

Gene Aliases

(None)

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.

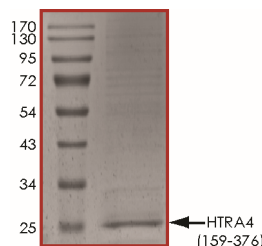
Scientific Background

HTRA1 is a serine protease with a conserved trypsin domain and a PDZ domain. As other related family members, HTRA4 may function as a secreted oligomeric chaperone protease to degrade misfolded secretory proteins. Htra4 may contribute to shallow trophoblast invasion in preeclampsia (1) and imply a potential application as a biomarker for this disorder (2).

References

1. Wang LJ. et.al: High-temperature requirement protein A4 (Htra4) suppresses the fusogenic activity of syncytin-1 and promotes trophoblast invasion. Mol Cell Biol. 2012 Sep;32(18):3707-17.
2. Inagaki A. et.al: Upregulation of Htra4 in the placentas of patients with severe pre-eclampsia. Placenta. 2012 Nov;33(11):919-26.

Purity



The purity of HTRA4 was determined to be **>80%** by densitometry. Approx. MW **25 kDa**.

HTRA4 Protein

Recombinant human protein expressed in E. coli cells

Catalog #	H534-31H
Lot #	Y972-4
Purity	>80%
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.