

HTRA3 Protein

Recombinant human protein expressed in *E. coli* cells

Catalog # H533-31H

Lot # Y935-5

Product Description

Recombinant human HTRA3 (132-353) protein was expressed in *E. coli* cells using an N-terminal His-tag. The protein accession number is [NM_053044](#).

Gene Aliases

PRSP; TASP

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.

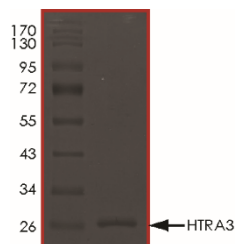
Scientific Background

High-temperature requirement factor A3 (HTRA3) is a member of the HTRA family of serine proteases that was initially identified in *E. coli* by two phenotypes of null mutants that were unable to grow at elevated temperatures. HTRA3 was discovered initially as a pregnancy-related serine protease (PRSP) identified by DDPCR (Differential Display Polymerase Chain Reaction) to be differentially expressed between implantation and interimplantation sites in the mouse uterus. A reduction of HTRA3 mRNA expression in ovarian cancer suggested that HTRA3 may act as a tumor suppressor.

References

1. Strauch KL, et al: An Escherichia coli mutation preventing degradation of abnormal periplasmic proteins. Proc Natl Acad Sci U S A. 85(5):1576-1580. 1988.
2. Nie GY, et al. A novel serine protease of the mammalian HtrA family is up-regulated in mouse uterus coinciding with placentation. Mol Hum Reprod. 9(5):279-290. 2003.
3. Bowden MA, et al: High-temperature requirement factor A3 (Htra3): a novel serine protease and its potential role in ovarian function and ovarian cancers. Mol Cell Endocrinol. 327(1-2):13-8. 2010.

Purity



The purity of HTRA3 was determined to be **>95%** by densitometry. Approx. MW **27 kDa**.

HTRA3 Protein

Recombinant human protein expressed in *E. coli* cells

Catalog #	H533-31H
Lot #	Y935-5
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.