

HTRA1 (161-379) Protein

Recombinant human protein expressed in E. coli cells

Catalog # H531-32H

Lot # T1039-4

Product Description

Recombinant human HTRA1 (161-379) protein was expressed in E. coli cells using an N-terminal His-tag. The protein accession number is [NM_002775](#).

Gene Aliases

ARMD7; CARASIL; HtrA; L56; ORF480; PRSS11

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.

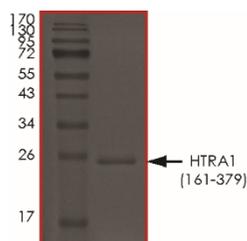
Scientific Background

HTRA1 is a serine protease. HTRA1 targets a variety of proteins (1 & 2), including fibronectin (induce synovial cells to up-regulate MMP1 and MMP3 production), proteoglycans (release soluble FGF-glycosaminoglycan complexes that promote the range and intensity of FGF signals in the extracellular space), IGF-binding proteins (regulates the availability of insulin-like growth factors) and TSC2 (leads to the activation of TSC2 downstream targets). HTRA1 relates to age-related macular degeneration, a multifactorial eye disease and the most common cause of irreversible vision loss (3).

References

- Hu S.J. et.al: Human HtrA, an evolutionarily conserved serine protease identified as a differentially expressed gene product in osteoarthritic cartilage. J. Biol. Chem. 273:34406-34412 (1998).
- Grau S. et.al: The role of human HtrA1 in arthritic disease. J. Biol. Chem. 281:6124-6129 (2006).
- Yang, Z. et.al: A variant of the HTRA1 gene increases susceptibility to age-related macular degeneration. Science 314: 992-993, 2006.

Purity



The purity of HTRA1 (161-379) was determined to be **>95%** by densitometry. Approx. MW **25 kDa**.

HTRA1 (161-379) Protein

Recombinant human protein expressed in E. coli cells

Catalog #	H531-32H
Lot #	T1039-4
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.