

ASK3 Protein

Recombinant human protein expressed in Sf9 cells

Catalog # M73-31G

Lot # K1802-3

Product Description

Recombinant human ASK3 (MAP3K15) (582-1169) was expressed by baculovirus in Sf9 insect cells using an N-terminal GST tag. This gene accession number is [NM_001001671](#).

Gene Aliases

MAP3K15; ASK3; bA723P2.3

Formulation

Recombinant protein stored in 50mM Tris-HCl, pH 7.5, 150mM NaCl, 10mM glutathione, 0.1mM EDTA, 0.25mM DTT, 0.1mM PMSF, 25% 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.

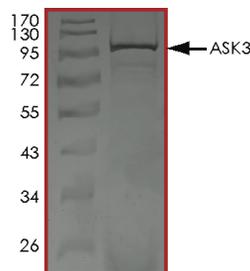
Scientific Background

ASK3 (MAP3K15) is a member of the mitogen-activated protein kinase (MAPK) family which functions in a protein kinase signal transduction cascade, where an activated MAPK kinase kinase (MAP3K) phosphorylates and activates a specific MAPK kinase (MAP2K), which then activates a specific MAPK. ASK3 plays an essential role in apoptotic cell death triggered by cellular stresses (1). ASK3 is also functions in a signal transduction pathway that is activated by various cell stresses and leads to apoptosis (2).

References

- Hartz, P. A.et.al: Personal Communication. Baltimore, Md. 7/26/2010.
- Kaji, T. et.al: ASK3, a novel member of the apoptosis signal-regulating kinase family, is essential for stress-induced cell death in HeLa cells.

Purity



The purity of ASK3 protein was determined to be **>85%** by densitometry.
Approx. MW **93 kDa**.

ASK3 Protein

Recombinant human protein expressed in Sf9 cells

Catalog #	M73-31G
Lot #	K1802-3
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.