

Mortalin Protein

Recombinant human protein expressed in Sf9 cells

Catalog # M54-31G

Lot # R255-1

Product Description

Recombinant human Mortalin (47-end) protein was expressed in Sf9 insect cells using an N-terminal GST tag. The gene accession number is [NM_004134](#).

Gene Aliases

HSPA9; CSA; GRP75; HSPA9B; MGC4500; MOT; mot-2; MOT2; MTHSP75; PBP74

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

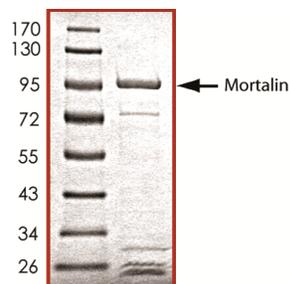
Mortalin is a mitochondrial chaperone and a member of the heat shock protein 70 family that is constitutively expressed in cells. Mortalin plays a central role in mitochondrial biogenesis through its capacity to direct the import of nuclear-encoded proteins into the mitochondria. Mortalin plays a role in the control of cell proliferation and elevated levels of mortalin have correlated with malignant transformation and poor cancer prognosis. Mortalin can support cancer cell resistance to complement-dependent cytotoxicity (1). Mortalin expression was decreased in the mitochondrial fraction of neurons from the substantia nigra of Parkinson disease patients (2).

References

1. Pilzer, D. et al: Mortalin inhibitors sensitize K562 leukemia cells to complement-dependent cytotoxicity. *Int J Cancer*. 2010 Mar 15;126(6):1428-35.
2. De Mena, L. et al: Mutational screening of the mortalin gene (HSPA9) in Parkinson's disease. *J Neural Transm*. 2009 Oct;116(10):1289-93.

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

Purity



The purity was determined to be **>80%** by densitometry. Approx. MW **97kDa**.

Mortalin Protein

Recombinant human protein expressed in Sf9 insect cells

Catalog Number	M54-31G
Specific Lot Number	R255-1
Purity	>80%
Concentration	0.2 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.