

AATK Protein

Recombinant protein expressed in Sf9 cells

Catalog # A02-35G

Lot # N356-1

Product Description

Recombinant mouse AATK (8-411) was expressed by baculovirus in Sf9 insect cells using an N-terminal GST tag. The gene accession number is [NM_007377](#).

Gene Aliases

AATYK; LMTK1; KIAA0641

Formulation

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

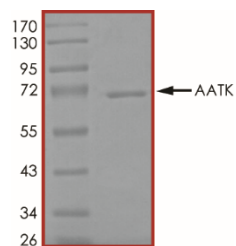
Scientific Background

AATK or apoptosis-associated tyrosine kinase contains a tyrosine kinase domain at the N-terminus and a proline-rich domain at the C-terminus. AATK is induced during apoptosis, and expression of this protein is a necessary pre-requisite for the induction of growth arrest and/or apoptosis of myeloid precursor cells (1). AATK is highly detected in brain, lung, kidney, and pancreas (2). AATK is also shown to produce neuronal differentiation in a neuroblastoma cell line.

References

- Gaozza, E. et.al: AATYK: a novel tyrosine kinase induced during growth arrest and apoptosis of myeloid cells. *Oncogene* 15: 3127-3135, 1997.
- Ishikawa, K. et.al: Prediction of the coding sequences of unidentified human genes. X. The complete sequences of 100 new cDNA clones from brain which can code for large proteins in vitro. *DNA Res.* 5: 169-176, 1998

Purity



The purity was determined to be **>95%** by densitometry. Approx. MW **71 kDa**.

AATK Protein

Recombinant protein expressed in Sf9 cells

Catalog #	A02-35G
Lot #	N356-1
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.