

MYT1 Protein

Full length recombinant protein expressed in Sf9 cells

Catalog # M67-34G

Lot # N209-2

Product Description

Recombinant full-length human MYT1 was expressed by baculovirus in Sf9 insect cells using an N-terminal GST tag. The gene accession number is [NM_004203](#).

Gene Aliases

PKMYT1, FLJ20093, DKFZp547K1610

Formulation

Recombinant protein stored in 50mM Tris-HCl, pH 7.5, 50mM 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

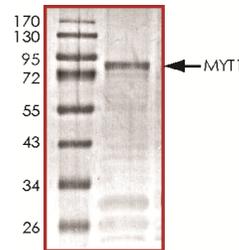
MYT1 is a member of a family of neural specific, zinc finger-containing DNA-binding proteins which binds to the promoter regions of proteolipid proteins of the central nervous system and plays a role in the developing nervous system. MYT1 of the oligodendrocyte lineage, along with a closely related CCHC zinc finger, is expressed in developing neurons in the mammalian central nervous system(1). The recombinant MYT1 fragments containing either the upstream 2 zinc fingers or the downstream 4 zinc fingers bound the same cis regulatory element in the PLP1 promoter in vitro and DNA binding required Zn(2+), but not other divalent cations (2).

References

1. Kim, J. G. et.al: Myelin transcription factor 1 (Myt1) of the oligodendrocyte lineage, along with a closely related CCHC zinc finger, is expressed in developing neurons in the mammalian central nervous system. *J. Neurosci. Res.* 50: 272-290, 1997.
2. Kim, J. G. et.al: Novel member of the zinc finger superfamily: a C(2)-HC finger that recognizes a glia-specific gene. *Molec. Cell. Biol.* 12: 5632-5639, 1992.

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 of MYT1 was determined to be **>80%** by densitometry. Approx. MW **93 kDa**.

MYT1 Protein

Full length recombinant protein expressed in Sf9 cells

Catalog Number	M67-34G
Specific Lot Number	N209-2
Purity	>80%
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.