

Anti-MYT1L

Rabbit Polyclonal Antibody

Catalog # M92-363R

Lot # O2121-51

Cited Applications

WB, ELISA, IF

Ideal working dilutions for each application should be empirically determined by the investigator.

Specificity

Recognizes the MYT1L protein

Cross Reactivity

Human, Mouse and Rat

Host/Isotype/Clone#

Rabbit, IgG

Immunogen

Rabbit polyclonal MYT1L antibody was raised against an 18 amino acid peptide near the carboxy terminus of human MYT1L

Formulation

PBS + 0.02% sodium azide

Stability

1yr at -20°C from date of shipment

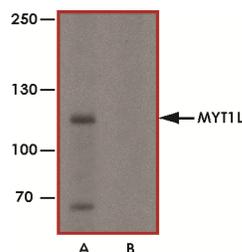
Scientific Background

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.

Sample Data



Western blot analysis of MYT1L in mouse brain tissue lysate with MYT1L antibody at 1 µg/ml in (A) the absence and (B) the presence of blocking peptide.

Anti-MYT1L

Rabbit Polyclonal Antibody

Catalog Number

M92-363R

Specific Lot Number

O2121-51

Purification

Affinity chromatography

Stability

1yr at -20°C from date of shipment

Storage & Shipping

Store product at -20°C. For optimal storage, aliquot antibody into smaller quantities after centrifugation and store at recommended temperature. For optimal performance, avoid repeated handling and multiple freeze/thaw cycles. Product shipped on ice packs.

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.