

MASH1 antibody (mAb)

Catalog Nos: 61271, 61272

RRID: AB_2793578

Clone: 24B7.2D11.1

Isotype: IgG1

Application(s): IHC, WB

Reactivity: Mouse, Rat

Quantities: 100 µg, 10 µg

Purification: Protein A Chromatography

Host: Mouse

Concentration: 1 µg/µl

Molecular Weight: 34 kDa

Background: The MASH1 (achaete-scute complex homolog 1 (Drosophila), ASCL1, ASH1, HASH1) gene encodes a protein that is a basic helix-loop-helix transcription factor that may play a role at early stages of development of specific neural lineages in most regions of the CNS, and of several lineages in the PNS. Essential for the generation of olfactory and autonomic neurons. Involved in the initiation of neuronal differentiation. Mediates transcriptional activation by binding to the E box motif.

Immunogen: This MASH1 antibody was raised against a full-length protein corresponding to rat MASH1.

Buffer: Purified IgG in PBS with 30% glycerol and 0.035% sodium azide. Sodium azide is highly toxic.

Application Notes:

Storage and Guarantee: Some products may be shipped at room temperature. This will not affect their stability or performance. Avoid repeated freeze/thaw cycles by aliquoting items into single-use fractions for storage at -20°C for up to 2 years. Keep all reagents on ice when not in storage. This product is guaranteed for 12 months from date of receipt.

This product is for research use only and is not for use in diagnostic procedures.