
MOBP Polyclonal Antibody

(Catalog # A73122)

Background

Predicted to enable actin binding activity and myosin binding activity. Predicted to be a structural constituent of myelin sheath. Predicted to be involved in nervous system development. Predicted to be located in mitochondrion. Predicted to be active in cortical actin cytoskeleton. Implicated in frontotemporal dementia.

Description

MOBP Polyclonal Antibody. Unconjugated. Raised in: Rabbit.

Formulation

Buffer: PBS with 0.01% thimerosal, 50% glycerol, pH7.3.

Specificity

Human, Mouse, Rat

Isotype

IgG

Uniprot ID

Q13875

Purification

Affinity Purification

Immunogen

Recombinant fusion protein containing a sequence corresponding to amino acids 1-70 of human MOBP (NP_891980.1).

Storage

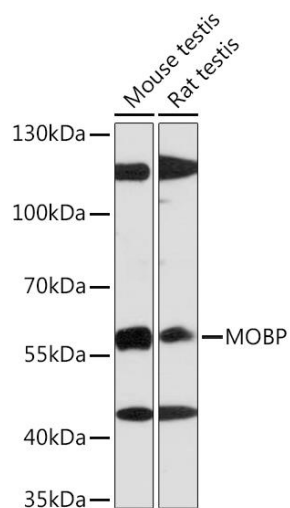
Shipped at 4°C. Upon receipt, store at -20°C. Avoid freeze / thaw cycles

Alternative Names

MOBP; myelin-associated oligodendrocyte basic protein

Application

WB; Recommended dilution: WB: 1:500-1:1000



Western blot analysis of extracts of various cell lines, using MOBP Polyclonal Antibody at 1:1000 dilution. Secondary antibody: HRP Goat Anti-Rabbit IgG (H+L) at 1:10000 dilution. Lysates/proteins: 25µg per lane. Blocking buffer: 3% nonfat dry milk in TBST.