
ZMYM3 Polyclonal Antibody

(Catalog # A72651)

Background

This gene is located on the X chromosome and is subject to X inactivation. It is highly conserved in vertebrates and most abundantly expressed in the brain. The encoded protein is a component of histone deacetylase-containing multiprotein complexes that function through modifying chromatin structure to keep genes silent. A chromosomal translocation (X;13) involving this gene is associated with X-linked mental retardation. Several alternatively spliced transcript variants have been found for this gene.

Description

ZMYM3 Polyclonal Antibody. Unconjugated. Raised in: Rabbit.

Formulation

Buffer: PBS with 0.02% sodium azide, 50% glycerol, pH7.3.

Specificity

Human

Isotype

IgG

Uniprot ID

Q14202

Purification

Affinity Purification

Immunogen

Recombinant fusion protein containing a sequence corresponding to amino acids 140-300 of human ZMYM3 (NP_001164634.1).

Storage

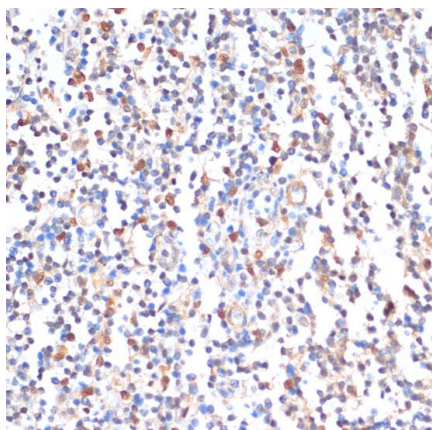
Shipped at 4°C. Upon receipt, store at -20°C. Avoid freeze / thaw cycles

Alternative Names

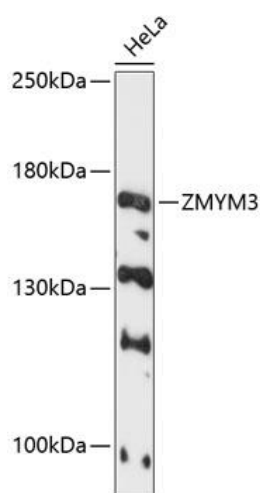
ZMYM3; DXS6673E; MYM; XFIM; ZNF198L2; ZNF261; zinc finger MYM-type protein 3

Application

WB, IHC; Recommended dilution: WB: 1:500-1:2000, IHC: 1:50-1:200



Immunohistochemistry of paraffin-embedded human tonsil using ZMYM3 Polyclonal Antibody at dilution of 1:100 (40x lens).



Western blot analysis of extracts of HeLa cells, using ZMYM3 Polyclonal Antibody at 1:3000 dilution.

Secondary Antibody: HRP Goat Anti-Rabbit IgG (H+L) at 1:10000 dilution.

Lysates/proteins: 25ug per lane.

Blocking buffer: 3% nonfat dry milk in TBST.

Exposure time: 90s.