
MBD3L1 Monoclonal Antibody [1C7]

[Catalog No. A-1106-100]

Background

MBD3L1, or methyl-CpG-binding domain protein 3-like 1, is a protein related to methyl-CpG-binding proteins but has not been shown to bind to methylated DNA because it lacks the methyl-CpG binding domain. Located in discrete areas within the nucleus, evidence suggests its expression is restricted to round spermatids. MBD3L1 is thought to play a role in male germ cell development during the postmeiotic stages.

Description

Mouse monoclonal antibody raised against a partial recombinant MBD3L1, clone 1C7.

Size

100 µg

Purification

Protein A purified

Immunogen

MBD3L1 (NP_660209, 1 a.a. ~ 110 a.a) partial recombinant protein with GST tag. MW of the GST tag alone is 26 KDa.

Cross Reactivity

Human

Isotype

IgG1 Kappa

Formulation

In 1x PBS, pH 7.2

Storage

Aliquot and store at -20°C or lower. Avoid repeated freezing and thawing. Multiple freeze/thaw cycles may result in decreased performance. Non-hazardous. No MSDS necessary.

Handling Recommendations

For maximum recovery of the products, centrifuge the vial prior to opening the cap.

Applications & Suggested Dilutions*

ELISA: 0.25-1 µg/ml

**The end user is responsible for determining optimal working dilutions.*

This product is for research purposes only. Not intended for use in diagnostic procedures.