

MBD5 Monoclonal Antibody [4A12-1B6]

[Catalog No. A-1105-100]

Background

MBD5, or methyl-CpG binding domain protein 5, is a part of the methyl-CpG-binding domain (MBD) family, which also includes MeCP2, MBD1, MBD2, MBD3, MBD4, CLLD8-like, and BAZ2A-like proteins. This protein contains a Pro-Trp-Trp-Pro motif (PWWP) domain as well as the MBD domain. The proteins of the PWWP domain play a role in cell division, cell growth, and cell differentiation. Evidence shows that diseases, such as mental retardation autosomal dominant type 1, can be caused by mutation to the MBD5 gene and some research indicates it may be involved in neurodevelopmental disorders and postnatal growth of mice.

Description

Mouse monoclonal antibody raised against a full length recombinant MBD5, clone 4A12-1B6.

Size

100 µg

Purification

Protein A purified

Immunogen

MBD5 (AAH14534, 1 a.a. ~ 291 a.a) full length 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.

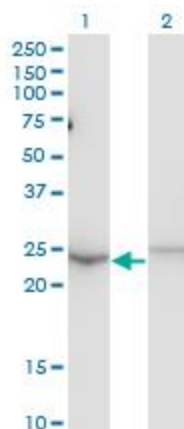
Applications & Suggested Dilutions*

Western Blot: 0.5-1 µg/ml

ELISA: 0.25-1 µg/ml

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

Data



Western Blot analysis of MBD5 expression in transfected 293T cell line by MBD5 Monoclonal Antibody, clone 4A12-1B6.

Lane 1: MBD5 transfected lysate (24.3 KDa).

Lane 2: Non-transfected lysate.

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