

GATAD2A Monoclonal Antibody [3F3]

[Catalog No. A-1109-100]

Background

GATAD2A (GATA zinc finger domain containing protein 2A), also known as Transcriptional repressor p66 alpha, is a transcriptional repressor that enhances repression mediated by MBD2 (methyl-CpG binding domain protein 2). For efficient repression by GATAD2A, the presence of GATAD2B is necessary. The transcriptional repressor p66-alpha protein binds MBD2 (methyl-CpG binding domain protein 2) and MBD3 (methyl-CpG binding domain protein 3) and requires interaction with MBD2 for targeting to the chromatin and for efficient repression. The protein also interacts with histones H2A, H2B, H3 and H4 and histone acetylation has been shown to reduce this interaction.

Description

Mouse monoclonal antibody raised against a partial recombinant GATAD2A, clone 3F3.

Size

100 µg

Purification

Protein A purified

Immunogen

Mouse monoclonal antibody raised against a partial recombinant GATAD2A.

Cross Reactivity

Human

Isotype

IgG2a 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.

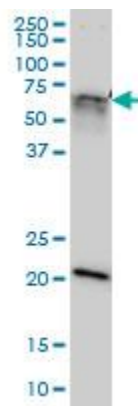
Applications & Suggested Dilutions*

Western Blot: 0.5-1 µg/ml

ELISA: 0.25-1 µg/ml

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

Data



GATAD2A Monoclonal Antibody, clone 3F3. Western Blot analysis of GATAD2A expression in HeLa NE.

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