
EIF2C4 Monoclonal Antibody [4G3]

[Catalog No. A-5102-100]

Background

EIF2C4 is a member of the Argonaute family of proteins that participate in RNA interference. This highly basic protein contains a PAS domain and a PIWI domain and is believed to play a role in short-interfering-RNA-mediated gene silencing. EIF2C4 is located on chromosome 1 in a cluster of similar family members including argonaute 3 and eukaryotic translation initiation factor 2C, 1.

Description

Mouse monoclonal antibody raised against a partial recombinant EIF2C4, clone 4G3.

Size

100 µg

Purification

Protein A purified

Immunogen

EIF2C4 (NP_060099, 1 a.a. ~ 101 a.a) partial recombinant protein with GST tag. MW of the GST tag alone is 26 KDa.

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*

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.