

FOXM1 Monoclonal Antibody [5G10]

[Catalog No. A-7179-100]

Description

Mouse monoclonal antibody raised against a partial recombinant FOXM1, clone 5G10.

Size

100 µg

Purification

Protein A purified

Immunogen

FOXM1 (NP_973731, 702 a.a. ~ 802 a.a) partial recombinant protein with GST tag. MW of the GST tag alone is 26 KDa.

Cross Reactivity

Human, Mouse, Rat

Isotype

IgG2b 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*

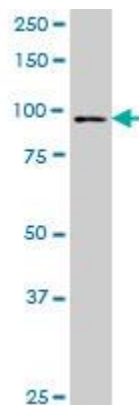
Western Blot: 0.5-1 µg/ml

ELISA: 0.25-1 µg/ml

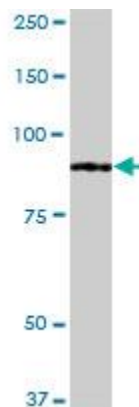
Immunofluorescence: 5 µg/ml

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

Data



FOXM1 Monoclonal Antibody, clone 5G10. Western Blot analysis of FOXM1 expression in PC-12.



FOXM1 Monoclonal Antibody, clone 5G10. Western Blot analysis of FOXM1 expression in NIH/3T3.

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