

RPS6KA3 Monoclonal Antibody [2G10]

[Catalog No. A-6148-100]

Description

Mouse monoclonal antibody raised against a partial recombinant RPS6KA3, clone 2G10.

Size

100 µg

Purification

Protein A purified

Immunogen

RPS6KA3 (NP_004577, 2 a.a. ~ 96 a.a) partial recombinant protein with GST tag. MW of the GST tag alone is 26 KDa.

Cross Reactivity

Human

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.

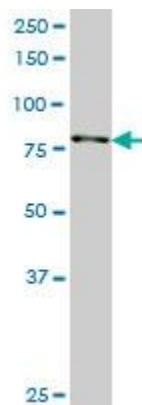
Applications & Suggested Dilutions*

Western Blot: 0.5-1 µg/ml

ELISA: 0.25-1 µg/ml

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

Data



RPS6KA3 Monoclonal Antibody, clone 2G10. Western Blot analysis of RPS6KA3 expression in HepG2.

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