

RPS6KA2 Monoclonal Antibody [1F6]

[Catalog No. A-6147-100]

Description

Mouse monoclonal antibody raised against a partial recombinant RPS6KA2, clone 1F6.

Size

100 µg

Purification

Protein A purified

Immunogen

RPS6KA2 (AAH02363, 631 a.a. ~ 733 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*

Western Blot: 0.5-1 µg/ml

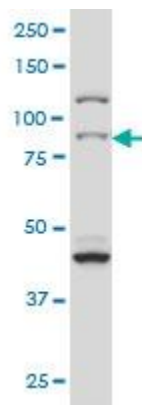
ELISA: 0.25-1 µg/ml

Immunohistochemistry (IHC): 2.5 µg/ml

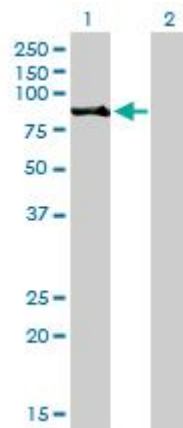
Immunofluorescence: 5 µg/ml

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

Data



RPS6KA2 Monoclonal Antibody, clone 1F6. Western Blot analysis of RPS6KA2 expression in HeLa NE.



Western Blot analysis of RPS6KA2 expression in transfected 293T cell line by RPS6KA2 Monoclonal Antibody, clone 1F6.

Lane 1: RPS6KA2 transfected lysate (83.2 KDa).
Lane 2: Non-transfected lysate.

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