

NEK10 Monoclonal Antibody [1C9]

[Catalog No. A-7141-100]

Description

Mouse monoclonal antibody raised against a partial recombinant NEK10, clone 1C9.

Size

100 µg

Purification

Protein A purified

Immunogen

NEK10 (NP_689747, 211 a.a. ~ 301 a.a) partial recombinant protein with GST tag. MW of the GST tag alone is 26 KDa.

Cross Reactivity

Human, Mouse

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*

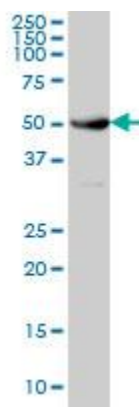
Western Blot: 0.5-1 µg/ml

ELISA: 0.25-1 µg/ml

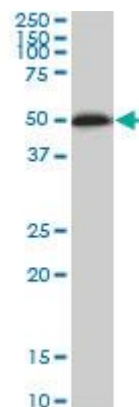
Immunofluorescence: 5 µg/ml

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

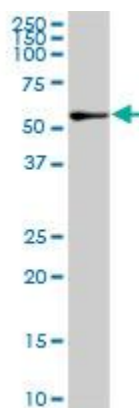
Data



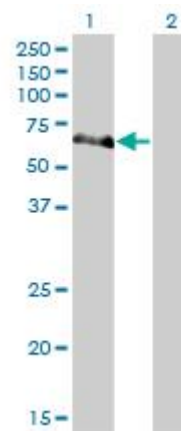
NEK10 Monoclonal Antibody, clone 1C9. Western Blot analysis of NEK10 expression in HeLa.



NEK10 Monoclonal Antibody, clone 1C9. Western Blot analysis of NEK10 expression in Raw 264.7.



NEK10 Monoclonal Antibody, clone 1C9. Western Blot analysis of NEK10 expression in NIH/3T3.



Western Blot analysis of NEK10 expression in transfected 293T cell line by NEK10 Monoclonal Antibody, clone 1C9.

Lane 1: NEK10 transfected lysate (53.1 KDa).
Lane 2: Non-transfected lysate.

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