

CDK4 Monoclonal Antibody [4F11]

[Catalog No. A-7153-100]

Description

Mouse monoclonal antibody raised against a partial recombinant CDK4, clone 4F11.

Size

100 µg

Purification

Protein A purified

Immunogen

CDK4 (AAH03644, 211 a.a. ~ 304 a.a) partial recombinant protein with GST tag. MW of the GST tag alone is 26 KDa.

Cross Reactivity

Human

Isotype

IgG1 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*

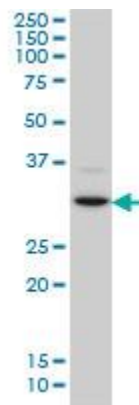
Western Blot: 0.5-1 µg/ml

ELISA: 0.25-1 µg/ml

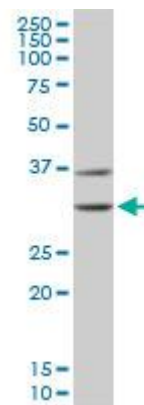
Immunofluorescence: 5 µg/ml

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

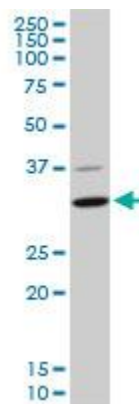
Data



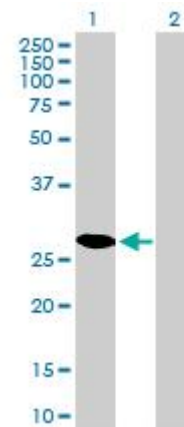
CDK4 Monoclonal Antibody, clone 4F11. Western Blot analysis of CDK4 expression in HeLa.



CDK4 Monoclonal Antibody, clone 4F11. Western Blot analysis of CDK4 expression in Raw 264.7.



CDK4 Monoclonal Antibody, clone 4F11. Western Blot analysis of CDK4 expression in NIH/3T3.



Western Blot analysis of CDK4 expression in transfected 293T cell line by CDK4 Monoclonal Antibody, clone 4F11.

Lane 1: CDK4 transfected lysate (33.7 KDa).
Lane 2: Non-transfected lysate.

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