

RNF168 Monoclonal Antibody [3E1]

[Catalog No. A-6143-100]

Description

Mouse monoclonal antibody raised against a partial recombinant RNF168, clone 3E1.

Size

100 µg

Purification

Protein A purified

Immunogen

RNF168 (NP_689830, 462 a.a. ~ 572 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

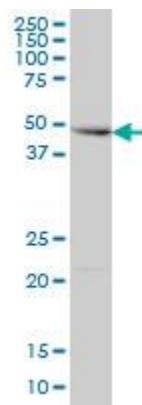
ELISA: 0.25-1 µg/ml

Immunohistochemistry (IHC): 2.5 µg/ml

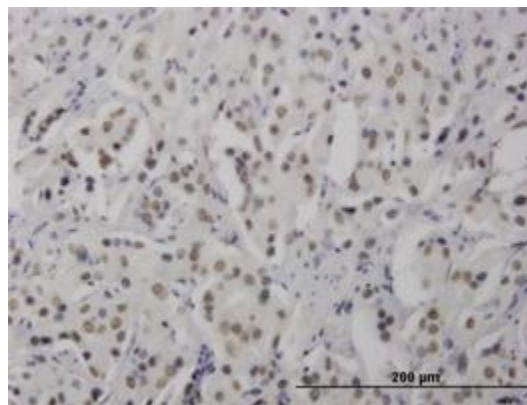
Immunofluorescence: 5 µg/ml

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

Data



RNF168 Monoclonal Antibody, clone 3E1. Western Blot analysis of RNF168 expression in HepG2.



Immunoperoxidase of monoclonal antibody to RNF168 on formalin-fixed paraffin-embedded human prostate cancer. [Antibody Concentration: 1.5 µg/ml]

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