

RNF138 Monoclonal Antibody [3G2]

[Catalog No. A-8133-200]

Description

Mouse monoclonal antibody raised against a partial recombinant RNF138, clone 3G2.

Size

200 µl

Purification

Protein A purified

Immunogen

RNF138 (-, 51 a.a. ~ 151 a.a) partial recombinant protein with GST tag. MW of the GST tag alone is 26 KDa.

Cross Reactivity

Human

Isotype

IgM Kappa

Formulation

In ascites fluid

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.

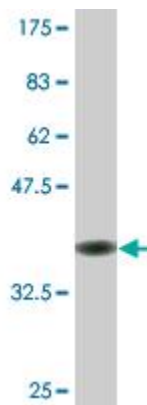
Applications & Suggested Dilutions*

Western Blot: 1: 500-1: 1000 dilution

ELISA: 1: 1000-1: 2000 dilution

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

Data



Western Blot detection against Immunogen (37.11 KDa).

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