

RAD17 Monoclonal Antibody [2G12]

[Catalog No. A-6134-100]

Description

Mouse monoclonal antibody raised against a partial recombinant RAD17, clone 2G12.

Size

100 µg

Purification

Protein A purified

Immunogen

RAD17 (AAH32304, 1 a.a. ~ 110 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.

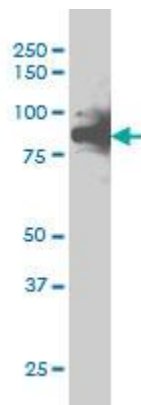
Applications & Suggested Dilutions*

Western Blot: 0.5-1 µg/ml

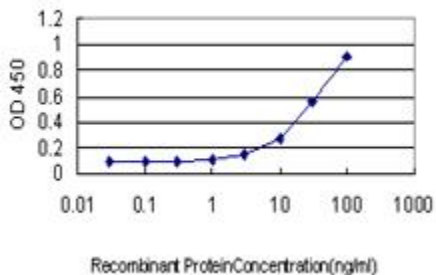
ELISA: 0.25-1 µg/ml

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

Data



RAD17 Monoclonal Antibody, clone 2G12. Western Blot analysis of RAD17 expression in HeLa NE.



Detection limit for recombinant GST tagged RAD17 is approximately 1 ng/ml as a capture antibody.

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