

RAD18 Monoclonal Antibody [3H7]

[Catalog No. A-6135-100]

Description

Mouse monoclonal antibody raised against a partial recombinant RAD18, clone 3H7.

Size

100 µg

Purification

Protein A purified

Immunogen

RAD18 (NP_064550, 332 a.a. ~ 431 a.a) partial recombinant protein with GST tag. MW of the GST tag alone is 26 KDa.

Cross Reactivity

Human, Rat

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

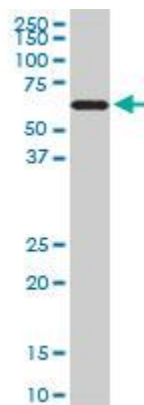
ELISA: 0.25-1 µg/ml

Immunohistochemistry (IHC): 2.5 µg/ml

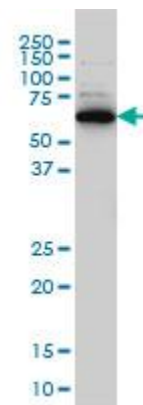
Immunofluorescence: 5 µg/ml

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

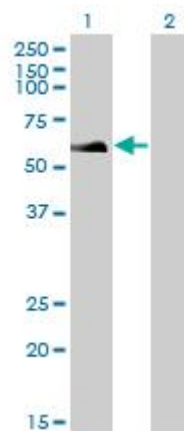
Data



RAD18 Monoclonal Antibody, clone 3H7. Western Blot analysis of RAD18 expression in PC-12.



RAD18 Monoclonal Antibody, clone 3H7. Western Blot analysis of RAD18 expression in HeLa NE.



Western Blot analysis of RAD18 expression in transfected 293T cell line by RAD18 monoclonal antibody, clone 3H7.

Lane 1: RAD18 transfected lysate (56.2 KDa).
Lane 2: Non-transfected lysate.

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