

## RAD23B Monoclonal Antibody [3H7]

[Catalog No. A-6136-100]

### Description

Mouse monoclonal antibody raised against a partial recombinant RAD23B, clone 3H7.

### Size

100 µg

### Purification

Protein A purified

### Immunogen

RAD23B (NP\_002865, 311 a.a. ~ 410 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\*

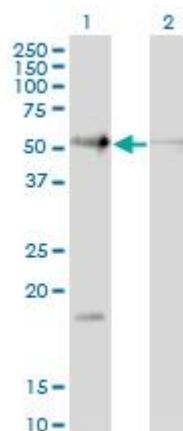
Western Blot: 0.5-1 µg/ml

ELISA: 0.25-1 µg/ml

Immunohistochemistry (IHC): 2.5 µg/ml

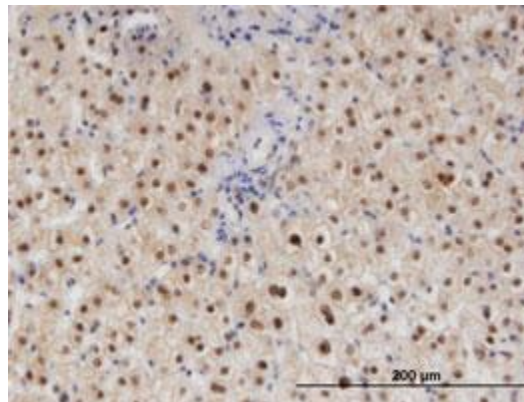
\*The end user is responsible for determining optimal working dilutions.

### Data



Western Blot analysis of RAD23B expression in transfected 293T cell line by RAD23B Monoclonal Antibody, clone 3H7.

Lane 1: RAD23B transfected lysate (43.2 KDa).  
Lane 2: Non-transfected lysate.



Immunoperoxidase of monoclonal antibody to RAD23B on formalin-fixed paraffin-embedded human liver. [Antibody Concentration: 3 µg/ml]

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