

## RPA3 Monoclonal Antibody [1F4]

[Catalog No. A-6188-100]

### Description

Mouse monoclonal antibody raised against a partial recombinant RPA3, clone 1F4.

### Size

100 µg

### Purification

Protein A purified

### Immunogen

RPA3 (AAH05264, 12 a.a. ~ 121 a.a) partial recombinant protein with GST tag. MW of the GST tag alone is 26 KDa.

### Cross Reactivity

Human

### Isotype

IgG1 Lambda

### 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\*

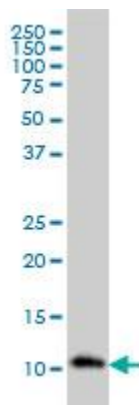
Western Blot: 0.5-1 µg/ml

ELISA: 0.25-1 µg/ml

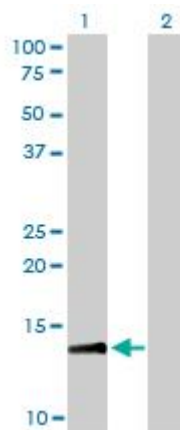
Immunohistochemistry (IHC): 2.5 µg/ml

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

### Data



RPA3 Monoclonal Antibody, clone 1F4. Western Blot analysis of RPA3 expression in HL-60.



Western Blot analysis of RPA3 expression in transfected 293T cell line by RPA3 Monoclonal Antibody, clone 1F4.

Lane 1: RPA3 transfected lysate (13.6 KDa).  
Lane 2: Non-transfected lysate.

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