

## EGF Monoclonal Antibody [2F1]

[Catalog No. A-7121-100]

### Description

Mouse monoclonal antibody raised against a partial recombinant EGF, clone 2F1.

### Size

100 µg

### Purification

Protein A purified

### Immunogen

EGF (NP\_001954, 926 a.a. ~ 1025 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.

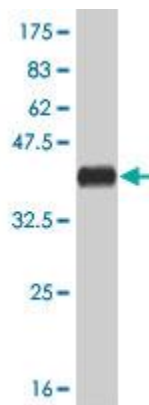
### Applications & Suggested Dilutions\*

Western Blot: 0.5-1 µg/ml

ELISA: 0.25-1 µg/ml

*\*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.*