

## TIPIN Monoclonal Antibody [1G11]

[Catalog No. A-6144-100]

### Description

Mouse monoclonal antibody raised against a full length recombinant TIPIN, clone 1G11.

### Size

100 µg

### Purification

Protein A purified

### Immunogen

TIPIN (AAH00870, 1 a.a. ~ 302 a.a) full length 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\*

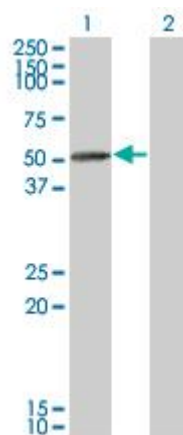
Western Blot: 0.5-1 µg/ml

ELISA: 0.25-1 µg/ml

Immunofluorescence: 5 µg/ml

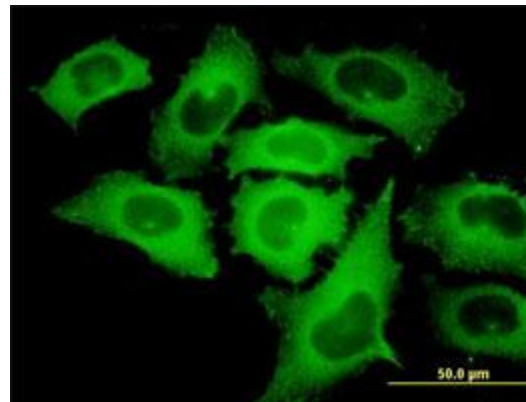
\*The end user is responsible for determining optimal working dilutions.

### Data



Western Blot analysis of TIPIN expression in transfected 293T cell line by TIPIN Monoclonal Antibody, clone 1G11.

Lane 1: TIPIN transfected lysate (34.496 KDa).  
Lane 2: Non-transfected lysate.



Immunofluorescence of monoclonal antibody to TIPIN on HeLa cell. [Antibody Concentration: 10 µg/ml]

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