

# SMURF1 Monoclonal Antibody [1D7]

[Catalog No. A-8131-100]

## Description

Mouse monoclonal antibody raised against a partial recombinant SMURF1, clone 1D7.

## Size

100 µg

## Purification

Protein A purified

## Immunogen

SMURF1 (NP\_065162, 165 a.a. ~ 269 a.a) partial recombinant protein with GST tag. MW of the GST tag alone is 26 KDa.

## Cross Reactivity

Human, Mouse

## Isotype

IgG2a 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

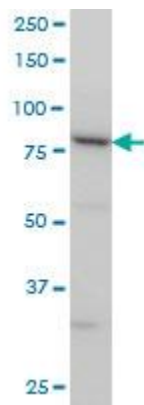
ELISA: 0.25-1 µg/ml

Immunohistochemistry (IHC): 2.5 µg/ml

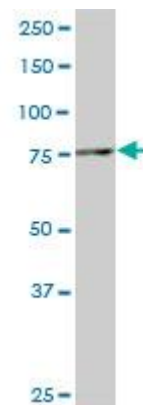
Immunofluorescence: 5 µg/ml

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

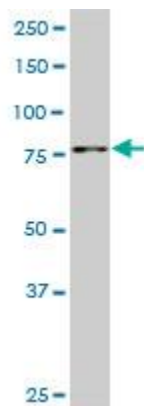
## Data



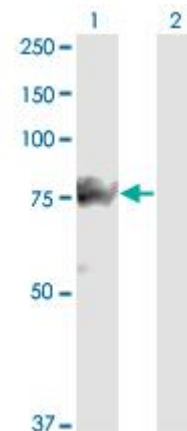
SMURF1 Monoclonal Antibody, clone 1D7. Western Blot analysis of SMURF1 expression in HeLa.



SMURF1 Monoclonal Antibody, clone 1D7. Western Blot analysis of SMURF1 expression in Raw 264.7.



SMURF1 Monoclonal Antibody, clone 1D7. Western Blot analysis of SMURF1 expression in NIH/3T3.



Western Blot analysis of SMURF1 expression in transfected 293T cell line by SMURF1 Monoclonal Antibody, clone 1D7.

Lane 1: SMURF1 transfected lysate (83.4 KDa).

Lane 2: Non-transfected lysate.

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