

---

## B-actin Monoclonal Antibody [A2-F6]

(Catalog No. A-0800)

---

### Background

Beta actin, or cytoplasmic actin, is a 43kDa highly conserved protein that is ubiquitously expressed in all eukaryotic cells. Actins, along with microtubules, comprise a major component of the cytoskeleton. B-Actin is a relatively stable cytoskeletal protein normally at a constant level in cells, regardless of experimental treatment or technical procedure. For this reason, measurement of Beta-actin is generally used as an internal control for experimental error.

### Description

Mouse monoclonal antibody to B-actin, clone A2-F6

### Specificity

Human, Mouse, Rat

### Isotype

IgG1

### Uniprot ID

P60709

### Purification

Affinity Purified

### Immunogen

Synthetic peptide (KLH-coupled)

### Formulation

Liquid. 1\*TBS (pH7.4), 0.5% BSA, 40% Glycerol. Preservative: 0.05% Sodium Azide.

### Storage

Store at 4°C after thawing. Aliquot store at -20°C (regular) or -80°C (long-term). Avoid repeated freeze / thaw cycles.

### Application

WB: 1:10,000

ICC: 1:100-1:200

IHC: 1:100-1:200

FC: 1:100-1:200

### Ordering Information

#### Products

B-actin Monoclonal Antibody [A2-F6]

#### Size

50 µg

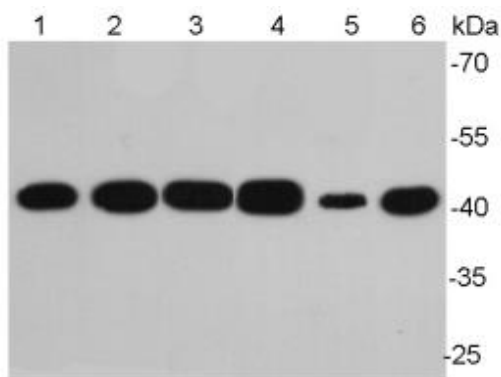
100 µg

#### Cat. No.

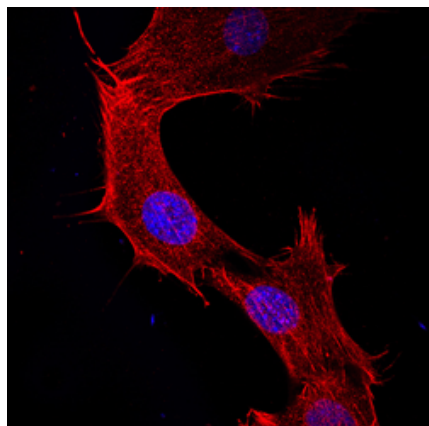
A-0800-050

A-0800-100

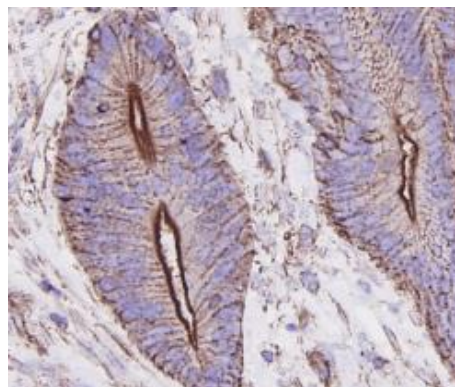
*This product is for research purposes only. Not intended for use in diagnostic procedures.*



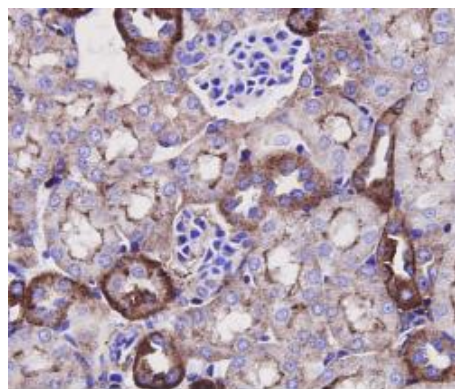
▲ Western blot analysis of Beta-actin on different cell lysates using B-actin Monoclonal Antibody [A2-F6] at 1/10000 dilution.  
Positive control:  
Lane 1: NIH/3T3  
Lane 2: PC12  
Lane 3: MCF-7  
Lane 4: HepG2  
Lane 5: Hela  
Lane 6: Mouse lung



▲ Immunofluorescent staining of NIH/3T3 cells using B-actin Monoclonal Antibody [A2-F6].



▲ Immunohistochemical analysis of paraffin-embedded human colon carcinoma tissue using B-actin Monoclonal Antibody [A2-F6]. Counter stained with hematoxylin.



▲ Immunohistochemical analysis of paraffin-embedded mouse kidney tissue using B-actin Monoclonal Antibody [A2-F6]. Counter stained with hematoxylin.

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