

## Anti-ADD1

Rabbit Polyclonal Antibody

**Catalog # A38-63R**

Lot # R027-4

### Cited Applications

- Western blot (1:1000 dilution)

*Ideal working dilutions for each application should be empirically determined by the investigator.*

### Specificity

Recognizes the ADD1 protein

### Cross Reactivity

- Western blot ADD1 from human and mouse cells

*ADD1 from other species may also be detectable*

### Host

Rabbit

### Immunogen

Synthetic peptide corresponding to amino acids 724-728 of human ADD1 (P35611)

### Formulation

PBS pH 7.4, 0.02% sodium azide, 50% glycerol

### Stability

Store at 4°C (add 0.1% NaN<sub>3</sub>) for several months, and at -20°C for longer periods. For optimal storage, aliquot antibody into smaller quantities after centrifugation and store at recommended temperature. For optimal performance, avoid repeated handling and multiple freeze/thaw cycles.

### Scientific Background

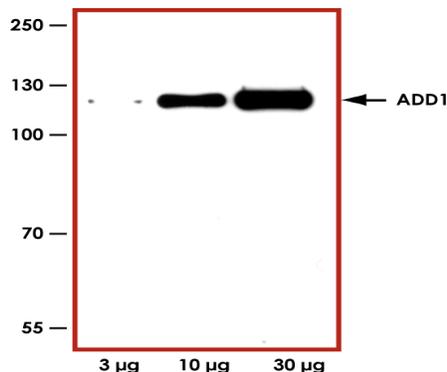
ADD1 or Adducin 1 is a heterodimeric protein associated with the erythrocyte membrane skeleton, which promotes binding of spectrin to actin, and is a substrate for PKC and PKA (1). Alpha- and gamma-adducins are ubiquitously expressed while beta-adducin is expressed at high levels in brain and hematopoietic tissues. ADD1 can completely block elongation and depolymerization at the fast-growing barbed ends of actin filaments in vitro and this suggests that adducin restrict actin filament length in erythrocytes (2).

### References

1. Gardner, K. et.al: A new erythrocyte membrane-associated protein with calmodulin binding activity: identification and purification. *J. Biol. Chem.* 261: 1339-1348, 1986.
2. Kuhlman, et.al: A new functions for adducin: calcium/calmodulin-regulated capping of the barbed ends of actin filaments. *J. Biol. Chem.* 271: 7986-7991, 1996.

To place your order, please contact us by phone 1-(604)-232-4600, fax 1-604-232-4601 or by email: [orders@signalchem.com](mailto:orders@signalchem.com)  
[www.signalchem.com](http://www.signalchem.com)

### Sample Data



Representative western blot with Anti-ADD1 (1:1000) using 3 µg, 10 µg and 30 µg of HEK-293T cell lysate.

## Anti-ADD1

Rabbit Polyclonal Antibody

Catalog Number

A38-63R

Specific Lot Number

R027-4

Purification

Affinity chromatography

Concentration

1.0 µg/µL

Stability

1yr at -20°C from date of shipment

Storage & Shipping

Store product at -20°C. For optimal storage, aliquot antibody into smaller quantities after centrifugation and store at recommended temperature. For optimal performance, avoid repeated handling and multiple freeze/thaw cycles. Product shipped on ice packs.