

---

## ARPC2 Polyclonal Antibody, Biotin Conjugated

(Catalog # A51680)

---

### Background

ARPC2 is one of seven subunits of the human Arp2/3 protein complex. The Arp2/3 protein complex has been implicated in the control of actin polymerization in cells and has been conserved through evolution. The exact role of ARPC2, the p34 subunit, has yet to be determined. Two alternatively spliced variants have been characterized to date. Additional alternatively spliced variants have been described but their full length nature has not been determined.

### Description

ARPC2 Polyclonal Antibody, Biotin Conjugated. Biotin. Raised in: Rabbit.

### Formulation

Liquid. 0.03% Proclin 300, 50% Glycerol, 0.01M PBS, PH 7.4.

### Specificity

Human

### Isotype

IgG

### Uniprot ID

O15144

### Purification

>95%, Protein G purified

### Immunogen

Recombinant Human Actin-related protein 2/3 complex subunit 2 protein (1-250AA)

### Storage

Shipped at 4°C. Upon delivery aliquot and store at -20°C (short-term) or -80°C (long-term). Avoid repeated freeze.

### Application

ELISA; Recommended dilution: ELISA 1:500-1:1000

---