

---

## PLEC Polyclonal Antibody, FITC Conjugated

(Catalog # A69885)

---

### Background

Interlinks intermediate filaments with microtubules and microfilaments and anchors intermediate filaments to desmosomes or hemidesmosomes. Could also bind muscle proteins such as actin to membrane complexes in muscle. May be involved not only in the filaments network, but also in the regulation of their dynamics. Structural component of muscle. Isoform 9 plays a major role in the maintenance of myofiber integrity.

### Description

PLEC Polyclonal Antibody, FITC Conjugated. FITC. Raised in: Rabbit.

### Formulation

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

### Specificity

Human

### Isotype

IgG

### Uniprot ID

Q15149

### Purification

>95%, Protein G purified

### Immunogen

Recombinant Human Plectin protein (1744-1929AA)

### Storage

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

### Alternative Names

Plectin, PCN, PLTN, Hemidesmosomal protein 1, HD1, Plectin-1, PLEC, PLEC1

### Application

ELISA; Recommended dilution: ELISA 1:100-1:500

---