

---

## PTPN3 Polyclonal Antibody, FITC Conjugated

(Catalog #A69669)

---

### Background

May act at junctions between the membrane and the cytoskeleton. Possesses tyrosine phosphatase activity.

### Description

PTPN3 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

P26045

### Purification

>95%, Protein G purified

### Immunogen

Recombinant Human Tyrosine-protein phosphatase non-receptor type 3 protein (365-505AA)

### Storage

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

### Alternative Names

Tyrosine-protein phosphatase non-receptor type 3, Protein-tyrosine phosphatase H1, PTP-H1, PTPN3, PTPH1

### Application

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

---