

---

## PTPRZ1 Polyclonal Antibody, FITC Conjugated

(Catalog # A56577)

---

### Background

Protein tyrosine phosphatase that negatively regulates oligodendrocyte precursor proliferation in the embryonic spinal cord. Required for normal differentiation of the precursor cells into mature, fully myelinating oligodendrocytes. May play a role in protecting oligodendrocytes against apoptosis. May play a role in the establishment of contextual memory, probably via the dephosphorylation of proteins that are part of important signaling cascades.

### Description

PTPRZ1 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

P23471

### Purification

>95%, Protein G purified

### Immunogen

Recombinant Human Receptor-type tyrosine-protein phosphatase zeta protein (36-300AA)

### Storage

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

### Alternative Names

Protein-tyrosine phosphatase receptor type Z polypeptide 1, Protein-tyrosine phosphatase receptor type Z polypeptide 2, R-PTP-zeta-2, HTPZP2, PTPRZ, PTPRZ2, PTPZ, PTPRZ1

### Application

ELISA (1:100-1:500)

---