
PRY Polyclonal Antibody, Biotin Conjugated

(Catalog #A63241)

Description

PRY 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

O14603

Purification

>95%, Protein G purified

Immunogen

Recombinant Human PTPN13-like protein, Y-linked protein (13-147AA)

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

PTPN13-like protein, Y-linked, Testis-specific PTP-BL-related Y protein, PRY, PRY1, PTPN13LY, AND, PRY2, PTPN13LY2, AND, PRYP3, AND, PRYP4

Application

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