

PTPRR Polyclonal Antibody

(Catalog # A71545)

Background

The protein encoded by this gene is a member of the protein tyrosine phosphatase (PTP) family. PTPs are known to be signaling molecules that regulate a variety of cellular processes including cell growth, differentiation, mitotic cycle, and oncogenic transformation. This PTP possesses an extracellular region, a single transmembrane region, and a single intracellular catalytic domain, and thus represents a receptor-type PTP. Silencing of this gene has been associated with colorectal cancer. Multiple transcript variants encoding different isoforms have been found for this gene. This gene shares a symbol (PTPRQ) with another gene, protein tyrosine phosphatase, receptor type, Q (GeneID 374462), which is also located on chromosome 12.

Description

PTPRR Polyclonal Antibody. Unconjugated. Raised in: Rabbit.

Formulation

Buffer: PBS with 0.02% sodium azide, 50% glycerol, pH7.3.

Specificity

Human, Mouse, Rat

Isotype

IgG

Uniprot ID

Q15256

Purification

Affinity Purification

Immunogen

Recombinant fusion protein containing a sequence corresponding to amino acids 240-400 of human PTPRR (NP_002840.2).

Storage

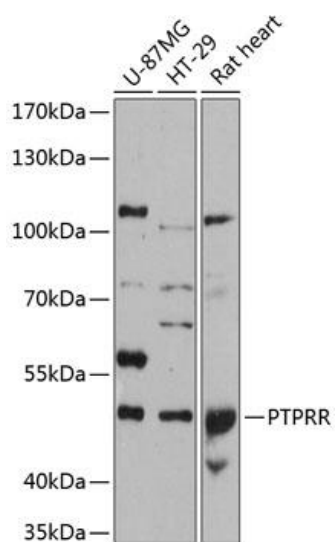
Shipped at 4°C. Upon receipt, store at -20°C. Avoid freeze / thaw cycles

Alternative Names

PTPRR; EC-PTP; PCPTP1; PTP-SL; PTPBR7; PTPRQ; receptor-type tyrosine-protein phosphatase R

Application

WB; Recommended dilution: WB: 1:500-1:2000



Western blot analysis of extracts of various cell lines, using PTPRR Polyclonal Antibody at 1:3000 dilution.
Secondary Antibody: HRP Goat Anti-Rabbit IgG (H+L) at 1:10000 dilution.
Lysates/proteins: 25ug per lane.
Blocking buffer: 3% nonfat dry milk in TBST.
Exposure time: 90s.