

## PTPRM (PTP-µ) Protein

Recombinant protein expressed in Sf9 cells

**Catalog # P58-35G**

Lot # C234-1

### Product Description

Recombinant human PTPRM (PTP-µ) (764-1452) was expressed by baculovirus in Sf9 insect cells using an N-terminal GST tag. The gene accession number is [BC051651](#).

### Gene Aliases

RPTPM, RPTPU, PTPRL1, hR-PTP $\mu$ , R-PTP-MU

### Formulation

Recombinant protein stored in 50mM Tris-HCl, pH 7.5, 150mM NaCl, 10mM glutathione, 0.1mM EDTA, 0.25mM DTT, 0.1mM PMSF, 25% glycerol.

### Storage and Stability

Store product at  $-70^{\circ}\text{C}$ . For optimal storage, aliquot target into smaller quantities after centrifugation and store at recommended temperature. For most favorable performance, avoid repeated handling and multiple freeze/thaw cycles.

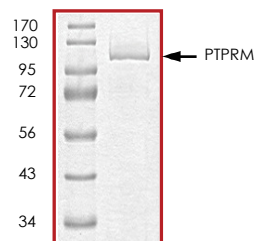
### Scientific Background

PTPRM is a member of the protein tyrosine phosphatase family and can participate in a variety of cellular processes including cell growth, differentiation, mitotic cycle, and oncogenic transformation (1). PTPRM has been shown to mediate cell-cell aggregation through the interaction with another molecule of PTPRM on an adjacent cell. PTPRM can interact with scaffolding protein RACK1/GNB2L1 and this interaction may be necessary for downstream signaling in response to cell-cell adhesion (2). PTPRM has been shown to be expressed in human pulmonary vascular endothelia where it directly binds to VE-cadherin and regulates both the tyrosine phosphorylation state of VE-cadherin and barrier integrity.

### References

- Phillips-Mason, P.J. et al: The receptor protein-tyrosine phosphatase PTPmu interacts with IQGAP1. *J. Biol. Chem.* 2006; 281(8):4903-10.
- Koop, E.A. et al: Receptor protein tyrosine phosphatase mu expression as a marker for endothelial cell heterogeneity; analysis of RPTPmu gene expression using LacZ knock-in mice. *Int. J. Dev. Biol.* 2003; 47(5):345-54

### Purity



The purity of PTPRM was determined to be **>90%** by densitometry. Approx. MW **110kDa**.

## PTPRM (PTP-µ) Protein

Recombinant protein expressed in Sf9 cells

Catalog Number	P58-35G
Specific Lot Number	C234-1
Purity	>90%
Concentration	0.2 µg/µl
Stability	1 yr at $-70^{\circ}\text{C}$ from date of shipment
Storage & Shipping	Store product at $-70^{\circ}\text{C}$ . For optimal storage, aliquot target into smaller quantities after centrifugation and store at recommended temperature. For most favorable performance, avoid repeated handling and multiple freeze/thaw cycles. Product shipped on dry ice.

To place your order, please contact us by phone 1-(604)-232-4600, fax 1-604-232-4601 or by email: [orders@signalchem.com](mailto:orders@signalchem.com)  
[www.signalchem.com](http://www.signalchem.com)

**FOR IN VITRO RESEARCH PURPOSES ONLY. NOT INTENDED FOR USE IN HUMAN OR ANIMALS.**