

PTP μ , cytoplasmic region (D1D2), Protein tyrosine phosphatase human, recombinant, *E. coli*

Cat. No.	Amount
PR-349	20 μ g

For *in vitro* use only
Quality guaranteed for 12 months
Store at -80°C

Avoid freeze / thaw cycles

Form

Liquid. Supplied as 0.4 mg/ml solution in 10 mM Tris-HCl, 10 mM Bis-Tris-HCl, pH 7.0.

Activity

1 nmol of PTP μ will hydrolyze 1.2 nmol of pNPP per minute at 37°C at pH 4.8.

Molecular weight

72.95 kDa.

Purity

>95% by SDS-PAGE.

Description

Protein Tyrosin Phosphatase, cytoplasmic region, containing catalytic domain D1 C-terminally linked to the non-catalytic domain D2.

PTP μ is an Ig superfamily receptor protein-tyrosine phosphatase which promotes cell-cell adhesion in a homophilic, Ca²⁺-independent manner and interacts with the cadherin-catenin complex. The intracellular region of PTP μ contains three domains: the juxtamembrane (JM, aa 765-923) and two phosphatase domains (D1 and D2). Only the D1 domain is catalitically active.

Selected References:

- Gebbink *et al.* (1993) Cell-cell adhesion mediated by a receptor-like protein tyrosine phosphatase. *J. Biol. Chem.* **268**:16101.
Brady-Kalnay *et al.* (1998) Dynamic interaction of PTP μ with multiple cadherins *in vivo*. *J. Cell Biol.* **141**:187.
Tonks *et al.* (1992) Protein tyrosine phosphatases: The problems of a growing family. *Cold Spring Harbor Symp. Quant. Biol.* **57**:87.