

PRL-3

human, recombinant, *E. coli*

Cat. No.	Amount
PR-330	20 µg

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

Avoid freeze / thaw cycles

Form

Liquid. Supplied in 50 mM Tris-HCl pH 7.5, 150 mM NaCl and 1 mM DTT.

Purity

>95% by SDS-PAGE

Description

Protein Tyrosine Phosphatase type IVA, member 3, isoform 1; whole sequence.

The proteins of regenerating liver (PRL-1, -2, and -3) are closely related protein tyrosine phosphatases (PTPs) having high sequence similarities to several dual-specificity PTPs. Elevated expression levels of PRL-3 have been reported in colon cancer cells that have metastasized to liver.

Selected References:

Matter *et al.* (2001) Role of PRL-3, a human muscle-specific tyrosine phosphatase, in angiotensin-II signaling. *Biochem. Biophys. Res. Commun.* **283**:1061.
Saha *et al.* (2001) A phosphatase associated with metastasis of colorectal cancer. *Science* **294**:1343.