

**PIM1, active, dephosphorylated**  
**PIM1<sup>His</sup>, active, dephosphorylated**  
**human, recombinant, *E. coli***

Cat. No.	Amount
PR-237	20 µg

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

**Avoid freeze / thaw cycles**

**Description**

Human PIM1 plays multiple roles in tumorigenesis. It promotes early transformation, cell proliferation and cell survival. In addition it may play a role in angiogenesis and vasculogenesis as a downstream effector of the VEGF-A/Flk1 pathway. PIM1 expression is correlated with tumor aggressiveness and is a marker for poor prognosis. PIM1 expression can be predictive of tumor outcome following chemotherapy and surgery and has been correlated with the enhanced metastatic potential of the tumor.

**Form**

Supplied in 50 mM Tris-HCl pH 7.5, 500 mM NaCl and 1 mM DTT.

**Activity**

>500,000 U/mg (1 Unit is defined as 1 picomole phosphate transferred to synthetic peptide KKRNRTLTV per min at 30 °C)

**Molecular Weight**

37.5 kDa

**Application**

Suitable for crystallization because it is purified by Ni-NTA agarose chromatography, ion exchange chromatography and gel filtration.

**Selected References:**

Shah *et al.* (2008) Potential roles for the PIM1 kinase in human cancer – A molecular and therapeutic appraisal. *Eur. J. Cancer.* **44**:2144.  
Bullok *et al.* (2005) Structure and substrate specificity of the Pim1 kinase. *J. Biol. Chem.* **280**:41675.