

PRODUCT DATA SHEET



DEP-1 protein tyrosine phosphatase (density-enhanced phosphatase-1, PTP-eta, CD148), catalytic domain (human, recombinant, N-terminal GST tag)

CATALOG NO.: SE-285

LOT NO.: temp

DESCRIPTION: catalytic domain, (a.a. 1024-1306) with N-terminal GST tag, MW=32.9 kDa, expressed in an *E. coli* expression system.

PURITY: >90% by SDS-PAGE.

SPECIFIC ACTIVITY: 114.0 U/μg. One unit will hydrolyze 1 nmol p-nitrophenyl phosphate per minute at pH 7.4 and 30°C. Assay buffer: 50 mM HEPES, pH 7.4, 100 mM NaCl, 2 mM EDTA, 3 mM dithiothreitol, 50 mM pNPP

USE : Remove phosphate specifically from tyrosine residues in proteins. Rate of dephosphorylation depends on DEP-1 substrate preference.

SUPPLIED AS: 1 mg/ml in 25 mM Tris-HCl, pH 8.0, 75 mM NaCl, 0.05% Tween-20, 50% glycerol, 2 mM EDTA, 1 mM dithiothreitol, 10 mM glutathione.

STORAGE: -20°C.

NOTE: When stored under the above conditions, this enzyme is stable at the concentration supplied, in its current storage buffer. Procedures such as dilution of the enzyme followed by refreezing, could lead to loss of activity.

REFERENCES:

1. A. Ostman, *et al. Proc. Natl. Acad. Sci (U.S.A.)* 1994 **91** 9680
2. L. Borges, *et al. Circ. Res.* 1996 **79** 570
3. M. de la Fuente-Garcia, *et al. Blood* 1998 **91** 2800
4. M. Kovalenko, *et al. J. Biol. Chem.* 2000 **275** 16219

FOR RESEARCH USE ONLY. NOT FOR DIAGNOSTIC OR THERAPEUTIC USE.