



Echelon Biosciences Inc.  
675 Arapeen Drive, Suite 302  
Salt Lake City, UT 84108  
Telephone 866-588-0455  
Fax 801-588-0497  
echelon@echelon-inc.com  
www.echelon-inc.com

## Technical Data Sheet

**For research use only**  
Not intended or approved for  
diagnostic or therapeutic use.

### **DEP1/CD148/PTPRJ, active**

(Recombinant human protein expressed in *E. coli*)

Catalog # **E-P017**  
Size: 20 µg  
Specific activity: 114,000 U/mg

#### **Product Description**

DEP1/PTP-eta/CD148/PTPRJ (density-enhanced phosphatase-1) is protein tyrosine phosphatase, receptor type, J. Recombinant human catalytic domain (amino acids 1024-1306) of DEP1 with N-terminal GST tag was expressed in an *E. coli* expression system. The GenBank Accession number is U10886. MW = 59 kDa.

#### Specific Activity

114,000 U/mg. 1 unit of DEP1 activity is equal to 1 nmol phosphate hydrolyzed from p-nitrophenyl phosphate substrate per minute at pH 7.4 and 30 °C.

#### Assay Condition

50 mM HEPES, pH 7.4, 2 mM EDTA, 3 mM DTT, 100 mM NaCl, 50 mM pNPP substrate, and 0.5 µg/ml recombinant DEP1.

#### Purity

> 95% by SDS-PAGE.

#### Formulation

Recombinant protein is in storage buffer (25 mM Tris-HCl, pH 8.0, 75 mM NaCl, 0.05% Tween-20, 2 mM EDTA, 3 mM DTT, and 50% glycerol).

#### Storage and Stability

Store product frozen at or below -70 °C. Enzyme will be stable for at least 6 months at -70 °C as undiluted stock. Freeze in working aliquots to avoid repeated thawing and freezing.

#### **Application**

DEP1 is useful for the study of enzyme kinetics, regulation, target substrate dephosphorylation, and inhibitor screening.

#### **References**

1. Massa, A. *et al.* (2004) *J. Biol. Chem.* **279**: 29004-12.
2. Kellie, S. *et al.* (2004) *J. Cell Sci.* **117**: 609-18.