



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.

MEG-2/PTPN9, active

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

Catalog # **E-P008**
Size: 20 µg
Specific activity: 2,000 U/mg

Product Description

MEG-2/PTP-MEG2/PTPN9 is protein tyrosine phosphatase, non-receptor type 9. Recombinant human catalytic domain (amino acids 285-593) of MEG-2 with N-terminal GST tag was expressed in an *E. coli* expression system. The GenBank Accession number is M83738. MW = 60 kDa.

Specific Activity

2,000 U/mg. 1 unit of MEG-2 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 10 µg/ml recombinant MEG-2.

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, 1 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

MEG-2 is useful for the study of enzyme kinetics, regulation, target substrate dephosphorylation, and inhibitor screening.

References

1. Qi, Y. *et al.* (2002) *J. Cell Biochem.* **86**: 79-89
2. Huynh, H. *et al.* (2003) *J. Immunol.* **171**: 6661-71