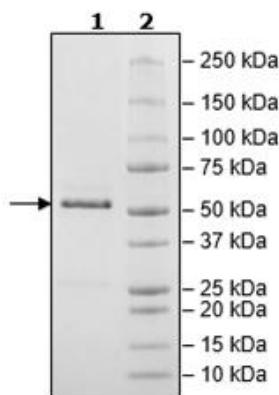


Product Information

Construct:	MEG-1 (GST-637-926(end))
Concentration:	0.32 mg/ml
Species:	Human
Formulated In:	40 mM Tris-HCl pH 8.0, 110 mM NaCl, 2.2 mM KCl, 20% glycerol, 3 mM DTT, 16 mM glutathione
Expression System:	<i>E. coli</i>
Format:	Aqueous buffer solution
Stability:	At least 6 months at -80°C. Avoid freeze/thaw cycles.
Storage:	-80°C
Genbank Accession:	NM_002830
MW:	60 kDa
Purity:	≥90%
Assay Conditions:	50 mM HEPES, pH 7.4, 2 mM EDTA, 3 mM DTT, 100 mM NaCl, 10 μM DIFMUP. Incubated for 5 minutes at room temperature with various amounts of MEG-1.
Applications:	Useful for the study of enzyme kinetics, screening inhibitors, and selectivity profiling.

Quality Control Data

4-20% SDS-Page Coomassie Staining



MEG-1 Activity

