

Product Information

Construct:	MGMT (His-2-207(end))
Concentration:	2.13 mg/ml
Species:	Human
Formulated In:	40 mM Tris-HCl, pH 8.0, 110 mM NaCl, 2.2 mM KCl, 0.04% Tween-20, 80 mM Imidazole, and 20% glycerol
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_002412
MW:	23 kDa
Purity:	85%

Quality Control Data

4-20% SDS-PAGE Coomassie Staining

