

Product Information

Construct:	MAOA (His-FLAG-2-497)
Concentration:	0.15 mg/ml
Species:	Human
Formulated In:	40 mM Tris-HCl, pH 8.0, 110 mM NaCl, 2.2 mM KCl, 0.04% Tween-20, 20% glycerol, and 100 µg/ml FLAG Peptide
Expression System:	Sf9
Format:	Aqueous buffer solution.
Stability:	At least 6 months at -80°C. Avoid freeze/thaw cycles.
Storage:	-80°C
Genbank Accession:	NM_000240
MW:	58 kDa
Purity:	≥90%

Quality Control Data

4-20% SDS-PAGE Coomassie Staining

