

Product Information

Construct:	ALOX15 (FLAG-2-662(end))
Concentration:	0.36 mg/ml
Species:	Human
Formulated In:	50 mM Tris-HCl, pH 8.0, 138 mM NaCl, 2.7 mM KCl, 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_001140
MW:	76 kDa
Purity:	≥90%

Quality Control Data

4-20% SDS-Page Coomassie Staining

