

Product Information

Description:	Recombinant human HER3 (Receptor tyrosine-protein kinase erbB-3), encompassing amino acids (667-1020) length. This construct contains an N-terminal His-tag. The recombinant protein was affinity purified.
Species:	Human
Construct:	HER3 (His-667-1020)
Concentration:	0.20 mg/ml
Expression System:	Sf9
Purity:	90%
Format:	Aqueous buffer solution.
Formulated In:	50 mM Sodium Phosphate, pH 7.5, 300 mM NaCl, 150 mM Imidazole, 0.25 mM DTT, 25% glycerol
MW:	46 kDa
Uniprot:	P21860
Stability:	At least 6 months at -80°C.
Storage:	-80°C
Instructions for Use:	Thaw on ice and gently mix prior to use. DO NOT VORTEX. Perform a quick spin before opening. Aliquot into small volumes and flash freeze for long term storage. Avoid multiple freeze/thaw cycles.

Quality Control Data

4-20% SDS-PAGE Coomassie Staining

