

## Product Information

<b>Description:</b>	Recombinant human HER3 (Receptor tyrosine-protein kinase erbB-3), encompassing amino acids (684-1020) length. This construct contains an N-terminal GST-tag and contains mutation of interest (E928G). The recombinant protein was affinity purified.
<b>Species:</b>	Human
<b>Construct:</b>	HER3 (E928G) (GST-684-1020)
<b>Mutation:</b>	E928G
<b>Concentration:</b>	0.20 mg/ml
<b>Expression System:</b>	Sf9
<b>Purity:</b>	90%
<b>Format:</b>	Aqueous buffer solution.
<b>Formulated In:</b>	50 mM Tris-HCl, pH 7.5, 150 mM NaCl, 10 mM glutathione, 0.1 mM EDTA, 0.25 mM DTT, 25% glycerol
<b>MW:</b>	65 kDa
<b>Uniprot:</b>	P21860
<b>Stability:</b>	At least 6 months at -80°C.
<b>Storage:</b>	-80°C
<b>Instructions for Use:</b>	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

