

## 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. The recombinant protein was affinity purified. |
| <b>Species:</b>              | Human   |
| <b>Construct:</b>            | HER3 (GST-684-1020)   |
| <b>Concentration:</b>        | 0.10 mg/ml  |
| <b>Expression System:</b>    | Sf9   |
| <b>Purity:</b>               | 85%   |
| <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>                   | 64 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

