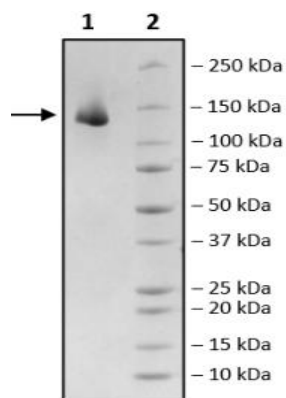


Product Information

Description:	Recombinant human HER3 (human epidermal growth factor receptor 3), encompassing amino acids 20-643 corresponding to the extracellular domain. This construct contains an Fc domain of IgG1 fused to the C-terminus, followed by a C-terminal Avi-Tag™. This protein was affinity purified.
Background:	HER3 (human epidermal growth factor receptor 3, also known as ERB3) is a transmembrane protein encoded by the ERBB3 gene. HER3 is broadly expressed in human tissues. However, increased expression of HER3 has been linked to a variety of solid tumors including ovarian, breast, colon, and gastric cancers and correlates with decreased overall survival in colorectal cancer patients. HER3 has the unique property of being inherently catalytically inactive but still being able to participate in ligand binding, forming heterodimers with nearby receptors of the HER family. The formation of these dimers initiates a cascade of downstream signaling steps critical to cell proliferation. The binding and subsequent phosphorylation of HER3 by HER2 or EGFR (epidermal growth factor receptor, also known as HER1) has been found to play a role in tumor growth and drug resistance. Due to its cell surface expression in a variety of cancers and correlation with decreased survival, HER3 is a prospective therapeutic target for antibody-drug conjugate (ADC) development.
Species:	Human
Construct:	HER3 (20-643-Fc(IgG1)-Avi)-(Biotin)
Concentration:	1.38 mg/ml
Expression System:	HEK293
Purity:	≥90%
Format:	Aqueous buffer solution.
Formulated In:	8 mM phosphate, pH 7.4, 110 mM NaCl, 2.2 mM KCl, and 20% glycerol
MW:	98 kDa + glycans
Glycosylation:	This protein runs at a higher MW by SDS-PAGE due to glycosylation.
Genbank Accession:	NM_001982.4
Label:	This protein is enzymatically biotinylated using Avi-Tag™ technology. Biotinylation is confirmed to be ≥90%.
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.
Applications:	Useful for SDS-PAGE and avidin-pull down assays.

Quality Control Data

4-20% SDS-PAGE Coomassie Staining



Biotin-Avidin Pulldown

