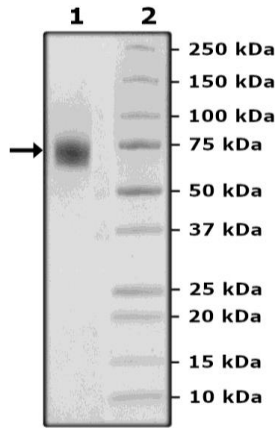


Product Information

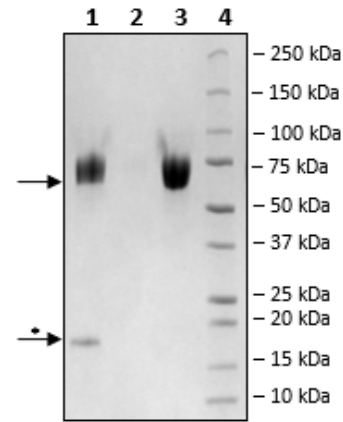
Description:	Recombinant human FGFR3 (fibroblast growth factor receptor 3, also known as CD333), encompassing amino acids 23-375 corresponding to the extracellular domain of the receptor. This construct contains a C-terminal Avi-Tag™ followed by a His-Tag (6xHis). This protein was affinity purified. HiP™ indicates a high purity protein (≥90% pure) and less than 10% aggregation as measured by gel filtration.
Background:	FGFR3, also known as fibroblast growth factor receptor 3 or CD333, is a transmembrane tyrosine kinase receptor which belongs to the fibroblast growth factor receptor family involved in osteogenesis and bone maintenance. It has a restricted pattern of expression and is found in the brain, kidneys, cartilage and intestine. Mutations in FGFR3 can result in achondroplasia, hypochondroplasia, Muenke syndrome, glioblastoma and urothelial carcinoma. Inhibitors of FGFR3 show promise for cancer therapy, with pemigatinib approved in 2020 for the treatment of metastatic cholangiocarcinoma.
Species:	Human
Construct:	FGFR3 (23-375-Avi-His)-(Biotin)
Concentration:	0.88 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:	41 kDa + glycans
Glycosylation:	This protein runs at a higher MW by SDS-PAGE due to glycosylation.
Aggregation:	<10%
Genbank Accession:	NM_000142.5
Label:	This protein is enzymatically biotinylated using Avi-Tag™ technology. Biotinylation was 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.

Quality Control Data

4-20% SDS-PAGE Coomassie Staining



Biotin-Avidin Pulldown



1. Beads
2. Flow thru
3. Control
4. Standards

* Avidin from beads.

Gel Filtration Curve

