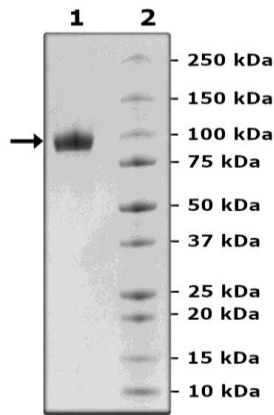


## Product Information

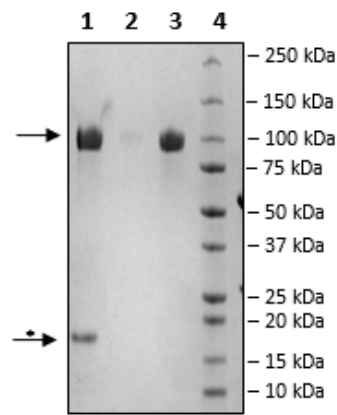
<b>Description:</b>	Recombinant human FGFR3 (fibroblast growth factor receptor 3, also known as CD333), encompassing amino acids 23-375 corresponding to the extracellular portion of the receptor. This construct contains a C-terminal Fc domain followed by an Avi-Tag™. This protein has a Factor Xa cleavage site prior to the Fc domain. This protein was affinity purified.
<b>Background:</b>	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.
<b>Species:</b>	Human
<b>Construct:</b>	FGFR3 (23-375-Fc(IgG1)-Avi)-(Biotin)
<b>Concentration:</b>	0.66 mg/ml
<b>Expression System:</b>	HEK293
<b>Purity:</b>	≥90%
<b>Format:</b>	Aqueous buffer solution.
<b>Formulated In:</b>	8 mM phosphate, pH 7.4, 110 mM NaCl, 2.2 mM KCl, and 20% glycerol
<b>MW:</b>	67 kDa + glycans
<b>Glycosylation:</b>	This protein runs at a higher MW by SDS-PAGE due to glycosylation.
<b>Genbank Accession:</b>	NM_000142.5
<b>Label:</b>	This protein is enzymatically biotinylated using Avi-Tag™ technology. Biotinylation was confirmed to be ≥90%.
<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



### Biotin-Avidin Pulldown



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

\* Avidin from beads.