

## Product Information

<b>Description:</b>	Recombinant human FGFR3 (fibroblast growth factor receptor 3), encompassing amino acids 23-375 corresponding to the extracellular domain. This construct contains a C-terminal Avi-Tag™, followed by an His-tag (6xHis). This protein was affinity purified.
<b>Background:</b>	FGFR3 (fibroblast growth factor receptor 3, also known as 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 FGFR3 inhibitor pemigatinib approved in 2020 for the treatment of metastatic cholangiocarcinoma.
<b>Species:</b>	Human
<b>Construct:</b>	FGFR3 (23-375-Avi-His)
<b>Concentration:</b>	2.33 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>	41 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>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.
<b>Applications:</b>	Useful for SDS-PAGE.

## Quality Control Data

### 4-20% SDS-PAGE Coomassie Staining

