

## Product Information

<b>Description:</b>	Recombinant human FGFR2C (Fibroblast Growth Factor Receptor 2C) protein corresponding to the extracellular domain, encompassing amino acids 22-377. This construct contains a C-terminal human IgG1 Fc domain, followed by an Avi-Tag™. This protein was affinity purified.
<b>Background:</b>	FGFR2 (Fibroblast Growth Factor Receptor 2) is a transmembrane receptor for FGF (fibroblast growth factor) with tyrosine kinase activity, involved in many pathways that play a significant role in cancer. Amplification or activation of FGFR2 has been reported in breast and gastric cancers, while FGFR2 mutations have been observed in endometrial and breast cancers. Mutations in FGFR2 are also associated with bone development disorders including Pfeiffer Syndrome and Crouzon Syndrome. The use of FGFR2 inhibitors and degraders are active areas of research and are promising approaches in cancer therapy.
<b>Species:</b>	Human
<b>Construct:</b>	FGFR2C (22-377-Fc(IgG1)-Avi)
<b>Concentration:</b>	1.32 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>	68 kDa + glycans
<b>Glycosylation:</b>	This protein runs at a higher MW by SDS-PAGE due to glycosylation.
<b>Genbank Accession:</b>	NM_000141.5
<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

