

Product Information

Description:	Human recombinant FGFR2C (Fibroblast growth factor receptor 2C, also known as CD332), corresponding to amino acids 22-377, the extracellular domain. This construct contains a C-terminal Avi-Tag™ followed by a His-Tag (6xHis). The recombinant protein was affinity purified.
Background:	Fibroblast Growth Factor Receptor 2 (FGFR2) belongs to a family of transmembrane receptors with tyrosine kinase activity playing a key role in cell growth, differentiation, and tissue repair. Receptor activation by FGF ligands triggers signaling pathways that ultimately regulate cell proliferation and differentiation. FGFR2 is essential for embryonic development, particularly in the formation of the skeletal system and the development of the central nervous system. It also plays a significant role in wound healing, tissue regeneration, and tissue maintenance in adults. Dysfunction of FGFR2 is associated with a range of disorders, including craniosynostosis and certain types of cancer, therefore, FGFR2 is a candidate therapeutic target in conditions where its function is dysregulated.
Species:	Human
Construct:	FGFR2C (22-377-Avi-His)
Concentration:	1.86 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:	43 kDa + glycans
Glycosylation:	This protein runs at a higher MW by SDS-PAGE due to glycosylation.
Genbank Accession:	NM_000141.5
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

