

FGF19, Avi-Tag, His-Tag, Biotin-Labeled Recombinant

Catalog: 101888

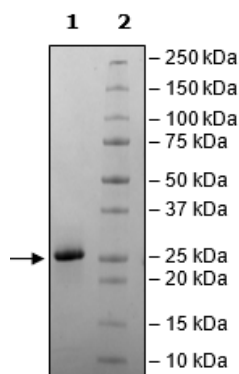
Lot: 230817

Product Information

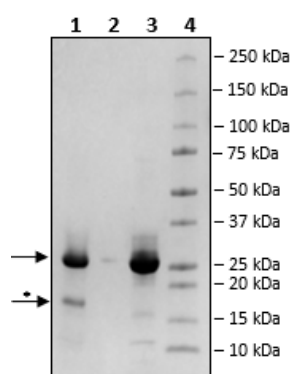
Description:	Recombinant human FGF19 (fibroblast growth factor 19), encompassing amino acids 25-216(end). This construct contains a C-terminal Avi-Tag™ followed by an His-tag (6x His). This protein was affinity purified.
Background:	FGF19, or fibroblast growth factor 19, is part of the FGF family of growth factors. It is involved in bile acid synthesis, where it functions as a hormone to regulate glucose and lipid metabolism. Defective levels of FGF19 can result in primary bile acid diarrhea, and FGF19 can be found in the liver of patients suffering from cholestasis. FGF19 overexpression has been found in cancer, such as prostate, bladder and lung cancer. Of note is its presence in almost half of the hepatocellular carcinomas. FGF19 is the ligand for FGFR4 (fibroblast growth factor receptor 4), and results in the activation of the receptor, acting like an oncogene. The development of specific inhibitors for FGFR4 activation by FGF19 may reveal as useful therapeutical tools for cancer treatment.
Species:	Human
Construct:	FGF19 (25-216(end)-Avi-His)-(Biotin)
Concentration:	1.30 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:	24 kDa
Genbank Accession:	NM_005117.3
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.
Applications:	Useful for SDS-PAGE and pulldown assays.

Quality Control Data

4-20% SDS-PAGE Coomassie Staining



Biotin-Avidin Pulldown



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

* Avidin from beads.