

VEGFR2 (KDR), Avi-Tag, His-Tag, Biotin-Labeled Recombinant

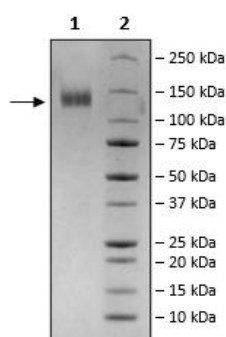
Catalog: 102133
Lot: 240306-2

Product Information

Description:	Recombinant human VEGFR2 (vascular endothelial growth factor receptor 2), also known as kinase insert domain receptor (KDR), encompassing amino acids 20-764 corresponding to the extracellular domain of the receptor. This construct contains a C-terminal Avi-Tag™ followed by an His-tag (6xHis). This protein was affinity purified.
Background:	The interaction between VEGF (Vascular Endothelial Growth Factor) and its receptor VEGFR is an important mediator of angiogenesis as well as endothelial cell survival and proliferation. VEGFR2 is a receptor tyrosine kinase that plays a major role in cellular responses driven by VEGF. Targeting the kinase activity of VEGFR2 has been an attractive tool in cancer therapy based on its anti-angiogenesis effects. In addition to the kinase inhibitors, humanized monoclonal antibodies interfering with VEGF/VEGFR2 interaction have been developed as anti-cancer treatments.
Species:	Human
Construct:	VEGFR2 (20-764-Avi-His)-(Biotin)
Concentration:	1.07 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:	87 kDa + glycans
Glycosylation:	This protein runs at a higher MW by SDS-PAGE due to glycosylation.
Genbank Accession:	NM_002253.4
Label:	This protein is enzymatically biotinylated using Avi-Tag™ technology. Biotinylation is 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 the SDS-PAGE and avidin-binding assays.

Quality Control Data

4-20% SDS-PAGE Coomassie Staining



Biotin-Avidin Pulldown

