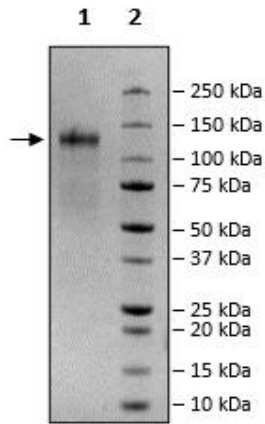


Product Information

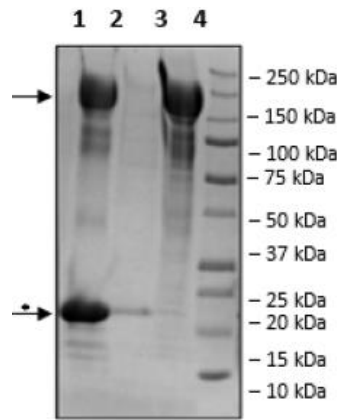
Description:	Recombinant human VEGFR3 (Vascular endothelial growth factor receptor 3), extracellular portion encompassing amino acids 25-776. This construct contains a C-terminal Avi-Tag™ followed by a His-tag (6x His). This protein was affinity purified.
Background:	VEGFR3 (also known as FLT4, fms related receptor tyrosine kinase 4) is one of three VEGFRs and has VEGFC and VEGFD as ligands. Its main roles are in angiogenesis, lymphangiogenesis, and vasculogenesis. VEGFR3 is a transmembrane receptor expressed as a homodimer, with tyrosine kinase activity. This protein can exist as three isoforms: full length, an isoform missing 65 amino acids on the C-terminus, and a third isoform missing a larger fraction of the C-terminus including the transmembrane domain. This third form is soluble and seems to function in sequestering VEGFC in the retina. VEGFR3 can form heterodimers with VEGFR2 and bind VEGFC, functioning in angiogenesis and hematopoiesis. Dysfunction of this protein can result in Milroy disease, congenital heart disease and cancer. The use of inhibitors targeting VEGFR3 can reduce angiogenesis and be beneficial in cancer therapy.
Species:	Human
Construct:	VEGFR3 (25-776-Avi-His)-(Biotin)
Concentration:	0.46 mg/ml
Expression System:	HEK293
Purity:	83%
Format:	Aqueous buffer solution.
Formulated In:	8 mM phosphate, pH 7.4, 110 mM NaCl, 2.2 mM KCl, and 20% glycerol
MW:	88 kDa + glycans
Glycosylation:	This protein runs at a higher MW by SDS-PAGE due to glycosylation.
Genbank Accession:	NM_182925
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 SDS-PAGE and avidin binding 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.