

Recombinant KDR / VEGFR2 (789-1356) protein

Catalog No: 81353, 81753

Expressed In: Baculovirus

Quantity: 20, 1000 µg

Concentration: 0.2 µg/µl

Source: Human

Buffer Contents: Recombinant KDR / VEGFR2 (789-1356) protein is supplied in 25 mM HEPES-NaOH pH 7.5, 300 mM NaCl, 10% glycerol, 0.04% Triton X-100, and 0.5 mM TCEP.

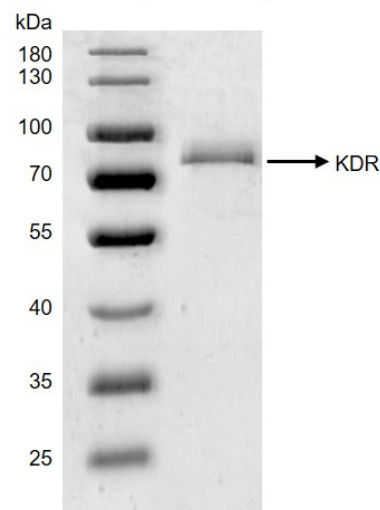
Background: KDR (Kinase Insert Domain Receptor) is a tyrosine-protein kinase that acts as a cell-surface receptor for VEGFA, VEGFC and VEGFD and plays an essential role in the regulation of angiogenesis, vascular development, vascular permeability, and embryonic hematopoiesis. KDR promotes proliferation, survival, migration and differentiation of endothelial cells and promotes reorganization of the actin cytoskeleton. Isoforms lacking a transmembrane domain, such as isoform 2 and isoform 3, may function as decoy receptors for VEGFA, VEGFC and/or VEGFD. Isoform 2 plays an important role as negative regulator of VEGFA- and VEGFC-mediated lymphangiogenesis by limiting the amount of free VEGFA and/or VEGFC and preventing their binding to FLT4. Modulates FLT1 and FLT4 signaling by forming heterodimers. Binding of vascular growth factors to isoform 1 leads to the activation of several signaling cascades. Diseases associated with KDR is infantile capillary hemangiomas.

Protein Details: Recombinant KDR / VEGFR2 (789-1356) protein that includes amino acids 789-1356 of human KDR protein (accession number NP_002244.1) with a N-terminal FLAG tag. The molecular weight of the protein is 64.4 kDa.

Application Notes: This product was manufactured as described in Protein Details. Where possible, Active Motif has developed functional or activity assays for recombinant proteins. Additional characterization such as enzyme kinetic activity assays, inhibitor screening or other biological activity assays may not have been performed for every product. All available data this product is shown.

Storage and Guarantee: Recombinant proteins in solution are temperature sensitive and must be stored at -80°C to prevent degradation. Avoid repeated freeze/thaw cycles and keep on ice when not in storage. This product is for research use only and is not for use in diagnostic procedures. This product is guaranteed for 6 months from date of arrival.

KDR (789-1356)



Recombinant KDR / VEGFR2 (789-1356) protein gel

10% SDS-PAGE with Coomassie staining

MW: 64.4 kDa

Purity: >90%