

Product Information

Description:	Recombinant human DLL4 (delta-like canonical Notch ligand 4), encompassing amino acids 27-529. This construct contains the Fc region of human IgG1 on the C-terminus, followed by an Avi-Tag™. This protein was affinity purified.
Background:	DLL4, or delta-like canonical Notch ligand 4, is part of the Notch signaling pathway and it is involved in normal tissue development and homeostasis. It is an agonist ligand and can bind to the four Notch receptors. It is expressed in the vascular system and plays a crucial role in angiogenesis. Loss of function of DLL4 results in VEGF upregulation, and the formation of immature vessels without a lumen. It has been shown that DLL4 inhibition can actually result in tumor regression by its role on VEGF upregulation. DLL4 is found in several tumor types, and it is found in tumor stem cells. Inhibition of DLL4 can thus contribute to cancer therapy via its several mechanisms of action.
Species:	Human
Construct:	DLL4 (27-529-Fc-Avi)
Concentration:	0.25 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:	84 kDa + glycans
Glycosylation:	This protein runs at a higher MW by SDS-PAGE due to glycosylation.
Genbank Accession:	NM_019074.4
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 to study the binding of DLL4 in ELISA and in cellular assays.

Quality Control Data

4-20% SDS-PAGE Coomassie Staining

