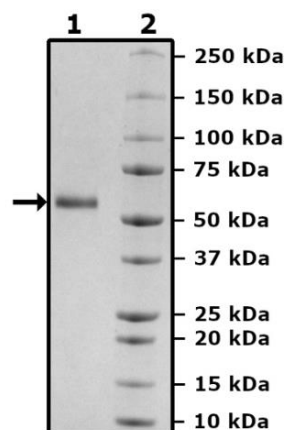


Product Information

Description:	Recombinant human DLL1 (delta like canonical Notch ligand 1), encompassing amino acids 18-540. This construct contains a C-terminal Avi-Tag™ followed by a His-tag (6xHis). The recombinant protein was affinity purified.
Background:	DLL1 is involved in cell-to-cell communication and in embryonic development, neurogenesis, and tissue homeostasis. It is a member of the Delta-Serrate-Lag2 (DSL) family of ligands that activate the Notch signaling pathway. DLL1 binds to the Notch receptors on neighboring cells, triggering signals that regulate cell fate determination, proliferation, and differentiation. Dysregulation of DLL1 has been implicated in several diseases.
Species:	Human
Construct:	DLL1 (18-540-Avi-His)-(Biotin)
Concentration:	0.22 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:	60 kDa + glycans
Glycosylation:	This protein runs at a higher MW by SDS-PAGE due to glycosylation.
Genbank Accession:	NM_005618.4
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 to study DLL1 binding in ELISA.

Quality Control Data

4-20% SDS-PAGE Coomassie Staining



Biotin-Avidin Pulldown

