

Product Information

Description:	Recombinant human Notch2 (neurogenic locus notch homolog protein 2), encompassing amino acids 26-530 (extracellular domain). This construct contains a C-terminal Fc domain from IgG1 followed by an Avi-tag™. This protein was affinity purified.
Background:	Notch2 (neurogenic locus notch homolog protein 2) is a type I transmembrane protein of the Notch family. It is involved in the development of the vascular network, brain, kidneys, liver, T and B lymphocytes. The protein undergoes cleavage in the trans-golgi network, forming a heterodimer. Binding of Notch2 to its ligands of the Delta/Serrate/LAG-2 family leads to its cleavage by ADAM (A disintegrin and metalloprotease) proteins. NICD (notch intracellular domain) is released and translocated to the nucleus, where it can form complexes with DNA-binding proteins and lead to transcription of target genes. Mutations in this protein can result in Hadju-Cheney syndrome. Overexpression of Notch2 is found in several cancer types, such as LSCC (laryngeal squamous cell carcinoma) and B cell malignancies. The use of inhibitors to target this protein is being studied with the goal of understanding better the role of this protein and potentially target it in cancer therapy.
Species:	Human
Construct:	Notch2 (26-530-Fc(IgG1)-Avi)
Concentration:	0.24 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:	83 kDa + glycans
Glycosylation:	This protein runs at a higher MW by SDS-PAGE due to glycosylation.
Genbank Accession:	NM_024408.4
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 for SDS-PAGE.

Quality Control Data

4-20% SDS-PAGE Coomassie Staining

