

Product Information

Description:	Recombinant human full length JosD1. The construct contains an N-terminal His-Tag. The protein was affinity purified and is active.
Background:	JOSD1 (Josephin-1) is a josephin domain-containing protein that has a deubiquitinase activity in vitro. Josephin domain-containing 1, often abbreviated as JOSD1, is a human gene that encodes a protein with a Josephin domain, which is a type of cysteine protease that can remove ubiquitin from other proteins. JOSD1 has been implicated in various studies to promote cancer cell growth and regulate membrane dynamics, cell motility, and endocytosis.
Species:	Human
Construct:	JosD1 (His-1-202(end))
Expression System:	<i>E. coli</i>
Purity:	≥90% by SDS-PAGE
Format:	Aqueous buffer solution.
Formulated In:	50 mM Tris pH 8.0, 150 mM NaCl, 10% glycerol
MW:	23 kDa
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.
Assay Conditions:	This enzyme has been validated using the Ub-CHOP2-Reporter Deubiquitination Assay Kit (LifeSensors #PR1101).
Applications:	Useful for the study of enzyme kinetics.

Quality Control Data

SDS-PAGE Coomassie Staining

