

Product Information

Description:	Recombinant human full length JosD2. The construct contains an N-terminal His-Tag. The protein was affinity purified and is active.
Background:	JOSD2 (Josephin-2) is a josephin domain-containing protein that has deubiquitinating activity <i>in vitro</i> . It cleaves K63-linked polyubiquitin chains, and with lesser efficiency K48-linked polyubiquitin chains.
Species:	Human
Construct:	JosD2 (His-1-188(end))
Expression System:	<i>E. coli</i>
Format:	Aqueous buffer solution.
MW:	22 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.