

Product Information

Description:	Recombinant human full length Praja1. The construct contains an N-terminal His-SUMO-Tag. The protein was affinity purified and is active.
Background:	Praja1 (along with Praja2, 53% identity) is a RING-H2 domain E3 (ubiquitin protein ligase) that is involved in the conjugation of ubiquitin to target substrates along with E1 and E2 enzymes. Substrates for Praja1 include Dlxin-1 (MAGE-D1, human) which regulates Dlx5 dependent transcriptional function (involved in osteoblast differentiation) and ELF and Smad3 involved in the TGF- β signaling pathway implicated in gastrointestinal cancers.
Species:	Human
Construct:	Praja1 (His-SUMO-1-643(end))
Expression System:	<i>E. coli</i>
Purity:	$\geq 90\%$ by SDS-PAGE
Format:	Aqueous buffer solution.
Uniprot:	Q8NG27
MW:	71 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:	Typical enzyme concentration of 100 nM - 5 mM is used for in vitro conjugation, depending on conditions.
Applications:	Useful for the study of enzyme kinetics.