

Praja1 (E3 ubiquitin protein ligase Praja1)

Cat. # UB303

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.

Alternate names: RNF70 (RING finger protein 70)

Product Information

Purity:	\geq 85% by SDS-PAGE
Molecular Weight:	71kDa
Quantity:	25 μ g
Physical State:	Liquid
Source:	Recombinant human
Tag:	His ₆ SUMO
Activity:	Typical enzyme concentration of 100 nM - 5 mM is used for in vitro conjugation, depending on conditions.
Storage:	-80° C. Avoid repeated freeze/thaw cycles

References

- 1) Yu, P., et al., PJA1, encoding a RING-H2 finger ubiquitin ligase, is a novel human X chromosome gene abundantly expressed in brain. *Genomics*, 2002. **79**:869-874.
- 2) Mishra, L., V. Katuri, and S. Evans, The role of PRAJA and ELF in TGF-beta signaling and gastric cancer. *Cancer Biol Ther*, 2005. **4**:694-699.
- 3) Saha, T., et al., RING finger-dependent ubiquitination by PRAJA is dependent on TGF-beta and potentially defines the functional status of the tumor suppressor ELF. *Oncogene*, 2006. **25**:693-705.
- 4) Sasaki, A., et al., A RING finger protein Praja1 regulates Dlx5-dependent transcription through its ubiquitin ligase activity for the Dlx/Msx-interacting MAGE/Necdin family protein, Dlxin-1. *J Biol Chem*, 2002. **277**:22541-22546.

All products are for research use only • Not intended for human or animal diagnostic or therapeutic uses
Copyright © 2009 LifeSensors, Inc. All Rights Reserved