

Praja1 (E3 ubiquitin protein ligase Praja1)

Cat. # UB303

Background

Praja1 (along with its homolog Praja2, which shares 53% identity) is a RING-H2 domain E3 ubiquitin protein ligase involved in the conjugation of ubiquitin to target substrates, in collaboration with E1 and E2 enzymes. Substrates for Praja1 include Dlxin-1 (also known as MAGE-D1 in humans), which regulates Dlx5-dependent transcriptional functions involved in osteoblast differentiation, as well as ELF and Smad3, which play roles in the TGF- β signaling pathway implicated in gastrointestinal cancers.

Alternate Names

Praja Ring Finger Ubiquitin Ligase 1, RING Finger Protein 70 (RNF70), Praja Ring Finger 1, Praja Ring Finger 1 E3 Ubiquitin Protein Ligase, RING-Type E3 Ubiquitin Transferase Praja-1

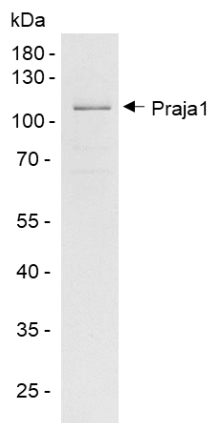
Application(s)

In vitro conjugation assay

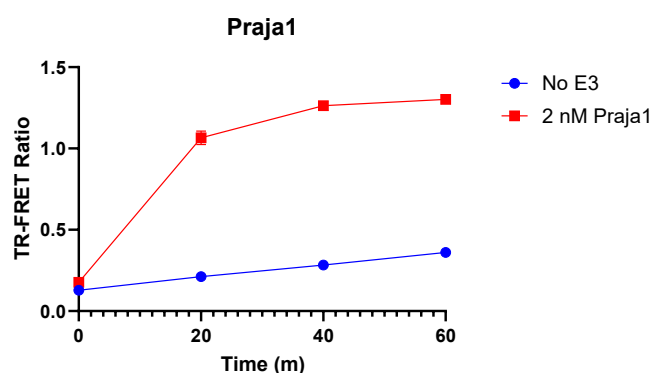
Product Specifications

| | |
|-------------------|---|
| Affinity tag | His6-SUMO |
| Purity | > 90% by SDS-PAGE |
| Molecular Weight | 71 kDa (without Tag), ~ 110 kDa with tags |
| Quantity | 25 μ g |
| Species | Human |
| Expression System | <i>E. coli</i> |
| Physical State | Liquid |
| Buffer | 50 mM Tris, 150 mM NaCl, 10 mM DTT, 10% glycerol |
| Activity | Typical enzyme concentration of 2-100 nM for in vitro conjugation, condition-dependent. |
| Storage | Store at -80°C. Avoid repeated freeze/thaw cycles |

Product QC



SDS-Page Analysis of purified Praja1. Two μ g of the protein was loaded on a 10-20% SDS-PAGE gel and stained with Coomassie brilliant blue.



Activity Assay of Praja1. 2 nM Praja1 was tested in a TR-FRET assay showing robust E3 ligase activity.

All products are for research use only • Not intended for human or animal diagnostic or therapeutic uses
Copyright © 2025 LifeSensors, Inc. All Rights Reserved

Praja1 (E3 ubiquitin protein ligase Praja1)

Cat. # UB303

References

1. Consalvi, S., et al., Nat Commun., 2017. 8:13956.
2. Watabe, K., et al., Neuropathology, 2022. 42(6):488-504.

All products are for research use only • Not intended for human or animal diagnostic or therapeutic uses
Copyright © 2025 LifeSensors, Inc. All Rights Reserved