

USP34core (Ubiquitin Specific Protease 34 catalytic domain)

Cat. # DB506

Background: USP34 regulates the Wnt signaling pathway by deubiquitinylation of AXIN1 and AXIN2. The catalytic core consists of residues 1892-2241. Deubiquitinylation leads to the stabilization and nuclear accumulation of AXIN1 and AXIN2[1]

Alternate names: FLJ43910, KIAA0570, KIAA0729

Product Information

Accession No.	Q70CQ2
Molecular Weight:	42 kDa
Quantity:	25 µg
Physical State:	Liquid, 25 mM Tris-HCl, pH 7.5, 150 mM NaCl, 5 mM DTT, 10% glycerol
Source:	Human Recombinant
Tag:	His ₆
Activity:	This enzyme is active in the Ub-CHOP assay.
Storage:	-80° C. Avoid repeated freeze/thaw cycles

References

1. Lui, T.T., Lacroix, C., Ahmed, S.M., Goldenberg, S.J., Leach, C.A., Daulat, A.M., and Angers, S., The ubiquitin specific protease USP34 regulates Axin stability and Wnt/beta-catenin signaling. *Mol Cell Biol.* **31**:2053-2065 (2011).

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