

BAP1 (BRCA1-associated protein 1)

Cat. # DB103

Background: BAP1 is the first large, nuclear-localized UCH to be identified. It is a protein of 90 kDa that binds to the BRCA RING finger and enhances BRCA1-mediated cell growth suppression¹.

Alternate names: BRCA1-associated protein 1, Cerebral protein 6, DKFZp686N04275, FLJ35406, FLJ37180, HUCEP-13, hucep-6, KIAA0272, Ubiquitin carboxyl-terminal hydrolase BAP1

Product Information

Molecular Weight:	82kDa
Quantity:	25µg
Physical State:	Liquid
Source:	Human Recombinant
Tag:	His6
Activity:	This enzyme cleaves Ub-AMC
Storage:	-80°C. Avoid repeated freeze/thaw cycles

References

- 1) David E. Jensen and Frank J. Rauscher III. BAP1, a Candidate Tumor Suppressor Protein That Interacts with BRCA1. *Ann N Y Acad Sci.* 1999. **886**:191-194.

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