

BAP1 (BRCA1-associated protein 1)

Cat. # DB103

Background

BAP1 plays a critical role in regulating various cellular processes, including DNA repair, cell cycle control, and gene expression. It is particularly known for its involvement in maintaining genomic stability by deubiquitinating histones and other proteins involved in chromatin remodeling. Mutations or loss of BAP1 function have been associated with a higher risk of developing certain cancers, including uveal melanoma and mesothelioma, underscoring its significance as a guardian of genome integrity.

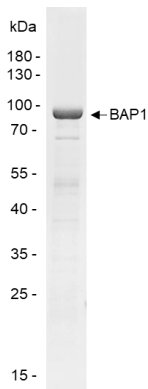
Alternate names

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

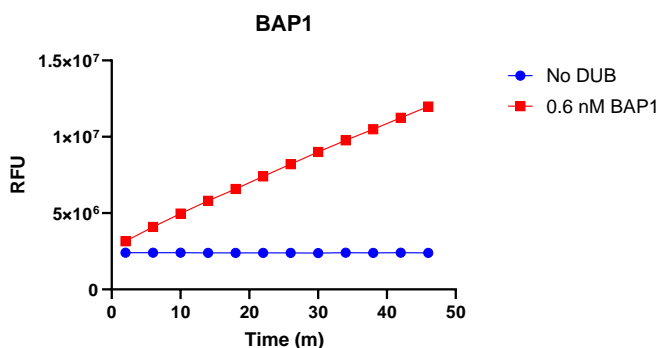
Product Information

Purity	≥ 85% by SDS-PAGE
Molecular Weight	82 kDa
Quantity	25
Physical State	Liquid
Source	Human Recombinant
Tag	His6
Activity	This enzyme cleaves Ub-AMC
Storage	-80°C. Avoid repeated freeze/thaw cycles

Product QC



SDS-Page Analysis of purified BAP1. Two μg of the enzyme was loaded on a 10-20% SDS-PAGE gel and stained with Coomassie brilliant blue.



Activity Assay of BAP1. 0.6 nM BAP1 was tested in a CHOP assay showing robust DUB activity.

References

1. Carbone, M., et al., *Cancer Discov.*, 2020. 10(8):1103-1120.
2. Masclef, L., et al., *Cell Death Differ.*, 2021. 28(2):606-625.

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