

BAP1 (human), (recombinant) (His- tag)

BAP1 is a BRCA1-associated ubiquitin C-terminal hydrolase. It shows homology to members of the ubiquitin C-terminal hydrolase (UCH) family and exhibits deubiquitylating activity *in vitro*.

Citations: 1

[View Online »](#)

Ordering Information

[Order Online »](#)

BML-UW9855-0050	50µg
-----------------	------

Manuals, SDS & CofA

[View Online »](#)

Handling & Storage

Long Term Storage -80°C

Shipping Dry Ice

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name BRCA1-associated protein 1, Ubiquitin C-terminal hydrolase BAP1

Application Notes **Uses:**
Deubiquitinylation assays using small molecule substrates (e.g. Ub-AMC) or ubiquitin-protein conjugates.
Studies of the tumor suppressor BRCA1, its interactions and associated proteins.

Formulation Liquid. In 50mM Na-phosphate, pH 8.0, 300mM NaCl.

MW 84.1 kDa

Purity ≥90% (SDS-PAGE)

Purity Detail Purified by Ni-affinity chromatography.

Source Produced in *E. coli*. Human BAP1 is fused at the N-terminus to a His-tag.

UniProt ID Q92560

Last modified: May 29, 2024



ENZO LIFE SCIENCES,
INC.
Phone: 800.942.0430
[info-
usa@enzolifesciences.com](mailto:info-usa@enzolifesciences.com)

European Sales Office
ENZO LIFE SCIENCES
(ELS) AG
Phone: +41 61 926 8989
[info-
eu@enzolifesciences.com](mailto:info-eu@enzolifesciences.com)

Belgium, The Netherlands
& Luxembourg
Phone: +32 3 466 0420
[info-
fr@enzolifesciences.com](mailto:info-fr@enzolifesciences.com)

France
Phone: +33 472 440 655
[info-
fr@enzolifesciences.com](mailto:info-fr@enzolifesciences.com)

Germany
Phone: +49 7621 5500 526
[info-
de@enzolifesciences.com](mailto:info-de@enzolifesciences.com)

UK & Ireland
Phone (UK customers):
0845 601 1488
Phone: +44 1392 825900
[info-
uk@enzolifesciences.com](mailto:info-uk@enzolifesciences.com)