BAP1 [GST-tagged]

Deconjugating enzyme: Deubiquitylase

Alternate Names: BRCA1 Associated Protein 1, BRCA1 Associated Protein-1 (Ubiquitin Carboxy-Terminal Hydrolase), Cerebral Protein 6, Ubiquitin Carboxy-Terminal Hydrolase 2, BRCA1-Associated Protein 1, Cerebral Protein-13

64-0052-050 Quantity: Cat. No. 50 µg Lot. No. 30421 Storage: -70°C

FOR RESEARCH USE ONLY NOT FOR USE IN HUMANS



CERTIFICATE OF ANALYSIS Page 1 of 2

Background

Deconjugating (DCEs) enzymes proteases that process ubiquitin ubiquitin-like gene products, reverse the modification of proteins by a single ubiquitin or ubiquitin-like protein (UBL) and remodel polyubiquitin (or poly-UBL) chains on target proteins (Reyes-Turcu et al. 2009). The deubiquitylating - or deubiquitinating enzymes (DUBs) represent the largest family of DCEs and regulate ubiquitin dependent signalling pathways. The activities of the DUBs include the generation of free ubiquitin from precursor molecules, the recycling of ubiquitin following substrate degradation to maintain cellular ubiquitin homeostasis and the removal of ubiquitin or ubiquitin-like proteins (UBL) modifications through chain editing to proteins from proteasomal degradation or to influence cell signalling events (Komander et al. 2009). There are two main classes of DUB, cysteine proteases and metalloproteases. BRCA1 Associated Protein 1 (BAP1) is a cysteine protease and member of the UCH family of ubiquitin Cterminal hydrolases. Cloning of human BAP1 gene was first described by Jensen et al. (1998). The nuclear DUB BAP1 is a tumour suppressor deleted and mutated in an increasing number of origins thereby of diverse cancers making BAP1 the most frequently widelv mutated **DUB-encoding** gene in cancer (Daou et al. 2015). BAP1 contains binding domains for BRCA1 (Breast cancer type 1) and BARD1 (BRCA1-associated **RING** domain protein 1), which form tumour heterodimeric suppressor complex, and HCFC1 (Host cell factor interacts with which histone-modifying complexes

Physical Characteristics

Protein Sequence: Please see page 2 Species: human

Source: E. coli Quantity: 50 µg

Concentration: 0.18 mg/ml

Formulation: 50 mM Tris/HCl pH 7.5,

0.1 mM EGTA, 150 mM NaCl, 0.1% ß-Mercaptoethanol, 270 mM Sucrose, 0.03% Brij-35

Molecular Weight: ~107 kDa

Purity: >80% by InstantBlue™ SDS-PAGE

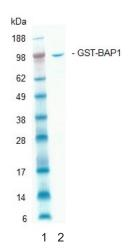
Stability/Storage: 12 months at -70°C;

aliquot as required

Quality Assurance

Purity:

4-12% gradient SDS-PAGE InstantBlue™ staining Lane 1: MW markers Lane 2: 1 µg GST-BAP1



Protein Identification:

Confirmed by mass spectrometry.

Deubiquitylase Enzyme Assay:

The activity of GST-BAP1 was validated by determining the increase in fluorescence measured as a result of the enzyme catalysed cleavage of the fluorogenic substrate Ubiquitin-Rhodamine110-Glycine generating Ubiquitin and Rhodamine110-Glycine. Incubation of the substrate in the presence or absence of GST-BAP1 was compared confirming the deubiquitylating activity of GST-BAP1.

Continued on page 2



Dundee, Scotland, UK

ORDERS / SALES SUPPORT

International: +1-617-245-0003 US Toll-Free: 1-888-4E1E2E3 (1-888-431-3233) Email: sales.support@ubiquigent.com

UK HQ and TECHNICAL SUPPORT International: +44 (0) 1382 381147 (9AM-5PM UTC) US/Canada: +1-617-245-0020

(9AM-5PM UTC) Email: tech.support@ubiquigent.com

Email services@ubiquigent.com for enquiries regarding compound profiling and/or custom assay development services.

© **Ubiquigent 2012**. Unless otherwise noted, Ubiquigent, Ubiquigent logo and all other trademarks are the property of Ubiquigent, Ltd.

Limited Terms of Use: For research use only. Not for use in humans or for diagnostics. Not for distribution or resale in any form, modification or derivative OR for use in providing services to a third party (e.g. screening or profiling) without the written permission of Ubiquigent, Ltd.

Lot-specific COA version tracker: v1.0.0

BAP1 [GST-tagged]

Deconjugating enzyme: Deubiquitylase

Alternate Names: BRCA1 Associated Protein 1, BRCA1 Associated Protein-1 (Ubiquitin Carboxy-Terminal Hydrolase), Cerebral Protein 6, Ubiquitin Carboxy-Terminal Hydrolase 2, BRCA1-Associated Protein 1, Cerebral Protein-13

Cat. No. 64-0052-050 Quantity: 50 µg Lot. No. 30421 Storage: -70°C

FOR RESEARCH USE ONLY NOT FOR USE IN HUMANS



CERTIFICATE OF ANALYSIS Page 2 of 2

Background

Continued from page 1

during cell division. BAP1 also interacts with ASXL1 (Putative Polycomb protein) to form the Polycomb group repressive de-ubiquitylase complex (PR-DUB), which is involved in stem pluripotency and other developmental (Harbour processes et al. 2010). BAP1 More recently, has been identified as a DUB for Krüppel-like zincfinger transcription factor (KLF5). a transcription factor that is highly expressed in certain types of breast cancers. KLF5 has been identified as an unstable protein that is ubiquitylated by WWP1 (NEDD4-like E3 ligase), SCFFBW7 SKP1-cullin-1-F-box complex that (a contains FBW7 as the F-box protein) and Smurf2 (SMAD Specific E3 ligase degraded. BAP1 promotes 2) and breast cancer cell proliferation and migration vitro and tumour growth and lung metastasis in vivo. The results from this study suggest that BAP1 KLF5 are potential therapeutic targets breast cancer for (Qin et al. 2015).

References:

Daou S, et al. (2015) The BAP1/ASXL2 histone H2A deubiquitinase complex regulates cell proliferation and Is disrupted in cancer. The Journal of biological chemistry 290:28643-28663.

Harbour JW, et al. (2010) Frequent mutation of BAP1 in metastasizing uveal melanomas. Science 330:1410-1413

Jensen DE, et al. (1998) BAP1: a novel ubiquitin hydrolase which binds to the BRCA1 RING finger and enhances BRCA1mediated cell growth suppression. Oncogene 16:1097-1112.

Komander D, et al. (2009) Breaking the chains: structure and function of the deubiquitinases. Nat Rev Mol Cell Biol 10:550-563.

Qin J, et al. (2015) BAP1 promotes breast cancer cell proliferation and metastasis by deubiquitinating KLF5. Nature communications 6:8471.

Reyes-Turcu FE et al. (2009) Regulation and cellular roles of ubiquitin-specific deubiquitinating enzymes. Annual review of biochemistry 78:363-397.

Physical Characteristics

Continued from page 1

Protein Sequence:

MSPILGYWKIKGLVQPTRLLLEYLEEKYEEH LYERDEGDKWRNKKFELGLEFPNLPYY IDGDVKLTQSMAIIRYIADKHNMLGGCPKER AEISMLEGAVLDIRYGVSRIAYSKDFETLKVD FLSKLPEMLKMFEDRLCHKTYLNGDHVTHP DFMLYDALDVVLYMDPMCLDAFPKLVCFK KRIEAIPQIDKYLKSSKYIAWPLQGWQATF GGGDHPPKSDLEVLFQGPLGSMNKGWLELES DPGLFTLLVEDFGVKGVQVEEIYDLQSKCQGPVY GFIFLFKWIEERRSRRKVSTLVDDTSVIDDDIVN NMFFAHQLIPNSCATHALLSVLLNCSSVDLGPTL SRMKDFTKGFSPESKGYAIGNAPELAKAHNSHAR PEPRHLPEKQNGLSAVRTMEAFHFVSYVPITGRL FELDGLKVYPIDHGPWGEDEEWTDKARRVIMERI GLATAGEPYHDIRFNLMAVVPDRRIKYEARLHVL KVNRQTVLEALQQLIRVTQPELIQTHKSQESQLP EESKSASNKSPLVLEANRAPAASEGNHTDGAEEA AGSCAQAPSHSPPNKPKLVVKPPGSSLNGVHPNP TPIVQRLPAFLDNHNYAKSPMQEEEDLAAGVGRS RVPVRPPQQYSDDEDDYEDDEEDDVQNTNSALRY KGKGTGKPGALSGSADGQLSVLQPNTINVLAEKL KESQKDLSIPLSIKTSSGAGSPAVAVPTHSQPSP TPSNESTDTASEIGSAFNSPLRSPIRSANPTRPS SPVTSHISKVLFGEDDSLLRVDCIRYNRAVRDLG PVTSTGLLHLAEDGVLSPLALTEGGKGSSPSTRP IQGSQGSSSPVEKEVVEATDSREKTGMVRPGEPL SGEKYSPKELLALLKCVEAEIANYEACLKEEVEK RKKFKIDDQRRTHNYDEFICTFISMLAQEGMLAN LVEQNISVRRRQGVSIGRLHKQRKPDRRKRSRPY KAKRQ

Tag (bold text): N-terminal GST Protease cleavage site: PreScission™ (LEVLFQ ▼ GP) BAP1 (regular text): Start bold italics (amino acid residues 1-729) Accession number: AAH01596



ORDERS / SALES SUPPORT

International: +1-617-245-0003

US Toll-Free: 1-888-4E1E2E3 (1-888-431-3233) Email: sales.support@ubiquigent.com

UK HQ and TECHNICAL SUPPORT

International: +44 (0) 1382 381147 (9AM-5PM UTC) US/Canada: +1-617-245-0020 (9AM-5PM UTC) Email: tech.support@ubiquigent.com

Email services@ubiquigent.com for enquiries regarding compound profiling and/or custom assay development services.

© Ubiquigent 2012. Unless otherwise noted, Ubiquigent, Ubiquigent logo and all other trademarks are the property of Ubiquigent, Ltd.

Limited Terms of Use: For research use only. Not for use in humans or for diagnostics. Not for distribution or resale in any form, modification or derivative OR for use in providing services to a third party (e.g. screening or profiling) without the written permission of Ubiquigent, Ltd.

Lot-specific COA version tracker: v1.0.0