

BAP1 monoclonal antibody

Other names: HUCEP-13, UCHL2, TPDS

Cat. No. C15200212

Type: Monoclonal

IgG isotype: IgG1 kappa

Source: Mouse

Lot #: 001

Size: 50 µg/50 µl

Concentration: 1 µg/µl

Specificity: Human, mouse: positive. Other species: not tested.

Purity: Protein G purified polyclonal antibody in PBS containing 0.05% azide.

Storage: Store at -20°C; for long storage, store at -80°C. Avoid multiple freeze-thaw cycles.

Precautions: This product is for research use only. Not for use in diagnostic or therapeutic procedures.

Description: Monoclonal antibody raised in mouse against human BAP1 (BRCA1 Associated Protein-1 (Ubiquitin Carboxy-Terminal Hydrolase)) using the full length recombinant protein.

Applications

	Suggested dilution	Results
Western blotting	1:1,000 - 1:2,000	Fig 1, 2
IP	1:400	Fig 2
IF	1:1,000	Fig 3

Target description

BAP1 (UniProt/Swiss-Prot entry Q92560) belongs to the ubiquitin C-terminal hydrolase subfamily of deubiquitinating enzymes that are involved in the removal of ubiquitin from proteins. It plays a key role in chromatin structure by mediating deubiquitination of histone H2A monoubiquitinated at 'Lys-119' (H2AK119ub). Recent studies have shown that BAP1 is a member of the polycomb-group proteins (PcG) of highly conserved transcriptional repressors required for long-term silencing of genes that regulate cell fate determination, stem cell pluripotency, and other developmental processes.

Results

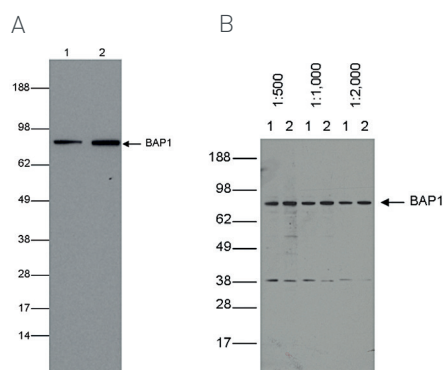


Figure 1. Western blot analysis using the Diagenode monoclonal antibody directed against BAP1

Figure 1A. Whole cell extracts from human NB4 cells (30 µg, lane 1) or mouse M1 cells (lane 2) were analysed by Western blot using the Diagenode antibody against BAP1 (Cat. No. C15200212), diluted 1:1,000 in PBS containing 10% milk. The position of the protein of interest (expected MW 80 kDa) is indicated on the right; the marker (in kDa) is shown on the left.

Figure 1B. Antibody titration on NB4 (1) and M1 (2) cells with the Diagenode BAP1 antibody.

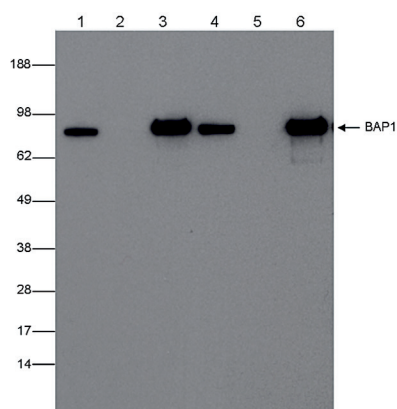


Figure 2. Immunoprecipitation using the Diagenode monoclonal antibody directed against BAP1

IP was performed on 1 mg RIPA cell lysate from human NB4 cells (lane 1-3) or mouse M1 cells (lane 4-6) using the Diagenode antibody against BAP1 (Cat. No. C15200212) diluted 1:400 (lane 3 and 6) or an IgG negative control (lane 2 and 5). The samples were analysed by Western blot analysis as described above. The input sample (30 µg RIPA lysate) was used as a positive control (lane 1 and 4).

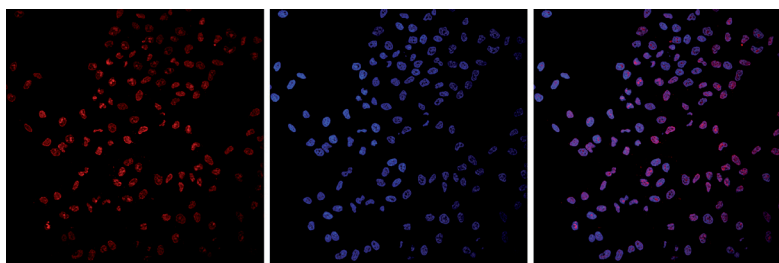


Figure 3. Immunofluorescence using the Diagenode monoclonal antibody directed against BAP1

HeLa cells were stained with the Diagenode antibody against BAP1 (Cat. No. C15200212) and with DAPI. Cells were fixed with 4% formaldehyde for 10' and blocked with PBS/TX-100 containing 1% BSA. The cells were immunofluorescently labelled with the BAP1 antibody (left) diluted 1:1,000 in blocking solution followed by an anti-mouse antibody conjugated to Alexa594. The middle panel shows staining of the nuclei with DAPI. A merge of the two stainings is shown on the right.

Diagenode sa. BELGIUM | EUROPE

LIEGE SCIENCE PARK
Rue Bois Saint-Jean, 3
4102 Seraing (Ougrée) - Belgium
Tel: +32 4 364 20 50
Fax: +32 4 364 20 51
orders@diagenode.com
info@diagenode.com

Diagenode Inc. USA | NORTH AMERICA

400 Morris Avenue, Suite 101
Denville, NJ 07834 - USA
Tel: +1 862 209-4680
Fax: +1 862 209-4681
orders.na@diagenode.com
info.na@diagenode.com

Last update: September, 5 2016