

UBXD7 (human; full length), pAb

Alternate Names: ABXN7, UBXN7, UBX domain protein 7, UBX domain containing 7

Cat. No. 68-0027-100
Lot. No. 30264

Quantity: 100 µg
Storage: -20°C

FOR RESEARCH USE ONLY

NOT FOR USE IN HUMANS

CERTIFICATE OF ANALYSIS

Page 1 of 2

This antibody was developed and validated by the Medical Research Council Protein Phosphorylation and Ubiquitylation Unit (University of Dundee, Dundee, UK).

Background

Ubiquitin signals are decoded in cells by at least 200 ubiquitin binding proteins, which interact with different types of polyubiquitin chains and ubiquitin-like modifiers. These interactions induce conformational changes that allow these proteins to transmit the ubiquitin signal to effector proteins (Dikic *et al.*, 2009). Cloning of the human UBX domain protein 7 (UBXD7) was first described by Seroussi *et al.* (1999). The UBX proteins bind ubiquitin conjugates as well as interact with E3 ubiquitin ligases. UBXD7 binds and interacts with several ubiquitin ligases including p97, an ATP-dependent chaperone involved in endoplasmic reticulum-associated degradation, and to the ubiquitin ligase CUL2/VHL and its substrate hypoxia-inducible factor 1α (HIF1α) (Alexandru *et al.*, 2008). UBXD7 has also been shown to directly interact with NEDDylated cullins and cause the accumulation of the Cullin2 substrate HIF1α (Bandua *et al.*, 2014).

Antibody Production:

Anti-UBXD7 (human) polyclonal antibody was raised in sheep against UBXD7 (residues 1-489 of human UBXD7). The antibodies were purified by the Medical Research Council Protein Phosphorylation and Ubiquitylation Unit (MRC-PPU, University of Dundee, Dundee, U.K.) by affinity purification of the anti-UBXD7 pAbs from the sheep serum using an antigen-

Physical Characteristics

Quantity: 100 µg

Concentration: to be provided on shipping

Source: sheep polyclonal antibody

Immunogen: human UBXD7 (residues 1-489) [GST-tagged]

Purification: affinity-purified using immobilized immunogen

Formulation: phosphate-buffered saline

Specificity: detects UBXD7 at ~55 kDa

Reactivity: human; other species not tested.

Stability/Storage: 12 months at -20°C; aliquot as required

Research Applications and Quality Assurance

Western Immunoblotting:
Dilute 1:3000

Immunoprecipitation:
Suitable for IP



Immunoprecipitation:

Immunoprecipitation of UBXD7 was performed on mouse brain lysate by capturing with 250 µg of anti-UBXD7 antibody bound to protein A agarose beads. UBXD7 was detected by Western blotting using (1:3000) of anti-UBXD7 antibody (Cat# 68-0027-100).

Continued on page 2



www.ubiquigent.com
Dundee, Scotland, UK

ORDERS / SALES SUPPORT

International: +1-617-245-0020
US Toll-Free: 1-888-4E1E2E3 (1-888-431-3233)
Email: sales.support@ubiquigent.com

UK HQ and TECHNICAL SUPPORT

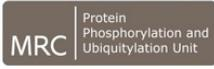
International: +44 (0) 1382 381147 (9AM-5PM UTC)
US/Canada: +1-617-245-0020 (9AM-5PM UTC)
Email: tech.support@ubiquigent.com

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

© Ubiquigent 2014. 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



UBXD7 (human; full length), pAb

Alternate Names: ABXN7, UBXN7, UBX domain protein 7, UBX domain containing 7

Cat. No. 68-0027-100
Lot. No. 30264

Quantity: 100 µg
Storage: -20°C

FOR RESEARCH USE ONLY

NOT FOR USE IN HUMANS

CERTIFICATE OF ANALYSIS

Page 2 of 2

Background

Continued from page 1

agarose column followed by depletion of any anti-GST pAbs using a GST-agarose column. Anti-UBXD7 (human) pAb was sourced by Ubiquigent directly from the MRC-PPU.

General References:

Alexandru G, Graumann J, Smith GT, Kolawa NJ, Fang R, Deshaies RJ (2008) UBXD7 binds multiple ubiquitin ligases and implicates p97 in H1-F1alpha turnover. *Cell* **5**, 804-16.

Bandau S, Knebel A, Gage ZO, Wood NT, Alexandru G (2012) UBXN7 docks on neddylated cullin complexes using its UIM motif and causes HIF1α accumulation. *BMC Biol* **10**, 36.

Dikic I, Wakatsuki S, Walters KJ (2009) Ubiquitin-binding domains - from structures to functions. *Nat Rev Mol Cell Biol* **10**, 659-671.

Seroussi E, Kedra D, Kost-Alimova M, Sandberg-Nordqvist AC, Fransson I, Jacobs JF, Fu Y, Pan HQ, Roe BA, Imreh S, Dumanski JP (1999) TOM1 genes map to human chromosome 22q13.1 and mouse chromosome 8C1 and encode proteins similar to the endosomal proteins HGS and STAM. *Genomics* **57**, 380-8.



www.ubiquigent.com
Dundee, Scotland, UK

ORDERS / SALES SUPPORT

International: +1-617-245-0020
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 2014. 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