

HOIL1 (human; full length), pAb

Alternate Names: Heme-oxidized IRP2 ubiquitin ligase 1, RanBP-type and C3HC4-type zinc finger-containing protein 1, HBV-associated factor 4, Hepatitis B virus X-associated protein 4, RING finger protein 54, Ubiquitin-conjugating enzyme 7-interacting protein 3

Cat. No. 68-0012-100
Lot. No. 30249

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

The linear ubiquitin chain assembly complex (LUBAC) mediates linear polyubiquitylation of proteins (Verhelst *et al.*, 2012) through ubiquitylation of the amino-terminal methionine of ubiquitin, repeated linear chain extension and attachment of such chains to the target substrate (Reiser *et al.*, 2012). It is an E3 ubiquitin ligase complex composed of a catalytic subunit HOIP (HOIL-1-interacting protein) and the two regulatory subunits HOIL-1 (heme-oxidized iron-regulatory protein 2 ubiquitin ligase-1) and SHARPIN (SHANK-associated RH domain-interacting protein) (Verhelst *et al.*, 2012; Tokunaga & Iwai, 2012). LUBAC plays an important role in TNF-induced NF-κB signalling (Haas *et al.*, 2009; Tokunaga *et al.*, 2009) and is involved in inflammatory responses, acquired and innate immunity, lymphocyte development, interferon production, the genotoxic stress response, and skeletal conditions. LUBAC has been implicated in various inflammatory, infectious and autoimmune diseases such as psoriasis-like dermatitis, rheumatoid arthritis, sepsis, and systemic lupus erythematosus (Tokunaga & Iwai, 2012). Various tumour tissues show enhanced SHARPIN expression which suggests a role for LUBAC in carcinogenesis (Jung *et al.*, 2010).

Continued on page 2

Physical Characteristics

Quantity: 100 µg

Formulation: phosphate-buffered saline

Concentration: to be provided on shipping

Specificity: detects HOIL1 at ~58 kDa

Source: sheep polyclonal antibody

Reactivity: human; other species not tested

Immunogen: human HOIL1 (residues 1-510) [His-tagged]

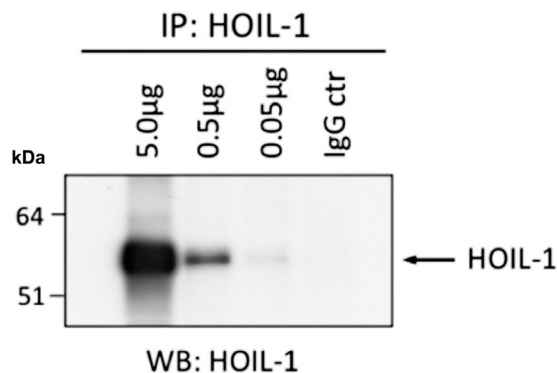
Stability/Storage: 12 months at -20°C; aliquot as required

Purification: affinity-purified using immobilized immunogen

Research Applications and Quality Assurance

Western Immunoblotting:
Use 1 µg/ml

Immunoprecipitation:
Use 5 µg/mg of cell extract



Immunoprecipitation Assay:

HOIL1 was immunoprecipitated from HeLa total cell extracts (1 mg) using various amounts of anti-HOIL1 antibody (Cat# 68-0012-100) or pre-immune serum (IgG). HOIL1 was subsequently detected by Western Blot using a commercially available anti-HOIL1 antibody.

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



HOIL1 (human; full length), pAb

Alternate Names: Heme-oxidized IRP2 ubiquitin ligase 1, RanBP-type and C3HC4-type zinc finger-containing protein 1, HBV-associated factor 4, Hepatitis B virus X-associated protein 4, RING finger protein 54, Ubiquitin-conjugating enzyme 7-interacting protein 3

Cat. No. 68-0012-100
Lot. No. 30249

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

Antibody Production:

Anti-HOIL1 (human) polyclonal antibody was raised in sheep against HOIL1 (residues 1-510 of human HOIL1). 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-HOIL1 pAbs from the sheep serum using a 6His-tagged-antigen-agarose column. Anti-HOIL1 (human) pAb was sourced by Ubiquigent directly from the MRC-PPU.

General References:

Haas TL, Emmerich CH, Gerlach B, Schmukle AC, Cordier SM *et al.* (2009) Recruitment of the linear ubiquitin chain assembly complex stabilizes the TNF-R1 signalling complex and is required for TNF-mediated gene induction. *Mol Cell* **36**, 831–844.

Jung JM, Kim B, Park Y, Cheon B *et al.* (2010) Newly identified tumor-associated role of human Sharpin. *Mol Cell Biochem* **340**, 161–167.

Rieser E, Cordier SM, Walczak H (2013) Linear ubiquitination: a newly discovered regulator of cell signalling. *Trends in Biochemical Sciences* **38**, 94–102.

Tokunaga F & Iwai K (2012) LUBAC, a novel ubiquitin ligase for linear ubiquitination, is crucial for inflammation and immune responses. *Microbes and Infection* **14**, 563–572.

Tokunaga F, Sakata S, Saeki Y, Satomi Y, Kirisako T *et al.* (2009) Involvement of linear polyubiquitylation of NEMO in NF-kappaB activation. *Nat Cell Biol* **11**, 123–132.

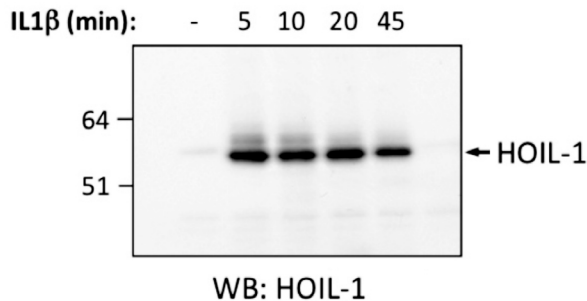
Verhelst K, Carpentier I, Kreike M, Meloni L, Verstrepen L, Kensche T, Dikic I, Beyaert R (2012) A20 inhibits LUBAC-mediated NF-κB activation by binding linear polyubiquitin chains via its zinc finger 7. *EMBO J* **31**, 3845–3855.

Application Reference:

Emmerich CH, Ordureau A, Strickson S, Arthur JSC, Pedriolo PGA, Komander D and Cohen P (2013) Activation of the canonical IKK complex by K63/M1-linked hybrid ubiquitin chains. *PNAS* **110**, 15247–52.

Research Applications and Quality Assurance

Continued from page 1



Western Blotting Analysis:

HEK293 IL-IR expressing cells were incubated with or without IL-1β for varying amounts of time, HOIL1 was precipitated from 3 mg cell lysates using immobilised NEMO (IKKγ); NEMO (Cat# 66-1002-050) captures linear and K63-linked ubiquitin chains. Western Blotting was carried out on eluted proteins using anti-HOIL1 antibody (Cat# 68-0012-100). The results show that NEMO captures HOIL1 from IL1β-stimulated and not unstimulated cells. HOIL1 was not captured when NEMO was replaced by the polyubiquitin-binding defective mutant NEMO D311N (data not shown; See Cat# 66-1013-050).



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