

TRIM65 (human; full length), pAb

Alternate Names: Tripartite Motif-Containing 65, 4732463G12Rik

Cat. No. 68-0024-100
Lot. No. 30261

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 enzymes of the ubiquitylation pathway play a pivotal role in a number of cellular processes including the regulated and targeted proteasome-dependent degradation of substrate proteins. Three classes of enzymes are involved in the process of ubiquitylation; activating enzymes (E1s), conjugating enzymes (E2s) and protein ligases (E3s). Tripartite Motif containing 65 (TRIM65) is a member of the E3 protein ligase family and cloning of the gene was first described by Strausberg *et al.* (2002). TRIM65, is a 517 amino acid protein that belongs to the TRIM/RING Finger-B Box-Coiled Coil (RBCC) family and is encoded by a gene located on human chromosome 17q25.1 (Jensen *et al.*, 2001; Fornage *et al.*, 2011). The TRIM family of proteins are characterized by a conserved TRIM domain that includes a RING finger domain, three zinc-binding domains, a coiled-coil region and a B box-type zinc Finger. Several of the TRIM proteins are upregulated by type I and type II interferons, which are essential for the development of pathogen-resistance (Nisole, *et al.*, 2005; Ozato *et al.*, 2008; McNab *et al.*, 2010).

Antibody Production:

Anti-TRIM65 (human) polyclonal antibody was raised in sheep against TRIM65 (residues 1-485)

Physical Characteristics

Quantity: 100 µg

Concentration: to be provided on shipping

Source: sheep polyclonal antibody

Immunogen: human TRIM65 (residues 1-485) [GST-tagged]

Purification: affinity-purified using immobilized immunogen

Formulation: phosphate-buffered saline

Specificity: detects TRIM65 at ~57 kDa

Reactivity: human; other species not tested.

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

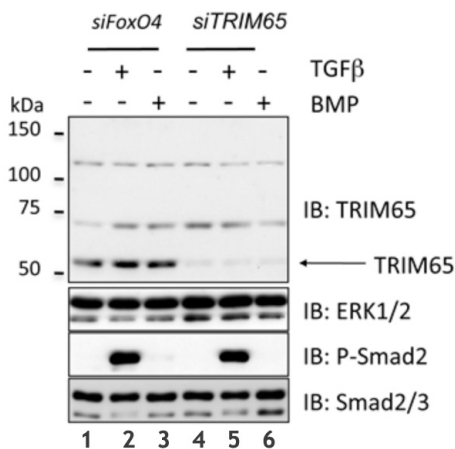
Research Applications and Quality Assurance

Western Immunoblotting:

Use 1.0 µg/ml

Immunoprecipitation:

Use 5.0 µg/mg of cell extract



Western Blotting Analysis:

By Western blotting TRIM65 was not detected in lysates from cells transfected with siTRIM65 (lanes 4-6) compared to those cells transfected with control siFoxO4 (lanes 1-3) where TRIM65 could be detected when probed with 0.1 µg/ml anti-TRIM65 antibody (Cat# 68-0024-100).

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

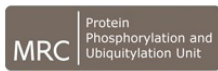
International: +44 (0) 1382 381147 (9AM-5PM UTC)
US/Canada: +1-617-245-0020 (9AM-5PM UTC)
Email: tech.support@ubiquigent.com

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



TRIM65 (human; full length), pAb

Alternate Names: Tripartite Motif-Containing 65, 4732463G12Rik

Cat. No. 68-0024-100
Lot. No. 30261

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

of human TRIM65). 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-TRIM65 pAbs from the sheep serum using an antigen-agarose column followed by depletion of any anti-GST pAbs using a GST-agarose column. Anti-TRIM65 (human) pAb was sourced by Ubiquigent directly from the MRC-PPU.

General References:

Fornage M, Debette S, Bis JC, Schmidt H, Ikram MA *et al.* (2011) Genome-wide association studies of cerebral white matter lesion burden: the CHARGE consortium. *Ann Neurol* **69**, 928-39.

Jensen K, Shiels C, Freemont PS (2001) PML protein isoforms and the RBCC/TRIM motif. *Oncogene* **20**, 7223-7233.

McNab FW, Rajsbaum R, Stoye JP, O'Garra A (2010) Tripartite-motif proteins and innate immune regulation. *Curr Opin Immunol* **23**, 46-56.

Nisole S, Stoye JP, Saib A (2005) TRIM family proteins: retroviral restriction and antiviral defence. *Nat Rev Microbiol* **3**, 799-808.

Ozato K, Shin DM, Chang TH, Morse HC (2008) TRIM family proteins and their emerging roles in innate immunity. *Nat Rev Immunol* **8**, 849-860.

Strausberg RL, Feingold EA, Grouse LH, Derge JG, Klausner RD (2002) Generation and initial analysis of more than 15,000 full-length human and mouse cDNA sequences. *PNAS* **99**, 16899-903.



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