



MYPT1 (human; residues 728 - 838), pAb

Alternate Names: Protein Phosphatase 1, Regulatory Subunit 12A, PPP1R12A, Myosin Phosphate target subunit 1

Cat. No. 68-0041-100
Lot. No. 30280

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

Protein ubiquitylation and protein phosphorylation are the two major mechanisms that regulate the functions of proteins in eukaryotic cells. However, these different post-translational modifications do not operate independently of one another, but are frequently interlinked to enable biological processes to be controlled in a more complex and sophisticated manner. Studying how protein phosphorylation events control the ubiquitin system and how ubiquitylation regulates protein phosphorylation has become a focal point of the study of cell regulation and human disease. The mammalian MYPT family consists of the products of five genes, denoted MYPT1, MYPT2, MBS85, MYPT3 and TIMAP (Grassie *et al.*, 2011). Myosin phosphatase (MP) activity, which regulates smooth muscle relaxation, is regulated by the phosphorylation of its regulatory subunit, myosin phosphatase targeting subunit 1 (MYPT1) (Cheng *et al.*, 2013). Cloning of human MYPT1 was first described by Takahashi *et al.* (1997). An example of the interplay between phosphorylation, ubiquitylation, and methylation, has been highlighted in a recent study showing that MYPT1 can be methylated *in vitro* and *in vivo* by histone lysine methyltransferase SETD7 and demethylated by histone demethylase LSD1. LSD1 silencing increased MYPT1 protein levels, decreasing the steady state level of phosphorylated retinoblastoma 1 (RB1; Ser

Continued on page 2

Physical Characteristics

Quantity: 100 µg

Concentration: to be provided on shipping

Source: sheep polyclonal antibody

Immunogen: human MYPT1 (residues 728-838) [GST-tagged]

Purification: affinity-purified using immobilized immunogen

Formulation: phosphate-buffered saline

Specificity: detects MYPT1 at ~ 115 kDa

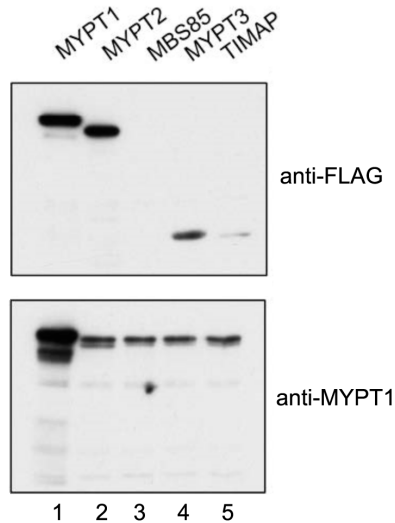
Reactivity: human; other species not tested

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

Research Applications and Quality Assurance

Western Immunoblotting: use 1 µg/ml

Immunoprecipitation: not tested



Western Blotting Analysis:

FLAG-tagged MYPT isoforms (Lanes 1-5) were overexpressed in HEK293 cells. By Western blotting using a commercially available anti-FLAG antibody, a band was observed in all lanes except Lane 3, where FLAG-MBS85 did not appear to express. By Western blotting the specific recognition of a band corresponding to MYPT1 was observed in all lanes when probed with the anti-MYPT1 antibody (Cat# 68-0041-100). No antibody cross reactivity was observed between the MYPT isoforms, this antibody recognises HEK293 endogenous MYPT1 as well as recombinant FLAG-tagged version.



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



MYPT1 (human; residues 728 - 838), pAb

Alternate Names: Protein Phosphatase 1, Regulatory Subunit 12A, PPP1R12A, Myosin Phosphate target subunit 1

Cat. No. 68-0041-100
Lot. No. 30280

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

807/811) and reducing E2F activity. Demethylation of MYPT1 has been shown to increase the ubiquitin proteasome pathway-dependent turnover of MYPT1. This study offers a novel cell cycle regulatory mechanism mediated by methylation/demethylation dynamics, and also reveals the significance of LSD1 overexpression in human carcinogenesis (Cho *et al.*, 2011).

Antibody Production:

Anti-MYPT1 (human) polyclonal antibody was raised in sheep against MYPT1 (residues 728-838 of human MYPT1). 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-MYPT1 pAbs from the sheep serum using an antigen-agarose column followed by depletion of any anti-GST antibodies using a GST-agarose column. Anti-MYPT1 (human) pAb was sourced by Ubiquigent directly from the MRC-PPU.

General References:

Cheng JC, Cheng HP, Tsai IC and Jiang MJ (2013) ROS-mediated downregulation of MYPT1 in smooth muscle cells: a potential mechanism for the aberrant contractility in atherosclerosis. *Lab Invest* **93**, 422-433.

Cho HS, Suzuki T, Dohmae N, Hayami S, Unoki M, Yoshimatsu M, *et al.* (2011) Demethylation of RB regulator MYPT1 by histone demethylase LSD1 promotes cell cycle progression in cancer cells. *Cancer Res* **71**, 655-660.

Grassie ME, Moffat LD, Walsh MP and MacDonald JA (2011) The myosin phosphatase targeting protein (MYPT) family: a regulated mechanism for achieving substrate specificity of the catalytic subunit of protein phosphatase type 1delta. *Arch Biochem Biophys* **510**, 147-159.

Takahashi N, Ito M, Tanaka J, Nakano T, Kaibuchi K, Odai H, *et al.* (1997) Localization of the gene coding for myosin phosphatase, target subunit 1 (MYPT1) to human chromosome 12q15-q21. *Genomics* **44**, 150-152.



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