



FAF1 (human; full length), pAb

Alternate Names: UBXD12, UBXN3A

Cat. No. 68-0011-100
Lot. No. 30248

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

Fas-associated factor 1 (FAF1) is a Fas-binding pro-apoptotic scaffolding protein that contains multiple protein interaction domains, including a FAS-interacting domain (FID), a death effector domain interacting domain (DEDID), and multi-ubiquitin-related domains (Menges *et al.*, 2009). FAF1 has been described as a component of the death-inducing signalling complex (DISC) in Fas-mediated apoptosis (De Zio *et al.*, 2008). Other DISC members include Fas-associated death domain protein (FADD) and caspase-8 (Ryu *et al.*, 2003). FAF1 has been shown to play an important role in normal development and neuronal cell survival, whereas FAF1 down regulation may contribute to multiple aspects of tumourigenesis. In particular, there is evidence implicating FAF1 as a tumour suppressor involved in the regulation of apoptosis and NF-κB activity, as well as in ubiquitylation and proteasomal degradation (Menges *et al.*, 2009). FAF1 is known to be associated with parkin, a key susceptibility protein for Parkinson's Disease (PD). Parkin acts as an E3 ubiquitin ligase to ubiquitylate FAF1 both *in vitro* and at the cellular level, identifying FAF1 as a direct substrate of parkin (Sul *et al.*, 2013).

Continued on page 2

Physical Characteristics

Quantity: 100 µg

Concentration: to be provided on shipping

Source: sheep polyclonal antibody

Immunogen: human FAF1 (residues 2-600) [His-tagged]

Purification: affinity-purified using immobilized immunogen

Formulation: phosphate-buffered saline

Specificity: detects FAF1 at ~74 kDa

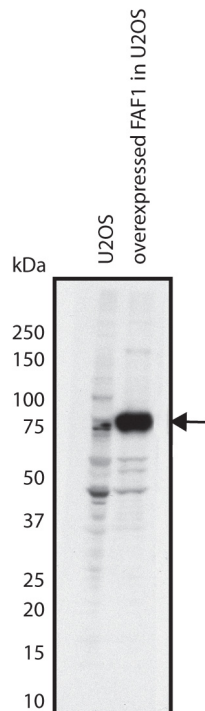
Reactivity: human; other species not tested

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

Research Applications and Quality Assurance

Western Immunoblotting:
Use 1 µg/ml

Immunoprecipitation:
Use 4 µg/mg of cell extract



Western Blotting Analysis:

Exogenously expressed FAF1 was detected in 40 µg U2OS total cell lysate by Western Blot using 1 µg/ml anti-FAF1 (Cat# 68-0011-100).



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



FAF1 (human; full length), pAb

Alternate Names: UBXD12, UBXN3A

Cat. No. 68-0011-100
Lot. No. 30248

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-FAF1 (human) polyclonal antibody was raised in sheep against FAF1 (residues 2-600 of human FAF1). 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-FAF1 pAbs from the sheep serum using a 6His-tagged-antigen-agarose column. Anti-FAF1 (human) pAb was sourced by Ubiquigent directly from the MRC-PPU.

General References:

De Zio D, Ferraro E, D'Amelio M, Simoni V, Bordini M, Soroldoni D, Berghella L, MeyerBI, Cecconi F (2008) Faf1 is expressed during neurodevelopment and is involved in Apaf1-dependent caspase-3 activation in proneural cells. *Cell Mol Life Sci* **65**, 1780-90.

Menges CW, Altomare DA, Testa JR (2009) FAS-associated factor 1 (FAF1): diverse functions and implications for oncogenesis. *Cell Cycle* **8**, 2528-34.

Ryu SW, Lee SJ, Park MY, Jun JI, Jung YK, Kim E (2003) Fas-associated factor 1, FAF1, is a member of Fas death-inducing signaling complex. *J Biol Chem* **278**, 24003-10.

Sul JW, Park MY, Juhee Shin J, Kim YR, Yoo SE, Kong YY, Kwon KS, Lee YH, Kim E (2013) Accumulation of the parkin substrate, FAF1, plays a key role in the dopaminergic neurodegeneration. *Hum Mol Genet* **22**, 1558-1573.



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