ATF6 monoclonal antibody (70B1413.1)

This antibody is covered by our Worry-Free Guarantee.

Citations: 30

View Online »

Ordering Information

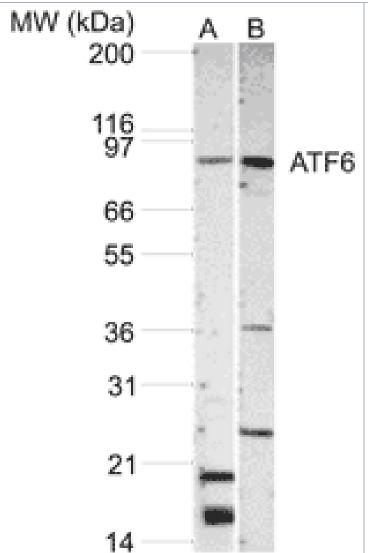
Order Online »

ALX-804-381-C100

100µg

Manuals, SDS & CofA

View Online »



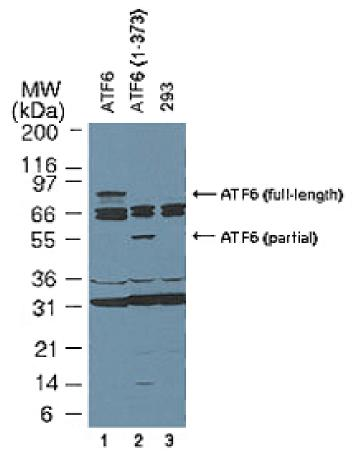
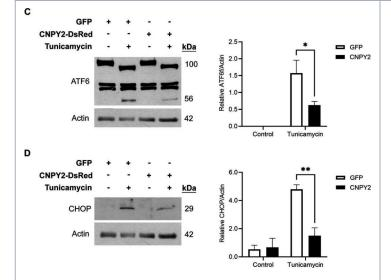


Figure 2: Western blot analysis of ATF6 using MAb to ATF6 (70B1413.1) (Prod. No. ALX-804-381). Method: Western blots were probed with 4μg/ml of MAb to ATF6 (70B1413.1), followed by an HRP-conjugated second step and visualized with PicoTect Western Blot Chemiluminescence Substrate. Film was exposed for 1 min. The top arrow corresponds to the ~90kDa form of ATF6 described as full-length in the literature. Lane 1: 293 cells transfected with full-length ATF6. Lane 2: 293 cells transfected with partial length ATF6 (aa 1-373). Lane 3: Untransfected 293 cells. The human full-length and partial length ATF6 plasmids are described in Luo and Lee (2002) [1].



CNPY2 regulates UPR signaling pathways. SH-SY5Y cells were transfected with control plasmid or with CNPY2-Myc-DDK (A,B) or CNPY2-dsRed (C,D) expressing plasmid for 24 h and cells were further treated with 2.5 µg/mL tunicamycin (Tun) for different times followed by immunoblotting as described above using β-actin (42 kDa) or GAPHD (37 kDa) as controls. Left, immunoblots, Right, quantification. Values are means \pm SD, n = 4–5. *p < 0.05 or **p < 0.01 for CNPY2 + Tun vs. Tun. (A,B) Tun was added for 6 h. The level of spliced XBP1 (XBP1s) was increased by Tun and was further elevated by CNPY2 expression (A). The level of phosphorylated eIF2α was increased by Tun but was not significantly affected by CNPY2 expression (B). (C) Tun was added for 16 h to study ATF6α signaling. The processing of ATF6α was induced by Tun as shown by the decrease in full length protein (100 kDa) and the appearance of the cleaved ATF6α fragment (56 kDa) in the immunoblots. CNPY2 expression reduced this ATF6α processing. (D) Tun was added for 24 h. The transcription factor CHOP (27 kDa) was induced by Tun and was decreased by CNPY2 expression.

Image collected and cropped by CiteAb under a CC-BY license from the following publication: CNPY2 protects against ER stress and is expressed by corticostriatal neurons together with CTIP2 in a mouse model of Huntington's disease. *Front Mol Neurosci* (2024)

Handling & Storage

Use/Stability Stable for 6 months when stored at +4°C.

Handling Avoid freeze/thaw cycles. After opening, prepare aliquots and store at -20°C.

Short Term Storage +4°C

Long Term Storage -20°C

Shipping Blue Ice

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name Activating transcription factor 6

Application ChIP, ICC, IHC (PS), IP, WB

Application Notes Detects bands of ~90kDa (full length ATF6) and ~50-

70kDa (cleaved forms of ATF6) by Western blot.

Clone 70B1413.1

Formulation Liquid. In PBS containing 0.05% BSA and 0.05% sodium

azide.

Host Mouse

Immunogen Partial human ATF6 (activating transcription factor 6)

containing aa 1-273.

lgG1

Purity Detail Protein G-affinity purified.

Recommendation Dilutions/Conditions Western Blot (1-5µg/ml)Suggested dilutions/conditions

may not be available for all applications. Optimal conditions

must be determined individually for each application.

Species Reactivity Hamster, Human, Mouse, Rabbit, Rat

P18850

Worry-free Guarantee

This antibody is covered by our Worry-Free Guarantee

Last modified: May 29, 2024



info-

ENZO LIFE SCIENCES, European Sales Office ENZO LIFE SCIENCES Phone: 800.942.0430 (ELS) AG Phone: +41 61 926 8989 usa@enzolifesciences.com eu@enzolifesciences.com Belgium, The Netherlands & Luxembourg Phone: +32 3 466 0420 infobe@enzolifesciences.com

Phone: +33 472 440 655 fr@enzolifesciences.com

Germany Phone: +49 7621 5500 526 de@enzolifesciences.com

UK & Ireland Phone (UK customers): 0845 601 1488 Phone: +44 1392 825900 info-uk@enzolifesciences.com