

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 »](#)

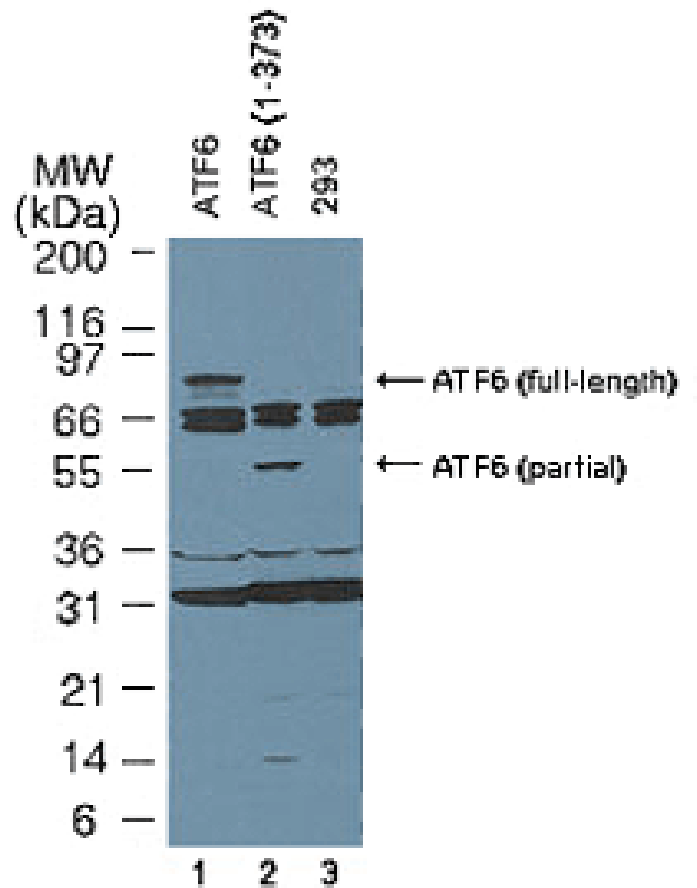
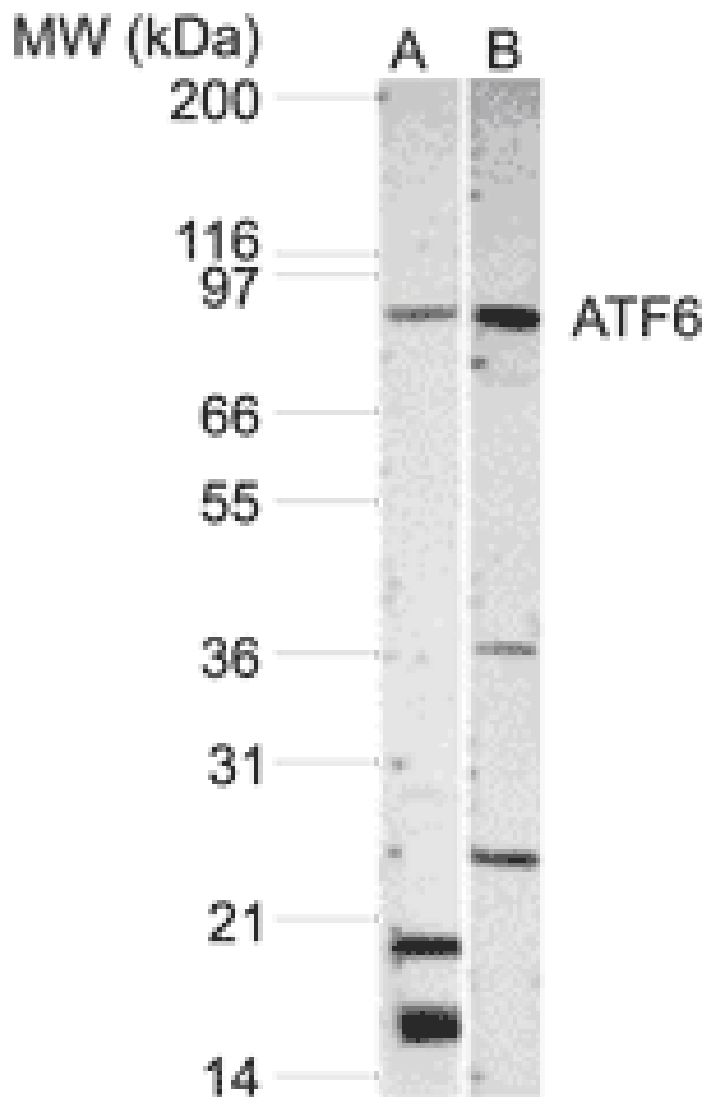
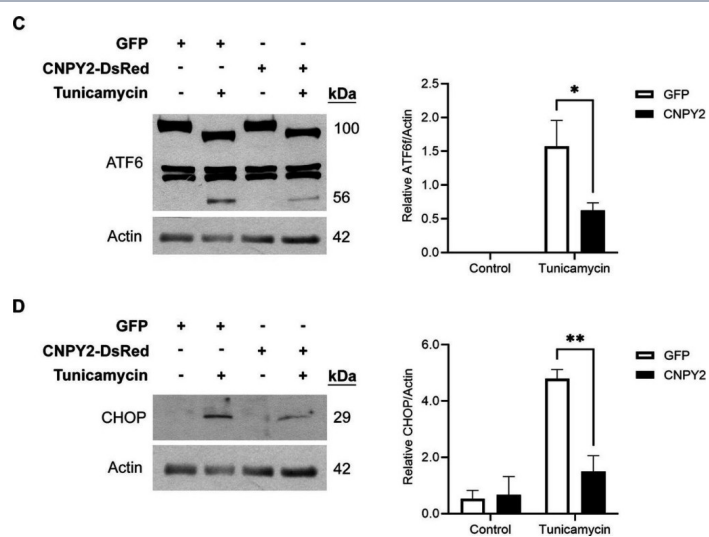


Figure 2: Western blot analysis of ATF6 using MAb to ATF6 (70B1413.1) (Prod. No. ALX-804-381). **Method:** Western blots were probed with 4µg/ml of MAb to ATF6 (70B1413.1), followed by an HRP-conjugated second step and visualized with PicoTect Western Blot Chemiluminescence Substrate. Film was exposed 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 μ g/mL tunicamycin (Tun) for different times followed by immunoblotting as described above using β -actin (42 kDa) or GAPDH (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 eIF2 α was increased by Tun but was not significantly affected by CNPY2 expression (B). (C) Tun was added for 16 h to study ATF6 α signaling. The processing of ATF6 α was induced by Tun as shown by the decrease in full length protein (100 kDa) and the appearance of the cleaved ATF6 α fragment (56 kDa) in the immunoblots. CNPY2 expression reduced this ATF6 α 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.
Isotype	IgG1
Purity Detail	Protein G-affinity purified.
Recommendation	Western Blot (1-5µg/ml)
Dilutions/Conditions	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



ENZO LIFE SCIENCES,
INC.
Phone: 800.942.0430
info-usa@enzolifesciences.com

P18850
ENZO LIFE SCIENCES
(ELS) AG
Phone: +41 61 926 8989
info-eu@enzolifesciences.com

Belgium, The Netherlands
& Luxembourg
Phone: +32 3 466 0420
info-fr@enzolifesciences.com

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

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

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

Worry-free Guarantee

This antibody is covered by our Worry-Free Guarantee.

For local distributors and detailed product information visit us online: enzo.com All product names, logos, and brands are the property of their respective owners.