

ATF6 monoclonal antibody (70B1413.1)

This antibody is covered by our [Worry-Free Guarantee](#).

Citations: 30

[View Online »](#)

Ordering Information

[Order Online »](#)

ALX-804-3810	100µg
--------------	-------

Manuals, SDS & CofA

[View Online »](#)

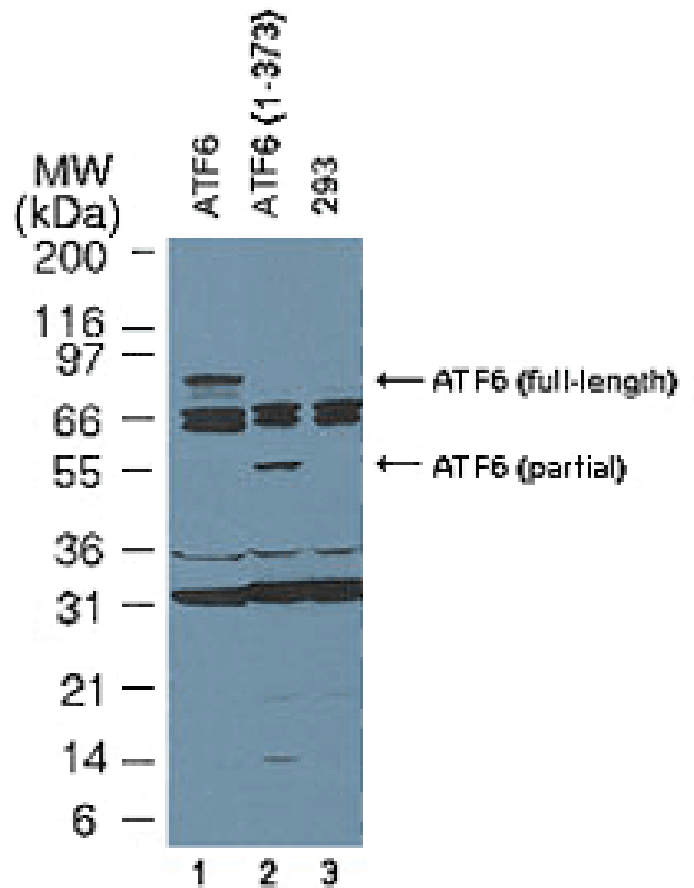
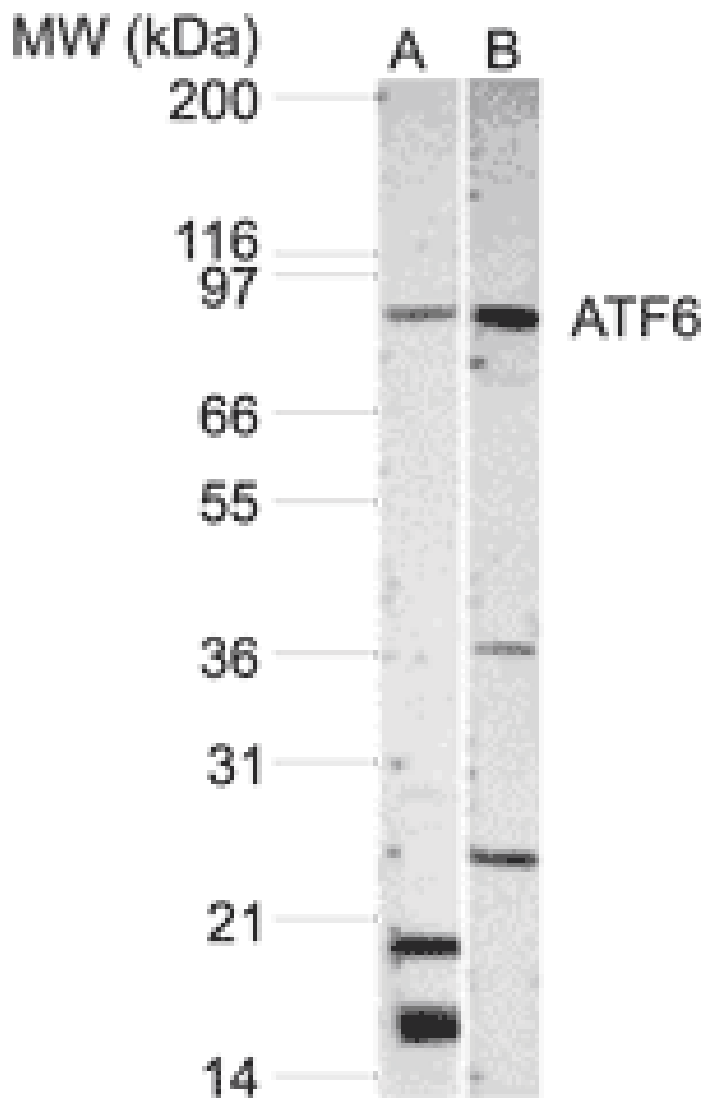
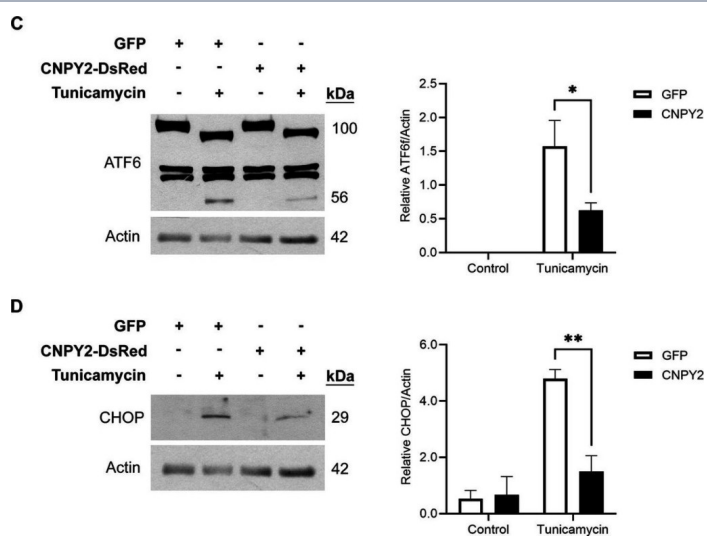


Figure 2: Western blot analysis of ATF6 using MAb to ATF6 (70B1413.1) (Prod. No. ALX-804-381). **Method:** Western blots were probed with 4 μ g/ml of MAb to ATF6 (70B1413.1), followed by an HRP-conjugated second step and visualized with PicoTect Western Blot Chemiluminescence Substrate. Film was exposed 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 $\mu\text{g}/\text{mL}$ 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 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

UniProt ID

P18850

Worry-free Guarantee

This antibody is covered by our [Worry-Free Guarantee](#)

Last modified: May 29, 2024



ENZO LIFE SCIENCES,
INC.
Phone: 800.942.0430
[info-
usa@enzolifesciences.com](mailto:info-usa@enzolifesciences.com)

European Sales Office
ENZO LIFE SCIENCES
(ELS) AG
Phone: +41 61 926 8989
[info-
eu@enzolifesciences.com](mailto:info-eu@enzolifesciences.com)

Belgium, The Netherlands
& Luxembourg
Phone: +32 3 466 0420
[info-
be@enzolifesciences.com](mailto:info-be@enzolifesciences.com)

France
Phone: +33 472 440 655
[info-
fr@enzolifesciences.com](mailto:info-fr@enzolifesciences.com)

Germany
Phone: +49 7621 5500 526
[info-
de@enzolifesciences.com](mailto:info-de@enzolifesciences.com)

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