

ATF6 polyclonal antibody

Disruptions of protein folding and maturation in the endoplasmic reticulum (ER) result in the activation of the unfolded protein response (UPR), an integrated cellular signaling pathway that transmits information from the ER lumen to the cytoplasm and nucleus. Activating transcription factor 6 (ATF6) as well as the ER-transmembrane protein kinases IRE1p and PERK are the major transducers of the UPR. ATF6 is an ER transmembrane protein that is normally bound to the ER chaperone Grp78, but upon ER stress is released from Grp78 and proteolytically cleaved to yield a cytosolic fragment which then migrates to the nucleus. Nuclear ATF6, together with the transcription factor XBP-1, activates transcription of UPR-responsive genes.

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

Citations: 2

[View Online »](#)

Ordering Information

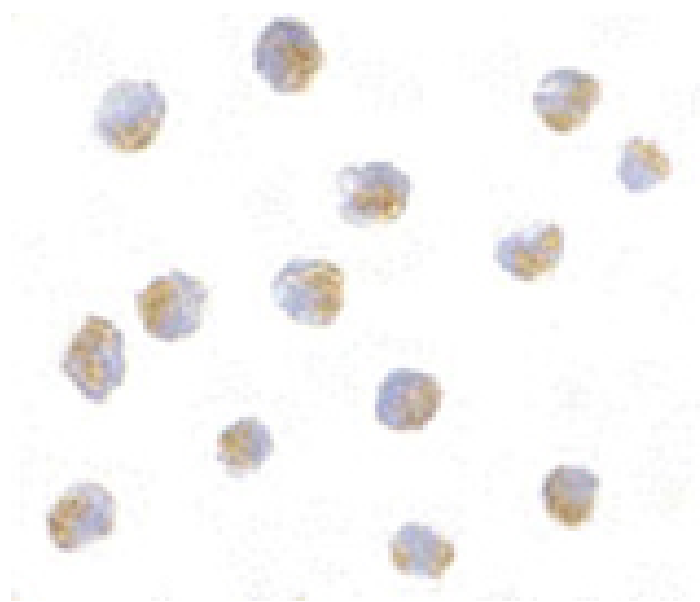
[Order Online »](#)

ADI-905-729-100

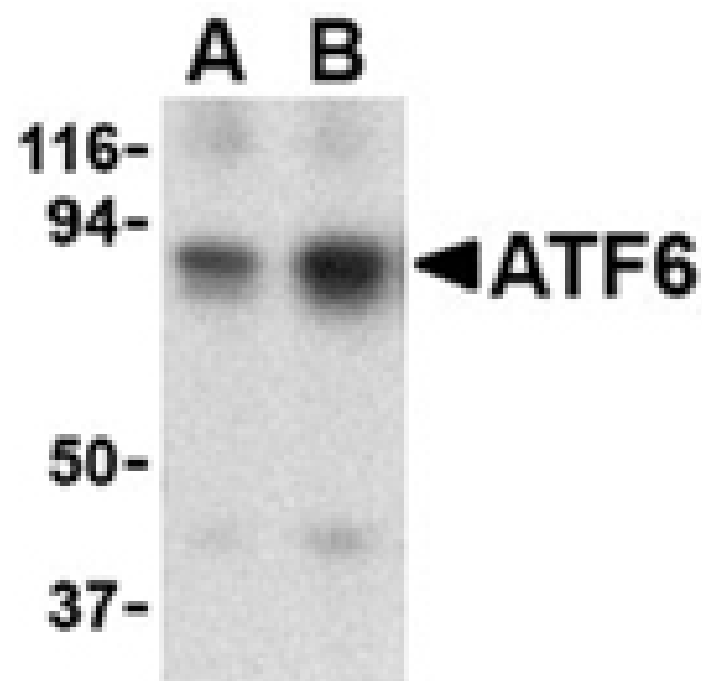
100µg

Manuals, SDS & CofA

[View Online »](#)



Immunocytochemistry analysis of ATF6 in A20 with ATF6 antibody at 10 μ g/ml.



Western blot analysis of ATF6 in MDA-MB-361 cell lysate with ATF6 antibody at (A) 0.5 and (B) 1 μ g/ml.

Handling & Storage

Long Term Storage -20°C

Shipping Blue Ice

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name Activating transcription factor 6

Application ICC, WB

Application Notes Detects a band of ~88kDa by Western blot.

Formulation Liquid. In PBS containing 0.02% sodium azide.

Host Rabbit

Immunogen Synthetic peptide corresponding to the sequence near the N-terminus of human ATF6.

Purity Detail Peptide affinity purified.

Recommendation Immunocytochemistry (10µg/ml)Western Blot (0.5-1.0µg/ml)
Dilutions/Conditions Suggested dilutions/conditions may not be available for all applications. Optimal conditions must be determined individually for each application.

Species Reactivity Human, Mouse

UniProt ID P18850

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



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

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

Belgium, The Netherlands
& Luxembourg
Phone: +32 3 466 0420
info-be@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