

XBP-1S monoclonal antibody (8F6.11)

X box binding protein 1 (XBP-1) is a key protein in the mammalian unfolded protein response (UPR) that protects the cell against the stress of misfolded proteins in the endoplasmic reticulum (ER). XBP-1 mRNA is spliced by IRE1p under ER stress, resulting in a 371 amino acid protein (XBP-1s) which is translocated to the nucleus where it binds to the regulatory elements of downstream genes. XBP-1 stimulates the production of ER stress proteins including the ER resident protein chaperones grp78 and grp94.

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

Citations: 1

[View Online »](#)

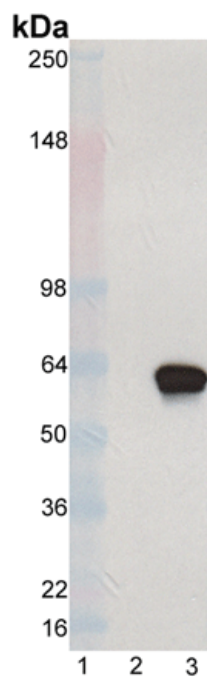
Ordering Information

[Order Online »](#)

ENZ-ABS255-0100	100µg
-----------------	-------

Manuals, SDS & CofA

[View Online »](#)



Western blot analysis of XBP-1S mAb (8F6.11) (Prod. No. ENZ-ABS255): Lane 1: MW markers, Lane 2: mock-transfected CHO cell, Lane 3: XBP-1S (spliced isoform)-transfected CHO cell. The observed molecular weight of XBP-1S is ~60 kDa (calculated MW: 40 kDa).

Handling & Storage

Handling	Avoid freeze/thaw cycles.
Long Term Storage	-20°C
Shipping	Blue Ice

Regulatory Status

RUO - Research Use Only

Product Details

Alternative Name	X-box binding protein 1, Tax-responsive element-binding protein 5, TREB5, XBP2
Application	ELISA, ICC, IP, WB
Application Notes	Detects a band of ~60kDa by Western blot.
Clone	8F6.11
Formulation	Liquid. In PBS containing 0.09% sodium azide.
Host	Mouse
Immunogen	Synthetic peptide corresponding to the sequence near the C-terminus of human XBP-1, spliced isoform XBP1(S).
Isotype	IgG1
Purity Detail	Protein G affinity purified.
Recommendation Dilutions/Conditions	Immunocytochemistry (1:100)Suggested dilutions/conditions may not be available for all applications.Optimal conditions must be determined individually for each application.
Source	Purified from concentrated hybridoma tissue culture supernatant.
Species Reactivity	Human, Mouse
UniProt ID	P17861

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

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