

# XBP-1 polyclonal antibody

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](#).

## Ordering Information

[Order Online »](#)

ENZ-ABS223-0100	100µg
-----------------	-------

## Manuals, SDS & CofA

[View Online »](#)

## 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** Flow Cytometry, ICC, WB

**Application Notes** Detects a band of ~60kDa by Western blot.

**Formulation** Liquid. In PBS containing 0.09% sodium azide.

**Host** Rabbit

**Immunogen** Synthetic peptide corresponding to the sequence near the C-terminus of human XBP-1, spliced isoform XBP1(S).

**Purity Detail** Peptide affinity purified.

**Recommendation Dilutions/Conditions** Flow Cytometry (1:50) 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 rabbit serum.

**Species Reactivity** Human, Mouse

**UniProt ID** P17861

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



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-  
fr@enzolifesciences.com](mailto:info-fr@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)