

# Bad monoclonal antibody (5E6)

Bad is an ~22 kDa pro-apoptotic protein that aids in the initiation of cell death. It exhibits sequence homology to Bcl-2, a protein that prevents apoptosis, within its BH1 and BH2 domains. Bad is phosphorylated at several sites including Ser112, Ser136; phosphorylation by kinases like Akt is known to alter the effects of Bad. Bad has been demonstrated to be phosphorylated at Ser170.

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

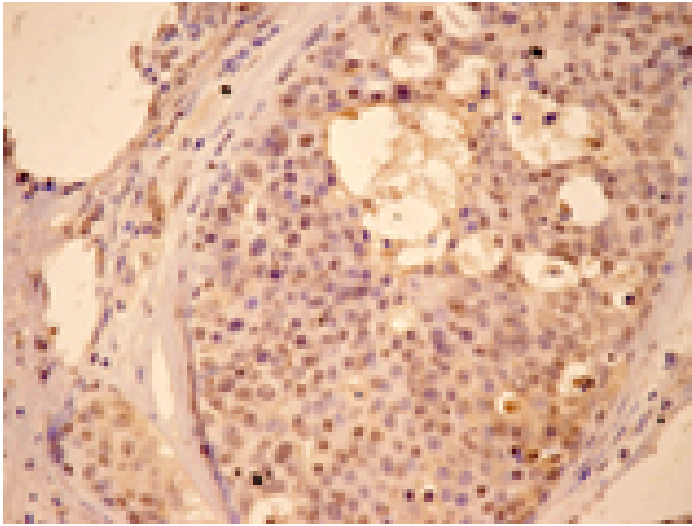
## Ordering Information

[Order Online »](#)

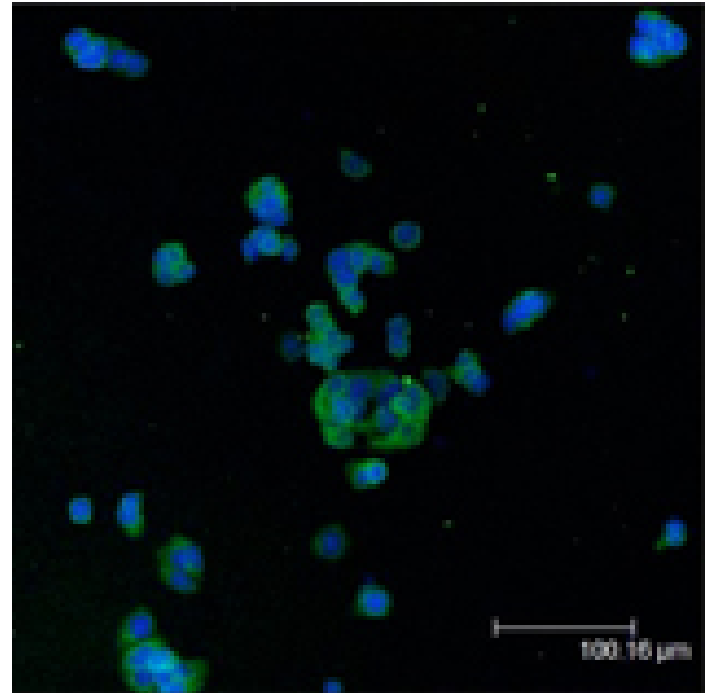
ADI-905-638-100	100µg
-----------------	-------

## Manuals, SDS & CofA

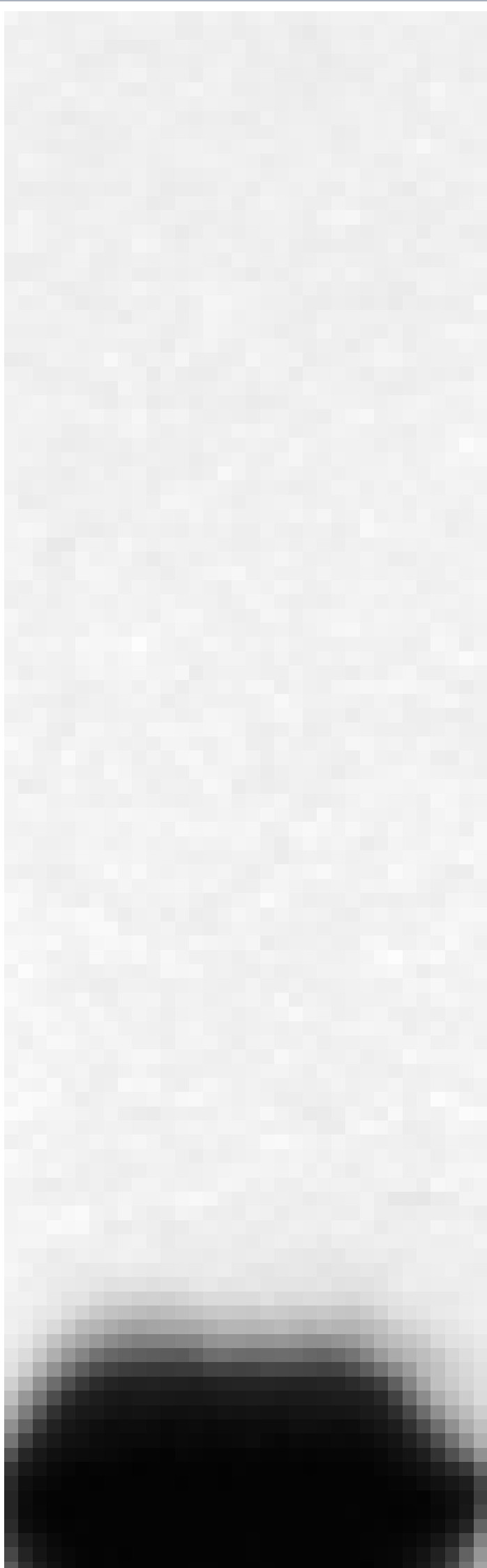
[View Online »](#)



Immunohistochemistry analysis of human breast cancer tissue with BAD mAb (5E6).



Immunofluorescent analysis (confocal) staining of MCF-7 cells using BAD mAb (5E6) (green); nuclei are stained in blue pseudocolor using DRAQ5.





# Handling & Storage

Use/Stability	Stable at -80°C up to 1 year, at 4°C up to 3 months.
Long Term Storage	-20°C
Shipping	Blue Ice

# Regulatory Status

RUO - Research Use Only

# Product Details

Alternative Name	BCL2 antagonist of cell death
Application	ICC, IHC, IP, WB
Application Notes	Detects a band of ~19kDa by Western blot.
Clone	5E6
Formulation	Lyophilized from 1ml of 2x PBS containing 0.09% sodium azide, PEG, and sucrose.
Host	Mouse
Immunogen	Recombinant human Bad.
Isotype	IgG1
Purity Detail	Thiophilic adsorption and size exclusion chromatography purified.
Recommendation Dilutions/Conditions	Immunohistochemistry (1:50)Immunoprecipitation (1-2µg/ml)Suggested dilutions/conditions may not be available for all applications.Optimal conditions must be determined individually for each application.
Reconstitution	Reconstitute with 1ml water (15 minutes at room temperature).
Species Reactivity	Human
UniProt ID	Q92934

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)