

BID monoclonal antibody (5C9)

BID is an approximately 22 kDa protein belonging to the Bcl-2 family. BID is a cytosolic protein that is cleaved by caspase-8 in response to death receptor stimulation. Proteolytic cleavage of BID results in the formation of a 15 kDa fragment, referred to as tBID. tBID migrates to the mitochondria where it aids in the release of cytochrome c and the subsequent activation of the apoptosome. tBID serves as a link between death receptor dependent and mitochondrial mediated apoptosis.

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

Ordering Information

[Order Online »](#)

ADI-AAM-141-E	100µg
---------------	-------

Manuals, SDS & CofA

[View Online »](#)

Handling & Storage

Long Term Storage -20°C

Shipping Blue Ice

Regulatory Status

RUO - Research Use Only

Product Details

Alternative Name	BH3 interacting domain death agonist
Application	WB
Application Notes	Detects a band of ~24kDa by Western blot.
Clone	5C9
Formulation	Liquid. In PBS containing 50% glycerol.
Host	Mouse
Immunogen	Synthetic peptide corresponding to a portion of human BID.
Isotype	IgG1
Purity Detail	Protein A 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	Human
UniProt ID	P55957

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
usa@enzolifesciences.com

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

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