

Malondialdehyde polyclonal antibody

MDA (Malondialdehyde) is a highly reactive three carbon dialdehyde produced as a byproduct of polyunsaturated fatty acid peroxidation and arachidonic acid metabolism. MDA readily combines with several functional groups on molecules including proteins, lipoproteins, and DNA. MDA modified proteins may show altered physico-chemical behavior and antigenicity. Antibodies to MDA will help to visualize the MDA-adducts.

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

Citations: 9

[View Online »](#)

Ordering Information

[Order Online »](#)

ALX-210-8790	100µl
--------------	-------

Manuals, SDS & CofA

[View Online »](#)

Handling & Storage

Use/Stability	Do not freeze, thaw or store diluted working solutions.
Handling	Avoid freeze/thaw cycles. After opening, prepare aliquots and store at -20°C.
Long Term Storage	-20°C
Shipping	Blue Ice

Regulatory Status

RUO - Research Use Only

Product Details

Alternative Name	MDA
Application	ELISA, IHC, WB
Formulation	Liquid. Contains 0.1% sodium azide.
Host	Rabbit
Immunogen	Malondialdehyde (MDA) conjugated to KLH.
Purity Detail	Unpurified
Recommendation Dilutions/Conditions	ELISA (1:1,000-1:10,000)Immunohistochemistry (1:500-1:1,000)Western blot (1:500-1:1,000)Suggested dilutions/conditions may not be available for all applications.Optimal conditions must be determined individually for each application.
Species Reactivity	Species independent
Specificity	Recognizes MDA adducts.
Worry-free Guarantee	This antibody is covered by our Worry-Free Guarantee