

# DMPO monoclonal antibody (1)

Oxidative stress resulting in the formation of biological radicals has been implicated in a number of human diseases, such as cancer and aging. There is, however, a paucity of reliable methods for *in vivo* or *ex vivo* detection of radical formation. Until now the only general technique that allowed for the detection of these highly reactive species was electron spin resonance (ESR) using spin traps. One of the most popular of these spin traps is 5,5-dimethyl-1-pyrroline N-oxide (DMPO). In the ESR method, radicals are trapped by DMPO, and the DMPO spin adduct signal is measured quantitatively by an ESR spectrometer. The DMPO, mAb (1) (Prod. No. ALX-803-340) specifically reacts with DMPO-protein and DMPO-DNA adducts and therefore highly simplifies the detection of radical formation compared to the ESR method.

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

Citations: 4

[View Online »](#)

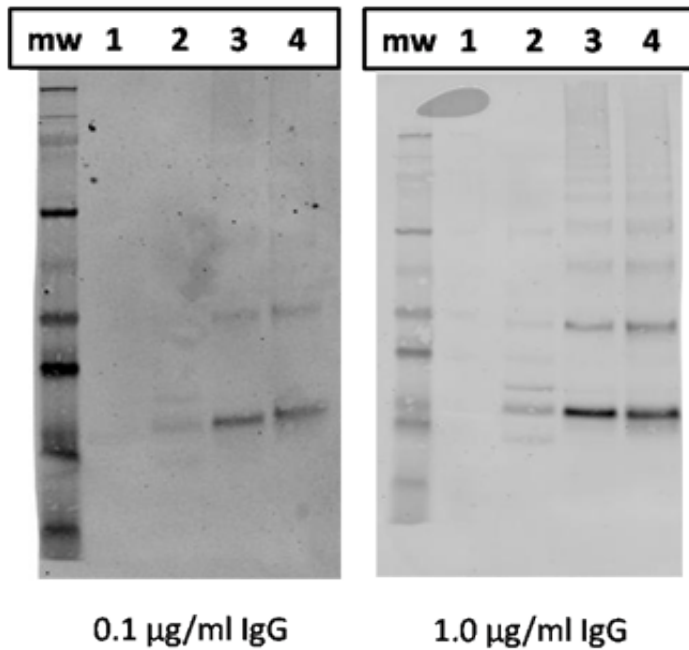
## Ordering Information

[Order Online »](#)

ALX-803-340-C100	100µg
------------------	-------

## Manuals, SDS & CofA

[View Online »](#)



**Figure:** Western Blot analysis using DMPO, mAb (1) (Prod. No. ALX-803-340).

**Method:** Samples are 2.5 µg each from reactions containing:

1= 10 µM Mb only

2= 10 µM Mb + 100 mM DMPO

3= 10 µM Mb + 100 µM H<sub>2</sub>O<sub>2</sub> + 100 mM DMPO

4= 10 µM Mb + 10 µM H<sub>2</sub>O<sub>2</sub> + 10 mM DMPO

Western conditions are:

-either 0.1 or 1.0 µg/ml anti-DMPO IgG for 1 hr at rt in 1XPBS/0.5% Tween

-wash 4X in 1XPBS/0.5% Tween

-1:10,000 anti-mouse IR800 for 1 hr at rt in 1XPBS/0.5% Tween

-wash 2X in 1XPBS/0.5% Tween, wash 2X in PBS

-rinse in H<sub>2</sub>O

-Scan on Licor Odyssey

## Handling & Storage

**Handling** Dilute only prior to immediate use.

**Long Term Storage** +4°C

**Shipping** Blue Ice

**Regulatory Status** RUO - Research Use Only

## Product Details

**Alternative Name** 5,5-dimethyl-1-pyrroline-N-oxide

**Application** ELISA, IP, WB

**Clone** 1

**Crossreactivity** Does not cross-react with non-adducted proteins or DNA. Does not cross-react with ovalbumin.

**Formulation** Liquid. In PBS, pH 7.4, containing 10% glycerol and 0.02% sodium azide.

**Host** Mouse

**Immunogen** DMPO-octanoic acid conjugated to ovalbumin.

**Isotype** IgG

**Purity Detail** Protein G-affinity purified.

**Recommendation** Western Blot (1-10ug/ml)ELISA (1-10ug/ml)Suggested dilutions/conditions may not be available for all applications.Optimal conditions must be determined individually for each application.

**Source** Hybridoma tissue culture supernatant.

**Species Reactivity** Species independent

**Specificity** Recognizes DMPO (Prod. No. ALX-430-090), DMPO-octanoic acid, DMPO-protein adducts and DMPO-DNA adducts.



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

**Worry-free Guarantee**

**This antibody is covered by our Worry-Free Guarantee.**