

Glutathione monoclonal antibody (D8)

Glutathione, a major intracellular non-protein thiol, protects against free radical damage by providing reducing equivalents for several key antioxidant enzymes. In addition, glutathione acts as a scavenger of hydroxyl radicals and nascent oxygen. Elevated glutathione levels are associated with a reduced rate of illness, higher levels of self-rated health, lower cholesterol, lower body mass index, and lower blood pressure among the elderly. Glutathione provides a primary defense system for the removal of oxidants in the brain. Studies have revealed a correlation between low glutathione levels and damage to neurons that manufacture dopamine, suggesting a link to Parkinson's disease.

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

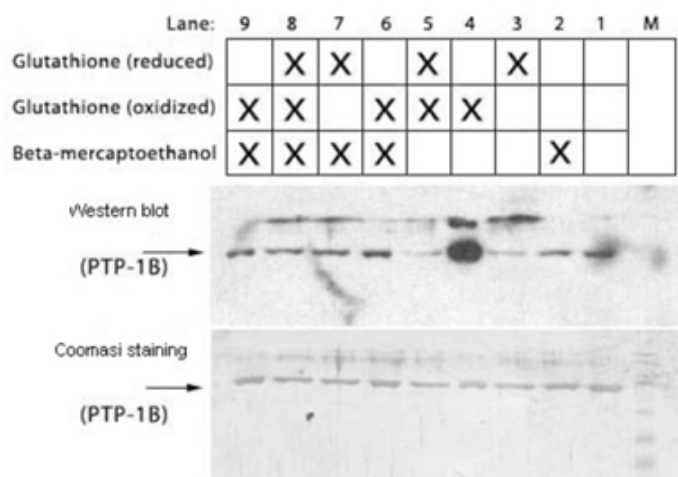
Ordering Information

[Order Online »](#)

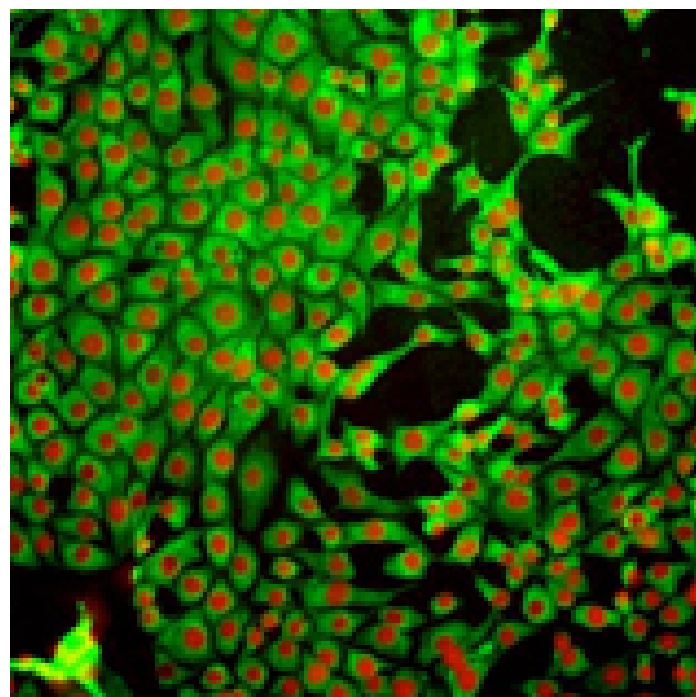
ADI-SPA-542-E	100µg
---------------	-------

Manuals, SDS & CofA

[View Online »](#)



Western blot analysis of Detection of protein-glutathione adducts on Western blot under non-reducing conditions. PTP-1B, phosphotyrosine phosphatase 1B.



Immunofluorescent analysis of HeLa cells with Glutathione mAb, clone D8, at 1:100 dilution. Red is DAPI nuclear stain.

Handling & Storage

Long Term Storage -20°C

Shipping Blue Ice

Regulatory Status

RUO - Research Use Only

Product Details

Application	ELISA, IP, WB
Clone	D8
Formulation	Liquid. In PBS, pH 7.2, containing 0.01% sodium azide.
Host	Mouse
Isotype	IgG2a
Purity Detail	Protein A affinity purified.
Recommendation Dilutions/Conditions	Western Blot (1:1,000, non-reducing)Suggested dilutions/conditions may not be available for all applications.Optimal conditions must be determined individually for each application.
Species Reactivity	Species independent
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-be@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)