

Glutathione peroxidase activity kit

The Glutathione Peroxidase activity kit is a convenient 96-well plate based assay for the higher throughput analysis of glutathione peroxidase activity in a variety of sample matrices in just 30 minutes.

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.

Citations: 28

[View Online »](#)

Ordering Information

[Order Online »](#)

ADI-900-158	5x96 wells
-------------	------------

Manuals, SDS & CofA

[View Online »](#)

- Validated in a variety of sample types
- Fast, accurate assay results in less than 30 minutes
- Convenient 96 well plate format for higher throughput.

Handling & Storage

Long Term Storage -20°C

Shipping Blue Ice

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name GPx

Application Activity assay, Colorimetric detection

Application Notes For the measurement of glutathione peroxidase in cell lysates, plasma, and tissue from any species.

Contents Microtiter plate, Assay buffer concentrate, Reaction mix, Glutathione reductase, GSSG

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)