

DNA damage ELISA kit

Rapid DNA damage ELISA kit for cancer, apoptosis and oxidative stress.

The DNA Damage EIA kit is a colorimetric competitive enzyme immunoassay kit with results in < 2.5 hours. This kit has been validated in-house in a variety of sample matrices and is supplied in a convenient 96 well plate format. Protocol (xls) available upon request.

Exposure of cells to oxidative and environmental stresses frequently results in the breakdown or oxidation of genomic DNA. Assays to evaluate the integrity of genomic DNA, or to assess the presence of oxidized DNA are frequently used as a means of verifying the onset of apoptosis or DNA damage.

Citations: 80

[View Online »](#)

Ordering Information

[Order Online »](#)

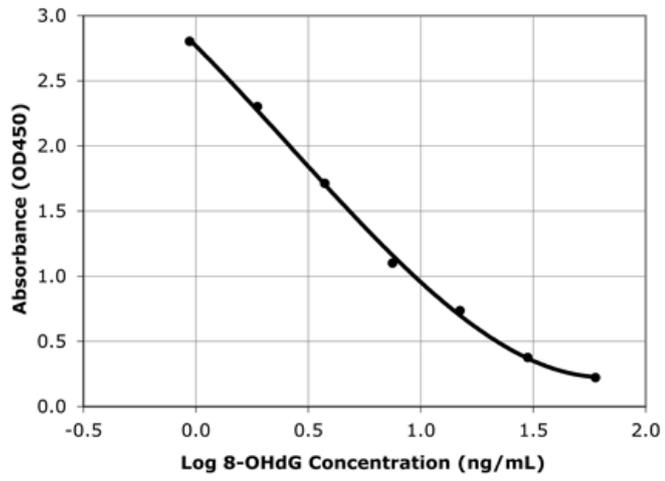
ADI-EKS-350

96 wells

Manuals, SDS & CofA

[View Online »](#)

- Highly sensitive measurements, detecting as low as 1 ng/ml
- Reliable and validated in-house in a variety of sample matrices
- Fully quantitative results that surpass semi-quantitative Western blot analysis
- No edge effect allowing you to assay more samples per kit



Standard curve of DNA damage ELISA kit

Handling & Storage

Use/Stability Store all components at 4°C, except standard at -20°C.

Shipping Dry Ice

Regulatory Status RUO - Research Use Only

Product Details

Application Colorimetric detection, ELISA

Application Notes For the measurement of DNA Damage (8-OHdG) in serum, saliva, and urine from any species. Cited sample types include plasma, seminal fluids, cell lysates, culture supernatants and DNA extracts.

Compatibility This product is compatible with the [Absorbance 96 Plate Reader](#).

Contents Microtiter plate, Standard, Sample diluent, Wash buffer concentrate, Antibody, Antibody diluent, Conjugate, Conjugate diluent, TMB Substrate, Stop solution 2

Crossreactivity The DNA Damage ELISA also detects 8-hydroxyguanosine (product of oxidative RNA damage) and 8-hydroxyguanine (product of oxidative DNA damage by hydroxyl radicals).

Sensitivity 0.59 ng/ml (range 0.94 – 60 ng/ml)

Species Reactivity Species independent

Wavelength 450 nm

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)