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 inhouse 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 »

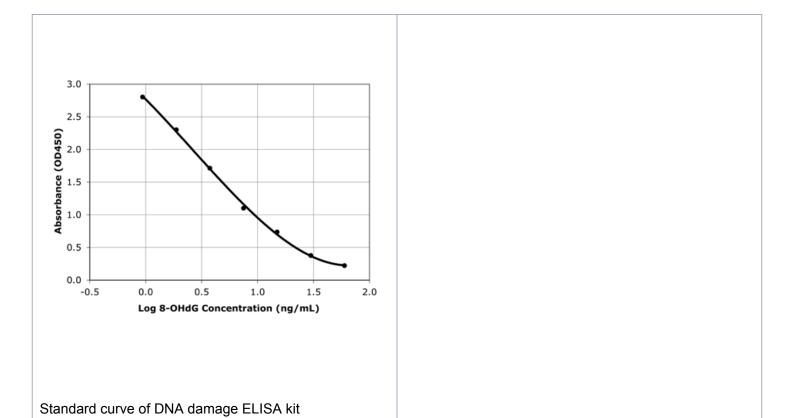
ADI-EKS-350

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



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 NotesFor 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.

Assay Time <2.5 hours

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 infousa@enzolifesciences.com European Sales Office ENZO LIFE SCIENCES (ELS) AG Phone: +41 61 926 8989 infoeu@enzolifesciences.com Belgium, The Netherlands & Luxembourg Phone: +32 3 466 0420 infobe@enzolifesciences.com

France
Phone: +33 472 440 655
infofr@enzolifesciences.com

Germany
Phone: +49 7621 5500 526
infode@enzolifesciences.com

UK & Ireland
Phone (UK customers):
0845 601 1488
Phone: +44 1392 825900
infouk@enzolifesciences.com