

OxiSelect™ UV-Induced DNA Damage ELISA Kit (CPD Quantitation)

Our OxiSelect™ UV-Induced DNA Damage ELISA Kit measures the formation of cyclobutane pyrimidine dimers (CPD) in DNA isolated from cells or tissues. Standards or unknown DNA samples are first heat denatured before adsorbed onto a 96-well DNA binding plate. The sample or standard are then probed with an anti-CPD antibody, followed by an HRP conjugated secondary antibody.

Ordering Information

[Order Online »](#)

CBI-STA-322	96 assays
-------------	-----------

Manuals, SDS & CofA

[View Online »](#)

- Measure CPD structures of DNA lesions induced by UV light
- Detect structures in isolated DNA using standard ELISA plate format

Handling & Storage

Use/Stability Upon receipt, store the CPD-DNA standard at -20°C and all other components at 4°C.

Shipping Ambient Temperature

Regulatory Status RUO - Research Use Only

Product Details

Application Colorimetric detection, ELISA

Technical Info / Product Notes For the Original Manufacturer's data sheet please [click here](#).



ENZO LIFE SCIENCES,
INC.
Phone: 800.942.0430
info-usa@enzolifesciences.com

European Sales Office
ENZO LIFE SCIENCES
(ELS) AG
Phone: +41 61 926 8989
info-eu@enzolifesciences.com

Belgium, The Netherlands
& Luxembourg
Phone: +32 3 466 0420
info-be@enzolifesciences.com

France
Phone: +33 472 440 655
info-fr@enzolifesciences.com

Germany
Phone: +49 7621 5500 526
info-de@enzolifesciences.com

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