

OxiSelect™ UV- Induced DNA Damage ELISA Combo Kit (CPD / 6-4PP Quantitation)

Our OxiSelect™ UV-Induced DNA Damage ELISA Combo Kit measures the formation of both CPD and 6-4PP structures 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-6-4PP or CPD antibody, followed by an HRP conjugated secondary antibody. CPD and 6-4PP structures are assayed on the same plate but in separate wells.

Ordering Information

[Order Online »](#)

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

Manuals, SDS & CofA

[View Online »](#)

- Combo kit measures both CPD and 6-4PP products on a single plate, but in separate wells
- Detects UV-damage structures from isolated DNA in standard ELISA plate format

Handling & Storage

Use/Stability Upon receipt, store the CPD-DNA and 6-4PP-DNA standards 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