

**NOTE: Revision to
Format and Assay Protocol**

Product Manual

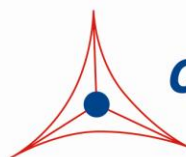
OxiSelect™ UV-Induced DNA Damage ELISA Kit (CPD / 6-4PP Quantitation)

Catalog Number

STA-322-C

96 assays

**FOR RESEARCH USE ONLY
Not for use in diagnostic procedures**



CELL BIOLABS, INC.
Creating Solutions for Life Science Research

Introduction

Absorption of ultraviolet (UV) light produces two predominant types of DNA damage, cyclobutane pyrimidine dimers (CPD) and pyrimidine (6-4) pyrimidone photoproducts (6-4PP) (Figure 1). The result is a transition of C to T and CC to TT, which are the most frequent mutations of p53 in both human and mouse skin cancers. UV damaged DNA is usually repaired by nucleotide excision repair (NER) or base excision repair (BER). After UV exposure, cells activate p53 and stall the cell cycle for repair. If the damage is too severe, the cell will trigger apoptosis to get rid of DNA damaged, potentially mutant cells.

Cell Biolabs' OxiSelect™ Oxidative UV-induced DNA Damage ELISA Combo Kit (CPD / 6-4PP Quantitation) is an enzyme immunoassay developed for rapid detection and quantitation of CPD or 6-4PP in any DNA samples. The quantity of CPD or 6-4PP in unknown sample is determined by comparing its absorbance respectively with that of a known CPD-DNA or 6-4PP-DNA standard curve. CPD and 6-4PP are assayed individually in separate wells, and each kit provides sufficient reagents to perform up to a total of 96 assays for CPD and 6-4PP combined, including standard curve and unknown samples. The 96 wells may be divided between CPD and 6-4PP assays at any ratio (e.g. 64 wells for CPD and 32 wells for 6-4PP, or 48 wells for each target).

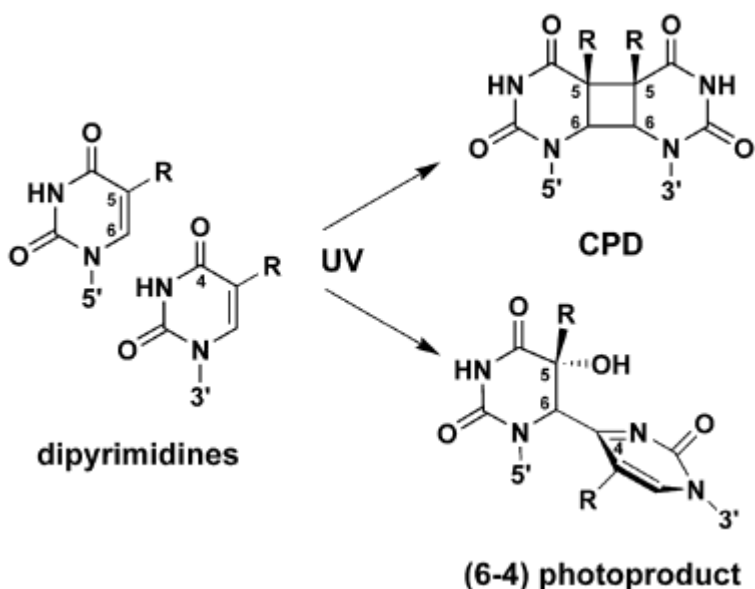


Figure 1: Structures of DNA lesions induced by UV Light

Assay Principle

CPD or 6-4PP-DNA standards or unknown DNA samples are first heat denatured before being adsorbed onto a 96-well DNA high-binding plate. The CPD or 6-4PPs present in the sample or standard are probed with an anti-CPD or anti-6-4PP antibody, followed by an HRP conjugated secondary antibody. The CPD or 6-4PP content in an unknown sample is determined by comparing with a standard curve that is prepared from predetermined CPD-DNA or 6-4PP-DNA standards.

Related Products

1. STA-320: OxiSelect™ Oxidative DNA Damage ELISA Kit (8-OHdG Quantitation)
2. STA-322: OxiSelect™ UV-induced DNA Damage ELISA Kit (CPD Quantitation)
3. STA-323: OxiSelect™ UV-induced DNA Damage ELISA Kit (6-4PP Quantitation)
4. STA-324: OxiSelect™ Oxidative DNA Damage Quantitation Kit (AP sites)
5. STA-325: OxiSelect™ Oxidative RNA Damage ELISA Kit (8-OHG Quantitation)
6. STA-326: OxiSelect™ Cellular UV-induced DNA Damage ELISA Kit (CPD)
7. STA-327: OxiSelect™ Cellular UV-induced DNA Damage Staining Kit (CPD)
8. STA-328: OxiSelect™ Cellular UV-induced DNA Damage ELISA Kit (6-4PP)
9. STA-329: OxiSelect™ Cellular UV-induced DNA Damage Staining Kit (6-4PP)

Kit Components

1. DNA High-Binding Plate (Part No. 232404): One 96-well strip plate.
2. DNA Binding Solution (Part No. 232405): One 6 mL bottle.
3. Anti-CPD Antibody (Part No. 232204): One 10 µL vial.
4. Anti-6-4PP Antibody (Part No. 232205): One 10 µL vial.
5. Secondary Antibody, HRP Conjugate (Part No. 10902): One 50 µL vial.
6. Assay Diluent (Part No. 310804): One 50 mL bottle.
7. 10X Wash Buffer (Part No. 310806): One 100 mL bottle.
8. Substrate Solution (Part No. 310807): One 12 mL amber bottle.
9. Stop Solution (Part. No. 310808): One 12 mL bottle.
10. CPD-DNA Standard (Part No. 232203): One 100 µL vial of 25 µg/mL CPD-DNA in 1X TE Buffer.
11. 6-4PP-DNA Standard (Part No. 232302): One 100 µL vial of 25 µg/mL 6-4PP-DNA in 1X TE Buffer.
12. Reduced DNA Standard (Part No. 232207): One 100 µL vial of 0.2 mg/mL reduced DNA in TE Buffer.

Materials Not Supplied

1. DNA samples such as cell or tissue genomic DNA
2. DNA Extraction Kit
3. Heating Block
4. PBS
5. 1X TE Buffer (10 mM Tris, pH 8.0, 1 mM EDTA)

6. 10 μL to 1000 μL adjustable single channel micropipettes with disposable tips
7. 50 μL to 300 μL adjustable multichannel micropipette with disposable tips
8. Multichannel micropipette reservoir
9. Microplate reader capable of reading at 450 nm (620 nm as optional reference wave length)

Storage

Upon receipt, aliquot and store the Reduced DNA, CPD-DNA, and 6-4PP-DNA Standards at -20°C to avoid multiple freeze/thaw cycles. Store all other components at 4°C .

Preparation of Reagents

- 1X Wash Buffer: Dilute the 10X Wash Buffer Concentrate to 1X with deionized water. Stir to homogeneity.
- Anti-CPD Antibody, Anti-6-4PP Antibody and Secondary Antibody: Immediately before use dilute the Anti-CPD or Anti-6-4PP Antibody 1:1000 and Secondary Antibody 1:1000 with Assay Diluent. Do not store diluted solutions.

Preparation of Standard Curve

1. Convert CPD-DNA (25 $\mu\text{g}/\text{mL}$), 6-4PP-DNA standard (25 $\mu\text{g}/\text{mL}$), and Reduced DNA (200 $\mu\text{g}/\text{mL}$) to single-stranded DNA by incubating the DNA at 95°C for 10 minutes and rapidly chilling on ice for 10 minutes.

Note: Aliquot and store denatured DNA at -20°C . Repeat the above denaturation step every time you prepare the CPD-DNA or 6-4PP-DNA standard.

2. Freshly prepare 4 $\mu\text{g}/\text{mL}$ of CPD-DNA or 6-4PP-DNA by diluting the denatured 25 $\mu\text{g}/\text{mL}$ stock in cold TE Buffer. Example: Add 8 μL to 42 μL of cold TE Buffer.
3. Freshly prepare 4 $\mu\text{g}/\text{mL}$ of Reduced DNA by diluting the denatured 200 $\mu\text{g}/\text{mL}$ stock in cold TE Buffer. Example: Add 40 μL to 1.96 mL of cold TE Buffer.
4. Prepare a series of CPD-DNA or 6-4PP-DNA standards according to Table 1.

Standard Tubes	4 $\mu\text{g}/\text{mL}$ Denatured CPD-DNA or 6-4PP-DNA (μL)	4 $\mu\text{g}/\text{mL}$ Denatured Reduced DNA (μL)	CPD-DNA or 6-4PP-DNA Conc. (ng/mL)
1	10	390	100
2	200 of tube #1	200	50
3	200 of tube #2	200	25
4	200 of tube #3	200	12.5
5	200 of tube #4	200	6.25
6	200 of tube #5	200	3.13
7	200 of tube #6	200	1.56
8	0	200	0

Table 1. Preparation of CPD-DNA or 6-4PP-DNA Standards

Assay Protocol

1. Extract DNA from cell or tissue samples using a commercial DNA Extraction kit or other desired method.
2. Convert DNA sample to single-stranded DNA by incubating the sample at 95°C for 10 minutes and rapidly chilling on ice for 10 minutes.
3. Dilute DNA samples to 4 µg/mL in cold TE Buffer.
Note: Samples with high concentrations of CPD or 6-4PP may be further diluted 2-4 fold in 4 µg/mL Reduced DNA. A titration may be performed to ensure the samples fall in the range of the standard curve.
4. Add 50 µL of unknown DNA samples, CPD-DNA, or 6-4PP-DNA standards to the wells of the DNA High-Binding plate.
5. Add 50 µL of DNA Binding Solution to each well. Mix well by pipetting and incubate at room temperature overnight on an orbital shaker. Each DNA sample including unknown and standard should be assayed in duplicate.
6. Remove the DNA solutions and wash twice with PBS. Blot plate on paper towels to remove excess fluid. Add 200 µL of Assay Diluent to each well and block for 1 hour at room temperature.
7. Remove the Assay Diluent. Blot plate on paper towels to remove excess fluid.
8. Add 100 µL of the diluted Anti-CPD or Anti-6-4PP Antibody to all wells and incubate for 1 hour at room temperature on an orbital shaker.
9. Wash 5 times with 250 µL of 1X Wash Buffer with thorough aspiration between each wash. After the last wash, empty wells and tap microwell strips on absorbent pad or paper towel to remove excess 1X Wash Buffer.
10. Add 100 µL of the diluted Secondary Antibody-HRP Conjugate to all wells and incubate for 1 hour at room temperature on an orbital shaker. Wash the strip wells 5 times according to step 9 above.
11. Warm Substrate Solution to room temperature. Add 100 µL of Substrate Solution to each well, including the blank wells. Incubate at room temperature on an orbital shaker. Actual incubation time may vary from 2-30 minutes.
Note: Watch plate carefully; if color changes rapidly, the reaction may need to be stopped sooner to prevent saturation.
12. Stop the enzyme reaction by adding 100 µL of Stop Solution to each well. Results should be read immediately (color will fade over time).
13. Read absorbance of each well on a microplate reader using 450 nm as the primary wave length. Use the Reduced DNA Standard as an absorbance blank.

Example of Results

The following figures demonstrate typical Oxidative UV-induced DNA Damage ELISA (CPD and 6-4PP Quantitation) results. One should use the data below for reference only. This data should not be used to interpret actual results.

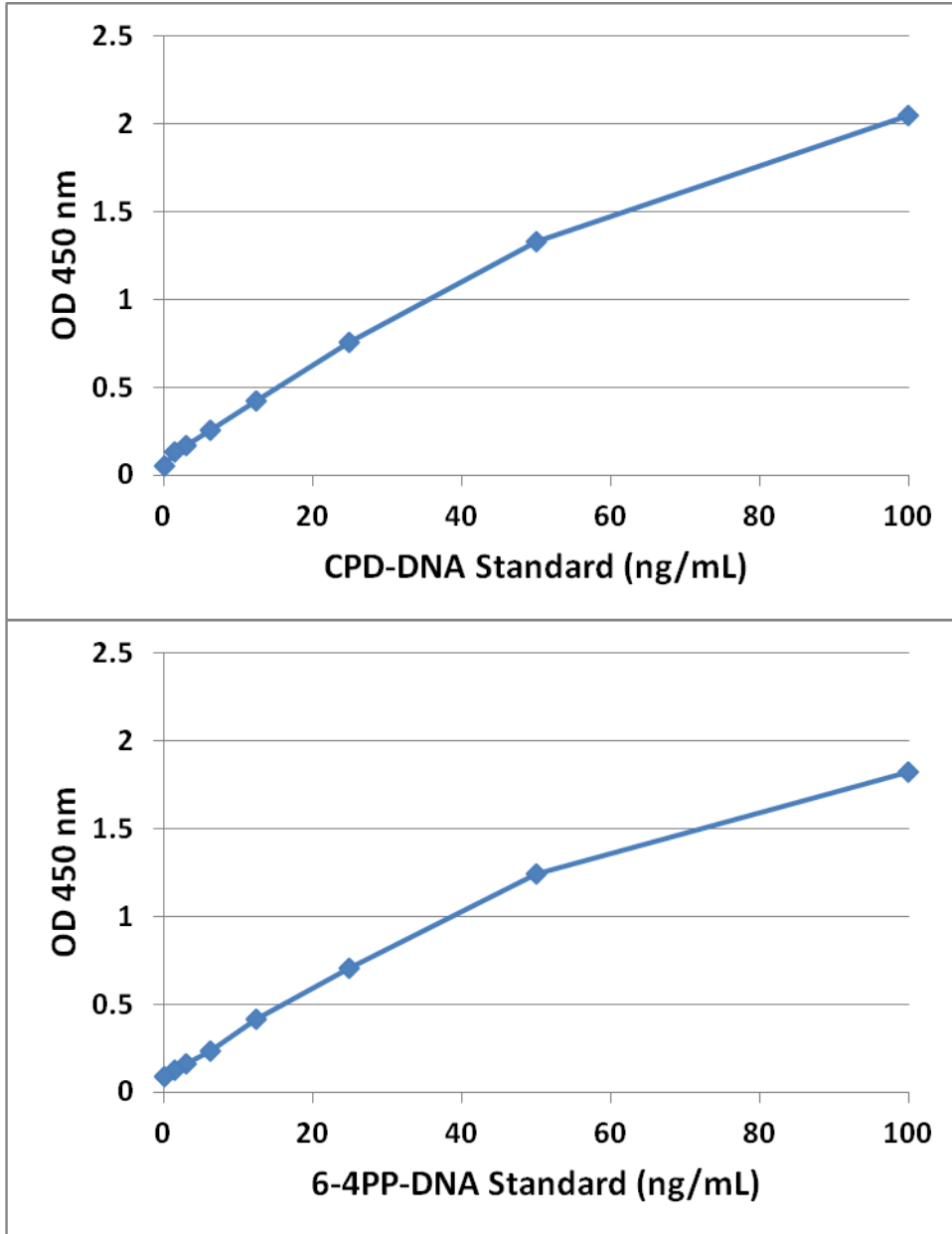


Figure 2: CPD-DNA and 6-4PP-DNA Standard Curves.

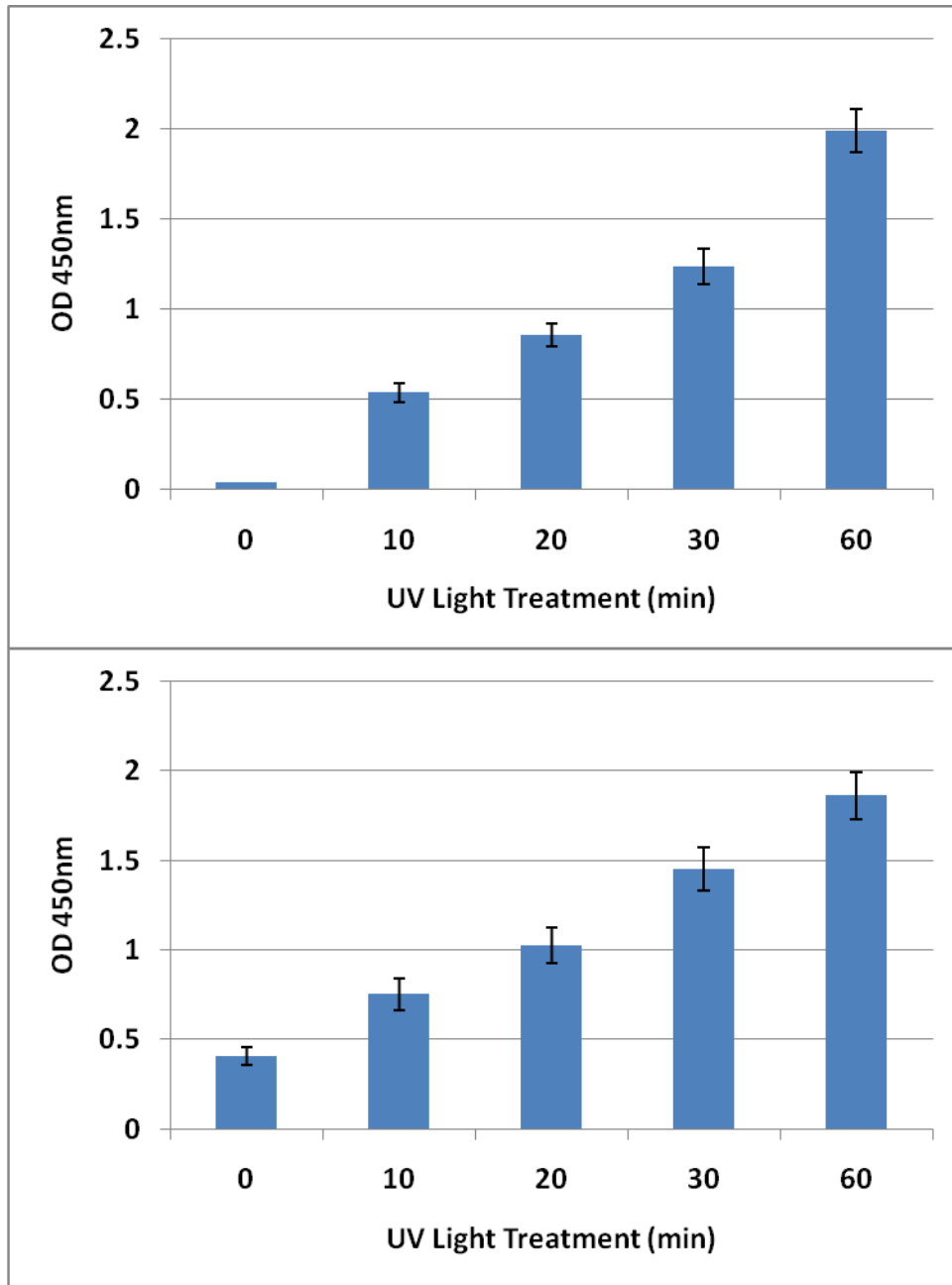


Figure 3: DNA Damage Induced by UV Light. 0.2 mg/mL Calf thymus DNA was exposed to UV light inside a cell culture hood for the time indicated. The CPD levels (**Upper panel**) and 6-4PP levels (**Lower panel**) in denatured DNA samples were determined as described in the Assay Protocol.

References

1. Lippke JA, Gordon LK, Brash DE, Haseltine WA. (1981) *Proc Natl Acad Sci U S A*. **78**:3388–3392.
2. Mitchell DL, Nairn RS. (1989) *Photochem Photobiol*. **49**:805–819.
3. Ananthaswamy HN, Loughlin SM, Cox P, Evans RL, Ullrich SE, Kripke ML. (1997) *Nat Med*. **3**:510–514.
4. Soehnge H, Ouhtit A, Ananthaswamy ON. (1997) *Front Biosci*. **2**:D538–D551.
5. el-Deiry WS, Tokino T, Velculescu VE, Levy DB, Parsons R, Trent JM, Lin D, Mercer WE, Kinzler KW, Vogelstein B. (1993) *Cell*. **75**:817–825.
6. Hermeking H, Lengauer C, Polyak K, He TC, Zhang L, Thiagalingam S, Kinzler KW, Vogelstein B. (1997) *Mol Cell*. **1**:3–11.
7. Hill LL, Ouhtit A, Loughlin SM, Kripke ML, Ananthaswamy HN, Owen-Schaub LB. (1999) *Science*. **285**:898–900.

Recent Product Citations

1. Luangpraditkun, K. et al. (2020). Photoprotective Potential of the Natural Artocarpin against In Vitro UVB-Induced Apoptosis. *Oxid Med Cell Longev*. doi: 10.1155/2020/1042451.
2. Helalat, S.H. et al. (2020). Investigating the efficacy of UVSE protein at repairing CPD and 6–4 pp DNA damages in human cells. *J Photochem Photobiol B*. doi: 10.1016/j.jphotobiol.2020.111843.
3. Wang, J. et al. (2020). Acetylation of XPF by TIP60 facilitates XPF-ERCC1 complex assembly and activation. *Nat Commun*. **11**(1):786. doi: 10.1038/s41467-020-14564-x.
4. Choong, P.F. et al. (2019). DNA repair efficiency associated with reprogrammed osteosarcoma cells. *Gene Reports*. 100409. doi: 10.1016/j.genrep.2019.100409.
5. Ben-Zvi, O. et al. (2019). Response of fluorescence morphs of the mesophotic coral *Euphyllia paradivisa* to ultra-violet radiation. *Sci Rep*. **9**(1):5245. doi: 10.1038/s41598-019-41710-3.
6. Tong, L. et al. (2018). The Mechanisms of Carnosol in Chemoprevention of Ultraviolet B-Light-Induced Non-Melanoma Skin Cancer Formation. *Sci Rep*. **8**(1):3574. doi: 10.1038/s41598-018-22029-x.
7. Fang, C. et al. (2016). Methyl-CpG binding domain protein acts to regulate the repair of cyclobutane pyrimidine dimers on rice DNA. *Sci Rep*. doi:10.1038/srep34569.
8. Gao, L. et al. (2015). The tomato DDI2, a PCNA ortholog, associating with DDB1-CUL4 complex is required for UV-damaged DNA repair and plant tolerance to UV stress. *Plant Science*. **235**:101-110.
9. Fujimori, N. et al. (2014). Plant DNA-damage repair/tolerant 100 protein repairs UV-B-induced DNA damage. *DNA Repair (Amst)*. **21**:171-176.

Warranty

These products are warranted to perform as described in their labeling and in Cell Biolabs literature when used in accordance with their instructions. THERE ARE NO WARRANTIES THAT EXTEND BEYOND THIS EXPRESSED WARRANTY AND CELL BIOLABS DISCLAIMS ANY IMPLIED WARRANTY OF MERCHANTABILITY OR WARRANTY OF FITNESS FOR PARTICULAR PURPOSE. CELL BIOLABS' sole obligation and purchaser's exclusive remedy for breach of this warranty shall be, at the option of CELL BIOLABS, to repair or replace the products. In no event shall CELL BIOLABS be liable for any proximate, incidental or consequential damages in connection with the products.

Contact Information

Cell Biolabs, Inc.
7758 Arjons Drive
San Diego, CA 92126
Worldwide: +1 858-271-6500
USA Toll-Free: 1-888-CBL-0505
E-mail: tech@cellbiolabs.com
www.cellbiolabs.com

©2011-2020: Cell Biolabs, Inc. - All rights reserved. No part of these works may be reproduced in any form without permissions in writing.