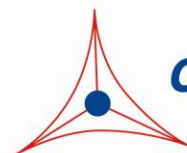

Product Manual

OxiSelect™ Comet Assay Kit (3-Well Slides)

Catalog Number

STA-351	75 assays
STA-351-5	5 x 75 assays

FOR RESEARCH USE ONLY
Not for use in diagnostic procedures



CELL BIOLABS, INC.
Creating Solutions for Life Science Research

Introduction

DNA damage, due to environmental factors and normal metabolic processes inside the cell, occurs at a rate of 1,000 to 1,000,000 molecular lesions per cell per day. While this counts for only a small part of the human genome's approximately 6 billion bases (3 billion base pairs), unrepaired lesions to critical genes can impede a cell's ability to carry out its function and appreciably increase the likelihood of cancer.

The comet assay, or single cell gel electrophoresis assay (SCGE), is a common technique for measurement of DNA damage in individual cells. Under an electrophoretic field, damaged cellular DNA (containing fragments and strand breaks) is separated from intact DNA, yielding a classic “comet tail” shape under the microscope. Extent of DNA damage is usually visually estimated by comet tail measurement; however, image analysis software is also available for measuring various parameters.

The OxiSelect™ Comet Assay is a fast and sensitive kit for the measurement of cellular DNA damage. Each kit provides sufficient reagents to perform up to 75 assays.

Assay Principle

The Cell Biolabs OxiSelect™ Comet Assay is a single cell gel electrophoresis assay (SCGE) for simple evaluation of cellular DNA damage. First, individual cells are mixed with molten agarose before application to the OxiSelect™ Comet Slide. These embedded cells are then treated with a lysis buffer and alkaline solution, which relaxes and denatures the DNA. Finally, the samples are electrophoresed in a horizontal chamber to separate intact DNA from damaged fragments. Following electrophoresis, the samples are dried, stained with a DNA dye, and visualized by epifluorescence microscopy. Under these conditions, the damaged DNA (containing cleavage and strand breaks) will migrate further than intact DNA and produce a “comet tail” shape (see Figure 1). Each kit provides sufficient reagents to perform up to 75 assays.

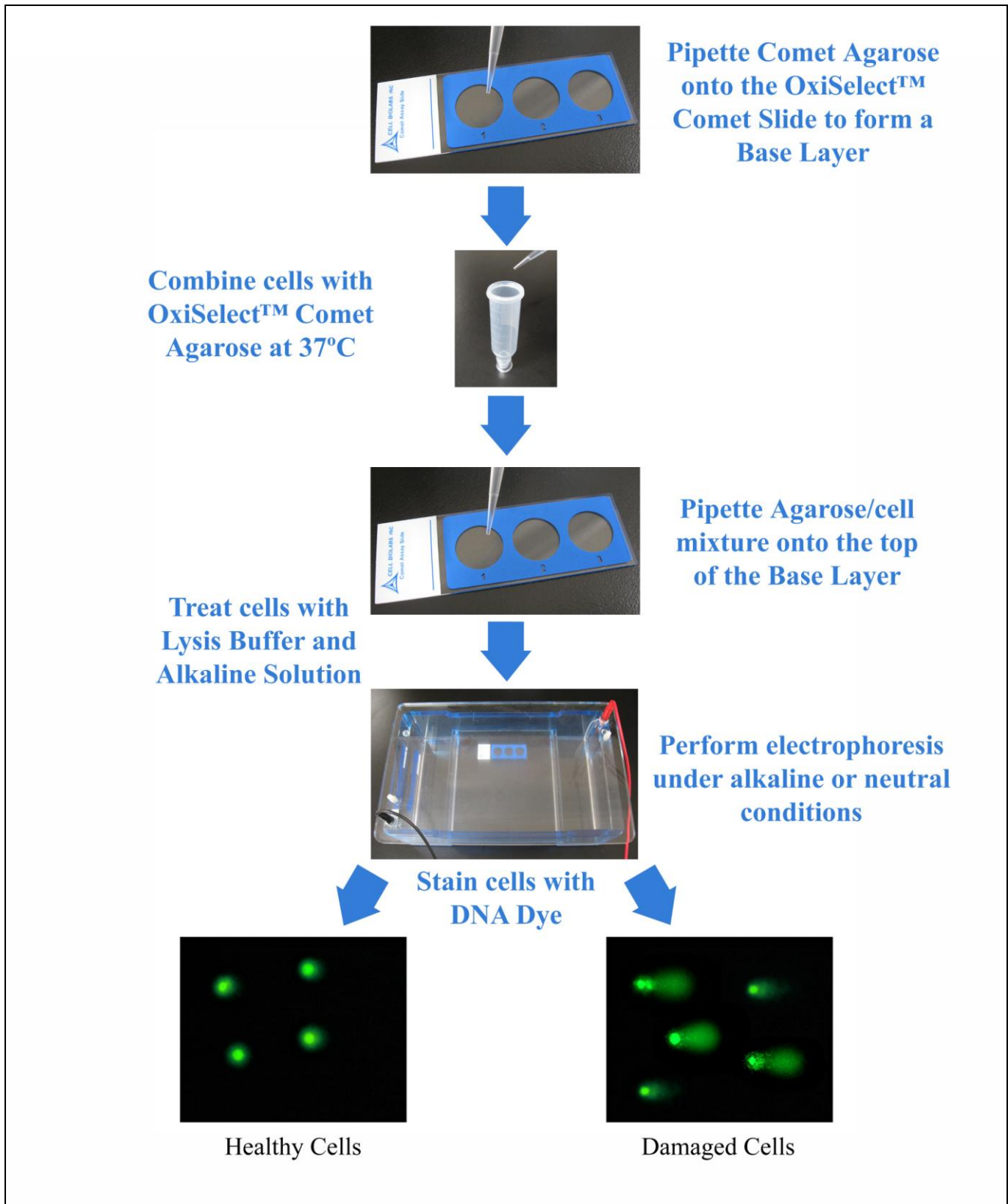


Figure 1: Comet Assay Principle

Related Products

1. STA-320: OxiSelect™ Oxidative DNA Damage ELISA Kit (8-OHdG Quantitation)
2. STA-321: OxiSelect™ DNA Double-Strand Break (DSB) Staining Kit
3. STA-324: OxiSelect™ Oxidative DNA Damage Quantitation Kit (AP sites)
4. STA-325: OxiSelect™ Oxidative RNA Damage ELISA Kit (8-OHG Quantitation)
5. STA-350: OxiSelect™ Comet Assay Kit (3-Well Slides), 15 Assays
6. STA-352: OxiSelect™ Comet Assay Slides (3-Well), 5 Slides
7. STA-353: OxiSelect™ Comet Assay Slides (3-Well), 25 Slides
8. STA-355: OxiSelect™ 96-Well Comet Assay Kit
9. STA-356: OxiSelect™ 96-Well Comet Assay Slide

Kit Components

1. OxiSelect™ 3-Well Comet Slides (Part No. STA-353): 25 slides.
2. OxiSelect™ Comet Agarose (Part No. 235002): One sterile 15 mL bottle.
3. Vista Green DNA Dye, 10000X (Part No. 235003): One 5 µL vial.
4. EDTA Solution, 500 mM (Part No. 235102): One 250 mL bottle.
5. 10X Lysis Solution (Part No. 235103): One 100 mL bottle.

Materials Not Supplied

1. NaCl powder
2. NaOH pellets
3. 10 N NaOH for pH adjustment
4. DMSO (optional)
5. 70% Ethanol
6. TE Buffer (10 mM Tris, pH 7.5, 1 mM EDTA)
7. PBS (without Mg²⁺ and Ca²⁺)
8. EDTA (disodium salt)
9. DI H₂O
10. 37°C and boiling water baths
11. Horizontal electrophoresis chamber
12. Adjustable single channel micropipettes with disposable tips
13. Epifluorescence microscope with FITC filter

Storage

Upon receipt, store the Vista Green DNA Dye at -20°C. Store all other kit components at room temperature.

Preparation of Reagents

- OxiSelect™ Comet Agarose: Heat the Comet Agarose bottle at 90-95°C in a water bath for 20 minutes, or until agarose liquefies. Transfer the bottle to a 37°C water bath for 20 minutes and maintain until needed.
- Vista Green DNA Dye: Prepare a 1X Vista Green DNA Staining Solution by diluting the provided stock 1:10000 in TE Buffer (10 mM Tris, pH 7.5, 1 mM EDTA). The solution can be stored at 4°C for up to 3 weeks, protected from light.
- Lysis Buffer: To prepare 100 mL of 1X Lysis Buffer

NaCl	14.6 g
EDTA Solution (provided)	20.0 mL
10X Lysis Solution (provided)	10.0 mL
DMSO	10.0 mL (optional for heme containing samples)
DI H ₂ O	Adjust volume to 90 mL

Mix thoroughly to dissolve NaCl. Slowly adjust the Lysis Buffer to pH 10.0 with 10 N NaOH, then QS to 100 mL with DI H₂O. Chill Lysis Buffer to 4 °C before use.

Note: Buffer will appear cloudy at room temperature, but will clear at 4°C. pH will also remain ~10.0.

- Alkaline Solution: To prepare 100 mL of Alkaline Solution

NaOH	1.2 g
EDTA Solution (provided)	0.2 mL
DI H ₂ O	Adjust volume to 100 mL

Mix thoroughly to dissolve NaOH. Chill Alkaline Solution to 4°C before use.

- Electrophoresis Running Solution: Choose the appropriate electrophoresis solution based on the desired running conditions and assay sensitivity. TBE is preferred for analysis of apoptosis and enables use of the tail length, rather than the tail moment, for data analysis. TBE electrophoresis will detect single-stranded and double-stranded DNA breaks, and may detect a few AP sites. Alkaline electrophoresis is more sensitive and will detect smaller amounts of DNA damage. Alkaline electrophoresis will detect single-stranded and double-stranded DNA breaks, the majority of AP sites, and alkali labile DNA adducts.

To prepare 1 L of Electrophoresis Solution

1. TBE Electrophoresis Solution

Tris Base	10.8 g
Boric Acid	5.5 g

EDTA (disodium salt)	0.93 g
DI H ₂ O	Adjust volume to 1 L

Mix thoroughly to dissolve solids. Chill TBE Running Solution to 4°C before use.

OR 2. Alkaline Electrophoresis Solution (300 mM NaOH, pH >13, 1 mM EDTA)

NaOH	12.0 g
EDTA Solution (provided)	2.0 mL
DI H ₂ O	Adjust volume to 1 L

Mix thoroughly to dissolve NaOH. Chill Alkaline Running Solution to 4°C before use.

Special Precautions

To avoid ultraviolet light damage to cell samples, perform the assay under low/dim light conditions.

Preparation of Samples and Slides

1. Prepare Lysis Buffer, Alkaline Solution, and Electrophoresis Running Solution (see Preparation of Reagents) prior to performing the assay. Chill all solutions to 4°C thoroughly.
2. Heat OxiSelect™ Comet Agarose to 90-95°C in a water bath for 20 minutes, or until agarose liquefies. Cool the agarose by transferring the bottle to a 37°C water bath for 20 minutes.
3. Add 75 µL of Comet Agarose per well onto the OxiSelect™ Comet Slide to create a Base Layer. **Ensure complete well coverage by spreading the solution over the well with the pipette tip.** Maintaining the slide horizontally, transfer the slide to 4°C for 15 minutes.
4. Prepare cell samples, including controls, as follows:
 - Suspension Cells: Centrifuge cells at 700 x g for 2 minutes and discard supernatant. Wash cell pellet once with ice-cold PBS (without Mg²⁺ and Ca²⁺), centrifuge, and discard the supernatant. Finally, resuspend the cells at 1 x 10⁵ cells/mL in ice-cold PBS (without Mg²⁺ and Ca²⁺).
 - Adherent Cells: Gently remove cells from flask/dish by scraping with a rubber policeman. Transfer cell suspension to a conical tube and centrifuge at 700 x g for 2 minutes, discarding the supernatant. Wash cell pellet once with ice-cold PBS (without Mg²⁺ and Ca²⁺), centrifuge, and discard the supernatant. Finally, resuspend the cells at 1 x 10⁵ cells/mL in ice-cold PBS (without Mg²⁺ and Ca²⁺).
 - Tissue Preparation: Using dissection scissors, mince a small piece of tissue in 1-2 mL of ice cold PBS containing 20 mM EDTA (without Mg²⁺ and Ca²⁺). Allow the tissue/cell suspension to stand for 5 minutes before transferring the supernatant to a centrifuge tube; avoid transferring debris. Centrifuge, discarding the supernatant, and then resuspend the cells at 1 x 10⁵ cells/mL in ice-cold PBS (without Mg²⁺ and Ca²⁺).
5. Combine cell samples with Comet Agarose (step 2) at 1:10 ratio (v/v), mix well by pipetting, and immediately transfer 75 µL/well onto the top of the Comet Agarose Base Layer (step 3). **Ensure complete well coverage by spreading the suspension very gently and carefully with the pipette tip, without disturbing the Base Layer.**

Note: For multiple samples, maintain suspensions at 37°C to avoid gelation. Titrates samples again just prior to slide addition.

6. Maintaining the slide horizontally, transfer the slide to 4°C in the dark for 15 minutes.
7. Carefully, transfer the slide to a small basin/container containing pre-chilled Lysis Buffer (~25 mL/slide). Immerse the slide in the buffer for 30-60 minutes at 4°C in the dark.
8. Carefully, aspirate the Lysis Buffer from the container and replace with pre-chilled Alkaline Solution (~25 mL/slide). Immerse the slide in the solution for 30 minutes at 4°C in the dark.

Assay Protocol

I. TBE Electrophoresis

1. Aspirate the Alkaline Solution from the container and replace with pre-chilled TBE Electrophoresis Solution. Immerse the slide for 5 minutes, and then repeat once more.
2. Maintaining the slide horizontally, carefully transfer the slide to a horizontal electrophoresis chamber. Fill the chamber with cold TBE Electrophoresis Solution until the buffer level covers the slide.
3. Apply voltage to the chamber for 10-15 minutes at 1 volt/cm (e.g. if the chamber electrodes are 35 cm apart, you would then apply 35 volts to the slide)
4. Maintaining the slide horizontally, carefully transfer the slide from the electrophoresis chamber to a clean, small basin/container containing pre-chilled DI H₂O (~25 mL/slide). Immerse the slide for 2 minutes, aspirate, and then repeat twice more.
5. Aspirate the final water rinse and replace with cold 70% Ethanol for 5 minutes.
6. Maintaining the slide horizontally, remove the slide from the 70% Ethanol and allow to air dry.
7. Once the agarose and slide is completely dry, add 100 µL/well of diluted Vista Green DNA Dye (see Preparation of Reagents). Incubate at room temperature for 15 minutes.
8. View slides by epifluorescence microscopy using a FITC filter.

II. Alkaline Electrophoresis

1. Maintaining the slide horizontally, carefully transfer the slide from the Alkaline Solution to a horizontal electrophoresis chamber. Fill the chamber with cold Alkaline Electrophoresis Solution until the buffer level covers the slide.
2. Apply voltage to the chamber for 15-30 minutes at 1 volt/cm (e.g. if the chamber electrodes are 35 cm apart, you would then apply 35 volts to the slide). Additionally, adjust the volume of Alkaline Electrophoresis Solution to produce a current setting of 300 mA.
3. Maintaining the slide horizontally, carefully transfer the slide from the electrophoresis chamber to a clean, small basin/container containing pre-chilled DI H₂O (~25 mL/slide). Immerse the slide for 2 minutes, aspirate, and then repeat twice more.
4. Aspirate the final water rinse and replace with cold 70% Ethanol for 5 minutes.
5. Maintaining the slide horizontally, remove the slide from the 70% Ethanol and allow to air dry.

6. Once the agarose and slide is completely dry, add 100 μL /well of diluted Vista Green DNA Dye (see Preparation of Reagents). Incubate at room temperature for 15 minutes.
7. View slides by epifluorescence microscopy using a FITC filter.

Example of Results

The following figures demonstrate typical OxiSelect™ Comet Assay Kit results. One should use the data below for reference only. This data should not be used to interpret actual results.

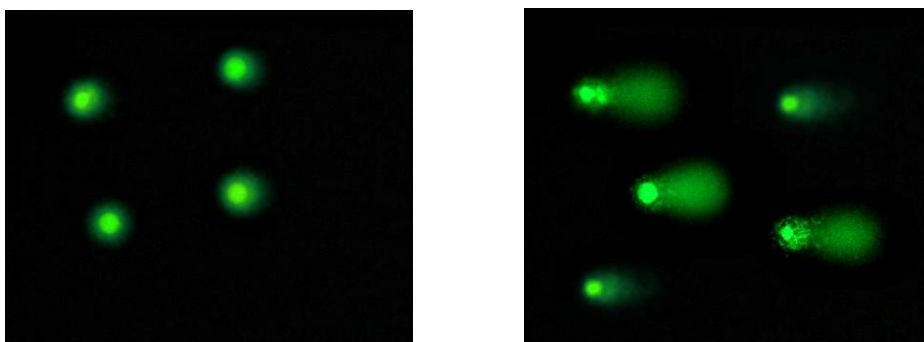


Figure 2. Etoposide Treatment of Jurkat Cells. Jurkat cells were untreated (left) or treated (right) with 20 μM Etoposide for 4 hours before performing Comet Assay (alkaline electrophoresis conditions, 33 V/300 mA for 15 minutes).

Calculation of Results

The DNA damage is quantified by measuring the displacement between the genetic material of the nucleus ('comet head') and the resulting 'tail'. Tail Moment and Tail DNA% are the two most common parameters to analyze Comet assay results. At least 50 -100 cells should be analyzed per sample. The Tail Moment has been suggested to be an appropriate index of induced DNA damage in considering both the migration of the genetic material as well as the relative amount of DNA in the tail.

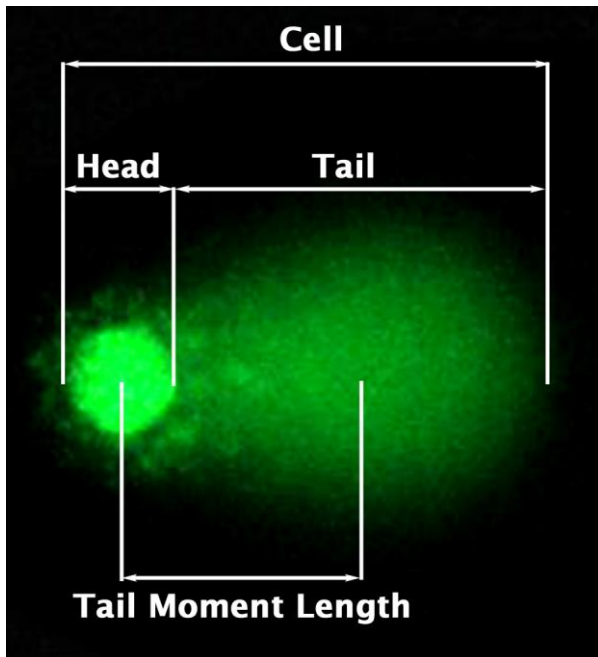


Figure 3: Typical Damaged DNA in Comet Assay.

$$\text{Tail DNA\%} = 100 \times \frac{\text{Tail DNA Intensity}}{\text{Cell DNA Intensity}}$$

Tail Moment can be measured using one of the following methods:

(a) Olive Tail Moment = Tail DNA% x Tail Moment Length*

(b) Extent Tail Moment = Tail DNA% x Length of Tail (see Figure 3)

A number of Comet Assay analysis software programs are commercially available, such as and Comet Assay IV (Perceptive Instruments) and CASPlab.

*Tail Moment Length is measured from the center of the head to the center of the tail (see Figure 3)

References

1. Ostling, O., and Johanson, K. J. (1984). *Biochem. Biophys. Res. Commun.* **123**, 291–298.
2. Singh, N. P., McCoy, M. T., Tice, R. R., and Schneider, E. L. (1988). *Exp. Cell. Res.* **175**, 184–191.
3. Olive, P. L., Banath, J. P., and Durand, R. E. (1990). *Radiat. Res.* **122**, 86–94.
4. De Boeck, M., Touil, N., De Visscher, G., Vande, P. A., and Kirsch-Volders, M. (2000). *Mutat. Res.* **469**, 181–197

Recent Product Citations

1. Siemionow, M. et al. (2020). Transplantation of Dystrophin Expressing Chimeric (DEC) Human Cells of Myoblast/MSC Origin Improves Function in Duchenne Muscular Dystrophy Model. *Stem Cells Dev.* doi: 10.1089/scd.2020.0161.
2. Lammert, C.R. et al. (2020). AIM2 inflammasome surveillance of DNA damage shapes neurodevelopment. *Nature.* doi: 10.1038/s41586-020-2174-3.
3. Shibayama, Y. et al. (2020). Aberrant (pro)renin receptor expression induces genomic instability in pancreatic ductal adenocarcinoma through upregulation of SMARCA5/SNF2H. *Commun Biol.* **3**(1):724. doi: 10.1038/s42003-020-01434-x.
4. Han, J. et al. (2020). Elevated CXorf67 Expression in PFA Ependymomas Suppresses DNA Repair and Sensitizes to PARP Inhibitors. *Cancer Cell.* doi: 10.1016/j.ccell.2020.10.009.
5. Hays, E. et al. (2020). The SWI/SNF ATPase BRG1 stimulates DNA end resection and homologous recombination by reducing nucleosome density at DNA double strand breaks and by promoting the recruitment of the CtIP nuclease. *Cell Cycle.* doi: 10.1080/15384101.2020.1831256.
6. Ibnu Rasid, E.N. et al. (2020). Effect of Dioscorea hispida var. Daemona (Roxb) Prain & Burkill on Oxidative Stress and DNA Damage in the Liver of Pregnant Rats. *Int J Biomed Sci.* **16**(3).

7. Le, B.V. et al. (2020). TGF β R-SMAD3 Signaling Induces Resistance to PARP Inhibitors in the Bone Marrow Microenvironment. *Cell Rep.* **33**(1):108221. doi: 10.1016/j.celrep.2020.108221.
8. Klotz-Noack, K. et al. (2020). SFPQ Depletion Is Synthetically Lethal with BRAFV600E in Colorectal Cancer Cells. *Cell Rep.* **32**(12):108184. doi: 10.1016/j.celrep.2020.108184.
9. Ito, S.S. et al. (2020). Inhibition of the ATR kinase enhances 5-FU sensitivity independently of non-homologous end-joining and homologous recombination repair pathways. *J Biol Chem.* doi: 10.1074/jbc.RA120.013726.
10. Klak, M. et al. (2020). Irradiation with 365 nm and 405 nm wavelength shows differences in DNA damage of swine pancreatic islets. *PLoS One.* **15**(6):e0235052. doi: 10.1371/journal.pone.0235052.
11. Khalil, A.M. et al. (2020). Association between Mobile Phone Using and DNA Damage of Epithelial Cells of the Oral Mucosa. *J Biotechnol Biomed.* **3**(2020): 50-66. doi: 10.26502/jbb.2642-91280027.
12. Wang, Y. et al. (2020). Targeting therapeutic vulnerabilities with PARP inhibition and radiation in IDH-mutant gliomas and cholangiocarcinomas. *Sci Adv.* doi: 10.1126/sciadv.aaz3221.
13. Fang, Y. et al. (2020). Epigenetic dysregulation of Mdr1b in the blood-testis barrier contributes to dyszoospermia in mice exposed to cadmium. *Ecotoxicol Environ Saf.* **190**:110142. doi: 10.1016/j.ecoenv.2019.110142.
14. Cupello, S. et al. (2019). Distinct roles of XRCC1 in genome integrity in *Xenopus* egg extracts. *Biochem J.* **476**(24):3791-3804. doi: 10.1042/BCJ20190798.
15. Naci, D. et al. (2019). Cell adhesion to collagen promotes leukemia resistance to doxorubicin by reducing DNA damage through the inhibition of Rac1 activation. *Sci Rep.* **9**(1):19455. doi: 10.1038/s41598-019-55934-w.
16. Lu, S. et al. (2019). Additive effects of a small molecular PCNA inhibitor PCNA-IIS and DNA damaging agents on growth inhibition and DNA damage in prostate and lung cancer cells. *PLoS One.* **14**(10):e0223894. doi: 10.1371/journal.pone.0223894.
17. Dedobbeleer, M. et al. (2019). MKP1 phosphatase is recruited by CXCL12 in Glioblastoma cells and plays a role in DNA Strand Breaks Repair. *Carcinogenesis.* doi: 10.1093/carcin/bgz151.
18. Zhou, J. et al (2019). Aberrantly Expressed Timeless Regulates Cell Proliferation and Cisplatin Efficacy in Cervical Cancer. *Hum Gene Ther.* doi: 10.1089/hum.2019.080.
19. Sarkar, R. et al. (2019). Rotavirus activates a noncanonical ATM-Chk2 branch of DNA damage response during infection to positively regulate viroplasm dynamics. *Cell Microbiol.* doi: 10.1111/cmi.13149.
20. Maner, J. et al. (2019). Hexachlorobenzene exerts genotoxic effects in a humpback whale cell line under stable exposure conditions. *RSC Adv.* **9**:39447-39457. doi: 10.1039/C9RA05352B.
21. Han, Y. et al. (2019). Exposure to waterborne nTiO₂ reduces fertilization success and increases polyspermy in a bivalve mollusc: A threat to population recruitment. *Environ Sci Technol.* doi: 10.1021/acs.est.9b03675.
22. Shansky, Y.D. et al. (2019). Human Platelet Lysate Sustains the Osteogenic/Adipogenic Differentiation Potential of Adipose-Derived Mesenchymal Stromal Cells and Maintains Their DNA Integrity in vitro. *Cells Tissues Organs.* doi: 10.1159/000502813.
23. Gothai, S. et al. (2019). In Vitro and In Vivo-Scientific Evaluation on Cytotoxicity and Genotoxicity of Traditional Medicinal Plant *Couroupita Guianensis* Aubl. Flower. *PhOL.* **2**:24-38.
24. Al Khateeb, W. et al. (2019). Growth, Yield and Genetic Integrity of Spinach and Chrysanthemum as Affected by Soil Supplementation with Dam Sediments Collected From King Talal and Al-Mujib Dams/ Jordan. *World Appl. Sci. J.* **37**(1):58-69. doi: 10.5829/idosi.wasj.2019.58.69.

25. Su, J. et al. (2019). Genomic Integrity Safeguards Self-Renewal in Embryonic Stem Cells. *Cell Rep.* **28**(6):1400-1409.e4. doi: 10.1016/j.celrep.2019.07.011.
26. Xia, Y. et al. (2019). Rescue of DNA damage after constricted migration reveals a mechano-regulated threshold for cell cycle. *J Cell Biol.* pii: jcb.201811100. doi: 10.1083/jcb.201811100.
27. Chesnokova, V. et al. (2019). Growth hormone induces colon DNA damage independent of IGF-1. *Endocrinology.* pii: en.2019-00132. doi: 10.1210/en.2019-00132.
28. Garzón, J. et al. (2019). Human RIF1-Protein Phosphatase 1 Prevents Degradation and Breakage of Nascent DNA on Replication Stalling. *Cell Rep.* **27**(9):2558-2566.e4. doi: 10.1016/j.celrep.2019.05.002.
29. Nakashima, A. et al. (2019). Autophagy is a new protective mechanism against the cytotoxicity of platinum nanoparticles in human trophoblasts. *Sci Rep.* **9**(1):5478. doi: 10.1038/s41598-019-41927-2.
30. Pillay, N. et al. (2019). DNA Replication Vulnerabilities Render Ovarian Cancer Cells Sensitive to Poly(ADP-Ribose) Glycohydrolase Inhibitors. *Cancer Cell.* **35**(3):519-533.e8. doi: 10.1016/j.ccell.2019.02.004.

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's 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

©2008-2021: Cell Biolabs, Inc. - All rights reserved. No part of these works may be reproduced in any form without permissions in writing.