



Trevigen Press Release

NEW PRODUCTS

Anti-8-oxo-dG (Clone 2E2)

DNA damage caused by oxidative radicals is associated with aging, inflammation, carcinogenesis, Parkinson's and Alzheimer's diseases. When exposed to oxidative radicals, 8-hydroxy-2'-deoxyguanosine (8-oxo-dG) is formed and can serve as a sensitive indicator of physiological and environmental damage to DNA. This mouse monoclonal antibody specifically binds to 8-hydroxy-2'- deoxyguanosine within DNA in H₂O₂-treated cells. It can be used to detect oxidative damage by ELISA and immunocytochemistry. Sufficient antibody is provided for approximately 50 slides, when a 1:250 dilution is used. Contact Trevigen, Inc. for details at Tel: 800-873-8443, Email: info@trevigen.com, or see our website at www.trevigen.com.

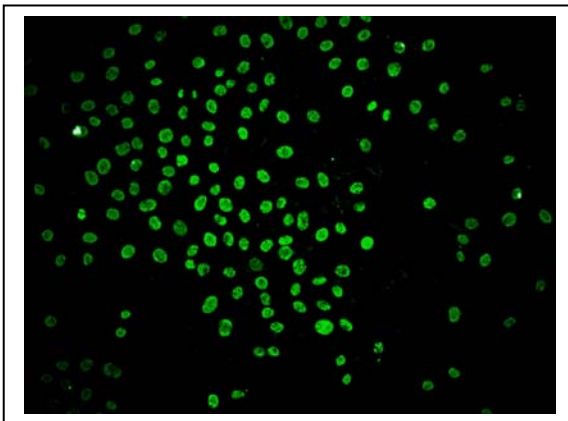


Figure A

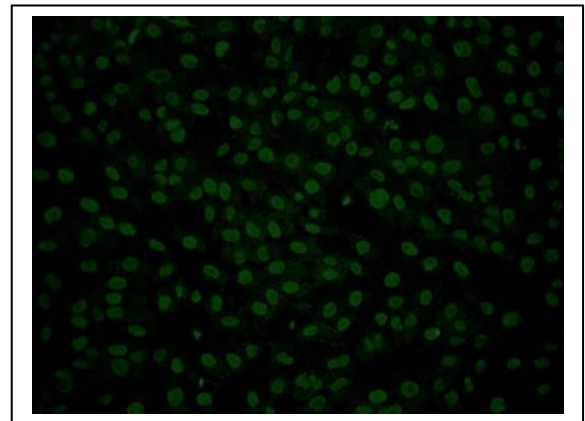


Figure B

H₂O₂ treated (A) and untreated (B) MCF-10A cells stained with 8-oxo-dG antibody (Cat# 4354-MC-050) according to the above protocol using AlexaFluor 488 conjugated anti-mouse antibody.

Trevigen, Inc.

8405 Helgerman Court, Gaithersburg, MD 20877 USA

Voice: 1-800-TREVIGEN (1-800-873-8443) • 301-216-2800 • Fax: 301-560-4973

Email: info@trevigen.com • <http://www.trevigen.com>