## **Comet SCGE assay kit**

Fast and simple electrophoresis method to measure DNA fragmentation

Exposure of cells to oxidative and environmental stresses frequently results in the breakdown or oxidation of genomic DNA. Assays to evaluate the integrity of genomic DNA, or to assess the presence of oxidized DNA are frequently used as a means of verifying the onset of apoptosis or DNA damage. The Comet SCGE Assay measures DNA damage by fluorescently detecting the integrity of DNA liberated from cells embedded in low melting point agarose. Upon electrophoresis, fragmented DNA produces a characteristic "comet" shaped tail as small DNA fragments migrate in the gel more rapidly than in-tact genomic DNA. The Comet SCGE Assay is a fast and simple electrophoresis method to detect and quantitate DNA fragmentation in cells associated with DNA damage and apoptosis. A unique nucleic acid stain, CYGREEN<sup>®</sup> Nucleic acid dye, provides improved sensitivity for DNA visualization compared to ethidium bromide.

Citations: 46

View Online »

**Ordering Information** 

Order Online »

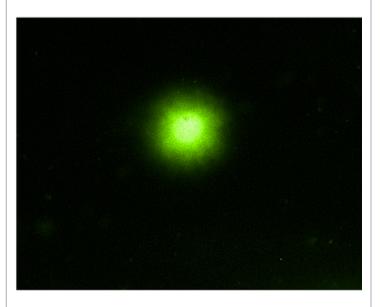
ADI-900-166

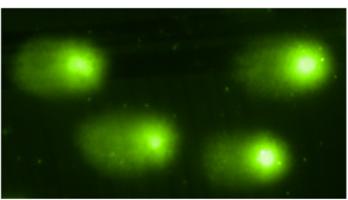
50 tests

Manuals, SDS & CofA

View Online »

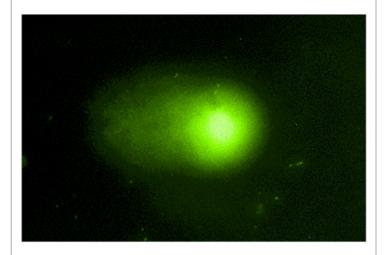
- Ready-to-use Comet Slides allow direct application of sample without pretreatment
- Shorter assay time allows for higher throughput sample analysis
- Hydrophobic barrier allows sample treatment with DNA repair enzymes
- Unique nucleic acid stain provides improved sensitivity for DNA visualization compared to ethidium bromide





Negative control. CaSki cells were untreated. Comet SCGE Assay was performed and slides were imaged using a FITC filter.

Positive control. CaSki cells were treated with 100mM hydrogen peroxide for 20 min at 4°C. Comet SCGE Assay was performed and slides were imaged using a FITC filter.



Positive control. CaSki cells were treated with 100mM hydrogen peroxide for 20 min at 4°C. Comet SCGE Assay was performed and slides were imaged using a FITC filter.

## **Handling & Storage**

**Use/Stability** Store CYGREEN® Nucleic Acid Dye at -20°C and all other reagents at +4°C.

Long Term Storage +4°C

**Shipping** Blue Ice

## Regulatory Status RUO - Research Use Only

## **Product Details**

**Application** Fluorescence microscopy

**Application Notes** For the measurement of DNA Damage, DNA Repair, and

Apoptosis in cells and tissue from any species via

electrophoresis.

Assay Time 3 hours

Contents Lysis solution, Comet LMAgarose, Comet slide, EDTA,

CYGREEN® Nucleic acid dye

Species Reactivity Species independent

**Technical Info / Product Notes**Please note that CYGREEN<sup>®</sup> Nucleic acid dye is now

included with the Comet SCGE assay kit. Please follow the

recommended staining procedures in the updated manual.

Last modified: February 19, 2025

European Sales Office