

## Product datasheet

# Comet Assay Kit (3-well slides) ab238544

★★★★★ [2 Abreviews](#) [29 References](#) [3 Images](#)

### Overview

**Product name** Comet Assay Kit (3-well slides)  
**Detection method** Fluorescent  
**Product overview** Comet Assay Kit (3-well slides) (ab238544) is a fast and sensitive kit for the measurement of cellular DNA damage.

The Comet Assay is a single cell gel electrophoresis assay (SCGE) for simple evaluation of cellular DNA damage. It is a convenient way to screen for general DNA damage, regardless of the source or nature of the damage. Kits include Comet Slides, reagents, and a fluorescent dye to visualize cells under an epifluorescence microscope.

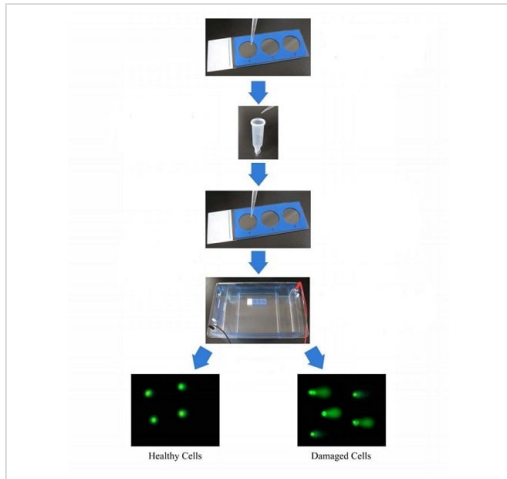
**Notes** Abcam has not and does not intend to apply for the REACH Authorisation of customers' uses of products that contain European Authorisation list (Annex XIV) substances. It is the responsibility of our customers to check the necessity of application of REACH Authorisation, and any other relevant authorisations, for their intended uses.

### Properties

**Storage instructions** Please refer to protocols.

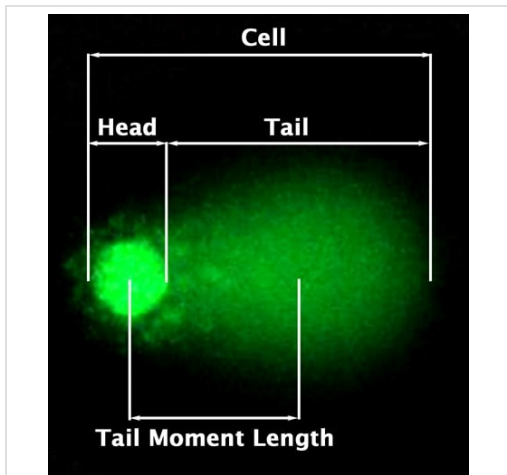
Components	15 tests	75 tests
10X Lysis Solution	1 x 20ml	1 x 100ml
3-Well Comet Slides	1 x 5 slides	1 x 25 slides
Comet Agarose	1 x 15ml	1 x 15ml
EDTA Solution	1 x 50ml	1 x 250ml
Vista Green DNA Dye, 10000X	1 x 5µl	1 x 5µl

### Images



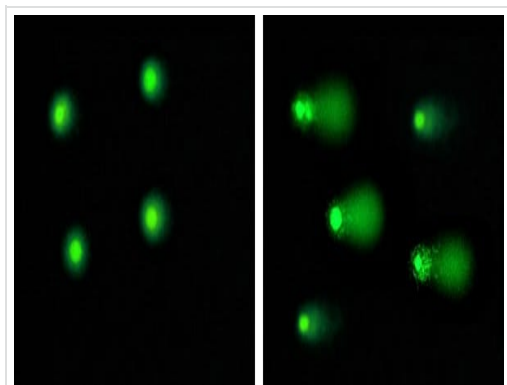
Comet Assay Principle

Comet Assay Principle



For Data Analysis please see Section 9 in the protocol booklet.

Epifluorescence Microscopy Visualization of DNA Damage.



Jurkat cells were untreated (left) or treated (right) with 20  $\mu$ M Etoposide for 4 hrs before performing Comet Assay (alkaline electrophoresis conditions, 33 V/300 mA for 15 mins).

Etoposide Treatment of Jurkat Cells.

**Please note:** All products are "FOR RESEARCH USE ONLY. NOT FOR USE IN DIAGNOSTIC PROCEDURES"

## **Our Abpromise to you: Quality guaranteed and expert technical support**

---

- Replacement or refund for products not performing as stated on the datasheet
- Valid for 12 months from date of delivery
- Response to your inquiry within 24 hours
  
- We provide support in Chinese, English, French, German, Japanese and Spanish
- Extensive multi-media technical resources to help you
- We investigate all quality concerns to ensure our products perform to the highest standards

If the product does not perform as described on this datasheet, we will offer a refund or replacement. For full details of the Abpromise, please visit <https://www.abcam.com/abpromise> or contact our technical team.

## **Terms and conditions**

---

- Guarantee only valid for products bought direct from Abcam or one of our authorized distributors