

Product datasheet

Comet Assay Slide (3-well) ab238546

1 Image

Overview

Product name	Comet Assay Slide (3-well)
Product overview	Comet Assay Slide (3-well) (ab238546) are a useful screening tool for various types of DNA damage. The slides are specially treated for adhesion of low-melting agarose used in the comet assay. These slides may be used in conjunction with reagents found in Comet Assay Kit (3-well slides) <a href="#">ab238544</a> or with your own comet assay reagents.

Properties

Storage instructions	Store at room temperature. Please refer to protocols.
----------------------	---

Components	5 slides	25 slides
3-Well Comet Slide	5 slides	25 slides

Images



3-well Comet Slide

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