

# TREVIGEN® Product Data

*For Research Use Only. Not For Use In Diagnostic Procedures*

---

## hOGG1 FLARE™ Module

**Catalog #:** 4130-100-FM

**Contents:**

<u>Component</u>	<u>Catalog #</u>	<u>Amount</u>	<u>Storage</u>
hOGG1	4100-100-01	100 units	-20°C
25X FLARE™ Buffer 1	3950-040-01	40 ml	Room Temp.
REC™ Dilution Buffer	3950-010-03	10 ml	-20°C
100X BSA Additive	3950-100-04	100 µl	-20°C

**Description:** Reagents for analysis of DNA damage in single cells using the Trevigen CometAssay®, FLARE™ Slides, and hOGG1.

**Assay & Analysis:** Refer to hOGG1 FLARE™ Assay Kit's (cat#: 4130-100-FK) Instructions for Use for further details.

Note: CometSlides and FLARE Slides are interchangeable in the FLARE Assay. The difference is in the sample number; the CometSlide™ holds 2 samples whereas the FLARE Slide holds 3.

**Lot Specific Data:** Contact Trevigen for lot specific activity of hOGG1.

## TREVIGEN®

8405 Helgerman Court, Gaithersburg, MD 20877 USA  
 Voice: 1-800-TREVIGEN (1-800-873-8443) • 301-216-2800  
 Fax: 301-560-4973 • e-mail: [info@trevigen.com](mailto:info@trevigen.com) • [www.trevigen.com](http://www.trevigen.com)

**Related Products:**

**FLARE™ Assay Kits:**

Catalog #	Description	Damage Recognized	Size
4040-100-FK	<b>Fpg Kit</b>	8-oxoguanine, DNA containing	<b>75 samples</b>
4040-100-FM		formamidopyrimidine moieties	<b>100 samples</b>
4055-100-FK	<b>T4 Endo V Kit</b>	Cis-syn isomers of cyclo-	<b>75 samples</b>
4055-100-FM		butane pyrimidine dimers	<b>100 samples</b>
4065-100-FK	<b>cv-PDG Kit</b>	Cis-syn and trans-syn isomers of	<b>75 samples</b>
4065-100-FM		cyclobutane pyrimidine dimers	<b>100 samples</b>
4130-100-FK	<b>hOGG1 Kit</b>	8-oxoguanine, DNA containing	<b>75 samples</b>
		formamidopyrimidine moieties	<b>100 samples</b>
4100-100-FK	<b>UVDE Kit</b>	Cyclobutane pyrimidine	<b>75 samples</b>
4100-100-FM		dimers, (6-4) photoproducts	<b>100 samples</b>
4045-010-FK	<b>Endonuclease III Kit</b>	Thymine Glycol, 5,6-dihydro-	<b>75 samples</b>
4045-010-FM		thymine, urea, 5-hydroxy-6-hydrothymine, 5,6-dihydro-uracil, alloxan, 5-hydroxy-6-hydrouracil, uracil glycol, 5-hydroxy-5-methylhydantoin, 5-hydroxycytosine, 5-hydroxy uracil, methyltaronyl urea, thymine ring saturated or fragmentation product	<b>100 samples</b>

**CometAssay® Kits:**

Catalog #	Description	Size
4250-050-ESK	CometAssay® Starter Kit	each
4250-050-ES	CometAssay® Electrophoresis System	1 system
4250-050-K	CometAssay®	50 samples
4252-040-K	CometAssay® Higher Throughput Kit	40 samples
4254-200-K	CometAssay® Silver Staining Kit	200 samples
4253-096-K	CometAssay® Kit 96 Wells	96 samples

**hOGG1 FLARE™  
Module**

Catalog #: 4130-100-FM  
Storage: -20 °C &  
Room Temp.

**TREVIGEN®**  
1-800-873-8443