

TREVIGEN® Product Data

For Research Use Only. Not For Use In Diagnostic Procedures

Fpg FLARE™ Module

Catalog #: 4040-100-FM

Contents:

<u>Component</u>	<u>Catalog #</u>	<u>Amount</u>	<u>Storage</u>
<i>E. coli</i> Fpg*	4040-010-01	500 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

*For optimal stability, freeze in working aliquots at -80°C to avoid repeated freeze-thawings.

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

Assay & Analysis: Refer to Fpg FLARE™ Assay Kit's (cat# 4040-100-FK) Instructions for Use for further details.

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

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

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 • www.trevigen.com

FLARE™ Assay Kits:

Catalog #	Description	Substrate	Size
4040-100-FK	Fpg Kit	8-oxoguanine	75 samples
4055-100-FK	T4-PDG Kit	Cis-syn isomers of cyclobutane pyrimidine dimers	75 samples
4065-100-FK	cv-PDG Kit	Cis-syn and trans-syn isomers of cyclobutane pyrimidine dimers	75 samples
4130-100-FK	hOGG1 Kit	8-oxoguanine, DNA containing formamidopyrimidine moieties	75 samples
4100-100-FK	UVDE Kit	Cyclobutane pyrimidine dimers, (6-4) photoproducts	75 samples
4045-01K-FK	Endonuclease III Kit	Thymine Glycol, 5,6-dihydrothymine 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, methyltartonyl urea, thymine ring saturated or fragmentation product	75 samples

Fpg FLARE™ Module
Catalog #: 4040-100-FM
Storage: -20 °C
 (Manual Defrost Freezer)
TREVIGEN®
 1-800-873-8443