

TREVIGEN® Product Data

For Research Use Only. Not For Use In Diagnostic Procedures

E. coli MutY DNA Glycosylase

Catalog #: 4000-500-EB

Contents: 4000-500-01 *E. coli* MutY DNA Glycosylase **Size:** 5000 Units
 3900-500-04 10X REC™ Buffer 4 1 ml

Description: *E. coli* MutY acts together with Fpg to prevent the potentially mutagenic consequences of 8-oxo-dG lesions. 8-oxo-dG lesions escaping repair by Fpg frequently pair with A during DNA replication, producing an 8-oxo-dG:A mispair. MutY removes the A from this base pair to initiate base excision repair. In the absence of MutY, DNA replication after an 8-oxo-dG:A mismatch results in thymine incorporation opposite the adenine in one of the daughter strands, creating a fixed mutation. MutY has an associated AP lyase activity.

Source: Purified from *E. coli* containing a recombinant plasmid harboring the *E. coli* MutY gene.

Unit Definition: One unit of enzyme cleaves 1 pmole of an oligonucleotide duplex containing an A/G mismatch in 1 hour at 37 °C. Only the strand with the A is cleaved.

Assay Conditions & Analysis: 4 pmoles of A/G mismatch oligonucleotide set with the A oligo end-labeled, 1X REC™ Buffer 4 (10 mM HEPES-KOH (pH 7.4), 100 mM KCl, 10 mM EDTA), and serial dilutions of enzyme in a 20 µl reaction volume are incubated for 1 hour at 37 °C. To complete cleavage of abasic site, fresh NaOH is added to final concentration of 166 mM then heated for 15 minutes at 95 °C. For analysis, 24 µl of 2X Loading Buffer (20 mM EDTA, 97% formamide, and 0.2% bromophenol blue) is added, the samples are heated at 95 °C for 5 min then fast cooled to 4 °C, and the cleavage products are resolved by 20% denaturing polyacrylamide gel electrophoresis. The bands are analyzed to quantify the cleavage products.

Storage Buffer: 20 mM Tris pH 7.5, 100 mM NaCl, 1 mM EDTA, 1 mM DTT, and 50% (v/v) glycerol.

Storage Conditions: Store at -20 °C in a manual defrost freezer. For long term storage, freeze in working aliquots at -80 °C. Avoid repeated freeze-thawings. Enzyme may be diluted in storage buffer containing 0.1 mg/ml BSA and stored at -20 °C for 2 weeks of experimental use.

References:

1. Lu, A. and I. Hsu. 1992. Detection of single DNA base mutations with mismatch repair enzymes. *Genomics* **14**:249-255.
2. Friedberg, E.C., G.C. Walker, and W. Siede. 1995. DNA Repair and Mutagenesis. American Society of Microbiology. Washington, D.C: ASM Press.
3. Hsu, I., W.E. Highsmith, J. Xu, and D. Kong. 1998. Mismatch cleavage detects base deletion in cystic fibrosis gene. *Biotechniques* **25**:692-696.

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

Related Products:

Catalog#	Description	Size
4020-100-EB	Human DNA Polymerase β	100 U
4025-100-EB	<i>E. coli</i> Uracil-N-Glycosylase (UNGase)	100 U
4040-100-EB	<i>E. coli</i> Formamidopyrimidine-DNA Glycosylase (Fpg)	500 U
4045-01K-EB	<i>E. coli</i> Endonuclease III (Thymine Glycol-DNA Glycosylase)	1000 U
4050-100-EB	<i>E. coli</i> Endonuclease IV (nfo protein)	100 U
4055-100-EB	T4 Endonuclease V (T4-Pyrimidine Dimer Glycosylase/T4-PDG)	10 ⁵ U
4060-01K-EB	<i>E. coli</i> Endonuclease VIII	1000 U
4065-100-EB	Chlorella Virus Pyrimidine Dimer Glycosylase (cv-PDG)	1000 U
4070-500-EB	Thermostable TDG Protein (Thymine DNA Glycosylase)	500 U
4100-100-EB	<i>S. pombe</i> UVDE	100 μ l
4110-01K-EB	Human Apurinic/Apyrimidinic Endonuclease (hAPE)	1000 U
4120-100-EB	Human FEN-1 (Flap Endonuclease)	100 U
4145-100-EB	<i>E. coli</i> Photolyase	2000 U
4130-100-EB	Human 8-oxoGuanine DNA Glycosylase (hOGG1)	100 U
4135-100-EB	Human Ku 70/80 Complex	250 U

Concentration:
Units:
Specific Activity:

E. coli MutY
DNA Glycosylase
Catalog #: 4000-500-EB
TREVIGEN[®]
1-800-873-8443