

TREVIGEN® Product Data

For Research Use Only. Not For Use In Diagnostic Procedures

Chlorella Virus Pyrimidine Dimer Glycosylase (cv-PDG)

Catalog #: 4065-100-EB

Contents: 4065-100-01 cv-PDG **Size:** 1000 Units
 3900-500-11 10X REC™ Buffer 11 1 ml

Description: cv-PDG is a DNA glycosylase of MW 16 kDa with an associated AP lyase activity that is equally specific for the *cis-syn* and *trans-syn* isomers of cyclobutane pyrimidine dimers induced by UV irradiation. The enzyme acts through formation of a protein-DNA imino intermediate. The associated AP lyase activity occurs by a β -elimination reaction, resulting in cleavage of the phosphodiester bond 3' to the AP site and formation of a 3' unsaturated aldehyde and 5' phosphate. This site is then incised on the 5' side by an AP endonuclease, generating an appropriate substrate for resynthesis and DNA repair (see Enzyme Activity Synopsis on reverse).

Source: Purified from *E. coli* containing a recombinant plasmid harboring the *Paramecium bursaria* Chlorella virus PDG gene.

Unit Definition: One Unit is the amount of enzyme required to completely relax 250 ng of a UV-irradiated supercoiled plasmid in 30 minutes at 37°C.

Substrate Specificity: cv-PDG recognizes *cis-syn* and *trans-syn* cyclobutane pyrimidine dimers and AP sites with equal efficiency. The enzyme cleaves AP sites on both double and single stranded DNA.

Assay Conditions & Analysis: 1X REC™ Buffer 11 (25 mM sodium phosphate (pH 6.8), 1 mM EDTA, 100 mM NaCl, 1 mM DTT, 0.1 mg/ml BSA), supercoiled plasmid (250 ng) irradiated with 100 J/m² UV light, and serial dilutions of enzyme in a 20 μ l reaction volume are incubated for 30 minutes at 37°C. For analysis, 5 μ l of 5X REC™ Loading buffer (cat# 4018-250: 20 mM EDTA, 20% Ficoll, and 0.2% bromophenol blue) are added and the supercoiled, linear, and open circle forms of the plasmid are resolved by 1% agarose gel electrophoresis. Bands are visualized by ethidium bromide staining.

Storage Buffer: 25 mM sodium phosphate (pH 6.8), 100 mM NaCl, 1 mM EDTA, 1 mM DTT, 0.1 mg/ml BSA.

Storage Conditions: Store at 4°C. Do not freeze enzyme. May be diluted in 1X REC™ Buffer 11 and used immediately. To preclude loss of activity due to adsorption to plastic or glass surfaces, include BSA at 100 μ g/ml in all buffers and assays.

References: see reverse.

TREVIGEN®

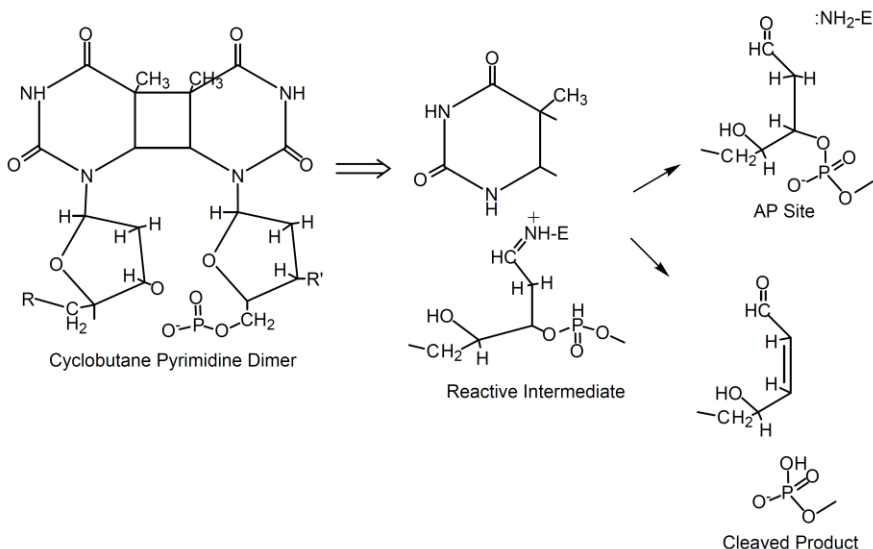
8405 Helgerman Court, Gaithersburg, MD 20877 USA

Voice: 1-800-TREVIGEN (1-800-873-8443) • 301-216-2800

Fax: 301-560-4973 • e-mail: info@trevigen.com • www.trevigen.com

References:

1. Garvish, J.F. and R.S. Lloyd. The catalytic mechanism of a pyrimidine dimer-specific glycosylase (pdg)/abasic lyase, chlorella virus-pdg. 1999. J Biol Chem **274**(14):9786-9794.
2. R.S. Lloyd. 1999. The initiation of DNA base excision repair of dipyrimidine photoproducts. Progress in Nuc Acid Res and Mol Biol **62**:155-175.
3. Garvish, J.F. and R.S. Lloyd. Active-site determination of a pyrimidine dimer glycosylase. 1999. J Mol Biol **295**(3):479-488.



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
4145-100-EB	<i>E. coli</i> Photolyase	2000 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
4125-100-EB	<i>E. coli</i> Mismatch Uracil DNA Glycosylase (Mug protein)	100 U
4130-100-EB	Human 8-oxoGuanine DNA Glycosylase (hOGG1)	100 U
4135-100-EB	Human Ku 70/80 Complex	250 U

**Chlorella Virus
Pyrimidine Dimer
Glycosylase (cv-PDG)**

Catalog#: 4065-100-EB

Storage: 4 °C

TREVIGEN®
1-800-873-8443