

TREVIGEN® Product Data

For Research Use Only. Not For Use In Diagnostic Procedures

Human 8-oxoGuanine DNA Glycosylase (hOGG1)

Catalog #: 4130-100-EB

Contents: 4130-100-01	hOGG1	Size: 100 Units
3900-500-06	10X REC™ Buffer 6	1 ml

Description: Reactive oxygen species generated from such things as ionizing radiation, cellular metabolism, and chemical genotoxins cause the DNA adducts 7,8-dihydro-8-oxo-guanine (8-oxoG) and 2,6-diamino-4-hydroxy-5-formamidopyrimidine (FaPy). Human 8-oxo-guanine DNA glycosylase (hOGG1) catalyzes the removal of the 8-oxoG and FaPy through cleavage of the DNA phosphodiester bond following Schiff base chemistry. hOGG1 does not recognize the C=O of 8-oxoG as expected, but rather recognizes a proton on N7 of the nucleotide. By mispairing with adenine during replication, 8-oxoG gives rise to G:C to T:A transversions, a frequent somatic mutation in human cancers. In contrast, a FaPy lesion leads to termination of replication and, therefore, is not considered a pre-mutagenic lesion.

Source: Purified from *E. coli* containing a recombinant plasmid harboring the α -hOGG1 gene (nuclear protein).

Unit Definition: One Unit cleaves 1 pmole of a labeled oligonucleotide probe containing 8-oxoG base paired with C within a duplex oligo.

Specificity: The catalytic activity of hOGG1 is dependent upon the base the 8-oxoG is paired with in the order of C>T>G, A. hOGG1 is also catalytically active when FaPy is paired with C. FaPy is only repaired when base paired to cytosine.

Assay Conditions and Analysis: 1X REC Buffer 6 (20 mM Tris-Cl (pH 8.0), 1 mM EDTA, 1 mM DTT, 100 µg/ml BSA), 4 pmoles of labeled 8-oxo-dG oligonucleotide annealed to the compliment oligonucleotide, and serial dilutions of enzyme in a reaction volume of 20 µl are incubated for 1 hour at 37°C. For analysis, 20 µl of 2X Loading Buffer (20 mM EDTA, 95% formamide, and 0.13% bromophenol blue) are added, the samples are heated to 95°C for 10 min then fast cooled to 4°C, and the cleavage products are resolved by 20% denaturing polyacrylamide gel electrophoresis, and percent cleavage quantified.

Storage Buffer: 20 mM Tris-Cl (pH 7.8), 1.0 mM DTT, 1 mM EDTA, 100 mM NaCl, 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 to avoid repeated freeze-thawing.

References:

1. Bruner, S.D., D.P.G. Norman, and G.L. Verdine. 2000. Structural basis for recognition and repair of the endogenous mutagen 8-oxoguanine in DNA. *Nature* **403**:859-866.
2. Boiteux, S. and J.P. Radicella. 2000. The human OGG1 gene: structure, functions, and its implications in the process of carcinogenesis. *Arch Biochem Biophys* **377**:1-8.

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
4090-500-EB	Mouse 3-mA DNA Glycosylase (Aag protein)	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
4070-500-EB	Thermostable TDG Protein	500 U
4125-100-EB	<i>E. coli</i> Mismatch Uracil DNA Glycosylase	100 U
4135-250-01	Human Ku 70/80	250 U

**Human 8-oxoGuanine
DNA Glycosylase
(hOGG1)**

Catalog#: 4130-100-EB

Storage: -20 °C

TREVIGEN®
1-800-873-8443