

TREVIGEN® Product Data

For Research Use Only. Not For Use In Diagnostic Procedures

Mouse 3-Methyladenine DNA Glycosylase Type II (Aag protein)

Catalog #: 4090-500-EB

Contents: 4090-100-01 Aag protein **Size:** 5 x 100 Units
 3900-500-09 10X REC™ Buffer 9 1 ml

Description: Mouse Aag is a 36 kDa constitutively expressed (1,000-2,000 copies/cell) protein. It acts on 3-methyladenine, 3-methylguanine, 7-methylguanine, hypoxanthine, and a number of other substrates.

Source: Purified from *E. coli* containing a recombinant plasmid encoding the mouse Aag protein.

Unit Definition: One Unit cleaves 1 pmole of a labeled oligonucleotide probe containing hypoxanthine within an oligonucleotide duplex in one hour at 37°C.

Specificity: Mouse Aag catalyzes the excision of the following forms of DNA damage: 3-methyladenine, 3-methylguanine, 7-methylguanine, hypoxanthine, and 1,N₆-ethenoadenine. It may also function on the following forms of DNA damage: 7- and 3-ethylpurines, 1-carboxyethyladenine, 7-carboxyethylguanine, O₂-methylpyrimidines, 7(2-ethoxyethyl)guanine, 7(2-hydroxyethyl)guanine, 7(2-chloroethyl)guanine, 1,2-bis(7-guanyl)ethane, 3-ethylthioethylpurines, N₂,3-ethenoguanine, N₂,3-ethanoguanine, 5-hydroxymethyluracil, 5-formyluracil, 3,N₄-ethenocytosine, 1,N₂-ethenoguanine, 3,N₂-ethenoguanine, chloroacetaldehyde cyclic adducts.

Assay Conditions: 1X REC™ Buffer 9 (10 mM HEPES-KOH (pH 7.4), 100 mM KCl, 1 mM EDTA, 1 mM EGTA, and 0.1 mM DTT), 4 pmole of labeled hypoxanthine 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. To complete cleavage of abasic site, fresh 1N 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, 95% formamide, and 0.13% bromophenol blue) are added, the samples are heated to 95°C for 10 minutes then fast cooled to 4°C, and the cleavage products are resolved by 20% denaturing polyacrylamide gel electrophoresis, and percent cleavage quantified.

Storage Buffer: 10 mM HEPES-KOH (pH 7.4), 100 mM KCl, 1 mM EDTA, 1 mM DTT, 0.2 mg/ml BSA, and 50% (v/v) glycerol.

Storage Conditions: Store at -20°C in a manual defrost freezer.

References:

1. Mattes, W.B., C.-S. Lee, J. Laval, and T.R. O'Conner. 1996. Excision of DNA adducts of nitrogen mustards by bacterial and mammalian 3-methyladenine-DNA glycosylases. *Carcinogenesis* 17:643-648.
2. Samson, L., B. Derfler, M. Boosalis, and K. Call. 1991. Cloning and characterization of a 3-methyladenine DNA glycosylase cDNA from human cells whose gene maps to chromosome 16. *Proc. Natl. Acad. Sci. USA* 88:9127-9131.

© 2013 Trevigen, Inc. All Rights Reserved. Trevigen is a registered trademark, and REC is a trademark of Trevigen, Inc.

8405 Helgerman Court, Gaithersburg, MD 20877 USA

TREVIGEN®

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 enzyme	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
4130-100-EB	Human 8-oxoGuanine DNA Glycosylase (hOGG1)	100 U
4135-250-01	Human Ku 70/80 Complex	250 U

Mouse 3-Methyladenine
DNA Glycosylase Type II
(Aag protein)

Catalog #: 4090-500-EB

TREVIGEN[®]
1-800-873-8443