

TREVIGEN® Product Data

For Research Use Only. Not For Use In Diagnostic Procedures

Human FEN-1 (Flap Endonuclease)

Catalog #: 4120-100-EB

| | | |
|------------------|---|------------------------|
| Contents: | 4120-100-01 Human FEN-1 | Size: 100 Units |
| | 4120-100-02 10X BSA Additive | 500 µl |
| | 3900-500-12 10X REC™ Reaction Buffer 12 | 1 ml |

Description: FEN-1 is a 50 kDa endonuclease/exonuclease that functions in mammalian replication and repair. The enzyme shows similarity to the yeast Rad2 and Rad13 genes. FEN-1 was identified as a necessary component of Okazaki fragment processing and functions on a number of branched DNA structures. It has been shown that FEN-1 specifically associates with PCNA and this binding stimulates FEN-1 up to 50 fold.

Source: Purified from *E. coli* containing a recombinant plasmid encoding the human FEN-1 protein.

Unit Definition: One Unit cleaves 10 pmole of a ³²P-labeled oligonucleotide flap complex in one hour at 30°C.

Specificity: FEN-1 catalyzes the nucleotide excision of DNA branch structures formed during replication of the lagging strand of DNA. The branch structures are formed due to the frequent mismatches that occur at the 5'-end of Okazaki fragments. FEN-1 is also involved in the endonucleolytic cleavage of branch structures that are formed during DNA double strand break end-joining. FEN-1 cleaves DNA flap structures one nucleotide distal to the elbow in the single-stranded region or one nucleotide proximal to the elbow in the double-stranded region, creating two possible cleavage products (see figure overleaf). FEN-1 also has a 5'-3' exonuclease activity specific for recessed 5' ends. This exonuclease activity may be involved in the removal of initiator RNA of mammalian Okazaki fragments.

Assay Conditions: The flap substrate is prepared by hybridizing 100 pmole Flap Oligonucleotide labeled with ³²P, 200 pmole Bridge Oligonucleotide, and 200 pmole Adjacent Oligonucleotide in 20 mM Tris (pH 7.4), 15 mM NaCl in a 100 µl volume at 99°C for 10 minutes and then cooling slowly to 4°C. 1X REC™ Reaction Buffer 12, 1X BSA Additive (0.1 mg/ml BSA, 5% glycerol), 8 µl of the above ³²P flap substrate, and serial dilutions of enzyme in a reaction volume of 20 µl are incubated for 1 hour at 30°C. For analysis, 5 µl of 5X REC™ Loading Buffer (Cat# 4018-250: 20 mM EDTA, 20% Ficoll, and 0.2% bromophenol blue) are added, and cleavage products are resolved by 20% denaturing polyacrylamide gel electrophoresis. The bands are cut out and radioactivity counted to quantify the cleavage products.

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

Storage Buffer: 50 mM Tris-HCl (pH 8.0), 50 mM NaCl, 1 mM DTT, 0.1 mg/ml BSA, and 50% (v/v) glycerol.

© 2010 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

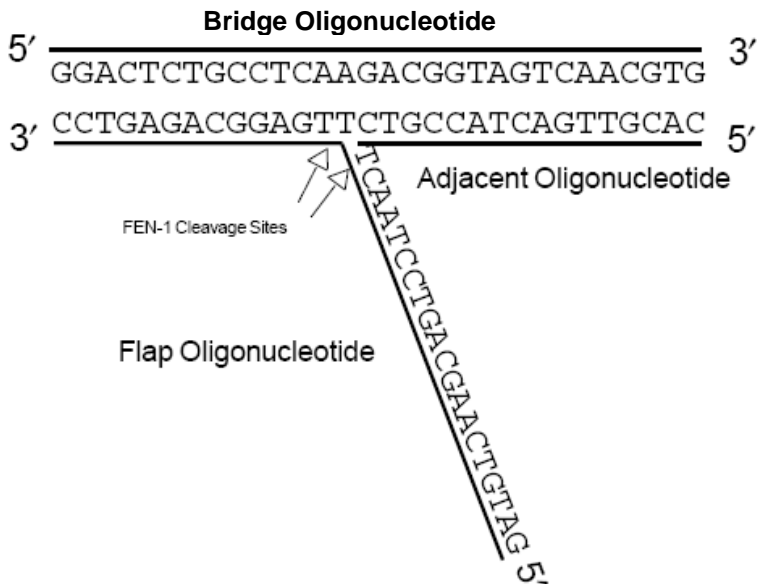
TREVIGEN®

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

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

References:

1. Harrington, J.J., and M.R. Lieber. 1994. The characterization of a mammalian DNA structure-specific endonuclease. *EMBO Journal* **13**:1235-1246.
2. Wu, X., J. Li, X. Li, C.-L. Hsieh, P.M.J. Burgers, and M.R. Lieber. 1996. Processing of branched DNA intermediates by a complex of human FEN-1 and PCNA. *Nucleic Acids Res.* **24**:2036-2043.



Notice: The supply of this Product by Trevigen conveys to you only a limited, non-transferable right to use this Product (or materials made from this Product) for purposes other than therapeutic and prophylactic applications and/or products, including, but not limited to drug screening or development. Should you desire to use this Product (or materials made from this Product) for therapeutic or prophylactic applications or products, please contact Athersys, Inc., Cleveland, Ohio, U.S.A. (www.athersys.com).

**Human Fen-1
(FLAP Endonuclease)**
Catalog #: 4120-100-EB
Storage: -20 °C
TREVIGEN®
1-800-873-8443