

# TREVIGEN® Product Data

For Research Use Only. Not For Use In Diagnostic Procedures

## Human DNA Polymerase $\beta$ Kit

Catalog#: 4020-100-K

### Contents:

Cat#	Description	Qty	Concentration
4020-100-01	Beta-polymerase	100 U	Lot specific
4020-050-02	Beta-pol Control DNA	10 $\mu$ l	0.5mg/ml
3900-200-08	10X REC™ Buffer 8	1 ml	see below*
4020-050-04	Aphidicolin	10 $\mu$ l	1 mM
4019-1	REC™ water	1 ml	N/A
4018-250	5X REC™ Loading buffer	250 $\mu$ l	see below**

\*10X REC™ Buffer 8 consists of 500 mM Tris-Cl (pH 8.8), 100 mM KCl, 100 mM MgCl<sub>2</sub>, 10 mM DTT, 10% Glycerol

\*\*20% Ficoll® 400; 20 mM EDTA, pH 8.0; 0.25% Bromophenol Blue.

**Description:** Human DNA Polymerase  $\beta$  is constitutively expressed in cells and functions by filling in gaps in DNA that are formed following base excision repair. The activity of DNA Polymerase  $\beta$  is not affected by aphidicolin, an inhibitor of DNA polymerases  $\alpha$ ,  $\delta$ , and  $\epsilon$ .

**Source:** Purified from *E. coli* containing a recombinant plasmid harboring the human DNA polymerase  $\beta$  gene.

**Unit Definition:** One Unit is the amount of enzyme required to catalyze the incorporation of 1 nmole of dNTP into an acid-insoluble form in 1 hour at 37°C.

**Specificity:** The enzyme can fill small gaps (up to 6 nucleotides) and nicks in DNA, catalyze DNA synthesis after nucleotide excision repair, and release 5'-terminal deoxyribose phosphate residues from incised AP sites.

**Assay Conditions:** 1X REC Buffer 8 (50 mM Tris-Cl (pH 8.8), 10 mM MgCl<sub>2</sub>, 10 mM KCl, 1.0 mM DTT, 1% glycerol), 50  $\mu$ M dCTP, 50  $\mu$ M dGTP, 50  $\mu$ M dATP, 50  $\mu$ M  $\alpha$ -<sup>32</sup>P-dTTP, and 100  $\mu$ g/ml of Activated DNA (Cat# 4667-50-06) in a reaction volume of 100  $\mu$ l are incubated for 5 min at 37°C.

**Storage Buffer:** 20 mM Tris-Cl (pH 7.8), 1.0 mM DTT, 1 mM EDTA, 100 mM NaCl, and 50% (v/v) glycerol.

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

© 2010 Trevigen, Inc. All Rights Reserved. Ficoll® 400 is a registered trademark of Pharmacia Biotech. Trevigen is a registered trademark of Trevigen, Inc. E6/8/10v1

## 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

### Application:

To use the  $\beta$ -polymerase enzyme, the following reaction can be set up with either radio-labeled  $\alpha$ -<sup>32</sup>P-dNTP or biotinylated nucleotides. The radiolabeled probe can be autoradiographed directly after electrophoresis, whereas the biotinylated nucleotide will need to be detected by Southern transfer of electrophoresed DNA, followed by Strep-HRP binding and colorimetric or chemiluminescent detection.

### Reaction:

Sample or Control DNA	1 $\mu$ l (0.5 $\mu$ g)
10X REC™ Buffer 8	1 $\mu$ l
$\beta$ -polymerase enzyme	2-5 units (diluted in 1X Buffer)
$\alpha$ - <sup>32</sup> P-dATP (3000 Ci/mMol)	10 pmol
10 mM each dCTP, dGTP, dTTP	1 $\mu$ l
REC™ water	<u>remainder</u>
Total	10 $\mu$ l

Let the reaction proceed at 37°C for one hour, then add 2  $\mu$ l of 5X REC™ Loading Buffer. Electrophorese the reaction products on a 0.8% TreviGel™ 5000/1X TAE gel until the blue dye has migrated 2/3 the distance of the gel, place onto a piece of 3MM paper, wrap in plastic wrap, (drying the gel down in a vacuum dryer is optional), then expose to autoradiography.

The treated control DNA will generate a ladder with the following molecular weights: 23.13 Kbp, 9.4 Kbp, 6.6 Kbp, 4.4 Kbp, 2.3 Kbp, 2.0 Kbp, 0.5 Kbp.

Aphidicolin inhibits DNA polymerase  $\alpha$ , but not  $\beta$  or  $\gamma$ . Therefore, it is useful to differentiate between different DNA polymerases.

### References:

1. Matsumoto, Y. and K. Kim. 1995. Excision of deoxyribose phosphate residues by DNA polymerase  $\beta$  during DNA repair. *Science* **269**:699-702.
2. Kunkel, T.A. and P.S. Alexander. 1986. The base substitution fidelity of eukaryotic DNA polymerases. *J Biol Chem* **261**:160-166.
3. Jenkins, T.M., J.K. Saxena, A. Kumar, S.H. Wilson, and E.J. Ackerman. 1992. DNA polymerase  $\beta$  and DNA synthesis in *Xenopus* oocytes and in a nuclear extract. *Science* **258**:475-478.
4. Vens C, E. Dahmen-Mooren, M. Venwijs-Janssen, W. Blyweert, L. Graversen, H. Bartelink, A.C. Begg. 2002. The role of DNA polymerase beta in determining sensitivity to ionizing radiation in human tumor cells. *Nucleic Acids Res.* **30**:2995-3004.
5. Bergoglio V, M.J. Pillaire, M. Lacroix-Triki, B. Raynaud-Messina, Y. Canitrot, A. Bieth, M. Gares, M. Wright, G. Delsol, L.A. Loeb, C. Cazaux, J.S. Hoffmann. 2002. Deregulated DNA polymerase beta induces chromosome instability and tumorigenesis. *Cancer Res.* **62**:3511-4.
6. Kedar P.S., S.J. Kim, A. Robertson, E. Hou, R. Prasad, J.K. Horton, S.H. Wilson. 2002. Direct interaction between mammalian DNA polymerase beta and proliferating cell nuclear antigen. *J Biol Chem.* **277**:31115-23.

Human DNA  
Polymerase  $\beta$  Kit  
Catalog#: 4020-100-K  
Storage: -20 °C  
**TREVIGEN®**  
1-800-873-8443