

Topoisomerase II, Alpha

Product number 78303

Relaxation of supercoiled DNA protocol

Topoisomerase II, Alpha alters the topological state of nucleic acids by passing an intact DNA helix through a transient break which generates a separate DNA helix. As a result of its double-stranded DNA passage mechanism, the enzyme can relax negatively or positively supercoiled DNA, as well as catenate/decatenate or knot/unknotted DNA molecules. Topoisomerase II, Alpha has an absolute requirement for divalent cation and ATP (or dATP).

Properties

Molecular weight: 340 kDa homodimer

Optimum pH: 7.9, rapidly loses activity below pH 7.5

Optimum temperature: 30°–37°C

Requirement for divalent cation: 5 mM Mg²⁺ is optimum for catalytic activity and 5 mM Ca²⁺ for DNA cleavage

Optimum ATP concentration: 0.5–1.0 mM

Optimum ionic strength: 100–170 mM NaCl/KCl

Inhibitors: N-ethylmaleimide, zinc, novobycin, coumermycin A₁, etoposide, amsacrine

10X Topoisomerase II Reaction Buffer

(PN 73592, included with enzyme):

100 mM Tris-HCl, pH 7.9, 500 mM NaCl, 500 mM KCl, 50 mM MgCl₂, 1 mM EDTA, 150 µg/ml BSA, 10 mM ATP

Topoisomerase II, Alpha and 10X Topoisomerase II Reaction Buffer (PN 73592) are functionally tested in the following protocol using 300 ng (0.1 pmol) pBR322 DNA at 30°C:

Relaxation of supercoiled DNA

- Mix

10X Topo II Reaction Buffer	2 µl
Supercoiled DNA	0.1–1.0 pmol
Topoisomerase II, Alpha	1–10 units
H ₂ O	to 20 µl
- Mix and incubate at 30°–37°C for 15 minutes.
- Stop the reaction by the addition of 3 µl of 7 mM EDTA or by the addition of SDS to 1%.

Documentation and support

Customer and technical support

Visit **thermofisher.com/support** for the latest in services and support, including:

- Worldwide contact telephone numbers
- Product support, including:
 - Product FAQs
 - Software, patches, and updates
- Order and web support
- Product documentation, including:
 - User guides, manuals, and protocols
 - Certificates of Analysis
 - Safety Data Sheets (SDSs; also known as MSDSs)

Note: For SDSs for reagents and chemicals from other manufacturers, contact the manufacturer.

Limited product warranty

Life Technologies Corporation and/or its affiliate(s) warrant their products as set forth in the Life Technologies' General Terms and Conditions of Sale found on Life Technologies' website at **www.thermofisher.com/us/en/home/global/terms-and-conditions.html**. If you have any questions, please contact Life Technologies at **thermofisher.com/support**.

The information in this guide is subject to change without notice.

DISCLAIMER: TO THE EXTENT ALLOWED BY LAW, LIFE TECHNOLOGIES AND/OR ITS AFFILIATE(S) WILL NOT BE LIABLE FOR SPECIAL, INCIDENTAL, INDIRECT, PUNITIVE, MULTIPLE, OR CONSEQUENTIAL DAMAGES IN CONNECTION WITH OR ARISING FROM THIS DOCUMENT, INCLUDING YOUR USE OF IT.

Important Licensing Information: These products may be covered by one or more Limited Use Label Licenses. By use of these products, you accept the terms and conditions of all applicable Limited Use Label Licenses.

Corporate entity: Life Technologies | Carlsbad, CA 92008 USA | Toll Free in USA 1.800.955.6288

©2017 Thermo Fisher Scientific Inc. All rights reserved. All trademarks are the property of Thermo Fisher Scientific and its subsidiaries unless otherwise specified. All other trademarks are properties of their respective owners.

For support visit **thermofisher.com/support** or email **techsupport@lifetech.com**

thermofisher.com

23 January 2017

ThermoFisher
SCIENTIFIC