

Supercoiled DNA Ladder

Cat. no. 15622-012	Size 25 µg	Conc. 0.25 µg/µL Store at -30°C to -10°C
Pub. Part no. 15622012.pps	MAN0001284	Rev. Date 11 Oct 2011

Description

The Supercoiled DNA Ladder contains eleven DNA plasmids for sizing supercoiled DNA from 2–16 kb. The bands may be visualized by ethidium bromide staining or by detection on Southern blots using a pBR322 β-lactamase (ampicillin-resistance) gene probe.

Storage Buffer

10 mM Tris-HCl (pH 8.0)

10 mM EDTA

Recommended Procedure

Apply approximately 0.1 µg of ladder per mm lane width. **Do not heat** before loading.

Notes

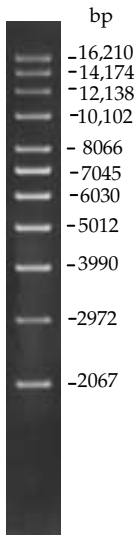
- Migration of supercoiled DNA plasmids in agarose gels can change depending on agarose concentration, buffer condition and electrophoretic field.
- For maximum accuracy, the intensity of the plasmid band to be sized should be about the same as the intensity of the bands of the supercoiled ladder. For example, apply 0.3 µg of supercoiled ladder (approximately 30 ng per band) next to 30 ng of the plasmid to be sized.

Certificate of Analysis

The Certificate of Analysis (CofA) is available at www.lifetechnologies.com/support, and it is searchable by product lot number, which is printed on each box.

Intended Use: For research use only.

Not intended for any animal or human therapeutic or diagnostic use.



0.2 µg/lane; 0.9% agarose gel stained with ethidium bromide

Limited Use Label License: Research Use Only

The purchase of this product conveys to the purchaser the limited, non-transferable right to use the purchased amount of the product only to perform internal research for the sole benefit of the purchaser. No right to resell this product or any of its components is conveyed expressly, by implication, or by estoppel. This product is for internal research purposes only and is not for use in commercial applications of any kind, including, without limitation, quality control and commercial services such as reporting the results of purchaser's activities for a fee or other form of consideration. For information on obtaining additional rights, please contact outlicensing@lifetech.com or Out Licensing, Life Technologies, 5791 Van Allen Way, Carlsbad, California 92008.

©2011 Life Technologies Corporation. All rights reserved. The trademarks mentioned herein are the property of Life Technologies Corporation or their respective owners.

For support visit www.lifetechnologies.com/support or
email techsupport@lifetech.com
www.lifetechnologies.com

