

50 bp DNA Ladder

Cat. no. 10416-014 **Size** 50 µg at 1 µg/µL **Store at** -30°C to -10°C

Doc. Part no. 10416014.pps **Pub. no.** MAN0000826 **Rev.** 2.0

Description

The 50 bp DNA Ladder consists of 16 blunt-ended fragments between 50 and 800 bp in multiples of 50 bp and an additional fragment at 2652 bp. The 350-bp band is approximately 2–3 times brighter than other ladder bands to provide internal orientation. This ladder is not designed for quantitation.

Storage Buffer

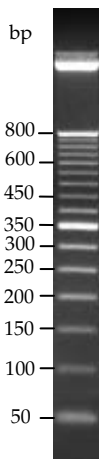
10 mM Tris-HCl (pH 7.5)
1 mM EDTA

Recommended Procedure

A final concentration of 20 mM NaCl is recommended for gel electrophoresis. Apply approximately 0.1 µg of ladder per mm lane width. *Do not heat* before loading.

Certificate of Analysis

The Certificate of Analysis provides detailed quality control and product qualification information for each product. Certificates of Analysis are available on our website. Go to **www.lifetechnologies.com/support** and search for the Certificate of Analysis by product lot number, which is printed on the box.



Example

- 50 bp DNA ladder loaded on a 2% agarose gel.
- 0.5 μ g/lane.
- Stained with ethidium bromide.

Note

The 50 bp DNA Ladder may be suitable for electrophoresis on native 6% polyacrylamide gels (Jordan, 1997).

References

Jordan, H. and Hartley, J. (1997) *FOCUS*[®] 19, 9.

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.lifetechnologies.com/termsandconditions. If you have any questions, please contact Life Technologies at www.lifetechnologies.com/support.

©2013 Life Technologies Corporation. All rights reserved. The trademarks mentioned herein are the property of Life Technologies Corporation and/or its affiliate(s) or their respective owners.

DISCLAIMER - LIFE TECHNOLOGIES CORPORATION AND/OR ITS AFFILIATE(S) DISCLAIM ALL WARRANTIES WITH RESPECT TO THIS DOCUMENT, EXPRESSED OR IMPLIED, INCLUDING BUT NOT LIMITED TO THOSE OF MERCHANTABILITY, FITNESS FOR A PARTICULAR PURPOSE, OR NON-INFRINGEMENT. TO THE EXTENT ALLOWED BY LAW, IN NO EVENT SHALL LIFE TECHNOLOGIES AND/OR ITS AFFILIATE(S) BE LIABLE, WHETHER IN CONTRACT, TORT, WARRANTY, OR UNDER ANY STATUTE OR ON ANY OTHER BASIS FOR SPECIAL, INCIDENTAL, INDIRECT, PUNITIVE, MULTIPLE OR CONSEQUENTIAL DAMAGES IN CONNECTION WITH OR ARISING FROM THIS DOCUMENT, INCLUDING BUT NOT LIMITED TO THE USE THEREOF.

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

lifetechnologies.com

25 January 2013

life
technologies™