

High DNA Mass Ladder

Cat. no. 10496-016 **Size** 200 μ L; 4 μ L per application **Store at** -30°C to -10°C

Doc. Part no. 10496016.pps

Pub. no. MAN0000844

Rev. 2.0

Description

The High DNA Mass Ladder is composed of an equimolar mixture of 6 DNA fragments of 10,000; 6000; 4000; 3000; 2000; and 1000 bp. Electrophoresis of 4 μ L of the High DNA Mass Ladder results in bands containing 200, 120, 80, 60, 40, and 20 ng of DNA, respectively.

Storage Buffer

10 mM Tris-HCl (pH 7.5)

1 mM EDTA

Recommended Procedure

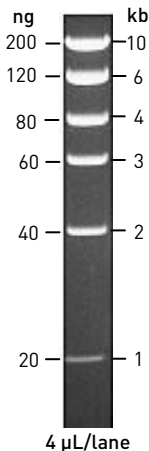
Mix 4 volumes of High DNA Mass Ladder with 1 volume of gel loading buffer (e.g. containing glycerol for density and bromophenol blue as a tracking dye). Apply an appropriate volume of ladder (see the table on page 2) to the gel loading dye to estimate the mass of unknown DNA samples. For reliable comparison of band intensities, the loading volume and buffer of the High DNA Mass Ladder and the experimental sample should be the same. Smaller volumes routinely give sharper bands. Do not heat the ladder before loading.

Note: The closer the size of the sample band relative to the band of comparable intensity in the High DNA Mass Ladder, the more accurate the mass estimation will be.

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.

For Research Use Only. Not for use in diagnostic procedures.



1% agarose gel stained
with ethidium bromide

Amount of DNA (ng) in each band			
Fragment size	2 µL	4 µL	8 µL
10,000 bp	100 ng	200 ng	400 ng
6000 bp	60 ng	120 ng	240 ng
4000 bp	40 ng	80 ng	160 ng
3000 bp	30 ng	60 ng	120 ng
2000 bp	20 ng	40 ng	80 ng
1000 bp	10 ng	20 ng	40 ng

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

22 August 2013

life
technologies™