

NoLimits 100 bp DNA Fragment

Catalog Number SM1441

Pub. No. MAN0013062 Rev. B.00



WARNING! Read the Safety Data Sheets (SDSs) and follow the handling instructions. Wear appropriate protective eyewear, clothing, and gloves. Safety Data Sheets (SDSs) are available from thermofisher.com/support.

Contents and storage

Contents	Amount	Storage
NoLimits 100 bp DNA Fragment	10 µg, 0.5 µg/µL	-25 °C to -15 °C

Description

Thermo Scientific™ NoLimits™ DNA Fragment is a chromatography-purified, exceptionally pure individual DNA fragment.

It is produced using specifically designed plasmid DNA purified by a patented technology. Plasmid DNA is digested with restriction endonucleases and the individual DNA fragment is chromatography-purified from the digestion mixture. The NoLimits™ DNA Fragment is suitable for gel and capillary electrophoresis, DNA quantification, HPLC analysis and other applications that require high purity and custom composition of DNA standards.

Convenient concentration of NoLimits™ DNA Fragments allows mixing of several individual fragments to produce a custom ladder of required range and composition.

Storage Buffer (TE buffer)

10 mM Tris-HCl (pH 7.6), 1 mM EDTA.

Important note: For NoLimits fragment dilution use TE buffer (10 mM Tris-HCl (pH 7.6), 1 mM EDTA) to avoid denaturation of the DNA fragment.

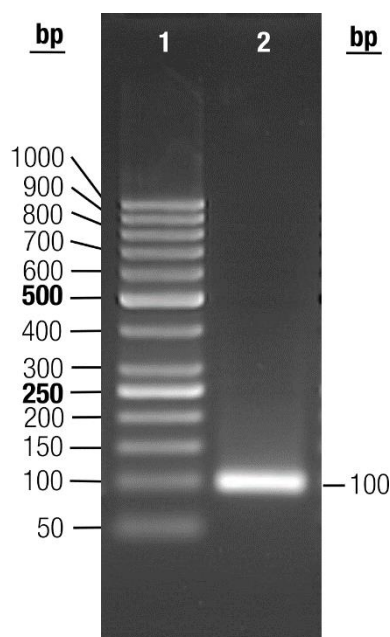


Fig.1. Analysis on agarose gel

1.1 % Agarose, 1X TAE, 7V/cm, 30 min

1 – Thermo Scientific GeneRuler 50 bp DNA Ladder – 0.5 µg

2 – NoLimits 100 bp DNA Fragment – 0.5 µg

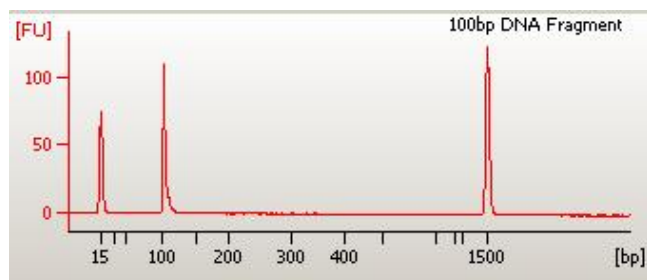


Fig.2. Analysis on Agilent Bioanalyzer 2100 and DNA 1000 LabChip® Kit

Collection of NoLimits DNA Fragments

DNA fragment size	Catalog #	Amount, µg	DNA fragment size	Catalog #	Amount, µg
10 bp	SM1391	10	850 bp	SM1661	10
15 bp	SM1381	10	900 bp	SM1491	10
20 bp	SM1401	10	1000 bp	SM1671	10
25 bp	SM1761	10	1200 bp	SM1681	10
35 bp	SM1411	10	1500 bp	SM1691	10
50 bp	SM1421	10	2000 bp	SM1701	10
75 bp	SM1431	10	2500 bp	SM1571	10
100 bp	SM1441	10	3000 bp	SM1711	10
150 bp	SM1601	10	3500 bp	SM1501	10
200 bp	SM1611	10	4000 bp	SM1721	10
250 bp	SM1451	10	5000 bp	SM1731	10
300 bp	SM1621	10	6000 bp	SM1511	10
400 bp	SM1631	10	7000 bp	SM1741	10
500 bp	SM1641	10	8000 bp	SM1521	10
600 bp	SM1461	10	10000 bp	SM1751	10
700 bp	SM1651	10	15000 bp	SM1531	10
750 bp	SM1471	10	17000 bp	SM1771	10
800 bp	SM1481	10	20000 bp	SM1541	10

See www.thermofisher.com for new additions to the collection.

Visit www.thermofisher.com for ordering the NoLimits Custom DNA Ladder Service, bulk quantities and custom formulations

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 at www.thermofisher.com/us/en/home/global/terms-and-conditions.html. If you have any questions, please contact Life Technologies at www.thermofisher.com/support.



Thermo Fisher Scientific Baltics UAB | V.A. Graiciuno 8, LT-02241 Vilnius, Lithuania
For descriptions of symbols on product labels or product documents, go to thermofisher.com/symbols-definition.

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

DISCLAIMER: TO THE EXTENT ALLOWED BY LAW, THERMO FISHER SCIENTIFIC INC. 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 this product, you accept the terms and conditions of all applicable Limited Use Label Licenses.

©2019 Thermo Fisher Scientific Inc. All rights reserved. All trademarks are the property of Thermo Fisher Scientific and its subsidiaries unless otherwise specified. LabChip is a registered trademark of Caliper Life Sciences, Inc.

thermofisher.com/support | thermofisher.com/askaquestion

thermofisher.com

15 October 2019

ThermoFisher
SCIENTIFIC