GeneRuler 1 kb Plus DNA Ladder

Catalog Number SM1331, SM1332

Pub. No. MAN0013047 **Rev.** D.00

 $\hat{}$

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

Cat. No.	Contents	Amount	Storage
SM1331	GeneRuler 1 kb Plus DNA Ladder	250 (5 x 50) μg (for 500 applications), 0.5 μg/μL	
	6X TriTrack DNA Loading Dye	2 x 1 mL	-25 °C to - 15 °C
SM1332	GeneRuler 1 kb Plus DNA Ladder	1250 (25 x 50) μg (for 2500 applications), 0.5 μg/μL	
	6X TriTrack DNA Loading Dye	10 x 1 mL	

Description

Thermo Scientific™ GeneRuler™ 1 kb Plus DNA Ladder is designed for sizing and quantification of DNA fragments in agarose gels. The ladder is composed of fifteen chromatography-purified individual DNA fragments (in base pairs): 20000, 10000, 7000, **5000**, 4000, 3000, 2000, **1500**, 1000, 700, **500**, 400, 300, 200, 75. It contains three reference bands (5000, 1500 and 500 bp) for easy orientation.

The ladder is dissolved in TE buffer.

Storage Buffer (TE buffer)

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

6X TriTrack DNA Loading Dye

10 mM Tris-HCl (pH 7.6), 0.03 % bromophenol blue, 0.03 % xylene cyanol FF, 0.15 % orange G, 60 % glycerol and 60 mM EDTA.

Protocol For Loading

Loading mixture for the 5 mm agarose gel lane*:

DNA ladder 1 μ L 6X TriTrack DNA Loading Dye 1 μ L Deionized water 4 μ L 6 μ L

Step 1: Mix gently Step 2: Load on the gel

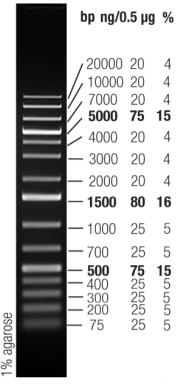
*For gels with other lane widths, the components of the mixture should be scaled either up or down. Use 0.2 μ L (0.1 μ g) of ladder per 1 mm of lane.



Recommendations

- Do not heat before loading.
- Dilute your DNA sample with the 6X TriTrack DNA Loading Dye (#R1161, supplied with the ladder): mix 1 volume of the dye solution with 5 volumes of the DNA sample;
- Load the same volumes of the DNA sample and the ladder;
- For quantification, adjust the concentration of the sample to equalize it approximately with the amount of DNA in the nearest band of the ladder.
- For DNA band visualization with SYBR™ Green and other intercalating dyes, do not add the dyes into the sample, use gel staining after electrophoresis or include dyes into agarose gel to avoid aberrant DNA migration.
- Important note: For DNA bands visualization with GelRed™ use gel staining after electrophoresis to avoid aberrant DNA migration.

GeneRuler 1 kb Plus DNA Ladder



0.5 µg/lane, 8 cm length gel, 1X TAE, 7 V/cm, 45 min

Note. Formation of diffused bands of small DNA fragments is a feature of agarose gel electrophoresis.

Limited product warranty

Life Technologies Corporation and/or it 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. GelRed is a registered trademark of Biotium Inc.

thermofisher.com/support | thermofisher.com/askaquestion

