

# GeneRuler DNA Ladder Mix, ready-to-use

Catalog Number SM0333, SM0334

Pub. No. MAN0013010 Rev. E00



**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](https://www.thermofisher.com/support).

## Contents and storage

Cat. No.	Contents	Amount	Storage
SM0333	GeneRuler DNA Ladder Mix, ready-to-use	250 µg (5 x 50) µg (for 500 applications), 0.1 µg/µL	at room temperature or at 4 °C for periods up to 24 months.*
	6X TriTrack DNA Loading Dye	2 x 1 mL	
SM0334	GeneRuler DNA Ladder Mix, ready-to-use	50 µg (for 100 applications), 0.1 µg/µL	
	6X TriTrack DNA Loading Dye	1 mL	

\*For longer periods store at -20 °C.

## Description

Thermo Scientific™ GeneRuler™ DNA Ladder Mix, ready-to-use, is designed for sizing and approximate quantification of wide range double-stranded DNA on agarose gel. The ladder is composed of 21 chromatography-purified DNA fragments (in base pairs): 10000, 8000, 6000, 5000, 4000, 3500, **3000**, 2500, 2000, 1500, 1200, **1000**, 900, 800, 700, 600, **500**, 400, 300, 200, and 100. It contains three reference bands (3000, 1000 and 500 bp) for easy orientation.

The ladder is ready to use – it is premixed with 6X TriTrack DNA Loading Dye for direct loading on gel.

## Storage and Loading Buffer

10 mM Tris-HCl (pH 7.6), 10 mM EDTA, 0.005 % bromophenol blue, 0.005 % xylene cyanol FF, 0.025 % orange G and 10 % glycerol.

## 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

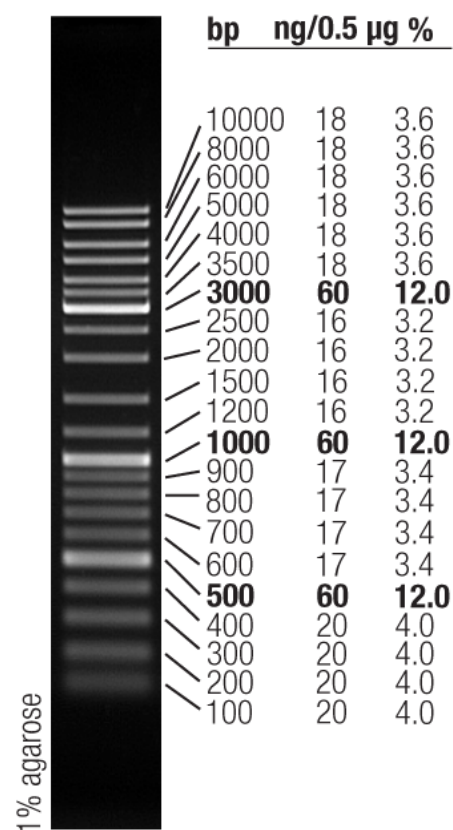
**Step 1:** Mix gently

**Step 2:** Load 1 µL per 1 mm gel 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 DNA 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<sup>™</sup> 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<sup>™</sup> use gel staining after electrophoresis to avoid aberrant DNA migration.

## GeneRuler DNA Ladder Mix, ready-to-use



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

Revision history: Pub. No. MAN0013015

Revision	Date	Description
E00	2025-04-08	Extended storage at room temperature to 24 months

## 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](http://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](http://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](http://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.

©2025 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](http://thermofisher.com/support) | [thermofisher.com/askaquestion](http://thermofisher.com/askaquestion)

[thermofisher.com](http://thermofisher.com)

08 April 2025

**ThermoFisher**  
SCIENTIFIC