

# pBR322 DNA

Catalog Number SD0041

Pub. No. MAN0012978 Rev. B.00



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

## Contents and storage

Contents	Amount	Storage
pBR322 DNA	100 µg, 0.5 µg/µL	-25 °C to -15 °C

## Description

Double-stranded closed circular medium copy plasmid, 4361 base pairs with a molecular weight of  $2.83 \times 10^6$  Da. Isolated from *E.coli* (*dam*<sup>+</sup>, *dcm*<sup>+</sup>).

## Features

- Purified by chromatography using proprietary patented technology.
- More than 90 % in the supercoiled form.

## Applications

- Cloning.
- Preparation of DNA molecular weight standards.

## Storage Buffer

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

## 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](https://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](https://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](https://www.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.

©2021 Thermo Fisher Scientific Inc. All rights reserved. All trademarks are the property of Thermo Fisher Scientific and its subsidiaries unless otherwise specified.

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

[thermofisher.com/support](https://www.thermofisher.com/support) | [thermofisher.com/askaquestion](https://www.thermofisher.com/askaquestion)  
[thermofisher.com](https://www.thermofisher.com)