

6X DNA Loading Dye

Catalog Number R0611

Pub. No. MAN0012965 Rev. C.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](https://www.thermofisher.com/support).

Contents and storage

Contents	Amount	Storage
6X DNA Loading Dye	5 x 1 mL	at room temperature or at 4 °C for periods up to 12 months. For longer periods, store at -20 °C.

Description

6X DNA Loading Dye is used for loading DNA markers and samples on agarose or polyacrylamide gels. It contains two dyes, bromophenol blue and xylene cyanol FF, for easy visual tracking of DNA migration during electrophoresis.

Composition

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

Note

- In 1 % agarose gels bromophenol blue comigrates with ~300 bp fragment and xylene cyanol FF – with ~4000 bp fragment.
- Add 1/6 volume of 6X DNA Loading Dye to your DNA sample.

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](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.

©2019 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)

ThermoFisher
SCIENTIFIC