

DNA Loading Buffer (6x) with TRIS and EDTA (Blue)

Ordering Information

[Order Online »](#)

ENZ-GEN321-LD15	1.5ml
-----------------	-------

Manuals, SDS & CofA

[View Online »](#)

Handling & Storage

Use/Stability	As indicated on product label or CoA when stored as recommended.
Handling	Spin vial briefly before opening, exercise care when dispensing DNA Loading Buffer with TRIS and EDTA (6x) (Blue) due to the intense staining of the contained tracking dyes.
Long Term Storage	+4°C
Shipping	Ambient Temperature

Regulatory Status

RUO - Research Use Only

Product Details

Formulation	Ready-to-use, sterile stock solution. 1.5ml Bromophenol Blue (0.5%), Xylene Cyanol FF (0.5%), Glycerol (40%), EDTA (Ethylenediaminetetraacetic acid) (40mM) and TRIS-HCl [Tris(hydroxymethyl)aminomethane hydrochloride] pH 7.5-8.0 (10mM) in sterile filtered double-distilled water (ddWater), without any additives.
-------------	---

Last modified: December 19, 2024



ENZO LIFE SCIENCES,
INC.
Phone: 800.942.0430
[info-
usa@enzolifesciences.com](mailto:info-usa@enzolifesciences.com)

European Sales Office
ENZO LIFE SCIENCES
(ELS) AG
Phone: +41 61 926 8989
[info-
eu@enzolifesciences.com](mailto:info-eu@enzolifesciences.com)

Belgium, The Netherlands
& Luxembourg
Phone: +32 3 466 0420
[info-
be@enzolifesciences.com](mailto:info-be@enzolifesciences.com)

France
Phone: +33 472 440 655
[info-
fr@enzolifesciences.com](mailto:info-fr@enzolifesciences.com)

Germany
Phone: +49 7621 5500 526
[info-
de@enzolifesciences.com](mailto:info-de@enzolifesciences.com)

UK & Ireland
Phone (UK customers):
0845 601 1488
Phone: +44 1392 825900
[info-
uk@enzolifesciences.com](mailto:info-uk@enzolifesciences.com)