

TRIS Buffer (1.5M)

Endotoxin-free (sterile)

Solubilisation and storage buffer for DNA and preparing other laboratory buffer reagents.

Ordering Information

[Order Online »](#)

ENZ-ACC180-0100	100ml
-----------------	-------

Manuals, SDS & CofA

[View Online »](#)

Handling & Storage

Use/Stability	As indicated on product label or CoA when stored as recommended.
Handling	Keep sterile.
Long Term Storage	+4°C
Shipping	Ambient Temperature

Regulatory Status

RUO - Research Use Only

Product Details

Endotoxin Content	<0.0002 EU/mg or <0.2 EU/ml (kinetic turbidimetric LAL method)
Formulation	Ready-to-use, sterile stock solution. Tris(hydroxymethyl)aminomethane hydrochloride (1.5M) in sterile filtered double-distilled water (ddWater), without any additives. pH value: 7.5-8.0

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)