

# TRIS Buffer (30mM) with NaCl (150mM) Endotoxin-free (sterile)

TRIS Buffer (30mM) with NaCl (150mM) Endotoxin-free (sterile) is a balanced salt solution used for a wide range of applications (cell culture, washing cells, transporting cells or tissue, diluting cells for counting, and preparing other laboratory reagents)

## Ordering Information

[Order Online »](#)

ENZ-ACC179-0100	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	<0.00002 EU/mg or <0.02 EU/ml (kinetic turbidimetric LAL method)
Formulation	Ready-to-use, sterile stock solution. TRIS [Tris(hydroxymethyl)aminomethane hydrochloride] Buffer (30mM) with NaCl [Sodium Chloride] (150mM) in sterile filtered double-distilled water (ddWater), without any additives. pH value: 7.0-7.5



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)