

HEPES Buffer (50mM) with NaCl (150mM) Endotoxin-free (sterile)

HEPES Buffer (50mM) with NaCl [Sodium Chloride] (150mM) Endotoxin-free (sterile) is a balanced salt solution used for a wide range of applications (cell culture, washing cells, transporting cells or tissue)

Ordering Information

[Order Online »](#)

ENZ-ACC181-0100	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

Formulation Ready-to-use, sterile stock solution. HEPES [4-(2-hydroxyethyl)-1-piperazineethanesulfonic acid] Buffer (50mM) with NaCl [Sodium Chloride] (150mM) in sterile filtered double-distilled water (ddWater), without any additives.
pH value: 7.2-7.6

Last modified: September 3, 2025



ENZO LIFE SCIENCES,
INC.
Phone: 800.942.0430
info-usa@enzolifesciences.com

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

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

France
Phone: +33 472 440 655
info-fr@enzolifesciences.com

Germany
Phone: +49 7621 5500 526
info-de@enzolifesciences.com

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