

# ddWater (endotoxin-free) (sterile)

Sterile, double distilled endotoxin-free water for use with TLRGRADE® or endotoxin-free grade reagents. Product has been subjected to multiple rounds of LPS removal by adsorption with activated charcoal.

Water is a ubiquitous solvent used throughout cell isolation and culture. Therefore, water contamination may affect experimental outcomes. The use of ultrapure water is advisable for culturing of cells sensitive to environmental contaminants, such as endotoxins. Cardiomyocyte isolation experiments showed that using water contaminated with bacteria and endotoxins, decreases cell viability, while using freshly produced ultrapure water dramatically increases it.

Citations: 1

[View Online »](#)

## Ordering Information

[Order Online »](#)

ALX-505-0080	1.5ml
--------------	-------

## Manuals, SDS & CofA

[View Online »](#)

## Handling & Storage

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

**Regulatory Status** RUO - Research Use Only

## Product Details

CAS	7732-18-5
Endotoxin Content	<0.00002 EU/mg or 0.02 EU/ml
MW	18.0

Last modified: October 18, 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)