

Wash buffer salts (3 packs)

Wash Buffer Salts can be used in tissue preparations, immunohistochemical staining procedures and in assays that require HRP-linked detection reagents. Each packet of salts yields 1 liter of buffered saline.

Ordering Information

[Order Online »](#)

ENZ-33802	3L
-----------	----

Manuals, SDS & CofA

[View Online »](#)

Handling & Storage

Use/Stability Store reconstituted buffer at +4°C.

Long Term Storage Ambient

Shipping Ambient Temperature

Regulatory Status RUO - Research Use Only

Product Details

Formulation Lyophilized.

Reconstitution Dissolve entire contents of packet in 1 liter of distilled or deionized water. The pH of the solution should be 7.1 ± 0.1 .

Last modified: May 29, 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)