

Streptavidin Blocker/Diluent

Ready-to-use blocking buffer and diluent, for use with SAVIEW® PLUS detection reagents

The Streptavidin Blocker/Diluent can be used to prevent non-specific binding in biotin-conjugated detection systems for IHC and ISH. It can also be used as a diluent for biotinylated antibodies.

This product is intended for qualitative immunohistochemistry and *in situ* hybridization with formalin-fixed paraffin-embedded tissue sections and frozen sections. The blocker/diluent may also be used with blood smears, cytosmears, and cell preparations.

Ordering Information

[Order Online »](#)

ENZ-ACC109-0100	100ml
-----------------	-------

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	Blue Ice

Regulatory Status

RUO - Research Use Only

Product Details

Technical Info / Product Notes

The Streptavidin Blocker/Diluent is a unique blocking solution used to prevent non-specific background staining. In addition to immunohistochemistry, the Streptavidin Blocker/Diluent can also be used for ISH.

The Streptavidin Blocker/Diluent can also be used as a diluent for SAVIEW[®] PLUS HRP reagent (ENZ-ACC102) and SAVIEW[®] PLUS AP reagent (ENZ-ACC111). It can also serve as a diluent for biotinylated antibodies.

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)