

Antibody Blocker/Diluent

Ready-to-use antibody blocking buffer and diluent, for use with POLYVIEW[®] PLUS detection reagents

The Antibody Blocker/Diluent can be used to block nonspecific staining due to antibodies during immunostaining. It can also be used as a diluent for the POLYVIEW[®] PLUS detection reagents or as a primary antibody diluent.

This product is intended for 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.

Citations: 2

[View Online »](#)

Ordering Information

[Order Online »](#)

ENZ-ACC108-0100	100ml
-----------------	-------

Manuals, SDS & CofA

[View Online »](#)

Handling & Storage

Long Term Storage +4°C

Shipping Blue Ice

Regulatory Status RUO - Research Use Only

Product Details

Technical Info / Product Notes

The Antibody Blocker/Diluent is a unique blocking solution used to reduce nonspecific background staining. It is a universal blocking agent which eliminates the need to match blocker to antibody species. In addition to immunohistochemistry, the blocker/diluent can also be used for *in situ* hybridization.

The Antibody Blocker/Diluent can also be used as a diluent for POLYVIEW[®] PLUS HRP (anti-rabbit) reagent (ENZ-ACC103), POLYVIEW[®] PLUS HRP (anti-mouse) reagent (ENZ-ACC104), and POLYVIEW[®] PLUS AP (anti-rabbit) reagent (ENZ-ACC110). It can also serve as a diluent for the primary antibody.



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)