

Avidin/Biotin Blocking Kit Set (50 ml)

Some tissue sections contain endogenous biotin, biotin binding proteins, lectins or non-specific binding substances. Tissues may also bind avidin, biotin, peroxidase biotin, peroxidase streptavidin in avidin biotin based Immunohistochemistry. These tissues will give high background in the absence of biotinylated secondary antibodies. It may be necessary to block the tissue with avidin, followed by biotin. These reagents are added before the primary antibody.

Ordering Information

[Order Online »](#)

ICT-AR-6585-02	50ml
----------------	------

Manuals, SDS & CofA

[View Online »](#)

Handling & Storage

Regulatory Status RUO - Research Use Only

Product Details

Contents Ready to use Reagent A (Avidin) 50 ml
Ready to use Reagent B (Biotin) 50 ml



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)