

HIGHDEF® blue IHC chromogen (AP)

High definition IHC chromogen producing brilliant blue color for AP-activated development.

HIGHDEF® blue IHC chromogen (AP) is intended to be used in conjunction with alkaline phosphatase (AP)-based immunostaining or in situ hybridization systems.

Application Note

[Three-color IHC Staining On Free-floating Sections From Human Brain](#)

Citations: 6

[View Online »](#)

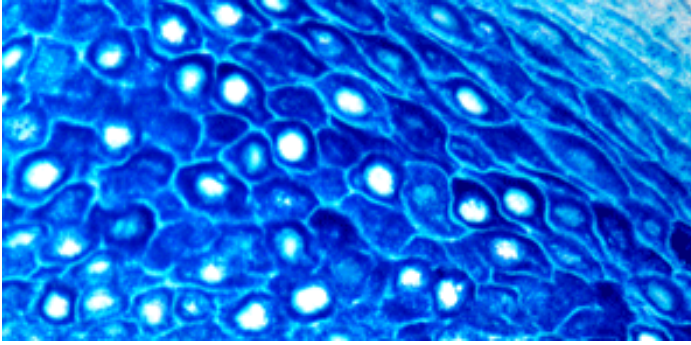
Ordering Information

[Order Online »](#)

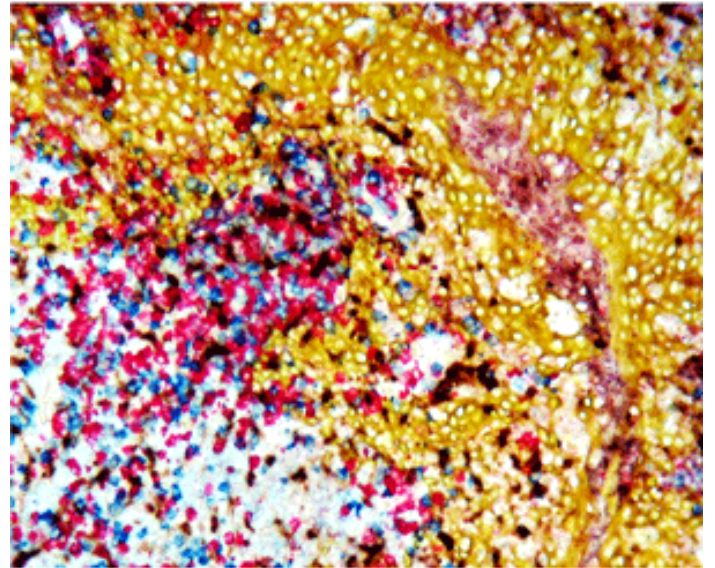
ADI-950-150-0030	30ml
------------------	------

Manuals, SDS & CofA

[View Online »](#)



Immunohistochemistry analysis: Formalin fixed paraffin embedded human tonsil stained with a high molecular weight CK antibody labeled with HIGHDEF[®] blue IHC chromogen (AP) (Prod. no. ADI-950-150) produces a distinct bright blue color.



Formalin-fixed paraffin-embedded human tonsil stained with pre-diluted mouse CD68 antibody labeled with HIGHDEF[®] IHC chromogen substrate (DAB, HC) (Prod. no. ADI-950-211, brown), pre-diluted rabbit lambda light-chain antibody labeled with HIGHDEF[®] blue IHC chromogen (AP) (Prod. no. ADI-950-150, blue), pre-diluted rabbit kappa light-chain antibody labeled with HIGHDEF[®] red IHC chromogen (AP, plus) (Prod. no. ADI-950-141, red), high molecular weight cytokeratin labeled with HIGHDEF[®] yellow IHC chromogen (HRP) (Prod. no. ADI-950-170, yellow), and methyl green nuclear counterstain.

Handling & Storage

Long Term Storage +4°C

Shipping Blue Ice

Regulatory Status RUO - Research Use Only

Product Details

Application IHC, ISH (in situ hybridization)

Contents HIGHDEF[®] blue IHC chromogen (AP) substrate buffer and chromogen, Empty mixing bottle

Technical Info / Product Notes

The HIGHDEF[®] blue IHC chromogen (AP) is a substrate-chromogen system designed to be used for either IHC or ISH when utilizing alkaline phosphatase. It has been formulated for increased stability of the working solution while producing a distinct bright blue color. It is insoluble in alcohol and xylene substitutes (both aliphatic hydrocarbon and citrus based); therefore sections can be dehydrated in alcohol, cleared in xylene substitute, and permanently mounted. However, we recommend air drying slides and then permanently mounting. This chromogen substrate system may be used for both automation and manual use.

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)