

HIGHDEF® yellow IHC chromogen (HRP)

Unique high definition yellow IHC chromogen for HRP-activated development.

HIGHDEF® yellow IHC chromogen (HRP) is intended to be used in conjunction with horseradish peroxidase (HRP)-based immunostaining or *in situ* hybridization (ISH) systems. With at least 1 day working stability, this substrate/chromogen system can be used in any automation system as well as manual use.

Application Note

[Three-color IHC Staining On Free-floating Sections From Human Brain Tumor Phenotyping and Immune Cell Regulation using Multiplex IHC](#)

Citations: 14

[View Online »](#)

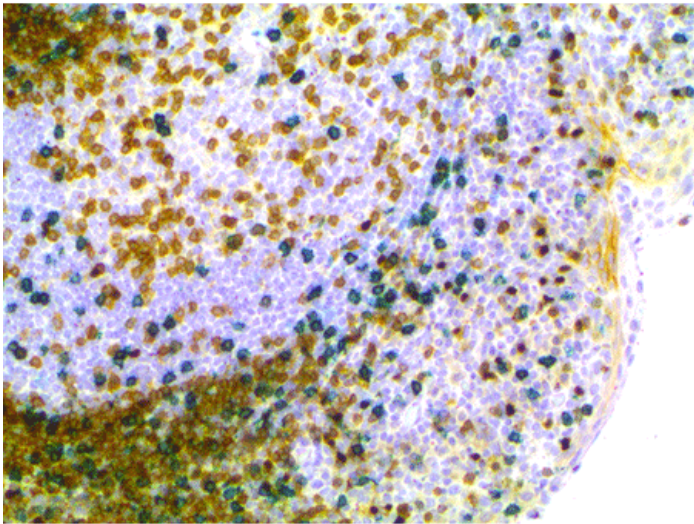
Ordering Information

[Order Online »](#)

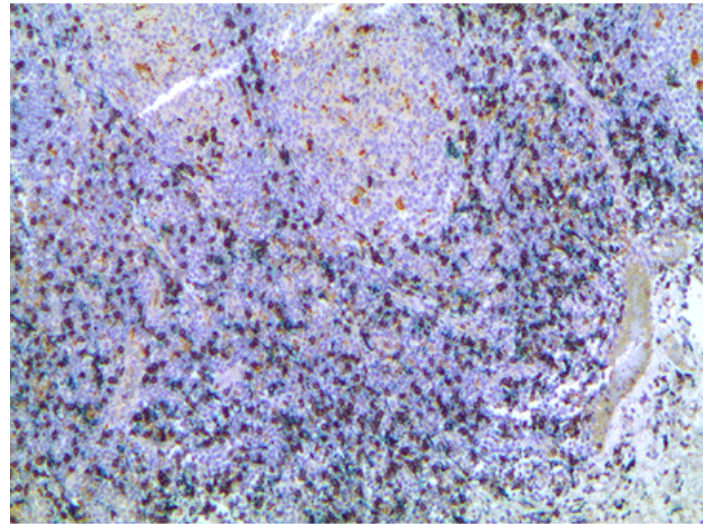
ADI-950-170-0030	30ml
------------------	------

Manuals, SDS & CofA

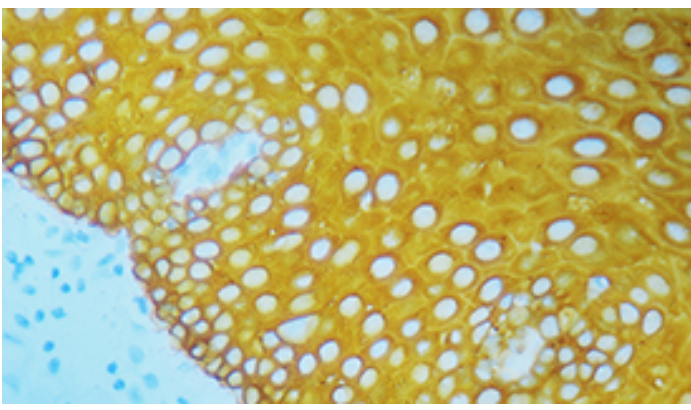
[View Online »](#)



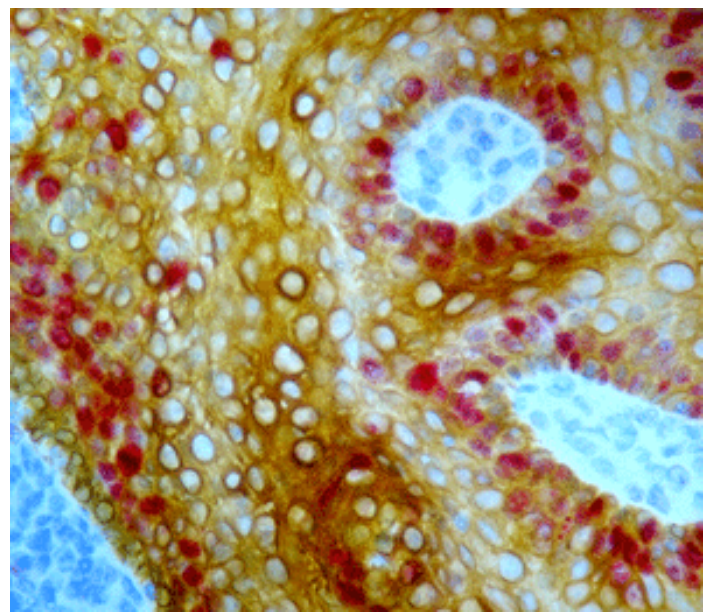
Multiplex IHC staining performed on formalin-fixed paraffin-embedded human tonsil stained with FOXP3 antibody labeled with HIGHDEF® black IHC chromogen (HRP) (Prod. Ref. ADI-950-171), CD3 antibody labeled with HIGHDEF® yellow IHC chromogen (HRP) (Prod. Ref. ADI-950-170), CD8 antibody labeled with HIGHDEF® green AP chromogen/substrate (Prod. Ref. ENZ-ACC130), and hematoxylin counterstain. Image used with permission from Plateforme de Pathologie Expérimentale et Translationnelle – PETRA, Unité AMMiCa, Gustave Roussy Cancer Campus, Villejuif, France.



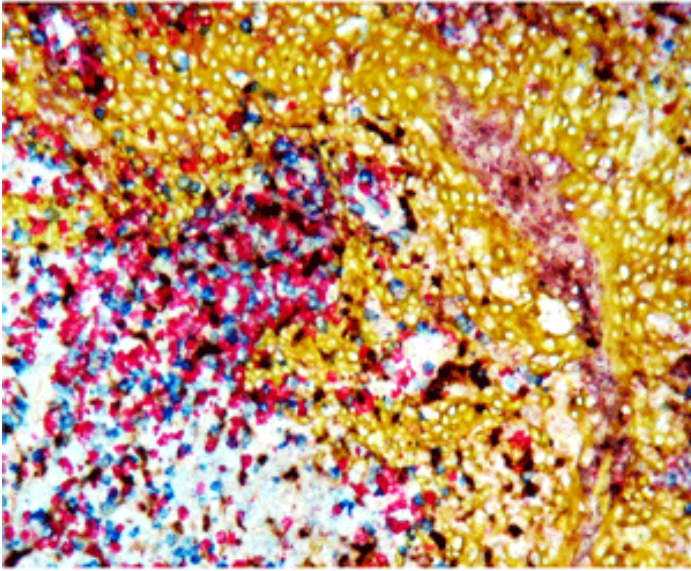
Multiplex IHC staining performed on formalin-fixed paraffin-embedded human tonsil stained with CD8 antibody labeled with HIGHDEF® black IHC chromogen (HRP) (Prod. Ref. ADI-950-171), CD68 antibody labeled with HIGHDEF® yellow IHC chromogen (HRP) (Prod. Ref. ADI-950-170), CD163 antibody labeled with HIGHDEF® green AP chromogen/substrate (Prod. Ref. ENZ-ACC130), and hematoxylin counterstain. Image courtesy from Plateforme de Pathologie Expérimentale et Translationnelle – PETRA, Unité AMMiCa, Gustave Roussy Cancer Campus, Villejuif, France.



Immunohistochemistry analysis: Formalin fixed paraffin embedded human tonsil stained with a high molecular weight CK antibody labeled with HIGHDEF® yellow IHC chromogen (HRP) (Prod. no. ADI-950-170) produces a distinct golden yellow color.



Immunohistochemical analysis: Formalin fixed paraffin embedded human tonsil double stained using MULTIVIEW® (mouse-HRP/rabbit-AP) IHC kit (Prod. no. ADI-950-100). Primary rabbit anti-Ki67 antibody was visualized with HIGHDEF® red IHC chromogen (AP) (Prod. no. ADI-950-140), while primary mouse anti-HMW CK antibody was visualized with HIGHDEF® yellow IHC chromogen (HRP) (Prod. no. ADI-950-170, sold separately), producing a distinct golden yellow color contrasting with the bright red.



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[®] yellow IHC chromogen (HRP) substrate buffer and chromogen, Empty mixing bottle

Technical Info / Product Notes

The HIGHDEF[®] yellow IHC chromogen (HRP) is a substrate-chromogen system designed to be used for either IHC or ISH when utilizing horseradish peroxidase. It produces a distinct bright yellow that can easily be distinguished from other stains.

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)