

HIGHDEF[®] black IHC chromogen (HRP)

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

HIGHDEF[®] black IHC chromogen (HRP) is intended to be used in conjunction with horseradish peroxidase (HRP)-based immunostaining or in situ hybridization 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

[Tumor Phenotyping and Immune Cell Regulation using Multiplex IHC](#)

Citations: 4

[View Online »](#)

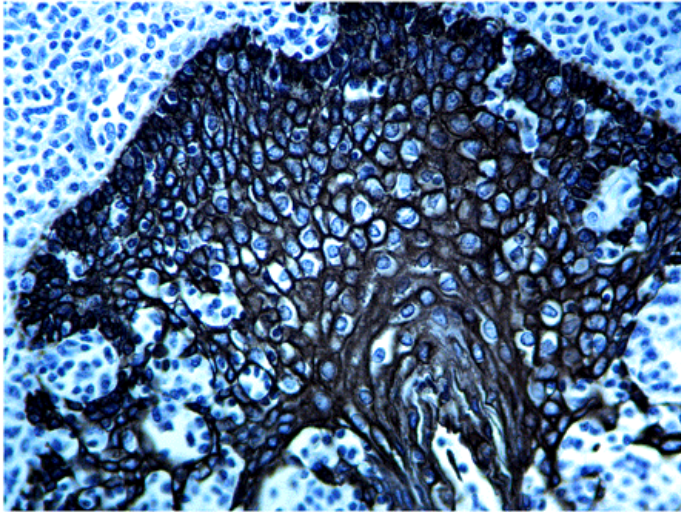
Ordering Information

[Order Online »](#)

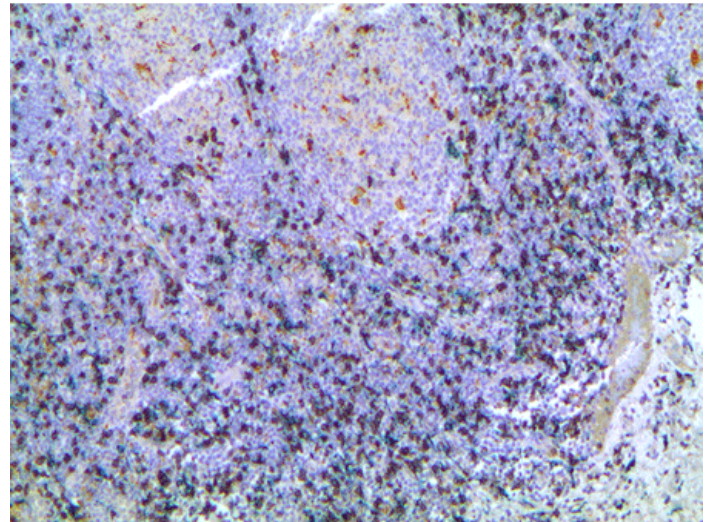
ADI-950-171-0030	30ml
------------------	------

Manuals, SDS & CofA

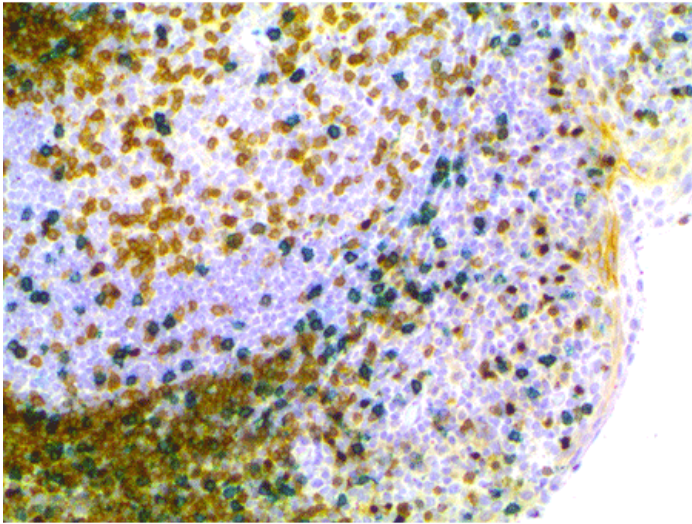
[View Online »](#)



Immunohistochemistry analysis: Formalin fixed paraffin embedded human tonsil stained with a high molecular weight cytokeratin antibody labeled and HIGHDEF® black IHC chromogen (HRP) (Prod. no. ADI-950-171) produces a sharply contrasting ebony black color.



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.



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 courtesy from Plateforme de Pathologie Expérimentale et Translationnelle – PETRA, Unité AMMICA, Gustave Roussy Cancer Campus, Villejuif, France.)

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

Technical Info / Product Notes

The HIGHDEF[®] black IHC chromogen (HRP) is a substrate-chromogen system designed to be used for either immunohistochemistry (IHC) or ISH when utilizing horseradish peroxidase. HIGHDEF[®] black IHC chromogen (HRP) produces a sharply contrasting ebony black color 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)