

HIGHDEF[®] blue IHC chromogen (HRP)

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

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

Application Note

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

Citations: 12

[View Online »](#)

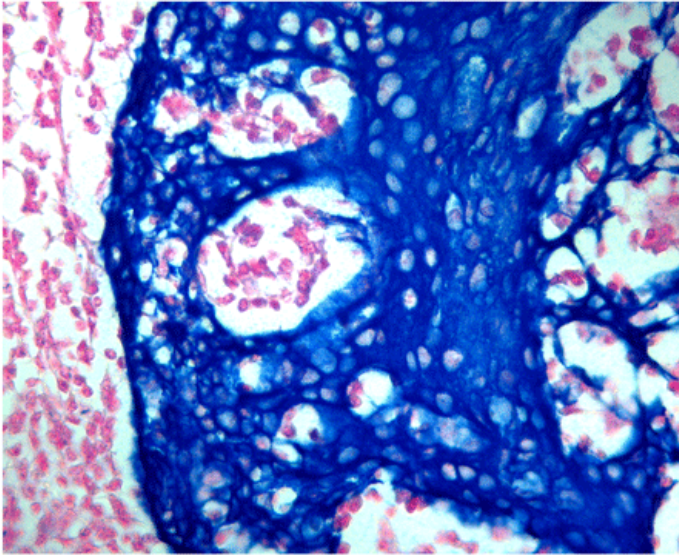
Ordering Information

[Order Online »](#)

ADI-950-151-0030	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[®] blue IHC chromogen (HRP) (Prod. no. ADI-950-151) produces a brilliant blue color.

Long Term Storage +4°C

Shipping Blue Ice

Product Details

Application	IHC, ISH (in situ hybridization)
--------------------	----------------------------------

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

Technical Info / Product Notes	The HIGHDEF[®] blue IHC chromogen (HRP) is a substrate-chromogen system designed to be used for either immunohistochemistry (IHC) or ISH when utilizing horseradish peroxidase. HIGHDEF[®] blue IHC chromogen (HRP) produces a sharply contrasting brilliant blue 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@enzolifesciences.com)