

HIGHDEF[®] Red IHC Chromogen (HRP)

The HIGHDEF[®] Red IHC Chromogen can be used in conjunction with peroxidase-based immunostaining systems. Peroxidase from the immunodetection system reacts with hydrogen peroxide substrate to degrade it, which reacts with AEC, precipitating it at positive sites yielding a reddish brown colored product.

Application Note

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

Citations: 4

[View Online »](#)

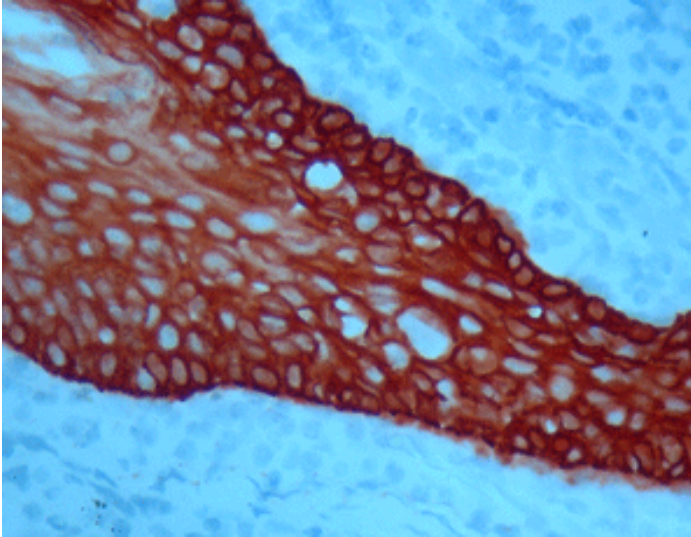
Ordering Information

[Order Online »](#)

ADI-950-210-0030	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[®] IHC substrate (AEC, stable) (Prod. no. ADI-950-210) produces a robust red color.

Handling & Storage

Handling Protect from light.

Long Term Storage +4°C

Shipping Blue Ice

Regulatory Status RUO - Research Use Only

Product Details

Appearance Amber solution.

Application IHC

Formulation Liquid. HIGHDEF[®] substrate (AEC, stable).

Technical Info / Product Notes

HIGHDEF[®] Red IHC Chromogen is a single, highly stable, AEC chromogen/substrate working solution. When used in conjunction with immunoperoxidase detection systems, AEC produces a red colored end product at positive sites, yielding strong contrast when combined with a blue hematoxylin counterstain. AEC has been well accepted among pathologists because it is less chemically hazardous than DAB. Specimens stained using HIGHDEF[®] Red IHC Chromogen cannot be dehydrated in ethanol and must be mounted in an aqueous-based mounting medium such as HIGHDEF[®] IHC mount (Prod. no. ADI-950-261).

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)