

# HIGHDEF<sup>®</sup> Red IHC Chromogen (HRP)

The HIGHDEF<sup>®</sup> 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 »](#)

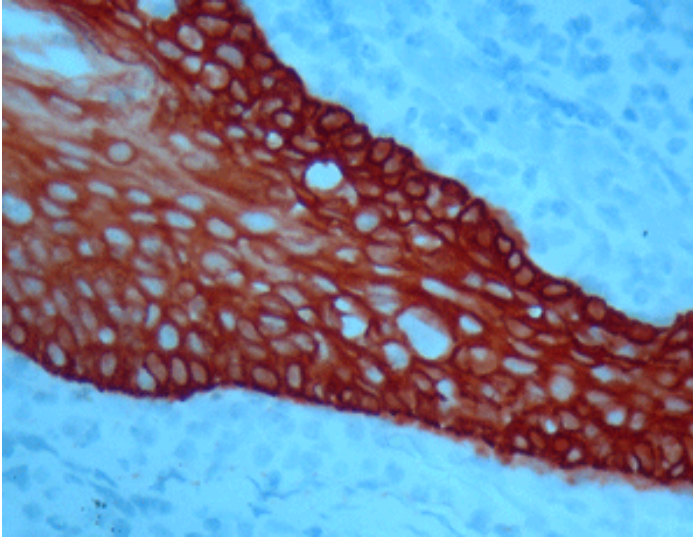
## Ordering Information

[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<sup>®</sup> 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<sup>®</sup> substrate (AEC, stable).

**Technical Info / Product Notes** HIGHDEF<sup>®</sup> 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<sup>®</sup> Red IHC Chromogen cannot be dehydrated in ethanol and must be mounted in an aqueous-based mounting medium such as HIGHDEF<sup>®</sup> IHC mount (Prod. no. ADI-950-261).



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)