

# HIGHDEF<sup>®</sup> red IHC chromogen (AP)

High definition IHC chromogen producing brilliant red color for AP-activated development.

HIGHDEF<sup>®</sup> red IHC chromogen (AP) is intended to be used in conjunction with alkaline phosphatase (AP)-based immunostaining or in situ hybridization systems.

Citations: 15

[View Online »](#)

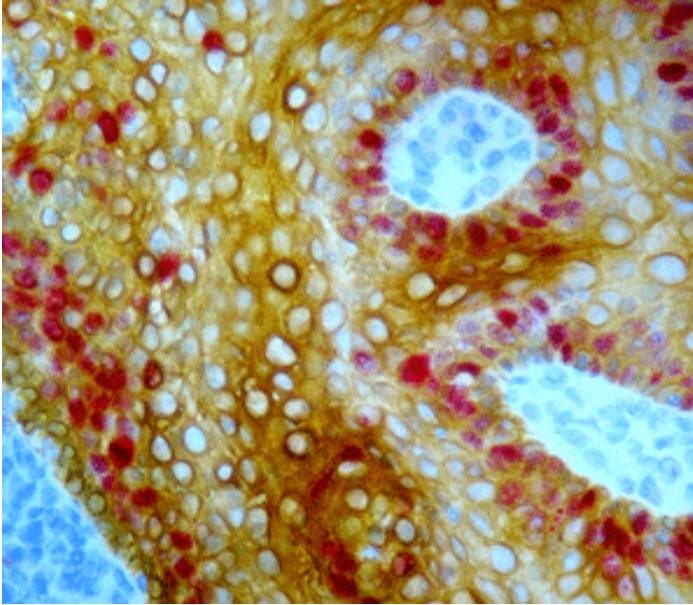
## Ordering Information

[Order Online »](#)

ADI-950-140-0030	30ml
------------------	------

## Manuals, SDS & CofA

[View Online »](#)



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

# 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<sup>®</sup> red IHC chromogen (AP) substrate buffer and chromogen, Empty mixing bottle

Technical Info / Product Notes                                    The HIGHDEF<sup>®</sup> red IHC chromogen (AP) is a substrate-chromogen system designed to be used for either IHC or ISH when using alkaline phosphatase detection. The system produces a brilliant dark red color and is insoluble in organic solvents; therefore, sections can be dehydrated in alcohol, cleared in xylene (or a xylene-substitute), and permanently mounted. This chromogen substrate system may be used for both automation and manual use.

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-  
fr@enzolifesciences.com](mailto:info-fr@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)