

HIGHDEF® IHC enhancer (DAB, plus)

The HIGHDEF® IHC enhancer (DAB, plus) is a DAB signal enhancer that can be used to increase intensity of immunostaining utilizing DAB as chromogen.

Citations: 1

[View Online »](#)

Ordering Information

[Order Online »](#)

ADI-950-182-0015	15ml
------------------	------

Manuals, SDS & CofA

[View Online »](#)

Handling & Storage

Long Term Storage Ambient

Shipping Ambient Temperature

Regulatory Status RUO - Research Use Only

Product Details

Appearance Amber solution

Application IHC

Formulation Liquid. 1X DAB chromogen enhancer.

Technical Info / Product Notes

One of the hallmarks of successful IHC is the strength of the signal generated during immunostaining. DAB, widely used as a chromogen for immunoperoxidase staining, has been well accepted by pathologists due to its superior performance compared to amino ethyl carbazol (AEC). Tissue specimens stained with DAB after only a brief incubation show increased staining intensity when the enhancer is applied. Increased sensitivity of the detection system allows the cost-effective option of diluting the primary antibody 5 to 10 fold. HIGHDEF[®] IHC enhancer (DAB, plus) increases the signal intensity without creating any background staining.

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)