

HIGHDEF® Green AP Chromogen/Substrate

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

HIGHDEF® Green AP Chromogen/Substrate is intended to be used in conjunction with alkaline phosphatase (AP)-based immunostaining or *in situ* hybridization systems.

Citations: 1

[View Online »](#)

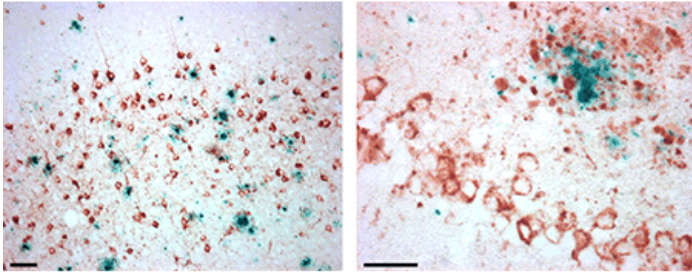
Ordering Information

[Order Online »](#)

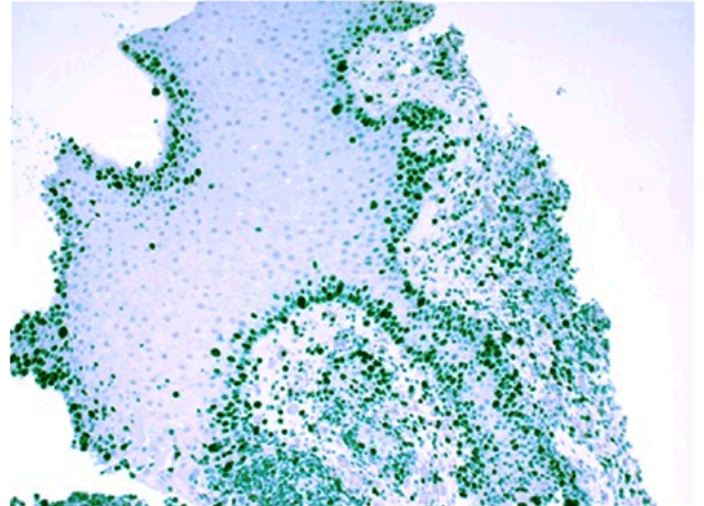
ENZ-ACC130-0030	30ml
-----------------	------

Manuals, SDS & CofA

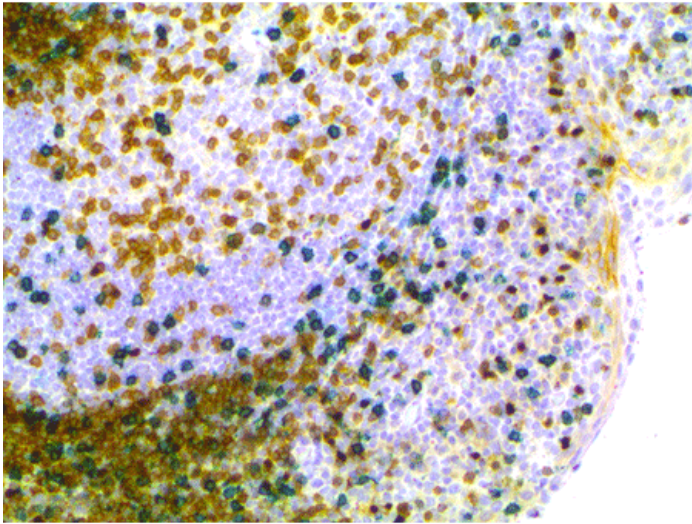
[View Online »](#)



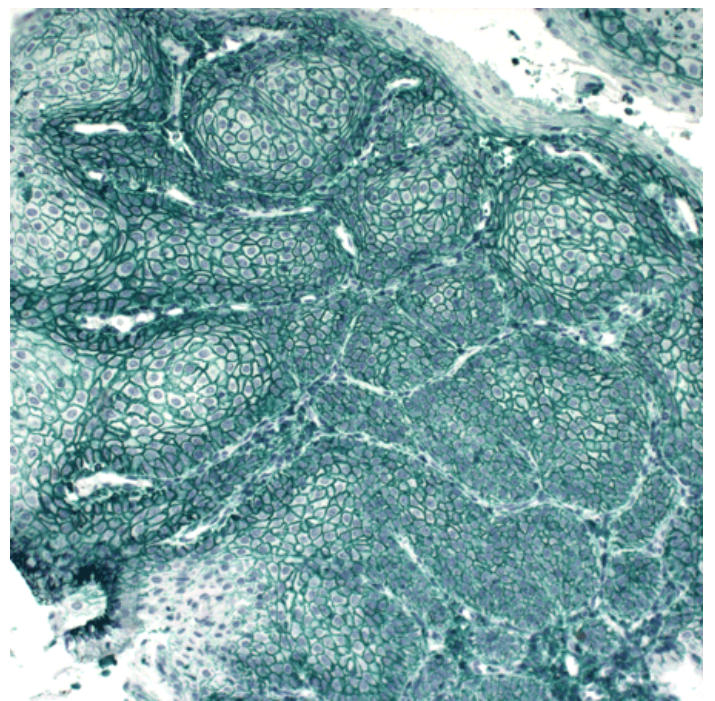
Dual IHC staining performed on formalin-fixed paraffin-embedded tissue section from human brain with Alzheimer's disease. The tissue sections were stained for alpha-amyloid with POLYVIEW® PLUS AP (anti-rabbit) reagent (Prod. No. ENZ-ACC110) and HIGHDEF® Green AP Chromogen/Substrate (Prod. No. ENZ-ACC130), and phospho-tau with POLYVIEW® PLUS HRP (anti-mouse) reagent (Prod. No. ENZ-ACC104) and HIGHDEF DAB Chromogen/Substrate Set (Prod. No. ENZ-ACC105). Picture courtesy of Dr. Karelle Leroy Laboratoire d'histologie, de neuroanatomie et de neuropathologie, Université Libre de Bruxelles, Bruxelles, Belgium.



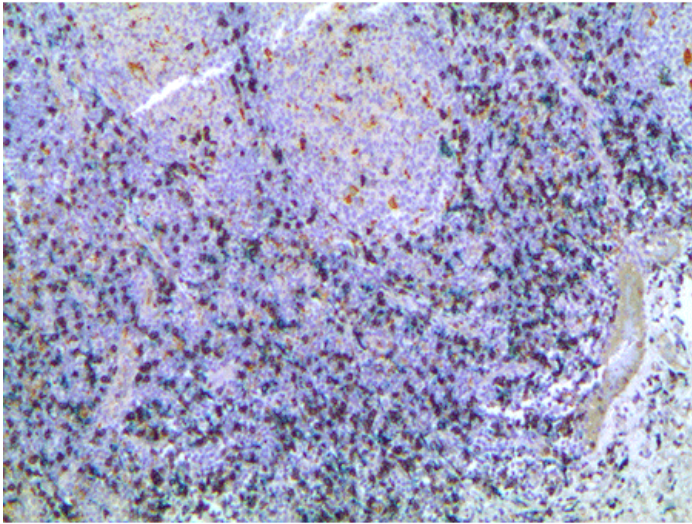
Cervical tissue incubated with a primary antibody to Ki-67, stained with POLYVIEW® PLUS (anti-Mouse) Reagent and HIGHDEF® Green AP Chromogen/Substrate



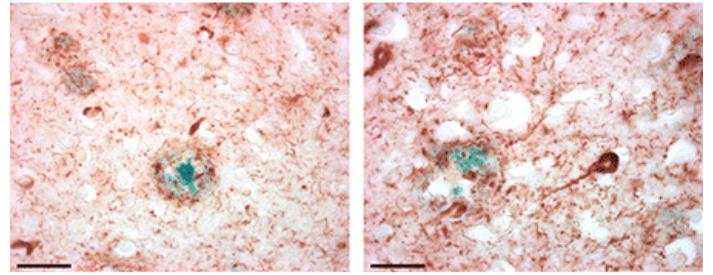
Multiplex IHC staining performed on formalin-fixed paraffin-embedded human tonsil stained with FOXP3 antibody labeled with HIGHDEF® black IHC chromogen (HRP) (Prod. Ref. ADI-950-171), CD3 antibody labeled with HIGHDEF® yellow IHC chromogen (HRP) (Prod. Ref. ADI-950-170), CD8 antibody labeled with HIGHDEF® green AP chromogen/substrate (Prod. Ref. ENZ-ACC130), and hematoxylin counterstain. Image courtesy from Plateforme de Pathologie Expérimentale et Translationnelle – PETRA, Unité AMMICA, Gustave Roussy Cancer Campus, Villejuif, France.



CD44 stained with SAVIEW® PLUS AP Reagent (ENZ-ACC111) and HIGHDEF® Green AP Chromogen/Substrate.



Multiplex IHC staining performed on formalin-fixed paraffin-embedded human tonsil stained with CD8 antibody labeled with HIGHDEF® black IHC chromogen (HRP) (Prod. Ref. ADI-950-171), CD68 antibody labeled with HIGHDEF® yellow IHC chromogen (HRP) (Prod. Ref. ADI-950-170), CD163 antibody labeled with HIGHDEF® green AP chromogen/substrate (Prod. Ref. ENZ-ACC130), and hematoxylin counterstain. Image used with permission from Plateforme de Pathologie Expérimentale et Translationnelle – PETRA, Unité AMMICA, Gustave Roussy Cancer Campus, Villejuif, France.



Dual IHC staining performed on formalin-fixed paraffin-embedded tissue section from transgenic mouse brain expressing human mutated APP and PS1. The tissue sections were stained for alpha-amyloid with POLYVIEW® PLUS AP (anti-rabbit) reagent (Prod. No. ENZ-ACC110) and HIGHDEF® Green AP Chromogen/Substrate (Prod. No. ENZ-ACC130), and phospho-tau with POLYVIEW® PLUS HRP (anti-mouse) reagent (Prod. No. ENZ-ACC104) and HIGHDEF DAB Chromogen/Substrate Set (Prod. No. ENZ-ACC105). Picture courtesy of Dr. Karelle Leroy Laboratoire d'histologie, de neuroanatomie et de neuropathologie, Université Libre de Bruxelles, Bruxelles, Belgium.

Handling & Storage

Use/Stability As indicated on product label or CoA when stored as recommended.

Long Term Storage +4°C

Shipping Blue Ice

Regulatory Status RUO - Research Use Only

Product Details

Application IHC, ISH (in situ hybridization)

Contents HIGHDEF[®] Green AP Chromogen, HIGHDEF[®] Green AP Substrate

Technical Info / Product Notes

The HIGHDEF[®] Green AP Chromogen/Substrate is a substrate-chromogen system designed to be used for either IHC or ISH when utilizing alkaline phosphatase. It has been modified to increase stability and staining intensity, producing a strong green color that is insoluble in alcohol and xylene substitutes (both aliphatic hydrocarbon and citrus based); therefore sections can be dehydrated in alcohol, cleared in xylene substitute, and permanently mounted. However, we recommend air drying slides and then permanently mounting. 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-
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)