

HIGHDEF[®] DAB

Chromogen/Substrate Set

Brown chromogen for use with HRP immunohistochemistry staining systems

The HIGHDEF[®] DAB Chromogen/Substrate Set can be used in conjunction with peroxidase-based immunostaining systems. Peroxidase from the antibody detection system reacts with hydrogen peroxide substrate to degrade it, which then reacts with DAB, precipitating it at positive sites and yielding a dark brown color.

3,3'-Diaminobenzidine (DAB), a widely used chromogen for immunoperoxidase staining, is well accepted among pathologists because of its increased sensitivity and decreased background when compared to amino ethyl carbazole (AEC). Specimens stained with DAB can be dehydrated, cleared, and mounted for permanent record keeping. HIGHDEF[®] DAB Chromogen/Substrate Set is more sensitive and stable than traditional working DAB solutions.

Citations: 14

[View Online »](#)

Ordering Information

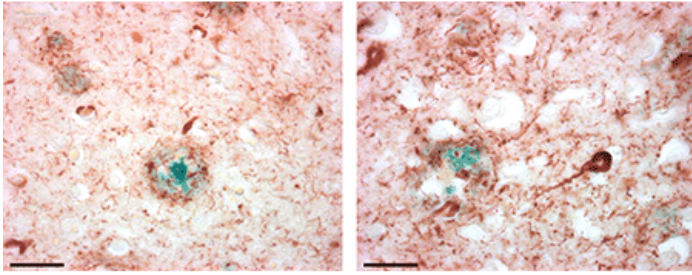
[Order Online »](#)

ENZ-ACC105-0200	200ml
-----------------	-------

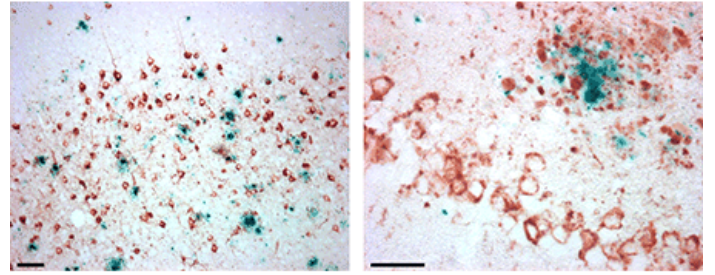
Manuals, SDS & CofA

[View Online »](#)

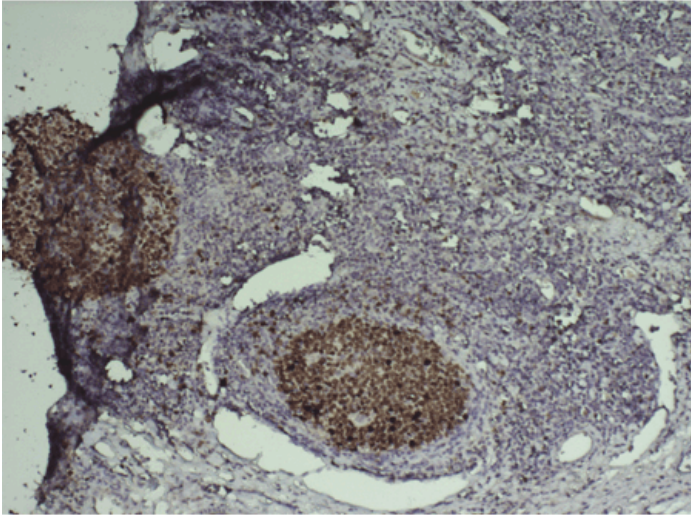
- Strong, clear color development
- Compatible with high-throughput systems



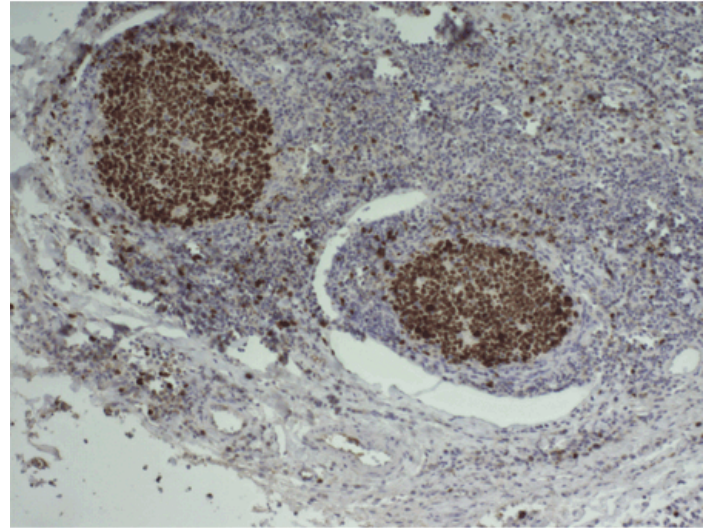
Dual IHC staining performed on formalin-fixed paraffin-embedded tissue section from transgenic mouse brain expressing human mutated APP and PS1. The tissue sections were stained for alpha-amyloid with POLYVIEW® PLUS AP (anti-rabbit) reagent (Prod. No. ENZ-ACC110) and HIGHDEF® Green AP Chromogen/Substrate (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.



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.



Cancerous tonsil tissue was incubated with an anti-Ki-67 primary antibody and then stained with POLYVIEW® PLUS HRP (anti-Rabbit) Reagent (Cat. No. ENZ-ACC104) and then color was developed using HIGHDEF® DAB Chromogen/Substrate Set.



Cancerous tonsil tissue was incubated with an anti-Ki-67 primary antibody and then stained with POLYVIEW® PLUS HRP (anti-Rabbit) Reagent (Cat. No. ENZ-ACC104) and then color was developed using HIGHDEF® DAB Chromogen/Substrate Set.

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[®] DAB Chromogen, HIGHDEF[®] DAB Substrate Buffer

Quantity Materials provided for 1,000 tests

Technical Info / Product Notes Interpretation of the results will be the sole responsibility of the user.

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)