

AEC peroxidase substrate kit

The AEC Peroxidase Substrate Kit contains reagents for visualization of peroxidase reactions using Aminoethylcarbazole (AEC). In the presence of H_2O_2 , peroxidases catalyze the oxidation of AEC into a brick red precipitate. The supplied reagents are sufficient for preparation of 300 ml of ready-to-use color reaction mixture.

Citations: 5

[View Online »](#)

Ordering Information

[Order Online »](#)

ENZ-43825	300ml
-----------	-------

Manuals, SDS & CofA

[View Online »](#)

Handling & Storage

Use/Stability Store the 8X Reaction Buffer/Substrate Reagent at +4°C.

Long Term Storage -20°C

Shipping Blue Ice

Regulatory Status RUO - Research Use Only

Product Details

Application Colorimetric detection

Contents

Aminoethylcarbazole, 6 ml 20 mg/ml in dimethylformamide

8X Reaction Buffer/Substrate Reagent, 40 ml Hydrogen peroxide in acetate buffer with stabilizers and enhancers

Mixing Bottle, 1 x 15 ml bottle for preparation of color reaction mixture

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)