

# 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: 4

[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)