

AEC Substrate Buffer (SIGNET)
Catalog Number: SIG-31030
Available Size: 2 x 50 mL



EC Rep: Cavendish Scott Ltd., P.O. Box 107, SG5 1FW, England.

Description: Substrate Buffer for use with AEC Chromogen, 2 x 50mL

(Formerly Signet Catalog No: 1030)

Intended Use: ** In Vitro Diagnostic (IVD) **

Form: Substrate Buffer

Specificity: Signet's AEC Substrate Buffer is a component of Signet's AEC Chromogen System which consists of a concentrated (50x) AEC chromogen and a substrate buffer. The AEC solution reacts to give a red staining with minimal background noise.

Uses: AEC Substrate Buffer is for use with AEC chromogen concentrate (Cat Nos: SIG-31023 and SIG-31027).

Suggested Working Dilution: The optimal working dilution should be determined for each specific assay condition.

Protocol:

1. Dilute AEC concentrate (Cat Nos: SIG-31023 & SIG-31027) 0.1:5 with Substrate Buffer
2. Apply mixture to tissue section
3. Incubate the slides for 5-10 min. at room temperature
4. Rinse tissue sections with deionized water
5. Counterstain and mount using aqueous mounting media
6. AEC induces a red precipitate with almost no background

Notes: AEC reaction product is soluble in alcohol. Do not use organic solvents after AEC staining.

Storage: Store this product between 2-8°C. AEC preparations should be stored in a manner which minimizes exposure to light.

Warranty/Conditions: Covance products may not be resold or modified for resale without prior written approval.

Related Products:

- AEC Chromogen Concentrate (SIGNET)
Catalog Number SIG-31023
- AEC Chromogen Concentrate (SIGNET)
Catalog Number SIG-31027

Product Revision Date: 5/1/2008