

**AEC Chromogen Concentrate (SIGNET)**  
**Catalog Number: SIG-31023**  
**Available Size: 2.0 mL**



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

**Description:** 50x AEC Chromogen, 2mL  
(Formerly Signet Catalog No: 1023)

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

**Form:** Chromogen Detection Reagent

**Specificity:** Signet's AEC Chromogen Systems consist of concentrated (50x) AEC chromogen and substrate buffer. The use of AEC in conjunction with peroxidase-based detection protocols results in the deposition of red chromogen deposits at reactions sites on tissue sections. The 2mL volume of 50x AEC chromogen provides enough material for 300 tests.

**Uses:** The AEC chromogen concentrate is to be diluted with the AEC chromogen substrate buffer (Cat No: SIG-31030).

**Suggested Working Dilution:**

**Protocol:**

1. Dilute AEC concentrate 0.1:5 with Substrate Buffer (Cat No: SIG-31030)
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 Substrate Buffer (SIGNET)  
Catalog Number SIG-31030
- AEC Chromogen Concentrate (SIGNET)  
Catalog Number SIG-31027
- AEC Chromogen Concentrate (SIGNET)  
Catalog Number SIG-31022

**Product Revision Date:** 5/1/2008