

**DAB Substrate Buffer with Stabilizer  
(SIGNET)**

**Catalog Number: SIG-31049**

**Available Size: 2 x 50 mL**

**Description:** DAB Substrate Buffer with Stabilizer, 2 x 50 mL  
(Formerly Signet Catalog No: 1049)

**Intended Use: \*\*Research Use Only (RUO)\*\***

This product is sold for laboratory research use only, not for human or in-vivo use.

**Form:** Chromogen Substrate Buffer

**Specificity:** Signet's DAB Substrate Buffer with Stabilizer consists of stable buffer. The DAB Substrate Buffer with Stabilizer in combination with DAB chromogen reacts with peroxidase to produce brown staining.

**Suggested Working Dilution:**

**Protocol:**

1. Add the DAB chromogen to the DAB substrate buffer as described below
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
6. DAB produces a brown precipitate with almost no background

**Notes:** Prepare the DAB solution by combining components as described in the following table.

Buffer	DAB Chromogen*
5 mL	+ 0.2 mL

25 mL + 1.0 mL

50 mL + 2.0 mL

\*Cat Nos: SIG-31041 & SIG-31043.

**Storage:** Store between 2-8°C.

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

**Related Products:**

- DAB Chromogen Concentrate (SIGNET)  
Catalog Number SIG-31041
- DAB Chromogen Concentrate (SIGNET)  
Catalog Number SIG-31043

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