

DAB Substrate Solution

Product Number: NEL938001EA

Description

This peroxidase substrate concentrate contains 3,3'-diaminobenzidine (DAB) at 25 mg/ml. 10 ml of liquid concentrate is provided. The substrate deposits a brown oxidation product in the presence of horseradish peroxidase (HRP) labeled reporters. It is useful for membrane blots, histochemistry, in situ hybridization and electron microscopy applications.

Storage Conditions and Stability

This product should be stored at 2-8°C. The concentrate is stable for a minimum of one year at 2-8°C. Discard solution if heavy precipitate develops. Nucleotides labeled with fluorophores should be protected from extended exposure to light. Minimizing freeze-thaw cycles and exposure to light are most critical factors to consider for long term usage.

Use of Reagent:

1. Warm to room temperature before use. Add 0.2 ml of DAB Substrate Solution to 10 ml of DAB Buffer (0.05M Tris-HCl, 0.1M imidazole pH 7.6 with HCl). To make 1 liter of DAB Buffer, add the following to 1 liter of distilled water:

6.06 g Tris
6.81 g Imidazole
pH with HCl to 7.6
2. Add 3.3 µl of 30% peroxide solution.
3. Mix thoroughly and use solution immediately.

Hazard Warning:

Contains hydrochloric acid (corrosive) and 3,3'-diaminobenzidine, a suspected carcinogen. Avoid contact with skin or clothing. Follow recommended disposal procedures. In case of accidental release, contain spill and clean-up with copious amounts of water.

The information provided in this document is for reference purposes only and may not be all-inclusive. Revvity, Inc., its subsidiaries, and/or affiliates (collectively, "Revvity") do not assume liability for the accuracy or completeness of the information contained herein. Users should exercise caution when handling materials as they may present unknown hazards. Revvity shall not be liable for any damages or losses resulting from handling or contact with the product, as Revvity cannot control actual methods, volumes, or conditions of use. Users are responsible for ensuring the product's suitability for their specific application. REVVITY EXPRESSLY DISCLAIMS ALL WARRANTIES, INCLUDING WARRANTIES OF MERCHANTABILITY OR FITNESS FOR A PARTICULAR PURPOSE, REGARDLESS OF WHETHER ORAL OR WRITTEN, EXPRESS OR IMPLIED, ALLEGEDLY ARISING FROM ANY USAGE OF ANY TRADE OR ANY COURSE OF DEALING, IN CONNECTION WITH THE USE OF INFORMATION CONTAINED HEREIN OR THE PRODUCT ITSELF

NEL938001EA-R REV01