

DAB Substrate Solution NEL938

Storage Conditions

This product should be stored at 2-8°C.

Product Description

This peroxidase substrate concentrate contains 3,3'-diaminobenzidine (DAB) at 25 mg/ml. 10 ml of liquid concentrate is provided.

Performance Characteristics

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.

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.

Stability

The concentrate is stable for a minimum of one year at 2-8°C. Discard solution if heavy precipitate develops.

Related Products

NEL300	4CN Plus Chromogenic Substrate	NEF812	Anti-Rabbit IgG (Goat), HRP-Labeled
NEL937	BCIP/NBT Substrate Solution	NEF822	Anti-Mouse IgG (Goat), HRP-Labeled
NEL750	Streptavidin-HRP Conjugate	NEL700	TSA Indirect For Chromogenic IHC
NEF710	Antifluorescein-HRP Conjugate	NEL761	BLAST [®] (Chromogenic) For Chromogenic Western
NEF802	Anti-Human IgG (Goat), HRP-Labeled		Blotting

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. Before using read Material Safety Data Sheet.