

ReadiUse™ NBT/BCIP Substrate Solution

Catalog number: 11631
Unit size: 50 mL

Component	Storage	Amount (Cat No. 11631)
ReadiUse™ NBT/BCIP Substrate Solution	Refrigerated (2-8 °C), Minimize light exposure	50 mL

OVERVIEW

ReadiUse™ NBT/BCIP Substrate Solution is a ready-to-use chromogenic substrate for alkaline phosphatase (AP)-based detection systems in immunohistochemistry (IHC), in situ hybridization (ISH), and blotting applications. Upon enzymatic reaction with AP, this stabilized formulation produces a permanent purple precipitate that is easily visualized by standard brightfield microscopy. The one-bottle format simplifies workflow by eliminating the need to mix multiple components. This solution supports variable incubation times, allowing both standard and high-sensitivity AP detection on fixed tissues and membrane-based platforms.

AT A GLANCE

1. Apply the ReadiUse™ NBT/BCIP Substrate Solution to the tissue sections
2. Incubate the tissue sections for 30 to 60 minutes
3. Rinse the tissue sections for 5 to 10 minutes
4. Apply the mounting medium to cover the section

Note: Before use, bring thw bottle to room temperature.

KEY PARAMETERS
Light microscope
SAMPLE EXPERIMENTAL PROTOCOL

1. After incubation with Alkaline Phosphatase (AP) detection system, rinse tissue sections with buffer, i.e TBS buffer (not provided).
2. Add enough volume of ReadiUse™ NBT/BCIP Substrate Solution (100 µL) to cover the tissue sections and incubate for 30 to 60 minutes in the dark at room temperature.
Note: Optimal development times should be determined by the end user. The incubation time of this substrate can be lengthened (up to 24 hours) providing significantly more sensitivity.
3. Wash slides in TBS buffer or water for 5 minutes.
4. For the permanent, non-aqueous mounting, dehydrate, clear and coverslip using non-aqueous mounting medium. For aqueous mounting medium, coverslip using an aqueous mounting medium such as FluoroQuest™ PLUS Antifade Mounting Medium (Cat# 20008).
5. Use the bright field to visualize the stain.

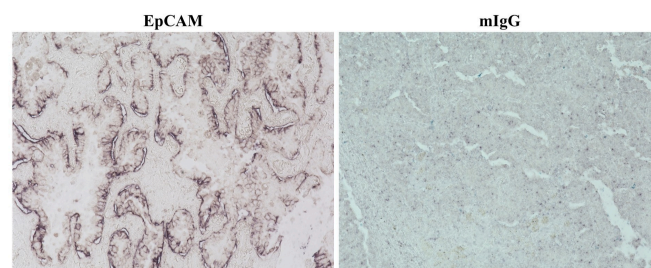
EXAMPLE DATA ANALYSIS AND FIGURES


Figure 1. Immunohistochemical detection of EpCAM in FFPE lung adenocarcinoma tissue. The tissue sections were incubated with EpCAM or Mouse IgG primary antibodies followed by ALP conjugated Goat anti-Mouse IgG and then developed with BCIP/NBT Alkaline Phosphatase (AP) substrate kit.

DISCLAIMER

AAT Bioquest provides high-quality reagents and materials for research use only. For proper handling of potentially hazardous chemicals, please consult the Safety Data Sheet (SDS) provided for the product. Chemical analysis and/or reverse engineering of any kit or its components is strictly prohibited without written permission from AAT Bioquest. Please call 408-733-1055 or email info@aatbio.com if you have any questions.