

Build a better assay with ELISA Solutions™ from ImmunoChemistry Technologies, LLC

BCIP/NBT 1-Component AP Membrane Substrate (SUBB)

A one-component formulation suitable for membrane assays using AP.

BCIP/NBT 1-Component AP Membrane Substrate (SUBB) is suitable for membrane applications using alkaline phosphatase (AP) as the conjugated detection enzyme. SUBB should not be used for microwell (ELISA) applications.

BCIP/NBT 1-Component AP Membrane Substrate is a one-component formulation containing 5-bromo-4-chloro-3-indolyl-phosphate (BCIP) and Nitroblue tetrazolium (NBT) substrate. SUBB reacts with alkaline phosphatase to yield an insoluble, dark blue reaction product. SUBB is supplied ready-to-use at 1X.

Allow BCIP/NBT 1-Component AP Membrane Substrate (SUBB) to equilibrate to room temperature (25°C) prior to use. After probing with the antibody and AP reagents, wash membrane thoroughly and transfer the membrane into a clean container. Add just enough SUBB to completely cover the membrane surface and incubate. The substrate will react with sites on the membrane containing alkaline phosphatase, producing an insoluble, permanent, dark blue reaction product.

For best results, monitor the substrate color development process until the target protein bands are visible. To stop the reaction, rinse the membrane with reagent quality water. If the reaction proceeds too long, there will be excessive background staining and diminished resolution of the target peptide or protein banding regions. If the color development is too rapid or intense, it is recommended to dilute the conjugate or shorten the conjugate incubation period.

SPECIFICATIONS

- Colorless to very light yellow liquid.
- 1X ready to use.

STORAGE

- 2-25°C.
- Protect from direct sunlight and UV sources.

SAFETY

- Warning! May cause an allergic skin reaction.
- See Safety Data Sheet (SDS) for safe handling and any precautions.
- SDS available at www.immunochemistry.com and by calling ICT.
- Not for use in diagnostic procedures.
- For research use only.

BCIP/NBT 1-COMPONENT AP MEMBRANE SUBSTRATE

Size Catalog#
100 mL 6281



HOW TO USE

1. Perform electro-blotting procedure.
2. Block membranes 4 hours - overnight.
3. Probe membranes with antibodies and AP conjugate.
4. Wash membranes after each antibody incubation step. Always transfer to a clean container for substrate development step.
5. Bring SUBB to room temperature; protect it from light.
6. Add just enough SUBB to cover the membrane surface.
7. Incubate SUBB at room temperature for several minutes.
8. Monitor the substrate color development to visualize the target peptide and protein bands.
9. Stop the color development reaction by transferring the membrane into diH₂O. Change the solution several times to ensure complete removal of all soluble BCIP/NBT components.
10. Analyze the data.

For more protocols, please visit www.immunochemistry.com.

Check out more ELISA
Solutions™ on our website
www.immunochemistry.com

