



Orders: orders@novusbio.com

Support: technical@novusbio.com

Web: www.novusbio.com

Protocols, Publications, Related Products, Reviews and more:

www.novusbio.com/NB900-78646

NB900-78646 Protocol

BCIP/NBT Alkaline Phosphatase Substrate Protocol (NB900-78646)

- 1.) Once tissue sections have been incubated with alkaline phosphatase, wash them with buffer thoroughly.
- 2.) Wipe the glass to remove excess of buffer and add enough drops of the BCIP/NBT solution to cover the tissue sections.
- 3.) Incubate for 10-30 minutes at room temperature. For the best results, look under the microscope for the signal development. Once desired signal to noise ratio is achieved, stop the reaction by washing the slides in the wash buffer.