



Orders: orders@novusbio.com

Support: technical@novusbio.com

Web: www.novusbio.com

Protocols, Publications, Related Products, Reviews and more:

www.novusbio.com/NB900-79773

NB900-79773 Protocol

AEC Chromogen/Substrate Kit Protocol (NB900-79773)

AEC Chromogen/Substrate Kit Protocol (NB900-79773):

Staining Procedure

1. Once sections have been incubated with peroxidase, wash sections with wash buffer.
2. Wipe slides to remove excess buffer. Add enough drops of AEC Chromogen/Substrate to cover tissue sections.
3. Incubate for 5-15 minutes at room temperature.
4. For best results, observe reaction under a microscope for signal development.
5. Once desired signal to noise ratio is achieved, stop reaction by washing slides in DI H₂O.