

ReadiUse™ TNB Blocking BufferCatalog number: 19999
Unit size: 100 mL

Component	Storage	Amount (Cat No. 19999)
ReadiUse™ TNB Blocking Buffer	Minimize light exposure, Refrigerated (2-8 °C)	100 mL

OVERVIEW

ReadiUse™ TNB Blocking Buffer is specifically designed to minimize background staining in applications such as tyramide signal amplification (TSA), power tyramide signal amplification (PSA), immunohistochemistry (IHC), in situ hybridization (ISH), and blotting. It contains a proprietary formulation in a buffered base, optimized to reduce non-specific interactions while preserving target signal intensity. This buffer is ready to use without further dilution and is compatible with peroxidase- and biotin-based labeling systems. It can be applied to both tissue sections and membrane formats to enhance signal clarity in high-sensitivity detection workflows. It can also be used as a dilution buffer for probes and antibodies.

SAMPLE EXPERIMENTAL PROTOCOL**For Blocking Step**

Incubate the tissue slides with enough volume to cover completely for 30 minutes at room temperature.

Note: Optimization of both incubation time and temperature is essential for best results.

For Dilutions

ReadiUse™ TNB Blocking Buffer can be used as dilution buffer for FISH probes, primary and secondary antibodies.

Note: The optimal working dilution of FISH probes and antibodies in ReadiUse™ TNB Blocking Buffer should be determined by the user for their specific experimental setup.

IMPORTANT NOTES

1. ReadiUse™ TNB Blocking Buffer is supplied in ready-to-use format. There is no need to dilute the buffer.
2. Bring the ReadiUse™ TNB Blocking Buffer to room temperature prior to use.

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.