

Alternative Block

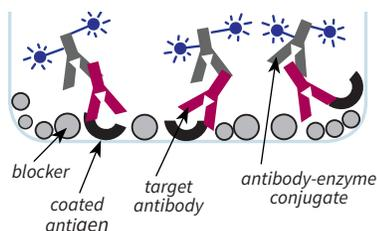
Reduces backgrounds without using proteins or detergents.

Alternative Block is a protein-free, detergent-free ELISA blocking buffer. This unique blocking buffer contains a heterogeneous mixture of proprietary blockers and synthetic stabilizers that block the uncoated regions of the plate without the use of conventional cross-reactive protein additives or detergents.

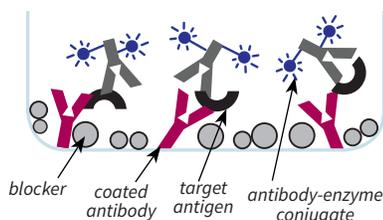
Alternative Block minimizes non-specific binding interactions and reduces background noise. It also stabilizes the coated protein during long-term storage by providing a microhydrated environment for improved retention of antigen epitope and antibody binding activity. Room temperature blocking of the plate and long-term refrigerated storage of dried plates are made possible by an antimicrobial component.

Alternative Block is suitable for use in most monoclonal and polyclonal antibody capture ELISA tests (also known as sandwich ELISAs) and peptide/protein antigen-down ELISAs. When preparing plates, the antibody or antigen is typically coated using 50-200 μ L of coating solution per well. After coating, plates are washed to remove unbound proteins and then blocked using a larger volume of blocking buffer than was used for coating, such as 300 μ L per well. This ensures that all uncoated regions inside the well are blocked. A 96-well plate blocked using this method will require 28.8 mL of blocking solution. Allow 10% extra blocking buffer to account for losses during pipetting.

Antigen-Down ELISA



Antibody Sandwich ELISA



ALTERNATIVE BLOCK

Size	Catalog #
100 mL	#6299
500 mL	#6300
1 L	#6301
10 L	#6302

INSTRUCTIONS:

1. Coat antibody or antigen onto the ELISA plate (use coating buffer catalog #645 or #6248).
2. Incubate covered plate 8-24 hours at room temperature.
3. Aspirate the coating solution.
4. Wash plate twice with ELISA Wash Buffer (catalog #652).
5. Block the uncoated regions of the ELISA plate by pipetting 300-400 μ L of blocking buffer into each well. Always use a greater volume of blocking buffer than was used for the coating solution.
6. Incubate 8-24 hours.
7. Aspirate the blocking buffer; do not wash.
8. Run the assay immediately, or dry the plate for long-term storage and seal in a foil bag (catalog #6288) with a desiccant pack (catalog #6289).

For more ELISA protocols and information, please visit www.immunochemistry.com.

SPECIFICATIONS:

- Clear liquid
- 1X ready to use
- pH 7.1-7.6

STORAGE:

- 24 months at 2-8°C
- 1 week at room temperature

SAFETY & USAGE:

- Contains $\leq 0.1\%$ sodium azide
- SDS available at immunochemistry.com
- Product intended for research use or for further manufacturing into in-vitro diagnostics reagents only.
- Not intended for use in human or therapeutics purposes.

BRIGHT MINDS, BRIGHT SOLUTIONS.

ImmunoChemistry Technologies, LLC gratefully acknowledges the significant contributions made by one of its founders, Brian W. Lee, Ph.D in the development of this product, including the creation and illustration of its strategy and protocol.



9401 James Avenue S., #155, Bloomington, MN 55431 USA

TOLL-FREE 1-800-829-3194 LOCAL 952-888-8788

FAX 952-888-8988

WWW.IMMUNOCHEMISTRY.COM