

# PhosphoBLOCKER™ Blocking Reagent

**CATALOG NUMBER:** AKR-104

**STORAGE:** Room Temperature

**QUANTITY AND CONCENTRATION:** 200 g dry blend; 5% concentration after reconstitution in 4L

## **Background**

Protein phosphorylation-dephosphorylation is one of the major signaling mechanisms for modulating the functional properties of proteins involved in gene expression, cell adhesion, cell cycle, cell proliferation, and differentiation. Proteins can be phosphorylated by protein kinases on specific serine, threonine, or tyrosine residues. The utilization of anti-phosphoprotein antibodies in western blotting has become a commonly used tool for signal transduction research. Unfortunately, low levels of endogenous phosphoprotein in various cell lysates often can not be detected, even with high concentrations of antibody and long exposure times. Most commercially available western blot blockers (e.g. dry milk, serum) are sufficient to block the unreacted sites on the membrane, reducing the amount of nonspecific antibody binding during the assay; however, they are not designed to preserve phosphoprotein antigens during blotting.

Cell Biolabs' PhosphoBLOCKER™ contains a proprietary formulation that provides several advantages over conventional blockers:

- Designed specifically for phosphoprotein blotting
- Enhances low level phosphoprotein signal without increasing background
- Premixed dry blend, easy to use

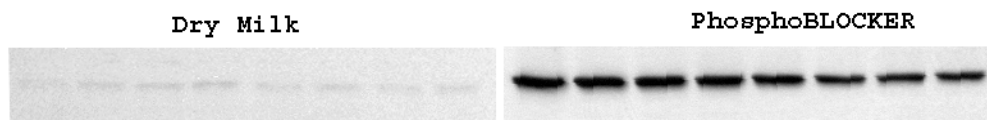
## **Methods**

Freshly prepare 5% PhosphoBLOCKER™ solution in TBST or PBST. Use the 5% PhosphoBLOCKER™ solution to block the blot. When probing the blot, use the 5% PhosphoBLOCKER™ solution to dilute primary and secondary antibodies.

*Note: Reconstituted PhosphoBLOCKER™ solution is only good for one week at 4°C.*

## **Example of results**

The following figures demonstrate typical titration results. One should use the data below for reference only. This data should not be used to interpret actual results.



**Figure 1.** Western Blot of Phospho-p38 in A549 cell lysate.

### **Recent Product Citations**

1. Martinez, P. et al. (2012). 53BP1 Deficiency Combined with Telomere Dysfunction Activates ATR-Dependent DNA Damage Response. *J.Cell.Biol.* **197**:283-300.
2. Kasuboski, J.M. et al. (2011). Zwint-1 is a Novel Aurora B Substrate Required for the Assembly of a Dynein-binding Platform on Kinetochores. *Mol. Biol. Cell.* **22**:3318-3330
3. Song, A. et al. (2010). A C. Elegans eIF4E-Family Member Upregulates Translation at Elevated Temperatures of MRNAs Encoding MSH-5 and Other Meiotic Crossover Proteins. *J. Cell Sci.* **123**:2228-2237.
4. Ishii, K. et al. (2010). Insect Cytokine Paralytic Peptide (PP) Induces Cellular and Humoral Immune Responses in the Silkworm *Bombyx mori*. *J. Biol. Chem.* **285**:28635-28642.
5. Ramakrishnan, R. et al. (2009). Characterization of Cdk9 T-loop Phosphorylation in Resting and Activated CD4+ T lymphocytes. *J. Leukoc. Biol.* **86**:1345-1350.

### **Warranty**

These products are warranted to perform as described in their labeling and in Cell Biolabs literature when used in accordance with their instructions. THERE ARE NO WARRANTIES THAT EXTEND BEYOND THIS EXPRESSED WARRANTY AND CELL BIOLABS DISCLAIMS ANY IMPLIED WARRANTY OF MERCHANTABILITY OR WARRANTY OF FITNESS FOR PARTICULAR PURPOSE. CELL BIOLABS's sole obligation and purchaser's exclusive remedy for breach of this warranty shall be, at the option of CELL BIOLABS, to repair or replace the products. In no event shall CELL BIOLABS be liable for any proximate, incidental or consequential damages in connection with the products.

***This product is for RESEARCH USE ONLY; not for use in diagnostic procedures.***

### **Contact Information**

Cell Biolabs, Inc.  
7758 Arjons Drive  
San Diego, CA 92126  
Worldwide: +1 858-271-6500  
USA Toll-Free: 1-888-CBL-0505  
E-mail: [tech@cellbiolabs.com](mailto:tech@cellbiolabs.com)  
[www.cellbiolabs.com](http://www.cellbiolabs.com)

©2004-2012: Cell Biolabs, Inc. - All rights reserved. No part of these works may be reproduced in any form without permissions in writing.