



## Product Data Sheet/Certificate of Analysis

# Bead Blocking Solution

<b>Compound:</b> Bead Blocking Solution		
<b>Storage:</b> Do not freeze! Upon receipt store at 4°C		
<b>Catalog Number:</b> S-4023		
<b>Lot Number:</b>	<b>Expiration:</b>	
<b>Components:</b> 5 mg/mL Hammarsten-grade casein in 25 mM Tris, 150 mM NaCl, pH 7.4 containing 600 ppm Kathon anti-microbial agent and 0.05% anti-foam A.	<b>Unit Size:</b>	
Test	Specification	Result
Appearance (color)	Colorless, Clear	Passed
Appearance (form)	Liquid	Passed
pH at 1X	7.4±.05	Passed

<b>QC Release Date:</b>	<b>Expiration Date:</b>
<b>Released By:</b>	

## Product Description

Bead Blocking Solution consists of biotin-free Hammarsten-grade casein (5 mg/mL) dissolved in Tris-buffered saline with a microbial preservative. During the manufacturing process this special block solution is filtered to produce a colloiddally stable reagent when stored at 4°C.

## Application

Bead Blocking Solution consists of a milk-based protein buffer solution specifically formulated to block excess non-specific protein binding sites on streptavidin magnetic beads. This solution rapidly blocks non-specific protein binding sites on NanoLink™ and MagnaLink™ Streptavidin without affecting biotin binding capacity. The use of Bead Blocking Solution in conjunction with ultra-sonication (5 minutes) also eliminates mild bead-to-bead aggregation and generates a microscopically mono-disperse bead population.

## Comments

**Do not freeze! Upon receipt store at 4°C.** Product is shipped at ambient temperature and contains an anti-microbial preservative. Not for internal or external use in humans. For research purposes only. Kathon® is a registered trademark of Rohm & Haas.

The products offered here are for research use only. Any commercial application will require a license from Solulink. The Solulink Conjugation System is patented and has multiple patents pending. Please contact Solulink for information regarding licensing information. Solulink products and methods may be covered by one or more of the following United States patents Nos. 6,686,461, 6,800,728, 7,102,024, 7,173,125, 7,462,689 and other pending patent applications. Information in this manual is subject to change without notice and does not constitute a commitment on the part of Solulink, Inc. It is supplied on an "as is" basis without any warranty of any kind, either explicit or implied. Information may be changed or updated in this manual at any time. This document may not be copied, transferred, reproduced, disclosed, or duplicated, in whole or in part, without the prior written consent of Solulink, Inc. This documentation is proprietary information and protected by the copyright laws of the United States and international treaties. The manufacturer of this documentation is Solulink, Inc.



The products offered here are for research use only. Any commercial application will require a license from Solulink. The Solulink Conjugation System is patented and has multiple patents pending. Please contact Solulink for information regarding licensing information. Solulink products and methods may be covered by one or more of the following United States patents Nos. 6,686,461, 6,800,728, 7,102,024, 7,173,125, 7,462,689 and other pending patent applications. Information in this manual is subject to change without notice and does not constitute a commitment on the part of Solulink, Inc. It is supplied on an "as is" basis without any warranty of any kind, either explicit or implied. Information may be changed or updated in this manual at any time. This document may not be copied, transferred, reproduced, disclosed, or duplicated, in whole or in part, without the prior written consent of Solulink, Inc. This documentation is proprietary information and protected by the copyright laws of the United States and international treaties. The manufacturer of this documentation is Solulink, Inc.

