

# Biotin-Catcher<sup>™</sup> Streptavidin Microwell Plates Data Sheet

## Description

The Biotin-Catcher streptavidin microwell plates are coated with highly purified recombinant streptavidin, a 60 kDa protein identified in *Streptomyces avidinii*. The purified streptavidin is immobilized to the surface of each well uniformly and in an oriented manner.

The Biotin-Catcher streptavidin plates are developed for capturing biotinylated proteins, peptides, oligonucleotides, and other biotin-conjugated molecules. Once bound to the well surface, these molecules can be used in various assays, such as DNA-protein interaction study, protein-protein interaction study, enzyme activity assays, immunoassays, etc.

- High affinity
- High specificity
- Uniformly coated
- Pre-blocked and ready to use

Cat. No.	Streptavidin Plate	Quantity	Detection Method		
			Colorimetric	Luminescent	Fluorescent
BCP-96C-05	96-well plate, clear	5 plates	$\checkmark$		$\checkmark$
BCP-96C-15	96-well plate, clear	15 plates	$\checkmark$		$\checkmark$
BCP-96B-05	96-well plate, black	5 plates	$\checkmark$	$\checkmark$	$\checkmark$
BCP-96B-15	96-well plate, black	15 plates	$\checkmark$	$\checkmark$	$\checkmark$
BCP-08C-05	8-well strip (12 strips per plate), clear	5 plates	$\checkmark$		$\checkmark$
BCP-08C-15	8-well strip (12 strips per plate), clear	15 plates	$\checkmark$		$\checkmark$
BCP-08B-05	8-well strip (12 strips per plate), black	5 plates	$\checkmark$	$\checkmark$	$\checkmark$
BCP-08B-15	8-well strip (12 strips per plate), black	15 plates		$\checkmark$	$\checkmark$

### Storage/Stability

The Biotin-Catcher<sup>™</sup> plates are stable for two years when stored at 2-8 °C in the sealed bag.

### **Coating and Blocking Volumes**

Each well of the plate was coated with 100  $\mu l$  of Streptavidin, and blocked with 300  $\mu l$  of 2% BSA.

### **Binding Capacity**

Each lot of the Biotin-Catcher plates is tested using biotinylated Avi-tagged EGFP. In PBST containing 0.1% Tween 20, over 5 pmoles of Avi-tagged EGFP (28kDa) can be captured in each well. For larger molecules the binding capacity may be less owing to steric hindrance.

#### **Binding and Washing Solutions**

Non-specific proteins or other molecules can be conveniently washed off by PBST or TBST containing 0.05-0.1%Tween 20. The pH should be 6.5-8.5.

The following reagents/chemicals may be included in the binding and washing solutions:

NaCl/KCI: 0.1-1 M Tween 20: 0.05-0.1% (w/v) EDTA: 0.5-5 mM BSA, casein, non-fat milk powder: 0.1-1% (w/v)

Although up to 1M of urea or 0.01% SDS can be tolerated by the coated streptavidin, they should be avoided in the binding and washing steps.

For more information, please contact us:

GeneCopoeia, Inc. 9620 Medical Center Drive, #101 Rockville, MD 20850 Tel: 301-762-0888 or 866-360-9531 Fax: 301-762-3888 support@genecopoeia.com www.genecopoeia.com

GeneCopoeia Products are for Research Use Only Trademarks: Biotin-Catcher™ (GeneCopoeia) Copyright © 2012 GeneCopoeia Inc. BCP-110912