

Pierce[®] Streptavidin Coated 384-Well Plates

15405 15406 15407

0880.2

Number	Description
--------	-------------

15405	Pierce Streptavidin Coated Plate (clear, 384-well), 5 each
-------	--

15406	Pierce Streptavidin Coated Plate (white, 384-well), 5 each
-------	--

15407	Pierce Streptavidin Coated Plate (black, 384-well), 5 each
-------	--

Blocking Buffer: These plates are supplied blocked with SuperBlock[®] Blocking Buffer

Binding Capacity: ~4pmol D-biotin/well

Coating Volume: 50µL

Blocking Volume: 100µL

Storage: Upon receipt store plates at 4°C in unopened pouches. Once opened, place unused plates in a resealable bag with desiccant and store at 4°C.

Introduction

The Thermo Scientific Pierce Streptavidin Coated Plates are made of polystyrene and are ideal for binding assays using biotinylated molecules. These plates are especially advantageous when direct adsorption to polystyrene plates denatures antibodies or the target molecule. Streptavidin has no carbohydrate groups and an isoelectric point of 5-6, resulting in low nonspecific interactions. The Pierce Streptavidin Coated Plates are available in clear for colorimetric assays, white for chemiluminescent assays, and black for fluorescent assays.

Example ELISA Protocol using Streptavidin Coated Plates

A. Materials Required

- Wash Buffer: Tris-buffered saline (25 mM Tris, 150 mM NaCl; pH 7.2; Product No. 28376), 0.1% BSA, 0.05% Tween[®]-20 Detergent; alternatively, use Thermo Scientific Blocker BSA (Product No. 37520) supplemented with 0.05% Tween-20 Detergent
- Biotinylated capture antibody adjusted to 10µg/mL, or other appropriate concentration, with Wash Buffer
- Antigen adjusted to appropriate concentration with Wash Buffer
- Primary antibody adjusted to appropriate concentration with Wash Buffer
- Enzyme-labeled secondary antibody adjusted to appropriate concentration with Wash Buffer
- Appropriate enzyme substrate: example substrates are the Thermo Scientific TMB Substrate Kit (Product No. 34021) for horseradish peroxidase and the Phosphatase Substrate Kit (Product No. 37620) for alkaline phosphatase

B. Method

1. Wash each well three times with 100µL of Wash Buffer. Add 50µL of the biotinylated capture antibody to each well and incubate for 2 hours at room temperature.
2. Wash each well three times with 100µL of Wash Buffer. Make a serial dilution of the antigen and add 50µL to each well. Incubate plate for 30 minutes at room temperature.
3. Wash each well three times with 100µL of Wash Buffer. Add 50µL of the primary antibody to each well and incubate plate for 30 minutes at room temperature.
4. Wash each well three times with 100µL of Wash Buffer. Add 50µL of the enzyme-labeled secondary antibody to each well. Incubate plate for 30 minutes at room temperature.

5. Wash each well three times with 100µL of Wash Buffer.
6. Follow the manufacturer's instructions for the specific detection system.

Related Thermo Scientific Products

37070	SuperSignal® ELISA Pico Chemiluminescent Substrate, 100mL, peroxidase substrate
15169	QuantaBlu™ Fluorogenic Peroxidase Substrate Kit
34028	1-Step™ Ultra TMB-ELISA, 250mL, colorimetric peroxidase substrate
37621	1-Step PNPP, 100mL, colorimetric phosphatase substrate
29339	Biotinylated Alkaline Phosphatase, 1mg
29139	Biotinylated Horseradish Peroxidase, 5mg
15075	Reagent Reservoirs, 200/pkg
15082	Microtube Racked System, 960 tubes
15036	Sealing Tape for 96-Well Plates, 100/pkg
15511	Pierce NeutrAvidin High Binding Capacity Coated Plates (clear, 384-well), 5 each
15512	Pierce NeutrAvidin High Binding Capacity Coated Plates (white, 384-well), 5 each
15513	Pierce NeutrAvidin High Binding Capacity Coated Plates (black, 384-well), 5 each

General References

- Denlinger, L.C., *et al.* (2001). Cutting Edge: The nucleotide receptor P2X7 contains multiple protein- and lipid-interaction motifs including a potential binding site for bacterial lipopolysaccharide. *J Immunol* **167**:1871-6.
- Ferre-Aubineau, V., *et al.* (1995). Colorimetric microtiter plate hybridization assay using monoclonal antibody for detection of an amplified human immunodeficiency virus target. *J Virol Meth* **55**:145-51.
- Hiller, Y. *et al.* (1987). Biotin binding to avidin. Oligosaccharide side chain not required for ligand association. *Biochem J* **248**:167-71.
- Holmstrom, K., *et al.* (1993). A highly sensitive and fast non-radioactive method for detection of polymerase chain reaction products. *Anal Biochem* **209**:278-283.
- Simon, M.D., *et al.* (2004). A phage display selection of engrailed homeodomain mutants and the importance of residue Q50. *Nucl Acid Res* **32**(12):3623-31.

Tween is a trademark of Croda International Plc.

This product ("Product") is warranted to operate or perform substantially in conformance with published Product specifications in effect at the time of sale, as set forth in the Product documentation, specifications and/or accompanying package inserts ("Documentation") and to be free from defects in material and workmanship. Unless otherwise expressly authorized in writing, Products are supplied for research use only. No claim of suitability for use in applications regulated by FDA is made. The warranty provided herein is valid only when used by properly trained individuals. Unless otherwise stated in the Documentation, this warranty is limited to one year from date of shipment when the Product is subjected to normal, proper and intended usage. This warranty does not extend to anyone other than the original purchaser of the Product ("Buyer").

No other warranties, express or implied, are granted, including without limitation, implied warranties of merchantability, fitness for any particular purpose, or non infringement. Buyer's exclusive remedy for non-conforming Products during the warranty period is limited to replacement of or refund for the non-conforming Product(s).

There is no obligation to replace Products as the result of (i) accident, disaster or event of force majeure, (ii) misuse, fault or negligence of or by Buyer, (iii) use of the Products in a manner for which they were not designed, or (iv) improper storage and handling of the Products.

Current product instructions are available at www.thermoscientific.com/pierce. For a faxed copy, call 800-874-3723 or contact your local distributor.

© 2011 Thermo Fisher Scientific Inc. All rights reserved. Unless otherwise indicated, all trademarks are property of Thermo Fisher Scientific Inc. and its subsidiaries. Printed in the USA.