

Pierce[®] Protein A, Protein G and Protein A/G Coated 96-Well Plates

0657.4

| Number | Description |
|--------|---|
| 15130 | Pierce Protein A Coated Plates (clear, 96-well), 5 each |
| 15132 | Pierce Protein A Coated Plates (clear, 8-well strips), 5 each |
| 15154 | Pierce Protein A Coated Plates (white, 96-well), 5 each |
| 15155 | Pierce Protein A Coated Plates (black, 96-well), 5 each Binding Capacity: ~4pmol rabbit IgG/well |
| 15131 | Pierce Protein G Coated Plates (clear, 96-well), 5 each |
| 15133 | Pierce Protein G Coated Plates (clear, 8-well strips), 5 each |
| 15156 | Pierce Protein G Coated Plates (white, 96-well), 5 each |
| 15157 | Pierce Protein G Coated Plates (black, 96-well), 5 each Binding Capacity: ~2pmol rabbit IgG/well |
| 15138 | Pierce Protein A/G Coated Plates (clear, 8-well strips), 5 each Binding Capacity: ~5pmol rabbit IgG/well |

Note: Plates are activated to 100µL and supplied pre-blocked with Thermo Scientific SuperBlock Blocking Buffer. Use these plates for single-antibody assays only. In multiple antibody assays, such as sandwich ELISAs, the first antibody cannot block all binding sites and subsequent antibodies will bind to the plate, resulting in false positives.

Storage: Upon receipt store plates at 4°C in unopened pouches. Place opened, unused plates in a resealable bag and store desiccated at 4°C. Plates are shipped at ambient temperature.

Introduction

Microplates coated with Proteins A, G and A/G specifically bind the Fc region of immunoglobulins resulting in antibody orientation such that the antigen binding site is directed away from the microplate and available for binding. These coated plates can increase antibody-binding capacity compared to coating antibodies directly to the plates.

These Thermo Scientific Pierce Coated Plates are useful for detecting antibodies using labeled antigens. The pre-coated plates can increase assay sensitivity and are ideal when a low amount of antibody is available. Antibodies bind to Protein A, Protein G or Protein A/G rapidly and dissociate with strong chaotropes and extreme pH. The coated plates are available in clear for colorimetric assays, white for chemiluminescent assays, and black for fluorescent assays.

The antibody being used determines whether to use Protein A, Protein G or Protein A/G coated plates. Protein G binds all subclasses of mouse, rabbit and goat IgG as well as most other commonly used species. Protein A binds strongly to rabbit antibodies and binds pig and guinea pig IgG better than Protein G; Protein A does not bind mouse IgG₁ well, and has weak binding of most rat IgG subclasses. Protein A/G is a secreted gene fusion product from a nonpathogenic form of *Bacillus* and contains four Fc-binding domains from Protein A and two from Protein G making it a more versatile tool. Please see the Pierce catalog website or for a list of binding characteristics of the different binding proteins.

Example Procedure for ELISA-based Applications

The following procedure is an example ELISA protocol. Specific applications and systems will require optimization. Do not use these plates for multiple antibody assays, such as sandwich ELISAs, because the first antibody will not block all binding sites and subsequent antibodies can still bind to the plate, resulting in false positives.

Note: For binding information concerning the various proteins, please see Tech Tip #34: Binding characteristics of Protein A, G, A/G and L from our website.

Additional Materials Required

- Capture antibody specific for the antigen of interest
- Wash Buffer: TBS (Product No. 28376) or PBS (Product No. 28374) with added 0.05% Tween®-20 Detergent
- Dilution Buffer: SuperBlock® Blocking Buffer (Product No. 37535) containing 0.05% Tween-20 Detergent (Product No. 28320) or other proteinaceous blocking solution
- Antigen labeled with HRP, alkaline phosphatase or biotin
- Enzyme substrate (see our catalog or website for a complete product listing)

Procedure

1. Rinse each well three times with 200µL of Wash Buffer.
2. Using the Dilution Buffer dilute antibody to ~1µg/mL and add 100µL to each well.
3. Incubate plate 30-60 minutes at room temperature. For best results, for all microplate incubations use a plate mixer that creates a vortex in each well.
4. Rinse each well three times with 200µL of Wash Buffer.
5. Add the labeled antigen to each well (~0.1µg/mL). Sample may be diluted in Dilution Buffer. Incubate at 37°C for 1 hour.
6. Rinse each well three times with 200µL of Wash Buffer.
7. If using a biotinylated antigen, add enzyme-labeled streptavidin or other biotin-binding protein and incubate at 37°C for 1 hour. Rinse each well three times with 200µL of Wash Buffer.
8. Detect the signal according to the instructions for the specific detection system being used.

Related Thermo Scientific Products

| | |
|--------------|---|
| 34028 | 1-Step Ultra TMB-ELISA, 250mL |
| 34022 | 1-Step Turbo TMB, 250mL |
| 37621 | 1-Step PNPP, 100mL |
| 15075 | Reagent Reservoirs, 200/pkg |
| 15036 | Sealing Tape for 96-Well Plates, 100/pkg |

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.

© 2012 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.