

Pierce[®] Nickel Coated Plates

15142 15242 15342 15442

0690.6

Number	Description
15142	Pierce Nickel Coated Plates (clear, 8-well strip), 5 plates/package
15242	Pierce Nickel Coated Plates (white, 96-well) 5 plates/package
15342	Pierce Nickel Coated Plates (black, 96-well), 5 plates/package
15442	Pierce Nickel Coated Plates (clear, 96-well), 5 plates/package

Activation level: 200 μ L

Binding Capacity: ~9pmol His-tagged protein (27kDa)/well

Note: Plates are supplied pre-blocked with bovine serum albumin.**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 Nickel Coated Plates are ideal for analyzing polyhistidine-tagged fusion proteins by ELISA-based methods. Proteins that contain a succession of several histidine residues at the amino or carboxyl terminus have a strong binding affinity for metal. Bacterial lysates containing polyhistidine-tagged fusion proteins can be added directly to the plates without the need for blocking. The clear, white or black plates can be used with colorimetric, chemiluminescent or fluorescent detection methods, respectively.

Important Product Information

- His-tagged protein binding to metal may be related to topology, microenvironment, reversible reactions, pH, ionic strength, and other factors. Protein folding may hinder binding of the histidine tag. Binding can occur in chaotropic agents such as 8M urea, 6M guanidine•HCl, or 3M thiocyanate. Nickel binding requires the presence of at least two histidine residues.
- A variety of solutes are compatible with metal interactions with macromolecules, including nonionic detergents, ethylene glycol and dimethylsulfoxide. Sodium chloride may enhance or decrease affinity.
- Avoid using solutions that contain metal chelators such as EDTA. Also, avoid reducing agents, such as mercaptoethanol. High concentrations of imidazole are commonly used to elute nickel-bound his-tagged proteins and, therefore, should be avoided for binding reactions.

General ELISA Procedure

A. Materials

- Cell lysate containing polyhistidine-tagged fusion protein
- Dilution Buffer for lysate: Tris-buffered saline (TBS, Product No. 28376) or phosphate-buffered saline (PBS, Product No. 28374)
- Wash Buffer: Dilution Buffer containing 0.05% Tween[®]-20 Detergent (Product No. 28320)
- Antibody to poly-histidine-tagged fusion protein
- Enzyme-conjugated or fluorescent-labeled secondary antibody
- Enzyme Substrate (e.g., for HRP use the Thermo Scientific TMB Substrate Kit, Product No. 34021)

B. Method

Note: Plates are supplied pre-blocked with bovine serum albumin.

1. Dilute lysate with Dilution Buffer. Add 100µL of diluted lysate per well and incubate with shaking for 1 hour at room temperature.
2. Wash wells three times using 200µL of Wash Buffer for each wash.
3. Add 100µL of primary antibody per well and incubate for 1 hour at room temperature.
4. Wash wells three times using 200µL of Wash Buffer for each wash.
5. Add 100µL per well of secondary antibody. Incubate for 1 hour at room temperature.
6. Wash wells three times using 200µL of Wash Buffer for each wash.
7. Develop and evaluate plate.

Related Thermo Scientific Products

37070	SuperSignal® ELISA Pico Chemiluminescent Substrate, 100mL
15169	QuantaBlu™ Fluorogenic Peroxidase Substrate Kit
34028	1-Step™ Ultra TMB-ELISA, 250mL
37621	1-Step PNPP, 100mL
15075	Reagent Reservoirs, 200/pkg
15082	Microtube Racked System, 960 tubes
15036	Sealing Tape for 96-Well Plates, 100/pkg
78248	B-PER® Bacterial Protein Extraction Reagent, 500mL
78410	Halt™ Protease Inhibitor Cocktail, EDTA-Free, 1mL
89968	HisPur™ Cobalt Spin Columns, 1mL
15165	HisProbe™-HRP, 2mg
78100	B-PER 6xHis Fusion Protein Purification Kit

General References

- Belew, M., and Porath, J. (1990). Immobilized metal ion affinity chromatography. Effect of solute structure, ligand density and salt concentration on the retention of peptides. *J Chrom* **516**:333-354.
- Hemdan, E., *et al.* (1989). Surface topography of histidine residues: a facile probe by immobilized metal ion affinity chromatography. *Proc Natl Acad Sci USA* **86**:1811-5.
- Porath, J. (1992). Immobilized metal ion affinity chromatography. *Protein Expr and Pur* **3**:263-281.

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.