Thermo Scientific Nunc Immobilizer Nickel-Chelate

Instruction protocol

Thermo Scientific Nunc Immobilizer Nickel-Chelate

The Thermo Scientific™ Nunc™ Immobilizer™ Nickel-Chelate plates are manufactured using a patented photochemical method¹ for covalent coupling of ligands to polymer surfaces.

A Nickel-Chelate complex is coupled via a spacer to the plates. The Nunc Immobilizer Nickel-Chelate Plates are designed for optimal binding of fusion proteins that have been tagged with an amino sequence containing six repetitive histidine (His) molecules.

Materials:

- Nunc Immobilizer Nickel-Chelate Plates
- · His-tagged biomolecule

Reagents:

- · Coupling buffer: 0.01 M KCI
- Washing buffer: PBST (Phosphate Buffered Saline pH 7.3 containing 0.05% (v/v) TWEEN® 20)

Coupling protocol for 96 well plate:

- 1. Pre-wash the plate with PBST (3 x 300 ul/well)
- 2. Prepare a solution of a His-tagged protein (0.01-1 µg/ml) in 0.01M KCl
- 3. Add the protein solution to the wells of the Nunc Immobilizer Nickel-Chelate Plate (100 µl/well)
- 4. Incubate the plate with gentle agitation (100 rpm) at 20-25°C for 2 hours, or at +4°C overnight
- 5. Aspirate the wells and wash with PBST (3 x 300 µl/well)
- 6. The plate is now ready for your assay

Coupling protocol for 384 well microplate:

- 1. Pre-wash the plate with PBST (3 x 100 ul/well)
- 2. Prepare a solution of a 6 x His-tagged protein (0.01-1 µg/ml) in 0.01M KCl
- 3. Add the protein solution to the wells of the Nunc Immobilizer Nickel-Chelate Plate (50 µl/well)
- 4. Incubate the plate with gentle agitation (100 rpm) at 20-25°C for 2 hours, or at +4°C overnight
- 5. Aspirate the wells and wash with PBST (3 x 100 ul/well)
- 6. The plate is now ready for your assay

Additional comments and recommendations

It is recommended to include small amounts of a non-ionic detergent like TWEEN $20\ (0.05\%\ (v/v))$ in washing buffers and in assay buffers, as this generally improves the signal to noise ratio of the assay.

Stability

Nunc Immobilizer Nickel-Chelate Plates can be stored at room temperature to their expiration date, which appears on the case label.

Trademarks and Patents:

TWEEN 20 is a registered trademark of ICI American Inc., U.S.A. Immobilizer is a trademark of Exigon A/S, Vedbaek, Denmark.

The product is produced under license from Exiqon A/S - EP 08 20483 and foreign applications and patents.

References:

Jensen, S. P., Rasmussen, S. E., Jakobsen, M. H. Photochemical Coupling of Peptides to Polystyrene MicroWell Plates. Innovations & Perspectives in Solid Phase Synthesis & Combinatorial Chemical Libraries, (1996), 419-422.

thermoscientific.com

© 2014 Thermo Fisher Scientific Inc. All rights reserved. All trademarks are the property of Thermo Fisher Scientific Inc. and its subsidiaries.

ANZ: Australia: 1300 735 292, New Zealand: 0800 933 966; Asia: China Toll-free: 800-810-5118 or 400-650-5118; India: +91 22 6716 2200, India Toll-free: 1 800 22 8374; Japan: +81-3-5826-1616; Other Asian countries: 65 68729717 Europe: Austria: +43 1 801 40 0; Belgium: +32 2 482 30 30; Denmark: +45 4631 2000; France: +33 2 2803 2180; Germany: +49 6184 90 6000, Germany Toll-free: 0800 1-536 376; Italy: +39 02 95059 554; Netherlands: +31 76 571 4440; Nordic/Baltic countries: +358 9 329 10200; Russia/CIS: +7 (812) 703 42 15; Spain/Portugal: +34 93 223 09 18; Switzerland: +41 44 454 12 22; UK/Ireland: +44 870 609 9203

North America: USA/Canada +1 585 586 8800; USA Toll-free: 800 625 4327

South America: USA sales support: +1 585 899 7198 Countries not listed: +49 6184 90 6000 or +33 2 2803 2000

