

Thermo Scientific Nunc Immobilizer Streptavidin

Instruction protocol

Thermo Scientific Nunc Immobilizer Streptavidin

The Thermo Scientific™ Nunc™ Immobilizer™ Streptavidin Plates/Strips are manufactured using a patented photochemical method¹ for covalent coupling of ligands to polymer surfaces. Streptavidin, a high affinity biotin-binding protein isolated from *Streptomyces avidinii*, is covalently coupled via a spacer to the plastic plates/strips. Streptavidin has a molecular weight of approx. 60,000 Dalton and has an isoelectric point (pI) close to 5. The Nunc Immobilizer Streptavidin Plates/Strips are designed and optimized for detection of various types of biotinylated biomolecules like biotinylated oligo-nucleotides, peptides and proteins.

Materials:

- Nunc Immobilizer Streptavidin Plates/Strips
- Biotinylated molecule of choice

Reagents:

- 5 x SSCT, pH 7.0 (5 x SSC (750 mM NaCl, and 75 mM Sodium Citrate) containing 0.05% (v/v) TWEEN® 20)
- 2 x SSCT, pH 7.0 (2 x SSC (300 mM NaCl, and 30 mM Sodium Citrate) containing 0.05% (v/v) TWEEN 20)
- PBST, pH 7.2 (Phosphate Buffered Saline containing 0.05% (v/v) TWEEN 20)

Note: Buffers should be used within one week from preparation.

Recommended coupling concentrations:

- Biotinylated oligonucleotides: We recommend using the following concentration range: 0.01 – 0.5 µM diluted in 5 x SSCT buffer pH 7.0
- Biotinylated peptides: We recommend using the following concentration range: 1 ng/ml – 1 µg/ml diluted in a PBST buffer pH 7.2
- Biotinylated proteins: We recommend using the following concentration range: 0.05 µg/ml – 5 µg/ml diluted in a PBST buffer pH 7.2

Coupling protocol for 96 well plate, and 8 well strips:

1. Pre-wash your plate with 3 x 300 µl/well PBST or 5 x SSCT buffer without any incubation step. This is done to ensure improved readouts and a very low coefficient of variation (CV%)
2. Prepare a solution of your biotinylated molecule in PBST buffer or 5 x SSCT (oligonucleotides)
3. Add the solution to the wells of a Nunc Immobilizer Streptavidin plate/strip (100 µl/well)
4. Incubate the plate with gentle agitation for 1 hour at room temperature (20° - 25°C)
5. Aspirate the wells and wash with PBST or 2 x SSCT (oligonucleotides) 3 x 300 µl/well
6. Your surface is now ready for assay application

Coupling protocol for 384 well plate:

1. Pre-wash your plate with 3 x 100 µl/well PBST or 5 x SSCT buffer without any incubation step. This is done to ensure improved readouts and a very low coefficient of variation (CV%)
2. Prepare a solution of your biotinylated molecule in PBST buffer or 5 x SSCT buffer (oligonucleotides)
3. Add the solution to the wells of a Nunc Immobilizer Streptavidin plate (50 µl/well)
4. Incubate the plate with gentle agitation for one hour at room temperature (20° - 25°C)
5. Aspirate the wells and wash with PBST or 2 x SSCT (oligonucleotides) 3 x 100 µl/well
6. Your surface is now ready for assay application

Specifications:

- Streptavidin coated area ~100 µl/well (96 well format)
- Stable when stored at room temperature to the 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 Exiqon A/S, Vedbaek, Denmark.

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

References:

- ¹ Koch, T., Jacobsen, N., Fensholdt, J., Boas, U., Fenger, M. Jakobsen, M. H., Photochemical Immobilization of Anthraquinone Conjugated Oligonucleotides and PCR Amplicons on Solid Surfaces. *Bioconjugate Chem.* 11 (2000), 474-483.

thermoscientific.com

© 2014 Thermo Fisher Scientific Inc. All rights reserved. All (other) 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

MNLSPIIMMOBSTREP 0614

Thermo
SCIENTIFIC

A Thermo Fisher Scientific Brand