

# Detection of His-tagged fusion proteins

## Key Words

Thermo Scientific™ Nunc™ Immobilizer™ Ni-Chelate plates, coated plates, ELISA, His-tagged protein, fusion proteins, competitor analysis, low detection limit, high signal/noise ratio.

## Goal

The goal of this application note is to detect His-tagged fusion proteins in Nunc Immobilizer Ni-Chelate plates compared to competitor plates.

The Nunc Immobilizer Ni-Chelate plates are designed for optimal binding of His-tagged fusion proteins without the need for purification prior to application.

Detection limit is very low. The functional Ni-Chelate surface is based on patented technology for covalent coupling of ligands to polymer surfaces<sup>1</sup> via a spacer.

Detection of histidine (His)-tagged fusion proteins on the Nunc Immobilizer Ni-Chelate plate compared with competitor plates.

## Recommended protocol for the detection of His-tagged fusion proteins

### Reagents and materials

- Nunc Immobilizer Ni-Chelate plate
- Coupling buffer: 0.01 M KCl
- Washing buffer: PBST (Phosphate Buffered Saline (pH 7.2) containing 0.05% (v/v) Tween 20)
- Chromogen/substrate: 3,3', 5,5' tetramethylbenzidine (TMB)
- Stop solution: Sulphuric acid 0.5 M

### Assay procedure 96 well transparent plates

- The plates are pre-washed three times with 300 µL/well PBST.
- Purified or crude His-tagged fusion protein (100 µL/well) is added to the Immobilizer Ni-Chelate plates in concentrations from 0.01 to 1 µg/mL. A control without His-tagged fusion protein is also included.



- The plates are incubated for two hours at room temperature with gentle agitation (100 rpm).
- The plates are washed three times with 300 µL/well PBST.
- Antibody against the fusion protein conjugated to horseradish peroxidase (HRP) (100 µL/well) diluted e.g. 1:5000 in PBST is added to the plates.
- The plates are incubated for one hour at room temperature with gentle agitation (100 rpm).
- The plates are washed three times with 300 µL/well PBST.
- TMB (100 µL/well) is added to the plates.
- The plates are incubated for 10 minutes at room temperature in the dark.

- The enzyme reaction is stopped with 0.5 M sulfuric acid (100  $\mu$ L/well).
- The absorbance is read at 450 nm using an ELISA reader.

### Application example

A comparison of the performance of the Immobilizer Ni-Chelate plate and other commercially available plates was performed.

Three different preparations (A, B, C) of a fusion protein b-2-microglobulin (b-2-m 12 kDa) tagged with six His residues were tested.

- A crude preparation of b-2-m tagged with a sequential His tag (crude His-b-2-m).
- A purified preparation of b-2-m tagged with a sequential His tag (purified His-b-2-m).
- A purified preparation of b-2-m tagged with a non-sequential His tag (purified HAT-b-2-m). The protein was tagged with a sequence of 20 amino acid, which included 6 histidines uniformly distributed throughout the sequence.

The following assay was conducted on the Immobilizer Ni-Chelate plate:

- The plates were pre-washed three times with 300  $\mu$ L/well PBST.
- A His-tagged fusion protein (A, B or C) was applied to the Immobilizer Ni-Chelate plates (100  $\mu$ L/well) in ten concentrations of a 1:2 dilution (500 ng–0.98 ng/mL).
- The plates were incubated for two hours at room temperature with gentle agitation (100 rpm).
- The plates were washed three times with 300  $\mu$ L/well PBST.

- Antibody against the fusion protein conjugated with HRP (HRP conjugated to anti-b-2-m) diluted in PBST was added to the plates (100  $\mu$ L/well). The antibody was diluted to an appropriate concentration, here 1:5000.

- The plates were incubated for one hour at room temperature with gentle agitation (100 rpm).
- The plates were washed three times with 300  $\mu$ L/well PBST.
- TMB (100  $\mu$ L/well) was added to the plates.
- The plates were incubated for 10 minutes at +20-25°C in the dark.
- The enzyme reaction was stopped with 0.5 M sulfuric acid (100  $\mu$ L/well).
- The absorbance was read at 450 nm.

The assays on the competitor plates were performed using the protocol supplied by the different competitors.

Figs. 1 and 2 illustrate the OD response of a 1:2 dilution curve of the same fusion protein (crude His-b-2-m or purified His-b-2-m) applied to a Immobilizer Ni-Chelate plate, as well as to plates from competitors A, B and C.

As seen in Figs. 1 and 2 the performance of the Immobilizer Ni-Chelate plate is superior to the performance of the competitor plates regardless of the degree of purification of the His-tagged fusion protein.

Even when testing an unpurified preparation of a cell lysate using the Immobilizer Ni-Chelate plates, very small concentrations of His-tagged fusion protein (0.5 ng per well) can be detected assuming an OD cut off value of 0.5.

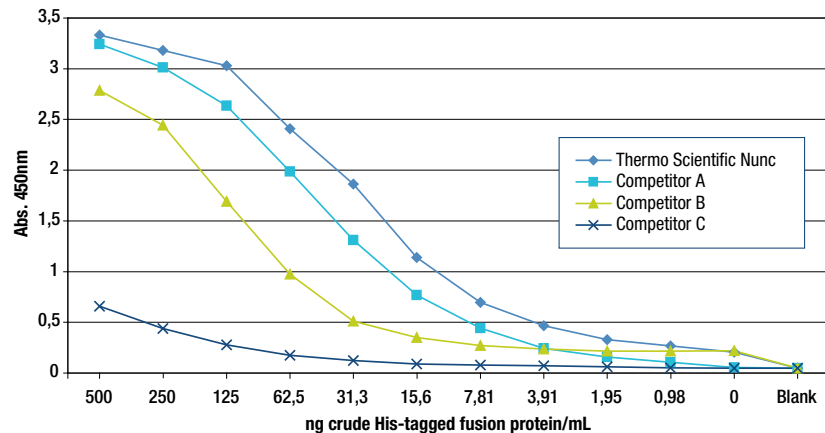


Fig. 1.

Detection of a His-tagged fusion protein from a crude cell lysate preparation (crude His- $\beta$ -2-m, preparation A). The His-tagged fusion protein was added in a 1:2 dilution to the plates. As a negative control, buffer without His-tagged fusion protein was added.

Fig. 3 illustrates a 1:2 dilution curve of the fusion protein tagged with a non-sequential His tag (purified HAT-b-2-m, preparation C). The purified HAT-b-2-m was applied to a Immobilizer Ni-Chelate plate, as well as to plates from competitors A, B and C. As seen in Fig. 3, the performance of the Immobilizer Ni-Chelate plate is superior to the performance of the competitor plates.

### Conclusion

As seen from Figs. 1 to 3, the Immobilizer Ni-Chelate plate shows a higher OD response when compared to the competitor plates. Also the calculated CV% is lower on the Immobilizer Ni-Chelate plate than on many competitor plates (data not shown).

This finding is true regardless of the histidine tagged fusion protein used (His in sequence or in a non-sequential distribution as well as a crude preparation of the His-tagged fusion protein).

Even when a fusion protein is tagged with a non-sequential His tag (HAT-b-2-m), the fusion protein binds equally well to the immobilizer Ni-Chelate plate. Using the competitor plates the HAT-b-2-m does not bind as well as the His-b-2-m.

### Acknowledgements

The His-tagged fusion proteins were kindly provided by Professor Søren Buus, Institute for Medical Microbiology and Immunology, Copenhagen University, Denmark.

### Credits

Assay and protocol for Ni-Chelate plate preparation was designed by Eva Jauho and Maria Nielsen.

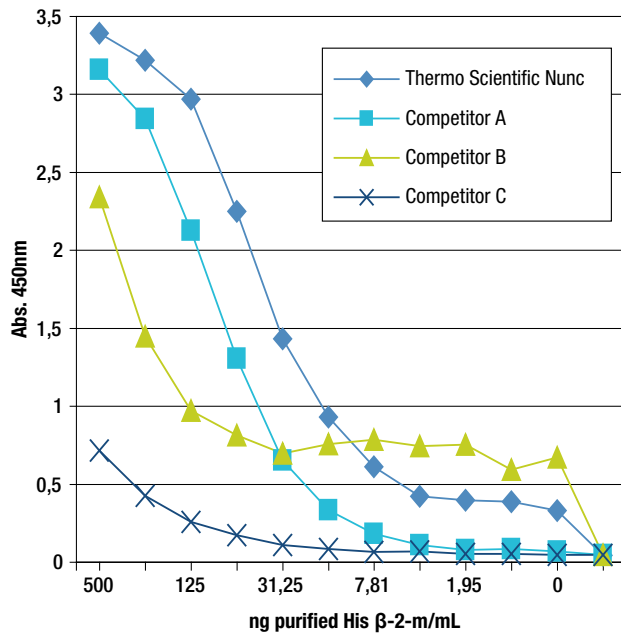


Fig. 2.

Detection of a purified His-tagged fusion protein (purified His-β-2-m, preparation B). The His-tagged fusion protein was applied in a 1:2 dilution to the plates. As a negative control, buffer without His-tagged fusion protein was applied.

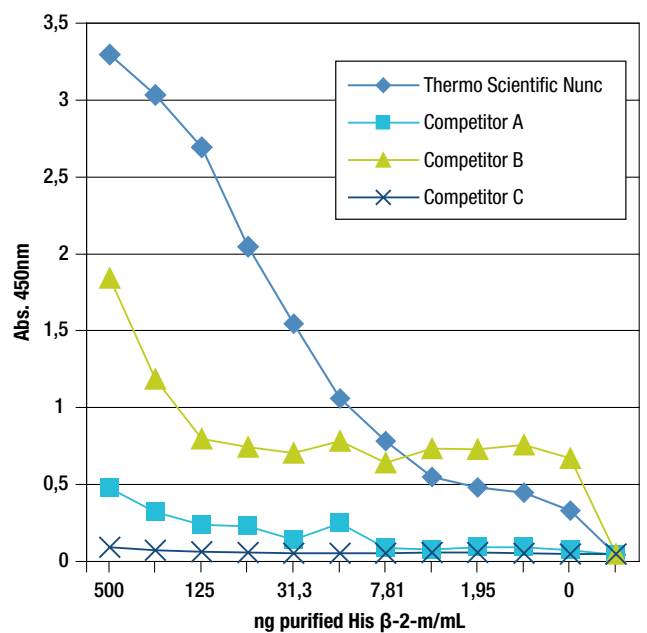


Fig. 3.

Detection of the fusion protein tagged with a non-sequential histidine tag (purified HAT-β-2-m, preparation C). The purified HAT-β-2-m was added in a 1:2 dilution to the plates. As a negative control, buffer without His-tagged fusion protein was added.

## References

- Jensen SP, Rasmussen SE, Jakobsen MH.  
Photochemical Coupling of Peptides to Polystyrene  
MicroWell Plates.  
Innovations & Perspectives in Solid Phase Synthesis &  
Combinatorial Chemical Libraries (1996), 419-422.
- Jacobsen MH, Koch T.  
Method of photochemical immobilization of ligands  
using quinones.  
WO 96/31557.

## References where covalent binding methods have been used

### Thermo Scientific™ CovaLink™

Krook M et al.  
Novel peptides binding to the Fc-portion of  
immunoglobulins obtained from a combinatorial phage  
display peptide library.  
J. Imm. Meth., 221:151-157.

Stadler S et al.  
Monoclonal anti-acid-labile subunit oligopeptide  
antibodies and their use in a two-site immunoassay for  
ALS measurement in humans.  
J. Imm. Meth. (2001), 252:73-82.

### Immobilizer

Davidson EJ et al.  
Human papillomavirus type 16 E2- and L1-specific  
serological and T-cell responses.  
J. Gen. Virol. (2003), 84:2089-2097.

Barend Bouma et al.  
Glycation Induces Formation of Amyloid Cross-Structure  
in Albumin.  
J. Biol. Chem. (2003), 278:41810-41819.

Nigel M. Page et al.  
Characterization of the endokinins: Human tachykinins  
with cardiovascular activity.  
PNAS (2003), 100:6245-6250.

Siao Ping Tsai et al.  
Nucleic acid capture assay, a new method for direct  
quantitation of nucleic acids.  
Nucleic Acids Research (2003), Vol. 31, No. 6, e25.

Irene Jauho E, Havsteen Jakobsen M.  
Rare cell isolation using antibodies covalently linked to  
slides: Application to fetal cells in maternal blood.  
Prenat. Diagn. (2003), 23 (11):898-900.

[thermoscientific.com/oemdiagnostics](http://thermoscientific.com/oemdiagnostics)

© 2014 Thermo Fisher Scientific Inc. All rights reserved. "Immobilizer" is a trademark of Exiqon A/S, Vedbaek, Denmark. The product is produced under license from Exiqon A/S - EP 0820483 and foreign applications and patents. "Tween" is a registered trademark of Uniqema Americas. 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

**Thermo**  
SCIENTIFIC

A Thermo Fisher Scientific Brand