# Custom-made Thermo Scientific Nunc Immobilizer for DNA Binding

Nana Jacobsen, M.Sc. and Jan Skouv, Ph.D.

#### **Key Words**

Thermo Scientific<sup>™</sup> Nunc<sup>™</sup> Immobilizer<sup>™</sup> DNA Surface, Thermo Scientific<sup>™</sup> Nunc<sup>™</sup> MicroWell<sup>™</sup> plate, high-dose hook effect, aminated PCR amplicon, covalent coupling, hybridization.

#### Goal

The goal of this application note is to show that the Immobilizer DNA plate is able to bind DNA using different concentrations of an aminated PCR amplicon. Futher to illustrate that the hybridization signal decreases to the background level if a too high concentration of amplicon is used in the assay.

> Nunc Immobilizer DNA MicroWell plates and strips for covalent immobilization of aminated DNA can be custom-made upon request.

> The production of the Nunc Immobilizer DNA surface introduces an ethylene glycol spacer and a stable electrophilic group that reacts with nucleophiles such as free amines. The spacer design and the density of electrophilic groups on this surface are optimized for detection of various types of molecules including aminated nucleic acids.



Fig. 1. Binding mechanism for aminated nucleic acids



#### **Coupling materials**

- Nunc Immobilizer DNA MicroWell plate
- 100 mm carbonate buffer, pH 9.6
- Aminated PCR amplicon
- SSC (1 x SSC is: 150 mm NaCl, 15 mm sodium citrate, pH 7.0)
- PBST is Phosphate Buffered Saline 1 (PBS) with 0.05% (v/v) Tween 20



#### **Coupling protocol**

- Prepare a solution of aminated DNA in 100 mm carbonate buffer, pH 9.6. It is recommended that the amount of aminated DNA is optimized, however for initial experiments we suggest: 1 nm aminated DNA PCR fragments (single stranded).
- Add the aminated DNA solution to the wells of the Immobilizer DNA plate (100 μL/well) (50 μL for 384).
- 3. Incubate the plate with gentle agitation at room temperature for two hours or overnight at +4°C.
- Aspirate the wells and wash with 3 x 300 μL 2 x SSC, 0.1% (v/v) Tween 20 (3 x 100 μL for 384).
- 5. The DNA surface is ready for use.

Detergents like Tween 20 effectively suppress covalent coupling of DNA and should consequently not be present in the coupling buffer. The use of competing nucleophiles like ethanolamine, lysine or tris (hydroxymethyl) amino methane (TRIS) should also be avoided in the coupling buffer.

The inclusion of small amounts of detergents like Tween 20 (0.05- 1% (v/v)) in subsequent wash and assay buffers, generally improves the signal to noise ratio of the assay. Other DNA concentrations, incubation times, temperatures, buffers or pH values than those recommended here can successfully be used.

#### **Application examples**

1. Detection of a PCR amplicon Using Nunc Immobilizer DNA for the detection of a PCR amplicon.

The DNA fragment to be detected was a 98 bp fragment from pUC19<sup>2,3</sup>. This 98 bp fragment was amplified using 5'- AAC AGC TAT GAC CAT

G-3' and 5'-GTA AAA CGA CGG CCA GT-3' as primers, pUC19 as template and a standard PCR kit. The fragment was amplified following the manufacturer's recommendations and incubating: 2 min. at 94°C; 30 cycles (94°C 1 min., 45°C 1 min., 72°C 2 min.); and 72°C 3 min. The yield was estimated by agarose gel electrophoresis.

#### **Protocol**

- Using the recommended coupling protocol, the capture sequence 5'-amine-AAC AGC TAT GAC CAT G-3' was covalently attached to the transparent Nunc Immobilizer DNA plate surface.
- Per well: 10 μL of the PCR reaction (approximately 40 ng) was dissolved in 2 x SSC, 0.1% (v/v) Tween 20. Boiled 5 min. and then placed on ice.
- 3. The specific detection probe (5'-biotin-ATG CCT GCA GGT CGA C-3') was added to the PCR/SSC mix. 0.5 pmol detection probe per μL, final vol. 100 μL. The mix was then added to the wells of the Nunc Immobilizer DNA plate and the PCR fragment was allowed to hybridise to the covalently attached capture sequence for 3 hours at 37°C (Fig. 2).
- The wells were aspirated and washed with 3 x 300 μL 2 x SSC, 0.1% (v/v) Tween 20.
- 5. A 1  $\mu$ g/mL solution of streptavidin/HRP in PBST was dispensed into the wells (100  $\mu$ L/ well), and the plate incubated for one hour.
- 6. The wells were aspirated and washed with 3 x 300  $\mu$ L PBST.
- 7. A solution of 6 mm orthophenylene- diamine (OPD), 4 mm  $H_2O_2$  in 100 mm citric acid buffer, pH 5.0 was added to the wells (100 µL/well) and left for color development.
- 8. After approximately 15 minutes, the enzyme reaction was stopped with  $H_2SO_4$ , 0.5 M (100 µL/ well) and the absorbance in this colorimetric assay was measured at 492 nm with an ELISA reader. The result are shown in Fig. 3. All incubations were carried out with gentle agitation at either room temperature or at 37°C when indicated.





Fig. 2. Detection of a PCR fragment by a target specific capture probe covalently linked to the DNA on an Nunc Immobilizer DNA plate using a biotinylated detection probe

Fig. 3. Detection of various amounts of the 98 bp PCR fragment from pUC19

#### Special application using PCR products

With the advent of polymerase chain reaction (PCR), ligase chain reaction (LCR 4), and similar techniques, double-stranded (ds) DNA fragments with a well defined DNA sequence can be prepared. In particular the efficient generations of dsDNA fragments by PCR have found numerous applications in diverse fields of biomedicine and molecular biology.

Oligos with an amino group attached to its 5'-end can be purchased from most commercial oligo suppliers. Including one such primer in a PCR (or LCR) reaction result in the synthesis of aminolabeled dsPCR fragment. Such fragments can easily be covalently linked to the surface of the Nunc Immobilizer DNA plates and strips and used for various applications.

### **Enzymatic activity**

A number of important enzymes, for instance restriction enzymes, kinases, phosphatases, polymerases, methylases etc. act on DNA. The possible relation between enzymatic activity and specific DNA sequences can conveniently be tested on DNA's covalently linked to the Immobilizer DNA surface. Analysis of DNA binding proteins We suggest that dsDNA's with various recognition sequences are generated and covalently attached to the Nunc Immobilizer DNA surface.

#### **Gene discovery**

A number of gene discovery methodologies (e.g. differential display 5) result in a large number of PCR fragments that have to be screened for the presence of a given consensus sequence. We suggest to attach such PCR fragments generated with one aminolabeled oligo to the surface of the Nunc Immobilizer DNA and screen for the presence of a particular DNA sequence as described below.

## Immobilization and detection of an aminolabeled PCR amplicon on the surface of Nunc Immobilizer DNA plate.

To illustrate various aspects of the performance of the Nunc Immobilizer DNA plates for detection of an amino PCR amplicon. The DNA fragment to be detected was a 630 bp fragment from human Nras 6,7,8. This 630 bp fragment was amplified using 5'-NH2-C6- spacer-CCA GCT CTC AGT AGT TTA GTA CA-3' (position 1427- 1449) and 5'-AAG TCA CAG ACG TAT CTC AGA C-3' (position 2035-2056) as primers, human Nras as template and a standard PCR kit. All oligos were purified by HPLC. The fragment was amplified following the manufacturer's recommendations and incubating 3 min. at 95°C; 30 cycles (55°C 2 min., 72°C 3 min., 95°C 1 min.); 55°C 2 min. and 72°C 3 min. The yield was estimated on a standard 1% agarose gel stained with ethidium bromide.

#### 2. Detection of amino PCR amplicon

- 1. Using the recommended coupling protocol described above for a 96 well plate, the NH2- Nras amplicon was covalently attached to the Nunc Immobilizer DNA plate. 10  $\mu$ L of the PCR reaction (approximately 120 ng) was diluted in 1:2 dilutions in 100 mm carbonate buffer, pH 9.6, and 100  $\mu$ L was dispensed per well. Incubation for two hours was allowed.
- After coupling the amplicon was denatured by 200 μL 0.4 M NaOH 0.25% (v/v) Tween 20 for 5 minutes and washed with 3 x 300 μL 2 x SSC, 0.1% (v/v) Tween 20. The Nras DNA was detected by hybridization with the specific detection probe 5'-TGT GTT TGT GCT GTG GAA GAA CCCbiotin- 3' (position 1549-1572). The probe was diluted in 2 x SSC, 0.1% (v/v) Tween 20 final concentration. 100 μL 0.5 μm probe was added per well of the Immobilizer DNA plate. The detection probe was allowed to hybridise to the covalently attached sequence for two hours at 37°C.
- The wells were aspirated and washed with 3 x 300 μL 2 x SSC, 0.1% (v/v) Tween 20.
- 4. A solution of streptavidin/ HRP in PBST (1  $\mu$ g/mL) was dispensed into the wells (100  $\mu$ L/ well), and the plate is incubated for one hour. The wells were aspirated and washed with 3 x 300  $\mu$ L PBST.
- 5. A solution of ortho-phenylenediamine (OPD), 6 mm and  $H_2O_2$ , 4 mm in 100 mm citric acid buffer, pH 5.0 was added to the wells (100 µL/well) and left for color development.

After approximately five minutes, the enzyme reaction was stopped with  $H_2SO_4$ , 0.5 M (100  $\mu$ L/well), and the absorbance was measured at 492 nm using an ELISA reader.

All incubations are carried out with gentle agitation at either room temperature or at 37°C when indicated.

# pplication Note

#### **Results**

The result of a typical experiment is shown in Fig. 4. The experiment indicates that at high concentrations of the amino amplicon the hybridization signal decreases to the background level. This effect appears somewhat similar to the 'high-dose hook effect' described for various immuno-assays <sup>9-11</sup> and emphasizes that an optimization of amino amplicon concentrations is necessary.



Fig. 4. Detection of various amounts of the 630 bp PCR fragments from Nras

### References

- Sambrook J, Fritsch EF, Maniatis T.Molecular Cloning: A Laboratory Manual. Cold Spring Harbour Laboratory Pres., NY (1989), 2'nd ed.
- 2. Yanisch-Perron C, Vieira J, Messing J. Improved M13 phage cloning vectors and host strains: Nucleotide sequences of the M13mp18 and pUC19 vectors. Gene (1985) 33:103-119.
- 3. pUC 19 accession no.: VB0026.
- 4. Wiedmann M, Wilson WJ, Czajka J, Luo J, Barany F, Batt CA. Ligase chain reaction (LCR) - overview and applications. PCR Methods Appl. (1994), 3:51-64.
- Liang P and Pardee AB. Differential display of eukaryotic messenger RNA by means of the polymerase chain reaction. Science (1992), 257:967-71.
- 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.
- Brown R and Hall I. Human N-ras: cDNA cloning and gene structure. Nucleic Acid Research (1985), 13:5255-5268.
- 8. Nras Accession no.: X02751.
- Rodbard D, Feldman Y, Jaffe ML, Miles LE. Kinetics of two-site immunoradiometric ('sandwich') assays— II: Studies on the nature of the-high-dose hook effect. Immunochemistry (1978), 15:77-82.
- Wolf BA, Garrett NC, Nahm MH. The "hook effect": high concentrations of prostate-specific antigen giving artifactually low values on one-step immunoassay. N Engl J Med. (1989), 320(26):1755-6.
- 11. Fernando SA, Wilson GS. Multiple epitope interaction in the twostep sandwich immunoassay. J Immunol Methods (1992), 151:67-86.

#### thermoscientific.com/oemdiagnostics

© 2015 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. Specifications, terms and pricing are subject to change. Not all products are available in all countries. Please consult your local sales representative for details.

Asia: Australia: 1300-735-292; New Zealand: 0800-933-966; China +86-21-6865-4588 or +86-10-8419-3588; China Toll-free: 800-810-5118 or 400-650-5118; Singapore +65-6872-9718; Japan: +81-3-5826-1616; Korea +82-2-2023-0640; Taiwan +886-2-87516655; India: +91-22-6680-3000 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/CIS countries: +358-10-329-2200; Russia: +7-(812)-703-42-15; Spain/Portugal: +34-93-223-09-18; Switzerland: +41-44-454-12-12; 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: +1-585-586-8800 Countries not listed: +49-6184-90-6000 or +33-2-2803-2000



TILSPNUNCTN56 0915