

# Thermo Scientific Nunc NucleoLink Procedure for Solid Phase PCR (DIAPOPS)

## Key Words

Thermo Scientific™ Nunc™ NucleoLink™ strips, covalent attachment, heat stable surface, binding of oligonucleotides, solid phase PCR, detection, hybridization.

## Goal

The goal of this application note is to give a detailed protocol for solid phase PCR followed by detection and hybridization in the same well (DIAPOPS).

The Nunc NucleoLink surface binds DNA covalently and heat stably.

The Nunc NucleoLink strips are V-shaped with flat bottoms and are compatible with thermal cyclers and plate readers. Nunc NucleoLink can therefore be used for solid phase PCR followed by detection by hybridization in the same well (DIAPOPS).

The strips are available in transparent colorimetric detection of DNA. This application note gives a detailed protocol for DIAPOPS.

## Preparation for and covalent immobilization of solid phase primer

1. The solid phase primer should be aminated or phosphorylated at the 5' end, and a linker of at least 10 T's (thymidine) should be added between the active primer sequence and the 5' end group.
2. Prepare a freshly made coating mix consisting of 100 nm solid phase primer and 10 mm EDC (1-ethyl-3-(3-dimethylaminopropyl)-carbodiimide) in 10 mm 1-methyl-imidazole (1-MeIm) (pH 7.0).
3. Add 100  $\mu$ L of this coating mix to each Nunc NucleoLink well. This gives a total of 10 pmol of the 5'-modified solid phase primer per well.
4. Seal the Nunc NucleoLink Strips (e.g. with Nunc Sealing Tape).
5. Incubate the Nunc NucleoLink Strips at 50°C for 4-24 hours.
6. Wash the empty Nunc NucleoLink wells three times,



soak for 5 minutes and wash three times, all with 100 mm Tris-HCl (pH 7.5), 150 mm NaCl and 0.1% Tween 20 at room temperature (RT). Empty the strips.

7. Wash once with deionized sterile water.
8. The empty, unsealed coated Nunc NucleoLink Strips can be stored at 4°C or below in a polyethylene bag.

## Amplification

1. BSA must be added to the PCR mix in a final concentration of 1 mg/mL which should not be exceeded as this may block the solid phase amplification. 0.1%-0.25% Tween 20 is also recommended.
2. The primers must be added to the PCR mix in a ratio of 1:8. We use 25 pmol per reaction of the liquid phase primer and 25/8 pmol per reaction of the primer used as the solid phase primer.
3. Add PCR mix to the wells (normally 20  $\mu$ L or 45  $\mu$ L).

4. Add DNA template to each well (the total reaction volume has been tested with both 25  $\mu$ L and 50  $\mu$ L).
5. Seal the Nunc NucleoLink Strips with Sealing Tape 8.
6. Place the strips in a thermal cycler block.
7. Place the silicone spacer plate on the tape sealed Nunc NucleoLink Strips.
8. Temperature cycle the strips with the temperatures and cycling parameters specific for the system.
9. Remove the Nunc NucleoLink Strips and empty them.
10. Wash the empty Nunc NucleoLink wells three times, soak for 5 minutes and wash three times, all with freshly made 0.2 M NaOH and 0.1% Tween 20 at RT.
11. Wash the empty Nunc NucleoLink wells three times, soak for 5 minutes and wash three times, all with 100 mm Tris-HCl (pH 7.5), 150 mm NaCl, and 0.1% Tween 20 at RT.

### Detection

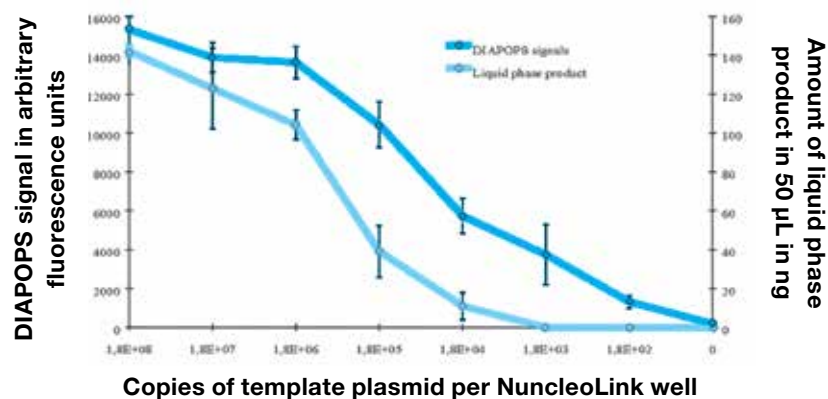
1. Add 100  $\mu$ L of 50-100 nm biotinylated hybridization probe diluted in 5 x SSC, 0.1% Tween 20 and 0.5% blocking reagent (BR) to each well.
2. Seal the Nunc NucleoLink Strips with sealing tape and incubate at 45-50°C for one to 20 hours.
3. Wash the empty Nunc NucleoLink wells three times at RT with 0.5 x SSC and 0.1% Tween 20.
4. Soak for 15 minutes at 50°C with 0.5 x SSC and 0.1% Tween 20.
5. Wash three times at RT with 0.5 x SSC and 0.1% Tween 20.
6. Detection of biotin label on the hybridized probe:  
When using alkaline phosphatase (AP): Add to each well 100  $\mu$ L AP conjugated streptavidin diluted 1:3000 (or as the producer suggests) in 100 mm Tris-HCl (pH 7.5), 150 mm NaCl, 0.1% Tween 20 and 0.5% BR.  
When using Horse Radish Peroxidase (HRP): Add to

each well 100  $\mu$ L HRP conjugated streptavidin diluted 1:5000 (or as the producer suggests) in 100 mm Tris-HCl (pH 7.5), 150 mm NaCl, 0.1% Tween 20 and 0.5% BR.

7. Seal the wells with sealing tape and incubate for one hour at 37°-50°C.
8. Wash the empty Nunc NucleoLink wells three times, soak for 5 minutes, and wash three times with 100 mm Tris-HCl (pH 7.5), 150 mm NaCl and 0.1% Tween 20 at RT.
9. Substrate incubation:  
When using 4-methylumbelliferyl phosphate: Add 100  $\mu$ L of 1 mm 4-MUP dissolved in 1 M diethanolamine (pH 9.8) and 1 mm MgCl<sub>2</sub> to each well. Incubate at 37°-50°C in the dark for 30-60 minutes. Add 50  $\mu$ L of 3 M K<sub>2</sub>HPO<sub>4</sub> to stop the hydrolyzation of 4-MUP.  
When using p-nitrophenyl phosphate: Add 100  $\mu$ L of 1 or 10 mg/mL pNPP in 1 M diethanolamine (pH 9.8) and 1 mm MgCl<sub>2</sub> to each well. Incubate at RT for 30 minutes (10 mg/mL) to 24 hours (1 mg/mL). Add 100  $\mu$ L of 1 M NaOH to stop hydrolyzation of pNPP.  
When using TMB: Add 100  $\mu$ L of the ready-to-use solution to each well. Incubate for 30 minutes at RT. Add 100  $\mu$ L 0.1 M H<sub>2</sub>SO<sub>4</sub> to stop the reaction. (Re-hybridization is not possible after addition of acid).
10. To measure the signal:  
Detection of hydrolyzed 4-MUP: Determine the signal in a fluorescence plate reader: Excitation 360 nm, emission 450 nm.  
Detection of hydrolyzed pNPP: Measure OD in a standard plate reader at 405 nm.  
Detection of TMB: Measure OD in a standard plate reader at 450 nm after the reaction has been stopped with acid (if the reaction is not stopped with acid, the color is blue and can be measured at 655 nm).

Fig. 1

**DIAPOPS signal and liquid phase PCR products comparison, n=3**

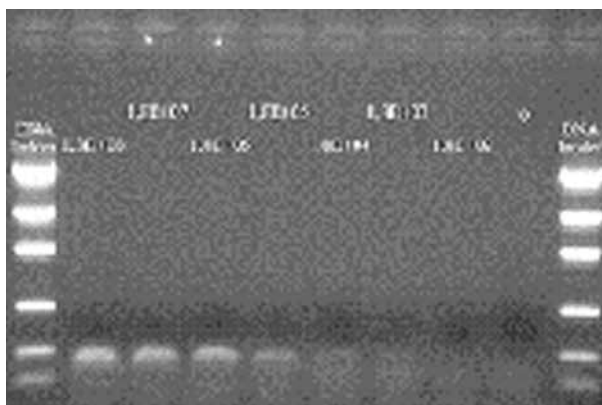


## Preparation for re-hybridization or storage

1. Remove substrate from wells after measuring the catalyzed substrate.
2. Wash the empty Nunc NucleoLink wells three times, soak for 5 minutes and wash three times with freshly made 0.2 M NaOH and 0.1% Tween 20 at RT.
3. Wash the empty Nunc NucleoLink wells three times, soak for 5 minutes, and wash three times with distilled water or 100 mM Tris-HCl (pH 7.5), 150 mM NaCl and 0.1% Tween 20 at RT.
4. The washed empty Nunc NucleoLink Strips can be stored at 4°C or below in an ethylene bag.

## Re-hybridization

1. No re-hydration of the Nunc NucleoLink Strips is necessary after storage. Commence with step 1 in the detection section and add hybridization solution directly to the empty dry wells.



Comparison of DIAPOPS and detection by gel electrophoresis. Using DIAPOPS the detection limit (1800 copies) was at least 1:10 improved compared to the gel (18000 copies).

## References

1. Rasmussen et al. *Biotechnology International* 1999, Vol. 2, 303-307.
2. Lundin BS, Karlsson MR, L. A. Svensson LA, Hanson LÅ, Dahlgren UIH, Telemo E. Active suppression in orally tolerized rats coincides with in situ transforming growth factor-beta (TGF-) expression in the draining lymph nodes. *Clinical & Experimental Immunology* 1999, Vol. 116(1), 181-187.
3. Niessen L, Klusmann J and Vogel RF. Quantitative estimation of *Fusarium graminearum* DNA using a solid phase PCR assay (DIAPOPS). *Journal of Food Mycology* 1998, 1:73-84.
4. Oroskar, Rasmussen SE, Rasmussen HN, Rasmussen SR, Sullivan BM and Johansson A. Detection of immobilized amplicons by ELISA-like techniques. *Clin Chem* 1996, 42:1547-1555.
5. Rasmussen SR, Rasmussen HB, Larsen MR, Hoff-Jorgensen R and Cano RJ. Combined polymerase chain reaction-hybridization microplate assay used to detect bovine leukemia virus and *Salmonella*. *Clin. Chem.* 1994, 40:200-205.

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

© 2014 Thermo Fisher Scientific Inc. All rights reserved. "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