

# Performance of Thermo Scientific Nunc Immuno C8 StarWell

The Nunc Immuno Module C8 StarWell, MaxiSorp or PolySorp surfaces, consists of Nunc MicroWells containing eight inner fins. As a result, the surface area is increased by a factor of approximately 1.5 when compared to standard, flat-bottom F-wells using reactant volumes ranging from 50 to 200  $\mu\text{L}$ .

In a solid phase assay, e.g. ELISA, a larger surface/volume ratio implies a faster adsorption of molecules from the liquid phase. Thus, the primary advantage of a larger surface/volume ratio is that assay time can be reduced.

Nunc StarWell incubation times may be reduced by a factor of approximately 2.25. According to adsorption kinetics modelling <sup>1</sup>, incubation times can be reduced by a factor equal to the square of the surface/volume increase factor, without reducing adsorption.

This Tech Note demonstrates Nunc StarWell vs. standard F-well performances for first layer, passive IgG adsorption according to the model <sup>1</sup>. The performance relationship holds for immobilization of any reactant in a solid phase bioassay sequence <sup>1</sup>.



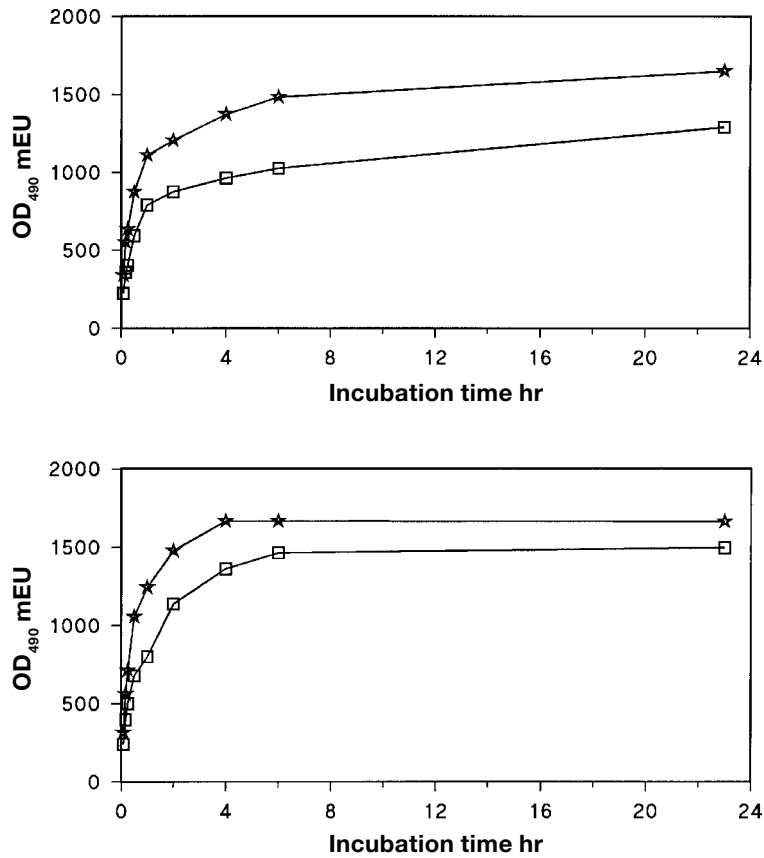


Fig. 1.

Data of IgG adsorption kinetics with Nunc MaxiSorp StarWell (★) and standard well (□) using IgG:HRP conjugate (Dako P 128) and IgG (Dako Z 181) diluted in carbonate buffer, pH 9.6.

In one case, 150  $\mu$ L/well of IgG:HRP conjugate was used in a dilution corresponding to 1% surface saturation of a standard well, i.e.  $F_{\square} = 0.01$  (left). In another case, 150  $\mu$ L/well of IgG + IgG:HRP (= 99+1) was used in a dilution corresponding to 100% standard well saturation, i.e.  $F_{\square} = 1$  (right).  $F$  generally denotes the ratio between the number of supplied molecules and the number of molecules that can be adsorbed. The MaxiSorp IgG adsorption capacity has been estimated to 650 ng/cm<sup>2</sup>, and at 150  $\mu$ L liquid volumes the Nunc StarWell and standard well surface/volume ratios are 1.5 x 8.3 and 8.3 cm<sup>-1</sup> respectively<sup>3</sup>.

After adsorption, the wells were washed 3 times with PBS + 0.2 M extra NaCl + 0.05% Triton X-100, pH 7.2. This was followed by substrate reaction with 150  $\mu$ L H<sub>2</sub>O<sub>2</sub>/OPD in phosphate-citrate buffer, pH 5.0, and addition of 110  $\mu$ L 2N H<sub>2</sub>SO<sub>4</sub> stopping solution, implying 260  $\mu$ L reading volumes.

Fig. 2.

Relationship between OD increase and liquid volume with Nunc StarWell due to the fins' elevation of liquid heights.

This relationship is empirically found to be approximately linear in a semilogarithmic plot. The dashed regression line represents the relationship with Nunc MaxiSorp, and the dotted line the relationship with Nunc PolySorp. The solid line and the equation represent the average data (○). In the hydrophobic Nunc PolySorp wells, the liquid stands higher than in the hydrophilic MaxiSorp wells due to the downward liquid surface curvature in Nunc MaxiSorp. This gives rise to a higher Nunc MaxiSorp elevation percentage. Actual percentages are determined by multiplying or dividing the average calculation by 1.15 for Nunc MaxiSorp or Nunc PolySorp respectively. In the present case, using Nunc MaxiSorp and reading volumes of 260  $\mu\text{L}$ , the elevation is approximately 7%. The Nunc StarWell readings must therefore be reduced by 7% to match the standard well readings for a comparative analysis of the respective adsorption performances (Figs. 3-5).

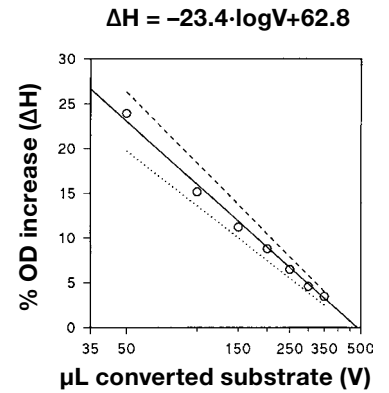


Fig. 3.

Simulations of the Fig. 1 data, adjusted according to Fig. 2, by the adsorption kinetics model <sup>1</sup>, showing correlation between real and theoretical kinetics.

At molecular supplies corresponding to  $F_{\square} = 1$  (Fig. 3A), the Nunc StarWell adsorption acceleration is much larger than at molecular supplies corresponding to  $F_{\square} = 0.01$  (Fig. 3B). The reason for this is explained in Fig. 4. As an example, the levels of 75% adsorption of supplied molecules (red horizontal lines) are used for comparison of the Nunc StarWell vs. standard well performances.

Between the lines of equal adsorption percentage (75%) and equal incubation time (red vertical lines), the StarWell curves exhibit regimes, elaborated in Fig. 5, where both time reduction and adsorption increase can be obtained.

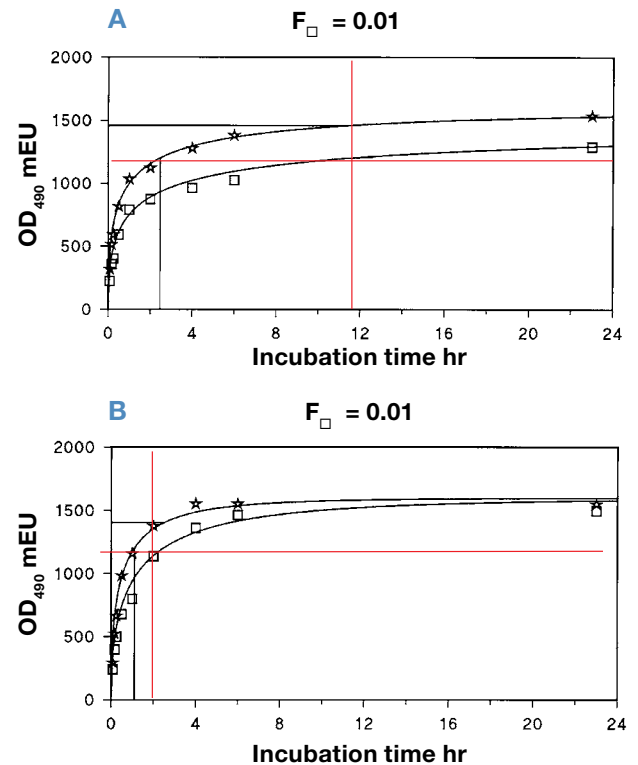


Fig. 4.

Model derived incubation times,  $T_{75}$ , for 75% adsorption of supplied molecules, as functions of  $F$ , in percent of the standard well  $T_{75}$  value at  $F = 1$ , with Nunc StarWell (lower curve) and standard well (upper curve).

The curves are almost constant for  $F$  values up to 0.1, because in this  $F$  regime, plenty of unoccupied sites are available on the surface. At  $F = 1$ , the curves assume maximum values, because the surfaces approximate saturation during adsorption; therefore, it will take a longer time for the last molecules to find the last unoccupied sites. In the present cases, where the molecular supplies correspond to  $F_{\square} = 0.01$  and 1, the corresponding Nunc StarWell values are 1.5 times smaller due to the 1.5 times larger surface area, i.e.  $F_{*} = 0.0067$  and 0.67, respectively. Therefore, a relatively larger Nunc StarWell adsorption acceleration is obtained at the high  $F$  values (dashed lines) than at the low  $F$  values (dotted lines) – in accordance with the data in Fig. 3. Thus, the claimed incubation time reduction factor of 2.25 is valid for  $F_{\square}$  values up to approximately 0.1. For  $F_{\square}$  values ranging from 0.1 to 1, the factor will be larger.

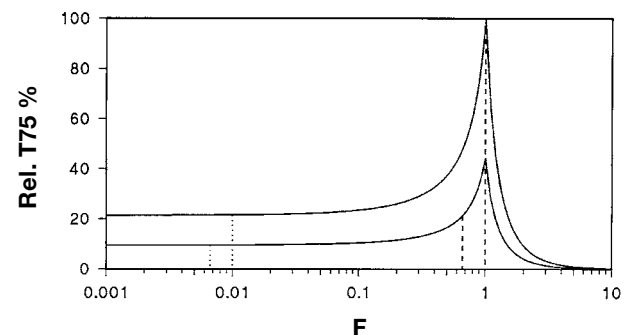
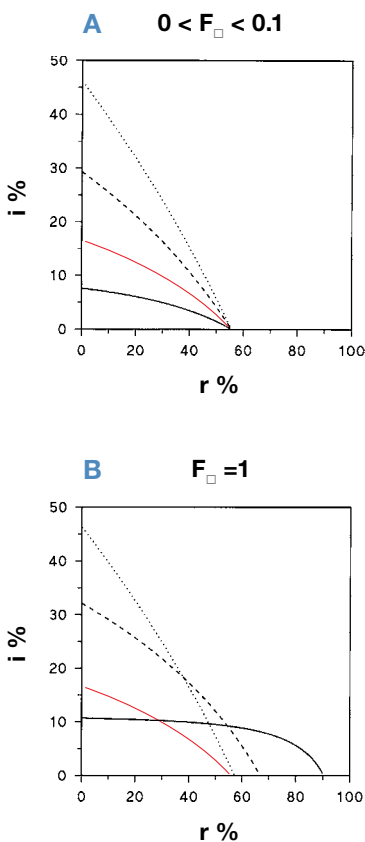


Fig. 5.

Model derived adsorption increase ( $i$ ) vs. incubation time reduction ( $r$ ) with Nunc StarWell compared to standard well for 10% (···), 50% (---), 75% (—) and 90% (—) adsorption of supplied molecules.

For  $F_{\square}$  values up to 0.1 (Fig. 5A) the maximum time reduction is approximately constant for any adsorption level. At  $F_{\square} = 1$  (Fig. 5B) the maximum time reduction increases with increasing adsorption level (Fig. 3). For  $F_{\square}$  values ranging from 0.1 to 1, a gradual transition from the left to the right diagram curves will occur.

Between the extremes of maximum  $r$  ( $i = 0$ ) and maximum  $i$  ( $r = 0$ ), intermediate  $r$  and  $i$  values can be obtained with StarWell.



## References

- Esser P. (1992).  
The surface/volume ratio in solid phase assays.  
Thermo Scientific Nunc Bulletin No. 10.
- Esser P. (1988).  
Principles in adsorption to polystyrene.  
Thermo Scientific Nunc Bulletin No. 6.
- Esser P. (1985).  
Adsorption geometry in Thermo Scientific Nunc  
Immuno products.  
Thermo Scientific Nunc Bulletin No. 1.

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

© 2014 Thermo Fisher Scientific Inc. All rights reserved. "Dako" is a registered trademark of Dako Denmark A/S. 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