

C8 Black Thermo Scientific Nunc LockWell FluoroNunc MaxiSorp and PolySorp for Fluorescence Detection

The new black Thermo Scientific Nunc LockWell format with MaxiSorp and PolySorp wells have a recommended maximum volume of 350 μL . These breakable strips have letters and notches on each well for easy identification of individual wells. The strips are suitable for all commonly used automated equipment. The dense pigmentation of black Nunc™ LockWell™ modules minimizes background fluorescence. MaxiSorp™ is optimized for binding IgG (antibodies) and PolySorp™ for binding more hydrophobic molecules.

In this study we compare the performance of the black LockWell modules with similar products from other large suppliers: Competitors A, B and B low volume (total volume of 205 μL /well). Plate uniformity and binding capability are determined by monitoring the reaction of immobilized horseradish peroxidase (HRP) from coating with a mixture of rabbit IgG and HRP conjugated rabbit anti-mouse IgG. Three of each plate type were tested on three different days.

When comparing high binding surfaces, the highest binding capability and lowest relative standard deviation was found to be LockWell MaxiSorp and Competitor A (using coating volumes of 150 μL /well) (Fig. 1).

Comparing the LockWell format to Competitor B low volume (using a coating volume of 100 μL /well) a very low relative standard deviation and high binding capability was found for the LockWell format. The binding capability of the LockWell MaxiSorp surface can easily be increased using a higher coating volume.

Assay

- Coating overnight at room temperature with antibody mixture (100 μL /well using low volume or 150 μL /well using standard 96 format).
- Wash three times with washing buffer.
- Addition of HPPA substrate (100 μL /well using low volume or 150 μL /well using standard 96 format).
- Addition of 50 μL /well sodium hydroxide after 14 minutes. The fluorescence was measured on EnVision 2101 using optimized fluorescence protocols with filter sets 340/405 and 485/535 nm.

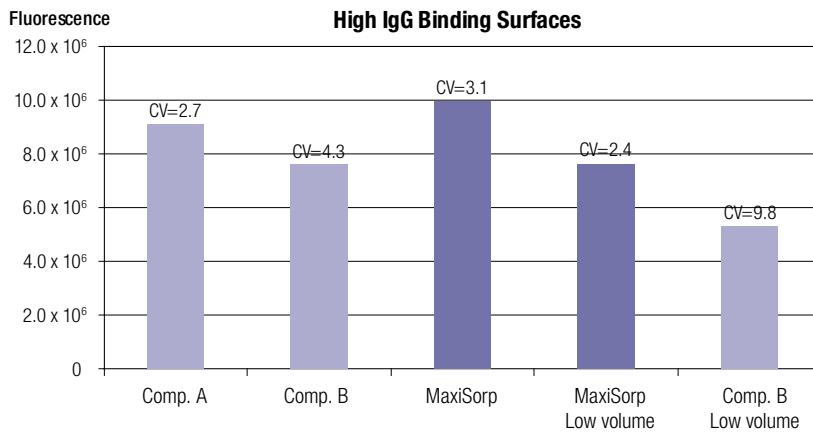


Fig. 1.

Fluorescence intensity (antibody binding capability) measured after performing antibody binding assay on different medium binding surfaces. Coating volume of Competitor A, B and MaxiSorp modules was 150 $\mu\text{L}/\text{well}$, and coating volume of PolySorp and competitor B low volume was 100 $\mu\text{L}/\text{well}$. Mean CV for tested module plates is shown above the respective binding capability column.

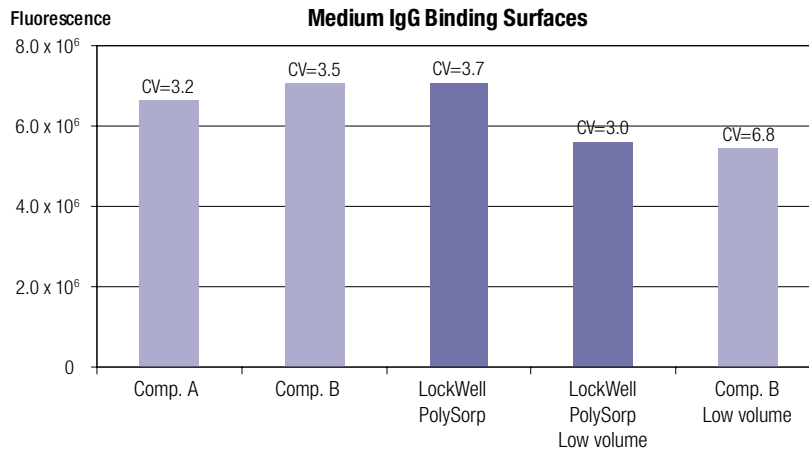


Fig. 2.

Fluorescence intensity (antibody binding capability) measured after performing antibody binding assay on different high medium surfaces. Coating volume of Competitor A, B and PolySorp modules was 150 $\mu\text{L}/\text{well}$, and coating volume of PolySorp and competitor B low volume was 100 $\mu\text{L}/\text{well}$. Mean CV for tested module plates is shown above each relevant binding capability column.

Reagents

Antibody mixture consisting of 65 ng/mL HRP conjugated rabbit anti-mouse IgG and 10 $\mu\text{g}/\text{mL}$ rabbit IgG, diluted in 0.05 M sodium carbonate buffer, pH 9.6.

Washing buffer: 0.15 M PBS, pH 7.2, with 0.05% detergent (Triton X for high binding surfaces and Tween 20 for medium binding surfaces).

Freshly prepared HPPA substrate: 2.5 mm 3-(p-hydroxyphenyl) propionic acid dissolved in 0.1 M TRIS buffer, pH 8.5, and 1.5 μL 30% hydrogen peroxide is added to 100 mL substrate.

Binding capability and relative standard deviations between medium binding surfaces was found to be comparable, using coating volumes of 150 $\mu\text{L}/\text{well}$ (Fig. 2). Comparing the LockWell format to Competitor B low volume, a very low relative standard deviation is

achieved using a coating volume at 100 $\mu\text{L}/\text{well}$. The binding capability when using the LockWell for low volume coatings is comparable to the Competitor B low volume. The binding capability of the LockWell PolySorp surface can easily be increased using a higher coating volume.

Fluorescence background measurements were performed using excitation and emission filter sets 485/535 and 340/405 nm. The data are not shown, as it was found relatively low for all products, regardless of the surface type and format.

Conclusion

Data show high binding capability and high uniformity using fluorescence detection on black C8 LockWell FluoroNunc™, demonstrated for different coating volumes by IgG binding assay on both the MaxiSorp and PolySorp surfaces.

thermoscientific.com/oemdiagnostics

© 2014 Thermo Fisher Scientific Inc. All rights reserved. "Triton" is a registered trademark of Dow Chemical Company; and "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