

Principles in Adsorption to Polystyrene

Peter Esser, M.Sc., Senior Scientist, Thermo Fisher Scientific

When considering the binding capacity of adsorbant plastic surfaces for biomacro-molecules, one must distinguish between the total amount of molecules that can be bound to the surface and the amount that can be bound and still remain biologically active. Both quantities are very much dependent on the nature of the molecules and the character of the surface.

Adsorption Forces

The adsorption of molecules to a polystyrene surface is due to inter-molecular attraction forces (van der Waals forces), to be distinguished from »true« chemical bonds, i.e. covalent bonds (through electron share) and ionic bonds (through stoichiometric charges of opposite signs) (see Fig. 1). Intermolecular attraction forces are based on intramolecular electric polarities of which two types can be distinguished: alternating polarities (AP) and stationary polarities (SP), i.e. dipoles.

AP arises when molecules approach each other, thereby creating disturbances in each other's electron clouds. This causes synchronously alternating polarities in the molecules, which may establish a bond between them, as illustrated in Fig. 2.

AP mediated binding is a common substance property, which is obviously the stronger, the larger the molecules implied. This is demonstrated by the fact that melting and boiling points increase with number of carbon atoms in the non-polar hydrocarbon series.

Indeed, it is due to this force that non-polar molecules at all aggregate into liquids and solids.

In addition to the AP attraction forces, molecules may possess SP (stationary polarity) through which they can bind to each other simply by bedding dipole against dipole, as illustrated in Fig. 1a.

Compared with SP, AP attraction decreases drastically with increasing distance between the molecules. Thus, AP attraction is inversely proportional to the seventh power of the distance, former has a much shorter range than the latter.

In general, van der Waals mediated bonds are about 100 times weaker than ionic and covalent bonds. However, among SP mediated bonds the hydrogen bond takes up an exceptional position because it is up to 10 times stronger than the others and because of its crucial importance for the properties of water and for the specific behaviors of bio-molecules.

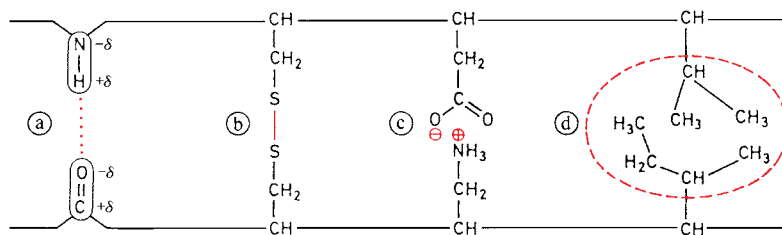


Fig. 1.

The four main types of possible bonds between macromolecules. »True« chemical bonds are represented by a covalent disulphide bond (b) and an ionic bond between a carboxyl ion and an amino ion (c). Van der Waals mediated bonds are represented by a hydrogen bond between two dipoles (a) and an alternating polarity bond between hydrocarbon residues protruding from the macromolecules' backbones (d), where the encircled area indicates a water-deprived zone. See text for further explanation.

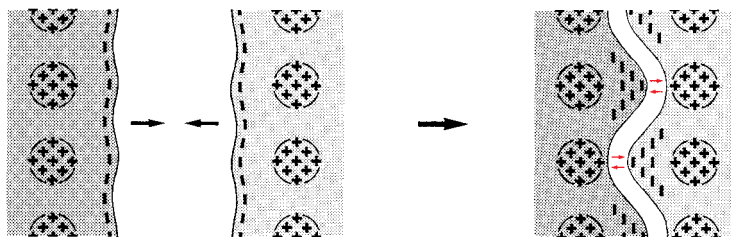


Fig. 2.

Schematic illustration of how synchronously alternating polarities (AP), created by reciprocal electron cloud disturbances in approaching molecules, can establish a bond between the molecules. Transient, minus-charged electron cloud condensations in one molecule will attract reciprocally exposed, plus charged nuclear regions in the other molecule.

Chemical groups, which can take part in hydrogen bonding, in particular $-OH$, $=O$, $-NH_2$, $=NH$, $\equiv N$, are called hydrophilic, as opposed to hydrophobic groups lacking this ability. Accordingly, hydrogen bonds may be called hydrophilic bonds, as opposed to AP mediated bonds, which are called hydrophobic bonds. The AP mediated attraction is also called hydrophobic interaction.

Adsorbing Surfaces

The Thermo Scientific Nunc Immuno program comprises different types of adsorbent polystyrene surfaces, including the Thermo Scientific Nunc PolySorp and MaxiSorp surfaces. While Nunc™ PolySorp™ predominantly presents hydrophobic groups, MaxiSorp™ has in addition many hydrophilic groups, which results in a fine patchwork of hydrophobic and hydrophilic binding sites. In aqueous medium, a repelling effect exists between the PolySorp surface and hydrophilic macromolecules (i.e. rich in hydrophilic groups), because these molecules will rather tend to intermingle with the water molecules (i.e. be dissolved) by the strong hydrogen bonds than bind to the surface by the weak hydrophobic bonds.

On a MaxiSorp surface, however, adsorption of hydrophilic macromolecules will be greatly facilitated, because not only can this surface compete with the water molecules for binding the macromolecules by hydrogen bonds, but the molecules can also be captured from a much longer distance by the long-range hydrogen bond

forces for establishment of both hydrogen bonds and eventually hydrophobic bonds (see Fig. 3).

On the other hand, hydrophobic macro-molecules (i.e. deficient in or lacking hydrophilic groups) can only be loosely adsorbed to MaxiSorp, because this surface tends to bind water molecules by hydrogen bonds, against which the macromolecules cannot compete and therefore exhibit poor ability for displacing water molecules and hydrophobic adsorption free from water pocket interruptions.

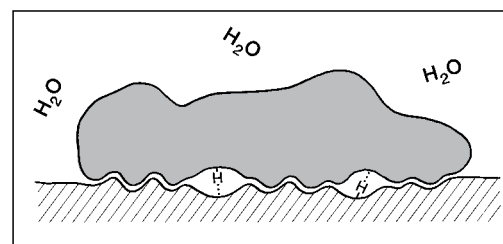


Fig. 3

Schematic illustration of how a hydrophilic macromolecule can be firmly adsorbed to MaxiSorp by »squeezing« out the water between the molecule and the surface through the combined action of hydrogen bond and AP bound forces. See text for further explanation.

Table 1

Theoretical PolySorp and MaxiSorp preferences for adsorption of various bio-macromolecules.

← PolySorp	MaxiSorp →
Proteins & Peptides*	
Lipoproteins	Glycoproteins
Compound lipids	Compound polyglycans
Lipids	Polyglycans

* Surface preference is dependent on predominance of hydrophobic or hydrophilic amino acid residues in the molecules.

On a PolySorp surface, however, no hindrance exists for stable hydrophobic adsorption of hydrophobic molecules, except that they may not be applicable in purely aqueous medium, wherefore addition of or substitution with detergent or organic solvents (e.g. ethanol or hexane) may be needed.

In summary, when no attention is paid to maintenance of specific activities, hydrophobic compounds bind preferably to PolySorp, and hydrophilic compounds preferably to MaxiSorp (see Table 1). However, with MaxiSorp binding events are more likely to occur, which means that adequate incubation conditions are more easy to establish – a fact that may extend the MaxiSorp application range well into the theoretical PolySorp domain. Nevertheless, considering the maintenance of the specific activities of the molecules (e.g. enzymatic, immunologic), which of course is crucial, the specific sites may well be obscured, impaired or destroyed through the binding to the surface.

Therefore, when constructing a solid phase assay, it is generally recommended to try adsorption to MaxiSorp first. If this does not work satisfactorily, it may be due to molecular malfunction, and one should then try PolySorp whereby specific activities may be maintained because of the different binding mechanism to this surface.

However with PolySorp, where molecules must come very close to the surface to establish hydrophobic bonds, one must anticipate more demanding incubation conditions, such as higher reactant concentration, longer duration, higher temperature, (more) agitation, to obtain adsorption efficiency comparable with MaxiSorp.

As mentioned above, van der Waals mediated bonds are relatively weak, wherefore they may be insufficient for stable binding when they are few in number, i.e. when the molecules are small. For binding of small molecules strong chemical bonds are needed. Ionic bonds would not do, because they normally dissociate in aqueous solution, leaving covalent bonds as the only possibility for direct, stable binding of small molecules. However, this difficulty may be overcome by using small molecules linked to (indifferent) carrier macromolecules. Small molecules

would in this context be e.g. peptides of less than 10 amino acids (corresponding to about 1500 dalton).

Geometric Estimation

Before making any experimental estimates of binding capacities on solid phase surfaces, it is worth making an estimate from geometric considerations of how many molecules can maximally be packed in one layer on a surface.

Taking immunoglobulin G antibody (IgG) as an example, and assuming that it is globular and packed in the densest monolayer packing (Fig. 4), the amount QGLOBE per cm^2 will be:

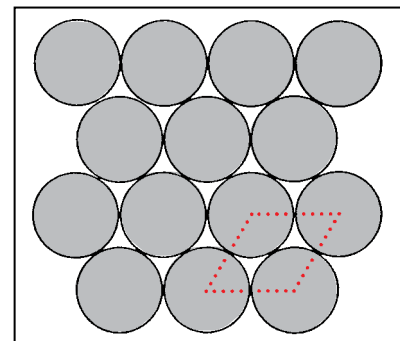


Fig. 4

The densest monolayer packing of globular molecules seen from above. The factor $2/\sqrt{3}$ in the text formulas for surface binding capacities originates in this non-quadratic pattern.

However, according to various sources, the IgG molecule is rather a lens shaped spheroid with a diameter, d , of about 15 nm and a thickness, t , of about 3 nm, as illustrated in Fig. 5.

Assuming the densest packing of these spheroids in »upright« or »lying« position (Fig. 6), the respective Q_{LENSE} values will accordingly be:

$$Q_{\text{LENSE}} = \frac{2}{\sqrt{3}} \cdot \frac{\text{MW}}{N} \cdot 10^9 \cdot \begin{cases} 1/dt = 650 \text{ ng/cm}^2 & \text{at upright position} \\ 1/d^2 = 130 \text{ ng/cm}^2 & \text{at lying position} \end{cases}$$

So, for geometrical reasons alone, the maximum amount of monolayer IgG that can be bound on a surface is 650 ng/cm². If an average is taken between the two Q_{LENSE} figures, the final estimate would be 400 ng/cm².

Assuming that molecular weight is proportional with volume, Q will not change considerably within wide molecular weight limits, other things being equal, because of the low power relationship between volume and profile area of a body. Fig. 7 illustrates the relationship between Q and molecular weight for globular molecules.

Experimental Estimation

Let us, as an example, stay with IgG, a glycoprotein with a structure shown schematically in Fig. 8.

On a MaxiSorp surface, one would expect an orientated adsorption in favor of exposing the antigen recognizing sites, because this surface would favor a binding through the hydrophilic carbohydrate moiety associated with the non-recognizing leg of the molecule.

On a PolySorp surface, on the other hand, one would expect an adsorption in favor of obscuring the antigen-recognizing sites, because of the repelling effect between this surface and the carbohydrate moiety.

To investigate the actual adsorption conditions, the following experiment was designed (see Fig. 9).

PolySorp and MaxiSorp MicroWell surfaces were coated with a dilution series of specific antibodies, starting with a concentration C well above saturation concentration, or with a corresponding dilution series mixed with unspecific antibodies to a constant total of C . The relative amount of specific antibody adsorbed in each case was determined by a sandwich ELISA for the antigen in question using excess antigen and excess HRP conjugated specific antibodies.

When it is assumed that equal signals mean equal amounts of specific antibody adsorbed in both dilution series, the quantity B/C is the fraction of the saturation concentration S formed by the maximum signal concentration A , i.e. $A = S \cdot B/C$, or:

$$S = \frac{A}{B} \cdot C$$

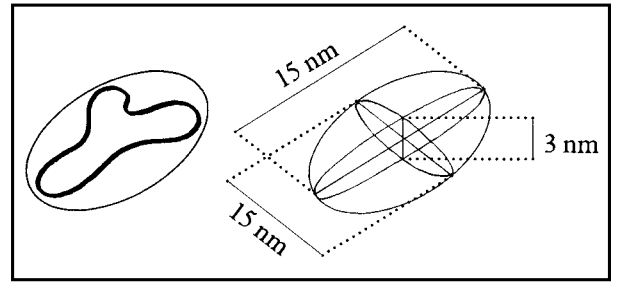


Fig. 5

The Y-shaped IgG antibody will approximately take up the volume of a lens shaped spheroid with a diameter of 15 nm and a thickness of 3 nm.

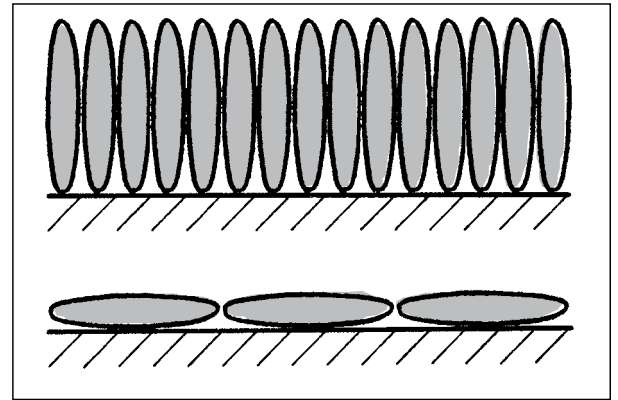


Fig. 6

Profiles of the densest IgG packings on a surface illustrating the density ratio of 5 to 1 between molecules packed in upright position (above) and in lying position (below).

$$Q_{\text{GLOBE}} = \frac{2}{\sqrt{3}} \cdot \frac{\text{MW}}{N} \cdot \frac{1}{(2r)^2} \cdot 10^9 \text{ ng/cm}^2 = 300 \text{ ng/cm}^2$$

where:

MW = molecular weight of IgG = 153,000 g v mole⁻¹

N = Avogadro's number = 6 · 10²³ mole⁻¹

r = Stokes radius of IgG = $\frac{R \cdot T_{20}}{6 \cdot \pi \cdot \eta_{20} \cdot D_{20} \cdot N}$ cm

R = gas constant = 8.3 · 10⁷ g · cm² · sec⁻² · °K⁻¹ · mole⁻¹

T₂₀ = room temperature (20°C) = 293°K

η_{20} = viscosity of water at 20°C = 1 · 10⁻² g · cm⁻¹ · sec⁻¹

D₂₀ = diff. coeff. of IgG ref. to water at 20°C = 4 · 10⁻⁷ cm² · sec⁻¹

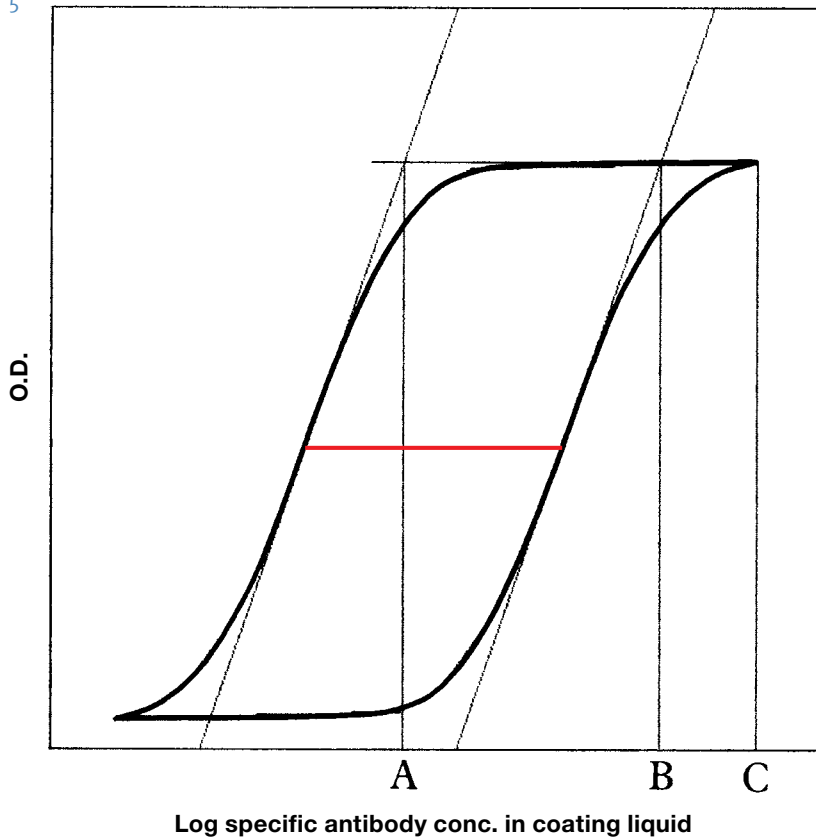


Fig. 9

Expected results from ELISA experiments with a dilution series of first layer specific IgG-antibodies, starting with a concentration C well above surface saturation concentration S (left sigmoid curve), or with a corresponding dilution series mixed with unspecific IgG to a constant total of C (right sigmoid curve). The ratio between S and C is A/B, which is represented by the distance between the two curves (red line). See text for further explanation.

In Fig. 10 are shown the results from experiments, designed as above, with four antibodies of different specificities. From the seemingly constant curve distances for MaxiSorp (MS) and PolySorp (PS) respectively, it is concluded that the adsorptions are independent of antibody specificity, and that they amount to the following quantities:

$$Q_{MS} = \frac{A_{MS}}{B_{MS}} \cdot C \cdot \frac{V}{F} \cdot 10^3 = 650 \text{ ng/cm}^2$$

$$Q_{PS} = \frac{A_{PS}}{B_{PS}} \cdot C \cdot \frac{V}{F} \cdot 10^3 = 220 \text{ ng/cm}^2$$

where:

$$A_{MS}/B_{MS} = 1/20$$

$$A_{PS}/B_{PS} = 1/60$$

$$C = \text{max. IgG conc.} = 100 \text{ } \mu\text{g/mL}$$

$$V = \text{reactant volume} = 0.2 \text{ mL}$$

$$F = \text{surface area} = 1.54 \text{ cm}^2$$

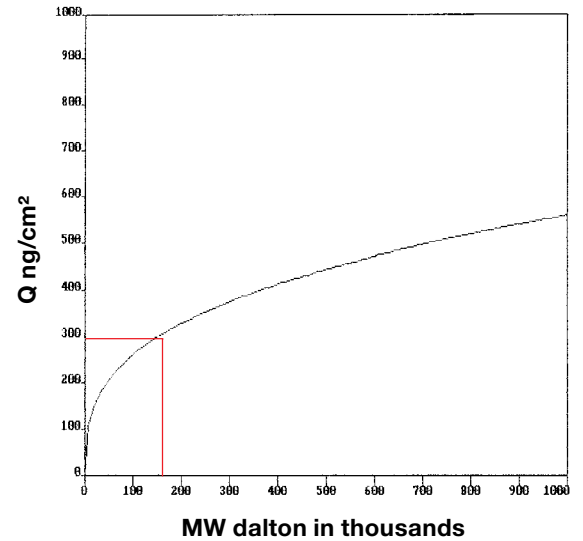


Fig. 7

Relationship between monolayer weight density (Q) and molecular weight (MW) of globular molecules illustrating that within a 10-factor MW range Q will roughly vary within only a 2-factor range. The curve is extrapolated on the basis of an idealized IgG molecule with an MW of 153,000 (red lines). See text for further explanation.

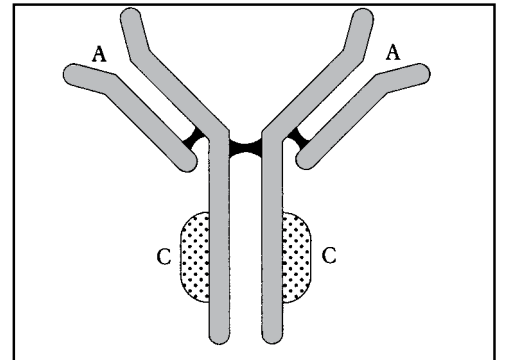


Fig. 8

Schematic illustration of the IgG antibody structure. Note the carbohydrate moiety (at C) associated with the leg opposite the antigen binding sites (at A) of the molecule.

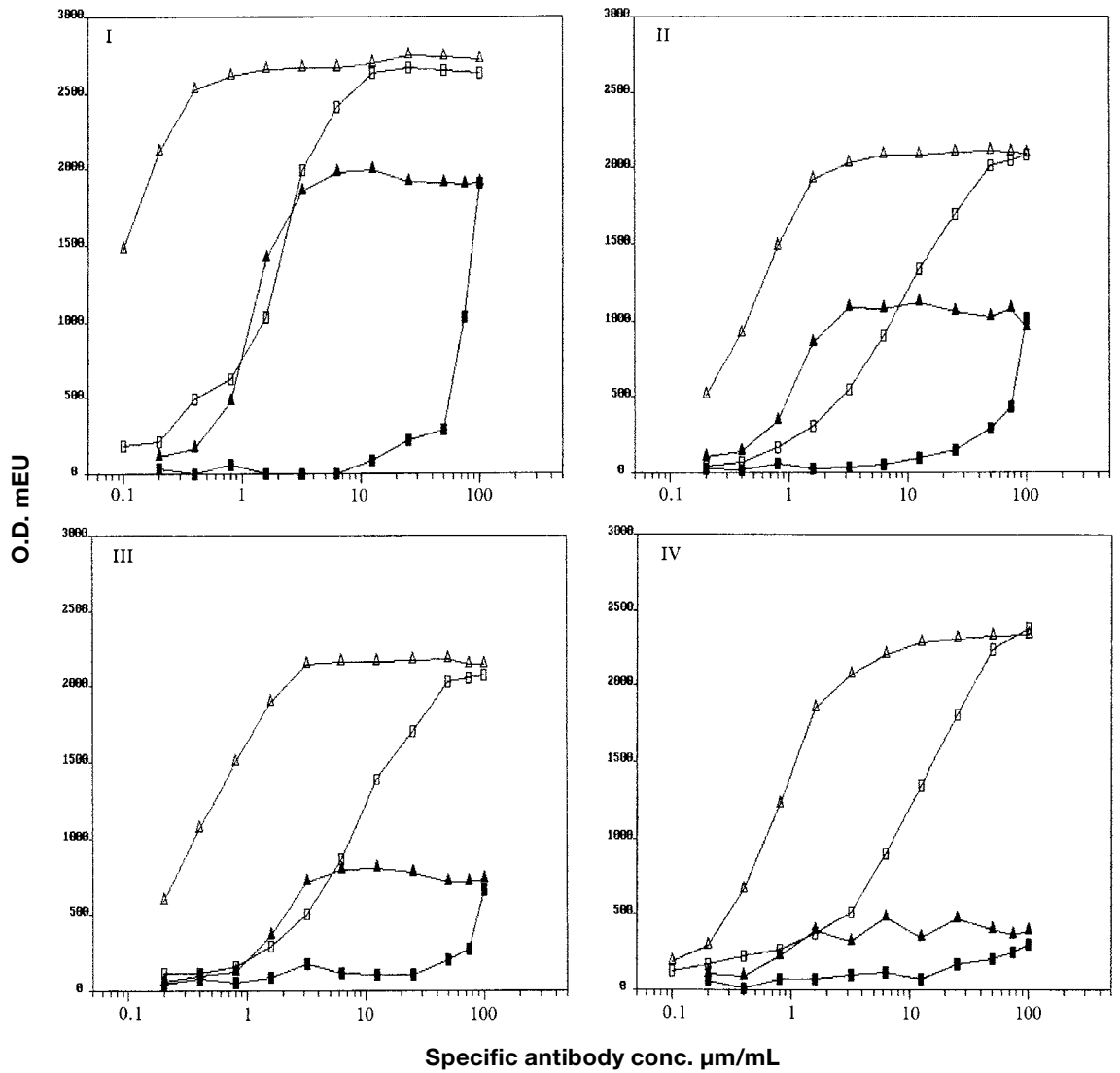


Fig. 10

Results from experiments prospected in Fig. 9 with four different antibody/antigen systems on MaxiSorp (open symbols) and on PolySorp (filled symbols).

I: ferritin antigen (MW 440,000);

II: fibronectin antigen (MW 450,000);

III: thyroglobulin antigen (MW 670,000);

IV: AFP (a – foetoprotein) antigen (MW 70,000).

Note the seemingly constant curve distances for MaxiSorp and PolySorp respectively, regardless of the system in question.

See text and Fig. 9 for further explanation.

Discussion

Whereas QMS is identical with the geometric maximum estimate for upright molecules, QPS is only one third hereof, which can be explained by assuming that on PolySorp upright and lying molecules are present in equal numbers, as illustrated in Fig. 11.

Because each IgG antibody can maximally bind two antigen molecules, this PolySorp decrease in number of adsorbed antibodies would have the greater effect, the smaller the antigen molecules are compared with the antibodies, as illustrated in Fig. 12.

This could, partly at least, explain the very low PolySorp signals for AFP, which has a molecular weight of less than half the weight of IgG, whereas the other antigens have 3-5 times the weight of IgG.

In addition, the variation of PolySorp-to-MaxiSorp maximum signal ratios may be due to differently obstructed affinities through the antibody adsorption to PolySorp, and/or due to different antibody-antigen affinities from one system to the other. In the ferritin system, the affinity seems relatively high since there is a long delay before maximum signal decline on MaxiSorp, which may be consistent with the high PolySorp maximum signal, i.e. the higher the affinity, the less it is obstructed by antibody adsorption to PolySorp.

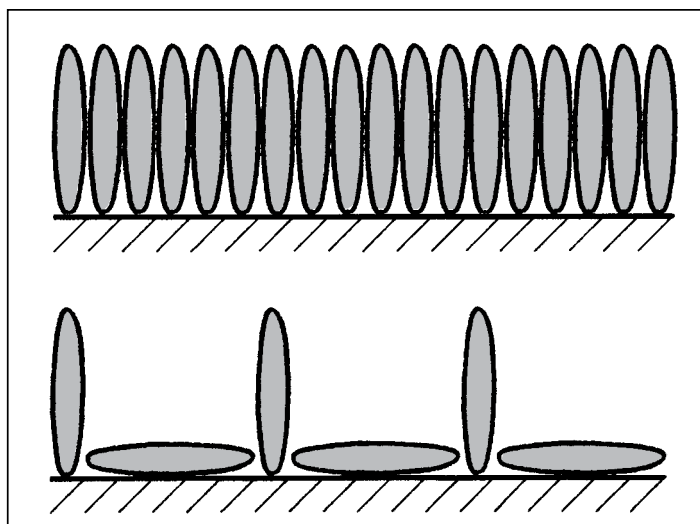


Fig. 11

Profiles of IgG adsorption patterns on MaxiSorp (above) and PolySorp (below) which can explain the experimentally found ratio of 3 to 1 between the densities on the respective surfaces.

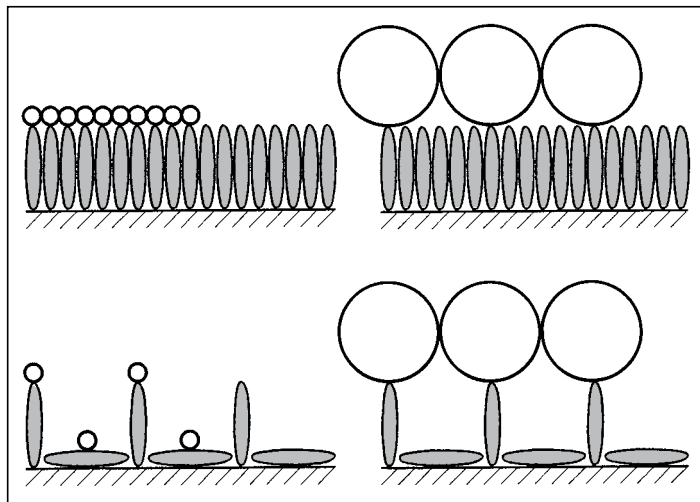


Fig. 12

Profiles of second layer (antigen) binding to antibody-coated MaxiSorp (above) and PolySorp (below) surfaces illustrating how the presumptive difference between the IgG adsorption patterns may imply a PolySorp decrease in bound amounts of small antigen molecules (left), but not of large antigen molecules (right). It should be noted that the third layer consisting of HRP conjugated antibodies would hardly influence the detection of this phenomenon, as HRP is a relatively small molecule (MW 40,000). See text for further explanation.

thermoscientific.com/oemdiagnostics

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