

Antibody Immobilization

Relative Affinity of immobilized Protein A for various species and subclasses of polyclonal and monoclonal IgGs^[1]

Relative avidity of Protein A for polyclonal antibodies

| Species / Subclass (Polyclonal) | Protein A |
|---------------------------------|-----------|
| Rabbit | ++++ |
| Cow | ++ |
| Horse | ++ |
| Goat | - |
| Guinea pig | ++++ |
| Sheep | +/- |
| Pig | +++ |
| Rat | +/- |
| Mouse | ++ |
| Chicken | - |
| Human IgG | ++++ |
| Human IgM | - |
| Human IgD | - |
| Human IgA | - |

++++ (strong binding) – (weak or no binding)

Relative affinity of Protein A for various isotypes

| Species / Subclass (Monoclonal) | Protein A |
|---------------------------------|-----------|
| Human | |
| IgG ₁ | ++++ |
| IgG ₂ | ++++ |
| IgG ₃ | - |
| IgG ₄ | ++++ |
| Mouse | |
| IgG ₁ | + |
| IgG _{2a} | ++++ |
| IgG _{2b} | +++ |
| IgG ₃ | ++ |
| Rat | |
| IgG ₁ | - |
| IgG _{2a} | - |
| IgG _{2b} | - |
| IgG _{2c} | + |

++++ (strong binding) – (weak or no binding)

Reference:

[1] Harlow and Lane (eds.) (1988) *Antibodies. A Laboratory Manual*. Cold Spring Harbor Laboratory, N.Y., 617-618.