

Immunology Tools to Enhance Your Research

BioIVT understands that the right reagent is required to properly sustain your cells. We offer reagents designed for your research needs.

Immunoglobulin Inhibiting Reagent

Product Numbers:

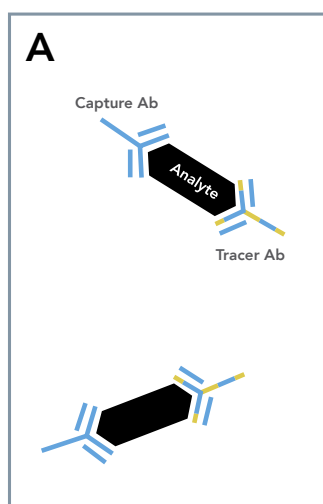
IIRLIQUID
IIRPOWDER

Immunoglobulin Inhibiting Reagent (IIR) is a formulation of immunoglobulins that specifically target against heterophilic antibodies to block their interference in immunoassays. IIR has an extremely high binding affinity to inhibit the action of human heterophilic antibody.

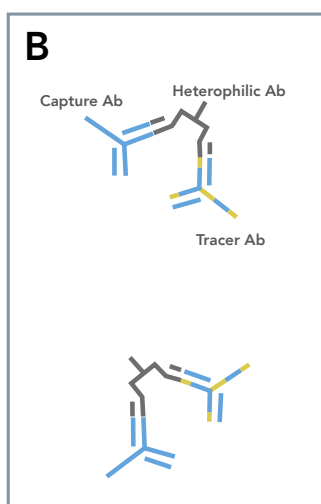
- Ready-to-use solution
- Widely cited (see website for detail)
- Available in 10mg, 1g, and custom sizes



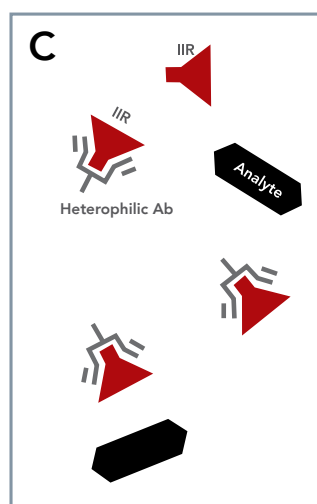
Binding of IIR



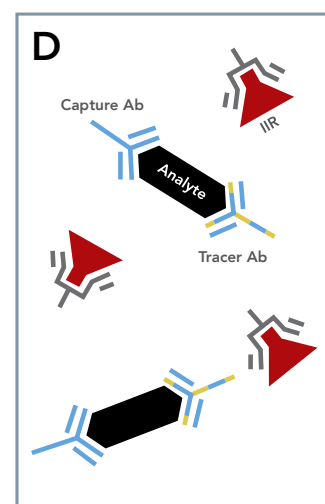
Normal Sandwich immunoassay where analyte concentration creates positive results



Presence of heterophilic antibody creates false positive results



Added immunoglobulin inhibiting reagent binds to heterophilic antibodies



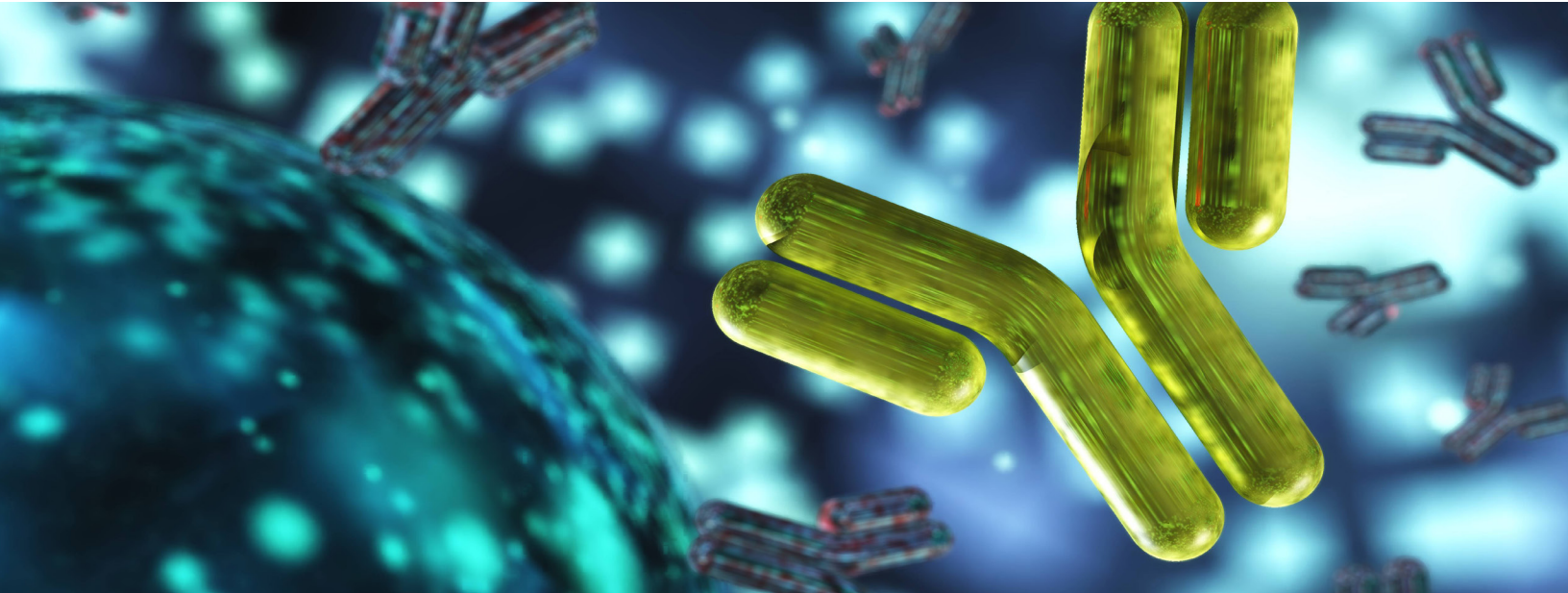
Bound heterophilic antibodies cannot cause a falsely elevated result

Rheumatoid Factors

Product Number:
RFACTOR

Rheumatoid Factor (RF) is typically IgM antibodies, but can also be IgG or IgA. RF is present in the serum of a majority of patients with rheumatoid arthritis. IgM Rheumatoid Factor is 20 mM Na Phosphate, with a pH of ~7. RF is mainly used as an interference control in immunoassays.

- Can be found in synovial fluid, pleural fluid, and pericardial fluid
- Ready-to-use solution
- Available in 1mL aliquots or custom sizes



Related Products



Human AB Serum



Rheumatoid Arthritis Biofluids



Human Systemic Lupus Erythematosus PBMCs

Other autoimmune disease related products available:

- Crohn's Disease
- Inflammatory Bowel Disease
- Osteoporosis Psoriasis
- Psoriasis
- Ulcerative Colitis
- Vitiligo and more!

For more tools for your immunology research visit BioIVT.com

BIOIVT

BioIVT.com

For Inquiries:

North America & Asia Pacific
T +1 516 483 1196 F +1 516 483 4683 E customerservice@bioivt.com

Europe, Middle East & Africa
T +44 (0) 1444 250010 F +44 (0) 1444 250066 E cseurope@bioivt.com