

AbX™

Goat Anti-Rabbit IgG Antibody

Catalog Number A009-10MG

Catalog Number A009-25MG



ARBOR
ASSAYS

FEATURES

- Affinity Purified Antibody to Rabbit IgG, Fc
- Extend primary antibody supplies
- Supplied as a PBS solution

INTRODUCTION

Solid phase immunoassay utilizing microtiter plates are the predominant assay format for the detection of low concentration small molecules in biological samples. The primary antibody to these small molecules serves two purposes; it confers specificity for the assay by selectively binding the small molecule, and, in binding it determines the sensitivity of the assay. Small molecule assays require a competitive immunoassay technique where the small molecule antigen binds to a limiting amount of the primary antibody in competition with a labeled reporter small molecule, such as HRP.

Primary antibodies can be coated on solid phases but the technique is difficult to reproducibly accomplish, as the dilution of the primary antibody has to be tightly controlled across all of the coated plates and the amounts of primary antibody required can be high. Secondary generic coated plates using goat anti-rabbit IgG antibody coating can be carried out easily and after being blocked and dried these plates are stable for many years. Typical primary antibody titers using coated generic plates can be 10 fold lower than directly coated primary antibodies.

- FORM:** Affinity purified against Rabbit IgG, Fc portion. In 10 mM sodium phosphate, 0.15M sodium chloride, 0.05% sodium azide, pH 7.2.
- CONCENTRATION:** Lot Specific C of A Supplied.
- STORAGE:** 4°C.
- IMMUNOGEN:** Highly purified Rabbit IgG Fc portion
- SPECIFICITY:** Based upon IEP this antibody reacts with heavy (γ) chains of rabbit IgG.
- CROSS REACTIVITY:** Based upon IEP, no reactivity is seen to non-immunoglobulin serum proteins, light chains on all rabbit immunoglobulins, or rabbit F(ab)₂ fragments.
- USES:** As a plate coating antibody for assays that utilize rabbit polyclonal antibodies as the primary antibody.
- SUGGESTED USE:** 10 μ g/mL in diluted Arbor Assays Coating Buffer, X108-10ML or X108-100ML.
- COUNTRY OF ORIGIN:** Goat serum was obtained from healthy animals of US origin and under the care of a registered veterinarian.