

AbX™

Donkey Anti-Sheep IgG Antibody



ARBOR
ASSAYS

Catalog Number A010-10MG Catalog Number A010-25MG
Catalog Number A010-50MG Catalog Number A010-100MG

FEATURES

- Affinity Purified Antibody to Sheep IgG, (H&L)
- 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 donkey anti-sheep 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 Sheep IgG 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 Sheep IgG whole molecule.
SPECIFICITY:	Based upon IEP this antibody reacts with all sheep IgG heavy (g) and light chains.
CROSS REACTIVITY:	Based upon IEP, no reactivity is seen to non-immunoglobulin serum proteins
USES:	As a plate coating antibody for assays that utilize sheep polyclonal antibodies as the primary antibody.
SUGGESTED USE:	10 µg/mL in diluted Arbor Assays Coating Buffer, X108-10ML or X108-100ML.
COUNTRY OF ORIGIN:	Donkey serum was obtained from healthy animals of US origin and under the care of a registered veterinarian.

FOR RESEARCH USE ONLY

info@ArborAssays.com • www.ArborAssays.com

11May20