

PRODUCT INFORMATION



Protein A/G-Coated 96-Well Strip Plates

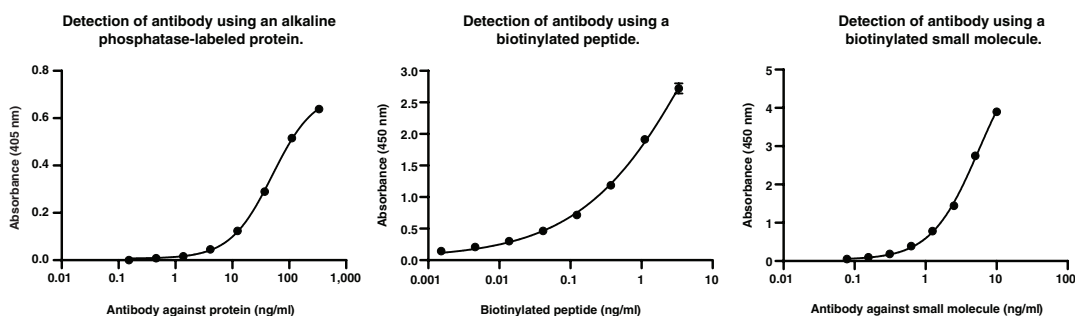
Item No. 33285

Storage: 4°C
Stability: ≥2 years

Description

Cayman's protein A/G-coated 96-well strip plates are high absorbance 96-well plates coated with a saturated concentration of highly pure protein A and protein G. The protein A/G-coated wells specifically bind the Fc region of IgGs resulting in an antibody orientation that will direct the antigen binding site away from the surface and offer maximum exposure of the binding site. These 96-well strip plates detect antibodies using labeled antigens, including small molecule, peptide, or protein antigens, and are ideal when a low amount of antibody is used. They are recommended for use in single antibody assays and are not suitable for multiple antibody assays, such as sandwich ELISAs. The protein A/G-coated 96-well strip plates provide an alternative to direct and passive adsorption methods of immobilizing antibodies that may render antibodies inactive. These 96-well strip plates have a high binding capacity (2-5 pmol antibody/well), low background, and minimal variation between wells (<7% well-to-well).

Image



The protein A/G-coated plates were incubated with antibody specific to a protein, peptide, or small molecule on a plate shaker for one hour at room temperature. The wells were rinsed to remove any unbound antibody, and the labeled antigen, alkaline phosphatase-labeled protein, biotinylated peptide, or biotinylated small molecule, was added to the wells and incubated for one hour at room temperature. The wells were rinsed to remove any unbound antigen, and the signal was detected according to the instructions for the specific detection system used.

WARNING

THIS PRODUCT IS FOR RESEARCH ONLY - NOT FOR HUMAN OR VETERINARY DIAGNOSTIC OR THERAPEUTIC USE.

SAFETY DATA

This material should be considered hazardous until further information becomes available. Do not ingest, inhale, get in eyes, on skin, or on clothing. Wash thoroughly after handling. Before use, the user must review the [complete](#) Safety Data Sheet, which has been sent via email to your institution.

WARRANTY AND LIMITATION OF REMEDY

Buyer agrees to purchase the material subject to Cayman's Terms and Conditions. Complete Terms and Conditions including Warranty and Limitation of Liability information can be found on our website.

Copyright Cayman Chemical Company, 05/01/2025

CAYMAN CHEMICAL

1180 EAST ELLSWORTH RD

ANN ARBOR, MI 48108 · USA

PHONE: [800] 364-9897

[734] 971-3335

FAX: [734] 971-3640

CUSTSERV@CAYMANCHEM.COM

WWW.CAYMANCHEM.COM