

PRODUCT INFORMATION



Anti-His-Coated 96-Well Strip Plate

Item No. 400701

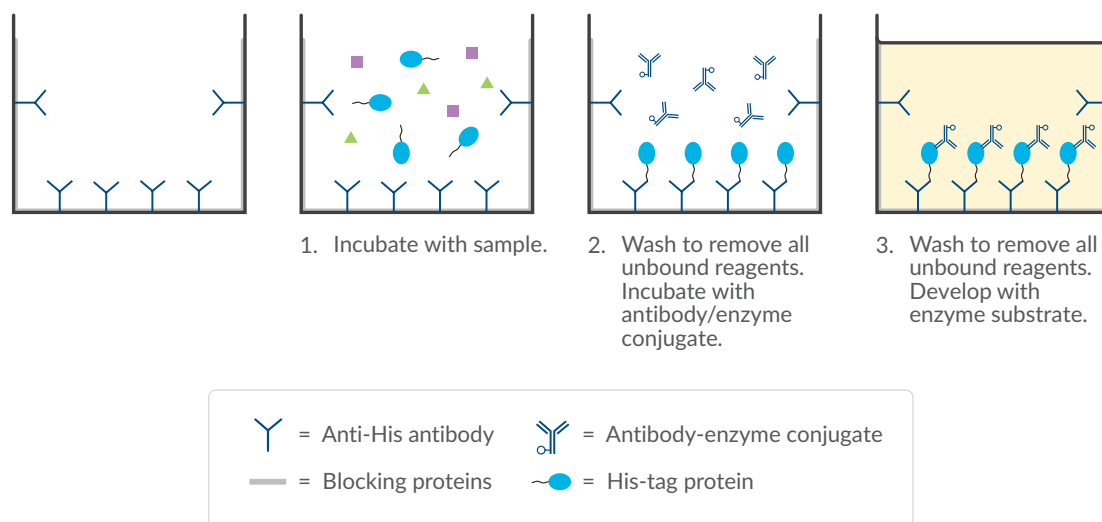
Storage: 4°C; ready to use as supplied

Description

Cayman's Anti-His-Coated 96-Well Strip Plate is coated with a His-tag recombinant monoclonal antibody and blocked with a proprietary formulation of proteins. The anti-His plate recognizes His tag-containing proteins with high specificity and binding capacity. This plate is versatile and adaptable for various applications, including but not limited to purification, quantifying His-tagged proteins in samples, studying protein-protein interactions, and monitoring of His-tag protein expression.

Suggested Applications

Figure 1. Quantification of His-tag proteins



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, 12/02/2025

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

PRODUCT INFORMATION

Figure 2. Assessment of protein-protein interactions

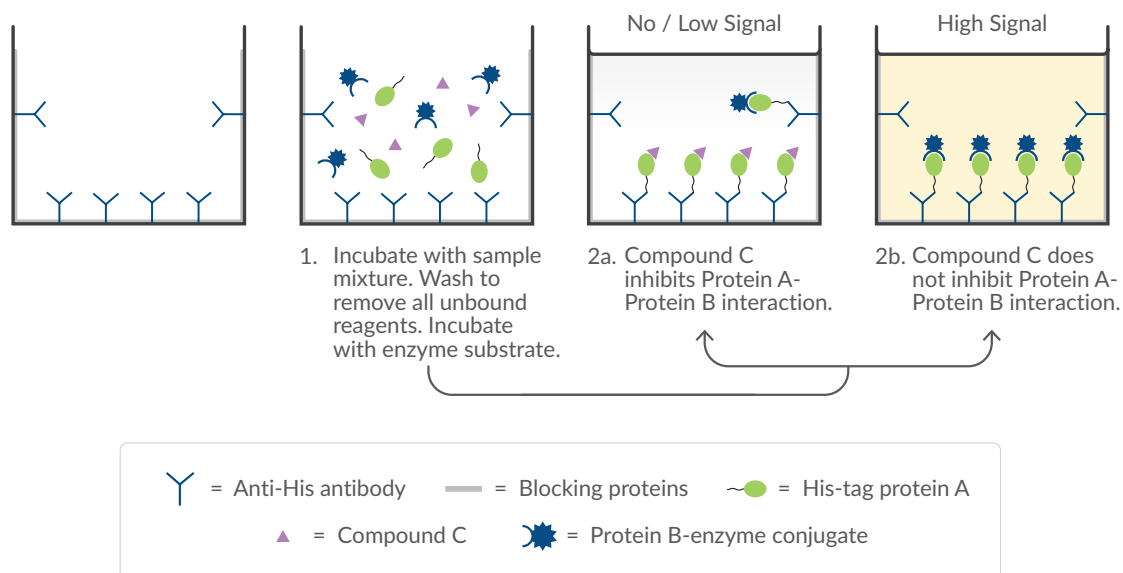
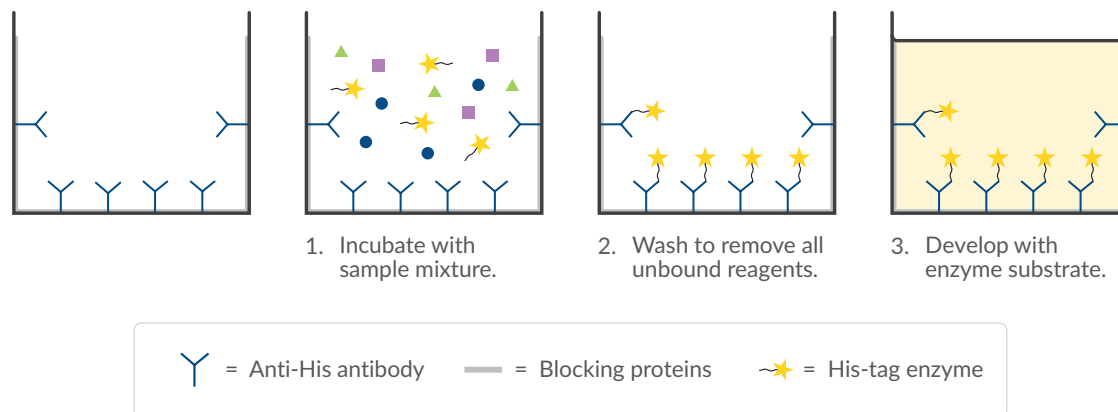


Figure 3. Assessment of His-tag enzyme expression



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