

# PRODUCT INFORMATION

Precoated (Mouse Anti-Rabbit IgG) ELISA 96-Well Strip Plate  
Item No. 400007



## Overview

---

**Contents:** This vial contains a 96-well plate pre-coated with Mouse Anti-Rabbit IgG.  
**Stability:**  $\geq 2$  years at 4°C

## Laboratory Procedures

---

This plate has been formulated and tested to work with Cayman's ELISA Kits. Detailed instructions for its use are contained in the respective Cayman ELISA Kit booklet.

## Precaution

---

This product may not perform as described if used with other assay reagents or protocols.

### 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, 12/08/2015

## 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