

PRODUCT INFORMATION

ELISA Blocking Buffer (1X)

Item No. 400829



Overview

Contents:	This vial contains 125 ml of ELISA Blocking Buffer (1X).
Storage:	4°C
Stability:	≥2 years
Application:	For use as a plate blocking buffer after coating for various ELISA applications.
Preparation:	Buffer is ready to use.

Suggested Protocol

1. Following plate coating at 100 µl/well, aspirate wells, then block with 200 µl of ELISA Blocking Buffer (1X).
2. Incubate the plate for 1 hour at room temperature.
3. Aspirate the ELISA Blocking Buffer (1X) from the wells. Plates can be used immediately or stored at 4°C.
4. For storage, plates must be dried for 1 hour at room temperature while inverted. If humidity levels are high, drying time may need to be extended. Store plates in well-sealed bags with a desiccant packet at 4°C.

NOTE: The amount of blocking buffer supplied is a sufficient volume for blocking five 96-well plates.

Precaution

The suggested protocol above is supplied as a guideline. Users may need to optimize the coating and blocking conditions based on specific applications.

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, 02/26/2026

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