

PRODUCT INFORMATION



Prostaglandin E2 Express AChE Tracer

Item No. 400140

Overview

Contents: This vial contains lyophilized Prostaglandin E₂ (PGE₂) Express Acetylcholinesterase (AChE) EIA Tracer (a covalent conjugate of PGE₂ and electric eel AChE (EC 3.1.1.7)).

Stability: ≥1 year at -20°C

Laboratory Procedures

Prostaglandin E2 Express AChE Tracer has been tested and formulated to work exclusively with Cayman's Prostaglandin E2 Express ELISA Kit (Item No. 500141). Please visit www.caymanchem.com/catalog/500141 for the kit protocol, procedures, and product handling.

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, 11/23/2015

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