

CERTIFICATE of ANALYSIS



EP4 Receptor (C-Term) Polyclonal PE Antibody

Item No. 10479 • Batch No. 0803808 • Expires 14MAY2027

Overview

Formulation:	100 µl of peptide affinity-purified polyclonal antibody conjugated to PE in PBS, pH 7.2, containing 50% glycerol with 0.02% sodium azide
Concentration:	1 mg/mL
Storage/stability:	-20°C/2yr
Applications:	FC and IF
Cross reactivity:	(-)EP1 receptor(-)EP2 receptor(-)EP3 receptor
Species reactivity:	(+)Human EP4 receptor(+)Mouse EP4 receptor(+)Ovine EP4 receptor(+)Rat EP4 receptor

Reviewed and approved by: Beth Hurst

Approved date: 27JUN2025

A handwritten signature in black ink that reads "Elizabeth Hurst".

WARNING

THIS PRODUCT IS FOR RESEARCH USE - NOT FOR HUMAN OR VETERINARY DIAGNOSTIC OR THERAPEUTIC USE. IT IS THE RESPONSIBILITY OF THE PURCHASER TO DETERMINE SUITABILITY FOR OTHER APPLICATIONS.

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, 10/12/2018

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