

CERTIFICATE of ANALYSIS



COX-2 Monoclonal Antibody (Clone 12C10)

Item No. 20198 • Batch No. 0824314 • Expires 15JUL2029

Overview

Formulation:	50 µg of protein G-purified IgG in PBS, pH 7.2, with 50% glycerol, 0.1% BSA, and 0.02% sodium azide
Concentration:	0.1 mg/mL
Storage/stability:	-20°C/3yr
Cross reactivity:	(-)COX-1 (all species)
Species reactivity:	(+)Human(+)Mouse(+)Ovine

Reviewed and approved by: Beth Hurst

Approved date: 20JUL2026

Beth 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