

PRODUCT INFORMATION



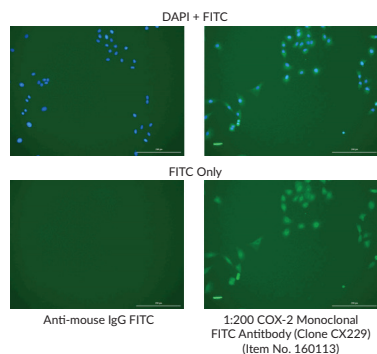
COX-2 Monoclonal FITC Antibody (Clone CX229)

Item No. 160113

Overview and Properties

Contents:	This vial contains 100 µg of FITC-labeled purified monoclonal antibody.
Synonyms:	Cyclooxygenase-2, PGHS-2, Prostaglandin H Synthase 2
Immunogen:	Synthetic peptide from the C-terminal region of human COX-2
Cross Reactivity:	(-) COX-1
Species Reactivity:	(+) Human, ovine; (-) Mouse, rabbit, rat; other species not tested
Uniprot No.:	P35354
Form:	Liquid
Storage:	-20°C (as supplied)
Stability:	≥1 year
Storage Buffer:	PBS, pH 7.2, with 50% glycerol, and 0.02% sodium azide
Clone:	CX229
Host:	Mouse
Isotype:	IgG1
Application:	Flow cytometry (FC), immunocytochemistry (ICC), and immunofluorescence (IF); the recommended starting concentration for FC is 1 µg antibody per 10 ⁶ cells and the recommended starting dilution for ICC is 1:50 and 1:200 for IF. Other applications were not tested, therefore optimal working concentration/dilution should be determined empirically.

Image



Immunofluorescence analysis of paraformaldehyde-fixed, A549 cells. After incubation with METTL16 Monoclonal Antibody (Item No. 29207), at a 1:200 dilution (or anti-mouse FITC as a negative control), followed by DAPI nuclear stain. Images show FITC alone or both fluorescence channels to highlight nuclear staining (where applicable).

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, 06/28/2023

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

PRODUCT INFORMATION



Description

This product is derived by labeling the COX-2 monoclonal antibody (Item No. 160112) with fluorescein.

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