

# PRODUCT INFORMATION



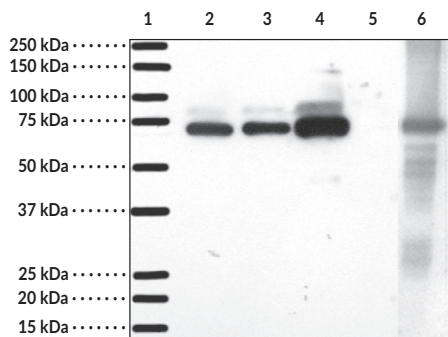
## COX-2 (mouse) Polyclonal Antiserum

Item No. 160116

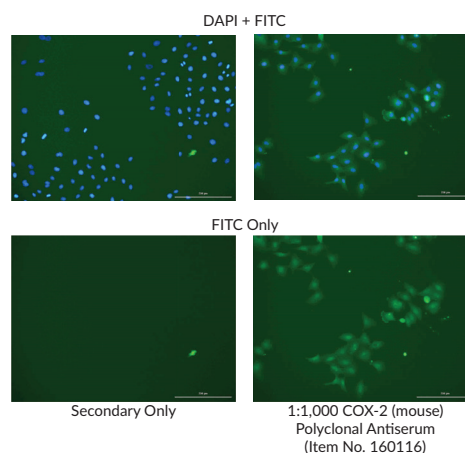
### Overview and Properties

**Contents:** This vial contains 100 µl of lyophilized antiserum.  
**Synonyms:** Cyclooxygenase 2, Prostaglandin H Synthase 2 (PGHS-2)  
**Immunogen:** Peptide from the C-terminal region of mouse COX-2  
**Cross Reactivity:** (+) COX-2; (-) COX-1  
**Species Reactivity:** (+) Human, mouse, ovine; other species not tested  
**Uniprot No.:** Q05769  
**Form:** Solid  
**Storage:** -20°C (as supplied)  
**Stability:** ≥1 year  
**Storage Buffer:** Polyclonal antiserum when reconstituted in 100 µl of double distilled water  
**Host:** Rabbit  
**Applications:** Immunofluorescence (IF), immunohistochemistry (IHC) and Western blot (WB); the recommended starting dilution for IHC is 1:300 and 1:1,000 for IF and WB. Other applications were not tested, therefore optimal working concentration/dilution should be determined empirically.

### Images



Lane 1: Precision Plus Protein Standard  
Lane 2: COX-2 (ovine) (0.05 µg)  
Lane 3: COX-2 (ovine) (0.1 µg)  
Lane 4: COX-2 (human recombinant) (5 µg)  
Lane 5: COX-1 (ovine) (0.1 µg)  
Lane 6: RAW264.7 microsomes (40 µg)



Immunofluorescence analysis of paraformaldehyde-fixed, A549 cells. After incubation with 1:1,000 COX-2 (mouse) Polyclonal Antiserum (Item No. 160116), at a 1:1,000 dilution (or negative control) cells were incubated with Goat Anti-Rabbit IgG FITC (Item No. 10006588), 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, 06/04/2021

**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

---

Cyclooxygenase 2 (COX-2) is a bifunctional enzyme that exhibits both COX and peroxidase activities and catalyzes the first step in the biosynthesis of prostaglandins, thromboxanes, and prostacyclins.<sup>1,2</sup> The COX component converts arachidonic acid to the hydroperoxy endoperoxide prostaglandin G<sub>2</sub> (PGG<sub>2</sub>; Item No. 17010), and the peroxidase component reduces the endoperoxide to the corresponding alcohol PGH<sub>2</sub> (Item No. 17020). COX2 expression is induced by a variety of stimuli, including phorbol esters, LPS, and cytokines and is responsible for the biosynthesis of PGs under acute inflammatory conditions.<sup>3,4</sup> Thus, COX-2 has been the focus of attention for nonsteroidal anti-inflammatory drug (NSAID) development. Cayman's COX-2 (mouse) Polyclonal Antiserum can be used for immunofluorescence (IF), immunohistochemistry (IHC), and Western blot (WB) applications. The antiserum recognizes a unique C-terminal region of COX-2 that is not present in COX-1, specifically detecting COX-2 at 70 kDa from human, mouse, and ovine samples.

## References

---

1. Nugteren, D.H. and Hazelhof, E. Isolation and properties of intermediates in prostaglandin biosynthesis. *Biochim. Biophys. Acta* **326(3)**, 448-461 (1973).
2. Hamberg, M., and Samuelsson, B. Detection and isolation of an endoperoxide intermediate in prostaglandin biosynthesis. *Proc. Natl. Acad. Sci. USA* **70(3)**, 899-903 (1973).
3. Kang, Y.-J., Mbonye, U.R., DeLong, C.J., *et al.* Regulation of intracellular cyclooxygenase levels by gene transcription and protein degradation. *Prog. Lipid Res.* **46(2)**, 108-25 (2007).
4. Blobaum, A.L. and Marnett, L.J. Structural and functional basis of cyclooxygenase inhibition. *J. Med. Chem.* **50(7)**, 1425-1441 (2007).

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