

Product Information



COX-2 (human) Western Ready Control

Catalog No. 10009624

Synonyms:	Prostaglandin H Synthase 2, Cyclooxygenase-2
Source:	Human recombinant N-terminal hexahistidine-tagged protein expressed in Sf21 cells
M_r:	73 kDa (as provided), 70 kDa (native)
Purity:	Whole cell lysate; a single band is detected on a western blot when probed with Cayman's COX-2 Monoclonal Antibody (Catalog No. 160112)
Supplied in:	Laemml buffer (2% SDS, 10% glycerol, 62.5 mM Tris-HCl, pH 6.8, 5% 2-mercaptoethanol, bromophenol blue)
Stability:	≥2 years at -20°C
Applications:	Positive control for western blotting

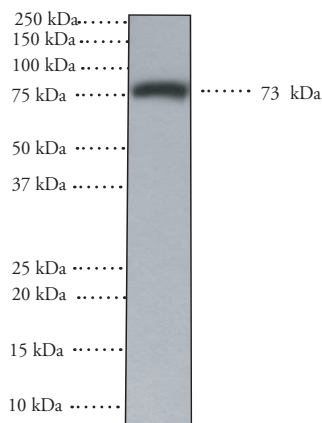
Cyclooxygenase-2 (COX-2) catalyzes the first step in the biosynthesis of PGs, thromboxanes, and prostacyclins: the conversion of arachidonic acid to PGH₂. Recent discoveries of the induction of the COX expression by a variety of stimuli such as phorbol esters, lipopolysaccharides, and cytokines led to the hypothesis that the inducible form of COX, COX-2, is responsible for the biosynthesis of prostaglandins under acute inflammatory conditions.¹ Thus, COX-2 has become the focus of attention for nonsteroidal anti-inflammatory drug (NSAID) development.

Laboratory Procedures

Western Ready Controls are designed to serve as positive controls on western blots. This COX-2 Western Ready Control is supplied pre-boiled in 55 µl of Laemml buffer. To use, thaw and load 10 µl of COX-2 Western Ready Control per well and follow a standard western blotting protocol. The amount needed per well will depend on the antibody and detection system used. Due to the presence of the 6X His-tag on the COX-2 provided in this control, the protein will migrate as a band ~3 kDa higher than the native, non-tagged protein.

Reference

1. Xie, W., Chipman, J.G., Robertson, D.L., *et al.* Expression of a mitogen-responsive gene encoding prostaglandin synthase is regulated by mRNA splicing. *Proc. Natl. Acad. Sci. USA* **88**, 2692-2696 (1991).



Western blot of human recombinant COX-2 probed with Cayman COX-2 Monoclonal Antibody (Catalog No. 160112)

Related Products

COX-2 (human recombinant) - Cat. No. 60122 • COX-2 (murine) Polyclonal Antibody - Cat. No. 160106 • COX-2 (human) Polyclonal Antibody - Cat. No. 160107 • COX-2 Monoclonal Antibody - Cat. No. 160112 • COX-2 (murine) Polyclonal Antiserum - Cat. No. 160116 • COX-2 (murine) Polyclonal Antibody - Affinity-Purified - Cat. No. 160126

WARNING: THIS PRODUCT IS NOT FOR HUMAN OR ANIMAL DISEASE DIAGNOSIS OR THERAPEUTIC DRUG USE.

MATERIAL SAFETY DATA

This material should be considered hazardous until information to the contrary becomes available. Do not ingest, swallow, or inhale. Do not get in eyes, on skin, or on clothing. Wash thoroughly after handling. This information contains some, but not all, of the information required for the safe and proper use of this material. Before use, the user must review the complete Material Safety Data Sheet, which has been sent under separate cover to the MSDS supervisor at your institution.

WARRANTY AND LIMITATION OF REMEDY

Cayman Chemical Company makes **no warranty or guarantee** of any kind, whether written or oral, expressed or implied, including without limitation, any warranty of fitness for a particular purpose, suitability and merchantability, which extends beyond the description of the chemicals hereof. Cayman **warrants only** to the original customer that the material will **meet our specifications at the time of delivery**.

Cayman will carry out its delivery obligations with due care and skill. Thus, in no event will Cayman have **any obligation or liability**, whether in tort (including negligence) or in contract, for any direct, indirect, incidental or consequential damages, even if Cayman is informed about their possible existence.

This limitation of liability does not apply in the case of intentional acts or negligence of Cayman, its directors or its employees. Buyer's **exclusive remedy** and Cayman's sole liability hereunder shall be limited to a **refund** of the purchase price, or at Cayman's option, the **replacement**, at no cost to Buyer, of all material that does not meet our specifications.

Said refund or replacement is conditioned on Buyer giving written notice to Cayman within thirty (30) days after arrival of the material at its destination. Failure of Buyer to give said notice within thirty (30) days shall constitute a waiver by Buyer of all claims hereunder with respect to said material.

For further details, please refer to our **Warranty and Limitation of Remedy located on our website and in our catalog**.

Copyright Cayman Chemical Company, 06/24/2008

Cayman Chemical

Mailing address

1180 E. Ellsworth Road
Ann Arbor, MI
48108 USA

Phone

(800) 364-9897
(734) 971-3335

Fax

(734) 971-3640

E-Mail

custserv@caymanchem.com

Web

www.caymanchem.com