

Product Information



COX-2 (murine) Polyclonal Antibody

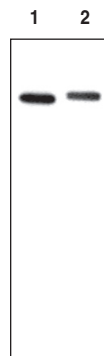
Item No. 160106 • Lot No. XXXX

Synonyms:	Cyclooxygenase 2 Polyclonal Antibody; Prostaglandin H Synthase 2 Polyclonal Antibody
Contents:	This vial contains (50-500 µg of peptide affinity-purified IgG, <i>lot specific</i>), lyophilized. Resuspend the IgG in 100-500 µl of purified water (<i>lot specific</i>) prior to use. Buffer will be 1X TBS.
Host:	Rabbit
Antigen:	Synthetic peptide from the C-terminal region of mouse/rat COX-2, conjugated to BSA (amino acids 570-598; DPQPTKTATINASASHSRLDDINPTVLIK). ¹
Cross-reactivity:	(+) Ovine, human, mouse, rabbit, monkey, guinea pig, and rat COX-2; (-) COX-1 (all species) Homology of peptide antigen with COX-2 from other species: Mouse DPQPTKTATINASASHSRLDDINPTVLIK Rat DPQPTKTATINASASHSRLDDINPTVLIK Human D P e l i K T v T I N A S s S r S g L D D I N P T V L I K Guinea pig D P Q l a K T v T I N A S A S H S R L e D I s P T V L I K Sheep D a h l T K T v T I N A S s S H S g L D D I N P T V L I K
Stability:	≥2 years at -20°C
Applications:	Western blot - 1:200. The antibody can also be used for immunohistochemistry (recommended starting dilution 1:300-1:600) and immunocytochemistry. The best working dilution must be determined empirically. ²⁻⁵
Concentration:	Varies by lot, from 0.1-1.0 mg/ml (50-500 µg/vial). Always 100 ml final working volume for western blotting.

COX catalyzes the first step in the biosynthesis of prostaglandins, thromboxanes, and prostacyclins: the conversion of arachidonic acid to prostaglandin H₂. Recent discoveries of the induction of COX 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.^{6,7} COX-2 is a 72 kDa protein which has been cloned from a variety of species including human, mouse, rat, and sheep.⁸⁻¹¹

References

1. Fletcher, B.S., Kujubu, D.A., Perrin, D.M., *et al. J. Biol. Chem.* **267**, 4338-4344 (1992).
2. Pomerantz, K.B., Summers, B., and Hajjar, D.P. *Biochemistry* **32**, 13624-13635 (1993).
3. Harris, R.C., McKanna, J.A., Akai, Y., *et al. J. Clin. Invest.* **94**, 2504-2510 (1994).
4. Langenbach, R., Morham, S.G., Tiano, H.F., *et al. Cell* **83**, 483-492 (1995).
5. Morham, S.G., Langenbach, R., Loftin, C.D., *et al. Cell* **83**, 473-482 (1995).
6. Xie, W., Chipman, J.G., Robertson, D.L., *et al. Proc. Nat. Acad. Sci USA* **88**, 2692-2696 (1991).
7. Smith, W.L., Garavito, R.M., and DeWitt, D.L. *J. Biol. Chem.* **271**, 33157-33160 (1996).
8. Hla, T. and Neilson, K. *Proc. Nat. Acad. Sci USA* **89**, 7384-7388 (1992).
9. Kujubu, D.A., Fletcher, B.S., Varnum, B.C., *et al. J. Biol. Chem.* **266**, 12866-12872 (1991).
10. Yamagata, K., Andreasson, K.I., Kaufmann, W.E., *et al. Neuron* **11**, 371-386 (1993).
11. Zhang, V., O'Sullivan, M., Hussain, H., *et al. Biochem. Biophys. Res. Commun.* **227**, 499-506 (1996).



Lane 1: Recombinant murine COX-2 (Cos-1 cell lysate)
Lane 2: Recombinant human COX-2 (Cos-1 cell lysate)

Related Products

COX-1 (ovine) - Item No. 60100 • COX-2 (ovine) - Item No. 60120 • COX Polyclonal Antibody - Item No. 160103 • COX-2 (human) Polyclonal Antibody - Item No. 160107 • COX-1 (ovine) Polyclonal Antiserum - Item No. 160108 • COX-1 (murine) Polyclonal Antibody - Item No. 160109 • COX-1 Monoclonal Antibody - Item No. 160110 • COX-1 Monoclonal FITC Antibody - Item No. 160111 • COX-2 Monoclonal Antibody - Item No. 160112 • COX-2 Monoclonal FITC Antibody - Item No. 160113 • COX-2 (murine) Polyclonal Antiserum - Item No. 160116 • COX-1 Monoclonal PE Antibody - Item No. 160120 • COX-2 Monoclonal PE Antibody - Item No. 160122 • COX-2 (murine) Blocking Peptide - Item No. 360106 • COX-2 (ovine) Electrophoresis Standard - Item No. 360120 • Goat Anti-Rabbit IgG HRP - Item No. 10004301

WARNING: THIS PRODUCT IS FOR LABORATORY RESEARCH ONLY: NOT FOR ADMINISTRATION TO HUMANS. NOT FOR HUMAN OR VETERINARY DIAGNOSTIC OR THERAPEUTIC 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 via email to 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, 08/31/2011

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