

# PRODUCT INFORMATION



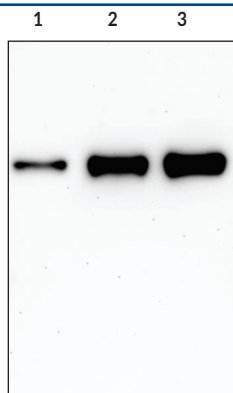
## COX-2 (ovine) Electrophoresis Standard

Item No. 360120

### Overview and Properties

<b>Contents:</b>	This vial contains 5 µg of purified COX-2
<b>Synonyms:</b>	Cyclooxygenase 2 (ovine) Electrophoresis Standard, Prostaglandin H Synthase 2 (ovine) Electrophoresis Standard
<b>Source:</b>	Isolated from sheep placenta. <sup>1</sup>
<b>MW:</b>	72 kDa/subunit
<b>Storage:</b>	-80°C (as supplied)
<b>Stability:</b>	≥1 year
<b>Purity:</b>	≥95%
<b>Storage Buffer:</b>	Laemmli buffer (2% SDS, 10% glycerol, 0.70 M Tris-HCl, pH 6.8, 1.5% 2-mercaptoethanol, 0.04 % bromophenol blue)
<b>Applications:</b>	This enzyme has been denatured and has no catalytic activity. It can be used as a standard for western blots and electrophoresis.

### Images



Lane 1: COX-2 Standard (10 ng)  
Lane 2: COX-2 Standard (50 ng)  
Lane 3: COX-2 Standard (100 ng)

Probed with COX-2 Polyclonal Antibody (Item No. 160126)

**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, 07/11/2019

**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 catalyzes the first step in the biosynthesis of prostaglandins, thromboxanes, and prostacyclins: the conversion of arachidonic acid to prostaglandin H<sub>2</sub>. Recent discoveries of the induction of cyclooxygenase by a variety of stimuli such as phorbol esters, lipopolysaccharides, and cytokines led to the hypothesis that the inducible form of cyclooxygenase, COX-2, is responsible for the biosynthesis of prostaglandins under acute inflammatory conditions.<sup>2,3</sup> COX-2 is a 72 kDa protein which has been cloned from a variety of species including human, mouse, rat, and sheep.<sup>4-7</sup>

## References

---

1. Johnson, J.L., Wimsatt, J., Buckel, S.D., *et al.* *Arch. Biochem. Biophys.* **324**, 26-34 (1995).
2. Xie, W., Chipman, J.G., Robertson, D.L., *et al.* *Proc. Nat. Acad. Sci USA* **88**, 2692-2696 (1991).
3. Smith, W.L., Garavito, R.M., and DeWitt, D.L. *J. Biol. Chem.* **271**, 33157-33160 (1996).
4. Hla, T. and Neilson, K. *Proc. Nat. Acad. Sci USA* **89**, 7384-7388 (1992).
5. Kujubu, D.A., Fletcher, B.S., Varnum, B.C., *et al.* *J. Biol. Chem.* **266**, 12866-12872 (1991).
6. Yamagata, K., Andreasson, K.I., Kaufmann, W.E., *et al.* *Neuron* **11**, 371-386 (1993).
7. Zhang, V., O'Sullivan, M., Hussain, H., *et al.* *Biochem. Biophys. Res. Commun.* **227**, 499-506 (1996).

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