

PRODUCT INFORMATION



COX-1 (mouse) Polyclonal Antibody

Item No. 160109

Overview and Properties

Contents:	This vial contains 500 µl of peptide affinity-purified antibody.
Synonyms:	Cyclooxygenase 1, PGHS-1, Prostaglandin H Synthase 1
Immunogen:	Peptide from the internal region of mouse COX-1
Cross Reactivity:	(-) COX-2
Species Reactivity:	(+) Mouse and ovine; (-) Human; other species not tested
Uniprot No.:	P22437
Form:	Liquid
Storage:	-20°C (as supplied)
Stability:	≥1 year
Storage Buffer:	PBS, pH 7.2, with 50% glycerol and 0.02% sodium azide
Host:	Rabbit
Application:	Western blot; the recommended starting dilution is 1:200. Other applications were not tested, therefore optimal working concentration/dilution should be determined empirically.

Image



Lane 1: Rat brain homogenate (30 µg)

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, 04/17/2019

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

COX-1 is responsible for the production of prostaglandins essential for the normal function of many organs including the stomach and kidney.¹ COX-1 is generally described as being constitutively expressed, although COX-1 expression is regulated developmentally and in response to a variety of other stimuli.²⁻⁵ Mouse COX-1 has a molecular weight of 70,000 and is expressed in nearly all tissues of the body.^{1,6}

References

1. Smith, W.L., Garavito, R.M., and DeWitt, D.L. Prostaglandin endoperoxide H synthases (cyclooxygenases)-1 and -2. *J. Biol. Chem.* **271**, 33157-33160 (1996).
2. Brannon, T.S., North, A.J., Wells, L.B., *et al.* Prostacyclin synthesis in ovine pulmonary artery is developmentally regulated by changes in cyclooxygenase-1 gene expression. *J. Clin. Invest.* **93**, 2230-2235 (1994).
3. Funk, C.D., Funk, L.B., Kennedy, M.E., *et al.* Human platelet/erythroleukemia cell prostaglandin G/H synthase: cDNA cloning, expression, and gene chromosomal assignment. *FASEB J.* **5**, 2304-2312 (1991).
4. Samet, J.M., Fasano, M.B., Fonteh, A.N., *et al.* Selective induction of prostaglandin G/H synthase I by stem cell factor and dexamethasone in mast cells. *J. Biol. Chem.* **270**, 8044-8049 (1995).
5. Hla, T. and Maciag, T. Cyclooxygenase gene expression is down-regulated by heparin-binding (acidic fibroblast) growth factor-1 in human endothelial cells. *J. Biol. Chem.* **266**, 24059-24063 (1991).
6. DeWitt, D.L., El-Harith, E.A., Kraemer, S. A., *et al.* The aspirin and heme-binding sites of ovine and murine prostaglandin endoperoxide synthases. *J. Biol. Chem.* **265**, 5192-5198 (1990).

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