

# PRODUCT INFORMATION



## Mouse Macrophage Microsomal Immunoblot Control

Item No. 10004910

### Overview and Properties

<b>Contents:</b>	This vial contains 500 µg of mouse macrophage microsomal protein.
<b>Synonym:</b>	Mouse Macrophage Microsomes Protein
<b>Form:</b>	Liquid
<b>Storage:</b>	-20°C (as supplied)
<b>Stability:</b>	≥3 years
<b>Storage Buffer:</b>	2% SDS, 10% glycerol, 0.70 M Tris-HCl pH 6.8, 1.5 % 2-mercaptoethanol, 0.04 % bromophenol blue
<b>Application:</b>	Western blot (WB); the recommended starting concentration is 10 µl/lane. Other applications were not tested, therefore optimal working concentration/dilution should be determined empirically.

### Description

Cyclooxygenase 2 (COX-2) is a 72 kDa protein that has been cloned from a variety of species including human, mouse, rat, and ovine.<sup>1-5</sup> This preparation of COX-2 microsomes was prepared by solubilization and dilution of the 100,000 x g pellet from the lysate of cytokine-stimulated mouse macrophage cultures. Cayman's preparation of mouse macrophage microsomes is a ready-to-use sample for SDS-PAGE. As such, it is an excellent positive control for the following Cayman Chemical polyclonal antibodies:

Item No.	Antibody Name
160106	COX-2 (mouse) Polyclonal Antibody (aa 570-598)
160126	COX-2 (mouse) Polyclonal Antibody (aa 584-598)
160116	COX-2 (mouse) Polyclonal Antiserum
101750	EP <sub>2</sub> Receptor Polyclonal Antibody

### References

1. Smith, W.L., Garavito, R.M., and DeWitt, D.L. Prostaglandin endoperoxide H synthases (cyclooxygenases)-1 and -2. *J. Biol. Chem.* **271**(52), 33157-33160 (1996).
2. Hla, T. and Neilson, K. Human cyclooxygenase-2 cDNA. *Proc. Natl. Acad. Sci. USA* **89**(16), 7384-7388 (1992).
3. Kujubu, D.A., Fletcher, B.S., Varnum, B.C., *et al.* TIS10, a phorbol ester tumor promoter-inducible mRNA from Swiss 3T3 cells, encodes a novel prostaglandin synthase/cyclooxygenase homologue. *J. Biol. Chem.* **266**(20), 12866-12872 (1991).
4. Yamagata, K., Andreasson, K.I., Kaufmann, W.E., *et al.* Expression of a mitogen-inducible cyclooxygenase in brain neurons: Regulation by synaptic activity and glucocorticoids. *Neuron* **11**(2), 371-386 (1993).
5. Zhang, V., O'Sullivan, M., Hussain, H., *et al.* Molecular cloning, functional expression, and selective regulation of ovine prostaglandin H synthase-2. *Biochem. Biophys. Res. Commun.* **227**(2), 499-506 (1996).

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, 10/11/2023

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