

# PRODUCT INFORMATION



## COX-2 (human) Blocking Peptide

Item No. 360107

### Overview and Properties

<b>Contents:</b>	This vial contains 200 µg of peptide.
<b>Synonyms:</b>	Cyclooxygenase 2, PGHS-2, Prostaglandin H Synthase 2
<b>Form:</b>	Lyophilized
<b>Storage:</b>	-20°C (as supplied)
<b>Storage Buffer:</b>	Peptide when reconstituted in 200 µl of PBS or distilled water
<b>Stability:</b>	≥3 years

### Procedures

This vial contains 200 µg of lyophilized peptide. The peptide is derived from the human COX-2 sequence. This blocking peptide can be used in conjunction with Cayman's COX-2 (human) Polyclonal Antibody (Item No. 160107) or the COX-2 (human) monoclonal antibodies (Item Nos. 160112 | 160113 | 160122) to block protein-antibody complex formation during immunochemical analysis for COX-2.

To block antibody/protein complex formation, the following procedure is recommended:

1. Mix the COX-2 (human) Polyclonal Antibody (Item No. 160107) and blocking peptide together in a 1:1 (v/v) ratio in a microfuge tube. For example, mix 20 µl of antibody and 20 µl of peptide.\*
2. Incubate for one hour at room temperature with occasional mixing prior to further dilution and application of the mixture to the immunoblot.
3. Dilute the mixture to the final working antibody concentration and apply to the slide or membrane as usual.

\*This is a recommended mixture. The minimum amount of peptide needed for complete blocking has not been precisely determined and may vary depending on the sample being analyzed. The amount of peptide required may need to be increased if sufficient blocking does not occur.

**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, 12/13/2023

**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