COX-2 polyclonal antibody

This antibody is covered by our Worry-Free Guarantee.

Ordering Information

Order Online »

BML-SA237-0025	25µg
BML-SA237-0100	100µg

Manuals, SDS & CofA

View Online »

Handling & Storage

Long Term Storage -80°C

Shipping Dry Ice

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name Prostaglandin H2 synthase, Cyclooxygenase-2

Application IHC, IP, WB

Application Notes Detects a band of ~68-70kDa by Western blot.

Crossreactivity Does not cross-react with COX-1.

Formulation Liquid. In PBS.

Host Rabbit

Immunogen Synthetic peptide corresponding to a portion of C-terminal

mouse COX-2.

Species Reactivity Human, Mouse, Sheep

UniProt ID Q05769

Worry-free Guarantee This antibody is covered by our Worry-Free Guarantee

.

Last modified: May 29, 2024

