

---

## COX-2 Polyclonal Antibody

(Cat. No. A-0641)

---

### Background

Cyclooxygenase (COX), also known as prostaglandin-endoperoxide synthase (PTGS), is a key enzyme in the conversion of arachidonic acid to Prostaglandin H<sub>2</sub>. Two isoforms that exist for this enzyme are COX-1 and COX-2. COX-1 is involved in the synthesis of prostaglandin and thromboxane in many cell types, including gastro-intestinal tract and blood platelets. COX-2 is an inducible enzyme that plays a significant role in the biosynthesis of prostaglandins (PGs) under acute inflammatory conditions. COX-2 can be induced by numerous stimuli and has been found to be overexpressed in adenocarcinomas and other tumors. COX-2 inhibitors have analgesic and anti-inflammatory function by blocking the transformation of arachidonic acid into prostaglandin H<sub>2</sub> selectively.

### Description

Rabbit polyclonal antibody to COX-2

### Formulation

Liquid. In PBS.

### Immunogen

Synthetic peptide corresponding to a portion of C-terminal mouse COX-2.

### Specificity

Human, Mouse, Sheep

### Purification

Serum

### Storage

Store at -80°C. Avoid multiple freeze/thaw cycles.

### Application

IHC, IP, WB

### Ordering Information

Products	Size	Cat. No.
COX-2 Polyclonal Antibody	25 µg	A-0641-025
	100 µg	A-0641-100

*This product is for research purposes only. Not intended for use in diagnostic procedures.*