

COX-2 (human), ELISA kit

Most sensitive COX-II immunoassay kit commercially available for inflammation research.

The COX-2 (human), EIA kit is a colorimetric immunometric enzyme immunoassay kit with results in 2 hours.

Cyclooxygenases-1 and 2 (COX-1 and COX-2) also known as Prostaglandin endoperoxide H synthases-1 and 2 are integral membrane proteins which catalyze the first step of prostaglandin, thromboxane, and prostacyclin production from arachidonic acid. The COX-1 and COX-2 enzymes are important pharmacologically as targets of nonsteroidal anti-inflammatory drugs (NSAIDs) including aspirin, ibuprofen and the COX-2 inhibitors. Inhibition of cyclooxygenases with NSAIDs acutely reduces inflammation, pain, and fever.

Citations: 5

[View Online »](#)

Ordering Information

[Order Online »](#)

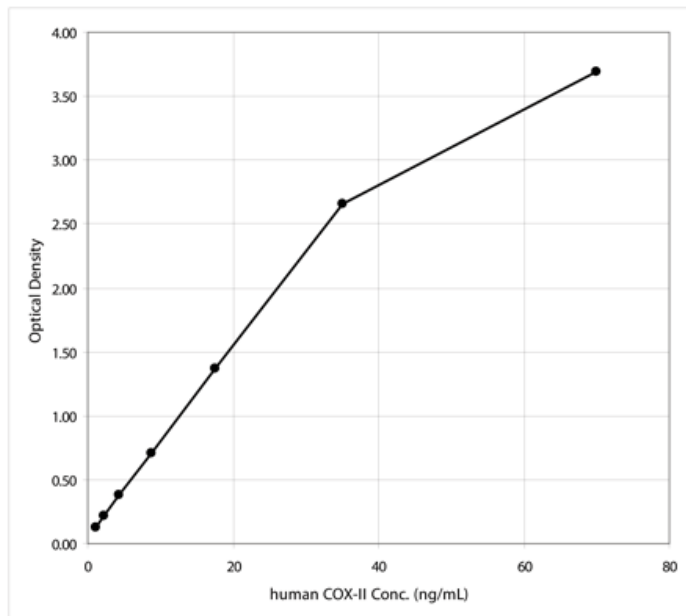
ADI-900-094

96 wells

Manuals, SDS & CofA

[View Online »](#)

- Easy-to-use - ready-to-use liquid color-coded reagents reduce errors
- Reliable - reproducible results day-after-day and lot-after-lot
- Fast - results in just 2 hours from up to 39 samples in duplicate



Handling & Storage

Long Term Storage +4°C

Shipping Blue Ice

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name Cyclooxygenase-2, Prostaglandin H synthase 2

Application Colorimetric detection, ELISA

Application Notes For the quantitative determination of human COX-2 in cell lysates.

Assay Time 2 hours

Compatibility This product is compatible with the [Absorbance 96 Plate Reader](#).

Contents Microtiter plate, Labeled antibody concentrate, Antibody diluent, Assay buffer, Wash buffer concentrate, Standard, TMB Substrate, Stop solution

Crossreactivity Human COX-2 (100%), human COX-1 (<0.1%)

Sensitivity 0.25 ng/ml (range 1.09 – 70 ng/ml)

Species Reactivity Human

UniProt ID P35354

Wavelength 450 nm



ENZO LIFE SCIENCES,
INC.
Phone: 800.942.0430
[info-
usa@enzolifesciences.com](mailto:info-usa@enzolifesciences.com)

European Sales Office
ENZO LIFE SCIENCES
(ELS) AG
Phone: +41 61 926 8989
[info-
eu@enzolifesciences.com](mailto:info-eu@enzolifesciences.com)

Belgium, The Netherlands
& Luxembourg
Phone: +32 3 466 0420
[info-
fr@enzolifesciences.com](mailto:info-fr@enzolifesciences.com)

France
Phone: +33 472 440 655
[info-
fr@enzolifesciences.com](mailto:info-fr@enzolifesciences.com)

Germany
Phone: +49 7621 5500 526
[info-
de@enzolifesciences.com](mailto:info-de@enzolifesciences.com)

UK & Ireland
Phone (UK customers):
0845 601 1488
Phone: +44 1392 825900
[info-
uk@enzolifesciences.com](mailto:info-uk@enzolifesciences.com)