

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

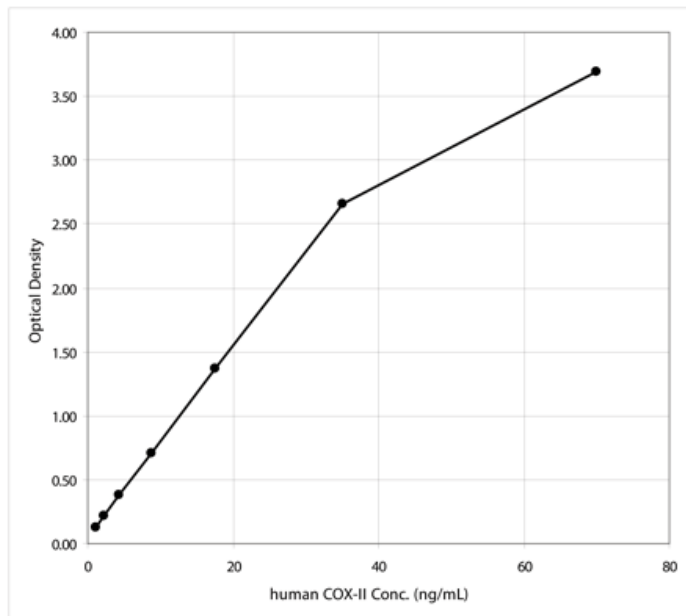
[Order Online »](#)

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

Last modified: May 29, 2024



ENZO LIFE SCIENCES,
INC.
Phone: 800.942.0430
info-usa@enzolifesciences.com

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

Belgium, The Netherlands
& Luxembourg
Phone: +32 3 466 0420
info-be@enzolifesciences.com

France
Phone: +33 472 440 655
info-fr@enzolifesciences.com

Germany
Phone: +49 7621 5500 526
info-de@enzolifesciences.com

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