

GRO/CINC-1 (rat), ELISA kit

The **GRO/CINC-1 (rat), ELISA kit** is a **colorimetric immunometric enzyme immunoassay kit**

The GRO/CINC-1 (rat), EIA kit is a colorimetric immunometric enzyme immunoassay kit with results in 2 hours. Absorbance is read at 450 nm. Assay up to 39 samples in duplicate per kit.

GRO/CINC-1 (Growth related oncogene/Cytokine-induced neutrophil chemoattractant-1) is a member of the α (CXC) subfamily of chemokines. It is the rat homologue to human interleukin-8 (IL-8). Three additional rat CXC chemokines (CINC-2a, CINC-2b, CINC-3/MIP-2) have been identified. CINC-3 appears to be the rat counterpart of human GROs.

Citations: 9

[View Online »](#)

Ordering Information

[Order Online »](#)

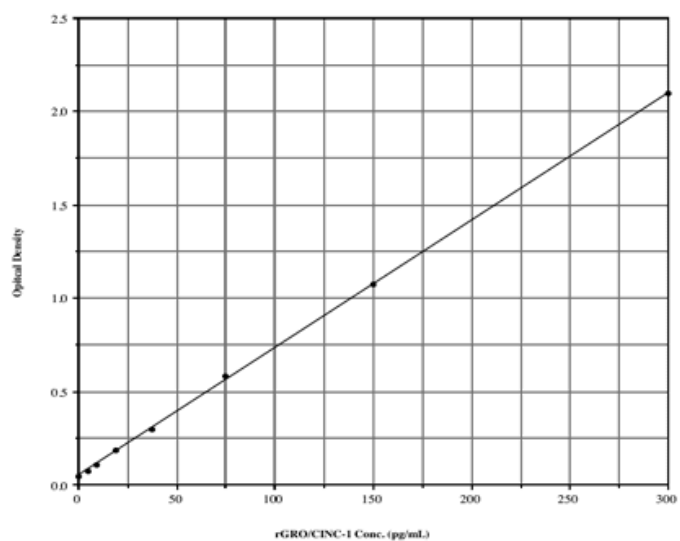
ADI-900-074

96 wells

Manuals, SDS & CofA

[View Online »](#)

- Fast - fully quantitative results in just 2 hours from up to 39 samples in duplicate
- Sensitive - measure < 2 pg/ml GRO/CINC-1
- Higher throughput - test 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	CXCL1
Application	Colorimetric detection, ELISA
Application Notes	For the quantitative determination of rat GRO/CINC-1 in culture supernatants, plasma, and serum.
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	Rat GRO/CINC-1 (100%) and <0.1%: IL-1β, MIP-1, GRO-CINC-3, GRO-CINC-2α, MCP-1, GRO-CINC-2β, Rantes , mouse GROβ, MIP-2
Sensitivity	1.99 pg/ml (range 4.7 – 300 pg/ml)
Species Reactivity	Rat
Technical Info / Product Notes	Cited samples: Cytokine ELISAs Cited Samples
UniProt ID	P14095
Wavelength	450 nm



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-fr@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

Last modified: May 29, 2024