

**Modified Reagent
Preparation and
New Storage
Conditions**



GRO/CINC-1 (rat), ELISA kit

Catalog No. ADI-900-074

96 Determination Kit

Table of Contents

Description	Page	2
Introduction		2
Precautions		2
Materials Supplied		3
Storage		3
Materials Needed but Not Supplied		3
Sample Handling		4
Procedural Notes		4
Reagent Preparation		5
Assay Procedure		6
Calculation of Results		7
Typical Results		7
Typical Standard Curve		8
Performance Characteristics		9
Sample Dilution Recommendations		11
References		11
Limited Warranty		12

FOR RESEARCH USE ONLY. NOT FOR USE IN DIAGNOSTIC PROCEDURES.

Description

The GRO/CINC-1 (rat), ELISA kit is a complete kit for the quantitative determination of rat GRO/CINC-1 in biological fluids. Please read the complete kit insert before performing this assay. The kit uses a polyclonal antibody to rat GRO/CINC-1 immobilized on a microtiter plate to bind the rat GRO/CINC-1 in the standards or sample. A recombinant rat GRO/CINC-1 Standard is provided in the kit. After a short incubation the excess sample or standard is washed out and a polyclonal antibody to rat GRO/CINC-1 labeled with the enzyme Horseradish peroxidase is added. This labeled antibody binds to the rat GRO/CINC-1 captured on the plate. After a short incubation the excess labeled antibody is washed out and substrate is added. The substrate reacts with the labeled antibody bound to the rat GRO/CINC-1 captured on the plate. After a short incubation, the enzyme reaction is stopped and the color generated is read at 450 nm. The measured optical density is directly proportional to the concentration of rat GRO/CINC-1 in either standards or samples. For further explanation of the principles and practice of immunoassays please see the excellent books by Chard¹ or Tijssen².

Introduction

Cytokine-induced Neutrophil Chemoattractant-1 (CINC-1) was originally purified from media conditioned by IL-1 β stimulated rat kidney epithelioid cells (NRK-52E)³⁻⁷. Watanabe's group at Toyoma Medical and Pharmaceutical University identified the amino acid sequence that encodes for rat CINC-1 in 1989. CINC-1 is a member of the alpha (CXC) subfamily of chemokines. Three additional rat CXC chemokines (CINC-2 α , CINC-2 β , CINC-3/MIP-2) have been identified. The protein sequence of CINC-1 is 63-67% identical to that of CINC-2 α , CINC-2 β , and CINC-3/MIP-2. In addition the Growth Related Oncogene (GRO), GRO α , GRO β and GRO γ shares 68%, 71% and 69% identity with CINC-1. It has been suggested that CINC-1 is the rat counter-part of human GROs.

Precautions

FOR RESEARCH USE ONLY. NOT FOR USE IN DIAGNOSTIC PROCEDURES.

1. Stop Solution is a 1N sulfuric acid solution. This solution is caustic; care should be taken in use.
2. The activity of the Horseradish peroxidase conjugate is affected by nucleophiles, such as azide, cyanide and hydroxylamine.
3. We test this kit's performance with a variety of samples, however it is possible that high levels of interfering substances may cause variation in assay results.
4. The rat GRO/CINC-1 Standard provided, Catalog No. 80-0676, should be handled with care, because of the known and unknown effects of GRO/CINC-1.

Materials Supplied

1. **rat GRO/CINC-1 Microtiter Plate, One Plate of 96 Wells, Catalog No. 80-0674**
A strip microtiter plate coated with rabbit antibody specific to rat GRO/CINC-1.
2. **rat GRO/CINC-1 Labeled Antibody Concentrate, 0.4 mL, Catalog No. 80-1381**
Rabbit antibody to rat GRO/CINC-1 conjugated to Horseradish peroxidase.
3. **Assay Buffer, 30 mL, Catalog No. 80-0170**
Phosphate buffered saline containing proteins and detergents.
4. **Labeled Antibody Diluent, 12 mL, Catalog No. 80-0182**
Phosphate buffered saline containing proteins and detergents.
5. **Wash Buffer Concentrate, 50 mL, Catalog No. 80-0171**
Phosphate buffered saline containing detergents.
6. **rat GRO/CINC-1 Standard, 2 each, Catalog No. 80-0676**
Two vials containing 300 pg each of recombinant rat GRO/CINC-1.
7. **TMB Substrate, 15 mL, Catalog No. 80-1342**
A solution of 3,3',5,5' tetramethyl benzidine (TMB) and hydrogen peroxide. Ready to use.
8. **Stop Solution, 12 mL, Catalog No. 80-0176**
A 1N solution of sulfuric acid in water. Keep tightly capped. Caution: **Caustic.**
9. **rat GRO/CINC-1 Assay Layout Sheet, 1 each, Catalog No. 30-0139**
10. **Plate Sealer, 2 each, Catalog No. 30-0012**

Storage

All components of this kit are stable at 4°C until the kit's expiration date.

Materials Needed but Not Supplied

1. Deionized or distilled water.
2. Precision pipets for volumes between 100 μ L and 1,000 μ L.
3. Disposable test tubes for dilution of samples and standards.
4. Repeater pipet for dispensing 100 μ L.
5. Disposable beakers for diluting buffer concentrates.
6. Graduated cylinders.
7. A 37°C incubator.
8. Adsorbent paper for blotting.
9. Microplate reader capable of reading at 450 nm, preferably with correction between 570 nm and 590 nm.
10. Graph paper for plotting the standard curve.

Sample Handling

The GRO/CINC-1 (rat), ELISA is compatible with rat GRO/CINC-1 (rat IL-8) samples in a wide range of matrices. Samples diluted sufficiently into Assay Buffer can be read directly from the standard curve. Please refer to the Sample Recovery recommendations on page 11 for details of suggested dilutions.

Culture fluids, serum and plasma are suitable for use in the assay. Samples containing visible precipitate should be clarified prior to use in the assay. Do not use grossly hemolyzed or lipemic specimens. Samples in the majority of tissue culture media, including those containing fetal bovine serum, can also be read in the assay if diluted into Assay Buffer. Users should only use standard curves generated in Assay Buffer to calculate concentrations of rat GRO/CINC-1.

Procedural Notes

1. Do not mix reagents from different kit lots or use reagents beyond the kit expiration date.
2. Allow all reagents to warm to room temperature for at least 30 minutes before opening.
3. Standards can be made up in either glass or plastic tubes.
4. Pre-rinse the pipet tip with reagent, use fresh pipet tips for each sample, standard and reagent.
5. Pipet standards and samples to the bottom of the wells.
6. Add the reagents to the side of the well to avoid contamination.
7. This kit uses plates with removable strips. Unused strips must be kept desiccated at 4 °C in the sealed bag provided. The strips should be used in the frame provided.
8. **Prior to addition of standard, antibody, and substrate, ensure that there is no residual wash buffer in these wells. Any remaining wash buffer may cause variation in assay results.**

Reagent Preparation

1. Wash Buffer

Prepare Wash Buffer by diluting 25 mL of the supplied concentrate with 975 mL of deionized water. This can be stored at 4°C until the kit expiration date, or for 3 months, whichever is earlier.

2. rat GRO/CINC-1 Standards

Add 500 µL of deionized water to the rat GRO/CINC-1 Standard. Let it sit at room temperature for 5 minutes. Mix it gently. This solution contains 600 pg/mL rat GRO/CINC-1.

Label seven 12 x 75 mm glass tubes #1 through 7. Pipet 230 µL of Assay Buffer into tubes #1 through #7. Add 230 µL of the 600 pg/mL standard to tube #1. Vortex. Add 230 µL of tube #1 to tube #2 and vortex thoroughly. Continue this for tubes #3 through #7.

The concentration of rat GRO/CINC-1 in tubes #1 through #7 will be 300, 150, 75, 37.5, 18.8, 9.38, and 4.69 pg/mL respectively. See rat GRO/CINC-1 Assay Layout Sheet for dilution details. STORE RECONSTITUTED STANDARD AT -20 °C OR BELOW; avoid repeated freeze/thaws.

3. Labeled Antibody Conjugate

Just before use, the rat GRO/CINC-1 Labeled Antibody Concentrate must be diluted 1:30 into the Labeled Antibody Diluent in a clean test tube and vortexed thoroughly. For example, if using one 8 well strip, dilute 30 µL of the Labeled Antibody Concentrate into 870 µL of the Labeled Antibody Diluent.

Assay Procedure

Bring all reagents to room temperature for at least 30 minutes prior to opening.

All standards and samples should be run in duplicate.

1. Refer to the Assay Layout Sheet to determine the number of wells to be used and put any remaining wells with the desiccant back into the pouch and seal the ziploc. Store unused wells at 4°C.
2. Pipet 100 µL of Assay Buffer into the S0 (0 pg/mL Standard) wells.
3. Pipet 100 µL of Standards #1 through #7 into the appropriate wells.
4. Pipet 100 µL of the Samples into the appropriate wells.
5. Tap the plate gently to mix the contents.
6. Seal the plate and incubate at 37°C for 1 hour.
7. Empty the contents of the wells and wash by adding 400 µL of wash solution to every well. Repeat the wash 6 more times for a total of **7 washes**. After the final wash, empty or aspirate the wells, and firmly tap the plate on a lint free paper towel to remove any remaining wash buffer.
8. Pipet 100 µL of the Labeled Antibody into each well, except the Blank.
9. Seal the plate and incubate at 37°C for 30 minutes.
10. Empty the contents of the wells and wash by adding 400 µL of wash solution to every well. Repeat the wash 8 more times for a total of **9 washes**. After the final wash, empty or aspirate the wells, and firmly tap the plate on a lint free paper towel to remove any remaining wash buffer.
11. Add 100 µL of the Substrate Solution to each well.
12. Incubate for 30 minutes at room temperature in the dark.
13. Add 100 µL of Stop Solution to each well.
14. Blank the plate reader against the Blank wells, read the optical density at 450 nm, preferably with correction between 570 and 590 nm. If the plate reader is not able to be blanked against the Blank wells, manually subtract the mean optical density of the Blank wells from all readings.

Calculation of Results

Several options are available for the calculation of the concentration of rat GRO/CINC-1 in the samples. We recommend that the data be handled by an immunoassay software package utilizing a 4 parameter logistic curve fitting program. If data reduction software is not readily available, the concentration of rat GRO/CINC-1 can be calculated as follows:

1. Calculate the average net Optical Density (OD) bound for each standard and sample by subtracting the average Blank OD from the average OD for each standard and sample.

$$\text{Average Net OD} = \text{Average OD} - \text{Average Blank OD}$$

2. Plot the Average Net OD for each standard versus rat GRO/CINC-1 concentration in each standard. Approximate a straight line through the points. The concentration of rat GRO/CINC-1 in the unknowns can be determined by interpolation.

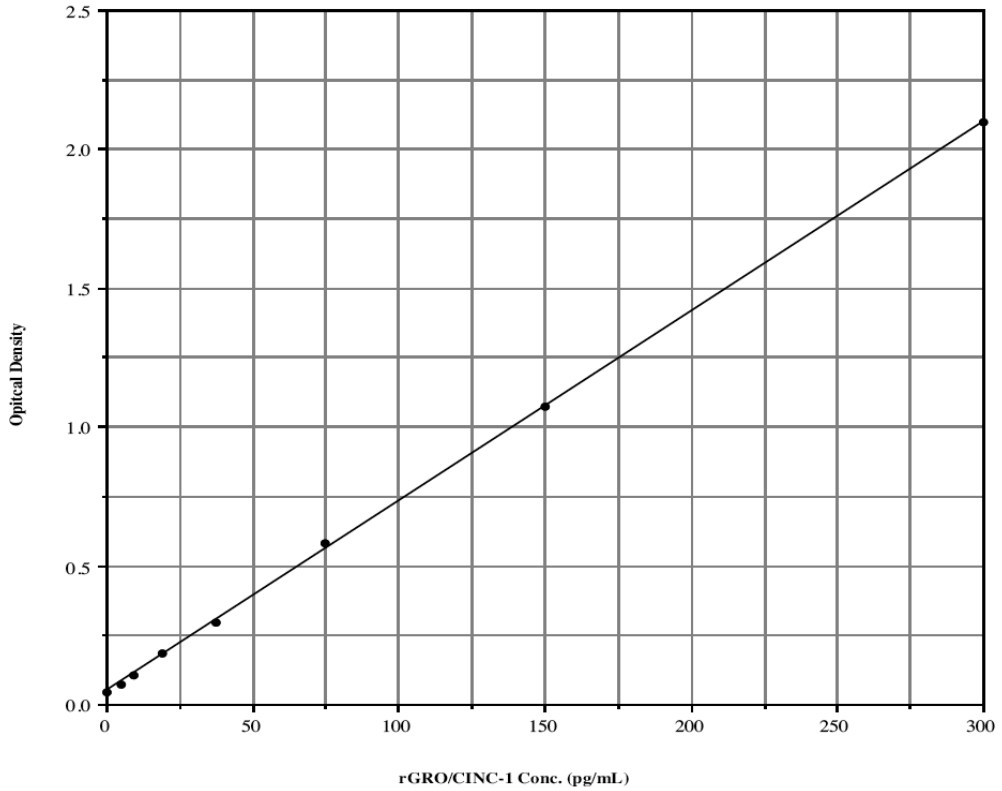
Typical Results

The results shown below are for illustration only and **should not** be used to calculate results from another assay.

<u>Sample</u>	<u>Average OD</u>	<u>Net OD</u>	rat GRO/CINC-1 (<u>pg/mL</u>)
Blank	(0.043)		
S0	0.087	0.044	0
S1	2.139	2.096	300
S2	1.117	1.074	150
S3	0.622	0.579	75
S4	0.338	0.295	37.5
S5	0.229	0.186	18.8
S6	0.152	0.109	9.38
S7	0.115	0.072	4.69

Typical Standard Curve

A typical standard curve is shown below. This curve **must not** be used to calculate rat GRO/CINC-1 concentrations; each user must run a standard curve for each assay.



Performance Characteristics

The following parameters for this kit were determined using the guidelines listed in the National Committee for Clinical Laboratory Standards (NCCLS) Evaluation Protocols⁸.

Sensitivity

Sensitivity was calculated by determining the average optical density bound for eight (8) wells run with 0 pg/mL Standard, and comparing to the average optical density for eight (8) wells run with Standard #7. The detection limit was determined as the concentration of rat GRO/CINC-1 measured at two (2) standard deviations from the 0 pg/mL Standard along the standard curve.

Average Optical Density for the S0 = 0.064 ± 0.007 (10.9%)

Average Optical Density for Standard #7 = 0.097 ± 0.007 (7.4%)

Delta Optical Density (4.69-0 pg/mL) = 0.033

2 SD's of the 0 pg/mL Standard = 2 x 0.007 = 0.014

Sensitivity = $\frac{0.014}{0.033} \times 4.69 \text{ pg/mL} = \mathbf{1.99 \text{ pg/mL}}$

Linearity

A sample containing 600 pg/mL rat GRO/CINC-1 was diluted 4 times 1:2 into Assay Buffer and measured in the assay. The data was plotted graphically as actual rat GRO/CINC-1 concentration versus measured rat GRO/CINC-1 concentration.

The line obtained had a slope of 1.085 and a correlation coefficient of 0.999.

Precision

Intra-assay precision was determined by taking samples containing low, medium and high concentrations of rat GRO/CINC-1 and running these samples multiple times (n=44) in the same assay. Inter-assay precision was determined by measuring three samples with low, medium and high concentrations of rat GRO/CINC-1 in multiple assays (n=24).

The precision numbers listed below represent the percent coefficient of variation for the concentrations of rat GRO/CINC-1 determined in these assays as calculated by a 4 parameter logistic curve fitting program.

	rat GRO/CINC-1 (pg/mL)	Intra-assay %CV	Inter-assay %CV
Low	7.75	7.6	
Medium	34.23	5.3	
High	136.46	4.4	
Low	8.17		5.1
Medium	35.34		4.7
High	133.20		2.9

Cross Reactivities

The cross reactivities for a number of related compounds was determined by dissolving the cross reactant in Assay Buffer. These samples were then measured in the rat GRO/CINC-1 assay, and the measured rat GRO/CINC-1 concentration calculated. The % cross reactivity was calculated by comparison with the actual concentration of cross reactant in the sample and expressed as a percentage.

<u>Compound</u>	<u>Cross Reactivity</u>
Rat GRO/CINC-1	100%
Rat GRO/CINC-2 α	<0.1%
Rat GRO/CINC-2 β	<0.1%
Rat GRO/CINC-3	<0.1%
Rat MCP-1	<0.1%
Rat Rantes	<0.1%
Rat MIP-1	<0.1%
Rat IL-1 β	<0.1%
Mouse GRO β /MIP-2	<0.1%

Sample Recoveries

Please refer to pages 4 and 5 for Sample Handling recommendations and Standard preparation.

Rat GRO/CINC-1 concentrations were measured in tissue culture media, rat serum and rat EDTA plasma. Rat GRO/CINC-1 was spiked into the undiluted samples of these media which were then diluted with the kit Assay Buffer and assayed in the kit. The following results were obtained:

<u>Sample</u>	<u>% Recovery*</u>	<u>Recommended Dilution*</u>
Tissue Culture Media	92.3	≥1:2
rat Serum	91.6	≥1:2
rat EDTA Plasma	96.6	≥1:4

* See Sample Handling instructions on page 4 for details.

References

1. T. Chard, "An Introduction to Radioimmunoassay & Related Techniques, 4th Ed.", (1990) Amsterdam: Elsevier.
2. P. Tijssen, "Practice & Theory of Enzyme Immunoassays", (1985) Amsterdam: Elsevier.
3. K. Watanabe, et al., J. Biol. Chem., (1989) 264: 19559-19563.
4. Y. Yamasaki, et al., Brain Research, (1997) 759: 103-111.
5. S. Hashimoto, et al., Am. J. Physiol., (1996) 270: L819-L828.
6. T. Tsuruma, et al., Transplant Proc., (1998) 30: 2644-2645.
7. H. Makita, et al., Am. J. Respir. Crit. Care Med., (1998) 158: 573-579.
8. National Committee for Clinical Laboratory Standards Evaluation Protocols, SC1, (1989) Villanova, PA: NCCLS.



USE FOR RESEARCH PURPOSES ONLY

Unless otherwise specified expressly on the packaging, all products sold hereunder are intended for and may be used for research purposes only and may not be used for food, drug, cosmetic or household use or for the diagnosis or treatment of human beings. Purchase does not include any right or license to use, develop or otherwise exploit these products commercially. Any commercial use, development or exploitation of these products or development using these products without the express written authorization of Enzo Life Sciences, Inc. is strictly prohibited. Buyer assumes all risk and liability for the use and/or results obtained by the use of the products covered by this invoice whether used singularly or in combination with other products.

LIMITED WARRANTY; DISCLAIMER OF WARRANTIES

These products are offered under a limited warranty. The products are guaranteed to meet all appropriate specifications described in the package insert at the time of shipment. Enzo Life Sciences' sole obligation is to replace the product to the extent of the purchasing price. All claims must be made to Enzo Life Sciences, Inc., within five (5) days of receipt of order. THIS WARRANTY IS EXPRESSLY IN LIEU OF ANY OTHER WARRANTIES OR LIABILITIES, EXPRESS OR IMPLIED, INCLUDING WARRANTIES OF MERCHANTABILITY, FITNESS FOR A PARTICULAR PURPOSE, AND NONINFRINGEMENT OF THE PATENT OR OTHER INTELLECTUAL PROPERTY RIGHTS OF OTHERS, AND ALL SUCH WARRANTIES (AND ANY OTHER WARRANTIES IMPLIED BY LAW) ARE EXPRESSLY DISCLAIMED.

TRADEMARKS AND PATENTS

Several Enzo Life Sciences products and product applications are covered by US and foreign patents and patents pending.

Global Headquarters

Enzo Life Sciences Inc.

10 Executive Blvd

Farmingdale, NY 11735

(p) 1-800-942-0430

(f) 1-631-694-7501

(e) info-usa@enzolifesciences.com

Enzo Life Sciences (ELS) AG

Industriestrasse 17, Postfach

CH-4415 Lause / Switzerland

(p) +41/0 61 926 89 89

(f) +41/0 61 926 89 79

(e) info-ch@enzolifesciences.com

