



## AlphaLISA<sup>®</sup> CXCL1/GRO-alpha Detection Kit

**Product number:** AL349C/F

Research Use Only. Not for use in diagnostic procedures.

### Product Information

- Application:** This kit is designed for the quantitative determination of human CXCL1/GRO- $\alpha$  in serum, buffered solution and cell culture supernatants using a homogeneous AlphaLISA assay (no wash steps).
- Sensitivity:** Lower Detection Limit (LDL): 0.5 pg/mL  
Lower Limit of Quantification (LLOQ): 1.8 pg/mL  
EC<sub>50</sub>: 4.4 ng/mL
- Dynamic range:** 0.5 – 30 000 pg/mL (Figure 1).

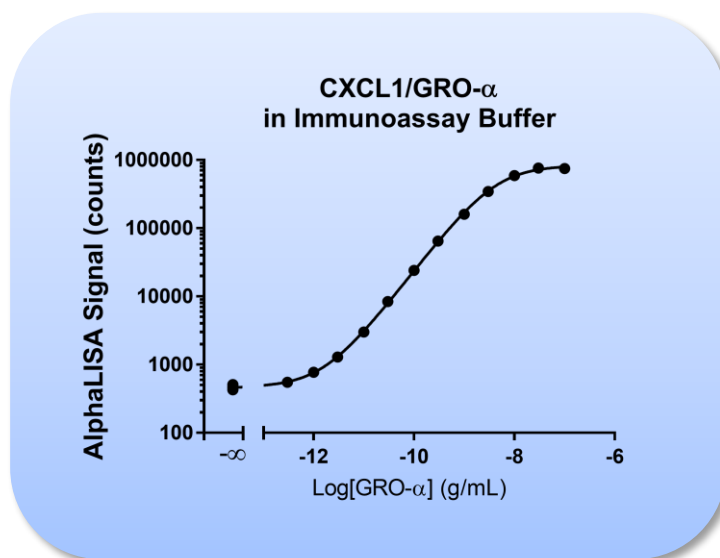


Figure 1. Typical sensitivity curves in AlphaLISA Immunoassay Buffer. The data was generated using a white Optiplate<sup>™</sup>384 microplate and the EnVision<sup>®</sup> Multilabel Plate Reader 2103 with Alpha option.

- Storage:** Store kit in the dark at +4°C. Store reconstituted analyte at -20°C.
- Stability:** This kit is stable for at least 12 months from the manufacturing date when stored in its original packaging and the recommended storage conditions.

## Analyte of Interest

C-X-C Motif Chemokine 1 (CXCL1) is an 11 kDa chemokine and is part of the intercrine alpha (chemokine CXC) family. The secreted protein is proteolytically processed at the N-terminus and the processed form is usually referred to as GRO- $\alpha$ . CXCL1 is expressed by macrophages, neutrophils and epithelial cells. Its synthesis is induced by a variety of inflammatory mediators such as PDGF, M-CSF, TNF $\alpha$ , LPS, and TLR 4. The major role of CXCL1 is its chemoattractant activity on neutrophils. CXCL1 plays a role in inflammation and exerts its effects on endothelial cells in an autocrine way. CXCL1 is implicated in spinal cord development, angiogenesis, inflammation, wound healing, and tumorigenesis. CXCL1 is overexpressed in invasive bladder cancer and is secreted by human melanoma cells

## Description of the AlphaLISA Assay

AlphaLISA technology allows the detection of molecules of interest in buffer, cell culture media, serum and plasma in a highly sensitive, quantitative, reproducible and user-friendly mode. In an AlphaLISA assay, a biotinylated anti-analyte antibody binds to the Streptavidin-coated Alpha Donor beads, while another anti-analyte antibody is conjugated to AlphaLISA Acceptor beads. In the presence of the analyte, the beads come into close proximity. The excitation of the Donor beads provokes the release of singlet oxygen molecules that triggers a cascade of energy transfer in the Acceptor beads, resulting in a sharp peak of light emission at 615 nm (Figure 2).

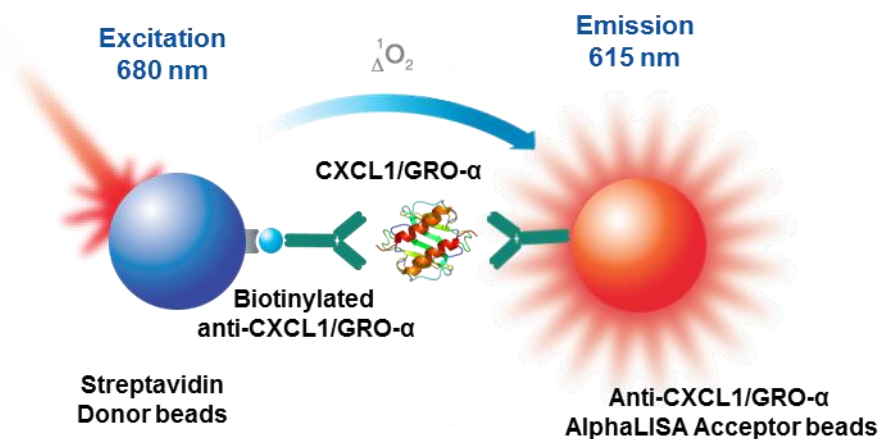


Figure 2. AlphaLISA Assay Principle.

## Precautions

- The Alpha Donor beads are light-sensitive. All the other assay reagents can be used under normal light conditions. All Alpha assays using the Donor beads should be performed under subdued laboratory lighting (< 100 lux). Green filters (LEE 090 filters (preferred) or Roscolux filters #389 from Rosco) can be applied to light fixtures.
- All blood components and biological materials should be handled as potentially hazardous.
- Some analytes are present in blood. Take precautionary measures to avoid contamination of the reagent solutions.
- The Biotinylated Anti-Analyte Antibody contains sodium azide. Contact with skin or inhalation should be avoided.

## Kit Contents

Kit components	AL349C (500 assay points <sup>***</sup> )	AL349F (5000 assay points <sup>***</sup> )
AlphaLISA Anti-CXCL1/GRO- $\alpha$ Acceptor beads stored in PBS, 0.05% Kathon, pH 7.2	50 $\mu$ L @ 5 mg/mL (1 brown tube, <u>white</u> cap)	500 $\mu$ L @ 5 mg/mL (1 brown tube, <u>white</u> cap)
Streptavidin (SA)-coated Donor beads stored in 25 mM HEPES, 100 mM NaCl, 0.05% Kathon, pH 7.4	200 $\mu$ L @ 5 mg/mL (1 brown tube, <u>black</u> cap)	2 X 1 mL @ 5 mg/mL (2 brown tubes, <u>black</u> caps)
Biotinylated Anti-CXCL1/GRO- $\alpha$ Antibody stored in PBS, 0.1% Tween-20, 0.05% NaN <sub>3</sub> , pH 7.4	50 $\mu$ L @ 500 nM (1 tube, <u>black</u> cap)	500 $\mu$ L @ 500 nM (1 tube, <u>black</u> cap)
Human CXCL1/GRO- $\alpha$ , analyte* lyophilized	0.1 $\mu$ g 1 tube, clear cap	0.1 $\mu$ g 1 tube, clear cap
AlphaLISA Immunoassay Buffer (10X) **	10 mL, 1 small bottle	100 mL, 1 large bottle

\* Reconstitute CXCL1/GRO- $\alpha$  in 100  $\mu$ L Milli-Q<sup>®</sup> grade H<sub>2</sub>O. The reconstituted analyte should be used within 60 minutes, if possible, or aliquoted into screw-capped polypropylene vials and stored at -20 °C for further experiments. Avoid repeated freezing and thawing. It has been demonstrated that reconstituted human CXCL1 / GRO- $\alpha$  is stable for at least 75 days at -20°C. One vial contains an amount of CXCL1/GRO- $\alpha$  sufficient for performing 10 standard curves. Additional vials can be ordered separately (cat # AL349S).

\*\* Contains 250 mM HEPES, pH 7.4, 1% Casein, 10 mg/mL Dextran-500, 5% Triton X-100 and 0.5% Kathon. Extra buffer can be ordered separately (cat # AL000C: 10 mL, cat # AL000F: 100 mL).  
Note: 10X buffer might be slightly yellow. However, this does not affect the assay results.

\*\*\* The number of assay points is based on an assay volume of 100  $\mu$ L in 96-well plates or 50  $\mu$ L in 384-well assay plates using the kit components at the recommended concentrations.

Sodium azide should **not** be added to the stock reagents. High concentrations of sodium azide (> 0.001 % final in the assay) might decrease the AlphaLISA signal.

### Specific additional required reagents and materials:

The following materials are recommended:

Item	Suggested source	Catalog #
TopSeal™-A Adhesive Sealing Film	Revvity Inc.	6050185
EnVision <sup>®</sup> -Alpha Reader	Revvity Inc.	-

## Recommendations

- The volume indicated on each tube is guaranteed for single pipetting. Multiple pipetting of the reagents may reduce the theoretical amount left in the tube. To minimize loss when pipetting beads, it is preferable not to pre-wet the tip.
- Centrifuge all tubes (including lyophilized analyte) before use to improve recovery of content (2000g, 10-15 sec). Re-suspend all reagents by vortexing before use.
- Use Milli-Q® grade H<sub>2</sub>O (18 MΩ•cm) to dilute 10X AlphaLISA Immunoassay Buffer to reconstitute the lyophilized analyte.
- When diluting the standard or samples, change tips between each standard or sample dilution. When loading reagents in the assay microplate, change tips between each standard or sample addition and after each set of reagents.
- When reagents are added to the microplate, make sure the liquids are at the bottom of the well.
- Small volumes may be prone to evaporation. It is recommended to cover microplates with TopSeal-A Adhesive Sealing Films to reduce evaporation during incubation. Microplates can be read with the TopSeal-A Film.
- The AlphaLISA signal is detected with an EnVision Multilabel Reader equipped with the Alpha option using the AlphaScreen standard settings (e.g. Total Measurement Time: 550 ms, Laser 680 nm Excitation Time: 180 ms, Mirror: D640as, Emission Filter: M570w, Center Wavelength 570 nm, Bandwidth 100 nm, Transmittance 75%).
- AlphaLISA signal will vary with temperature and incubation time. For consistent results, identical incubation times and temperature should be used for each plate.
- The standard curves shown in this technical data sheet are provided for information only. A standard curve must be generated for each experiment. The standard curve should be performed in the Immunoassay buffer for serum and/or plasma samples.

## Assay Procedure

IMPORTANT: PLEASE READ THE RECOMMENDATIONS BELOW BEFORE USE

- The manual described below is an example for generating one standard curve in a 50 µL final assay volume (48 wells, triplicate determinations). The manuals also include testing samples in 452 wells. If a different amount of samples are tested, the volumes of all reagents have to be adjusted accordingly, as shown in the table below. These calculations do not include excess reagent to account for losses during transfer of solutions or dead volumes.
- The standard dilution manual is provided for information only. As needed, the number of replicates or the range of concentrations covered can be modified.
- Use of four background points in triplicate (12 wells) is recommended when LDL/LLOQ is calculated. One background point in triplicate (3 wells) can be used when LDL/LLOQ is not calculated.

Format	# of data points	Volume					Plate recommendation
		Final	Sample	AlphaLISA beads	Biotin Antibody	Donor beads	
AL349C	250	100 µL	10 µL	10 µL	10 µL	70 µL	White OptiPlate-96 (cat # 6005290) White ½ AreaPlate-96 (cat # 6005560)
	500	50 µL	5 µL	5 µL	5 µL	35 µL	White ½ AreaPlate-96 (cat # 6005560) White OptiPlate-384 (cat # 6007290) Light gray AlphaPlate™-384 (cat # 6005350)
	1 250	20 µL	2 µL	2 µL	2 µL	14 µL	Light gray AlphaPlate-384 (cat # 6005350) ProxiPlate™-384 Plus (cat # 6008280) White OptiPlate-384 (cat # 6007290)
	2 500	10 µL	1 µL	1 µL	1 µL	7 µL	Light gray AlphaPlate-1536 (cat # 6004350)
AL349F	5 000	50 µL	5 µL	5 µL	5 µL	35 µL	White ½ AreaPlate-96 (cat # 6005560) White OptiPlate-384 (cat # 6007290) Light gray AlphaPlate-384 (cat # 6005350)
	12 500	20 µL	2 µL	2 µL	2 µL	14 µL	Light gray AlphaPlate-384 (cat # 6005350) ProxiPlate-384 Plus (cat # 6008280) White OptiPlate-384 (cat # 6007290)
	25 000	10 µL	1 µL	1 µL	1 µL	7 µL	Light gray AlphaPlate-1536 (cat # 6004350)

The manual (3 incubation steps) described below is for 500 assay points including one standard curve (48 wells) and samples (452 wells).

If a different amount of samples are tested, the volumes of all reagents have to be adjusted accordingly.

1) Preparation of 1X AlphaLISA Immunoassay Buffer:

Add 3 mL of 10X AlphaLISA Immunoassay Buffer to 27 mL H<sub>2</sub>O.

2) Preparation of Anti-CXCL1/GRO- $\alpha$  analyte standard dilutions:

- a. Reconstitute lyophilized CXCL1/GRO- $\alpha$  (0.1  $\mu$ g) in 100  $\mu$ L H<sub>2</sub>O.
- b. Prepare standard dilutions as follows in 1X AlphaLISA Immunoassay Buffer (change tip between each standard dilution):

Tube	Vol. of CXCL1/GRO- $\alpha$ ( $\mu$ L)	Vol. of diluent ( $\mu$ L) *	[CXCL1/GRO- $\alpha$ ] in standard curve	
			(g/mL in 5 $\mu$ L)	(pg/mL in 5 $\mu$ L)
A	10 $\mu$ L of reconstituted CXCL1/GRO- $\alpha$	90	1.00E-07	100 000
B	60 $\mu$ L of tube A	140	3.00E-08	30 000
C	60 $\mu$ L of tube B	120	1.00E-08	10 000
D	60 $\mu$ L of tube C	140	3.00E-09	3 000
E	60 $\mu$ L of tube D	120	1.00E-09	1 000
F	60 $\mu$ L of tube E	140	3.00E-10	300
G	60 $\mu$ L of tube F	120	1.00E-10	100
H	60 $\mu$ L of tube G	140	3.00E-11	30
I	60 $\mu$ L of tube H	120	1.00E-11	10
J	60 $\mu$ L of tube I	140	3.00E-12	3
K	60 $\mu$ L of tube J	120	1.00E-12	1
L	60 $\mu$ L of tube K	140	3.00E-13	0.3
M ** (background)	0	100	0	0
N ** (background)	0	100	0	0
O ** (background)	0	100	0	0
P ** (background)	0	100	0	0

\* Dilute standards in diluent (e.g. 1X AlphaLISA Immunoassay Buffer).  
At low concentrations of analyte, a significant amount of analyte can bind to the vial. Therefore, load the analyte standard dilutions in the assay microplate within 60 minutes of preparation.

\*\* Four background points in triplicate (12 wells) are used when LDL is calculated. If LDL does not need to be calculated, one background point in triplicate can be used (3 wells).

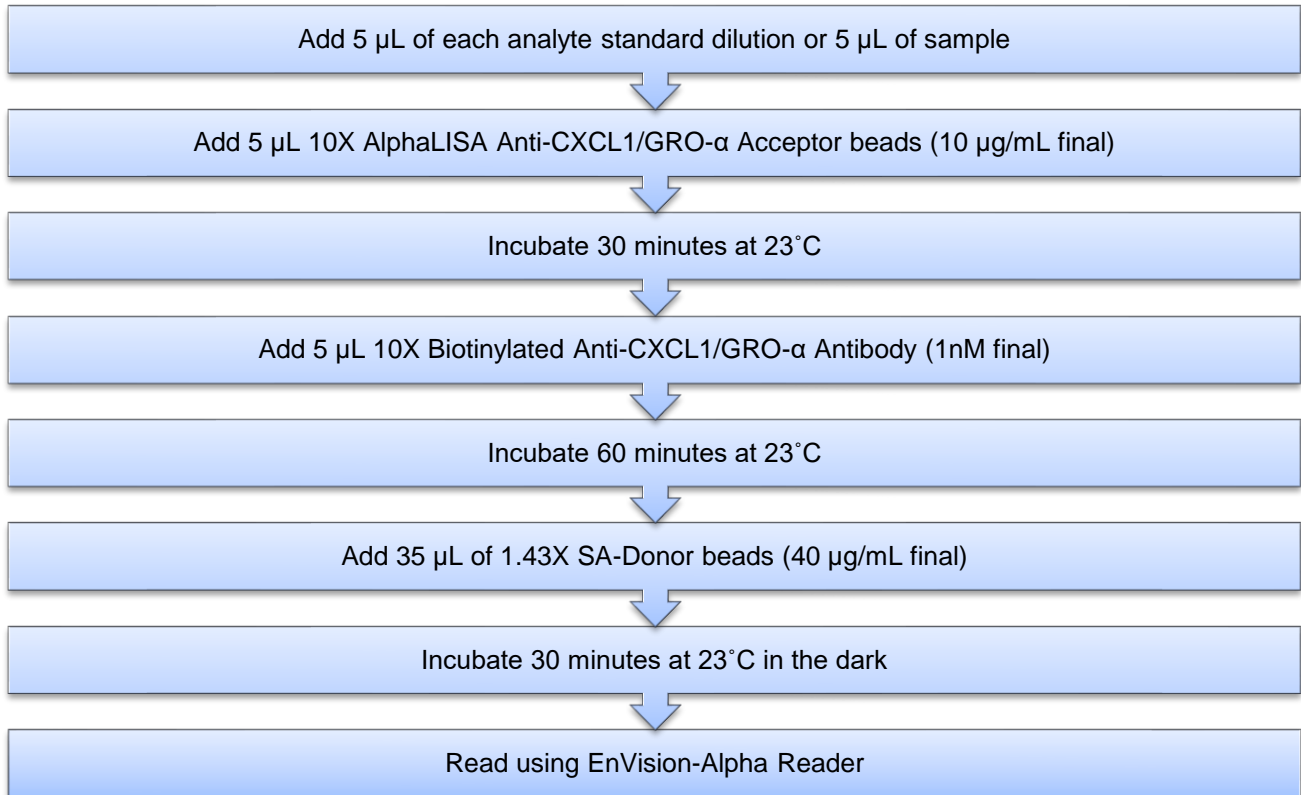
3) Preparation of 10X AlphaLISA Anti-CXCL1/GRO- $\alpha$  antibody Acceptor beads (100  $\mu$ g/mL):

- a. Prepare just before use.
- b. Add 50  $\mu$ L of 5 mg/mL AlphaLISA Anti-CXCL1/GRO- $\alpha$  antibody Acceptor beads to 2450  $\mu$ L of 1X AlphaLISA Immunoassay Buffer.

4) Preparation of 10X Biotinylated Anti- CXCL1/GRO- $\alpha$  antibody (10 nM):

- a. Prepare just before use.
- b. Add 50  $\mu$ L of 500 nM Biotinylated Anti- CXCL1/GRO- $\alpha$  antibody to 2450  $\mu$ L of 1X AlphaLISA Immunoassay Buffer.

- 5) Preparation of 1.43X Streptavidin (SA) Donor beads (57.2 µg/mL): Keep the beads under subdued laboratory lighting.
- Prepare just before use
  - Add 200 µL of 5 mg/mL SA-Donor beads to 17 300 µL of 1X AlphaLISA Immunoassay Buffer.
- 6) In a white Optiplate (384 wells):



## Data Analysis

- Calculate the average count value for the background wells.
- Generate a standard curve by plotting the AlphaLISA counts versus the concentration of analyte. A log scale can be used for either or both axes. No additional data transformation is required.
- Analyze data according to a nonlinear regression using the 4-parameter logistic equation (sigmoidal dose-response curve with variable slope) and a  $1/Y^2$  data weighting (the values at maximal concentrations of analyte after the hook point should be removed for correct analysis).
- The LDL is calculated by interpolating the average background counts (12 wells without analyte) + 3 x standard deviation value (average background counts + (3xSD)) on the standard curve.
- The LLOQ as measured here is calculated by interpolating the average background counts (12 wells without analyte) + 10 x standard deviation value (average background counts + (10xSD)) on the standard curve. Alternatively, the true LLOQ can be determined by spiking known concentrations of analyte in the matrix and measuring the percent recovery, and then determining the minimal amount of spiked analyte that can be quantified within a given limit (usually +/- 20% or 30% of the real concentration).
- Read from the standard curve the concentration of analyte contained in the samples.
- If samples have been diluted, the concentration read from the standard curve must be multiplied by the dilution factor.

## Assay Performance Characteristics

AlphaLISA assay performance described below was determined using the 3 step manual.

- Assay Sensitivity:

The LDL was calculated as described above. The values correspond to the lowest concentration of analyte that can be detected in a volume of 5  $\mu$ L using the recommended assay conditions.

LDL (pg/mL)	Buffer/Media	# of experiments
0.5	AlphaLISA Immunoassay Buffer	9
1.6	Fetal Bovine Serum	6
0.7	DMEM with 10% FBS	6
0.9	RPMI with 10% FBS	6

\*Note that LDL can be decreased (i.e. sensitivity increased) by increasing the volume of analyte in the assay (e.g. use 10  $\mu$ L of analyte in a final assay volume of 50  $\mu$ L).

- Assay Precision:

The following assay precision data were calculated from the three independent assays using two different kit lots. In each lot, the analytes were prepared in AlphaLISA Immunoassay Buffer, or RPMI. Each assay consisted of one standard curve comprising 12 data points (each in triplicate) and 12 background wells (no analytes). The assays were performed in 384-well format using AlphaLISA Immunoassay Buffer.

- Intra-assay precision:

The intra-assay precision was determined using a total of 7 independent determinations in triplicate, shown as CV%.

Human CXCL1/GRO- $\alpha$	Immunoassay Buffer	DMEM with 10% FBS	RPMI with 10% FBS	FBS
CV%	6	6	7	7

- Inter-assay precision:

The inter-assay precision was determined using a total of 7 independent determinations with 21 measurements. Shown as CV%.

CXCL1/GRO- $\alpha$	Immunoassay Buffer	DMEM with 10% FBS	RPMI with 10% FBS	FBS
CV%	10	9	11	11



- Spike Recovery:

Four known concentrations of analyte were spiked in Immunoassay Buffer and cell culture media containing 10% FBS. All samples, including non-spiked Immunoassay Buffers and culture media were measured in the assay. The average recovery from three independent measurements is reported.

Spiked Human CXCL1/GRO- $\alpha$ (ng/mL)	% Recovery			
	Immunoassay Buffer	DMEM with 10% FBS	RPMI with 10% FBS	FBS
3	94	94	104	106
1	95	88	95	93
0.3	89	89	102	104

- Specificity:

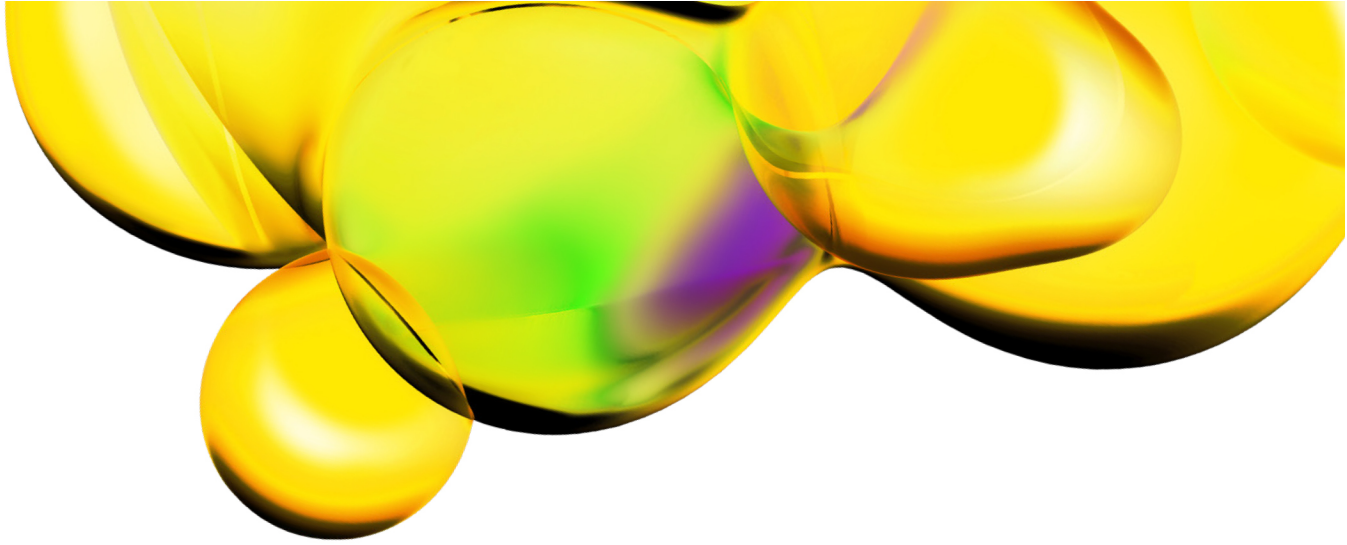
Cross-reactivity of the CXCL1/GRO- $\alpha$  AlphaLISA Kit was tested using the following proteins at 30 ng/mL in AlphaLISA Immunoassay Buffer.

Protein	% Cross-reactivity
Mouse CXCL1/GRO- $\alpha$	<0.1
Rat CXCL1/GRO- $\alpha$	<0.1
Human CXCL2/GRO- $\beta$	19
Human CXCL3/GRO- $\gamma$	33
Human CXCL9/MIG	<0.1

## Troubleshooting Guide

You will find detailed recommendations for common situations you might encounter with your AlphaLISA Assay kit at: [www.revivity.com](http://www.revivity.com)

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



The information provided in this document is for reference purposes only and may not be all-inclusive. Revvity, Inc., its subsidiaries, and/or affiliates (collectively, "Revvity") do not assume liability for the accuracy or completeness of the information contained herein. Users should exercise caution when handling materials as they may present unknown hazards. Revvity shall not be liable for any damages or losses resulting from handling or contact with the product, as Revvity cannot control actual methods, volumes, or conditions of use. Users are responsible for ensuring the product's suitability for their specific application. REVVITY EXPRESSLY DISCLAIMS ALL WARRANTIES, INCLUDING WARRANTIES OF MERCHANTABILITY OR FITNESS FOR A PARTICULAR PURPOSE, REGARDLESS OF WHETHER ORAL OR WRITTEN, EXPRESS OR IMPLIED, ALLEGEDLY ARISING FROM ANY USAGE OF ANY TRADE OR ANY COURSE OF DEALING, IN CONNECTION WITH THE USE OF INFORMATION CONTAINED HEREIN OR THE PRODUCT ITSELF

[www.revvity.com](http://www.revvity.com)

revvity

**Revvity, Inc.**  
940 Winter Street  
Waltham, MA 02451 USA  
[www.revvity.com](http://www.revvity.com)

For a complete listing of our global offices, visit [www.revvity.com](http://www.revvity.com)  
Copyright ©2023, Revvity, Inc. All rights reserved.