

MSD[®] MULTI-SPOT Assay System

Human I-309 Kit

1-Plate Kit	K151NXD-1
5-Plate Kit	K151NXD-2
25-Plate Kit	K151NXD-4



MSD Biomarker Assays

Human I-309 Kit

This package insert must be read in its entirety before using this product.

FOR RESEARCH USE ONLY.

NOT FOR USE IN DIAGNOSTIC PROCEDURES.

MESO SCALE DISCOVERY[®]

A division of Meso Scale Diagnostics, LLC.

1601 Research Boulevard

Rockville, MD 20850-3173 USA

www.mesoscale.com

MESO SCALE DISCOVERY, MESO SCALE DIAGNOSTICS, DISCOVERY WORKBENCH, MULTI-ARRAY, MULTI-SPOT, QUICKPLEX, SECTOR PR, SECTOR, SECTOR HTS, SULFO-TAG, www.mesoscale.com, SMALL SPOT (design), 96 WELL 1, 4, 7, & 10-SPOT (designs), 384 WELL 1 & 4-SPOT (designs), MSD, MSD (design), and SPOT THE DIFFERENCE are trademarks and/or service marks of Meso Scale Diagnostics, LLC.

© 2012 Meso Scale Diagnostics, LLC. All rights reserved.

Table of Contents

Introduction	4
Principle of the Assay	4
Reagents Supplied	5
Required Material and Equipment (not supplied)	5
Safety	5
Reagent Preparation	6
Protocol	8
Curve Fitting.....	8
Typical Data.....	9
Sensitivity	9
Assay Components	9
References.....	10
Summary Protocol	11
Plate Diagrams	13

Ordering Information

MSD Customer Service

Phone: 1-301-947-2085
Fax: 1-301-990-2776
Email: CustomerService@mesoscale.com

MSD Scientific Support

Phone: 1-301-947-2025
Fax: 1-240-632-2219 attn: Scientific Support
Email: ScientificSupport@mesoscale.com

Introduction

Chemokines are the largest cytokine family; they exert effects by binding to G-protein-coupled receptors. I-309/CCL1 is a human chemokine which is structurally related to other pro-inflammatory cytokines sharing three conserved C-terminal basic amino acids.¹ I-309 is a selective ligand for the CCR8 receptor; receptor binding modulates recruitment of monocytes, natural killer cells, immature B cells, and dendritic cells.²⁻⁵ In addition to pro-inflammatory functions, I-309 modulates apoptosis and angiogenesis⁶ and protects murine thymic lymphoma cell lines against dexamethasone-induced apoptosis.²

Principle of the Assay

MSD biomarker assays provide a rapid and convenient method for measuring the levels of protein targets within a single, small-volume sample. Human I-309 is a sandwich immunoassay (Figure 1). MSD provides a plate pre-coated with capture antibodies. The user adds the sample and a solution containing detection antibodies conjugated with electrochemiluminescent labels (MSD SULFO-TAG™) over the course of one or more incubation periods. Analytes in the sample bind to capture antibodies immobilized on the working electrode surface; recruitment of the detection antibodies by the bound analytes completes the sandwich. The user adds an MSD buffer that provides the appropriate chemical environment for electrochemiluminescence and loads the plate into a SECTOR® Imager where a voltage applied to the plate electrodes causes the captured labels to emit light. The instrument measures the intensity of emitted light to provide a quantitative measure of analytes in the sample.

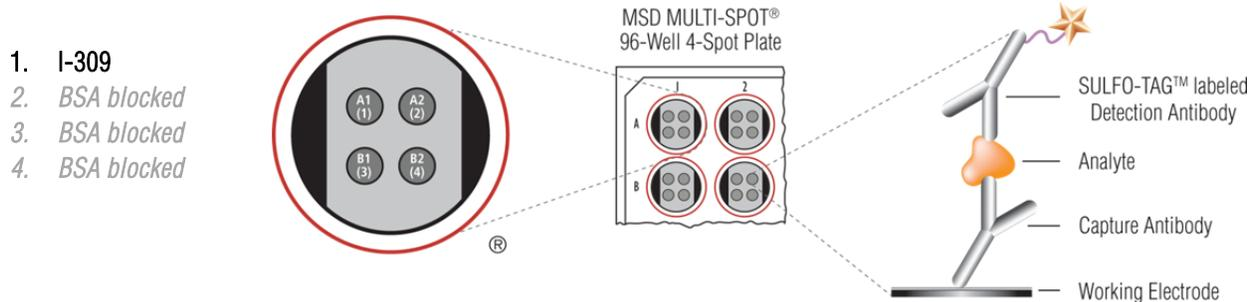


Figure 1. Spot diagram showing placement of analyte capture antibodies. The numbering convention for the different spots is maintained in the software visualization tools, on the plate packaging, and in the data files. A unique bar code label on each plate allows complete traceability back to MSD manufacturing records.

Reagents Supplied

Product Description	Storage	Quantity per Kit		
		K151NXD-1	K151NXD-2	K151NXD-4
MULTI-SPOT 96-Well, 4-Spot Human I-309 Plate N451NXA	2–8°C	1 plate	5 plates	25 plates
SULFO-TAG Anti-hu I-309 Antibody ¹ (50X)	2–8°C	1 vial (75 µL)	1 vial (375 µL)	5 vials (375 µL ea)
Human I-309 Calibrator (20X)	≤-70°C	1 vial (60 µL)	5 vials (60 µL ea)	25 vials (60 µL ea)
Diluent 12 R50JA-4 (10 mL), R50JA-3 (50 mL)	≤-10°C	1 bottle (10 mL)	1 bottle (50 mL)	5 bottles (50 mL ea)
Diluent 3 R51BA-4 (5 mL), R51BA-5 (25 mL)	≤-10°C	1 bottle (5 mL)	1 bottle (25 mL ea)	5 bottles (25 mL ea)
Blocker A Kit (Blocker A [dry] in 250 mL bottle and 50 mL bottle of 5X Phosphate Buffer) R93AA-2 (250 mL)	RT	1 kit (250 mL)	1 kit (250 mL)	5 kits (250 mL ea)
Read Buffer T (4X) R92TC-3 (50 mL)	RT	1 bottle (50 mL)	1 bottle (50 mL)	5 bottles (50 mL ea)

Required Material and Equipment (not supplied)

- Appropriately sized tubes for reagent preparation
- Microcentrifuge tubes for preparing serial dilutions
- Phosphate-buffered saline plus 0.05% Tween-20 (PBS-T) for plate washing
- Liquid handling equipment for desired throughput, capable of dispensing 10 to 150 µL/well into a 96-well microtiter plate
- Plate washing equipment: automated plate washer or multichannel pipette
- Adhesive plate seals
- Microtiter plate shaker
- Deionized water

Safety

Use safe laboratory practices and wear gloves, safety glasses, and lab coats when handling kit components. Handle and dispose of all hazardous samples properly in accordance with local, state, and federal guidelines.

Additional safety information is available in the product Material Safety Data Sheet, which can be obtained from MSD Customer Service.

¹ SULFO-TAG–conjugated detection antibodies should be stored in the dark.

Reagent Preparation

Bring all reagents to room temperature. Thaw the stock calibrator on ice.

Important: Upon first thaw, separate Diluent 3 and Diluent 12 into aliquots appropriate for the size of your needs before refreezing.

Prepare Blocker A Solution

Follow the Blocker A instructions included in the kit.

Prepare Standards

MSD supplies calibrator for the Human I-309 Kit at 20-fold higher concentration than the recommended highest standard. We recommend a 7-point standard curve with 3-fold serial dilution steps and a zero calibrator blank. Signals from the blank should be excluded when generating the curve. Thaw the stock calibrator and keep on ice, then add to diluent at room temperature to make the standard curve solutions.

Standard	Human I-309 (pg/mL)	Dilution Factor
Stock Calibrator	5000	
STD-01	250	20
STD-02	83.3	3
STD-03	27.8	3
STD-04	9.26	3
STD-05	3.09	3
STD-06	1.03	3
STD-07	0.343	3
STD-08	0	n/a

To prepare 7 standard solutions plus a zero calibrator blank for up to 3 replicates:

- 1) Prepare the highest standard by adding 15 μ L of stock calibrator to 285 μ L of Diluent 12. Mix well.
- 2) Prepare the next standard by transferring 100 μ L of the highest standard to 200 μ L of Diluent 12. Mix well. Repeat 3-fold serial dilutions 5 additional times to generate 7 standards.
- 3) Use Diluent 12 as the blank.

Dilute Samples

For human serum and plasma samples, MSD recommends 2-fold dilution in Diluent 12; however, you may need to adjust the dilution factor for the sample set under investigation. To dilute sample 2-fold, add 75 μ L of sample to 75 μ L of Diluent 12.

Prepare Detection Antibody Solution

MSD provides detection antibody as a 50X stock solution. The working detection antibody solution is 1X.

For 1 plate, combine:

- 60 μ L of 50X SULFO-TAG Anti-hu I-309 Antibody
- 2940 μ L of Diluent 3

Prepare Read Buffer

MSD provides Read Buffer T as a 4X stock solution. The working solution is 2X.

For 1 plate, combine:

- 10 mL of Read Buffer T (4X)
- 10 mL of deionized water

You may prepare diluted read buffer in advance and store it at room temperature in a tightly sealed container.

Prepare MSD Plate

MSD plates are pre-coated with capture antibodies (Figure 1) and exposed to a proprietary stabilizing treatment to ensure the integrity and stability of the immobilized antibodies. Plates can be used as delivered; no additional preparation (e.g., pre-wetting) is required.

Protocol

- Add Blocker A Solution:** Add 150 μL of Blocker A solution to each well. Seal the plate with an adhesive plate seal and incubate for 1 hour with vigorous shaking (300–1000 rpm) at room temperature.
- Wash and Add Sample:** Wash the plate 3 times with 300 μL /well of PBS-T. Add 50 μL of sample (standards, controls, or unknowns) per well. Seal the plate with an adhesive plate seal and incubate for 2 hours with vigorous shaking (300–1000 rpm) at room temperature. You may prepare detection antibody solution during incubation.
- Wash and Add Detection Antibody Solution:** Wash the plate 3 times with 300 μL /well of PBS-T. Add 25 μL of detection antibody solution to each well. Seal the plate with an adhesive plate seal and incubate for 1 hour with vigorous shaking (300–1000 rpm) at room temperature. You may prepare diluted read buffer during incubation.
- Wash and Read:** Wash the plate 3 times with 300 μL /well of PBS-T. Add 150 μL of 2X Read Buffer T to each well. Analyze the plate on the SECTOR Imager. No incubation in read buffer is required before reading the plate.

Notes

Shaking the plate typically accelerates capture at the working electrode.

You may keep excess diluted read buffer in a tightly sealed container at room temperature for later use.

Bubbles introduced when adding read buffer will interfere with imaging of the plate and produce unreliable data. Use reverse pipetting technique to avoid creating bubbles.

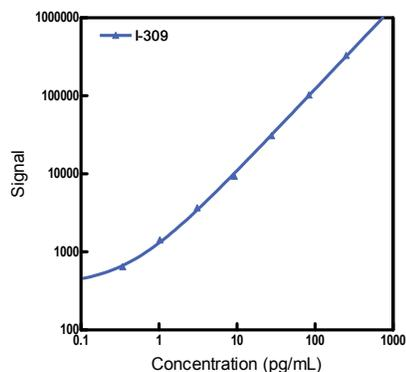
Due to the varying nature of each research application, you should assess assay stability before allowing plates to sit with read buffer for extended periods.

Curve Fitting

MSD DISCOVERY WORKBENCH[®] software uses least-squares fitting algorithms to generate the standard curve that will be used to calculate the concentration of analyte in the samples. The assays have a wide dynamic range (3–4 logs) that allows accurate quantification without the need for dilution in many cases. By default, the software uses a 4-parameter logistic model (or sigmoidal dose-response) and includes a $1/Y^2$ weighting function. The weighting function is important because it provides a better fit of data over a wide dynamic range, particularly at the low end of the standard curve.

Typical Data

The following standard curve graph illustrates the dynamic range of the assay. Actual signals will vary. Best quantification of unknown samples will be achieved by generating a standard curve for each plate using a minimum of 2 replicates of the standards.



I-309		
Conc. (pg/mL)	Average Signal	%CV
0	208	6.6
0.343	647	11.9
1.03	1424	6.5
3.09	3677	7.3
9.26	9314	3.6
27.8	30 691	6.0
83.3	102 529	3.8
250	330 631	7.1

Sensitivity

The lower limit of detection (LLOD) is a calculated concentration based on a signal 2.5 standard deviations above the background (zero calibrator blank). The LLOD shown below was calculated based on 4 runs.

I-309	
Average LLOD (pg/mL)	0.431
LLOD Range (pg/mL)	0.408–0.461

Assay Components

Calibrator

The assay calibrator uses recombinant I-309 protein, residues 24–96, expressed in *E. coli*.

Antibodies

Analyte	Source Species	
	MSD Capture Antibody	MSD Detection Antibody
I-309	Mouse monoclonal	Goat Polyclonal

References

1. Miller MD, Krangel MS. The human cytokine I-309 is a monocyte chemoattractant. *Proc Natl Acad Sci USA*. 1992 Apr 1;89(7):2950-4.
2. Zingoni A, et al. The chemokine receptor CCR8 is preferentially expressed in Th2 but not Th1 cells. *J Immunol*. 1998 Jul 15;161(2):547-51.
3. Tiffany HL, et al. Identification of CCR8: a human monocyte and thymus receptor for the CC chemokine I-309. *J Exp Med*. 1997 Jul 7;186(1):165-70.
4. Roos RS, et al. Identification of CCR8, the receptor for the human CC chemokine I-309. *J Biol Chem*. 1997 Jul 11;272(28):17251-4.
5. Goya I, et al. Identification of CCR8 as the specific receptor for the human beta-chemokine I-309 cloning and molecular characterization of murine CCR8 as the receptor for TCA-3. *J Immunol*. 1998 Feb 15;160(4):1975-81.
6. Denis C, et al. C-Terminal Clipping of Chemokine CCL1/I-309 Enhances CCR8-Mediated Intracellular Calcium Release and Anti-Apoptotic Activity. *PLoS One*. 2012;7(3):e34199. Epub 2012 Mar 27

Summary Protocol

Human I-309 Kit

*MSD provides this summary protocol for your convenience.
Please read the entire detailed protocol prior to performing
the Human I-309 assay.*

Sample and Reagent Preparation

Bring all reagents to room temperature and thaw the calibrator on ice.

Prepare Blocker A solution.

Prepare standard solutions using the supplied calibrator:

- Dilute the stock calibrator 20-fold in Diluent 12.
- Perform a series of 3-fold dilution steps and prepare a zero calibrator blank.

Dilute samples 2-fold in Diluent 12 before adding to the plate.

Prepare detection antibody solution by diluting stock detection antibody 50-fold in Diluent 3.

Prepare 2X Read Buffer T by diluting stock 4X Read Buffer T 2-fold with deionized water.

Step 1: Add Blocker A Solution

Add 150 μ L/well of Blocker A solution.

Incubate at room temperature with vigorous shaking (300–1000 rpm) for 1 hour.

Step 2: Wash and Add Sample

Wash plate 3 times with 300 μ L/well of PBS-T.

Add 50 μ L/well of sample (standards, controls, or unknowns).

Incubate at room temperature with vigorous shaking (300–1000 rpm) for 2 hours.

Step 3: Wash and Add Detection Antibody Solution

Wash plate 3 times with 300 μ L/well of PBS-T.

Add 25 μ L/well of 1X detection antibody solution.

Incubate at room temperature with vigorous shaking (300–1000 rpm) for 1 hour.

Step 4: Wash and Read Plate

Wash plate 3 times with 300 μ L/well of PBS-T.

Add 150 μ L/well of 2X Read Buffer T.

Analyze plate on SECTOR Imager.

Plate Diagrams

