



Human Adiponectin

www.mesoscale.com®

Ordering Information

MSD Customer Service
 Phone: 1-240-314-2795
 : 1-301-990-2776
 Email: CustomerService@mesoscale.com

Scientific Support

Phone: 1-240-314-2798
 Email: ScientificSupport@mesoscale.com

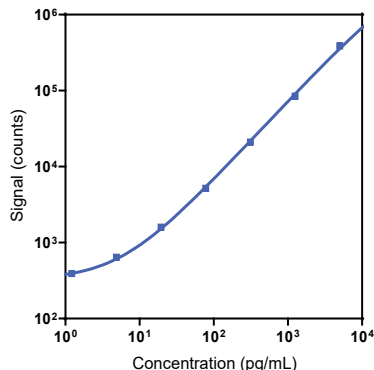
Company Address

MESO SCALE DISCOVERY®
 A division of
 Meso Scale Diagnostics, LLC.
 1601 Research Boulevard
 Rockville, MD 20850-3173 USA

| Product Options | Catalog Number | Description |
|-----------------|--|---|
| Multiplex | K151AGM, K251AGM | U-PLEX Biomarker Group 3 (human) |
| | K151R9K-1/-2/-4 | U-PLEX Human Adiponectin Assay with SECTOR™ plates |
| Singleplex | K151R9K-21/-22/-24 | U-PLEX Human Adiponectin Assay with QuickPlex® plates |
| | K251R9K-2/-4 | U-PLEX Human Adiponectin Assay with 384-well plates |
| Antibody Set | B21R9-2/-3 | U-PLEX Human Adiponectin Antibody Set |
| Protocol | U-PLEX Product Inserts are available at www.mesoscale.com . | |

The U-PLEX® platform was designed to provide ultimate flexibility for detection of biomarkers in a wide variety of sample types. This datasheet provides the representative performance of the U-PLEX Human Adiponectin Assay tested on U-PLEX 96-well SECTOR plates run as a multiplex. The data do not represent the product specifications. Under your experimental conditions, the assay may perform differently from the representative data. U-PLEX assays are offered in either singleplex or multiplex; both are available in 96- or 384-well plates. See a U-PLEX product insert for instrument compatibility.

Representative Calibration Curve and Sensitivity



| Assay | Median LLOD (pg/mL) | LLOD Range (pg/mL) |
|-------------|---------------------|--------------------|
| Adiponectin | 1.5 | 1.3-1.7 |

The Calibrator curve was fitted with a 4-parameter logistic model with a 1/Y² weighting. The lower limit of detection (LLOD) is a calculated concentration corresponding to the signal 2.5 standard deviations above the background (zero Calibrator).

Precision

| Control | Average Conc. (pg/mL) | Average intra-run Conc. (%CV) | Inter-run Conc. (%CV) |
|---------|-----------------------|-------------------------------|-----------------------|
| High | 179 | 4.1 | 9.0 |
| Mid | 148 | 3.4 | 8.7 |
| Low | 109 | 3.4 | 9.4 |

Controls were made by spiking Calibrator into assay diluent at 3 levels within the quantitative range of the assay. Average intra-run concentration %CV is the average %CV of the control replicates within an individual run. Inter-run concentration %CV is the variability of controls across multiple runs.

For Research Use Only.
 Not for use in diagnostic procedures.

MSD[®] U-PLEX Human Adiponectin

Tested Samples

| Sample Type | Serum (N=10) | EDTA Plasma (N=10) | Citrate Plasma (N=5) |
|----------------|--------------|--------------------|----------------------|
| Median (µg/mL) | 20 | 16 | 22 |
| Range (µg/mL) | 10–45 | 9.2–24 | 12–45 |
| % Detected | 100 | 100 | 100 |

Normal serum, EDTA plasma, and citrate plasma samples were diluted 200,000-fold prior to the assay.

Parallelism

| Serum | | | EDTA Plasma | | |
|---------------|--------------------|------------------|---------------|--------------------|------------------|
| Fold Dilution | Average % Recovery | % Recovery Range | Fold Dilution | Average % Recovery | % Recovery Range |
| 100,000 | 99 | 95–102 | 100,000 | 98 | 94–102 |
| 400,000 | 103 | 99–109 | 400,000 | 102 | 99–106 |
| 800,000 | 102 | 97–109 | 800,000 | 98 | 89–112 |

Normal human serum and EDTA plasma were tested at different dilutions. Percent recovery at each dilution level was normalized to the concentration of 200,000-fold diluted samples.

% Recovery = (measured concentration / expected concentration) x 100

Specificity

To assess specificity, the Adiponectin Antibody Set was tested individually against a larger panel of analytes for nonspecific binding (A2M, Adiponectin, ApoA1, ApoC3, CA1, Clusterin, Complement C9, Complement factor D, CRP, Cystatin C, DPPIV, Factor VII, ICAM-1, NGAL/LCN2, RBP4, SAA, Serpin A1, SHBG, sTfR-1, VCAM-1, vWF). Nonspecific binding was less than 0.5%.

% Nonspecificity = (nonspecific signal / specific signal) x 100

The Adiponectin capture antibody cross reacts with the CA1 calibrator 3.3%.

Diluent Compatibility

The data included in this document have been collected with Assay Diluent 12 and Antibody Diluent 11. MSD offers a range of assay and antibody diluents for separate purchase. Depending on your assay needs, other diluents may be tested.

Assay Components

Calibrator: Adiponectin is included in Calibrator 26. The human Adiponectin Calibrator is a full-length recombinant protein expressed in a mouse cell line.

Antibodies: The U-PLEX Human Adiponectin Assay uses a mouse monoclonal antibody for capture and a mouse monoclonal antibody for detection.

Assay generation: A

Note: This datasheet contains representative assay performance data. In custom multiplex formats, the assay may perform differently from the representative data shown.

Note: MSD recommends that samples be diluted 200,000-fold prior to analysis in this assay.

