



www.mesoscale.com

Ordering Information

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

Scientific Support

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

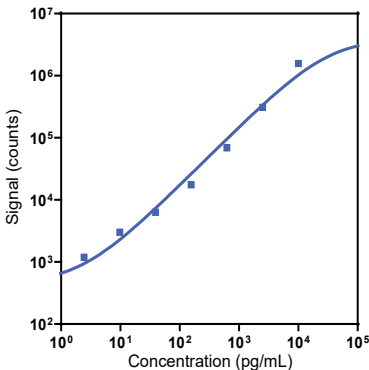
Company Address

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) |
| Singleplex | K151T9K-1/-2/-4 | U-PLEX Human ApoC3 Assay with SECTOR™ plates |
| | K151T9K-21/-22/-24 | U-PLEX Human ApoC3 Assay with QuickPlex Ultra™ plates |
| | K251T9K-2/-4 | U-PLEX Human ApoC3 Assay with 384-well plates |
| Antibody Set | B21T9-2/-3 | U-PLEX Human ApoC3 Antibody Set |
| Protocol | U-PLEX Product Inserts are available at www.mesoscale.com | |

The MESO SCALE DISCOVERY® 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 ApoC3 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) |
|-------|---------------------|--------------------|
| ApoC3 | 1.1 | 0.84-1.3 |

The Calibrator curve was fitted with a 4-parameter logistic model with a $1/Y^2$ 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 | 992 | 4.5 | 16.9 |
| Mid | 138 | 4.2 | 40.3 |
| Low | 23 | 4.9 | 55.8 |

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 ApoC3

Tested Samples

| Sample Type | Serum (N=10) | EDTA Plasma (N=10) | Citrate Plasma (N=5) |
|----------------|-----------------|-----------------------|-------------------------|
| Median (µg/mL) | 535 | 311 | 82 |
| Range (µg/mL) | 112-1,510 | 112-824 | 40-307 |
| % 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 | 94 | 86-102 | 100,000 | 92 | 85-95 |
| 400,000 | 107 | 102-111 | 400,000 | 99 | 93-106 |
| 800,000 | 106 | 97-112 | 800,000 | 102 | 94-110 |

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 ApoC3 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

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: ApoC3 is included in Calibrator 26. The human ApoC3 Calibrator is a native protein purified from human plasma.

Antibodies: The MESO SCALE DISCOVERY® U-PLEX Human ApoC3 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.

