

Mouse NGAL/LCN2



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

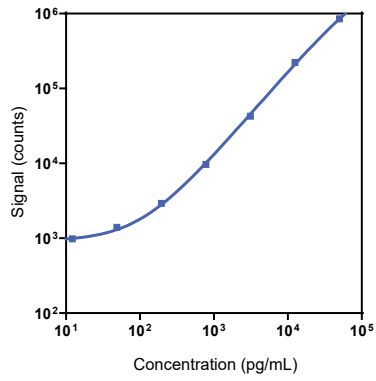
Company Address

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

Product Options	Catalog Number	Description
Multiplex	K15069M, K25069M	U-PLEX Biomarker Group 1 (mouse)
Singleplex	K152Z1K-1/-2/-4	U-PLEX Mouse NGAL/LCN2 Assay with SECTOR [™] plates
	K152Z1K-21/-22/-24	U-PLEX Mouse NGAL/LCN2 Assay with QuickPlex [®] plates
	K252Z1K-2/-4	U-PLEX Mouse NGAL/LCN2 Assay with 384-well plates
Antibody Set	B22Z1-2/-3	U-PLEX Mouse NGAL/LCN2 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 Mouse NGAL/LCN2 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)
NGAL/LCN2	24	21-27

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 2.5X the standard deviations above the background (zero Calibrator).

Precision

Control	Average Conc. (pg/mL)	Average Intra-run Conc. (%CV)	Inter-run Conc. (%CV)
High	6,230	3.1	9.2
Mid	1,460	2.9	10.2
Low	418	4.1	12.2

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 Mouse NGAL/LCN2

Tested Samples

Sample Type	Serum (N=6)	EDTA Plasma (N=6)
Median (pg/mL)	107,000	146,000
Range (pg/mL)	64,900-184,000	109,000-162,000
% Detected	100	100

Normal serum and plasma samples were diluted 2-fold prior to the assay.

Dilution Linearity

Serum			EDTA Plasma		
Fold Dilution	Average % Recovery	% Recovery Range	Fold Dilution	Average % Recovery	% Recovery Range
2	149	138-165	2	159	146-175
4	184	150-225	4	222	187-265
8	193	102-243	8	266	201-319

Normal mouse serum and EDTA plasma were spiked with Calibrator and tested at different dilutions. Undiluted samples were tested to determine the expected concentration of the analyte. Samples may benefit from additional dilution with assay diluent to reduce matrix effects.

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

Spike Recovery

Spike Level	Serum		EDTA Plasma	
	Average % Recovery	% Recovery Range	Average % Recovery	% Recovery Range
High	93	92-97	94	82-101
Mid	98	96-102	99	97-102
Low	98	97-101	97	95-100

Normal serum and plasma were spiked with Calibrator at 3 levels. Undiluted samples were tested to determine the expected concentration of the analyte. Samples may benefit from additional dilution with assay diluent to reduce matrix effects.

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

Specificity

To assess specificity, the NGAL/LCN2 Antibody Set was tested individually against a larger panel of analytes for nonspecific binding (6CKine/Ccl21, BAFF, BCA-1/BLC, CD40, Eotaxin, EPO, GM-CSF, IFN- α , IFN- β , IFN- γ , IL-1 β , IL-2, IL-4, IL-5, IL-6, IL-9, IL-10, IL-12/IL-23p40, IL-12p70, IL-13, IL-15, IL-16, IL-17A, IL-17A/F, IL-17C, IL-17E/IL-25, IL-17F, IL-21, IL-22, IL-23, IL-27p28/IL-30, IL-31, IL-33, IP-10, KC/GRO, MCP-1, MCP-5/Ccl12, MDC, MIP-1 α , MIP-1 β , MIP-2, MIP-3 α , MMP-9 (total), NGAL/LCN2, RANTES, SDF-1 α , TARC, TNF-RI, TNF- α , VEGF-A). 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 41 and Antibody Diluent 45. 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: NGAL/LCN2 is included in Calibrator 17. The mouse NGAL/LCN2 Calibrator is a full-length recombinant protein expressed in a mouse cell line.

Antibodies: The U-PLEX Mouse NGAL/LCN2 Assay uses a rat monoclonal antibody for capture and a rat 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.

