



Human Factor VII

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

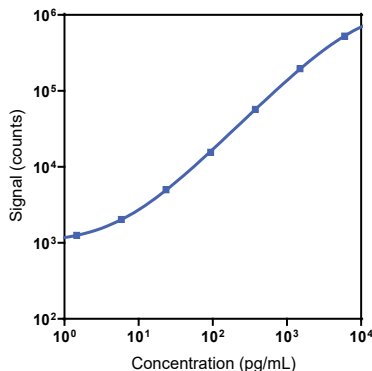
Company Address

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)
	K151N9K-1/-2/-4	U-PLEX Human Factor VII Assay with SECTOR™ plates
Singleplex	K151N9K-21/-22/-24	U-PLEX Human Factor VII Assay with QuickPlex® plates
	K251N9K-2/-4	U-PLEX Human Factor VII Assay with 384-well plates
Antibody Set	B21N9-2/-3	U-PLEX Human Factor VII 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 Factor VII 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)
Factor VII	0.60	0.48-0.63

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	857	3.6	9.0
Mid	199	3.0	5.3
Low	45	5.6	10.6

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 Factor VII

Tested Samples

Sample Type	Serum (N=10)	EDTA Plasma (N=10)	Citrate Plasma (N=5)
Median (µg/mL)	0.10	0.10	0.13
Range (µg/mL)	0.05-0.21	0.07-0.13	0.09-0.14
% Detected	100	100	100

Normal serum, EDTA plasma, and citrate plasma samples were diluted 4,000-fold prior to the assay.

Parallelism

Serum			EDTA Plasma		
Fold Dilution	Average % Recovery	% Recovery Range	Fold Dilution	Average % Recovery	% Recovery Range
2,000	96	93-103	2,000	107	96-120
8,000	94	88-104	8,000	105	97-114
16,000	94	87-102	16,000	94	90-98

Normal human serum and EDTA plasma were tested at different dilutions. Percent recovery at each dilution level was normalized to the concentration of 4,000-fold diluted samples.

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

Specificity

To assess specificity, the Factor VII 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: Factor VII is included in Calibrator 25. The human Factor VII Calibrator is a full-length recombinant protein expressed in a human cell line.

Antibodies: The U-PLEX Human Factor VII Assay uses a sheep polyclonal antibody for capture and a goat polyclonal 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 4,000-fold prior to analysis in this assay.

