

MSD[®] Total GLP-1 (ver. 2) Assay Kit



DATA SHEET

For the quantitative determination of Total GLP-1 in human, mouse, rat serum and plasma

- Alzheimer's Disease
- BioProcess
- Cardiac
- Cell Signaling
- Clinical Immunology
- Cytokines
- Hypoxia
- Immunogenicity
- Inflammation
- Metabolic**
- Oncology
- Toxicology
- Vascular

MSD Technology Overview

MESO SCALE DISCOVERY'S MULTI-ARRAY[®] Technology is a multiplex immunoassay system that enables the measurement of biomarkers utilizing the next generation of electrochemiluminescent detection. In an MSD assay, specific capture antibodies for the analytes are coated in arrays in each well of a 96-well carbon electrode plate surface. The detection system uses patented SULFO-TAG[™] labels that emit light upon electrochemical stimulation initiated at the electrode surfaces of the MULTI-ARRAY and MULTI-SPOT[®] plates. MSD assays have low background, are highly stable and are non-radioactive.

MSD assays require minimal sample volume compared to traditional ELISA. With an MSD assay, ten different biomarkers can be analyzed simultaneously with typical sample volumes less than 25 μ L. These assays have high sensitivity, up to five logs of linear dynamic range, and excellent performance in complex biological matrices. Combined, these advantages enable the measurement of native levels of biomarkers in normal and diseased samples without multiple dilutions.

This datasheet outlines the performance of MSD Total GLP-1 (ver. 2) assay.

Glucagon-like peptide-1 (GLP-1), a post-translational product of proglucagon, is a 3.5 kD protein hormone produced in intestinal L cells and plays a key role in the promotion of glucose-dependent insulin secretion and insulin biosynthesis. In addition, GLP-1 works in concert with insulin to inhibit glucose secretion and thus lower overall blood glucose levels. Through the activation of different physiological systems, it plays roles in gastric emptying upon nutrient intake and in the regulation of short-term feeding behavior. Upon release, its action is mediated through a single G-protein-coupled receptor. The cleaved peptides, commonly referred to as GLP-1 (7-36) amide and GLP-1 (7-37) are the biologically active forms of GLP-1. In vivo, these active isoforms are rapidly cleaved by dipeptidyl peptidase IV (DPP IV). The primary amino acid sequence for GLP-1 is conserved among mammalian species, i.e. human, mouse, rat, monkey, canine, etc. MSD offers a comprehensive array of GLP-1 assays that measure active, total and amidated isoforms of the GLP-1 protein using detection antibodies that recognize the amino acids in the C-terminus region of the peptide.

Assay Sensitivity

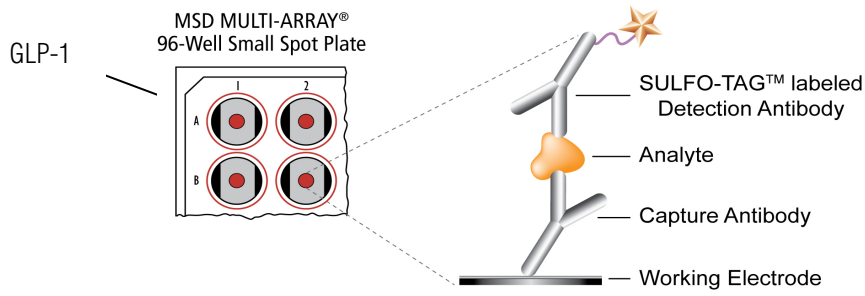
Note: 1 pmol/L = 3.297pg/mL

Total GLP-1	
LLOD (pg/mL)	0.98

The lower limit of detection (LLOD) is the calculated concentration of the signal that is 2.5 standard deviations over the zero Calibrator.

The MSD Total GLP-1 (ver. 2) assay is available in the singleplex format on MSD 96-well Small Spot plate. This assay detects all isoforms of GLP-1 using 25 μ L sample volume.

For a complete list of products, please check our website at www.mesoscale.com.



Catalog Numbers

Total GLP-1 (ver. 2) Kit	
Kit size	
1 plate	K150JVC-1
5 plates	K150JVC-2
25 plates	K150JVC-4

Ordering information

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

FOR RESEARCH USE ONLY. NOT FOR USE IN DIAGNOSTIC PROCEDURES.



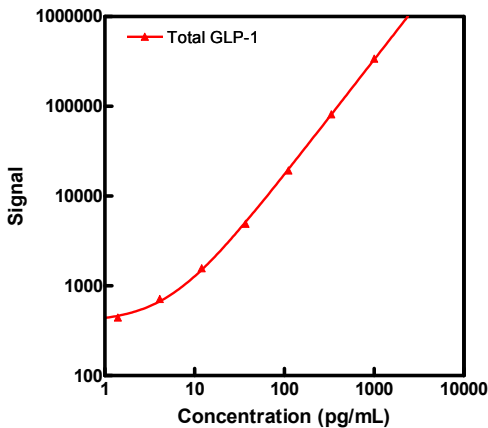
MSD Human Metabolic Assays

Example Protocol

1. Add Blocking solution, incubate for 1 hour at RT.
2. Wash. Add Assay Diluent + Calibrator / Sample, incubate for 2 hours at RT.
3. Wash. Add Detection Antibody, incubate for 1 hour at RT.
4. Wash. Add Read Buffer T, read.

Typical Standard Curve:

The following standard curves demonstrate the wide dynamic range (3-4 logs) of the MSD Total GLP-1 (ver. 2) assay. This allows for accurate quantification of many biological samples without the need for dilution.



Total GLP-1		
Conc. (pg/mL)	Average Signal	%CV
0	328	5.8
1.4	445	5.1
4.1	714	3.8
12	1570	4.6
37	4913	7.6
111	19309	4.7
333	81466	3.1
1000	338128	1.1

Spike Recovery:

Serum, EDTA plasma, and heparin plasma samples from human, mouse and rat were spiked with the Calibrators at multiple values throughout the range of the assay. Measured analyte represents average spike recovery in multiple pooled serum and plasma samples. Results of spike-recovery may vary based on the individual samples.

% Recovery = measured /expected x 100

Sample	Human			% Recovery
	Spike Conc. (pg/mL)	Measured Conc. (pg/mL)	Measured Conc. %CV	
Serum	0	16	7.3	
	10	27	1.0	105
	100	112	1.2	97
	800	817	0.0	100
EDTA Plasma	0	25	0.1	
	10	34	0.7	98
	100	119	2.8	96
	800	835	2.1	101
Heparin Plasma	0	6.1	3.9	
	10	14	4.2	86
	100	125	3.8	118
	800	644	3.3	80

MSD Human Metabolic Assays

Sample	Mouse				Rat			
	Spike Conc.	Measured Conc.	Measured Conc.	% Recovery	Spike Conc.	Measured Conc.	Measured Conc.	% Recovery
	(pg/mL)	(pg/mL)	%CV		(pg/mL)	(pg/mL)	%CV	
Serum	0	45	0.4		0	13	1.6	
	10	56	5.5	101	10	24	2.0	107
	100	142	1.7	98	100	130	2.6	115
	800	896	3.6	106	800	874	0.2	108
EDTA Plasma	0	58	1.7		0	1.4	27.9	
	10	71	0.1	104	10	10	6.0	90
	100	172	1.5	108	100	70	6.3	69
	800	1005	2.2	117	800	647	8.9	81
Heparin Plasma	0	45	2.0		0	2.4	40.4	
	10	58	2.3	105	10	11	22.6	88
	100	157	2.4	108	100	68	10.6	66
	800	937	0.3	111	800	637	8.9	79

Linearity:

Linearity was measured by spiking Calibrator levels in pooled serum, EDTA and heparin plasma samples from human, mouse and rat followed by subsequent dilution. Percent recovery is calculated as the measured concentration divided by the concentration of the previous dilution (expected).

% Recovery = measured x dilution factor / expected x 100

Sample	Fold Dilution	Human		
		Conc. (pg/mL)	Conc. %CV	% Recovery
Serum	1	112	2.3	
	2	56	1.2	100
	4	27	0.6	98
	8	15	0.6	111
EDTA Plasma	1	129	10.9	
	2	66	2.5	102
	4	33	1.5	102
	8	18	2.4	107
Heparin Plasma	1	91	4.6	
	2	51	0.9	112
	4	29	3.5	115
	8	17	6.5	114



MSD Human Metabolic Assays

Sample	Fold Dilution	Mouse			Rat		
		Conc. (unit)	Conc. %CV	% Recovery	Conc. (unit)	Conc. %CV	% Recovery
Serum	1	88	0.0		85	8.3	
	2	41	1.2	94	43	7.5	100
	4	20	0.3	100	18	4.1	82
	8	10	3.7	102	10	4.8	113
EDTA Plasma	1	87	6.3		56	1.4	
	2	42	4.9	96	26	2.2	94
	4	20	0.5	97	11	2.3	86
	8	11	0.8	108	5	1.1	83
Heparin Plasma	1	73	5.0		37	6.0	
	2	36	4.8	97	20	1.6	109
	4	18	0.4	99	11	2.4	104
	8	10	0.4	109	5	0.8	87

Cross Reactivity:

The cross-reactivity shown below is calculated based on signal generated using different GLP-1 isoforms.

Total GLP-1	
Form	Cross-Reactivity
GLP-1 (7-36) amide	100%
GLP-1 (9-36) amide	38%
GLP-1 (1-36) amide	25%
GLP-1 (7-37)	34%
GLP-1 (1-37)	15%

References using MSD platform for the measurement of GLP-1:

1. Read, P.A., Khan, F.Z., Heck, P.M., Hoole, S.P., Dutka, D.P. (2010) *DPP-4 Inhibition by Sitagliptin Improves the Myocardial Response to Dobutamine Stress and Mitigates Stunning in a Pilot Study of Patients with Coronary Artery Disease*. Circ Cardiovasc Imaging. 2010 Jan 14 [Epub ahead of print]
2. Sauve, M., Ban, K., Momen, M.A., Zhou, Y.Q., Henkelman, R.M., Husain, M., Drucker, D.J. (2010) *Genetic deletion or pharmacological inhibition of dipeptidyl peptidase-4 improves cardiovascular outcomes following myocardial infarction in mice*. Diabetes. Vol. 59(4):1063-73
3. Fujita, Y., Wideman, R.D., Speck, M., Asadi, A., King, D.S., Webber, T.D., Haneda, M., Kieffer, T.J. (2009) *Incretin release from gut is acutely enhanced by sugar but not by sweeteners in vivo*. Am J Physiol Endocrinol Metab. Vol. 296(3):E473-9
4. Lauffer, L.M., Iakoubov, R., Brubaker, P.L. (2009) *GPR119 is essential for oleylethanolamide-induced glucagon-like peptide-1 secretion from the intestinal enteroendocrine L-cell*. Diabetes. Vol. 58(5):1058-66

Company Address:

Meso Scale Discovery
A division of Meso Scale Diagnostics, LLC.
9238 Gaither Road
Gaithersburg, MD 20877 USA
Website: www.mesoscale.com

MESO SCALE DISCOVERY, MESO SCALE DIAGNOSTICS, WWW.MESOSCALE.COM, MSD, MSD (DESIGN), DISCOVERY WORKBENCH, QUICKPLEX, MULTI-ARRAY, MULTI-SPOT, SULFO-TAG, SECTOR, SECTOR HTS and SECTOR PR are trademarks and/or service marks of Meso Scale Diagnostics, LLC.
© 2010 Meso Scale Diagnostics, LLC. All rights reserved.

FOR RESEARCH USE ONLY. NOT FOR USE IN DIAGNOSTIC PROCEDURES.