

# MSD<sup>®</sup> Mouse/Rat C-peptide Kit

For quantitative determination in mouse and rat serum and plasma



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

## Catalog Numbers

Mouse/Rat C-peptide Kit	
Kit Size	Catalog #
1 plate	K1500ID-1
5 plates	K1500ID-2
25 plates	K1500ID-4

## Ordering Information

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

## Scientific Support

Phone: 1-301-947-2025  
Email: ScientificSupport@mesoscale.com

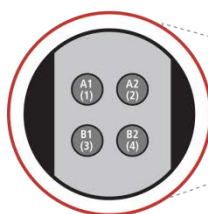
## Company Address

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

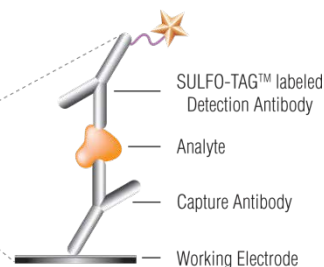
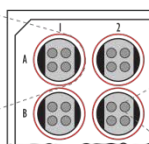
[www.mesoscale.com](http://www.mesoscale.com)<sup>®</sup>

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

1. **C-peptide**
2. *BSA blocked*
3. *BSA blocked*
4. *BSA blocked*



MSD MULTI-SPOT<sup>®</sup>  
96-Well 4-Spot Plate

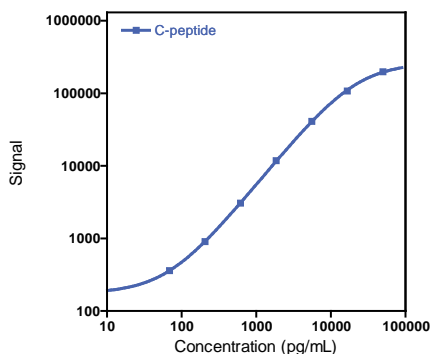


Connecting peptide, or C-peptide, is a 31 amino acid protein constituent of proinsulin produced by pancreatic  $\beta$ -cells in the Islets of Langerhans. C-peptide connects the A-chain and the B-chain of the proinsulin polypeptide and facilitates proper processing of proinsulin in the endoplasmic reticulum. Proinsulin is cleaved by  $\text{Ca}^{2+}$ -dependent endopeptidases; equimolar quantities of active insulin and C-peptide are stored in secretory granules, released into the portal vein, enter circulation, and are excreted by the kidneys.<sup>1</sup> Circulating C-peptide binds multiple cell types, including endothelial cells, fibroblasts, neuronal cells, and renal tubular cells. Bound C-peptide activates  $\text{Ca}^{2+}$ -dependent signaling pathways, including MAPK,  $\text{PLC}\gamma$ , and PKC, resulting in both increased eNOS and  $\text{Na}^+\text{K}^+\text{ATPase}$  activity and upregulation of transcription factors.<sup>2</sup>

Plasma levels of C-peptide have traditionally been used to distinguish between Type 1 and Type 2 diabetes.<sup>3</sup> Type 1 diabetics fail to produce proinsulin so that C-peptide levels are decreased, while Type II diabetics produce increased quantities of proinsulin so that C-peptide levels are increased. C-peptide is an active biomolecule of interest in the study of hormone-dependent conditions, including gastromas associated with multiple endocrine neoplasms syndrome, polycystic ovarian syndrome, metabolic syndrome X, and sub-sets of renal disease.<sup>4</sup> The MSD Mouse/Rat C-peptide assay is available on 96-well, 4-spot plates. Representative data from the assay is presented below.

## Assay Sensitivity

The following standard curve illustrates the dynamic range of the Mouse/Rat C-peptide assay.



C-peptide	
Average LLOD (pg/mL)	25
LLOD Range (pg/mL)	18–32

The lower limit of detection (LLOD) is a calculated concentration based on a signal 2.5 standard deviations above the background (zero calibrator blank). The LLOD shown above was calculated based on 3 runs.

## MSD Advantage

- **Multiplexing:** Multiple analytes can be measured in one well using typical sample volumes of 25  $\mu\text{L}$  or less without compromising speed or performance
- **Large dynamic range:** Linear range of up to five logs enables the measurement of native levels of biomarkers in normal and diseased samples without multiple dilutions
- **Minimal background:** The stimulation mechanism (electricity) is decoupled from the response (light signal), minimizing matrix interference
- **Simple protocols:** Only labels bound near the electrode surface are excited, enabling assays with fewer washes
- **Flexibility:** Labels are stable, non-radioactive, and conveniently conjugated to biological molecules
- **High sensitivity and precision:** Multiple rounds of label excitation and emission enhance light levels and improve sensitivity

For a complete list of products, please visit our website at [www.mesoscale.com](http://www.mesoscale.com).

# MSD Metabolic Assays

## Tested Samples

Normal mouse plasma and rat serum samples were collected and tested with the Mouse/Rat C-peptide Kit. Median and range of concentrations for each sample set are displayed below.

Sample Type	Statistic	C-peptide
Mouse EDTA plasma	Median (pg/mL)	335
	Range (pg/mL)	284–391
	Number of Samples	4
	Samples in Quantitative Range	4
Rat serum	Median (pg/mL)	209
	Range (pg/mL)	108–693
	Number of Samples	3
	Samples in Quantitative Range	3

## References

1. Steiner DF, et al. Insulin biosynthesis: evidence for a precursor. *Science*. 1967 Aug 11;157(3789):697-700.
2. Hills CE, et al. Intracellular signaling by C-peptide. *Exp Diabetes Res*. 2008;2008:635158.
3. Vasic C, et al. C-peptide: a new mediator of atherosclerosis in diabetes. *Mediators Inflamm*. 2012;2012:858692.
4. Wahren J. C-peptide: new findings and therapeutic implications in diabetes. *Clin Physiol Funct Imaging*. 2004 Jul;24(4):180-9.

MESO SCALE DISCOVERY, MESO SCALE DIAGNOSTICS, DISCOVERY WORKBENCH, MULTI-ARRAY, MULTI-SPOT, QUICKPLEX, SECTOR PR, SECTOR, SECTOR HTS, SULFO-TAG, www.mesoscale.com, SMALL SPOT (design), 96 WELL 1, 4, 7, & 10-SPOT (designs), 384 WELL 1 & 4-SPOT (designs), MSD, MSD (design), and V-PLEX, STREPTAVIDIN GOLD, SPOT THE DIFFERENCE are trademarks and/or service marks of Meso Scale Diagnostics, LLC. © 2013 Meso Scale Diagnostics, LLC. All rights reserved.

