

Product Number:	100199
Lot Number:	502666
Expiration:	31-Aug-2021
Platform(s):	HD-1/HD-X & SR-X



Data below represents results generated on the Simoa™ HD-1 Analyzer.

Component	Part Number	Lot Number
Bead Reagent	101045	035606
Detector Reagent	101047	035108
SBG Reagent	101057	035613
Sample Diluent	101055	024420
Calibrator Diluent	101051	031908
Calibrator Concentrate	101049	006306
SBG Diluent (SR-X)†	103158	N/A ¹
RGP Reagent	103159	N/A ¹

¹ Reagents are not Kit Lot Specific

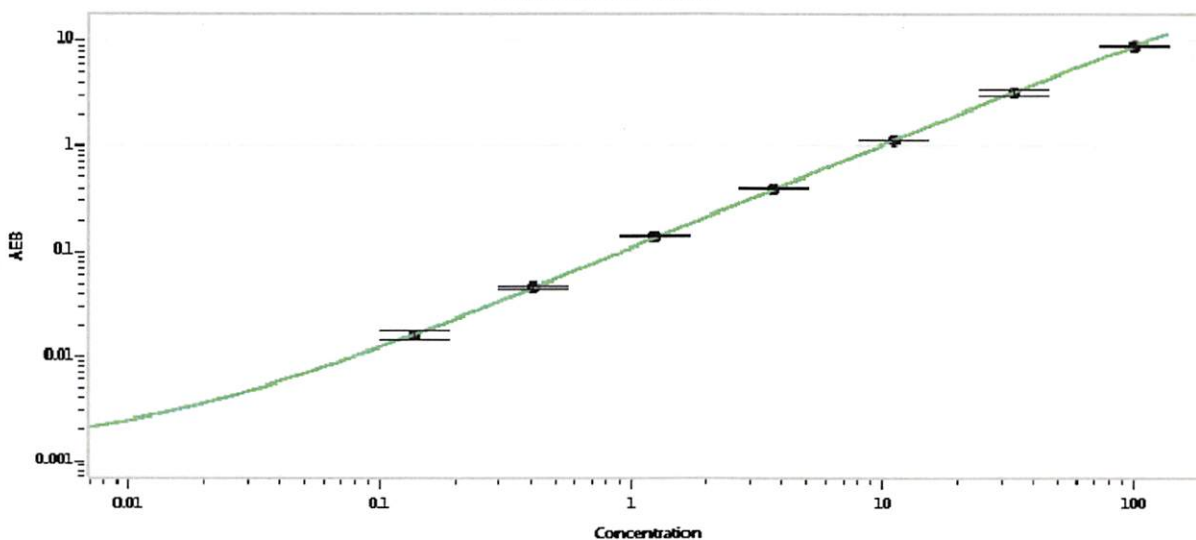
† Specific for use on the SR-X platform - refer to the Simoa™ C-Peptide Advantage SR-X Kit Instructions (KI-0109).

Release Materials*			
	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	MF-032017-04	1.77	1.32-1.97
Control 2	MF-032017-04	55.0	41.9-62.9
Panel 1	818703	72.2	67.3-111
Panel 2	MF-060118-02	1534	1481-2310
Panel 3	MF-060118-03	2711	2168-3264

* Ranges shown are generated internally for new lot release only. Customer should generate their own control ranges.

Calibrator Concentrate Value Assignment (pg/mL)	9338
---	------

Calibration Curve



Calibrator Levels (pg/mL)	
A	0.000
B	0.137
C	0.412
D	1.23
E	3.70
F	11.1
G	33.3
H	100

Review/Approval

Junpil Park
Name

QC Supervisor
Title

Signature/Date
30 Dec 2020