

Product Number:	100199
Lot Number:	502272
Expiration:	12-Dec-2020
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	005512
Detector Reagent	101047	005506
SBG Reagent	101057	006609
Sample Diluent	101055	004911
Calibrator Diluent	101051	005207
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

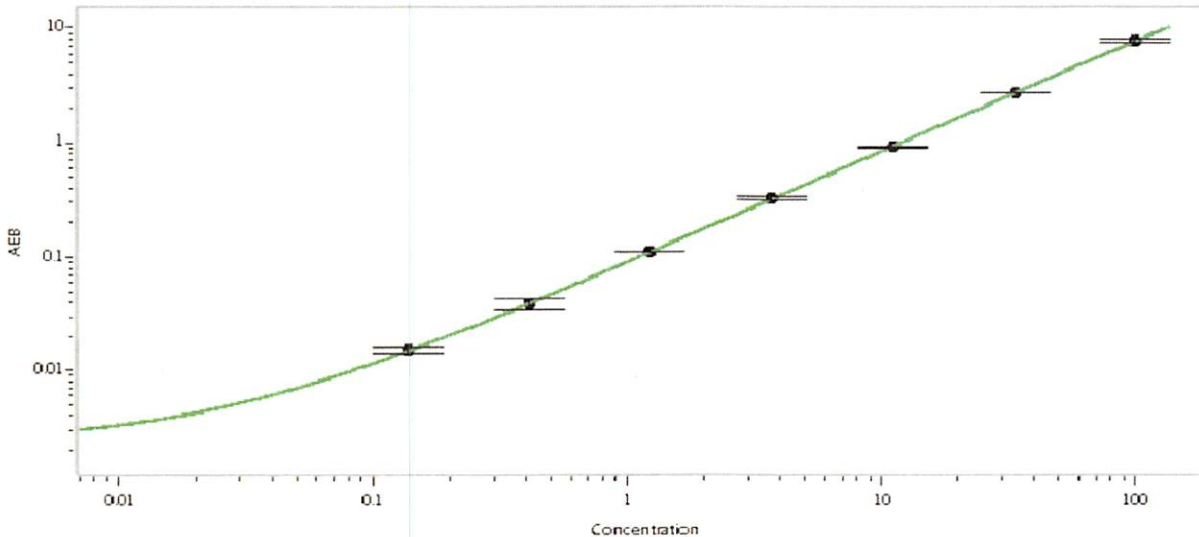
† Specific for use on the SR-X platform - refer to the Simoa™ C-Peptide Advantage SR-X Kit Instructions (KI-0109).

Release Materials*			
	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	MF-032017-04	1.63	1.32-1.97
Control 2	MF-032017-04	48.6	41.9-62.9
Panel 1	818704	107	91.6-137
Panel 2	MF-060118-02	2040	1481-2310
Panel 3	MF-060118-03	2522	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

Patrizia Stadler Assoc. Director, QC

Name

Title

P. Stadler

11 MAR 2020

Signature/Date