

Product Number:	100199
Lot Number:	501860
Expiration:	21-Dec-2019
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	916303
Detector Reagent	101047	916405
SBG Reagent	101057	913406
Sample Diluent	101055	835512
Calibrator Diluent	101051	912016
Calibrator Concentrate	101049	81302
SBG Diluent (SR-X)†	103158	N/A ¹
RGP Reagent	103159	N/A ¹

¹ Reagents are not Kit Lot Specific

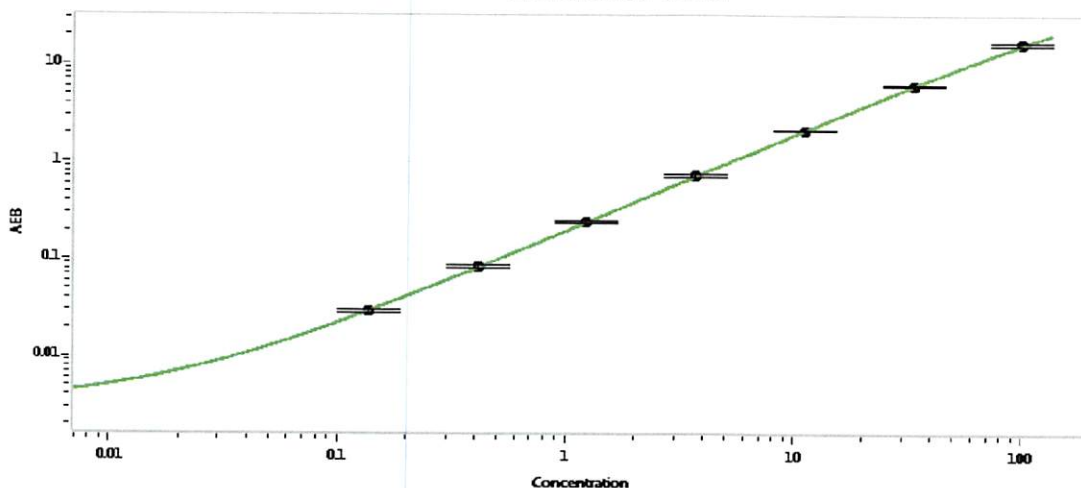
† Specific for use on the SR-X platform - refer to the Simoa™ C-Peptide Advantage SR-X Kit Instructions (KI-0109).

Release Materials*			
	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	MF-032017-04	1.78	1.32-1.97
Control 2	MF-032017-04	50.7	42.4-63.5
Panel 1	818704	110	91.6-137
Panel 2	MF-060118-02	1625	1458-2290
Panel 3	MF-060118-03	2443	2193-3290

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

Calibrator Concentrate Value Assignment (pg/mL)	1832
---	------

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 24 JUN 2019
Signature/Date