

Product Number:	100199
Lot Number:	501672
Expiration:	9-Jul-2019
Platform(s):	HD-1 & SR-X



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

Component	Part Number	Lot Number
Bead Reagent	101045	821411
Detector Reagent	101047	821505
SBG Reagent	101057	821317
Sample Diluent	101055	819024
Calibrator Diluent	101051	819815
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

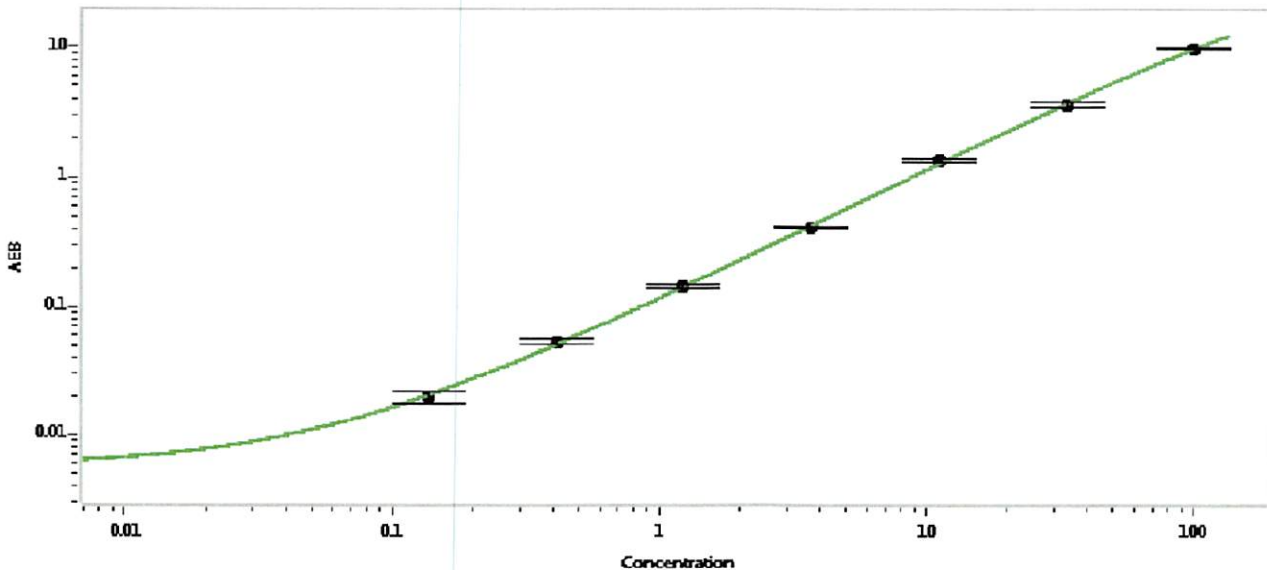
† Specific for use on the SR-X platform - refer to the Simoa™ C-Peptide Advantage SR-X Kit Instructions (KI-0109).

Release Materials*			
	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	MF-032017-04	1.60	1.30-1.95
Control 2	MF-032017-04	49.2	42.7-64.0
Panel 1	818701	0.867	0.354-1.32
Panel 2	MF-060118-02	1850	1534-2301
Panel 3	MF-060118-03	2824	2230-3345

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

Calibrator Concentrate Value Assignment (pg/mL)	1,832
---	-------

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: Patrizia Stadler, Title: Assoc. Director, QC, Signature: p. Stadler, Date: 21 Feb 2019