

Product Number:	100199
Lot Number:	503178
Expiration:	3-Mar-2022
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	128201
Detector Reagent	101047	115910
SBG Reagent	101057	116707
Sample Diluent	101055	115919
Calibrator Diluent	101051	111728
Calibrator Concentrate	101049	006306
SBG Diluent (SR-X) †	103158	N/A <sup>1</sup>
RGP Reagent	103159	N/A <sup>1</sup>

<sup>1</sup> Reagents are not Kit Lot Specific

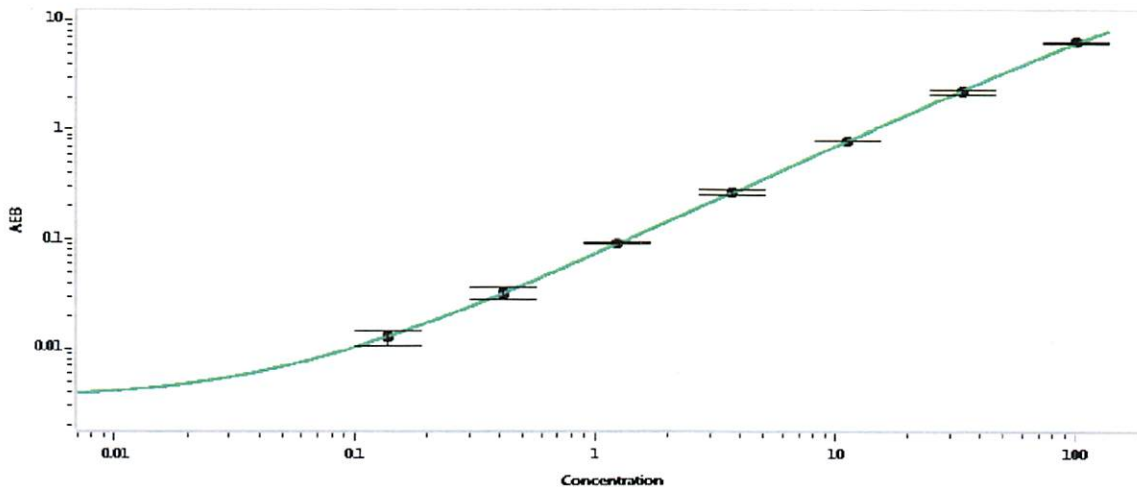
† Specific for use on the SR-X platform - refer to the Simoa™ C-Peptide Advantage SR-X Kit Instructions (KI-0109).

Release Materials*			
	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	006306	57.1	45.0 - 75.0
Control 2	006306	1.93	1.50 - 2.50
Panel 1	102505	1800	1209 - 2076
Panel 2	102505	691	618 - 1042
Panel 3	102505	674	596 - 894

\* 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 19 Oct 2021