

Product Number:	100199
Lot Number:	503810
Expiration:	28-Nov-2023
Platform(s):	HD-1/HD-X & SR-X

Component	Part Number	Lot Number
Bead Reagent	101045	304111
Detector Reagent	101047	304405
SBG Reagent	101057	304406
Sample Diluent	101055	304407
Calibrator Diluent	101051	303217
Calibrator Concentrate	101049	133212
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).



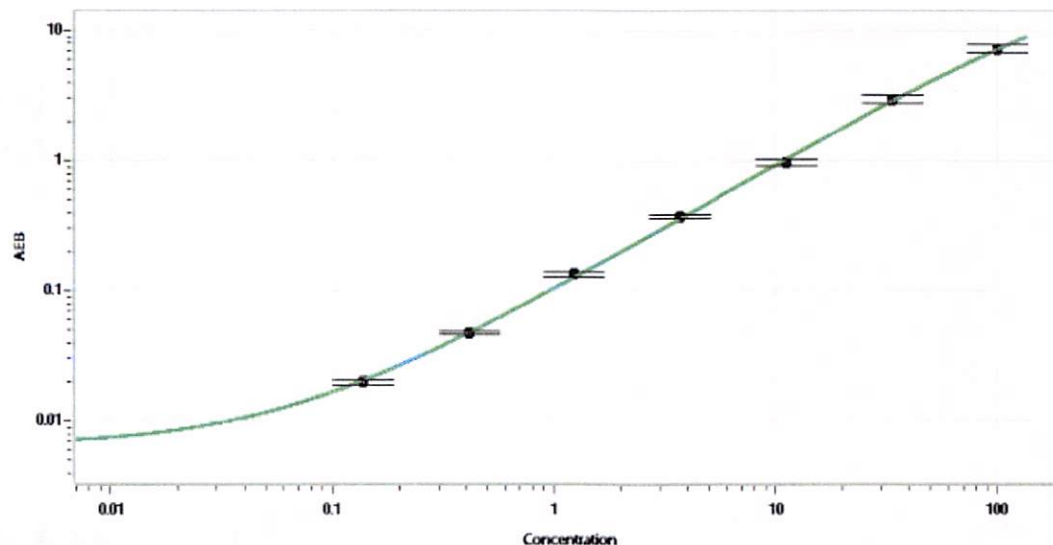
Data below represents results generated on the Simoa® HD-X.

Release Materials*			
	Lot	Result (pg/mL)	Spec. Range (pg/mL)
Control 1	119615	1.75	1.00 - 2.46
Control 2	119615	49.5	36.5 - 58.6
Panel 1	320804	13.7	10.8 - 16.9
Panel 2	320804	1195	937 - 1527
Panel 3	320804	2863	2226 - 3694

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

Calibrator Concentrate Value Assignment (pg/mL)	7027
---	------

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: Sonia Khillan Name, QC Manager Title, Sonia Khillan 22 AUG 2023 Signature/Date