

Product Number:	100199
Lot Number:	503953
Expiration:	23-Jun-2024
Platform(s):	HD-1/HD-X & SR-X



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

Component	Part Number	Lot Number	Control Lot Number
Bead Reagent	101045	327505	N/A
Detector Reagent	101047	327804	N/A
SBG Reagent	101057	327805	N/A
Sample Diluent	101055	317410	317410A
Calibrator Diluent	101051	325613	N/A
Calibrator Concentrate	101049	324007	N/A
SBG Diluent (SR-X) †	103158	N/A ¹	N/A ¹
RGP Reagent	103159	N/A ¹	N/A ¹

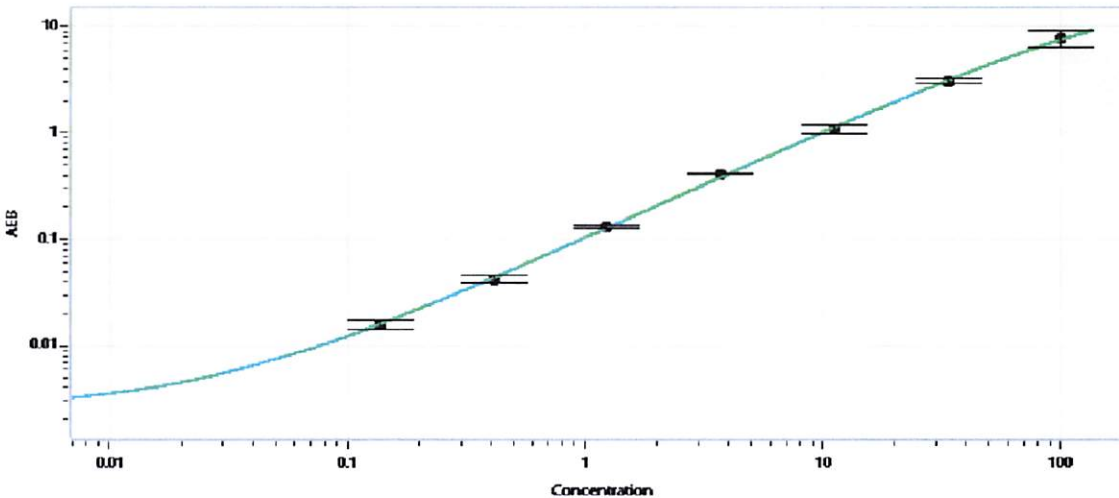
¹ Reagents are not Kit Lot Specific

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

Release Materials*				* Ranges shown are generated internally for new lot release only. Customer should generate their own control ranges.
	Lot	Result (pg/mL)	Spec. Range (pg/mL)	
Control 1	119615	1.72	1.07 - 2.39	
Control 2	119615	45.4	36.9 - 57.3	
Panel 1	320804	12.2	10.2 - 16.8	
Panel 2	320804	1110	912 - 1492	
Panel 3	320804	2993	2298 - 3638	

Calibrator Concentrate Value Assignment (pg/mL)	1491
---	------

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 12 OCT 2023
Signature/Date