

Product Number:	100199
Lot Number:	504188
Expiration:	8-Nov-2024
Platform(s):	SR-X and HD-1/HD-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	327805A
Sample Diluent	101055	317410	317410A
Calibrator Diluent	101051	325613	N/A
Calibrator Concentrate	101049	324007	N/A
SBG Diluent (SRX)†	103158	N/A ¹	N/A ¹
RGP Reagent	103159	N/A ¹	N/A ¹

¹ Reagents are not Kit Lot Specific

† Specific to use on the SRX platform.

Refer to the Simoa® C-Peptide SRX Kit Instructions (KI-0109).

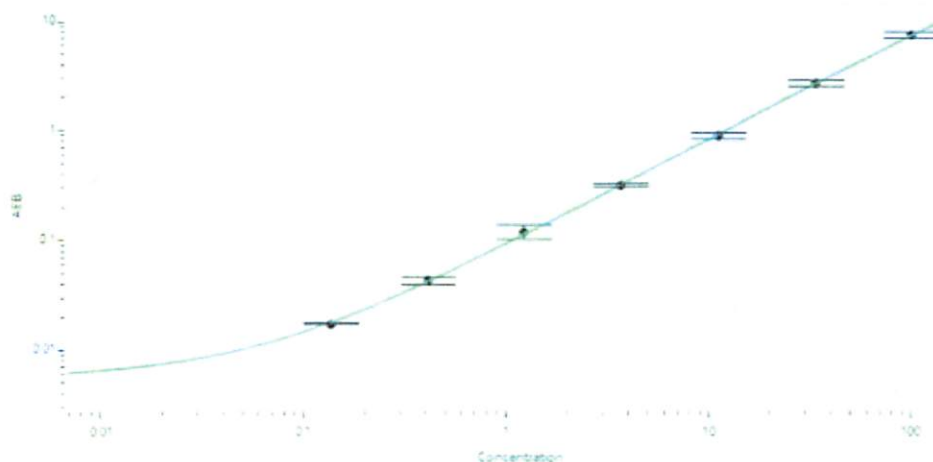
Release Materials*			
	Lot	Result (pg/mL)	Spec. Range (pg/mL)
Control 1	324007	56.6	45.0 - 75.0
Control 2	324007	2.33	1.50 - 2.50
Panel 1	320804	13.6	10.3 - 16.7
Panel 2	320804	1259	933 - 1491
Panel 3	320804	2965	2287 - 3647

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

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

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

Cody Carlson
Name

QC Manager
Title

Cody Carlson 22 MAR 2024
Signature/Date