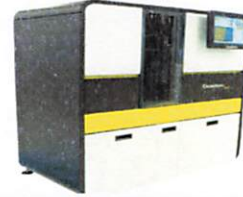


<b>Product Number:</b>	100199
<b>Lot Number:</b>	504170
<b>Expiration:</b>	23-Jun-2024
<b>Platform(s):</b>	SR-X and HD-1/HD-X



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

Component	Part Number	Lot Number	Control 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 (SR-X)†	103158	N/A <sup>1</sup>	N/A <sup>1</sup>
RGP Reagent	103159	N/A <sup>1</sup>	N/A <sup>1</sup>

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

† Specific to use on the SRX platform.

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

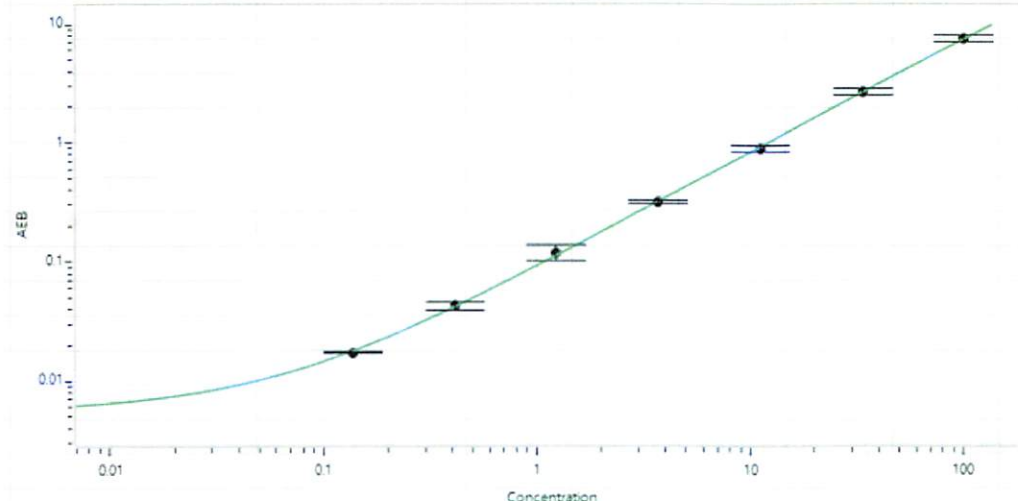
Release Materials*			
	Lot	Result (pg/mL)	Spec. Range (pg/mL)
Control 1	324007	56.6	45.0 - 75.0
Control 2	324007	2.33	1.50 - 2.50
Panel 1	320804	13.6	10.3 - 16.7
Panel 2	320804	1259	933 - 1491
Panel 3	320804	2965	2287 - 3647

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

<b>Calibrator Concentrate Value Assignment (pg/mL)</b>	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

*[Signature]* 14 MAR 2024  
Signature/Date