

<b>Product Number:</b>	104428
<b>Lot Number:</b>	504269
<b>Expiration:</b>	02-Nov-2025
<b>Platform(s):</b>	HD-X



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

Component	Part Number	Lot Number
Bead Reagent	104409	413808
Detector Reagent	104411	414111
SBG Reagent	104413	414112
Sample Diluent	104415	413816
Calibrator Diluent	104418	412334
Control Diluent	104476	412335
Calibrator Concentrate	104483	330609
Control 1 Concentrate	104426	330611
Control 2 Concentrate	104427	330612
RGP 2.0 Reagent	104321	N/A <sup>1</sup>
Activation Buffer Reagent	104324	N/A <sup>1</sup>

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

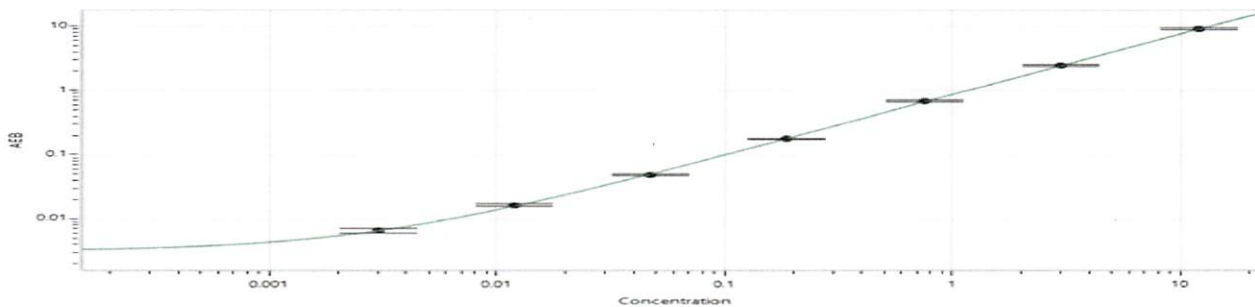
Release Materials*			
	Lot	Result (pg/mL)	Spec. Range (pg/mL)
Control 1 Concentrate	330611	0.179	0.125 - 0.233
Control 2 Concentrate	330612	5.53	3.85 - 7.15
Panel 1	328501	0.014	0.010 - 0.030
Panel 2	328501	0.532	0.364 - 0.676
Panel 3	328501	2.26	1.59 - 2.96
Panel 4	328501	5.93	3.63 - 6.73

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

<b>Calibrator Concentrate Reconstitution Volume (uL)</b>	756
<b>Control 1 Concentrate Reconstitution Volume (uL)</b>	756
<b>Control 2 Concentrate Reconstitution Volume (uL)</b>	756

Calibrator Levels (pg/mL)	
A	0.000
B	0.003
C	0.012
D	0.047
E	0.188
F	0.750
G	3.00
H	12.0

Calibration Curve



Review/Approval

Cody Carlson

QC Manager

Name

Title

*[Signature]* 18 JUN 2024  
Signature/Date