

<b>Product Number:</b>	104691
<b>Lot Number:</b>	504093
<b>Expiration:</b>	07Jan2025
<b>Platform(s):</b>	HD-X



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

Component	Part Number	Lot Number	Control Lot Number
Bead Reagent	104752	334515	N/A
Detector Reagent	104675	334711	N/A
SBG Reagent	104677	334710	N/A
Plasma Sample Diluent	104719	329811	N/A
CSF Sample Diluent	104720	329624	N/A
Calibrator Diluent	104678	318811	318811A
Control Diluent	104718	318812	318812A
Calibrator Concentrate	104687	402909	N/A
Control 1 Concentrate	104689	402910	N/A
Control 2 Concentrate	104690	402911	N/A
RGP 2.0 Reagent	104321	N/A <sup>1</sup>	N/A <sup>1</sup>
Activation Buffer Reagent	104324	N/A <sup>1</sup>	N/A <sup>1</sup>

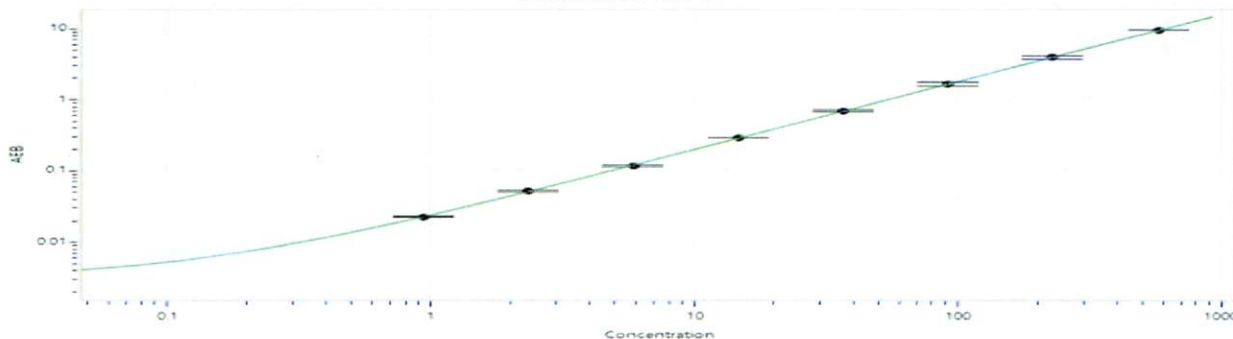
<sup>1</sup> Reagents are not Kit Lot Specific

Release Materials*			
	Lot	Result (pg/mL)	Spec. Range (pg/mL)
Control 1 Concentrate	402910	96.0	60.5 - 112
Control 2 Concentrate	402911	1163	796 - 1478
Panel 1	327811	61.2	28.8 - 77.6
Panel 2	327811	1605	744 - 1965

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

<b>Calibrator Concentrate Reconstitution Volume (uL)</b>	1144
<b>Control 1 Concentrate Reconstitution Volume (uL)</b>	1144
<b>Control 2 Concentrate Reconstitution Volume (uL)</b>	1144

Calibration Curve



Calibrator Levels (pg/mL)

A	0.000
B	0.940
C	2.34
D	5.86
E	14.7
F	36.6
G	91.2
H	229
I	572

Review/Approval

Cody Carlson

QC Manager

*Cody Carlson* 21 MAR 2024

Name

Title

Signature/Date