

<b>Product Number:</b>	104797
<b>Lot Number:</b>	504381
<b>Expiration:</b>	28-Feb-2026
<b>Platform(s):</b>	HD-X and SR-X



Component	Part Number	Lot Number
Bead Reagent	104778	424704
Detector Reagent	104780	424801
SBG Reagent	104782	424802
Plasma Sample Diluent	104783	424309
CSF Sample Diluent	104784	424211
Calibrator Diluent	104786	424715
Control Diluent	104787	424716
Calibrator Concentrate	104793	409518
Control 1 Concentrate	104795	409519
Control 2 Concentrate	104796	409520
RGP 2.0 Reagent	104321	N/A <sup>1</sup>
Activation Buffer Reagent	104324	N/A <sup>1</sup>

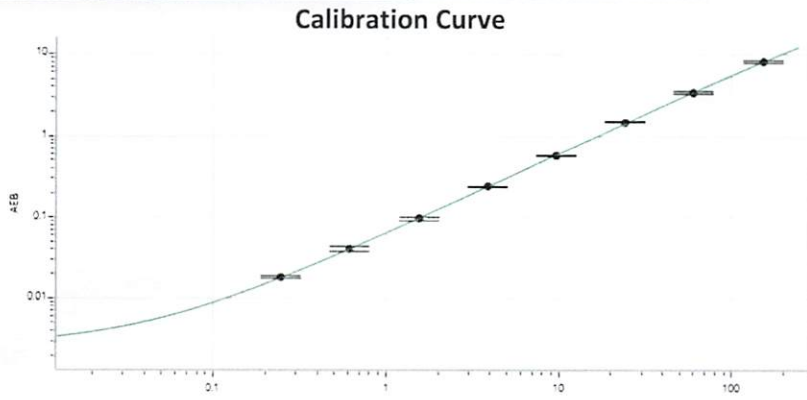
Data below represents results generated on the Simoa® HD-X Analyzer.

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

Release Materials*			* Ranges shown are generated internally for new lot release only. Customer should generate their own control ranges.
	Lot	Result (pg/mL)	
Control 1 Concentrate	409513	8.47	5.77 - 10.7
Control 2 Concentrate	409514	344	231 - 438

<b>Calibrator Concentrate Reconstitution Volume (uL)</b>	1579
<b>Control 1 Concentrate Reconstitution Volume (uL)</b>	1579
<b>Control 2 Concentrate Reconstitution Volume (uL)</b>	1579

Calibrator Levels (pg/mL)	
A	0.000
B	0.246
C	0.614
D	1.54
E	3.84
F	9.60
G	24.0
H	60.0
I	150



Review/Approval

Lynne Allen SR. mgr Doc Control Lynne Allen 15/NOV/2024

Name

Title

Signature/Date