

<b>Product Number:</b>	105796
<b>Lot Number:</b>	504628
<b>Expiration:</b>	22-Nov-2026
<b>Platform(s):</b>	HD-X



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

Component	Part Number	Lot Number
Bead Reagent	105772	515701
Detector Reagent	105774	516706
SBG Reagent	105776	516707
Sample Diluent	105777	514206
Calibrator Diluent	105779	516308
Control Diluent	105780	516309
Calibrator Concentrate	105792	507910
Control 1 Concentrate	105794	507911
Control 2 Concentrate	105795	507912
RGP 2.0 Reagent	104321	N/A <sup>1</sup>
Activation Buffer Reagent	104324	N/A <sup>1</sup>

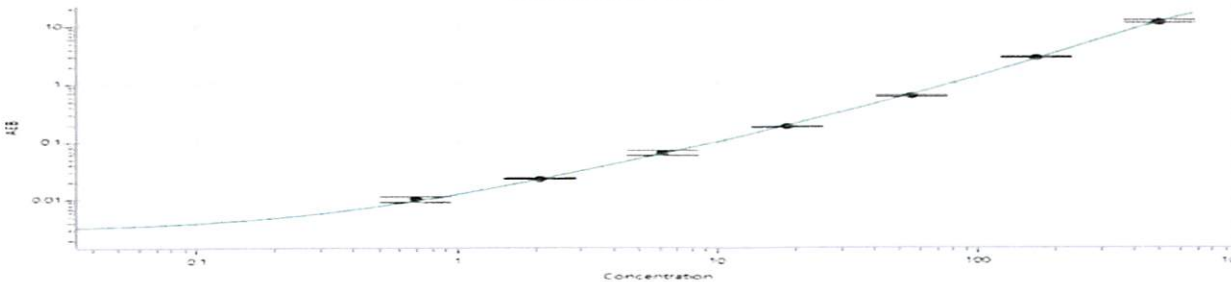
<sup>1</sup> Reagents are not Kit Lot Specific

Release Materials*			
	Lot	Result (pg/mL)	Spec. Range (pg/mL)
Control 1 Concentrate	507911	157	91.5 - 233
Control 2 Concentrate	507912	7421	5009 - 10863

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

<b>Calibrator Concentrate Reconstitution Volume (uL)</b>	1303
<b>Control 1 Concentrate Reconstitution Volume (uL)</b>	500
<b>Control 2 Concentrate Reconstitution Volume (uL)</b>	500

Calibration Curve



Calibrator Levels (pg/mL)	
A	0.000
B	0.686
C	2.06
D	6.17
E	18.5
F	55.6
G	167
H	500

Review/Approval

*Dhruvi Patel* QA Specialist II, QA *[Signature]* / 19 Aug 2025  
 Name Title Signature/Date