

<b>Product Number:</b>	104588
<b>Lot Number:</b>	504483
<b>Expiration:</b>	22-May-2026
<b>Platform(s):</b>	HD-X



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

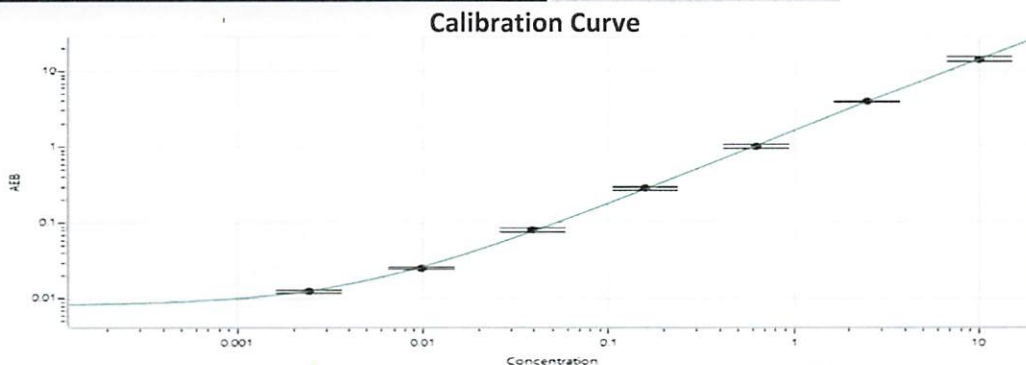
Component	Part Number	Lot Number
Bead Reagent	104572	432703
Detector Reagent	104573	433102
SBG Reagent	104574	433101
Sample Diluent Reagent	104575	434506
Calibrator Diluent Reagent	104577	432713
Control Diluent Reagent	104578	432714
Calibrator Concentrate	104584	329610
Control 1 Concentrate	104586	329612
Control 2 Concentrate	104587	329613
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*			* Ranges shown are generated internally for new lot release only. Customer should generate their own control ranges.
	Lot	Result (pg/mL)	
Control 1 Concentrate	329612	0.095	0.072 - 0.133
Control 2 Concentrate	329613	2.70	1.79 - 3.32

<b>Calibrator Concentrate Reconstitution Volume (uL)</b>	873
<b>Control 1 Concentrate Reconstitution Volume (uL)</b>	873
<b>Control 2 Concentrate Reconstitution Volume (uL)</b>	873

Calibrator Levels (pg/mL)	
A	0.000
B	0.00244
C	0.00977
D	0.0391
E	0.156
F	0.625
G	2.50
H	10.0



Review/Approval

*Dhruvi Patel*

*Doc Control I, QA*

*DP 23 Dec 2024*

Name

Title

Signature/Date