

<b>Product Number:</b>	104570
<b>Lot Number:</b>	504451
<b>Expiration:</b>	11-May-2026
<b>Platform(s):</b>	HD-X and SR-X



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

Component	Part Number	Lot Number
Bead Reagent	104546	431601
Detector Reagent	104548	431801
SBG Reagent	104550	431802
Sample Diluent	104552	433113
Calibrator Diluent Reagent	104745	433114
Control Diluent Reagent	104554	433115
Calibrator Concentrate	104566	405210
Control 1 Concentrate	104568	405211
Control 2 Concentrate	104569	405212
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

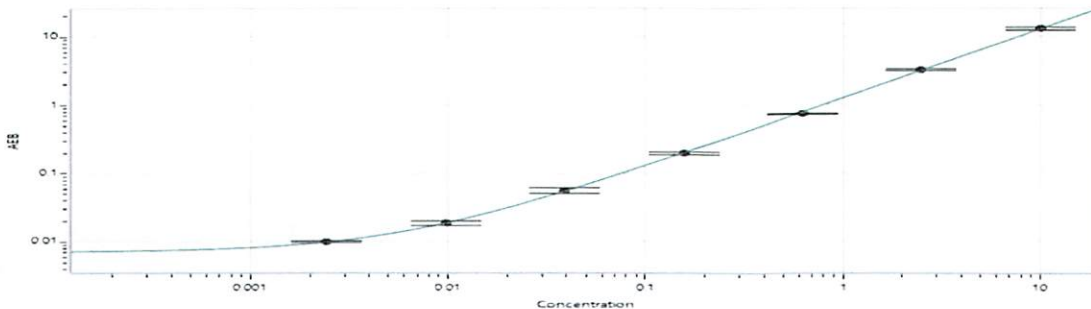
Release Materials*			
	Lot	Result (pg/mL)	Spec. Range (pg/mL)
Control 1 Concentrate	405211	0.768	0.496 - 0.922
Control 2 Concentrate	405212	7.71	5.11 - 9.49

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

<b>Calibrator Concentrate Reconstitution Volume (uL)</b>	1035
<b>Control 1 Concentrate Reconstitution Volume (uL)</b>	1035
<b>Control 2 Concentrate Reconstitution Volume (uL)</b>	1035

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

Calibration Curve



Review/Approval

*Dipesh Patel*  
Name

*Doc control IIRA*  
Title

*[Signature]* 26 DEC 2024  
Signature/Date