

Product Number:	104904
Lot Number:	504411
Expiration:	30-Mar-2026
Platform(s):	HD-X



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

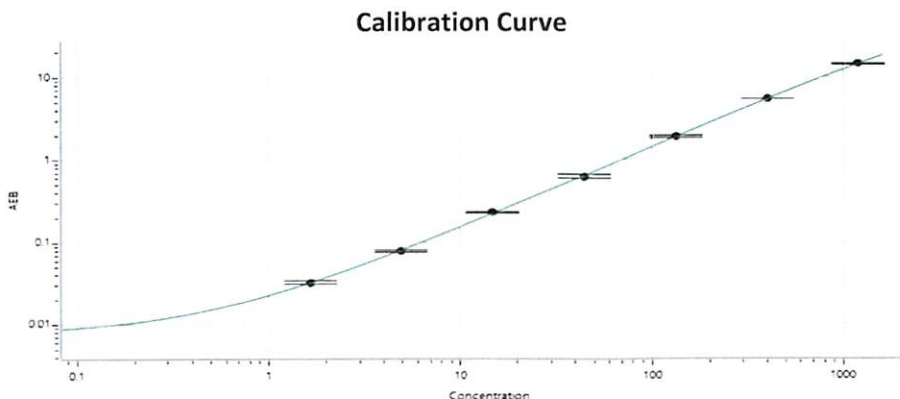
Component	Part Number	Lot Number
Bead Reagent	104880	428109
Detector Reagent	104882	428204
SBG Reagent	104884	428205
Sample Diluent	104885	427411
Calibrator Diluent	104887	427412
Control Diluent	104888	427413
Calibrator Concentrate	104900	411616
Control 1 Concentrate	104902	411618
Control 2 Concentrate	104903	411619
RGP 2.0 Reagent	104321	N/A ¹
Activation Buffer Reagent	104324	N/A ¹

¹ Reagents are not Kit Lot Specific

Release Materials*			
	Lot	Result (pg/mL)	Spec. Range (pg/mL)
Control 1 Concentrate	411618	27.0	19.6 - 36.4
Control 2 Concentrate	411619	1055	829 - 1539

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

Calibrator Concentrate Reconstitution Volume (uL)	488
Control 1 Concentrate Reconstitution Volume (uL)	488
Control 2 Concentrate Reconstitution Volume (uL)	488



Calibrator Levels (pg/mL)	
A	0.000
B	1.65
C	4.94
D	14.8
E	44.4
F	133
G	400
H	1200

Review/Approval

Dipesh Patel
Name

Doc control IT QA
Title

at 12 NOV 2024
Signature/Date