

Product Number:	104364
Lot Number:	504534
Expiration:	7-May-2027
Platform(s):	HD-X and SR-X

Component	Part Number	Lot Number	Control Lot Number
Bead Reagent	104352	517103	N/A
Detector Reagent	104353	506607	N/A
SBG Reagent	104354	506610	506610A
Sample Diluent Reagent	104355	506605	N/A
Calibrator Diluent Reagent	104358	506916	N/A
Control Diluent Reagent	104359	506917	N/A
Calibrator Concentrate	104361	412713	N/A
Control 1 Concentrate	104362	412714	N/A
Control 2 Concentrate	104363	412715	N/A
RGP 2.0 Reagent	104321	N/A ¹	N/A ¹
Activation Buffer Reagent	104324	N/A ¹	N/A ¹

¹ Reagents are not Kit Lot Specific



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

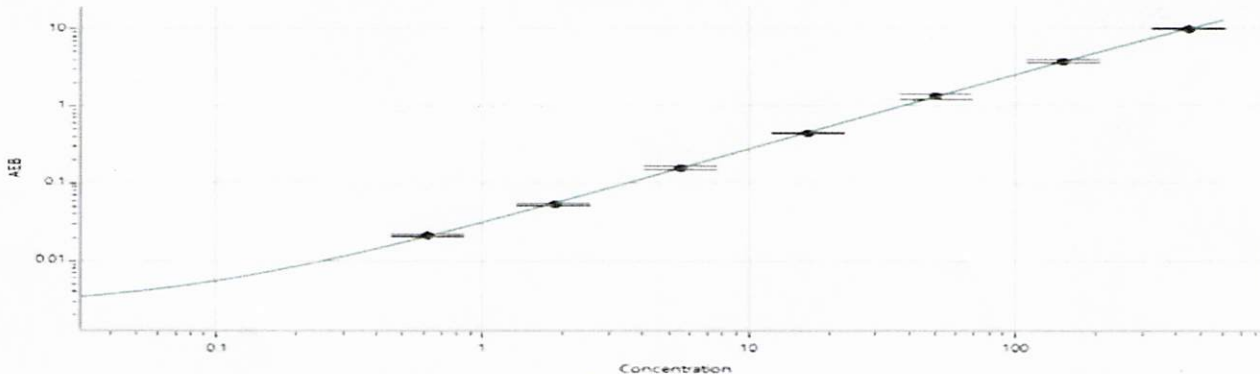
Release Materials*			
	Lot	Result (pg/mL)	Spec. Range (pg/mL)
Control 1	412714	4.14	3.56 - 6.61
Control 2	412715	344	265 - 499

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

Calibrator Concentrate Reconstitution Volume (uL)	1388
Control 1 Concentrate Reconstitution Volume (uL)	1388
Control 2 Concentrate Reconstitution Volume (uL)	1388

Calibrator Levels (pg/mL)	
A	0.000
B	0.617
C	1.85
D	5.56
E	16.7
F	50.0
G	150
H	450

Calibration Curve



Review/Approval

Dilish Patel
Name

Sr. Dir. Control QA
Title

27 Jun 2025
Signature/Date