

Product Number:	104512
Lot Number:	536615
Expiration:	15-Apr-2027
Platform(s):	HD-X and SR-X

Component	Part Number	Lot Number
Bead Reagent	104485	529603
Detector Reagent	104487	530008
SBG Reagent	104489	530009
Sample Diluent	104491	528820
Calibrator Diluent	104494	528821
Control Diluent	104495	528822
Calibrator Concentrate	104508	330613
Control 1 Concentrate	104510	330615
Control 2 Concentrate	104511	330616
RGP 2.0 Reagent	104321	N/A ¹
Activation Buffer Reagent	104324	N/A ¹

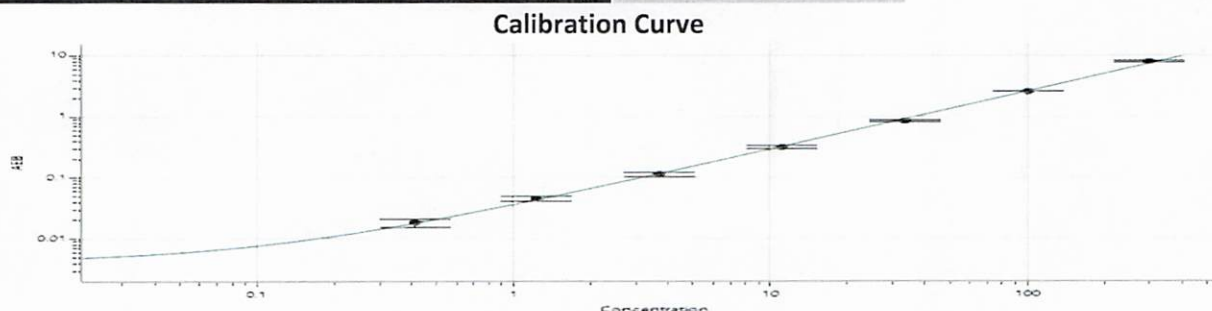
¹ Reagents are not Kit Lot Specific

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

Release Materials*				* Ranges shown are generated internally for new lot release only. Customer should generate their own control ranges.
	Lot	Result (pg/mL)	Spec. Range (pg/mL)	
Control 1 Concentrate	330615	2.81	2.56 - 4.82	
Control 2 Concentrate	330616	263	196 - 365	

Calibrator Concentrate Reconstitution Volume (uL)	749
Control 1 Concentrate Reconstitution Volume (uL)	749
Control 2 Concentrate Reconstitution Volume (uL)	749

**Calibrator Levels
(pg/mL)**



A	0.000
B	0.412
C	1.23
D	3.70
E	11.1
F	33.3
G	100
H	300

Review/Approval

Dhvani Patel QA Specialist II [Signature] / 25 Nov 2025

Name	Title	Signature/Date
Dhvani Patel	QA Specialist II	[Signature] / 25 Nov 2025