

Product Number:	104364
Lot Number:	504598
Expiration:	24-Apr-2027
Platform(s):	HD-X and SR-X

Component	Part Number	Lot Number
Bead Reagent	104352	505501
Detector Reagent	104353	511102
SBG Reagent	104354	505506
Sample Diluent Reagent	104355	505504
Calibrator Diluent Reagent	104358	505607
Control Diluent Reagent	104359	505608
Calibrator Concentrate	104361	412713
Control 1 Concentrate	104362	412714
Control 2 Concentrate	104363	412715
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.

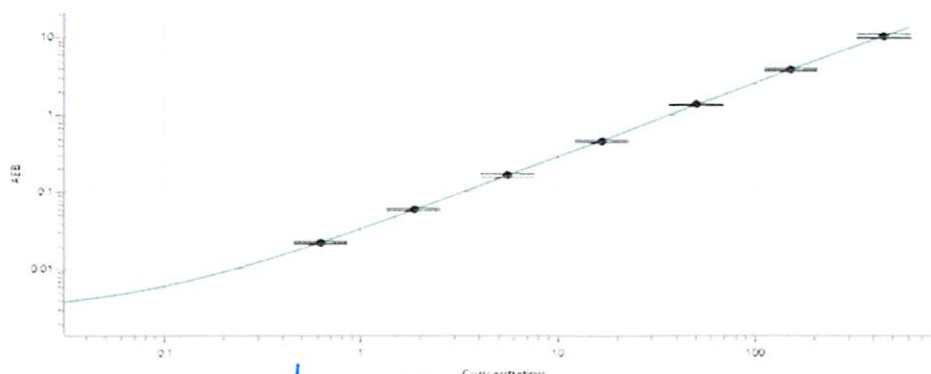
Release Materials*			
	Lot	Result (pg/mL)	Spec. Range (pg/mL)
Control 1	412714	3.68	3.11 - 5.78
Control 2	412715	315	246 - 457

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

Calibrator Concentrate Reconstitution Volume (uL)	1201
Control 1 Concentrate Reconstitution Volume (uL)	1201
Control 2 Concentrate Reconstitution Volume (uL)	1201

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

Dhruvi Patel Name
DOC Control II, QA Title

DP 12 May 2025 Signature/Date