

Product Number:	102336
Lot Number:	504559
Expiration:	4-Sep-2026
Platform(s):	SR-X and HD-1/HD-X



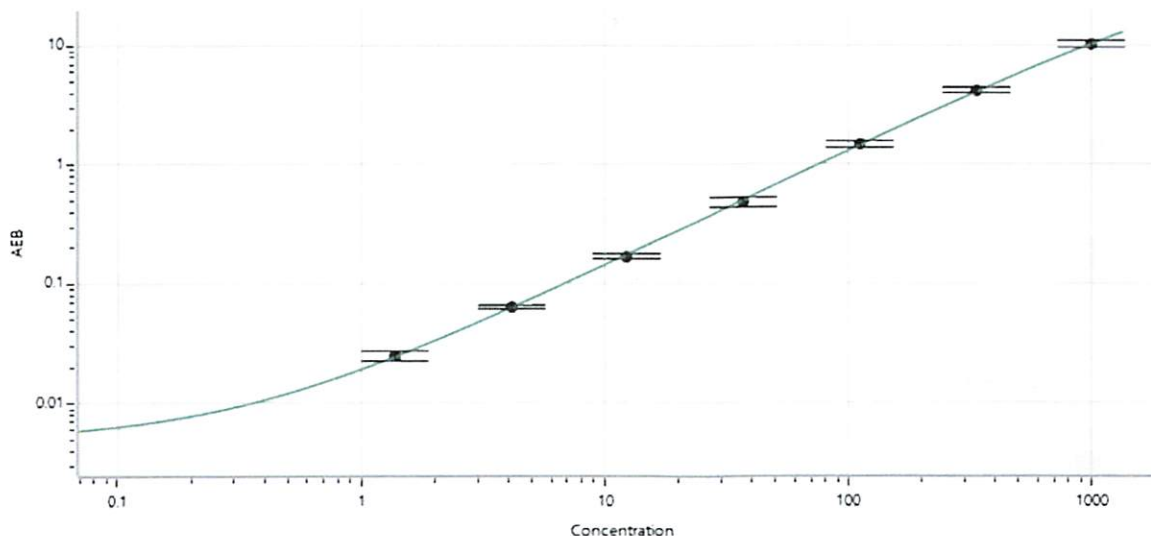
Data below represents results generated on the Simoa® HD-X Analyzer.

Component	Part Number	Lot Number
Bead Stock	102330	507604
Detector Stock	102331	507602
Calibrator Concentrate	102335	507307
Calibrator Diluent	102332	507612
Sample Diluent	102333	507614

Generic Components - Included With Kit But Not Lot Number Specific	
Discovery Bead Diluent	101954
Discovery Detector Diluent	101955
Discovery SBG Diluent	101956
Discovery SBG Stock	103297
RGP Reagent	103159

Calibrator Concentrate Target Value (ng/mL)	20
---	----

Calibration Curve



Calibrator Levels (pg/mL)	
A	0.000
B	1.37
C	4.12
D	12.3
E	37.0
F	111
G	333
H	1000

Review/Approval

Dipesh Patel Sr. Doc control-QA CT / 24 MAR 2025
Name Title Signature/Date