

Product Number:	102336
Lot Number:	504261
Expiration:	14-Dec-2024
Platform(s):	SR-X and HD-1/HD-X

Component	Part Number	Lot Number	Control Lot Number
Bead Stock	102330	317304	317304A
Detector Stock	102331	317404	317404A
Calibrator Concentrate	102335	317707	317707A
Calibrator Diluent	102332	317407	317407A
Sample Diluent	102333	317706	317706A

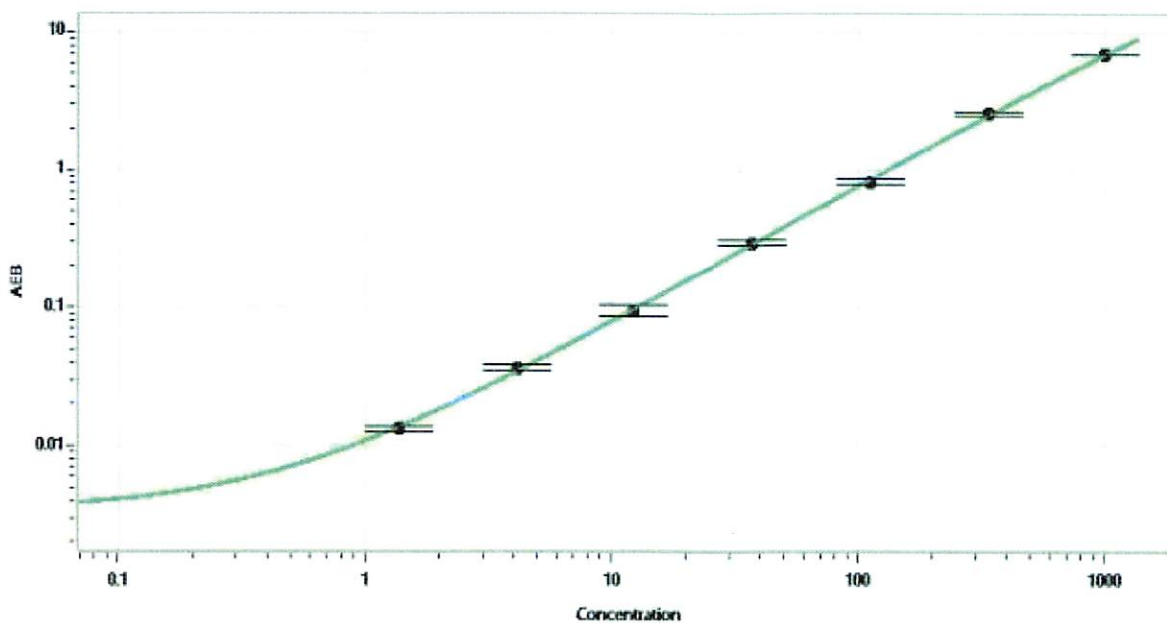


Data below represents results generated on the Simoa® HD-X Analyzer.

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

Calibrator Concentrate Target Value	20
--	----

Calibration Curve



Calibrator Levels (pg/mL)

A	0.000
B	1.37
C	4.11
D	12.3
E	37.0
F	111
G	333
H	1000

Review/Approval

ROBERT SOMOGYI

Name

QC MANAGER

Title

RT S 09 MAY 2024

Signature/Date