

<b>Product Number:</b>	102336
<b>Lot Number:</b>	504192
<b>Expiration:</b>	26-Mar-2025
<b>Platform(s):</b>	SR-X and HD-1/HD-X

Component	Part Number	Lot Number
Bead Stock	102330	331306
Detector Stock	102331	326905
Calibrator Concentrate	102335	331504
Calibrator Diluent	102332	331506
Sample Diluent	102333	331414

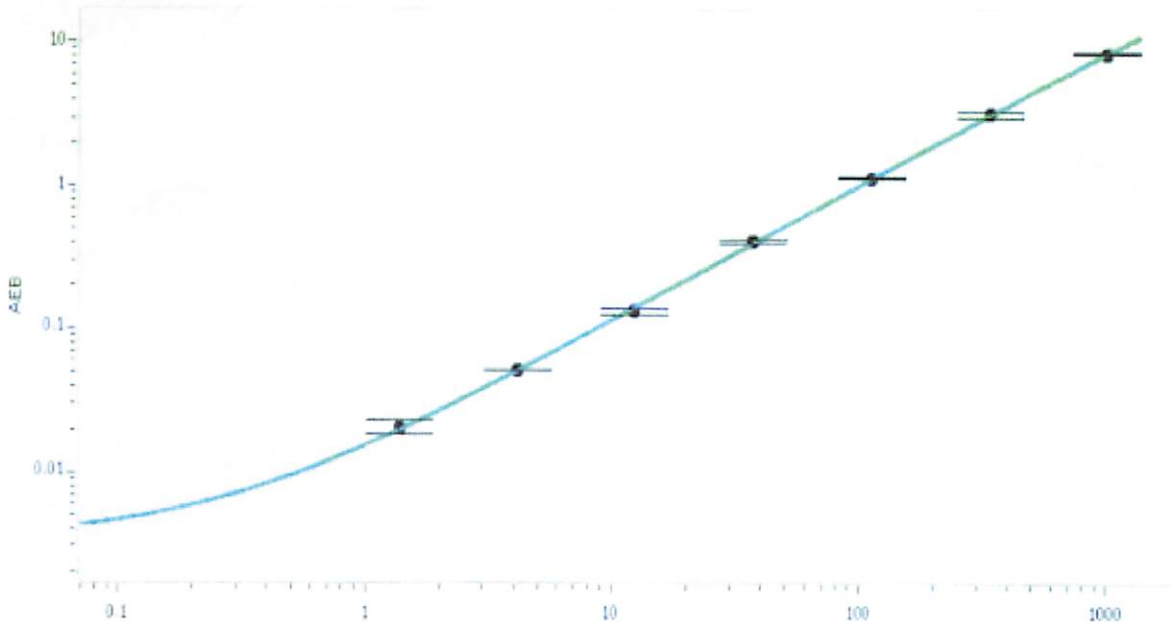
Generic Components - Included With Kit But Not Lot Number Specific	
Discovery Bead Diluent	101954
Discovery Detector Diluent	101955
Discovery SBG Diluent	101956
Discovery SBG Stock	103297
RGP Reagent	103159



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

<b>Calibrator Concentrate Target Value (ng/mL)</b>	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

Cody Carlson

QC Manager

Name

Title

*Cody Carlson* 27 MAR 2024  
Signature/Date