

<b>Product Number:</b>	102039
<b>Lot Number:</b>	504144
<b>Expiration:</b>	10-Aug-2025
<b>Platform(s):</b>	SR-X and HD-1/HD-X

Component	Part Number	Lot Number
Bead Stock	102028	406810
Detector Stock	102029	407208
Calibrator Concentrate	102032	407909
Calibrator Diluent	102030	404110
Sample Diluent	103402	406816

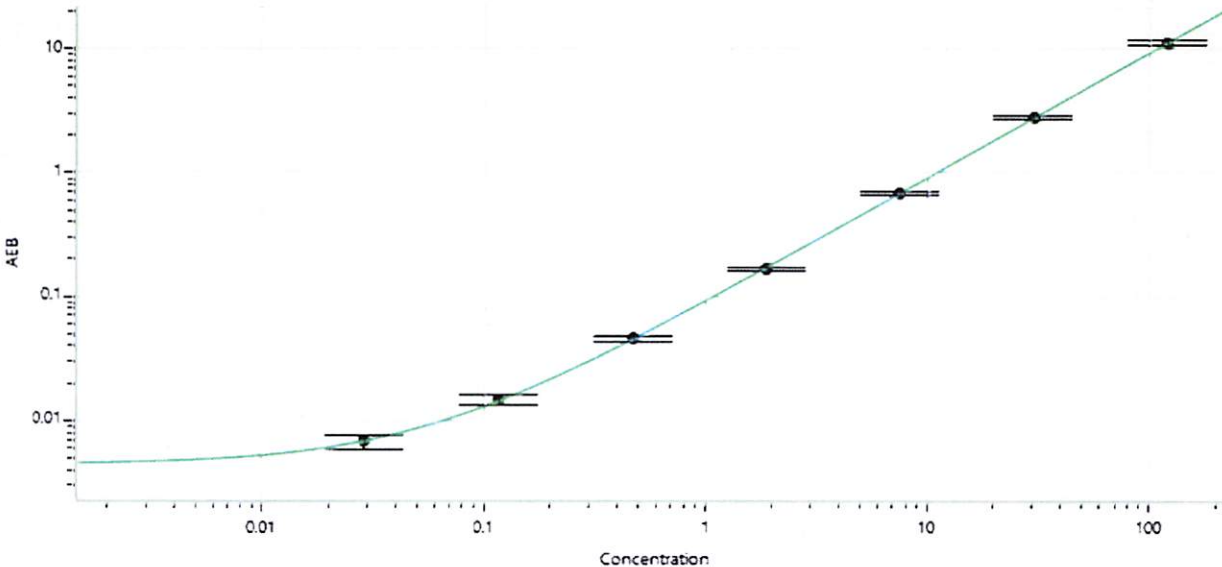
Generic Components - Included With Kit But Not Lot Number Specific		
Discovery Bead Diluent	101954	N/A
Discovery Detector Diluent	101955	N/A
Discovery SBG Diluent	101956	N/A
Discovery SBG Stock	103297	N/A
RGP Reagent	103159	N/A



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

<b>Calibrator Concentrate Target Value (ng/mL)</b>	2
--	---

**Calibration Curve**



**Calibrator Levels (pg/mL)**

A	0.000
B	0.029
C	0.117
D	0.469
E	1.88
F	7.50
G	30.0
H	120

Review/Approval

Cody Carlson

QC Manager

Name

Title

*Cody Carlson* 27 MAR 2024  
Signature/Date