

<b>Product Number:</b>	102336
<b>Lot Number:</b>	504298
<b>Expiration:</b>	2-Mar-2025
<b>Platform(s):</b>	SR-X and HD-1/HD-X

Component	Part Number	Lot Number
Bead Stock	102330	326406
Detector Stock	102331	326905
Calibrator Concentrate	102335	326601
Calibrator Diluent	102332	324511
Sample Diluent	102333	326809

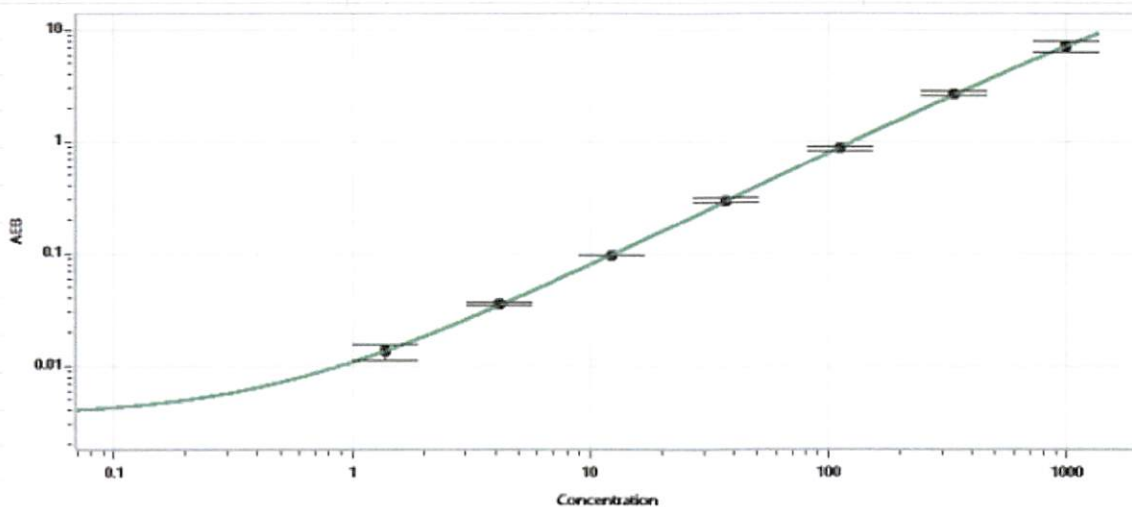


*Data below represents results generated on the Simoa® HD-X Analyzer.*

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

<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

*Dipesh Patel* Name      *Doc control II QA* Title      *DT 16 Oct 2024* Signature/Date