

<b>Product Number:</b>	102860
<b>Lot Number:</b>	504251
<b>Expiration:</b>	17-Oct-2025
<b>Platform(s):</b>	SR-X and HD-1/HD-X



Component	Part Number	Lot Number
Bead Reagent	102852	412002
Detector Reagent	102854	412110
SBG Reagent	102855	412204
Sample Diluent	102856	412309
Calibrator Diluent	102859	412310
Calibrator Concentrate	102857	413405
RGP Reagent	103159	N/A <sup>1</sup>

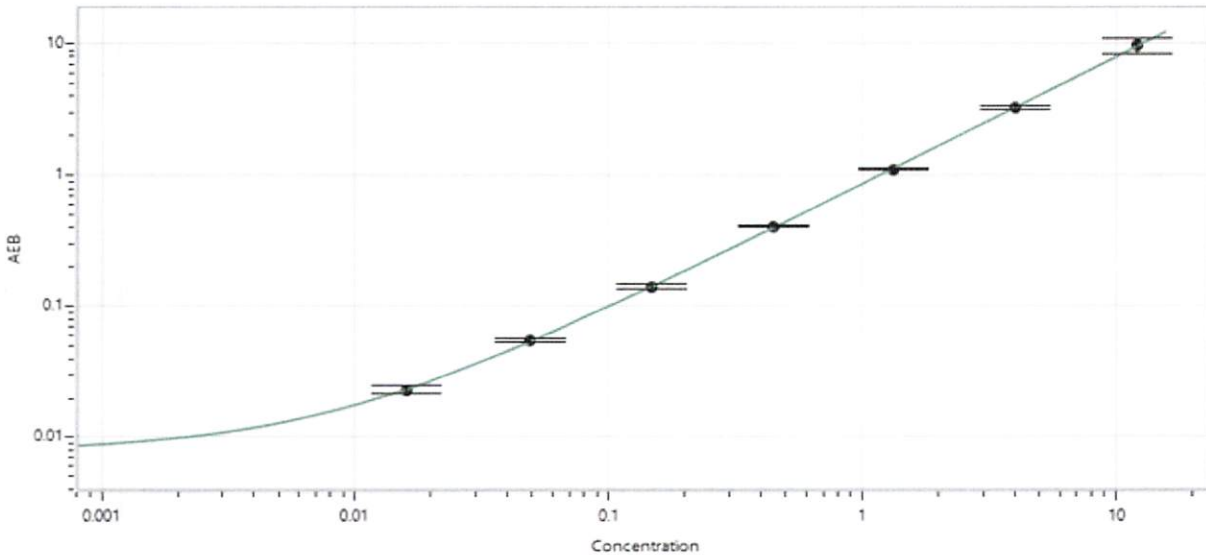
<sup>1</sup> Reagents are not Kit Lot Specific

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

Release Materials*			
	Lot	Result (pg/mL)	Spec. Range (pg/mL)
Control 1	413405	0.222	0.180 - 0.300
Control 2	413405	4.57	3.60 - 6.00
Panel 1	231306	0.253	0.223 - 0.335
Panel 2	234108	3.03	2.40 - 3.75
Panel 3	234108	10.0	8.02 - 13.5

\* Ranges shown are generated internally for new lot release only. Customer should generate their own control ranges.

<b>Calibrator Concentrate Value Assignment (pg/mL)</b>	305
--	-----



Calibrator Levels (pg/mL)	
A	0.000
B	0.016
C	0.049
D	0.148
E	0.444
F	1.33
G	4.00
H	12.0

Review/Approval: Cody Carlson QC Manager  
 Name: Cody Carlson Title: QC Manager Signature/Date: [Signature] 06 JUN 2024