

<b>Product Number:</b>	103464
<b>Lot Number:</b>	504146
<b>Expiration:</b>	19-Jul-2025
<b>Platform(s):</b>	SR-X and HD-1/HD-X



Component	Part Number	Lot Number
Bead Reagent	103452	403808
Detector Reagent	103454	406801
SBG Reagent	103456	406802
Sample Diluent	103458	403814
Calibrator Diluent	103460	403815
Calibrator Concentrate	103459	405307
RGP Reagent	103159	N/A <sup>1</sup>

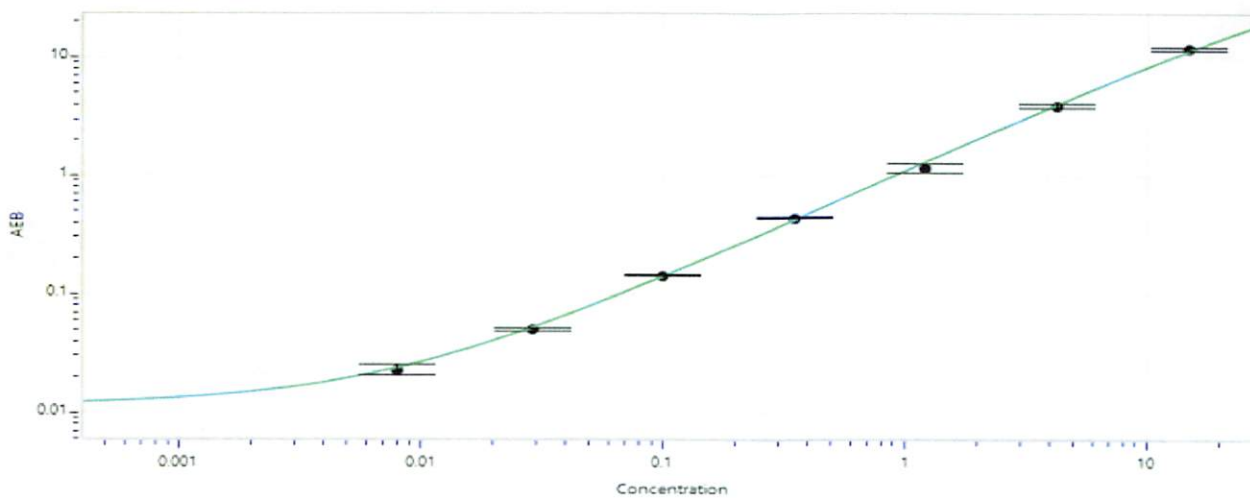
<sup>1</sup> Reagents are not Kit Lot Specific

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

Release Materials*			
	Lot	Result (pg/mL)	Spec. Range (pg/mL)
Control 1	027502	0.211	0.134 - 0.249
Control 2	027502	10.9	7.77 - 11.7
Panel 1	106413	0.098	0.050 - 0.148
Panel 2	106413	0.040	0.019 - 0.076
Panel 3	106413	4.12	3.40 - 5.79
Panel 4	106413	20.5	10.6 - 24.9

\* 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>	356
--	-----



**Calibrator Levels (pg/mL)**

A	0.000
B	0.008
C	0.029
D	0.100
E	0.350
F	1.22
G	4.29
H	15.0

Review/Approval

*Cody Carlson*  
Name Title

*[Signature]* 19 MAR 2024  
Signature/Date