

Product Number:	101599
Lot Number:	504168
Expiration:	13-Oct-2024
Platform(s):	SR-X and HD-1/HD-X



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

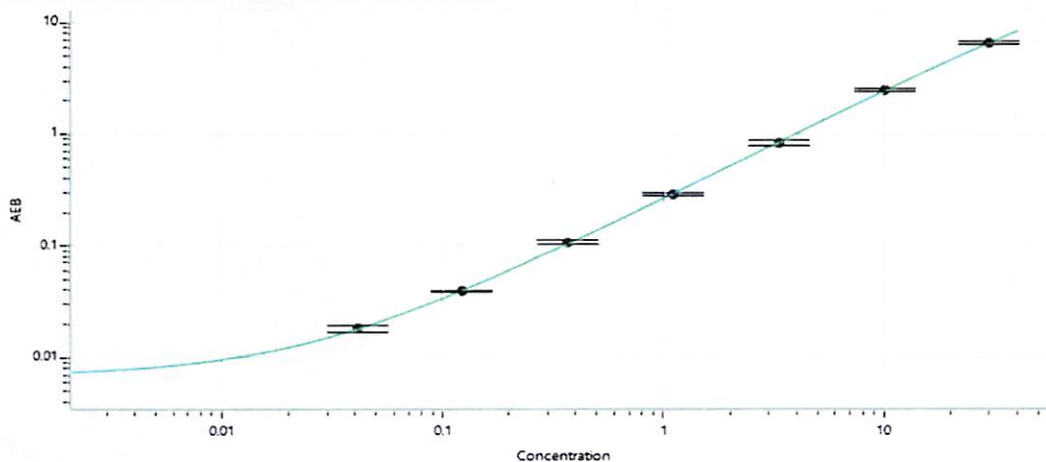
Component	Part Number	Lot Number	Control Lot Number
Bead Reagent	101601	328606	N/A
Detector Reagent	102012	330318	N/A
SBG Reagent	101602	329102	329102A
Sample Diluent	103886	329213	N/A
Calibrator Diluent	102014	328616	N/A
Calibrator Concentrate	100778	329107	N/A
RGP Reagent	103159	N/A ¹	N/A ¹

¹ Reagents are not Kit Lot Specific

Release Materials*			
	Lot	Result (pg/mL)	Spec. Range (pg/mL)
Control 1	209305	0.508	0.432 - 0.749
Control 2	209305	28.2	9.37 - 50.1
Panel 1	321205	2.05	1.53 - 2.29
Panel 2	321205	0.328	0.252 - 0.431
Panel 3	319303	9.21	5.30 - 10.9
Panel 4	319303	29.7	21.2 - 34.4

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

Calibrator Concentrate Value Assignment (pg/mL)	418
--	-----



Calibrator Levels (pg/mL)	
A	0.000
B	0.041
C	0.123
D	0.370
E	1.11
F	3.33
G	10.0
H	30.0

Review/Approval

Cody Carlson QC Manager *Cody Carlson* 08 MAR 2024
 Name Title Signature/Date