

Product Number:	104111
Lot Number:	503537
Expiration:	19-Jul-2023
Platform(s):	HD-1 / HD-X & SR-X



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

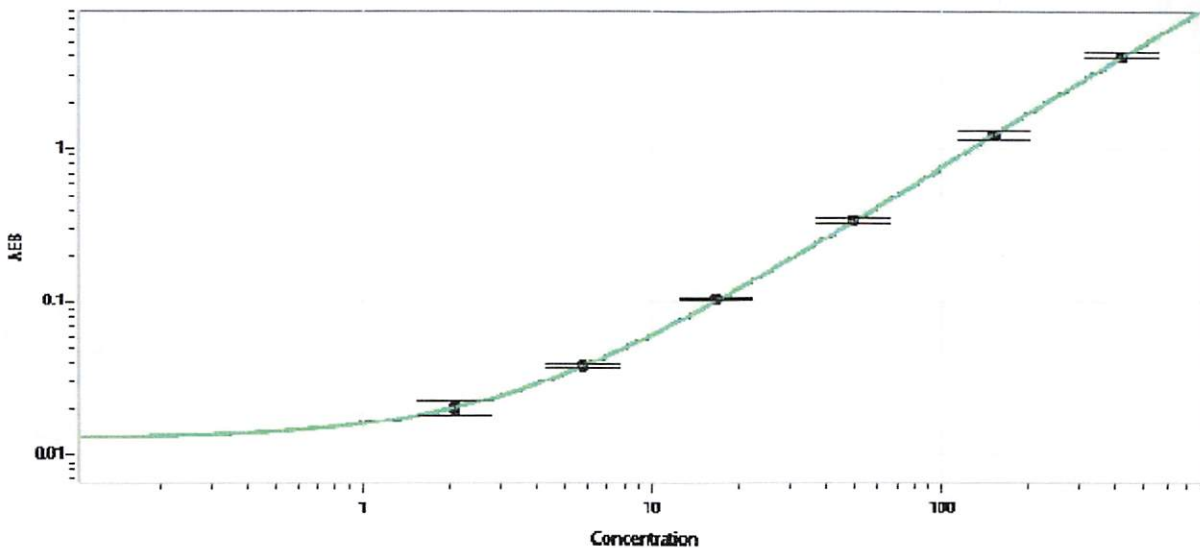
Component	Part Number	Lot Number
Bead Reagent	104098	220002
Detector Reagent	104099	220003
SBG Reagent	104100	220705
Sample Diluent	104101	220005
Reference Calibrator	104106	222002
Control 1	104107	222003
Control 2	104108	222003
RGP Reagent	103159	N/A ¹

¹ RGP is not Kit Lot Specific

Release Materials*			
	Lot	Result (pg/mL)	Spec. Range (pg/mL)
Control 1	222003	55.4	41.5 - 62.3
Control 2	222003	746	539 - 809
Panel 1	214602	23.8	18.7 - 28.0
Panel 2	214602	25.5	18.5 - 28.8
Panel 3	214602	721	544 - 817
Panel 4	214602	242	177 - 265

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

Calibration Curve



Calibrator Levels (pg/mL)	
A	0.00
B	2.07
C	5.76
D	16.7
E	50.0
F	153
G	423

Review/Approval: Sonia Khillan QC Manager Sonia Khillan / 06 OCT 2022
 Name Title Signature/Date