

Product Number:	103186
Lot Number:	503448
Expiration:	14-Apr-2023
Platform(s):	HD-1 / HD-X



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

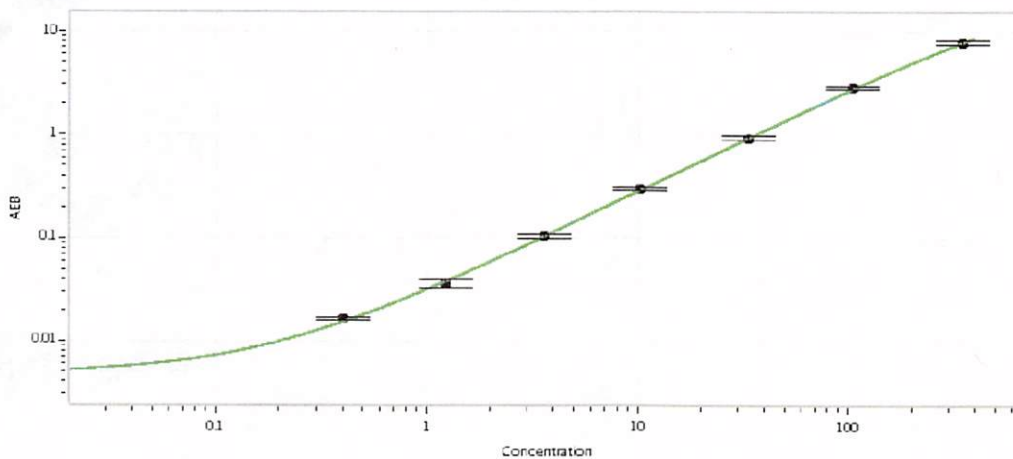
Component	Part Number	Lot Number
Bead Reagent	102246	210408
Detector Reagent	102248	210502
SBG Reagent	102250	210503
Sample Diluent	102252	210504
Calibrators	102255	216601
Control 1	102256	216602
Control 2	102257	216602
RGP Reagent	103159	N/A ¹

¹ RGP is not Kit Lot Specific

Release Materials*			
	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	216602	3.48	2.67-4.16
Control 2	216602	136	102-173
Panel 1	134208	7.42	3.95-9.43
Panel 2	134208	4.89	1.86-7.15
Panel 3	212511	78.8	67.6-116
Panel 4	212511	232	148-361

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

Calibration Curve



Calibrator Levels (pg/mL)	
A	0.000
B	0.400
C	1.24
D	3.63
E	10.2
F	33.5
G	105
H	348

Review/Approval Matt Smart Director, QC [Signature] 30 Jun 2022
Name Title Signature/Date