

Product Number:	103180
Lot Number:	502317
Expiration:	13-Mar-2021
Platform(s):	HD-1/HD-x



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

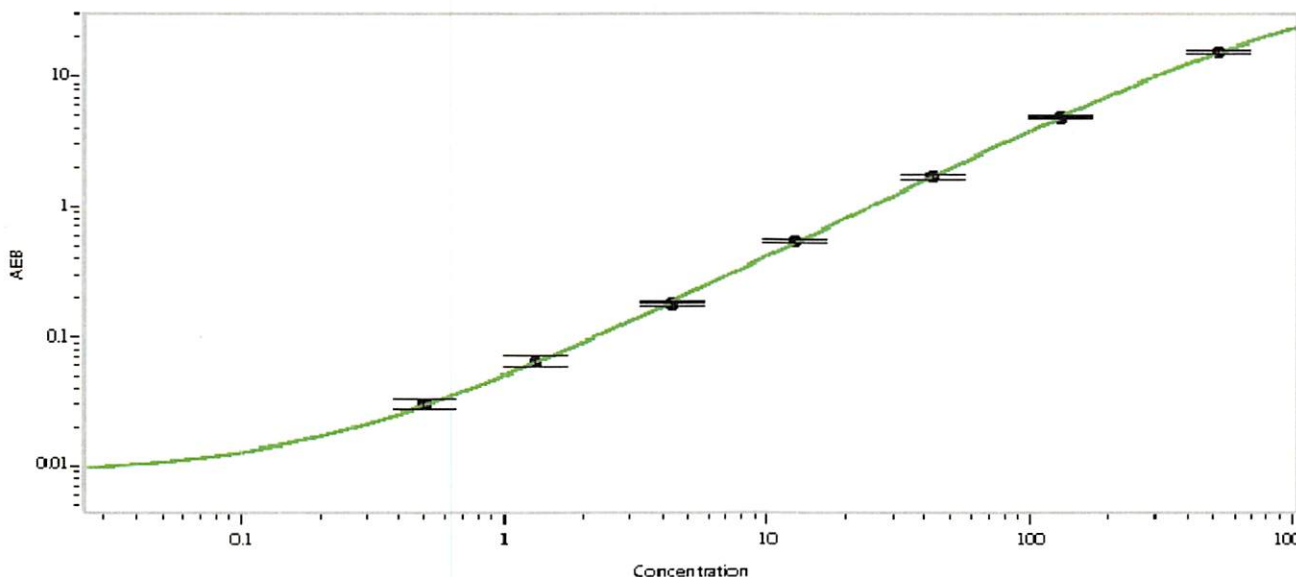
Component	Part Number	Lot Number
Bead Reagent	102246	010001
Detector Reagent	102248	009701
SBG Reagent	102250	009707
Sample Diluent	102252	009702
Calibrators	102255	924701
Controls	102256 - 102257	924803
RGP Reagent	103159	N/A ¹

¹ RGP is not Kit Lot Specific

Release Materials*			
	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	924803	3.36	2.50-4.49
Control 2	924803	136	116-173
Panel 1	922601	17.3	14.1-21.1
Panel 2	922602	5.96	4.94-7.41
Panel 3	922603	39.4	28.4-42.6
Panel 4	922604	158	129-204

* 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	0.503
C	1.33
D	4.32
E	12.8
F	42.3
G	130
H	521

Review/Approval: Patrizia Stadler Name, Assoc. Director, QC Title, P. Stadler Signature, 14 Apr 2020 Date