

Product Number:	103180
Lot Number:	502164
Expiration:	8-Nov-2020
Platform(s):	HD-1/HD-x



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

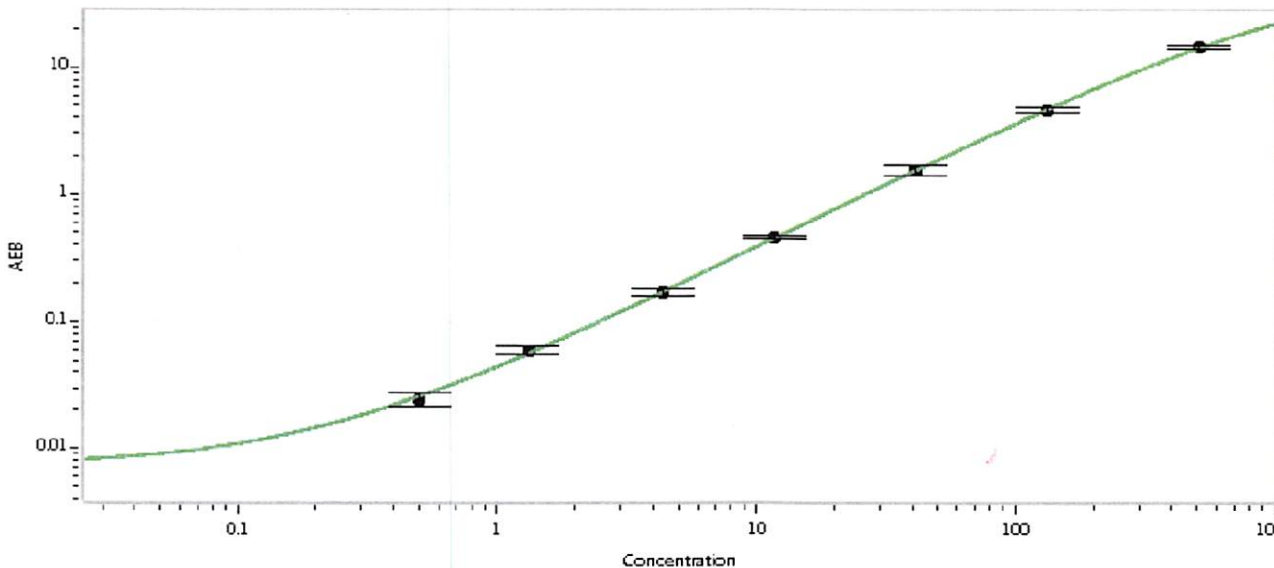
Component	Part Number	Lot Number
Bead Reagent	102246	933607
Detector Reagent	102248	932706
SBG Reagent	102250	931201
Sample Diluent	102252	931203
Calibrators	102255	933601
Controls	102256 - 102257	933901
RGP Reagent	103159	N/A <sup>1</sup>

<sup>1</sup> RGP is not Kit Lot Specific

Release Materials*			
	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	933901	4.10	2.37-4.99
Control 2	933901	147	116-181
Panel 1	922601	18.0	14.1-21.1
Panel 2	922602	6.48	4.94-7.41
Panel 3	922603	40.3	28.4-42.6
Panel 4	922604	169	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	0.501
C	1.32
D	4.32
E	11.8
F	41.1
G	131
H	506

Review/Approval: Patrizia Stadler Assoc. Director, QC p. stadler 12 Dec 2019  
 Name Title Signature/Date