

Product Number:	103186
Lot Number:	502157
Expiration:	10-Dec-2020
Platform(s):	HD-1/HD-x



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

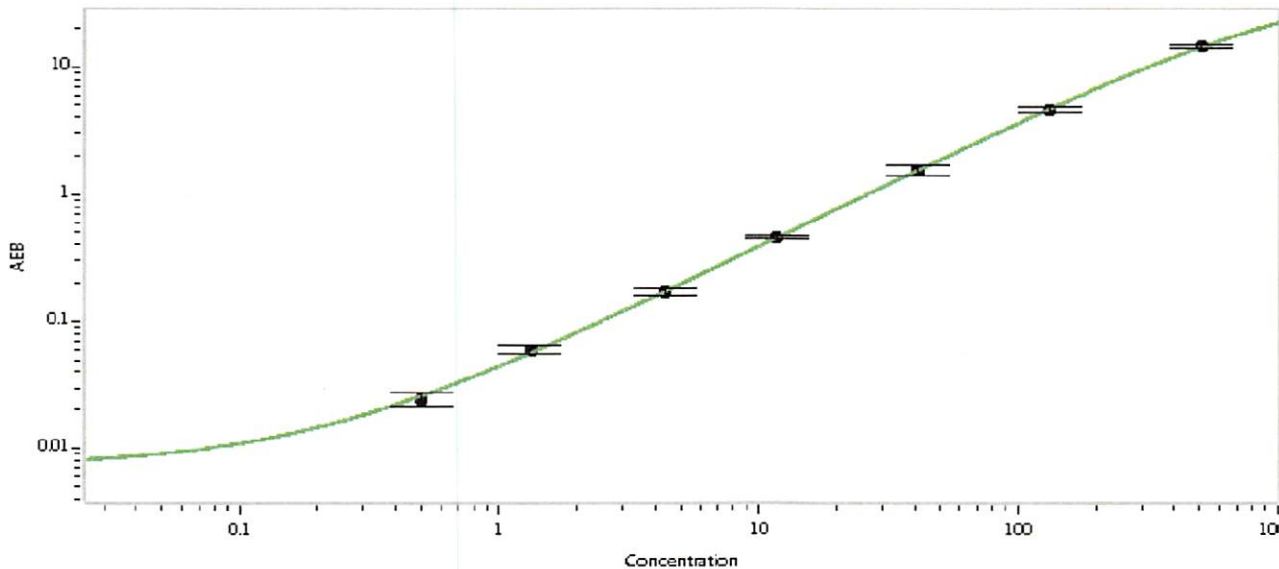
Component	Part Number	Lot Number
Bead Reagent	102246	934507
Detector Reagent	102248	934405
SBG Reagent	102250	934404
Sample Diluent	102252	934403
Calibrators	102255	933601
Controls	102256 - 102257	933901
RGP Reagent	103159	N/A ¹

¹ RGP is not Kit Lot Specific

Release Materials*			
	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	933901	3.94	2.56-4.80
Control 2	933901	154	121-181
Panel 1	922601	18.6	14.1-21.1
Panel 2	922602	6.50	4.94-7.41
Panel 3	922603	38.1	28.4-42.6
Panel 4	922604	166	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	<i>p. Stadler</i>
	Name	Title	Signature/Date
			23 Dec 2019