

Product Number:	103186
Lot Number:	501992
Expiration:	1-Aug-2020
Platform(s):	HD-1/HD-x & SR-X



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

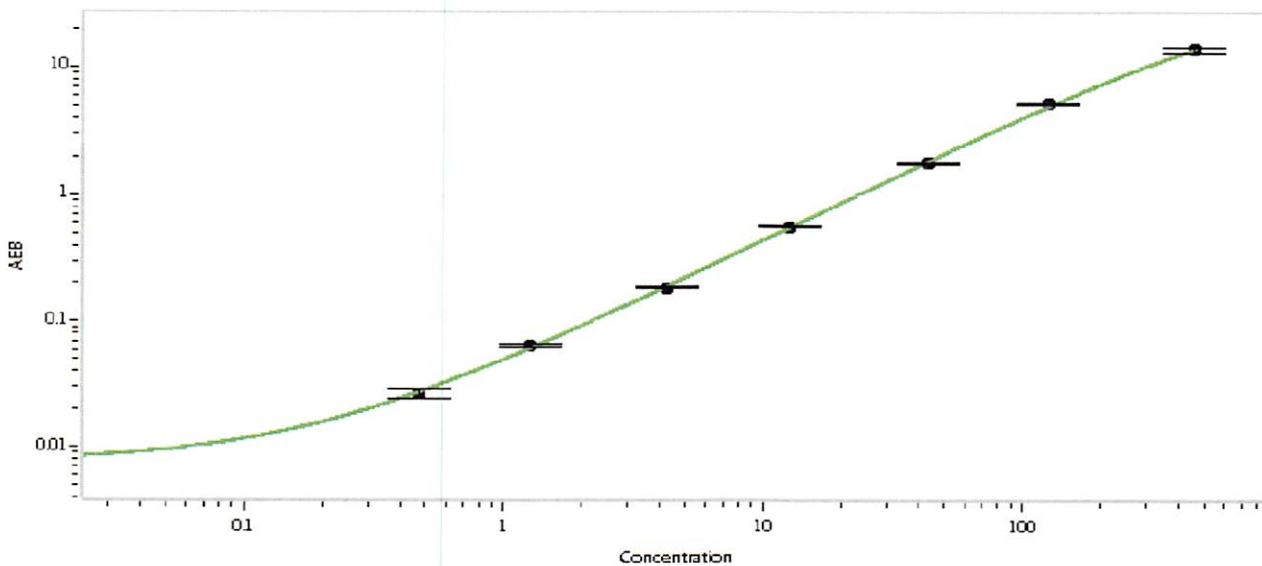
Component	Part Number	Lot Number
Bead Reagent	102246	921706
Detector Reagent	102248	921701
SBG Reagent	102250	921313
Sample Diluent	102252	921702
Calibrators	102255	922701
Controls	102256 - 102257	922702
RGP Reagent	103159	N/A ¹

¹ RGP is not Kit Lot Specific

Release Materials*			
	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	922702	3.91	2.95-5.07
Control 2	922702	148	108-198
Panel 1a	922601	16.7	13.8-20.8
Panel 1b	922602	5.38	4.90-7.36
Panel 2	922603	35.6	27.7-41.6
Panel 3	922604	168	128-206

* 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	0.471
C	1.26
D	4.23
E	12.5
F	42.8
G	126
H	453

Review/Approval: Daniel Ritchie, Senior QC Analyst
 Name: Daniel Ritchie, Title: Senior QC Analyst, Signature/Date: Daniel Ritchie / 23 Aug 2019
 *Signing on behalf of Patrizia Stadler