

Product Number:	103186
Lot Number:	501844
Expiration:	16-Feb-2020
Platform(s):	HD-1 & SR-X

Component	Part Number	Lot Number
Bead Reagent	102246	906503
Detector Reagent	102248	904703
SBG Reagent	102250	904702
Sample Diluent	102252	904701
Reference Calibrators	102255	913501
Controls	102256 - 102257	913601
RGP Reagent	103159	N/A ¹

¹ RGP is not Kit Lot Specific

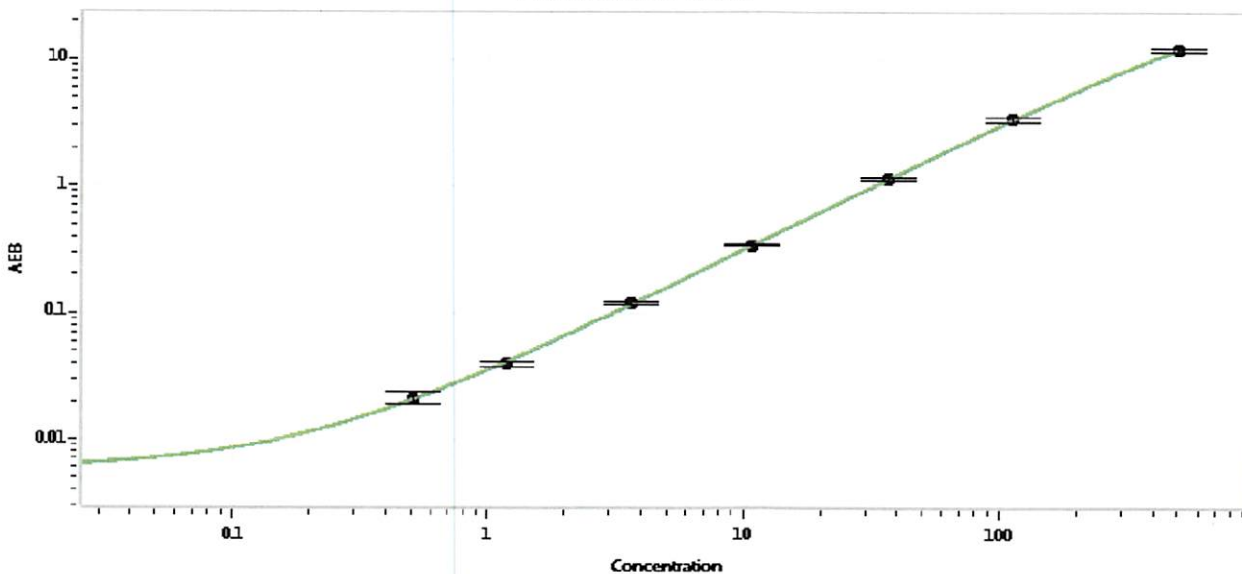


Data below represents results generated on the Simoa™ HD-1 Analyzer.

Release Materials*			
	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	913601	4.70	3.56-5.35
Control 2	913601	196	145-218
Panel 1-A	823301	4.95	4.08-6.68
Panel 1-B	823302	6.19	4.90-7.61
Panel 2	823601	80.4	59.3-101
Panel 3	823602	313	236-402

* 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	0.509
C	1.18
D	3.63
E	10.8
F	36.7
G	112
H	493

Review/Approval: Patrizia Stadler QC Assoc. Director P. Stooler 22/04/2019
Name Title Signature/Date