

Product Number:	101664
Lot Number:	502094
Expiration:	19-Apr-2020
Platform(s):	HD-1/HD-X & SR-X

Component	Part Number	Lot Number
Bead Reagent	101666	910905
Detector Reagent	101665	911703
SBG Reagent	101667	930203
Sample Diluent	101668	930201
Calibrator Diluent	101669	929803
Calibrator Concentrate	103217	912801
RGP Reagent	103159	N/A ¹

¹ RGP is not Kit Lot Specific

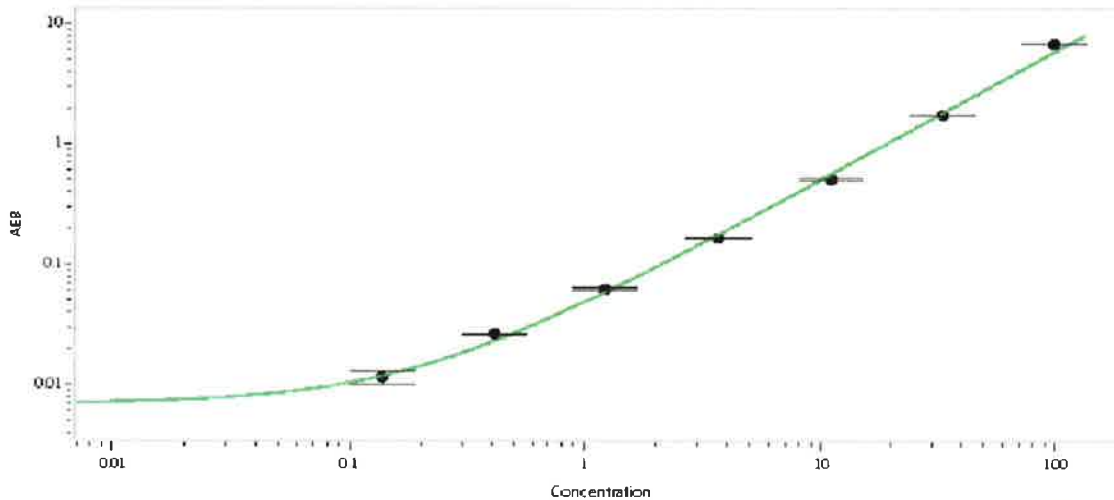


Data below represents results generated on the Simoa™ HD-1 Analyzer.

Release Materials*				* Ranges shown are generated internally for new lot release only. Customer should generate their own control ranges.
	Lot	Result (pg/mL)	Mean Range (pg/mL)	
Control 1	52520	0.915	0.740-1.33	
Control 2	52520	31.9	27.1-40.6	
Panel 1	934301	4.65	2.56-6.16	
Panel 2	934302	18.8	13.6-23.5	
Panel 3	934701	92.0	74.8-112	

Calibrator Concentrate Value Assignment (pg/mL)	1749
---	------

Calibration Curve



Calibrator Levels (pg/mL)	
A	0.000
B	0.137
C	0.412
D	1.23
E	3.70
F	11.1
G	33.3
H	100

Review/Approval

Patrizia Stadler Assoc. Director, QC

P. Stadler

2 Jan 2020

Name

Title

Signature/Date