

Product Number:	103562
Lot Number:	502152
Expiration:	29-Oct-2020
Platform(s):	HD-1/HD-x & SR-X



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

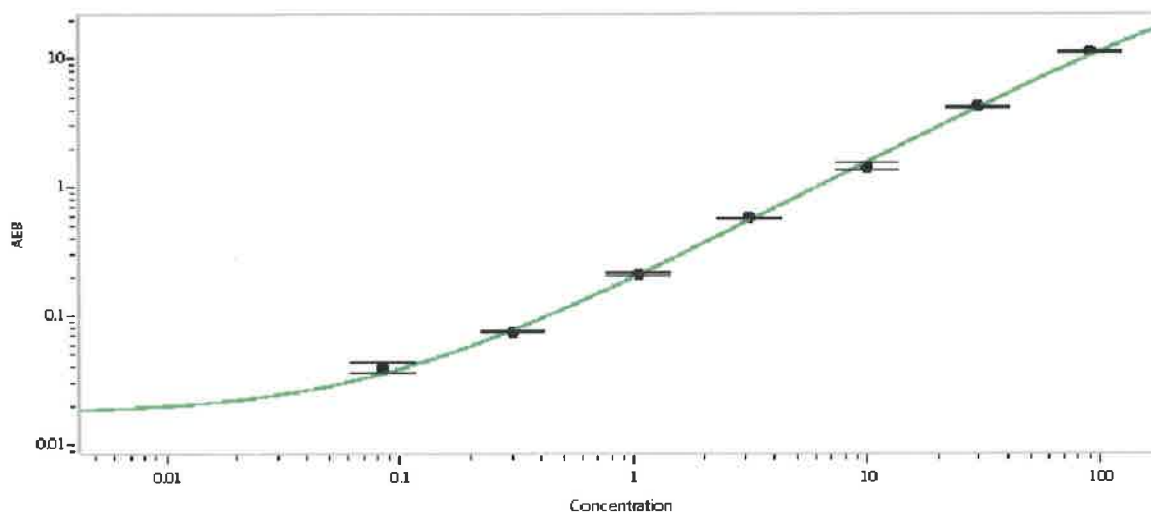
Component	Part Number	Lot Number
Bead Reagent	101554	934411
Detector Reagent	101553	933113
SBG Reagent	101555	933114
Sample Diluent	101556	933116
Calibrators	101559	935401
Controls	101560-101561	935402
SBG Diluent (SR-X)†	103158	N/A ¹
RGP Reagent	103159	N/A ¹

¹ RGP is not Kit Lot Specific

Release Materials*			
	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	935402	1.13	0.686-1.71
Control 2	935402	59.1	47.2-70.8
Panel 1	916901	0.322	0.089-1.59
Panel 2	916902	0.849	0.538-2.13
Panel 3	916903	25.5	15.5-33.4
Panel 4	916904	91.1	63.0-102

* 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	0.084
C	0.303
D	1.05
E	3.13
F	9.96
G	29.7
H	90.7

Review/Approval

Patrizia Stadler Assoc. Director, QC
Name Title

P. Stadler 14 Jan 2020
Signature/Date