

Product Number:	101622
Lot Number:	502077
Expiration:	7-Oct-2020
Platform(s):	HD-1/HD-x & SR-X



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

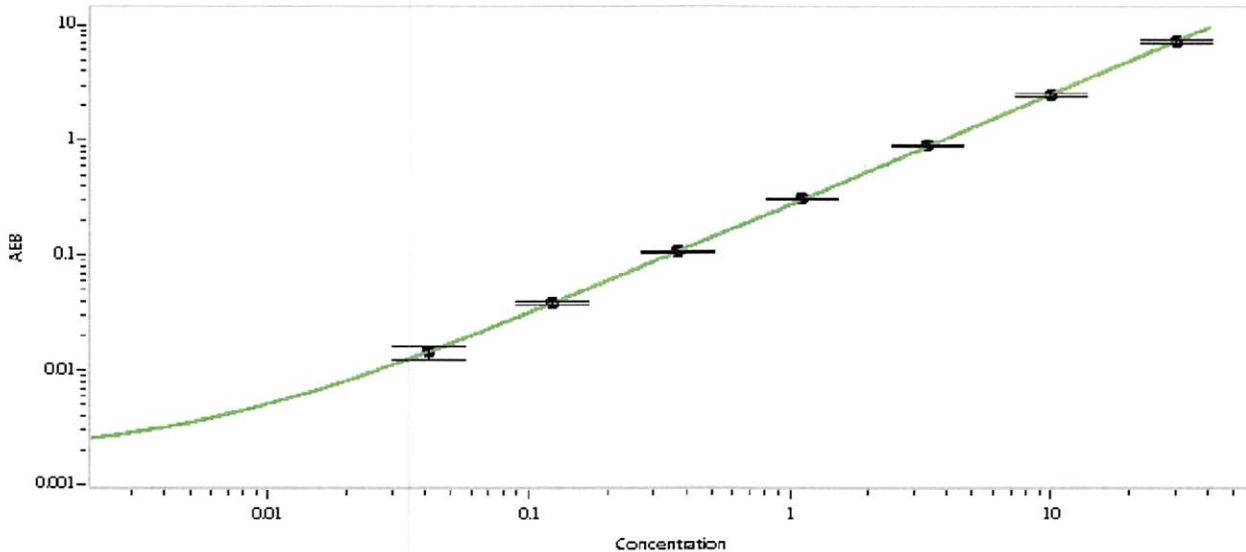
Component	Part Number	Lot Number
Bead Reagent	101624	928906
Detector Reagent	101623	928809
SBG Reagent	101625	928802
Sample Diluent	103203	928021
Calibrator Diluent	101627	928013
Calibrator Concentrate	100652	929802
RGP Reagent	103159	N/A ¹

¹ RGP is not Kit Lot Specific

	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	901702	1.16	0.934-1.40	
Control 2	901702	20.5	15.1-22.7	
Panel 1	MF-050918-01	60.1	41.6-67.5	
Panel 2	MF-050918-02	8.24	5.03-8.84	
Panel 3	MF-050918-03	1.18	0.644-1.24	

Calibrator Concentrate Value Assignment (pg/mL)	673
---	-----

Calibration Curve



Calibrator Levels (pg/mL)	
A	0.000
B	0.041
C	0.123
D	0.370
E	1.11
F	3.33
G	10.0
H	30.0

Review/Approval

Patrizia Stadler Assoc. Director, QC
Name Title

P. Stadler 26 Nov 2019
Signature/Date