

Product Number:	102215
Lot Number:	501601
Expiration:	10-Sep-2019
Platform(s):	HD-1 & SR-X



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

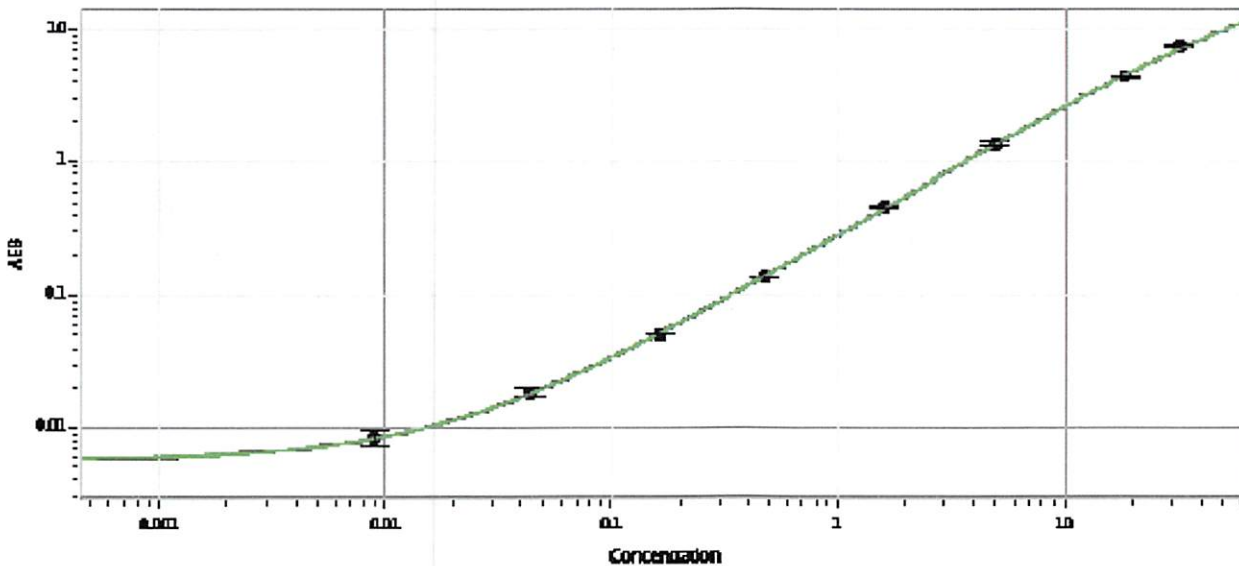
Component	Part Number	Lot Number
Bead Reagent	102211	826307
Detector Reagent	102212	901704
SBG Reagent	102213	825309
Assay Diluent	102214	825408
Calibrators	101310	829501
Controls	100702-100703	829502
RGP Reagent	103159	N/A ¹

¹ RGP is not Kit Lot Specific

Release Materials*			
	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	829502	0.167	0.118-0.206
Control 2	829502	19.0	14.2-22.4
Panel 1	MF-092217-01	0.291	0.177-0.353
Panel 2	MF-092217-02	2.89	1.99-3.51
Panel 3	MF-101117-01	19.0	12.8-20.8

* 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 0.009
C 0.044
D 0.165
E 0.472
F 1.60
G 4.97
H 18.8
I 32.5

Review/Approval

Patrizia Stadler

Associate Director, QC

P. Stadler

28 Jun 2019

Name

Title

Signature/Date