

Product Number:	103714
Lot Number:	502582
Expiration:	13-Apr-2021
Platform(s):	HD-1 / HD-X & SR-X



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

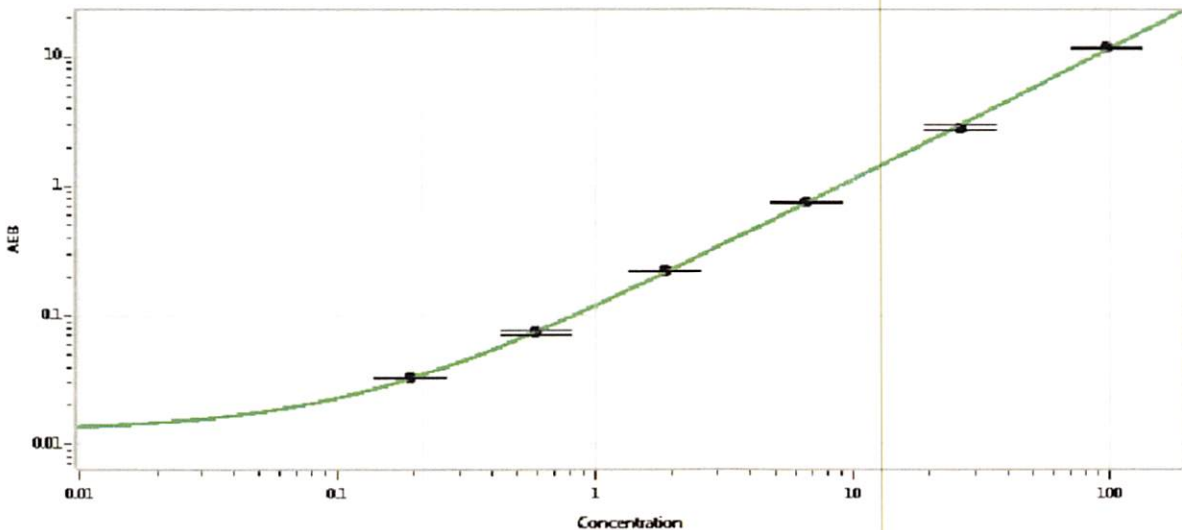
Component	Part Number	Lot Number
Bead Reagent	103698	028710
Detector Reagent	103700	028906
SBG Reagent	103702	028904
Sample Diluent	103704	028903
Reference Calibrator	103711	029302
Control 1	103712	029303
Control 2	103713	029303
RGP Reagent	103159	N/A ¹

¹ RGP is not Kit Lot Specific

Release Materials*			
	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	029303	3.81	2.43-4.64
Control 2	029303	60.0	36.2-67.5
Panel 1	013903	0.801	0.530-0.934
Panel 2	013903	1.84	1.28-2.07
Panel 3	013903	2.69	1.61-4.72

* 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	0.193
C	0.589
D	1.87
E	6.57
F	26.0
G	96.0

Review/Approval

Shivani Goel
Name

QC Manager
Title

Shivani Goel 29 Oct 2020
Signature/Date