

Product Number:	103714
Lot Number:	502508
Expiration:	6-Jan-2021
Platform(s):	HD-1 / HD-X & SR-X



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

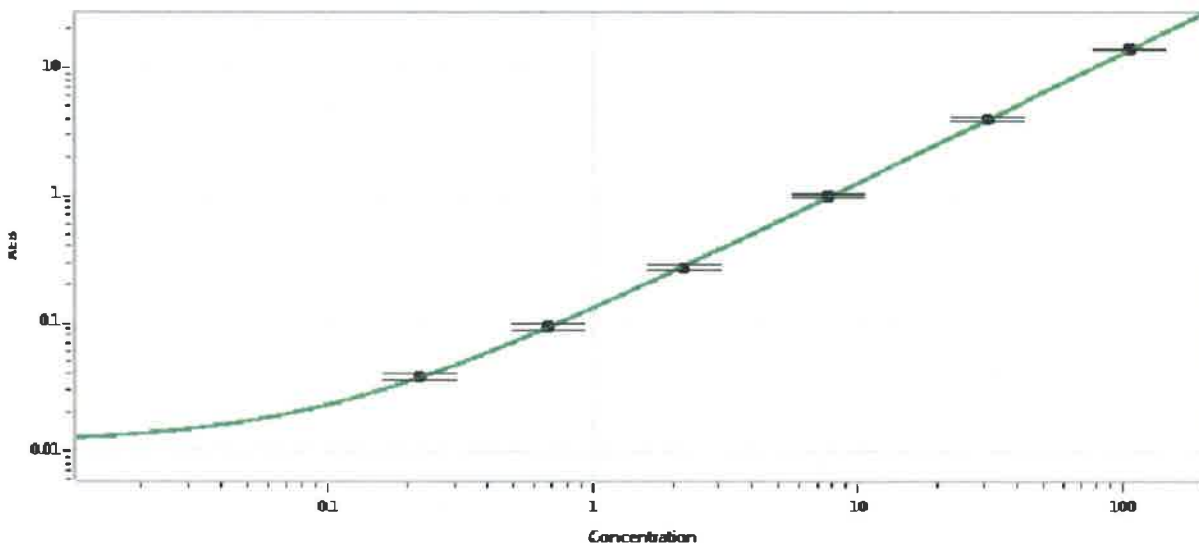
Component	Part Number	Lot Number
Bead Reagent	103698	018804
Detector Reagent	103700	018902
SBG Reagent	103702	018803
Sample Diluent	103704	019705
Reference Calibrator	103711	023201
Control 1	103712	021901
Control 2	103713	021901
RGP Reagent	103159	N/A ¹

¹ RGP is not Kit Lot Specific

Release Materials*			
	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	021901	3.82	2.98-4.47
Control 2	021901	62.2	48.8-75.1
Panel 1	013903	0.769	0.575-0.916
Panel 2	013903	1.68	1.35-2.04
Panel 3	013903	2.68	1.73-4.89

* 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	0.222
C	0.678
D	2.19
E	7.77
F	30.9
G	106

Review/Approval

Shivani Goel
Name

QC Manager
Title

Shivani Goel 08 Sep 2020
Signature/Date

