

Product Number:	101622
Lot Number:	502570
Expiration:	3-Aug-2021
Platform(s):	HD-1 / HD-X & SR-X

Component	Part Number	Lot Number
Bead Reagent	101624	021704
Detector Reagent	101623	021804
SBG Reagent	101625	028108
Sample Diluent	103732	028106
Calibrator Diluent	101627	024716
Calibrator Concentrate	100652	005601
RGP Reagent	103159	N/A ¹

¹ RGP is not Kit Lot Specific



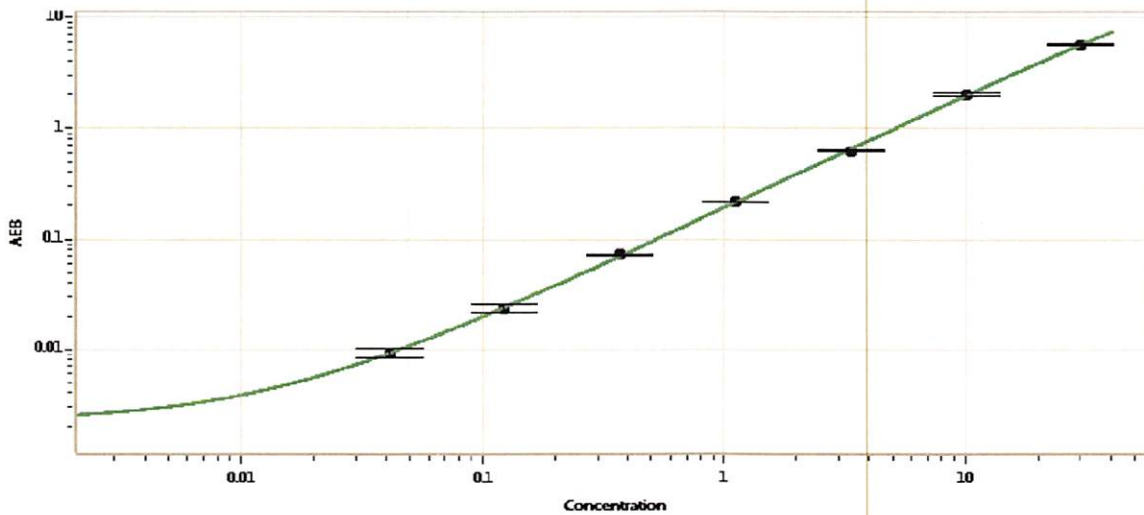
Data below represents results generated on the Simoa™ HD-1 Analyzer.

Release Materials*			
	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	901702	1.29	0.944-1.42
Control 2	901702	18.2	15.4-23.1
Panel 1	029301	52.4	40.2-60.3
Panel 2	029401	5.53	4.33-6.50
Panel 3	029301	1.60	1.21-1.81

* Ranges shown are generated internally for new lot release only. Customer should generate their own control ranges.

Calibrator Concentrate Value Assignment (pg/mL)	645
---	-----

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

Shivani Goel

QC Manager

Shivani Goel 30 Oct 2020

Name

Title

Signature/Date