

Product Number:	101580
Lot Number:	503142
Expiration:	6-Aug-2022
Platform(s):	HD-1/HD-x & SR-X

Component	Part Number	Lot Number
Bead Reagent	101582	125807
Detector Reagent	101581	126001
SBG Reagent	101583	126002
Sample Diluent	103735	123309
Calibrator Diluent	101585	126004
Calibrator Concentrate	100655	126005
RGP Reagent	103159	N/A <sup>1</sup>

<sup>1</sup> RGP is not Kit Lot Specific

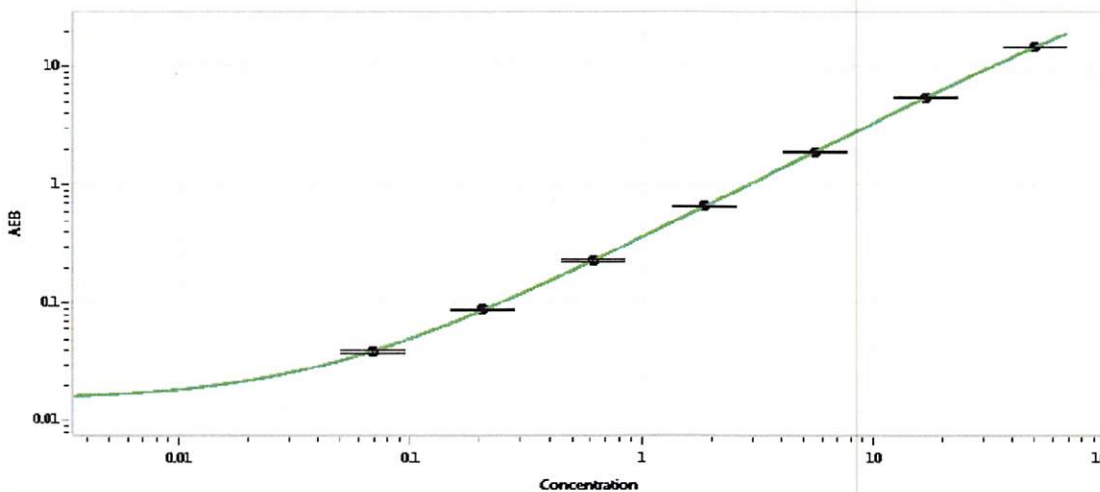


Data below represents results generated on the Simoa™ HD-1 Analyzer.

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	024603	1.33	0.594 - 2.32	
Control 2	024603	23.6	13.9 - 36.9	
Panel 2	115317	2.33	1.37 - 3.90	
Panel 3	116301	27.3	14.8 - 39.9	

Calibrator Concentrate Value Assignment (pg/mL)	1120
---	------

Calibration Curve



Calibrator Levels (pg/mL)	
A	0.000
B	0.069
C	0.206
D	0.617
E	1.85
F	5.56
G	16.7
H	50.0

Review/Approval

Junpil Park  
Name

QC Supervisor  
Title

*[Signature]* 23 Sep 2021  
Signature/Date