

Product Number:	103575
Lot Number:	503115
Expiration:	10-Jun-2022
Platform(s):	HD-1 / HD-X & SR-X

Component	Part Number	Lot Number
Bead Reagent	103564	123305
Detector Reagent	103566	123407
SBG Reagent	103568	123408
Sample Diluent	103569	117923
Calibrator Diluent	103571	120411
Calibrator Concentrate	103570	123503
SR-X SBG Diluent	103158	N/A ¹
RGP Reagent	103159	N/A ¹

¹ Generic items are not Kit Lot Specific



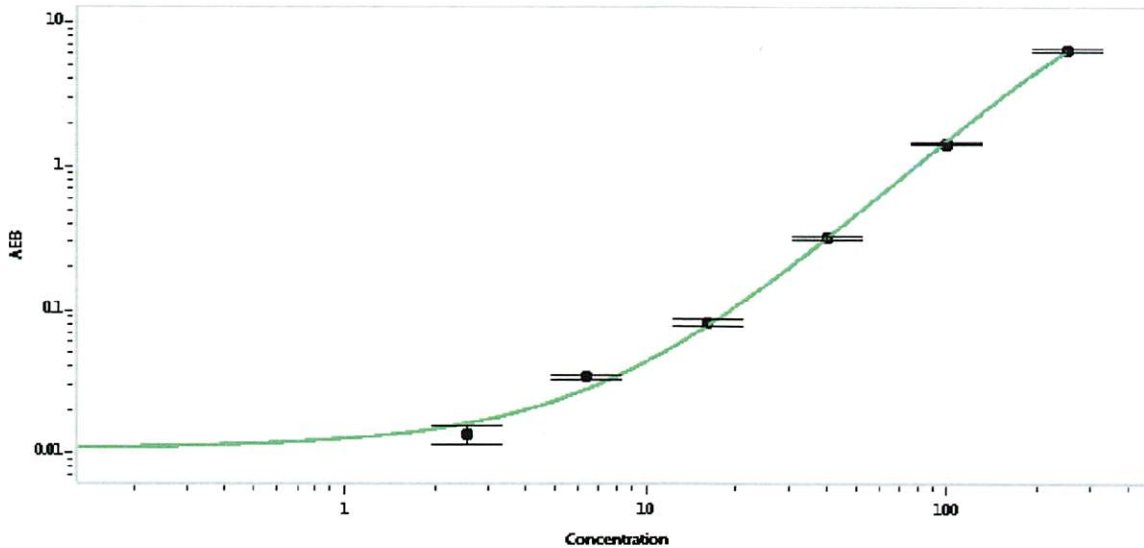
Data below represents results generated on the Simoa™ HD-1 Analyzer.

Release Materials*			
	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	023306	49.3	39.2 - 58.8
Control 2	023306	432	333 - 500
Panel 1	020204	59.0	47.2 - 70.8
Panel 2	020204	72.1	58.4 - 87.7
Panel 3	102806	536	390 - 586

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

Calibrator Concentrate Value Assignment (pg/mL)	4669
---	------

Calibration Curve



Calibrator Levels (pg/mL)	
A	0.000
B	2.56
C	6.40
D	16.0
E	40.0
F	100
G	250

Review/Approval

Junpil Park
Name

QC Supervisor
Title

[Signature] 02 Sep 2021
Signature/Date