

Product Number:	103575
Lot Number:	502783
Expiration:	26-Jan-2022
Platform(s):	HD-1 / HD-X & SR-X

Component	Part Number	Lot Number
Bead Reagent	103564	105803
Detector Reagent	103566	107010
SBG Reagent	103568	106002
Sample Diluent	103569	105515
Calibrator Diluent	103571	105615
Calibrator Concentrate	103570	106802
SR-X SBG Diluent	103158	N/A ¹
RGP Reagent	103159	N/A ¹

¹ Generic items are not Kit Lot Specific



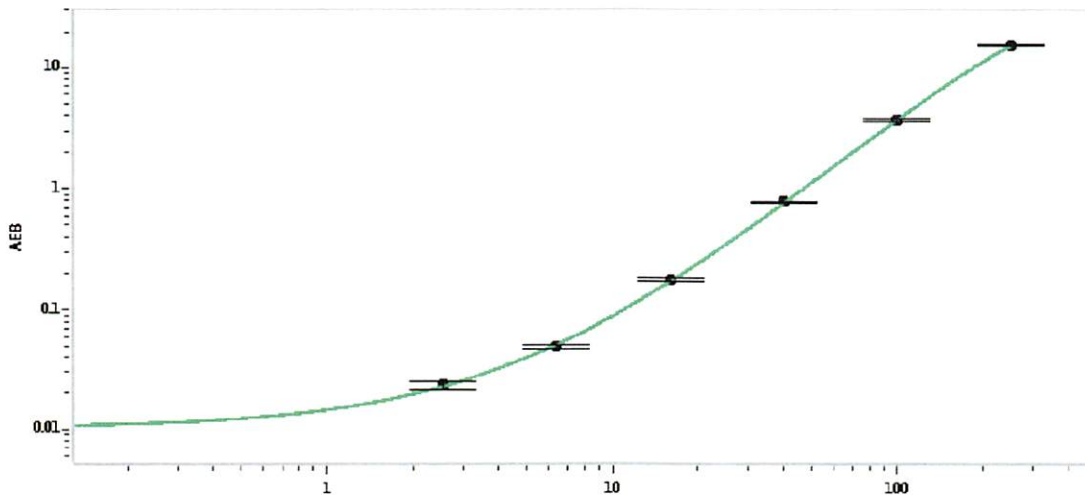
Data below represents results generated on the Simoa™ HD-1 Analyzer.

Release Materials*			
	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	023306	58.1	39.1-58.7
Control 2	023306	449	329-494
Panel 1	020204	64.8	47.2-70.7
Panel 2	020204	85.3	58.7-88.0
Panel 3	102806	484	378-566

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

Calibrator Concentrate Value Assignment (pg/mL)	4185
---	------

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

08 Apr 2021
Date

Signature/Date