

Product Number:	103575
Lot Number:	503385
Expiration:	5-Dec-2022
Platform(s):	HD-1 / HD-X & SR-X

Component	Part Number	Lot Number
Bead Reagent	103564	207501
Detector Reagent	103566	207607
SBG Reagent	103568	207608
Sample Diluent	103569	133940
Calibrator Diluent	103571	204615
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



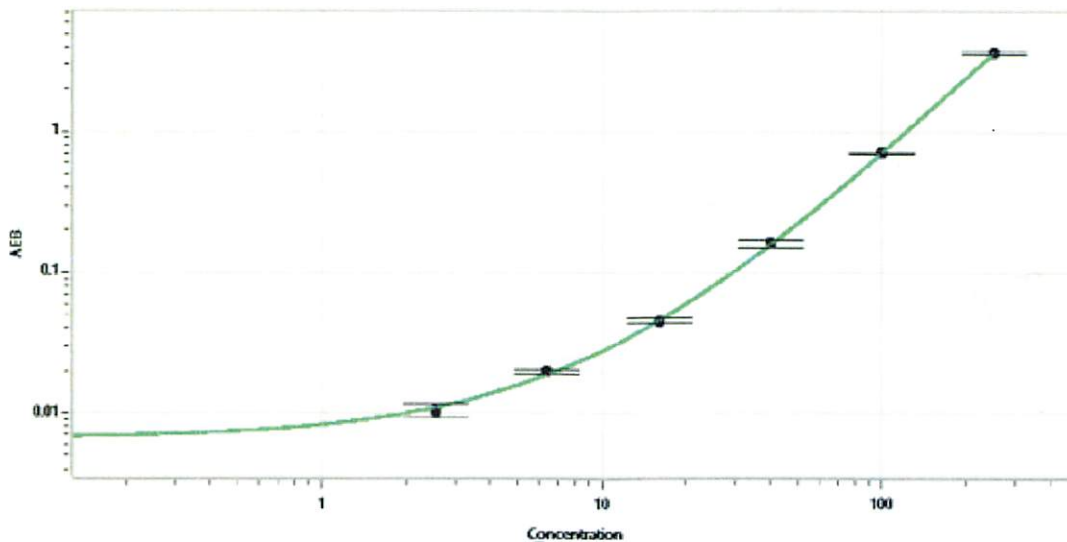
Data below represents results generated on the Simoa® HD-1 / HD-X & SR-X Analyzer.

Release Materials*			
	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	023306	56.8	40.4-60.7
Control 2	023306	494	344-521
Panel 1	020204	65.5	48.2-72.3
Panel 2	020204	80.9	59.5-89.3
Panel 3	102806	613	363-663

* 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: Shivani Goel QC Manager Shivani Goel 11 Apr 2022
Name Title Signature/Date