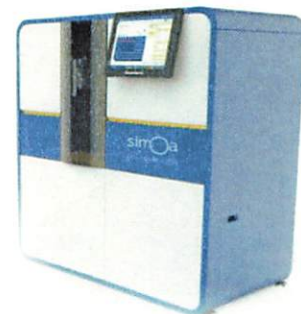


Product Number:	103714
Lot Number:	503057
Expiration:	15-May-2022
Platform(s):	HD-1 / HD-X & SR-X



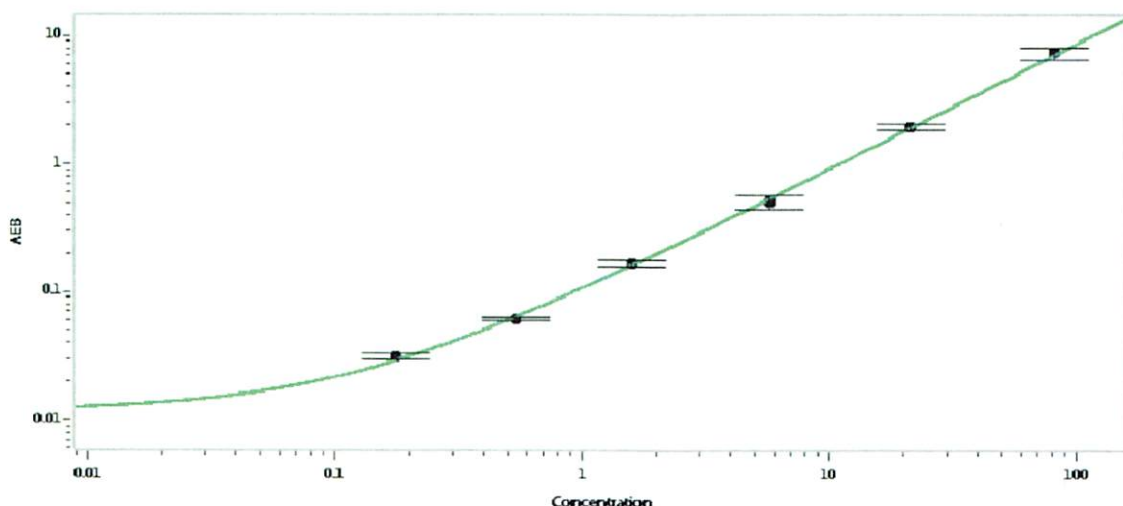
Data below represents results generated on the Simoa™ HD-1 Analyzer.

Component	Part Number	Lot Number
Bead Reagent	103698	119601
Detector Reagent	103700	119703
SBG Reagent	103702	119704
Sample Diluent	103704	119705
Reference Calibrator	103711	120006
Control 1	103712	120007
Control 2	103713	120007
RGP Reagent	103159	N/A ¹

¹ RGP is not Kit Lot Specific

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	120007	4.06	3.08 - 4.62	
Control 2	120007	152	91.9 - 172	
Panel 1	013903	0.779	0.258 - 1.11	
Panel 2	013903	1.72	0.840 - 2.44	
Panel 3	013903	2.74	1.79 - 4.31	

Calibration Curve



Calibrator Levels (pg/mL)	
A	0.000
B	0.177
C	0.540
D	1.59
E	5.76
F	21.4
G	81.7

Review/Approval

Shivani Goel
Name

QC

Title

Shivani Goel 10 Mar 2022
Signature/Date