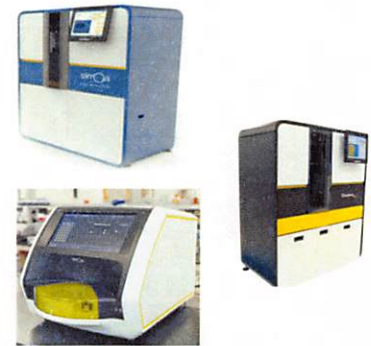


Product Number:	101552
Lot Number:	503307
Expiration:	19-Feb-2023
Platform(s):	HD-1/HD-x & SR-X



Component	Part Number	Lot Number
Bead Reagent	101554	205402
Detector Reagent	101553	205405
SBG Reagent	101555	205406
Sample Diluent	101556	205407
Calibrators	101559	206805
Control 1	101560	206806
Control 2	101561	206806
SBG Diluent (SR-X) †	103158	N/A <sup>1</sup>
RGP Reagent	103159	N/A <sup>1</sup>

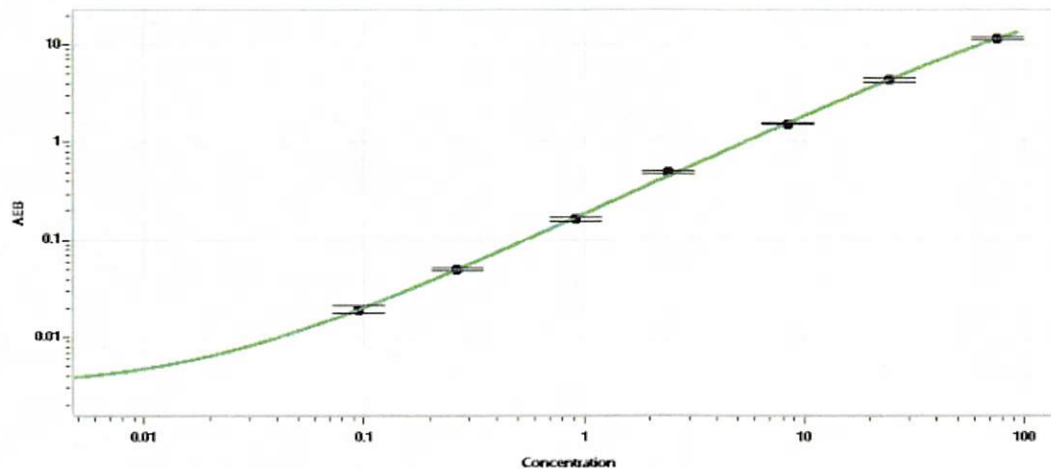
Data below represents results generated on the Simoa® HD-X / HD-1 & SR-X Analyzer.

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

Release Materials*			
	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	206806	0.979	0.731-1.10
Control 2	206806	47.3	39.7-59.6
Panel 1	117205	4.41	3.49-5.23
Panel 2	117205	2.83	1.89-4.80
Panel 3	117205	21.7	17.8-26.7
Panel 4	119612	34.8	23.3-41.8

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

Calibration Curve



Calibrator Levels (pg/mL)

A	0.000
B	0.095
C	0.266
D	0.916
E	2.43
F	8.39
G	24.3
H	75.3

Review/Approval

Shivani Goel  
Name

QC Manager  
Title

Shivani Goel  
Signature/Date 13Apr2022