

Product Number:	102215
Lot Number:	503353
Expiration:	15-Feb-2023
Platform(s):	HD-1/HD-x & SR-X



Component	Part Number	Lot Number
Bead Reagent	102211	204808
Detector Reagent	102212	204903
SBG Reagent	102213	204904
Assay Diluent	102214	204905
Calibrators	101310	116506
Control 1	100702	116507
Control 2	100703	116507
RGP Reagent	103159	N/A ¹

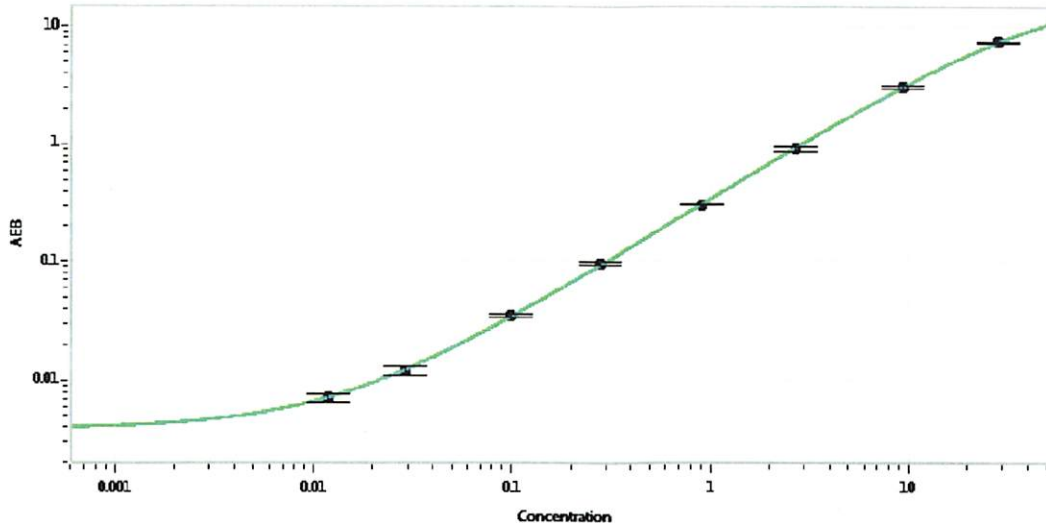
Data below represents results generated on the Simoa® HD-1/ HD-X & SR-X Analyzer.

¹ RGP is not Kit Lot Specific

Release Materials*			
	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	116507	0.097	0.077-0.115
Control 2	116507	9.20	7.30-10.9
Panel 1	019601	0.204	0.127-0.243
Panel 2	019601	1.94	1.56-2.34
Panel 3	019601	11.1	9.44-14.2

* 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	0.012
C	0.029
D	0.099
E	0.280
F	0.903
G	2.67
H	9.35
I	28.2

Review/Approval Shivani Goel QC Manager Shivani Goel 23 Mar 2022
 Name Title Signature/Date