

Product Number:	102215
Lot Number:	502464
Expiration:	27-May-2021
Platform(s):	HD-1/HD-x & SR-X



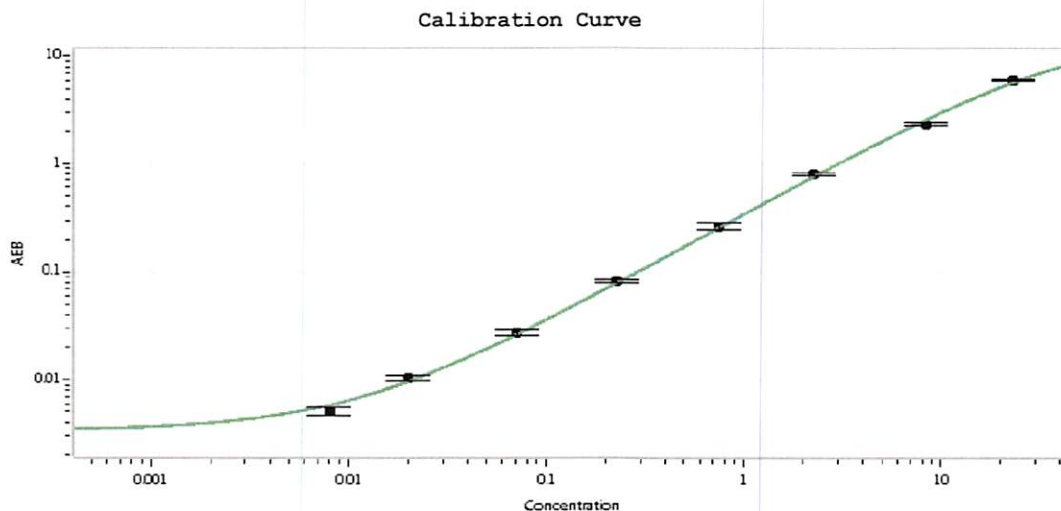
Data below represents results generated on the Simoa™ HD-1 Analyzer.

Component	Part Number	Lot Number
Bead Reagent	102211	019205
Detector Reagent	102212	019201
SBG Reagent	102213	019501
Sample Diluent	102214	019502
Calibrators	101310	004401
Control 1	100702	004402
Control 2	100703	004402
RGP Reagent	103159	N/A ¹

¹ RGP is not Kit Lot Specific

Release Materials*			
	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	004402	0.076	0.058-0.087
Control 2	004402	7.56	6.17-9.26
Panel 1	019601	0.191	0.142-0.212
Panel 2	019601	2.06	1.51-2.27
Panel 3	019601	11.6	9.43-14.1

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



Calibrator Levels (pg/mL)

A	0.000
B	0.008
C	0.020
D	0.071
E	0.225
F	0.754
G	2.28
H	8.33
I	22.8

Review/Approval	Shivani Goel	QC Manager	Shivani Goel 03 Aug 2020
	Name	Title	Signature/Date