

Product Number:	101235
Lot Number:	503103
Expiration:	5-Aug-2022
Platform(s):	HD-1/HD-X



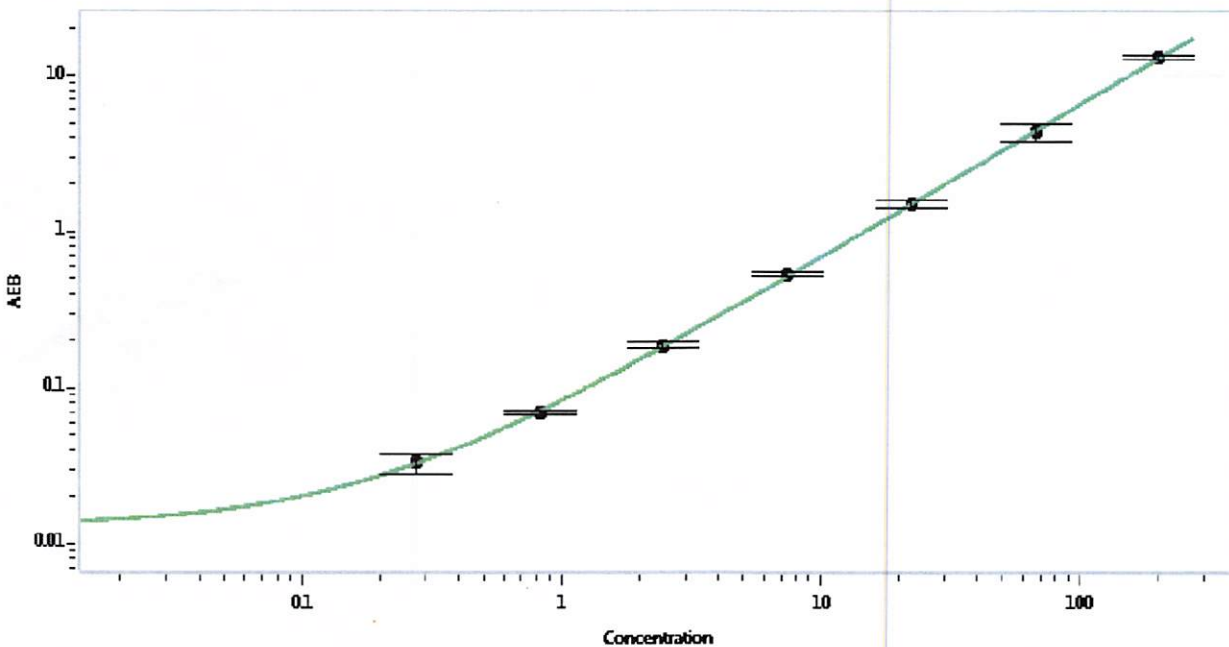
Data below represents results generated on the Simoa™ HD-1 Analyzer.

Component	Part Number	Lot Number
Bead Reagent	101222	122606
Detector Reagent	101224	122701
SBG Reagent	101234	122702
Sample Diluent	101232	122703
Calibrator Diluent	101228	121715
Calibrator Concentrate	101226	122704
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	122704	15.7	11.3 - 18.8
Control 2	122704	396	300 - 500
Panel 1	MF-081516-01	5.81	5.14 - 10.6
Panel 2	MF-081516-02	55.4	47.5 - 95.2
Panel 3	MF-081516-03	2.53	1.14 - 5.12

* 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	0.274
C	0.823
D	2.47
E	7.41
F	22.2
G	66.7
H	200

Review/Approval

Junpil Park QC Supervisor *[Signature]* 27 Aug 2021
 Name Title Signature/Date