

Product Number:	103186
Lot Number:	502907
Expiration:	23-Apr-2022
Platform(s):	HD-1/ HD-X



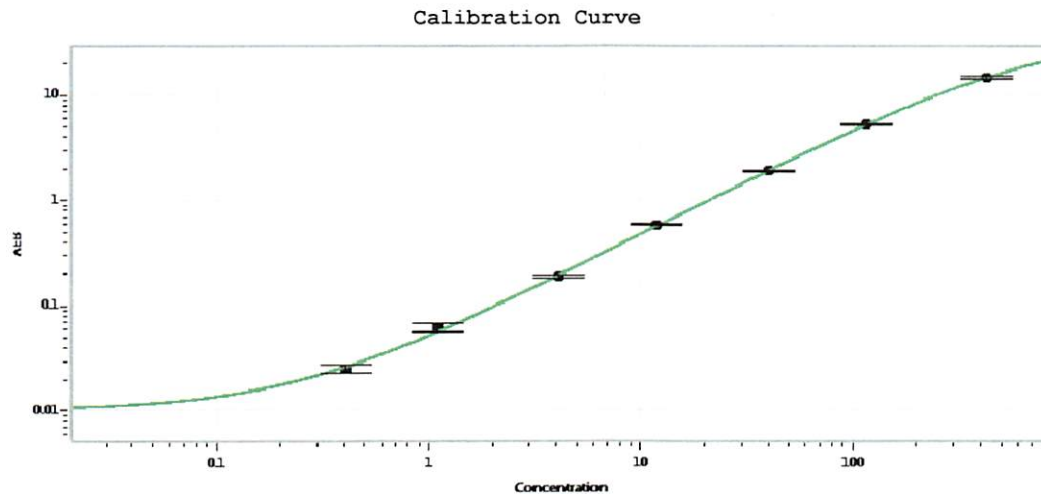
Data below represents results generated on the Simoa HD-1® Analyzer.

Component	Part Number	Lot Number
Bead Reagent	102246	111804
Detector Reagent	102248	111715
SBG Reagent	102250	111406
Sample Diluent	102252	111405
Calibrators	102255	112301
Control 1	102256	112302
Control 2	102257	112302
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	112302	3.69	2.77-4.15
Control 2	112302	203	167-251
Panel 1	008501	13.3	8.59-20.5
Panel 2	008501	10.4	7.39-15.1
Panel 3	008501	58.8	43.8-91.8
Panel 4	009401	191	144-253

* 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 0.411
C 1.10
D 4.12
E 11.8
F 40.6
G 117
H 427

Review/Approval

Junpil Park
Name

QC Supervisor
Title


Signature/Date 19 May 2021