

Product Number:	103186
Lot Number:	503396
Expiration:	2-Dec-2022
Platform(s):	HD-1 / HD-X



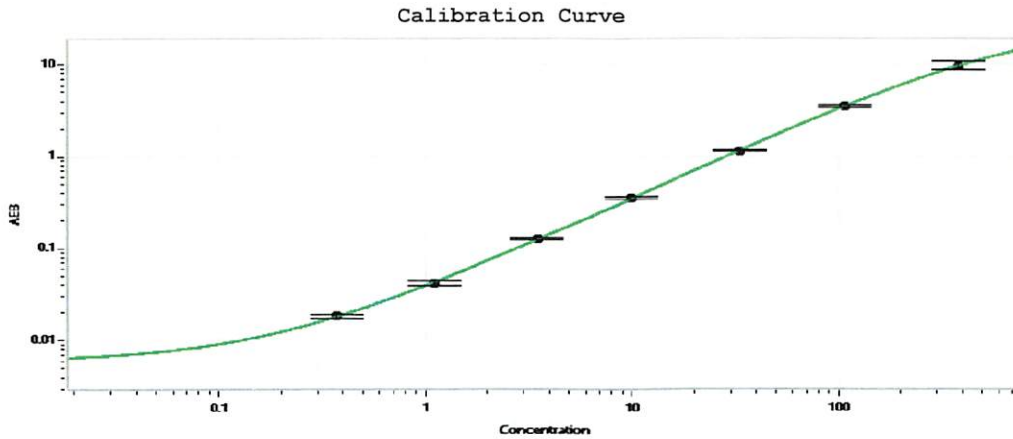
Data below represents results generated on the Simoa HD-1® / HD-X® Analyzer.

Component	Part Number	Lot Number
Bead Reagent	102246	133607
Detector Reagent	102248	134604
SBG Reagent	102250	134709
Sample Diluent	102252	134708
Calibrators	102255	206605
Control 1	102256	206606
Control 2	102257	206606
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	206606	3.32	2.61-3.92
Control 2	206606	197	156-245
Panel 1	008501	12.4	8.59-20.5
Panel 2	008501	10.1	7.39-15.1
Panel 3	008501	61.2	43.8-91.8
Panel 4	009401	183	128-302

* 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 0.371
C 1.11
D 3.50
E 10.0
F 33.5
G 107
H 382

Review/Approval

Richa Shrestha QC Supervisor
Name Title

[Signature] 19APR2022
Signature/Date