

Product Number:	103186
Lot Number:	503706
Expiration:	3-Aug-2023
Platform(s):	HD-1 / HD-X



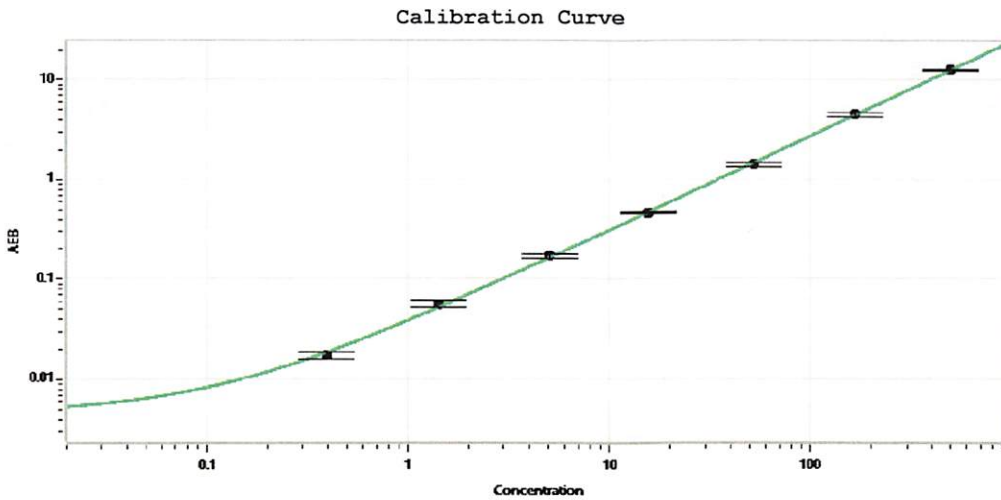
Data below represents results generated on the Simoa HD-X® Analyzer.

Component	Part Number	Lot Number
Bead Reagent	102246	227107
Detector Reagent	102248	227218
SBG Reagent	102250	229301
Sample Diluent	102252	227220
Calibrators	102255	227605
Control 1	102256	227606
Control 2	102257	227606
RGP Reagent	103159	N/A ¹

¹ RGP is not Kit Lot Specific

Release Materials*			
	Lot	Result (pg/mL)	Spec. Range (pg/mL)
Control 1	227606	10.7	8.95-13.4
Control 2	227606	518	386-602
Panel 1	224402	8.69	7.80-12.5
Panel 2	224402	6.80	5.71-9.38
Panel 3	224402	64.8	60.3-91.5
Panel 4	212511	268	166-325

* 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	0.398
C	1.41
D	5.04
E	15.5
F	52.3
G	166
H	493

Review/Approval: Sonia Khilkar Name: Sonia Khilkar Title: QC Manager Signature/Date: [Signature] / 07 NOV 2022