

Product Number:	104073
Lot Number:	503744
Expiration:	21-Nov-2023
Platform(s):	HD-1 / HD-X / SR-X



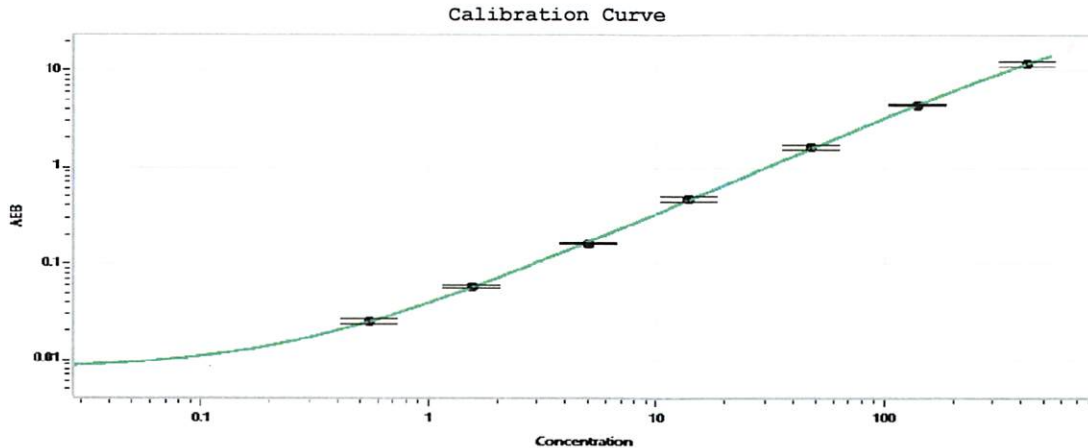
Data below represents results generated on the Simoa HD-X® Analyzer.

Component	Part Number	Lot Number
Bead Reagent	104067	232501
Detector Reagent	104068	232602
SBG Reagent	104069	232604
Sample Diluent	104071	232605
Calibrators	102255	234306
Control 1	102256	234307
Control 2	102257	234307
RGP Reagent	103159	N/A <sup>1</sup>

<sup>1</sup> RGP is not Kit Lot Specific

Release Materials*			
	Lot	Result (pg/mL)	Spec. Range (pg/mL)
Control 1	234307	5.16	4.21 - 6.32
Control 2	234307	220	178 - 267
Panel 1	224402	11.1	8.93 - 13.4
Panel 2	224402	7.99	6.52 - 9.78
Panel 3	224402	79.5	63.9 - 95.8
Panel 4	212511	266	175 - 376

\* 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	0.548
C	1.55
D	5.07
E	13.9
F	48.0
G	140
H	421

Review/Approval Sonia Khilkan QC manager Sonia Khilkan / 31 JAN 2023  
 Name Title Signature/Date