

Product Number:	103714
Lot Number:	503703
Expiration:	22-Aug-2023
Platform(s):	HD-1 / HD-X & SR-X



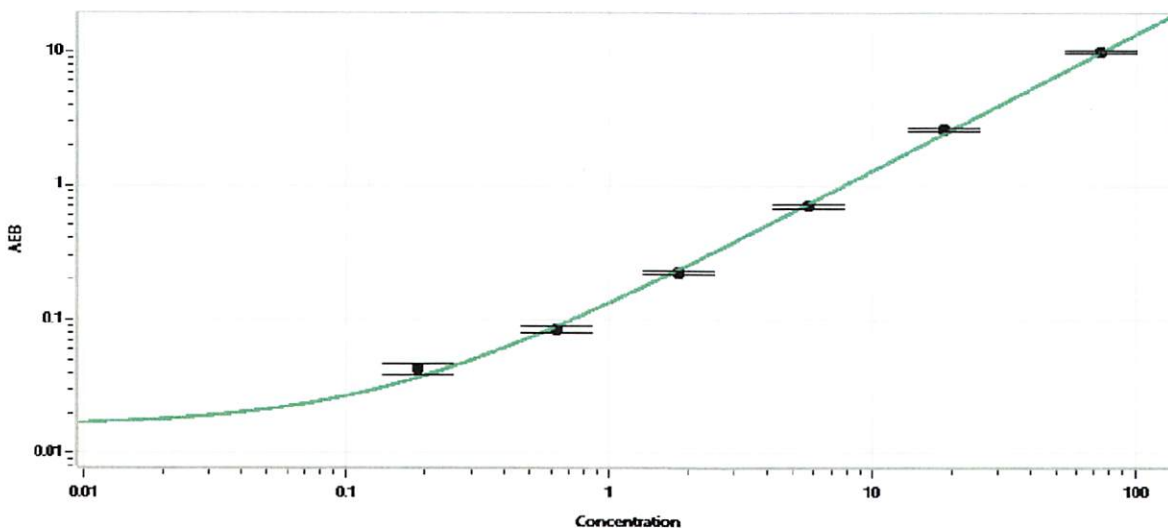
Data below represents results generated on the Simoa® HD-X Analyzer.

Component	Part Number	Lot Number
Bead Reagent	103698	227609
Detector Reagent	103700	227707
SBG Reagent	103702	227708
Sample Diluent	103704	227802
Reference Calibrator	103711	228603
Control 1	103712	228604
Control 2	103713	228604
RGP Reagent	103159	N/A ¹

¹ RGP is not Kit Lot Specific

Release Materials*				* Ranges shown are generated internally for new lot release only. Customer should generate their own control ranges.
	Lot	Result (pg/mL)	Spec. Range (pg/mL)	
Control 1	228604	3.94	3.27-4.90	
Control 2	228604	119	88.5-157	
Panel 1	205508	0.926	0.423-1.21	
Panel 2	205508	2.88	1.76-3.31	
Panel 3	211609	23.0	16.6-27.7	

Calibration Curve



Calibrator Levels (pg/mL)	
A	0.000
B	0.188
C	0.629
D	1.82
E	5.73
F	18.6
G	73.4

Review/Approval

Sonia Khillan
Name

QC Manager
Title

Sonia Khillan / 25 OCT 2022
Signature/Date