

Product Number:	103714
Lot Number:	503403
Expiration:	21-Mar-2023
Platform(s):	HD-1 / HD-X & SR-X



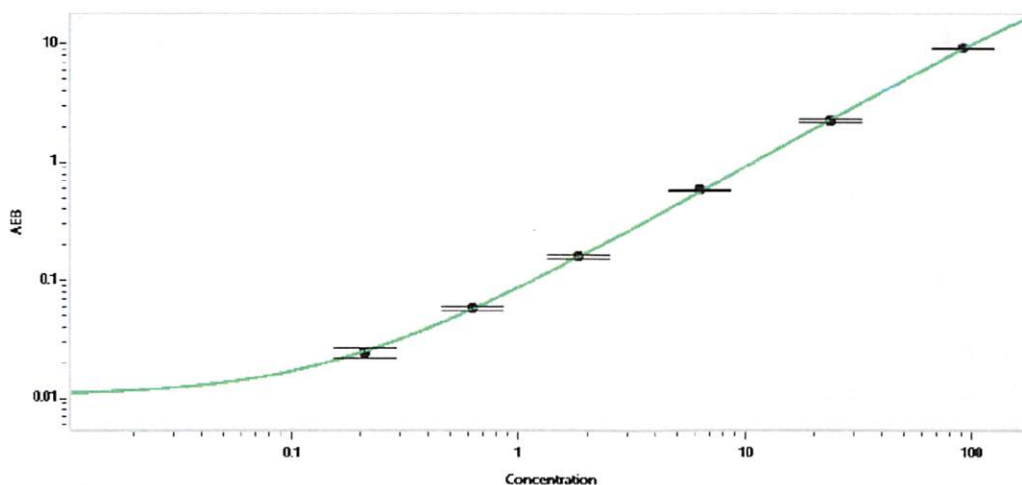
Data below represents results generated on the Simoa® HD-1 / HD-X & SR-X Analyzer.

Component	Part Number	Lot Number
Bead Reagent	103698	208907
Detector Reagent	103700	208007
SBG Reagent	103702	208006
Sample Diluent	103704	208102
Reference Calibrator	103711	208003
Control 1	103712	208004
Control 2	103713	208004
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)	Mean Range (pg/mL)	
Control 1	208004	4.12	3.21-4.82	
Control 2	208004	109.9	91.4-137	
Panel 1	205508	0.848	0.383-1.21	
Panel 2	205508	2.62	1.70-3.27	
Panel 3	211609	21.47	16.0-27.9	

Calibration Curve



Calibrator Levels (pg/mL)

A	0.000
B	0.210
C	0.626
D	1.83
E	6.31
F	23.8
G	91.2

Review/Approval

Matt Smart
Name

QC Director
Title

M. Smart 22 Jun 2022
Signature/Date