

Product Number:	103714
Lot Number:	503356
Expiration:	15-Feb-2023
Platform(s):	HD-1 / HD-X & SR-X

Component	Part Number	Lot Number
Bead Reagent	103698	206104
Detector Reagent	103700	206105
SBG Reagent	103702	206106
Sample Diluent	103704	206107
Reference Calibrator	103711	207204
Control 1	103712	207205
Control 2	103713	207205
RGP Reagent	103159	N/A ¹

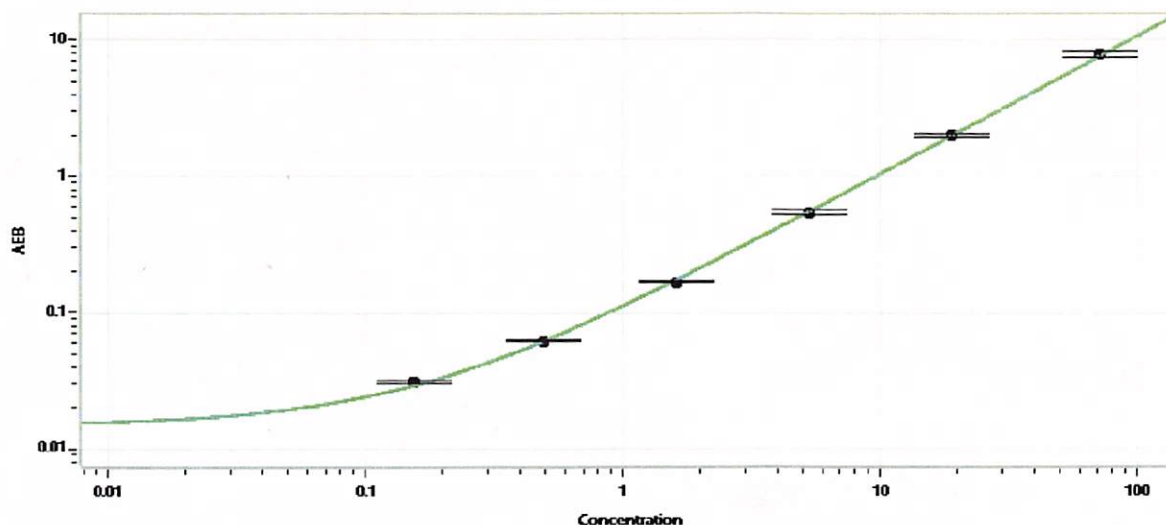
¹ RGP is not Kit Lot Specific



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

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	207205	2.99	2.32-3.47	
Control 2	207205	100	79.3-119	
Panel 1	013903	0.770	0.258-1.11	
Panel 2	013903	1.78	0.841-2.44	
Panel 3	013903	2.54	1.79-4.31	

Calibration Curve



Calibrator Levels (pg/mL)

A	0.000
B	0.155
C	0.488
D	1.61
E	5.32
F	19.1
G	72.3

Review/Approval

Shivani Goel
Name

QC Manager
Title

Shivani Goel 23 Mar 2022
Signature/Date