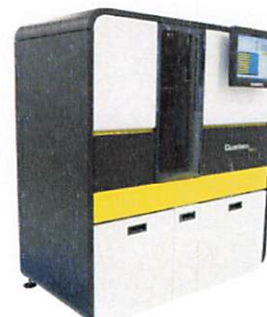


Product Number:	103714
Lot Number:	503732
Expiration:	3-Oct-2023
Platform(s):	HD-1 / HD-X & SR-X

Component	Part Number	Lot Number
Bead Reagent	103698	227609
Detector Reagent	103700	231308
SBG Reagent	103702	231309
Sample Diluent	103704	231310
Reference Calibrator	103711	228603
Control 1	103712	228604
Control 2	103713	228604
RGP Reagent	103159	N/A ¹

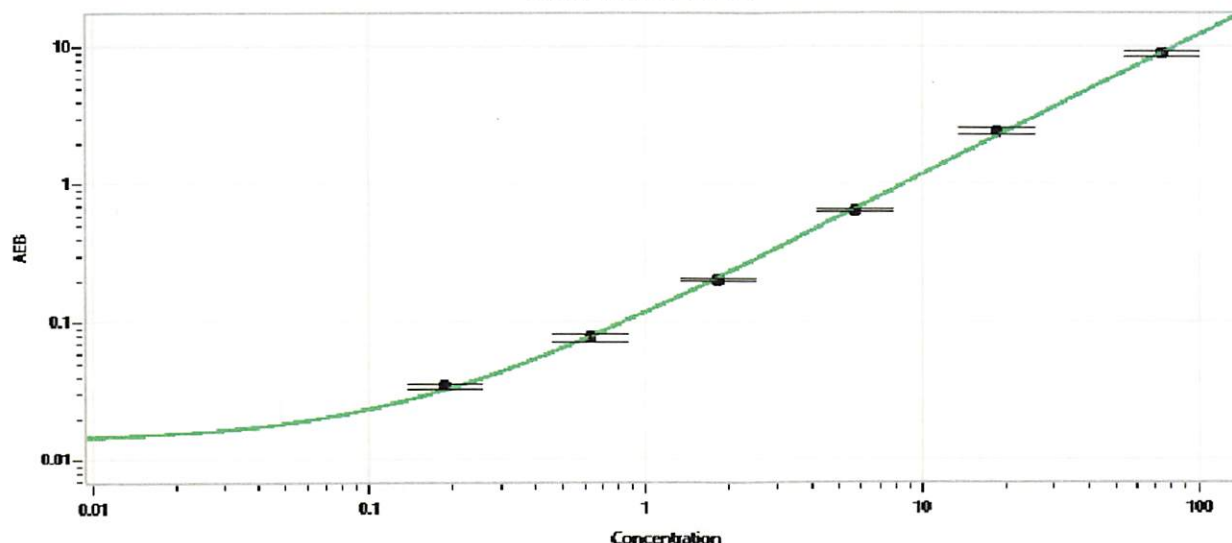
¹ RGP is not Kit Lot Specific



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

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	4.01	3.25 - 4.88	
Control 2	228604	112	87.5 - 154	
Panel 1	205508	0.857	0.423 - 1.21	
Panel 2	205508	2.84	1.76 - 3.31	
Panel 3	211609	21.5	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

Title

Sonia Khillan / 18 Nov 2022
Signature/Date