

Product Number:	101580
Lot Number:	503457
Expiration:	29-Apr-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	101582	211908
Detector Reagent	101581	232708
SBG Reagent	101583	212002
Sample Diluent	103735	230415
Calibrator Diluent	101585	212006
Calibrator Concentrate	100655	212004
RGP Reagent	103159	N/A ¹

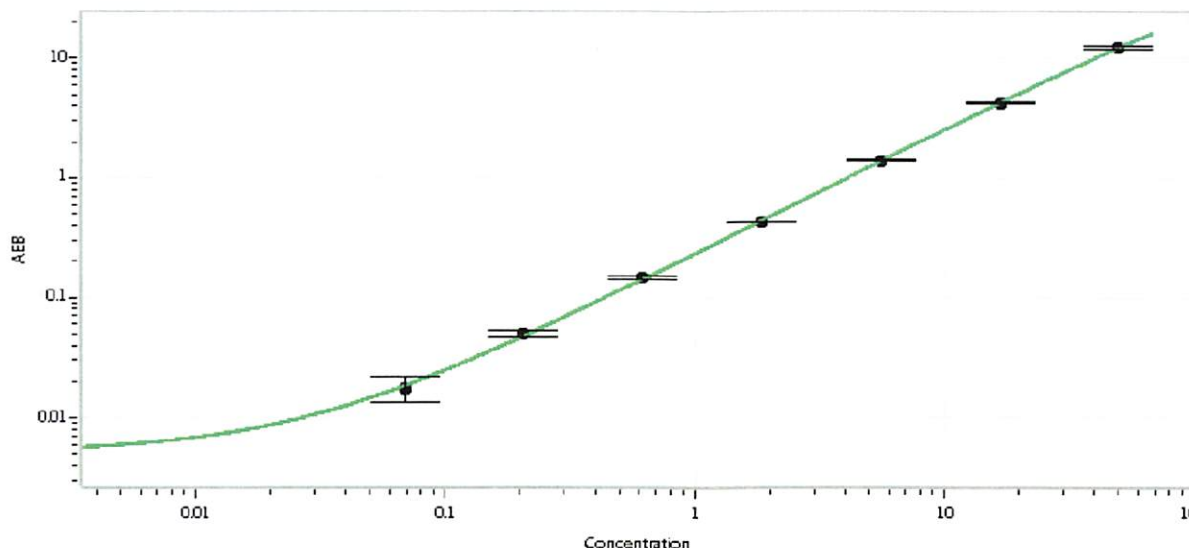
¹ RGP is not Kit Lot Specific

Release Materials*			
	Lot	Result (pg/mL)	Spec. Range (pg/mL)
Control 1	206502	2.85	2.23-3.36
Control 2	206502	45.5	35.8-56.5
Panel 2	115317	2.19	1.51-3.73
Panel 3	116301	27.4	15.5-39.3

* Ranges shown are generated internally for new lot release only. Customer should generate their own control ranges.

Calibrator Concentrate Value Assignment (pg/mL)	1065
---	------

Calibration Curve



Calibrator Levels (pg/mL)	
A	0.000
B	0.069
C	0.206
D	0.617
E	1.85
F	5.56
G	16.7
H	50.0

Review/Approval: Sonia Khillan AC Manager Sonia Khillan / 06 DEC 2022

Name: Sonia Khillan Title: AC Manager Signature/Date: Sonia Khillan / 06 DEC 2022