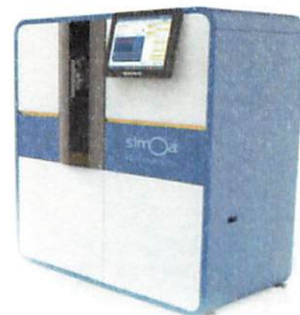


Product Number:	102732
Lot Number:	503734
Expiration:	26-Oct-2023
Platform(s):	HD-1/HD-X & SR-X



Data below represents results generated on the Simoa® Analyzer.

Component	Part Number	Lot Number
Bead Reagent	102722	229908
Detector Reagent	102724	230111
SBG Reagent	102726	230110
Sample Diluent	102728	229912
Calibrator Diluent	102731	230619
Calibrator Concentrate	100658	204302
RGP Reagent	103159	N/A ¹

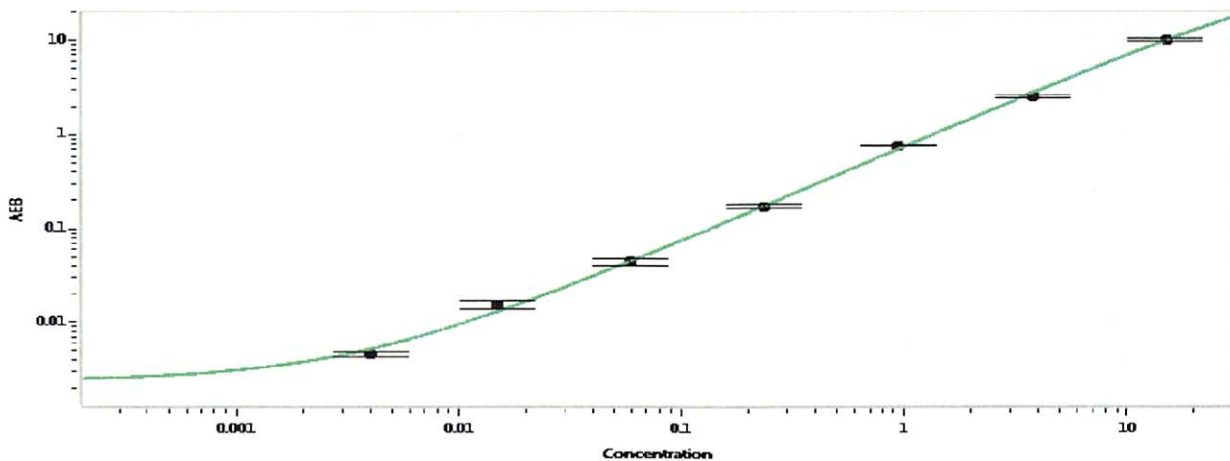
¹ RGP is not Kit Lot Specific

Release Materials*			
	Lot	Result (pg/mL)	Spec. Range (pg/mL)
Control 1	114108	0.107	0.067 - 0.165
Control 2	114108	2.93	2.01 - 3.94
Panel 2	112714	0.035	0.011 - 0.054
Panel 3	112714	1.53	0.832 - 2.29
Panel 4	112714	5.15	2.50 - 8.63

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

Calibrator Concentrate Value Assignment (pg/mL)	375
---	-----

Calibration Curve



Calibrator Levels (pg/mL)	
A	0.000
B	0.004
C	0.015
D	0.059
E	0.234
F	0.938
G	3.75
H	15.0

Review/Approval

Sonia Khillan
Name

QC
Title

Sonia Khillan 08 FEB 2023
Signature/Date