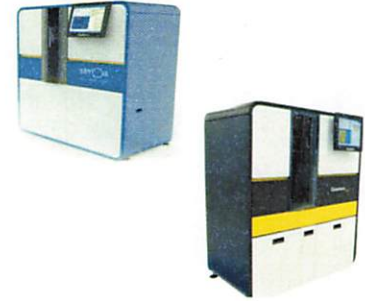


Product Number:	101635
Lot Number:	503379
Expiration:	27-Feb-2023
Platform(s):	HD-1/ HD-X



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

Component	Part Number	Lot Number
Bead Reagent	101637	206804
Detector Reagent	101636	206803
SBG Reagent	101638	206904
Sample Diluent	101639	206905
Calibrator Diluent	101640	206906
Calibrator Concentrate	100661	206907
RGP Reagent	103159	N/A ¹

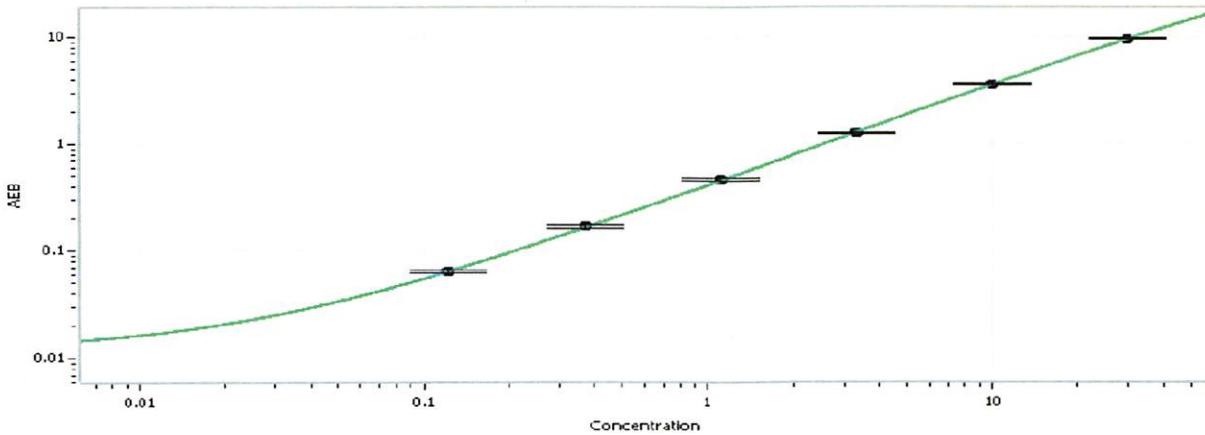
¹ RGP is not Kit Lot Specific

Release Materials*			
	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	129010	0.493	0.389-0.605
Control 2	129010	50.2	38.9-59.7
Panel 1	035301	0.315	0.206-0.492
Panel 2	035301	3.59	2.65-4.99
Panel 3	035301	44.0	28.3-55.8

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

Calibrator Concentrate Value Assignment (pg/mL)	2568
---	------

Calibration Curve



Calibrator Levels (pg/mL)	
A	0.000
B	0.123
C	0.370
D	1.11
E	3.33
F	10.0
G	30.0

Review/Approval

Shivani Goel
Name

QC Manager
Title

Shivani Goel 12 Apr 2020
Signature/Date