

Product Number:	101635
Lot Number:	503323
Expiration:	24-Jan-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	202406
Detector Reagent	101636	203305
SBG Reagent	101638	202804
Sample Diluent	101639	202805
Calibrator Diluent	101640	202806
Calibrator Concentrate	100661	127604
RGP Reagent	103159	N/A ¹

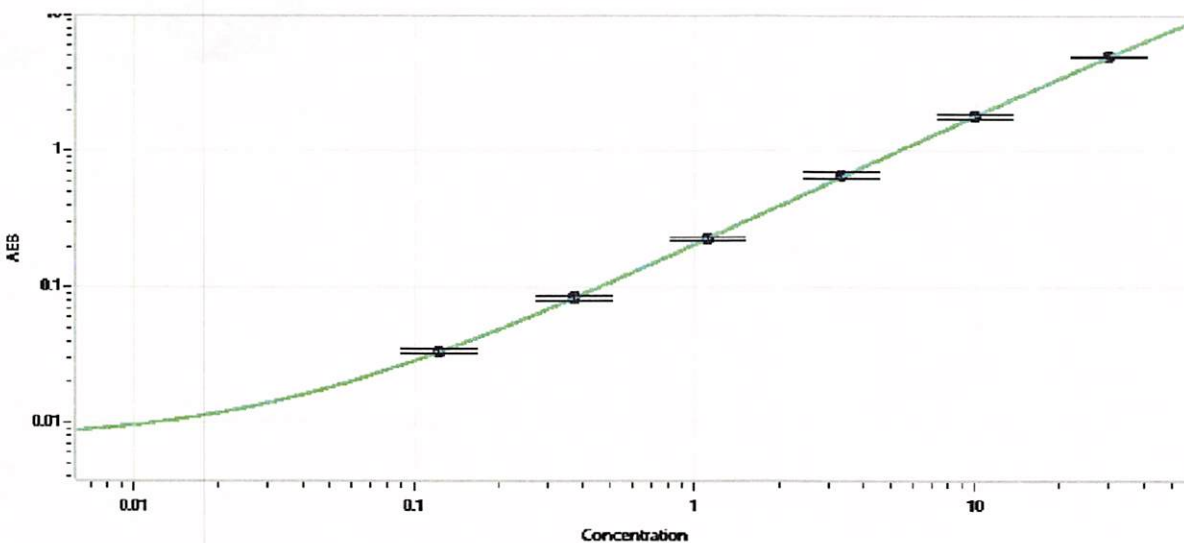
¹ RGP is not Kit Lot Specific

Release Materials*			
	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	129010	0.469	0.379-0.617
Control 2	129010	46.1	37.6-60.7
Panel 1	035301	0.343	0.209-0.503
Panel 2	035301	3.04	2.62-5.10
Panel 3	035301	31.1	26.9-56.5

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

Calibrator Concentrate Value Assignment (pg/mL)	1974
---	------

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 28 Mar 2022
Signature/Date