

Product Number:	101132
Lot Number:	502594
Expiration:	2-Jun-2021
Platform(s):	HD-1/HD-X



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

Component	Part Number	Lot Number
Bead Reagent	101119	029509
Detector Reagent	101121	029706
SBG Reagent	101131	029707
Sample Diluent	103319	015435
Calibrator Diluent	101125	029712
Calibrator Concentrate	101123	016712
RGP Reagent	103159	N/A ¹

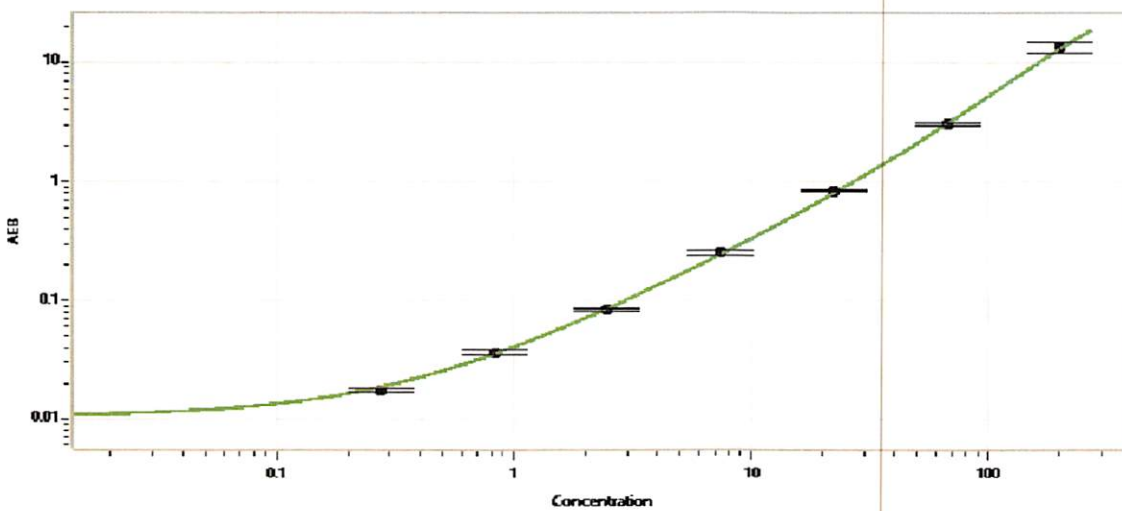
¹ RGP is not Kit Lot Specific

Release Materials*			
	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	920702	2.51	1.66-3.21
Control 2	920702	258	122-349
Panel 1	830204	89.1	78.7-118
Panel 2	830205	77.2	68.7-103
Panel 3	830206	213	193-289
Panel 4	830207	281	227-393

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

Calibrator Concentrate Value Assignment (pg/mL)	8099
---	------

Calibration Curve



Calibrator Levels (pg/mL)	
A	0.000
B	0.274
C	0.823
D	2.47
E	7.41
F	22.2
G	66.7
H	200

Review/Approval

Shivani Goel
Name

QC Manager
Title

Shivani Goel 19 Nov 2020
Signature/Date