

Product Number:	102329
Lot Number:	502273
Expiration:	29-Jan-2021
Platform(s):	HD-1/HD-X



Data below represents results generated on the Simoa™ HD-1 Analyzer.

Component	Part Number	Lot Number
Bead Reagent	102322	005706
Detector Reagent	102324	005513
SBG Reagent	102325	005602
Sample Diluent	102326	005218
Calibrator Diluent	102327	005217
Calibrator Concentrate	100849	006603
RGP Reagent	103159	N/A ¹

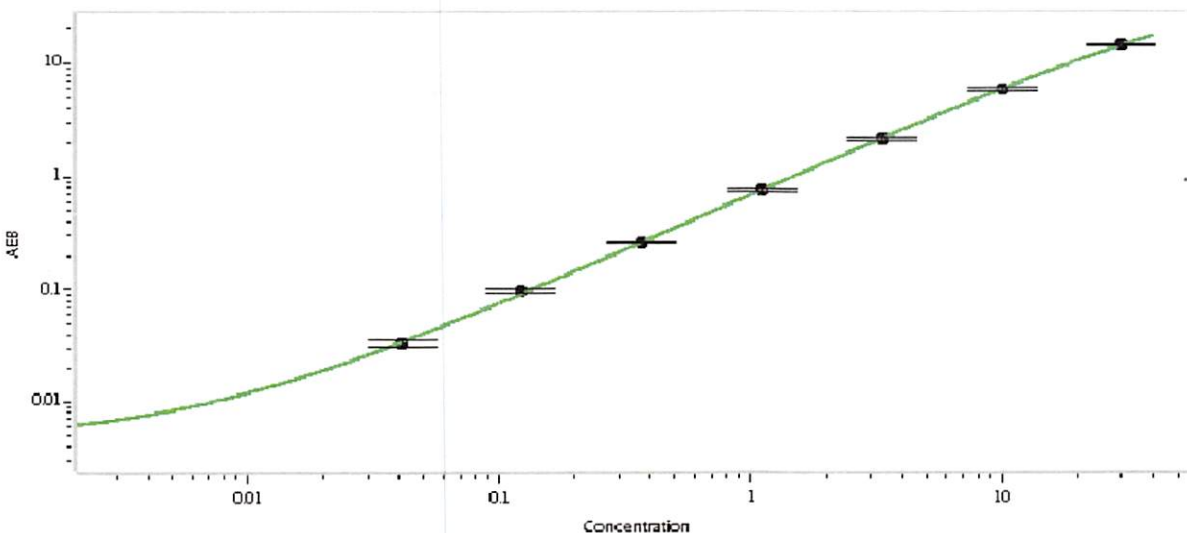
¹ RGP is not Kit Lot Specific

Release Materials*			
	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	MF-092716-01	0.208	0.139-0.285
Control 2	MF-092716-01	27.0	19.4-37.3
Panel 1	MF-092116-03	0.096	0.065-0.135
Panel 2	MF-092116-04	0.084	0.065-0.135
Panel 3	MF-092116-05	16.8	10.5-25.3

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

Calibrator Concentrate Value Assignment (pg/mL)	4586
---	------

Calibration Curve



Calibrator Levels (pg/mL)

A	0.000
B	0.041
C	0.123
D	0.370
E	1.11
F	3.33
G	10.0
H	30.0

Review/Approval

Patrizia Stadler Assoc. Director, QC

P. Stadler

12/17/2020

Name

Title

Signature/Date