

Product Number:	101132
Lot Number:	502429
Expiration:	29-Apr-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	101119	016704
Detector Reagent	101121	016707
SBG Reagent	101131	016708
Sample Diluent	103319	015416
Calibrator Diluent	101125	015113
Calibrator Concentrate	101123	016712
RGP Reagent	103159	N/A ¹

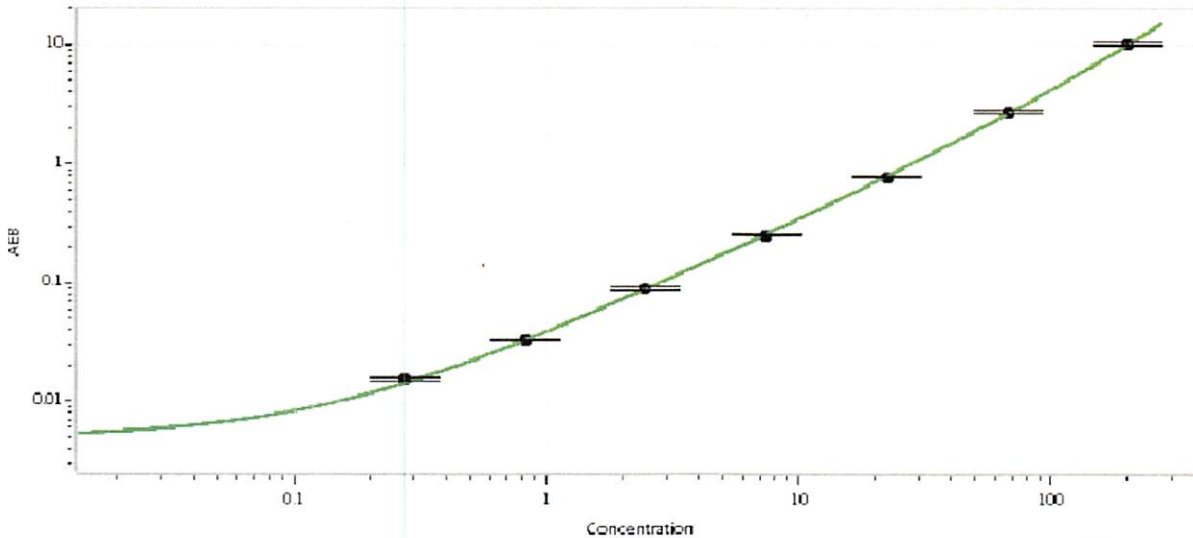
¹ RGP is not Kit Lot Specific

Release Materials*			
	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	920702	2.61	1.62-3.21
Control 2	920702	288	108-351
Panel 1	830204	104	79.6-119
Panel 2	830205	89.7	69.6-104
Panel 3	830206	269	195-293
Panel 4	830207	376	233-393

* 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	Patrizia Stadler Assoc. Director, QC	<i>p. stadler</i>	08 Jul 2020
	Name	Title	Signature/Date