

Product Number:	101154
Lot Number:	502398
Expiration:	21-Feb-2021
Platform(s):	HD-1/HDX



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

Component	Part Number	Lot Number
Bead Reagent	101141	014707
Detector Reagent	101143	014702
SBG Reagent	101153	014701
Sample Diluent	103320	005233
Calibrator Diluent	101147	012512
Calibrator Concentrate	101145	004202
RGP Reagent	103159	N/A ¹

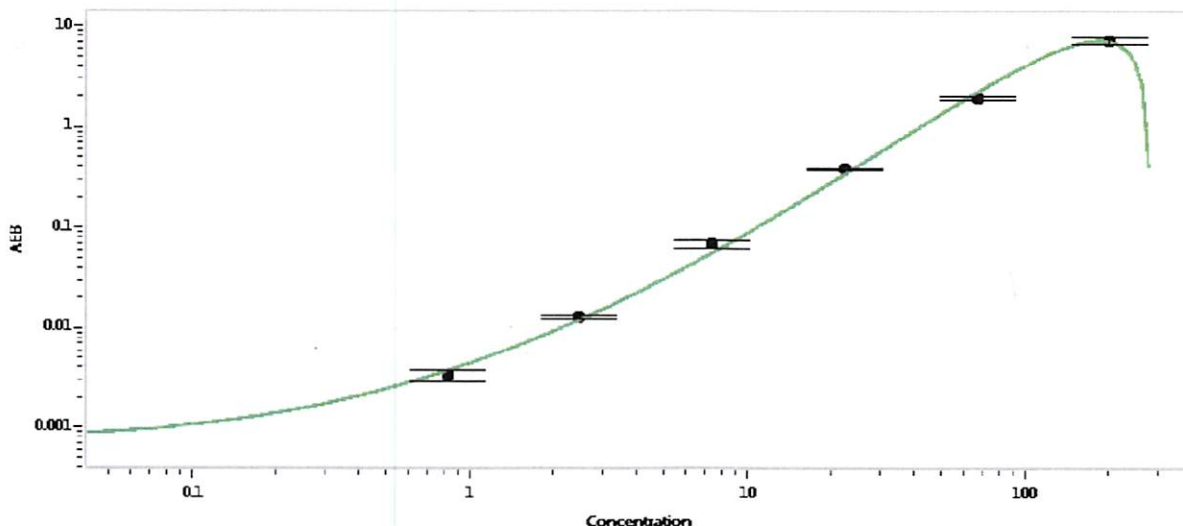
¹ RGP is not Kit Lot Specific

Release Materials*			
	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	918402	49.0	31.7-50.4
Control 2	918402	398	251-461
Panel 1	909901	104	66.3-112
Panel 2	909902	93.2	58.0-102
Panel 3	909903	317	161-367

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

Calibrator Concentrate Value Assignment (pg/mL)	4130
---	------

Calibration Curve



Calibrator Levels (pg/mL)

A	0.000
B	0.823
C	2.47
D	7.41
E	22.2
F	66.7
G	200

Review/Approval: Patrizia Stadler Assoc. Director, QC *p.stadler* 03 Jun 2020
Name Title Signature/Date