

Product Number:	102732
Lot Number:	502282
Expiration:	12-Feb-2021
Platform(s):	HD-1/HD-X & SR-X



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

Component	Part Number	Lot Number
Bead Reagent	102722	007201
Detector Reagent	102724	006403
SBG Reagent	102726	006404
Sample Diluent	102728	006412
Calibrator Diluent	102731	006411
Calibrator Concentrate	100658	007202
RGP Reagent	103159	N/A ¹

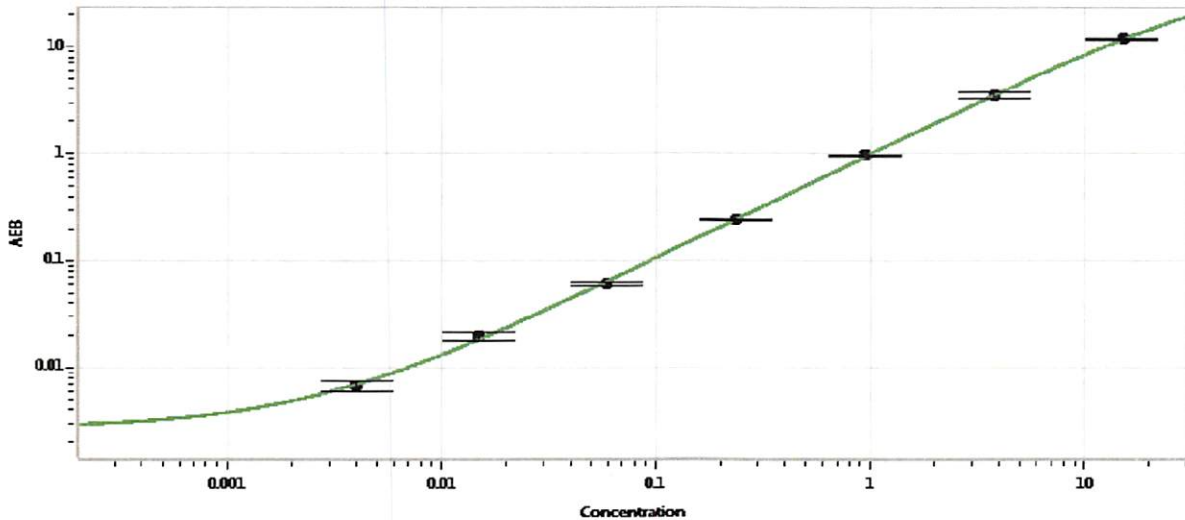
¹ RGP is not Kit Lot Specific

Release Materials*			
	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	909502	0.117	0.096-0.157
Control 2	909502	3.33	2.92-4.53
Panel 1	911401	0.033	0.013-0.051
Panel 3	911403	1.40	1.22-2.02
Panel 4	911404	4.53	3.94-7.52

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

Calibrator Concentrate Value Assignment (pg/mL)	229
---	-----

Calibration Curve



Calibrator Levels (pg/mL)

A	0.000
B	0.004
C	0.015
D	0.059
E	0.234
F	0.938
G	3.75
H	15.0

Review/Approval	Patrizia Stadler	Assoc. Director, QC	<i>P. Stadler</i>	25 Mar 2020
	Name	Title	Signature/Date	