

Product Number:	102732
Lot Number:	503022
Expiration:	22-May-2022
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	117910
Detector Reagent	102724	118018
SBG Reagent	102726	117916
Sample Diluent	102728	114311
Calibrator Diluent	102731	114204
Calibrator Concentrate	100658	118008
RGP Reagent	103159	N/A ¹

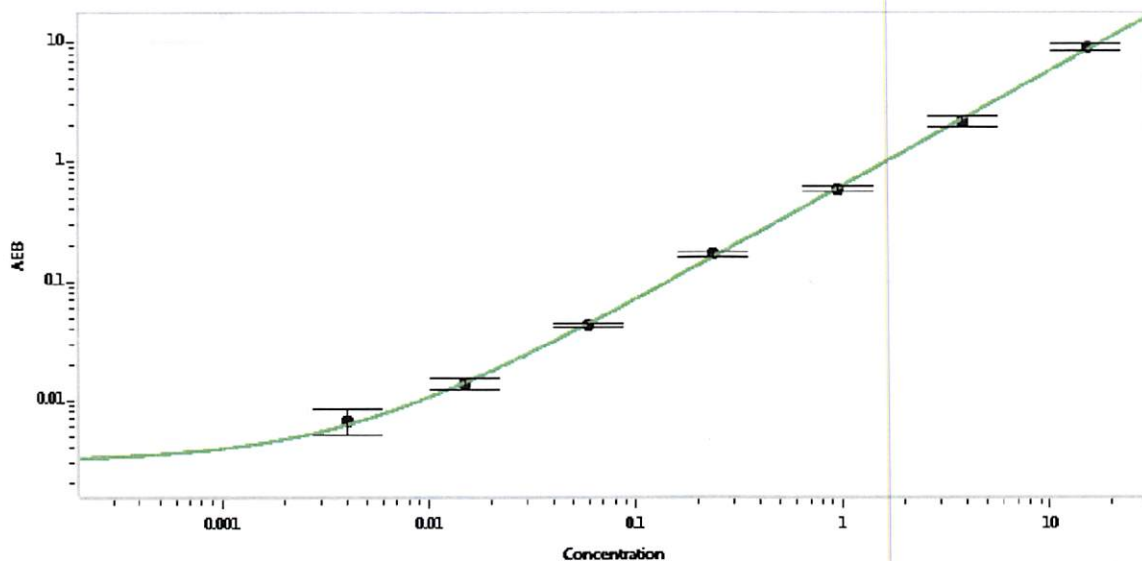
¹ RGP is not Kit Lot Specific

Release Materials*			
	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	114108	0.076	0.067-0.159
Control 2	114108	2.57	1.91-3.81
Panel 2	112714	0.028	0.010-0.049
Panel 3	112714	1.27	1.02-1.87
Panel 4	112714	4.59	2.79-7.70

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

Calibrator Concentrate Value Assignment (pg/mL)	352
-------------------------------------------------	-----

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

Shivani Goel

QC Manager

Shivani Goel 14 Jul 2021

Name

Title

Signature/Date