

Product Number:	102329
Lot Number:	502488
Expiration:	2-Jun-2021
Platform(s):	HD-1/HD-X

Component	Part Number	Lot Number
Bead Reagent	102322	020503
Detector Reagent	102324	020501
SBG Reagent	102325	022001
Sample Diluent	102326	015420
Calibrator Diluent	102327	020420
Calibrator Concentrate	100849	006603
RGP Reagent	103159	N/A ¹

¹ RGP is not Kit Lot Specific

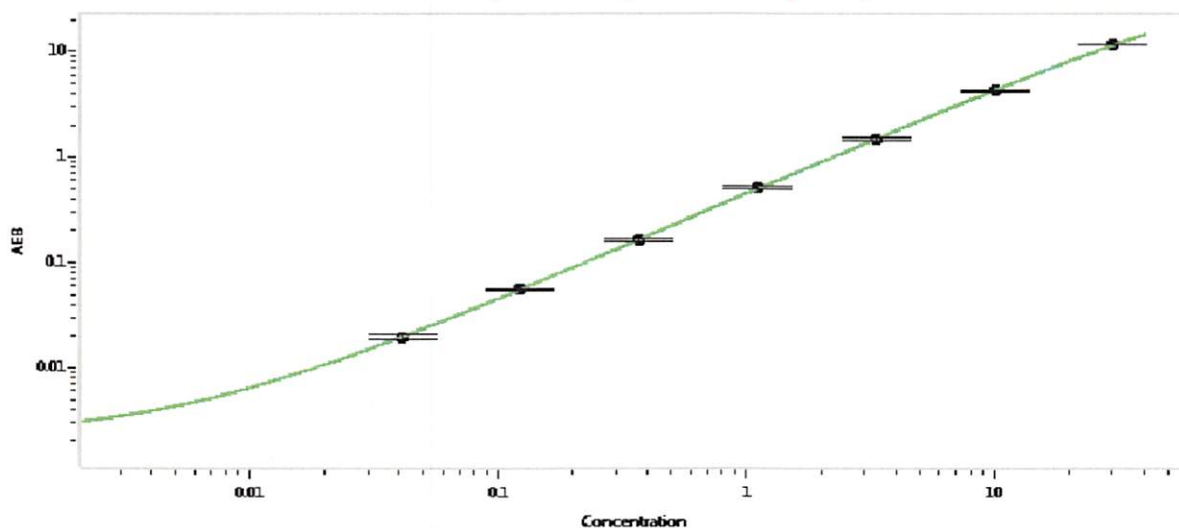


Data below represents results generated on the Simoa™ HD-1 Analyzer.

Release Materials*				* Ranges shown are generated internally for new lot release only. Customer should generate their own control ranges.
	Lot	Result (pg/mL)	Mean Range (pg/mL)	
Control 1	023702	0.193	0.127-0.211	
Control 2	023702	23.9	18.4-30.0	
Panel 1	027607	0.098	0.032-0.136	
Panel 2	027607	0.136	0.064-0.171	
Panel 3	027607	13.4	8.06-15.5	

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

Junpil Park
Name

QC Supervisor
Title

[Signature] 15 Jan 2021
Signature/Date