

Product Number:	102732
Lot Number:	502565
Expiration:	2-Oct-2021
Platform(s):	HD-1/HD-X & SR-X

Component	Part Number	Lot Number
Bead Reagent	102722	027907
Detector Reagent	102724	031126
SBG Reagent	102726	031111
Sample Diluent	102728	031202
Calibrator Diluent	102731	027613
Calibrator Concentrate	100658	013605
RGP Reagent	103159	N/A <sup>1</sup>

<sup>1</sup> RGP is not Kit Lot Specific



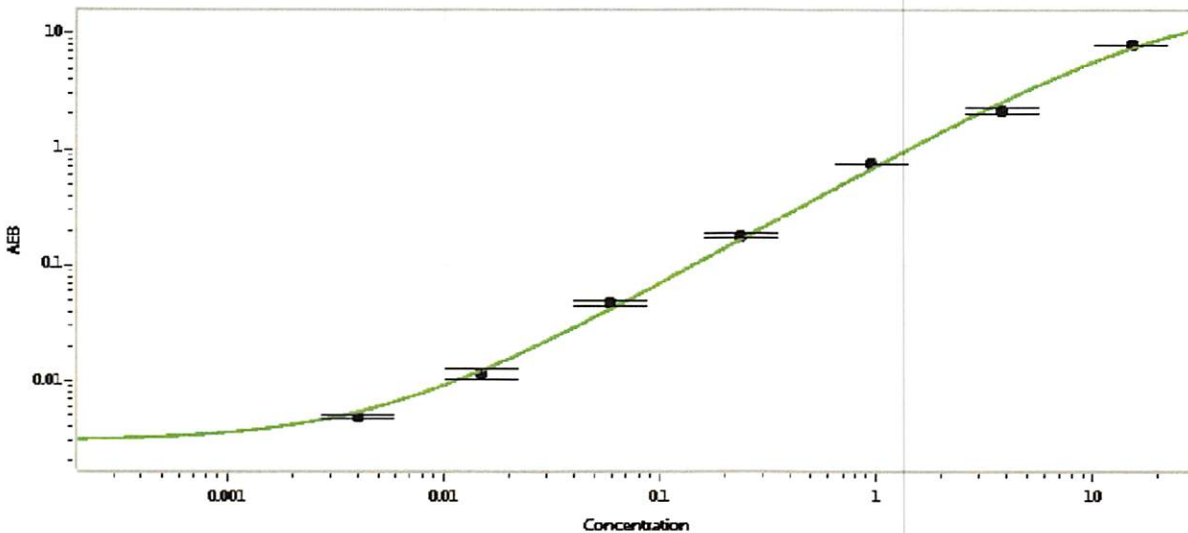
Data below represents results generated on the Simoa™ HD-1 Analyzer.

Release Materials*			
	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	909502	0.158	0.093-0.167
Control 2	909502	2.98	2.83-4.48
Panel 1	911401	0.041	0.017-0.050
Panel 3	911403	1.65	1.18-2.02
Panel 4	911404	4.20	3.80-7.25

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

Calibrator Concentrate Value Assignment (pg/mL)	497
---	-----

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  
Name

QC Manager  
Title

Shivani Goel 24 Nov 2020  
Signature/Date