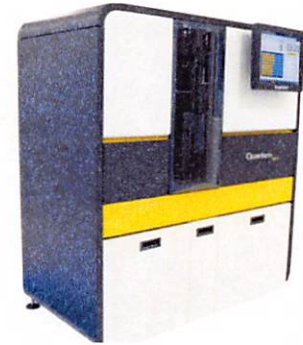


Product Number:	101552
Lot Number:	503264
Expiration:	26-Sep-2022
Platform(s):	HD-1/HD-x & SR-X

Component	Part Number	Lot Number
Bead Reagent	101554	126905
Detector Reagent	101553	134411
SBG Reagent	101555	126906
Sample Diluent	101556	126907
Calibrators	101559	127104
Control 1	101560	127105
Control 2	101561	127105
SBG Diluent (SR-X)†	103158	N/A <sup>1</sup>
RGP Reagent	103159	N/A <sup>1</sup>

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

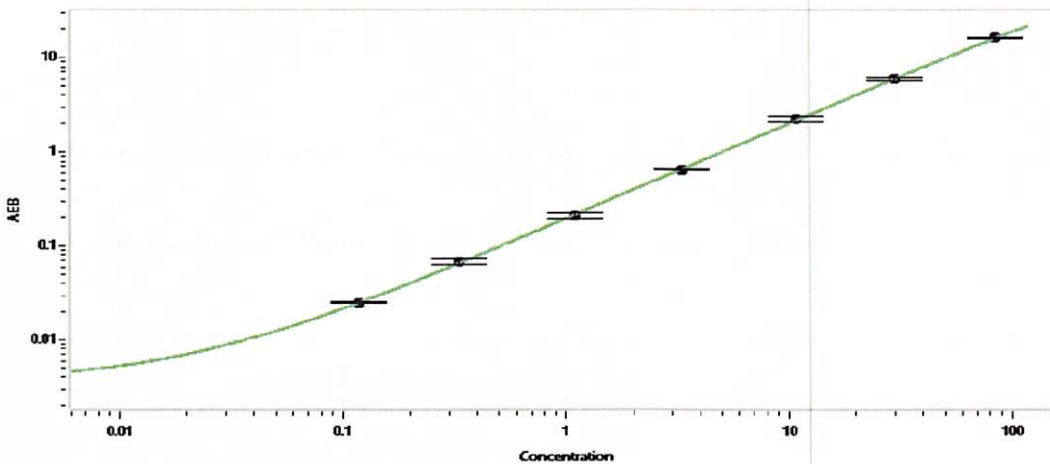


Data below represents results generated on the Simoa™ HD-X Analyzer.

Release Materials*			
	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	127105	1.20	0.784-1.52
Control 2	127105	63.7	47.8-71.7
Panel 1	117205	4.85	3.48 - 5.22
Panel 2	117205	3.27	2.00 - 4.90
Panel 3	117205	23.3	17.9 -26.8
Panel 4	119612	35.5	22.6 - 41.6

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

Calibration Curve



Calibrator Levels (pg/mL)	
A	0.000
B	0.118
C	0.331
D	1.10
E	3.27
F	10.6
G	29.4
H	83.7

Review/Approval

Wenkai Tsai  
Name

QC Supervisor  
Title

*[Signature]*  
Signature/Date

13 Dec 2021