

Product Number:	103789
Lot Number:	502707
Expiration:	25-Jul-2021
Platform(s):	HD-X



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

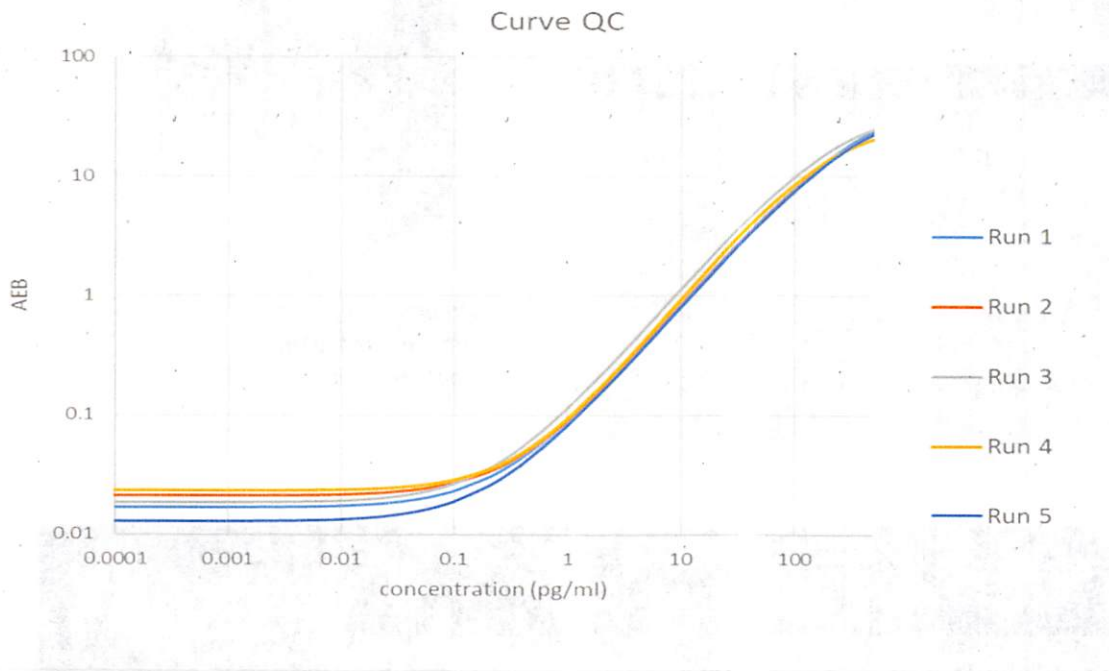
Component	Part Number	Lot Number
Bead Reagent	103776	102513
Detector Reagent	103778	102519
SBG Reagent	103780	102608
Sample Diluent	103782	102525
Calibrators	103786	102603
Control 1	103787	102604
Control 2	103788	102604
Control 3	103809	102604
RGP Reagent	103844	N/A <sup>1</sup>

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

Release Materials*				* Ranges shown are generated internally for new lot release only. Customer should generate their own control ranges.
Sample	Lot	Result (pg/mL)	Mean Range (pg/mL)	
Control 1	102604	5.09	4.04-6.29	
Control 2	102604	108	86.6-136	
Control 3	102604	0.137	< 3.00	
Sample	Lot	Bridging Result	Kit Release Result	Pass/ Fail
Panel 1	104301	Negative	Negative	PASS
Panel 2	104302	Negative	Negative	PASS
Panel 3	104303	Negative	Negative	PASS
Panel 4	104304	Negative	Negative	PASS
Panel 5	104305	Negative	Negative	PASS
Panel 6	104306	Positive	Positive	PASS
Panel 7	104307	Positive	Positive	PASS
Panel 8	104308	Positive	Positive	PASS
Panel 9	104309	Positive	Positive	PASS
Panel 10	104310	Positive	Positive	PASS
Panel 11	104311	Positive	Positive	PASS
Panel 12	104312	Positive	Positive	PASS
Panel 13	104313	Positive	Positive	PASS
Panel 14	104314	Positive	Positive	PASS

<b>Review/Approval</b>	Junpil Park	QC Supervisor	22 Feb 2021
	Name	Title	Signature/Date

Calibration Curve



Calibrator Levels  
(ng/mL)

<b>A</b>	0.000
<b>B</b>	0.204
<b>C</b>	0.517
<b>D</b>	1.61
<b>E</b>	4.74
<b>F</b>	13.4
<b>G</b>	40.2
<b>H</b>	147