

<b>Product Number:</b>	103789
<b>Lot Number:</b>	503131
<b>Expiration:</b>	13-Mar-2022
<b>Platform(s):</b>	HD-X



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

Component	Part Number	Lot Number
Bead Reagent	103776	125601
Detector Reagent	103778	125602
SBG Reagent	103780	125606
Sample Diluent	103782	125607
Calibrators	103786	126204
Control 1	103787	126205
Control 2	103788	126205
Control 3	103809	126206
RGP Reagent	103844	N/A <sup>1</sup>

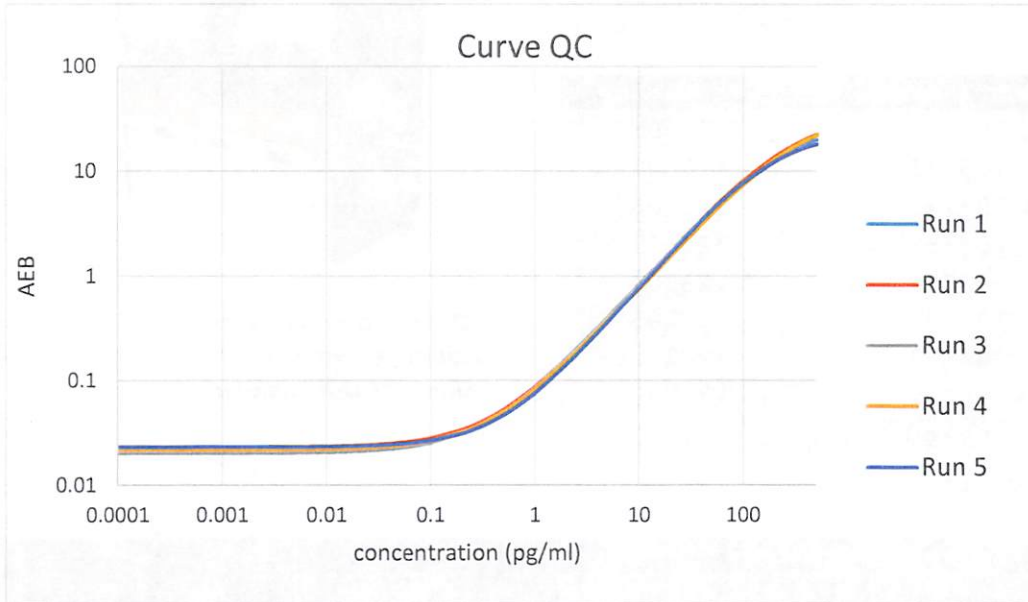
<sup>1</sup> RGP is not Kit Lot Specific

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	126205	7.75	6.21-9.31	
Control 2	126205	147	107-172	
Control 3	126206	0.017	< 3.00	
Sample	Lot	Bridging Result	Kit Release Result	Pass/ Fail
Panel 1	202101	Negative	Negative	PASS
Panel 2	202102	Negative	Negative	PASS
Panel 3	202103	Negative	Negative	PASS
Panel 4	202104	Negative	Negative	PASS
Panel 5	202105	Negative	Negative	PASS
Panel 6	202106	Positive	Positive	PASS
Panel 7	202107	Positive	Positive	PASS
Panel 8	202108	Positive	Positive	PASS
Panel 9	202109	Positive	Positive	PASS
Panel 10	202110	Positive	Positive	PASS
Panel 11	202111	Positive	Positive	PASS
Panel 12	202112	Positive	Positive	PASS
Panel 13	202113	Positive	Positive	PASS
Panel 14	202114	Positive	Positive	PASS

<b>Review/Approval</b>	Junpil Park	QC Supervisor		08 Oct 2021
	<b>Name</b>	<b>Title</b>	<b>Signature/Date</b>	



**Calibration Curve**



**Calibrator Levels  
(pg/mL)**

<b>A</b>	0.000
<b>B</b>	0.301
<b>C</b>	0.762
<b>D</b>	2.38
<b>E</b>	6.98
<b>F</b>	17.4
<b>G</b>	52.0
<b>H</b>	191