

Product Number:	103789
Lot Number:	502623
Expiration:	16-May-2021
Platform(s):	HD-X



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

Component	Part Number	Lot Number
Bead Reagent	103776	036314
Detector Reagent	103778	032108
SBG Reagent	103780	032105
Sample Diluent	103782	032102
Calibrators	103786	035306
Control 1	103787	032204
Control 2	103788	032204
Control 3	103809	032304
RGP Reagent	103844	N/A ¹

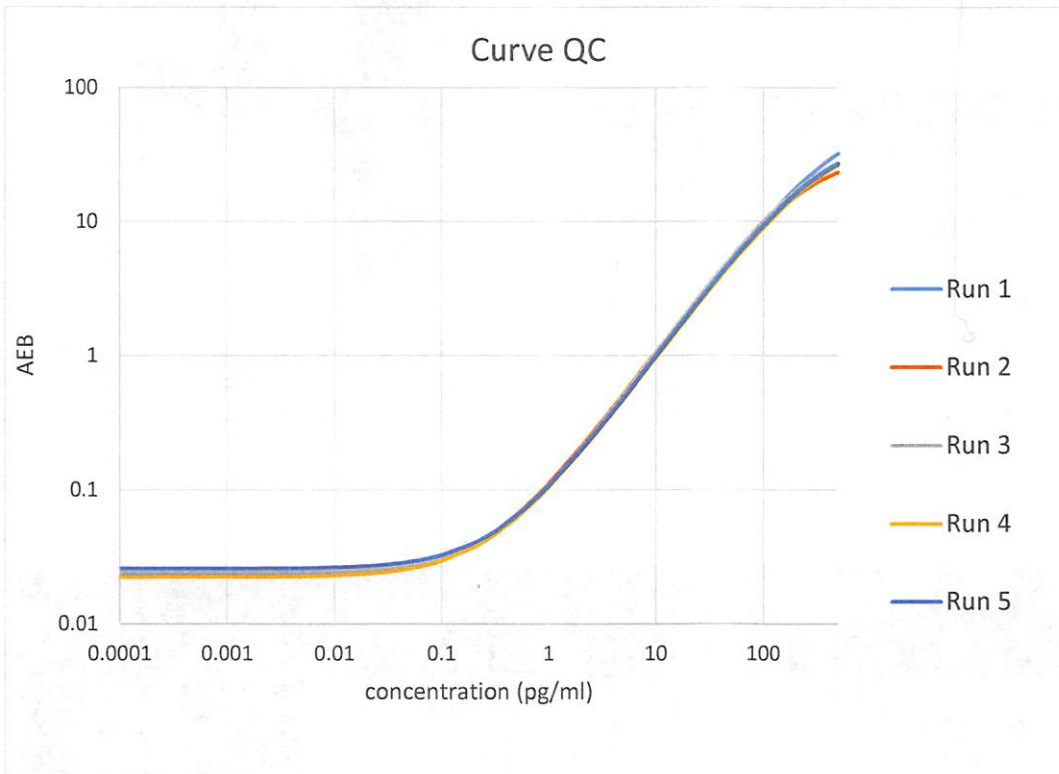
¹ RGP is not Kit Lot Specific

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	032204	3.34	1.71-4.23	
Control 2	032204	74.1	57.2-111	
Control 3	032304	0.175	< 3.00	

Sample	Lot	Bridging Result	Kit Release Result	Pass/ Fail
Panel 1	023202	Negative	Negative	Pass
Panel 2	023202	Negative	Negative	Pass
Panel 3	023202	Negative	Negative	Pass
Panel 4	023202	Negative	Negative	Pass
Panel 5	023202	Negative	Negative	Pass
Panel 6	023202	Positive	Positive	Pass
Panel 7	023202	Positive	Positive	Pass
Panel 8	023202	Positive	Positive	Pass
Panel 9	023202	Positive	Positive	Pass
Panel 10	023202	Positive	Positive	Pass
Panel 11	023202	Positive	Positive	Pass
Panel 12	023202	Positive	Positive	Pass

Review/Approval	Shivani Goel	QC Manager	Shivani Goel 02 Feb 2021
	Name	Title	Signature/Date

Calibration Curve



**Calibrator Levels
(ng/mL)**

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