

Product Number:	103769
Lot Number:	502437
Expiration:	22-Dec-2020
Platform(s):	HD-X

Component	Part Number	Lot Number
Bead Reagent	103758	019106
Detector Reagent	103760	017703
SBG Reagent	103762	017701
Sample Diluent	103763	017405
Calibrators	103766	017602
Control 1	103767	017603
Control 2	103768	017603
Control 3	103773	017603
RGP Reagent	103159	N/A ¹

¹ RGP is not Kit Lot Specific



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

Release Materials*				* Ranges shown are generated internally for new lot release only. Customer should generate their own control ranges.
Sample	Lot	Result (ng/mL)	Mean Range (ng/mL)	
Control 1	017603	417	234-584	
Control 2	017603	3830	2553-4367	
Control 3	017603	18330	16129-24193	
Sample	Lot	Result (ug/mL)		Range (ug/mL)
Panel 2	016209	0.835		0.706-1.07
Panel 3	016209	1.17		0.963-1.44
Panel 4	016209	0.964		0.759-1.14
Panel 6	016209	2.11		1.62-2.43
Panel 7	016209	5.95		4.41-7.38
Panel 8	016209	10.0		6.16-11.2
Panel 9	016209	30.0		15.6-43.7
Panel 10	016209	5.13		3.60-5.47
Control 5	013323	29.4		20.5-38.8
Control 6	013324	6.38		3.81-7.55
Panel 11	016209	0.089		Reported Result
Panel 12	016209	0.033		Reported Result
Panel 13	016209	0.064		Reported Result
Panel 14	016209	0.067		Reported Result
Panel 15	016209	0.101		Reported Result
Panel 16	016209	0.020		Reported Result
Panel 17	016209	0.039		Reported Result
Panel 18	016209	0.308		Reported Result
Panel 19	016209	0.094		Reported Result
Panel 20	016209	0.189		Reported Result

Review/Approval

Shivani Goel

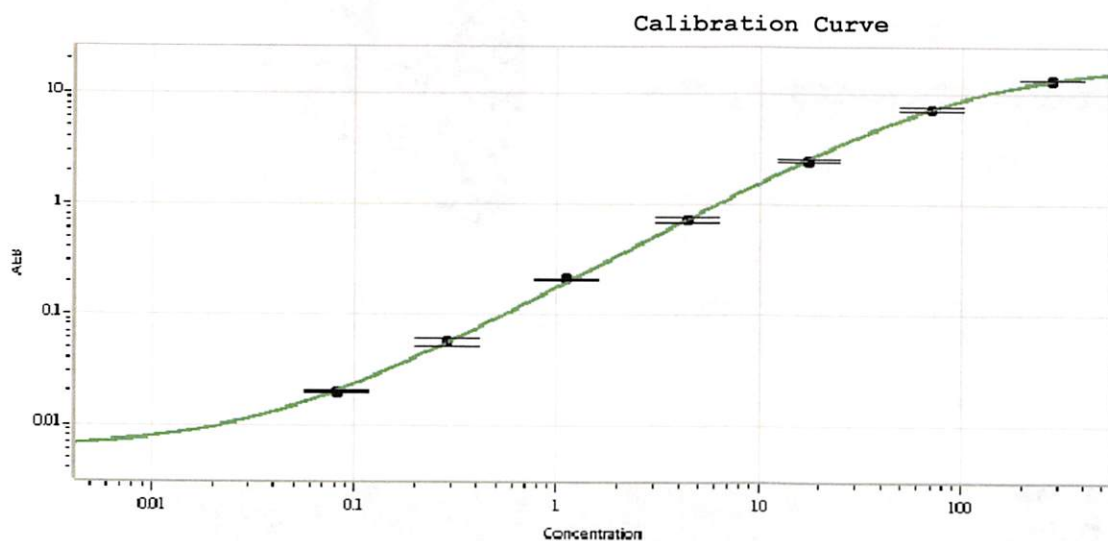
QC Manager

Name

Title

Shivani Goel 13Oct2020

Signature/Date



**Calibrator Levels
(ng/mL)**

A	0.000
B	0.082
C	0.287
D	1.11
E	4.39
F	17.5
G	69.9
H	280