

Product Number:	103769
Lot Number:	502579
Expiration:	6-Oct-2021
Platform(s):	HD-X
Product Number (Bulk):	103845
Lot Number (Bulk):	502605



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

Component	Part Number	Lot Number
Bead Reagent	103758	028007
Detector Reagent	103760	028114
SBG Reagent	103762	028117
Sample Diluent	103763	028702
Calibrators	103766	029501
Control 1	103767	029502
Control 2	103768	029502
Control 3	103773	029502
RGP Reagent	103159	N/A ¹

¹ RGP is not Kit Lot Specific

Release Materials*			
Sample	Lot	Result (ng/mL)	Mean Range (ng/mL)
Control 1	029502	631	424-679
Control 2	029502	4008	2694-4042
Control 3	029502	17751	11939-21841
Sample	Lot	Result (ug/mL)	Range (ug/mL)
Panel 1	013323	27.6	17.9-38.6
Panel 2	016209	0.985	0.700-1.07
Panel 3	016209	1.40	0.928-1.50
Panel 4	016209	1.13	0.730-1.20
Panel 5	013324	5.40	3.65-7.61
Panel 6	016209	2.28	1.63-2.50
Panel 7	016209	6.77	4.40-7.65
Panel 8	016209	11.4	5.58-12.5
Panel 9	016209	27.0	16.1-42.5
Panel 10	016209	5.22	3.56-5.70
Panel 11	016209	0.058	Reported Result
Panel 12	016209	0.023	Reported Result
Panel 13	016209	0.068	Reported Result
Panel 14	016209	0.059	Reported Result
Panel 15	016209	0.107	Reported Result
Panel 16	016209	0.012	Reported Result
Panel 17	016209	0.040	Reported Result
Panel 18	016209	0.289	Reported Result
Panel 19	016209	0.110	Reported Result
Panel 20	016209	0.186	Reported Result

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

Review/Approval

Shivani Goel

QC Manager

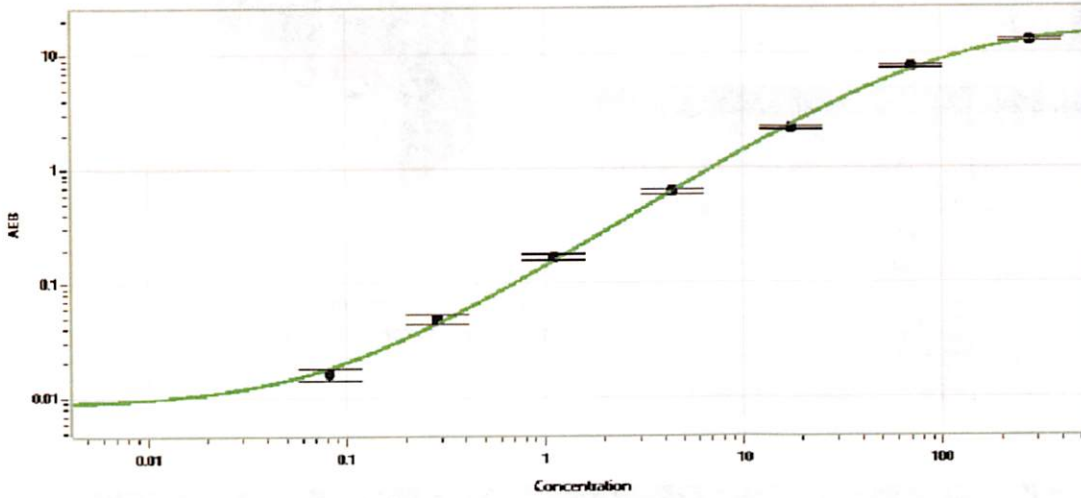
Shivani Goel 24 Nov 2020

Name

Title

Signature/Date

Calibration Curve



**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