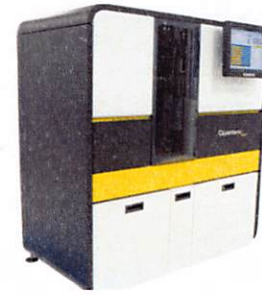


Product Number:	103624
Lot Number:	503456
Expiration:	19-Jul-2023
Platform(s):	HD-1/HD-X and SR-X



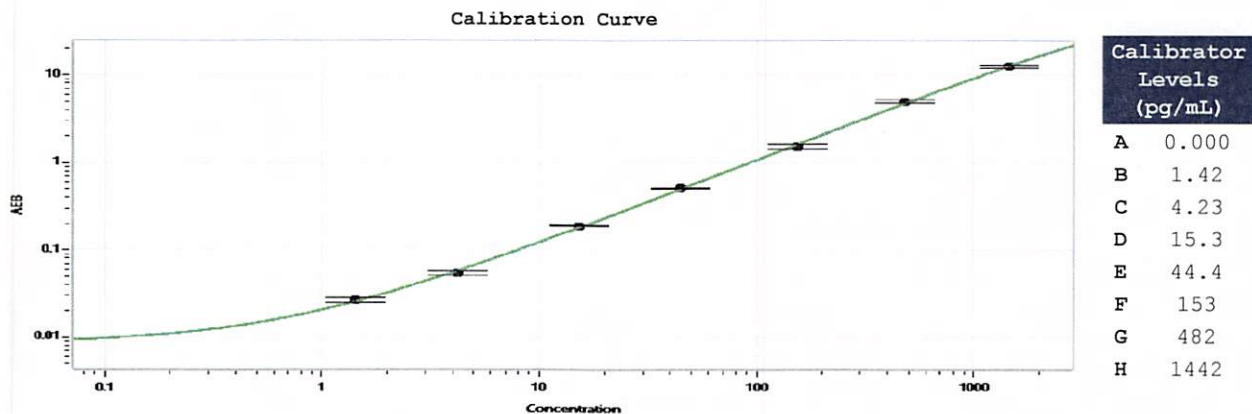
Data below represents results generated on the Simoa® HD-X Analyzer.

Component	Part Number	Lot Number
Bead Reagent	103617	233305
Detector Reagent	103619	221702
SBG Reagent	103620	221703
Sample Diluent	103621	220018
Calibrators	103893	220006
Control 1	103894	220007
Control 2	103895	220007
RGP Reagent	103159	N/A ¹

¹ Reagents are not Kit Lot Specific.

Release Materials			
	Lot	Result (pg/mL)	Spec. Range (pg/mL)
Control 1	220007	25.4	17.0 - 28.5
Control 2	220007	1286	946 - 1418
Panel 1	211003	85.0	42.6 - 99.6
Panel 2	211003	85.4	52.3 - 92.9
Panel 3	218001	438	224 - 480
Panel 4	218001	2569	1865 - 2797

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



Review/Approval

Matt Koulouris
Name

QC Supervisor
Title

M. Koulouris 25 Jan 2023
Signature/Date