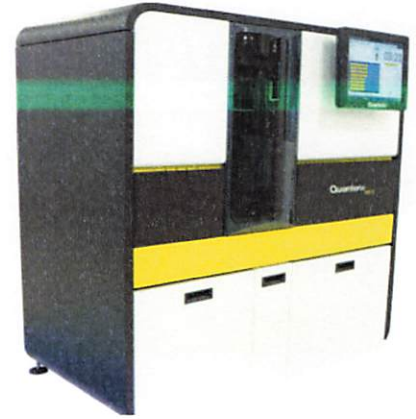


Product Number:	103337
Lot Number:	504248
Expiration:	4-Apr-2025
Platform(s):	SR-X and HD-1/HD-X



Component	Part Number	Lot Number
Bead Reagent	103334	328411
Detector Reagent	101091	328503
SBG Reagent	101101	328504
Sample Diluent	103336	328505
Calibrator Diluent	101095	327712
Calibrator Concentrate	101093	328513
RGP Reagent	103159	N/A ¹

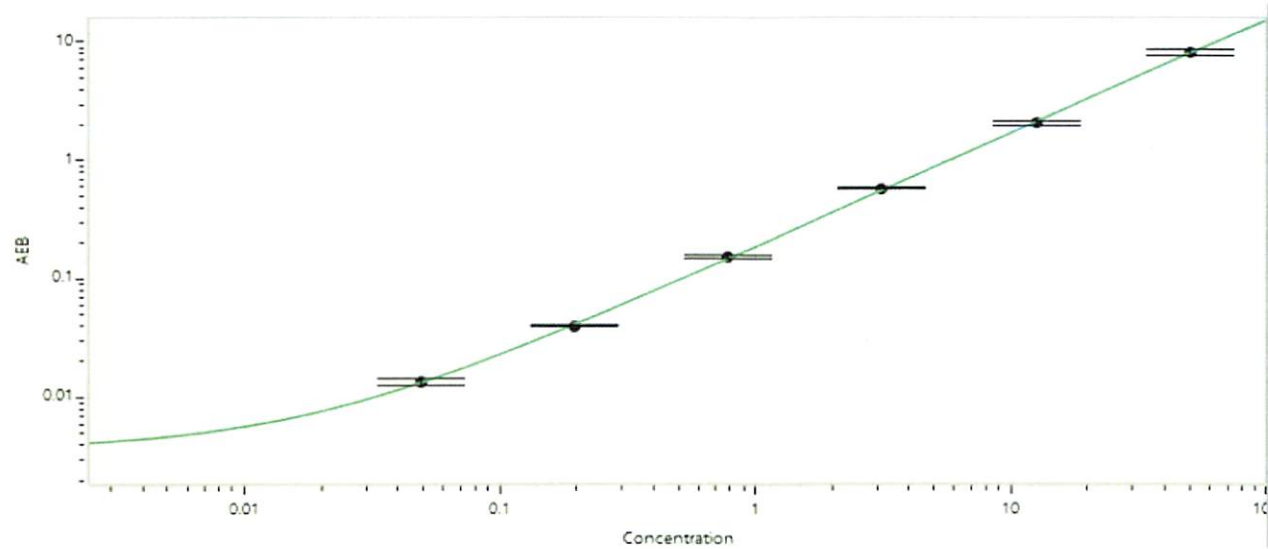
¹ Reagents are not Kit Lot Specific

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

Release Materials*			
	Lot	Result (pg/mL)	Spec. Range (pg/mL)
Control 1	328513	0.786	0.750 - 1.25
Control 2	328513	22.6	15.0 - 25.0
Panel 1	211906	1.94	1.26 - 2.41
Panel 2	211906	3.40	2.37 - 4.51
Panel 3	211906	33.6	21.0 - 45.2

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

Calibrator Concentrate Value Assignment (pg/mL)	2,499
---	-------



Calibrator Levels (pg/mL)	
A	0.000
B	0.049
C	0.195
D	0.781
E	3.13
F	12.5
G	50.0

Review/Approval

ROBERT SOMOGYI
Name

QC MANAGER
Title

[Signature]
Signature/Date
30 APR 2024