

Product Number:	103806
Lot Number:	503876
Expiration:	2-Feb-2024
Bulk Kit	
Product Number:	103872
Lot Number:	503761
Platform(s):	HD-X



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

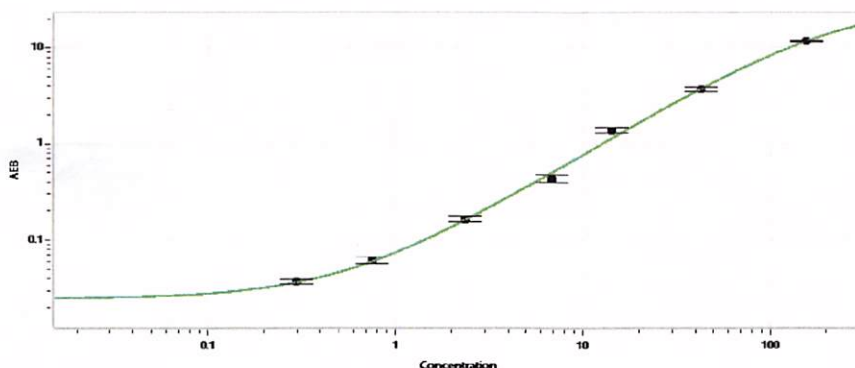
Component	Part Number	Lot Number
Bead Reagent	103794	303308
Detector Reagent	103796	303707
SBG Reagent	103798	303710
Sample Diluent	103800	303908
Calibrators	103803	305502
Control 1	103804	305503
Control 2	103805	305503
RGP Reagent	103159	N/A ¹

¹ RGP is not Kit Lot Specific

Release Materials*			
Sample	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	305503	8.18	5.66-9.17
Control 2	305503	116	93.6-140
Panel 1	129117	7.73	6.66-12.0
Panel 2	021111	47.6	34.9-91.4
Panel 3	021111	92.9	64.6-208

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

Calibration Curve



Calibrator Levels (pg/mL)

A	0.000
B	0.298
C	0.752
D	2.35
E	6.89
F	14.2
G	42.6
H	156

Review/Approval

Sonia Khillan QC Manager
Name Title

Soni D. 1/6 30 MAY 2023
Signature/Date