

Product Number:	103806
Lot Number:	503539
Expiration:	27-Jan-2023
Bulk Kit	
Product Number:	103872
Lot Number:	503552
Platform(s):	HD-X

Component	Part Number	Lot Number
Bead Reagent	103794	202711
Detector Reagent	103796	203312
SBG Reagent	103798	202712
Sample Diluent	103800	202713
Calibrators	103803	203015
Control 1	103804	203016
Control 2	103805	203016
RGP Reagent	103159	N/A ¹

¹ RGP is not Kit Lot Specific

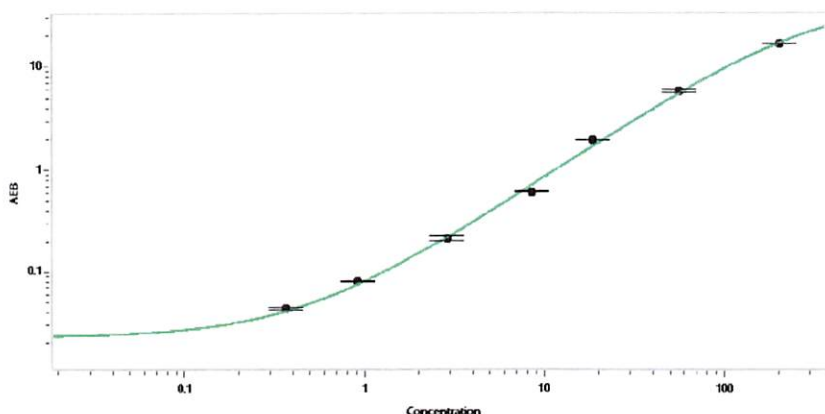


Data below represents results generated on the Simoa® HD-X Analyzer.

Release Materials*			
Sample	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	203016	6.81	6.50-9.75
Control 2	203016	116	116-173
Panel 1	021305	3.38	2.55-4.27
Panel 2	021111	53.2	42.2-90.4
Panel 3	021111	116	89.1-202

* 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	0.365
C	0.923
D	2.88
E	8.45
F	18.6
G	55.6
H	204

Review/Approval

Sonia Khillan

QC Manager

Name

Title

Signature/Date

Sonia Khillan / 21 JUL 2022