

Product Number:	103575
Lot Number:	503521
Expiration:	3-May-2023
Platform(s):	HD-1 / HD-X & SR-X

Component	Part Number	Lot Number
Bead Reagent	103564	217210
Detector Reagent	103566	217406
SBG Reagent	103568	217205
Sample Diluent	103569	217214
Calibrator Diluent	103571	214213
Calibrator Concentrate	103570	216901
SR-X SBG Diluent	103158	N/A <sup>1</sup>
RGP Reagent	103159	N/A <sup>1</sup>

<sup>1</sup> Generic items are not Kit Lot Specific



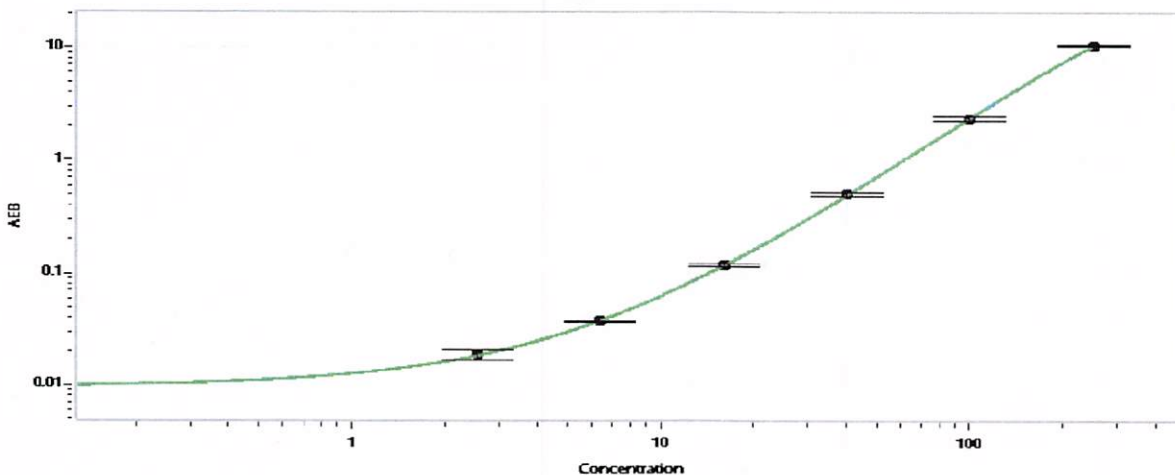
Data below represents results generated on the Simoa® HD-X Analyzer.

Release Materials*			
	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	216901	48.0	37.5-62.5
Control 2	216901	367	300-500
Panel 1	020204	63.8	48.2-72.3
Panel 2	020204	79.3	59.5-89.3
Panel 3	102806	479	363-663

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

Calibrator Concentrate Value Assignment (pg/mL)	4171
---	------

Calibration Curve



Calibrator Levels (pg/mL)	
A	0.000
B	2.56
C	6.40
D	16.0
E	40.0
F	100
G	250

Review/Approval

*Sonia Khillan*  
Name

*QC Manager*  
Title

*Sonia Khillan / 20 JUL 2022*  
Signature/Date