

Product Number:	103575
Lot Number:	503835
Expiration:	10-Mar-2024
Platform(s):	HD-1 / HD-X & SR-X



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

Component	Part Number	Lot Number
Bead Reagent	103564	308108
Detector Reagent	103566	308207
SBG Reagent	103568	308208
Sample Diluent	103569	308210
Calibrator Diluent	103571	306908
Calibrator Concentrate	103570	309702
SR-X SBG Diluent	103158	N/A <sup>1</sup>
RGP Reagent	103159	N/A <sup>1</sup>

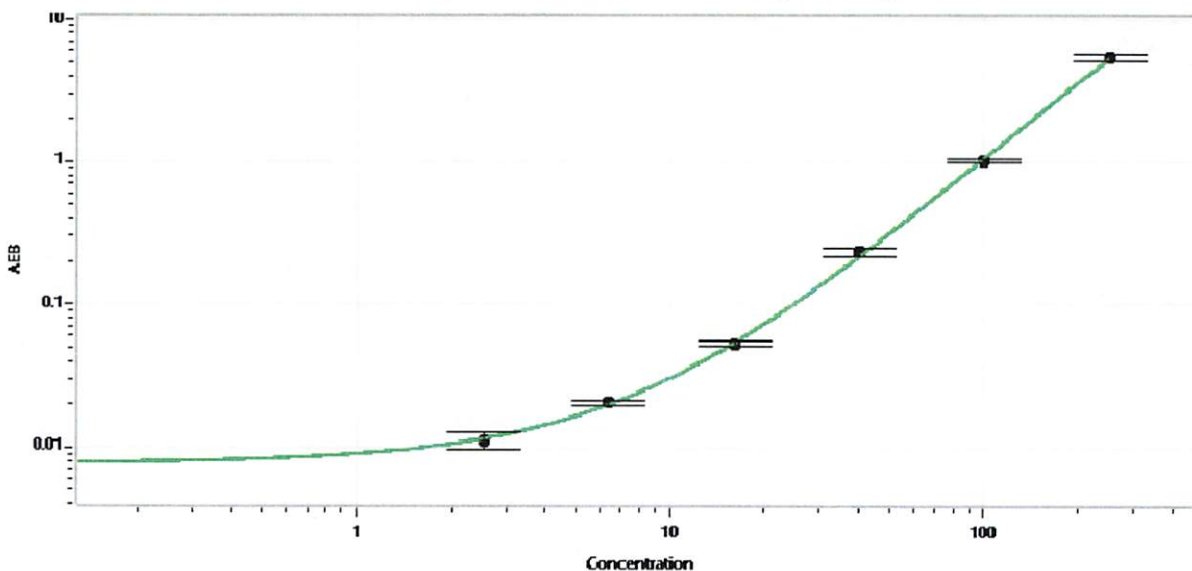
<sup>1</sup> Generic items are not Kit Lot Specific

Release Materials*			
	Lot	Result (pg/mL)	Spec. Range (pg/mL)
Control 1	23306	56.0	41.2 - 61.7
Control 2	23306	415.1	344 - 515
Panel 1	313805	41.3	30.9 - 46.3
Panel 3	313805	537.3	413 - 619

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

Calibrator Concentrate Value Assignment (pg/mL)	4195
---	------

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

ROBERT SOMOSI, MANAGER, QC  
Name Title

[Signature] 06 JUN 2023  
Signature/Date