

Product Number:	101552
Lot Number:	503724
Expiration:	21-Oct-2023
Platform(s):	HD-1/HD-x & SR-X



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

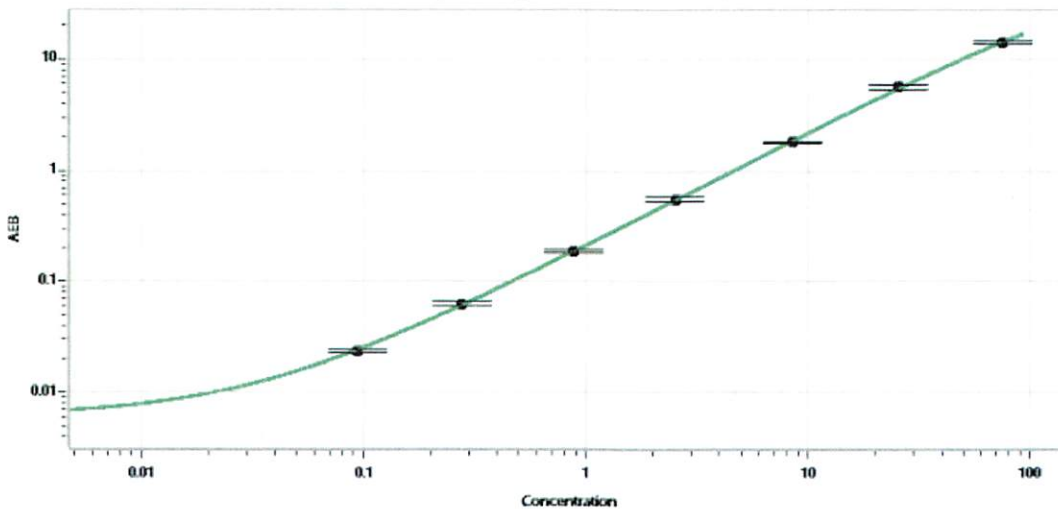
Component	Part Number	Lot Number
Bead Reagent	101554	229402
Detector Reagent	101553	231202
SBG Reagent	101555	229807
Sample Diluent	101556	230413
Calibrators	101559	235311
Control 1	101560	235312
Control 2	101561	235312
SBG Diluent (SR-X)†	103158	N/A <sup>1</sup>
RGP Reagent	103159	N/A <sup>1</sup>

<sup>1</sup> RGP is not Kit Lot Specific

Release Materials*			
	Lot	Result (pg/mL)	Spec. Range (pg/mL)
Control 1	235312	0.983	0.610 - 1.13
Control 2	235312	50.3	40.8 - 61.2
Panel 1	232607	0.822	0.491 - 1.52
Panel 2	232607	0.757	0.516 - 1.53
Panel 3	234908	20.1	16.7 - 25.1
Panel 4	234908	29.9	25.6 - 38.5

\* 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	0.094
C	0.276
D	0.884
E	2.53
F	8.55
G	25.5
H	74.6

Review/Approval

*Sonia Khillan*  
Name

QC  
Title

*Sonia Khillan / 17 JAN 2023*  
Signature/Date