

<b>Product Number:</b>	Individual: 103806; Bulk: 103872
<b>Lot Number:</b>	Individual: 504228; Bulk: 504295
<b>Expiration Date:</b>	24-Feb-25
<b>Platform(s):</b>	HD-X



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

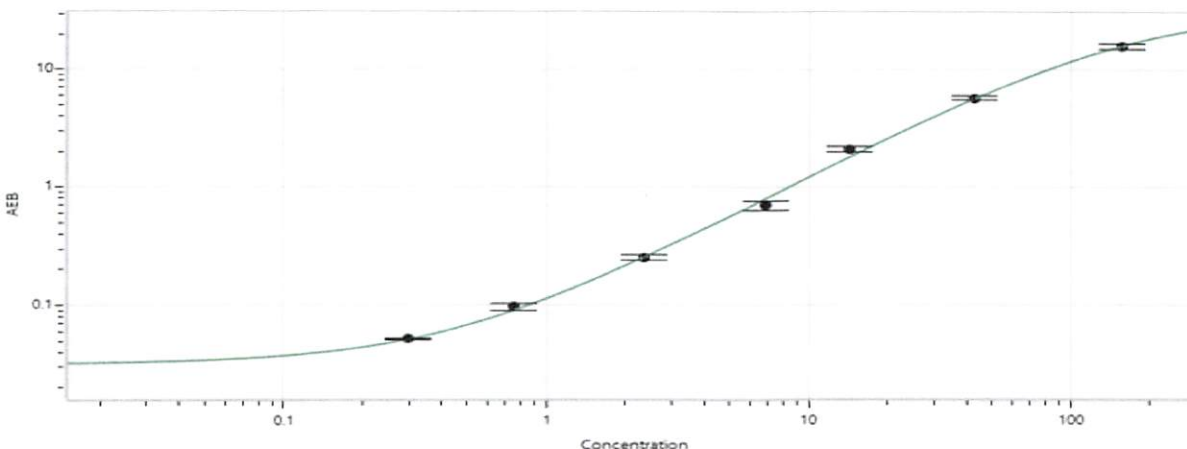
Component	Part Number	Lot Number
Bead Reagent	103794	411401
Detector Reagent	103796	411402
SBG Reagent	103798	411403
Sample Diluent	103800	411404
Calibrators	103803	305502
Control 1	103804	305503
Control 2	103805	305503
RGP Reagent	103159	N/A <sup>1</sup>

<sup>1</sup> Reagents are not Kit Lot Specific

Release Materials*			
	Lot	Result (pg/mL)	Spec. Range (pg/mL)
Control 1	305503	6.93	5.66 - 9.17
Control 2	305503	119	93.6 - 140
Panel 1	414901	4.52	4.12 - 6.18
Panel 2	412010	47.7	45.3 - 68.0
Panel 3	414901	105	92.8 - 139

\* 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

Cody Carlson

QC Manager

*[Signature]* 24 JUN 2024

Name

Title

Signature/Date