

Product Number:	100196
Lot Number:	503906
Expiration:	18-Jul-2024
Platform(s):	HD-1/HD-X



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

Component	Part Number	Lot Number	Control Number
Bead Reagent	100391	324306	N/A
Detector Reagent	100315	324805	N/A
SBG Reagent	100316	324803	N/A
Sample Diluent	100317	319916	319916A
Calibrator Diluent	100665	324009	N/A
Calibrator Concentrate	100664	320709	N/A
RGP Reagent	103159	N/A ¹	N/A ¹

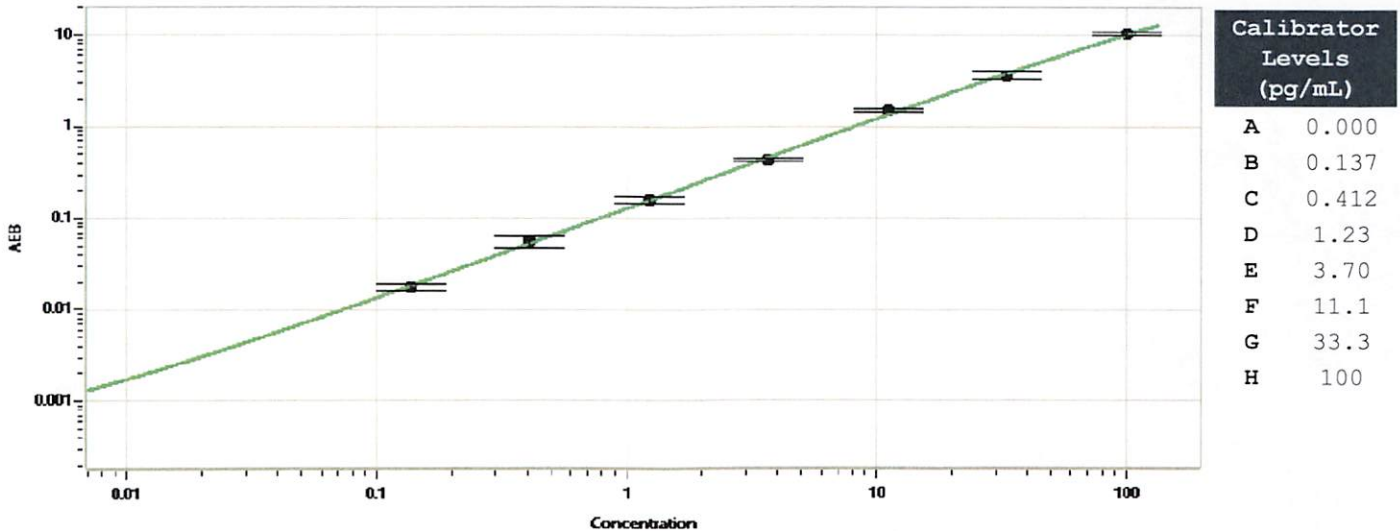
¹ RGP is not Kit Lot Specific

Release Materials*			
	Lot	Result (pg/mL)	Spec. Range (pg/mL)
Control 1	123410	8.30	5.34 - 13.4
Control 2	123410	83.7	51.9 - 143
Panel 1	029002	1.61	1.19 - 2.82
Panel 2	029002	7.42	5.13 - 14.0
Panel 3	029002	81.1	46.8 - 134

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

Calibrator Concentrate Value Assignment (pg/mL)	1910
---	------

Calibration Curve



Review/Approval

ROBERT SOMOGYI
Name

QC MANAGER
Title

[Signature] 14SEP2023
Signature/Date