

Product Number:	103186
Lot Number:	502076
Expiration:	8-Oct-2020
Platform(s):	HD-1/HD-x



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

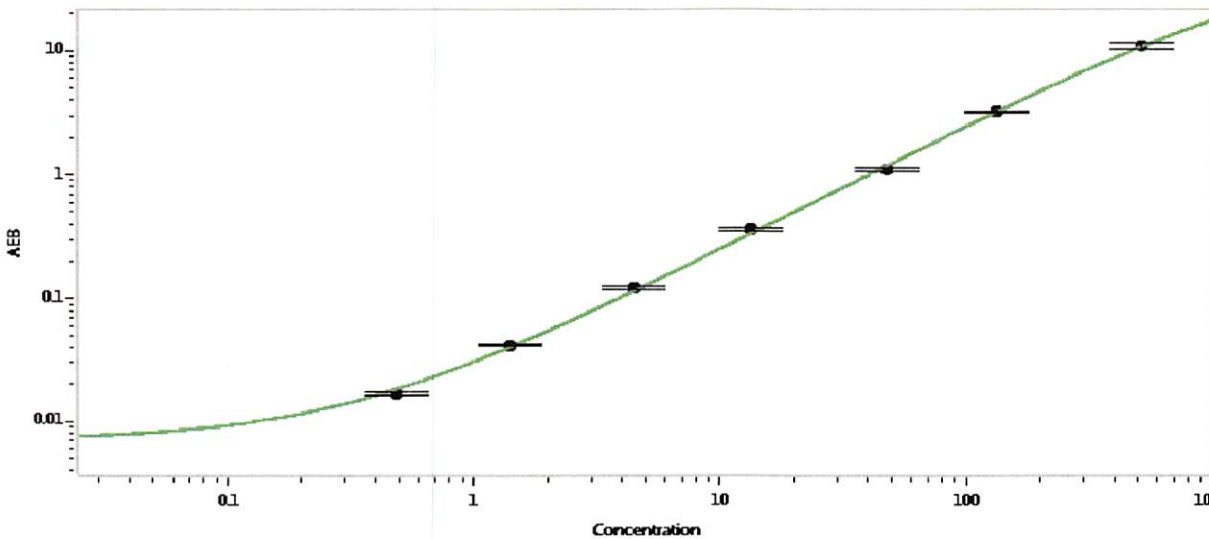
Component	Part Number	Lot Number
Bead Reagent	102246	928701
Detector Reagent	102248	928805
SBG Reagent	102250	928704
Sample Diluent	102252	929010
Calibrators	102255	929601
Controls	102256 - 102257	929703
RGP Reagent	103159	N/A ¹

¹ RGP is not Kit Lot Specific

Release Materials*			
	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	929703	4.10	3.05-4.58
Control 2	929703	178	130-194
Panel 1a	922601	21.0	14.1-21.1
Panel 1b	922602	6.80	4.94-7.41
Panel 2	922603	42.2	28.4-42.6
Panel 3	922604	174	129-204

* 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	0.488
C	1.40
D	4.47
E	13.4
F	47.5
G	133
H	514

Review/Approval: Patrizia Stadler Name, Assoc. Director, QC Title, P. Stadler Signature, 06 NOV 2019 Date