

Product Number:	103186
Lot Number:	502495
Expiration:	6-Apr-2021
Platform(s):	HD-1/ HD-X & SR-X

Component	Part Number	Lot Number
Bead Reagent	102246	022302
Detector Reagent	102248	022404
SBG Reagent	102250	022403
Sample Diluent	102252	009702
Calibrators	102255	022401
Control 1	102256	022402
Control 2	102257	022402
RGP Reagent	103159	N/A ¹

¹ RGP is not Kit Lot Specific

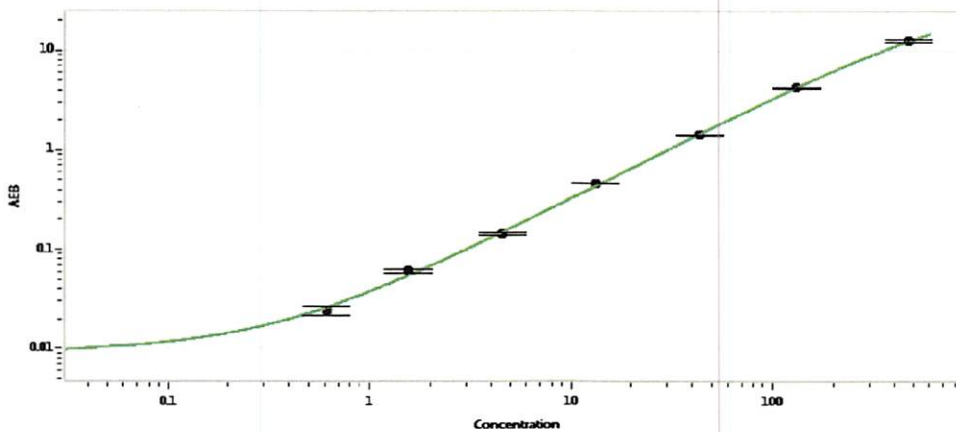


Data below represents results generated on the Simoa HD-X® Analyzer.

Release Materials*			
	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	022402	5.05	3.33-5.59
Control 2	022402	179	137-205
Panel 1	922601	18.7	14.1-21.1
Panel 2	922602	7.05	4.94-7.41
Panel 3	922603	42.6	28.4-42.6
Panel 4	922604	178	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	0.609
C	1.56
D	4.56
E	13.1
F	43.1
G	131
H	468

Review/Approval: Shivani Goel QC Manager Shivani Goel 04 Sep 2020.
Name Title Signature/Date