

Product Number:	103180
Lot Number:	502547
Expiration:	14-Sep-2021
Platform(s):	HD-1/ HD-X & SR-X

Component	Part Number	Lot Number
Bead Reagent	102246	026804
Detector Reagent	102248	026904
SBG Reagent	102250	026908
Sample Diluent	102252	026905
Calibrators	102255	933601
Control 1	102256	933901
Control 2	102257	933901
RGP Reagent	103159	N/A ¹

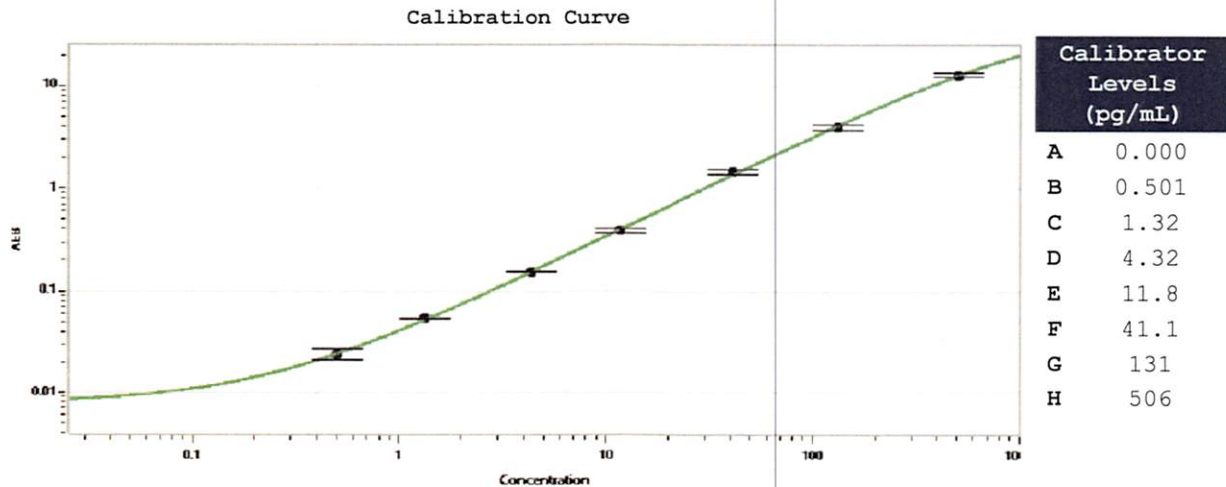
¹ RGP is not Kit Lot Specific



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

Release Materials*			
	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	933901	3.27	2.60-4.70
Control 2	933901	151	122-183
Panel 1	008501	13.4	10.5-15.7
Panel 2	008501	9.88	8.22-12.3
Panel 3	008501	61.6	49.1-73.6
Panel 4	009401	214	151-234

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



Review/Approval

Patrizia Stadler Assoc. Director, QC

P. Stadler 29 Sep 2020

Name

Title

Signature/Date