

Product Number:	103377
Lot Number:	502345
Expiration:	11-Jan-2021
Platform(s):	HD-1/HD-X and SR-X

Component	Part Number	Lot Number
Bead Reagent	103364	010905
Detector Reagent	103366	009501
SBG Reagent	103368	009302
Sample Diluent	103370	009305
Calibrators	103374	901102
Controls	103375/103376	901103/901801
SBG Diluent (SRX)†	103158	N/A ¹
RGP Reagent	103159	N/A ¹

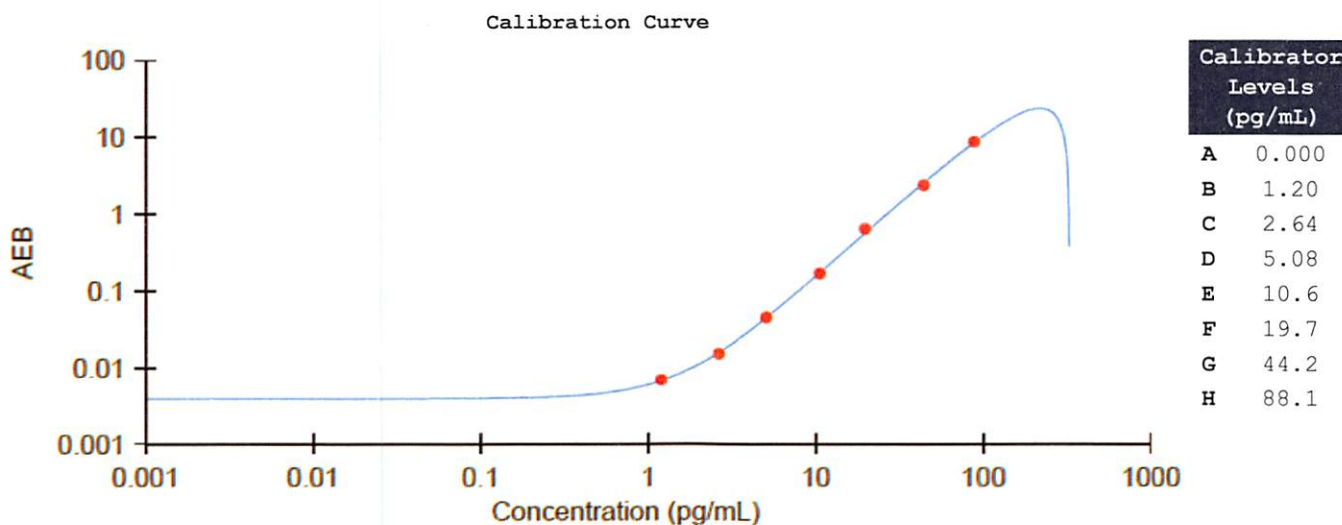
¹ Reagents are not Kit Lot Specific.

† Specific to use on the SRX platform - refer to the SRX Kit Instructions (KI-0045).



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

Release Materials				Ranges shown are generated internally for new lot release only. Customer should generate their own control ranges.
	Lot	Result (pg/mL)	Mean Range (pg/mL)	
Control 1	901103	21.4	16.1-24.2	
Control 2	901801	159	125-188	
Panel 1	900701	40.2	29.7-44.5	
Panel 2	900702	48.7	35.6-53.4	
Panel 3	900703	27.1	20.7-31.1	



Review/Approval

Shivani Goel
Name

QC Manager
Title

Shivani Goel 30Apr2020
Signature/Date