

Product Number:	103377
Lot Number:	502668
Expiration:	20-Nov-2021
Platform(s):	HD-1/HD-X and SR-X



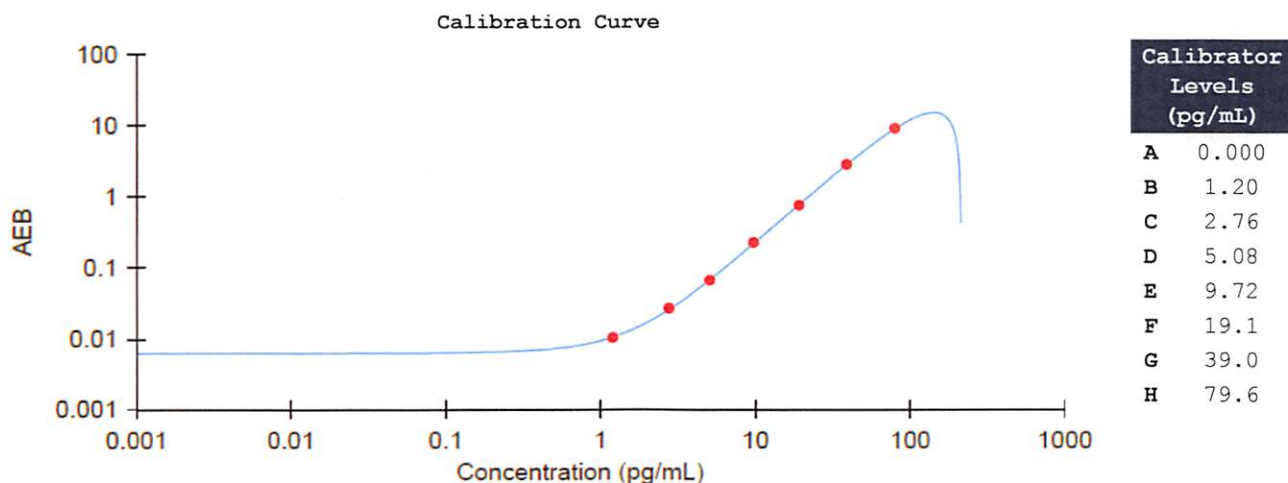
Data below represents results generated on the Simoa™ SR-X Analyzer.

Component	Part Number	Lot Number
Bead Reagent	103364	035302
Detector Reagent	103366	035608
SBG Reagent	103368	035607
Sample Diluent	103370	035709
Calibrators	103374	035701
Control 1	103375	035702
Control 2	103376	035702
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).

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	035702	18.1	14.6-21.9	
Control 2	035702	154	122-182	
Panel 1	900701	36.5	29.6-44.5	
Panel 2	900702	43.8	35.5-53.3	
Panel 3	900703	26.8	20.8-31.2	



Review/Approval

Shivani Goel
Name

QC Manager
Title

Shivani Goel 15 Jan 2021
Signature/Date