

Product Number: 103377
Lot Number: 501754
Expiration: 6-Mar-2020
Platform(s): HD-1 and SR-X

Component	Part Number	Lot Number
Bead Reagent	103364	906513
Detector Reagent	103366	906701
SBG Reagent	103368	906802
Sample Diluent	103370	906602
Calibrators	103374	909401
Controls	103375/103376	909402
SBG Diluent (SRX) †	103158	N/A ¹
RGP Reagent	103159	N/A ¹

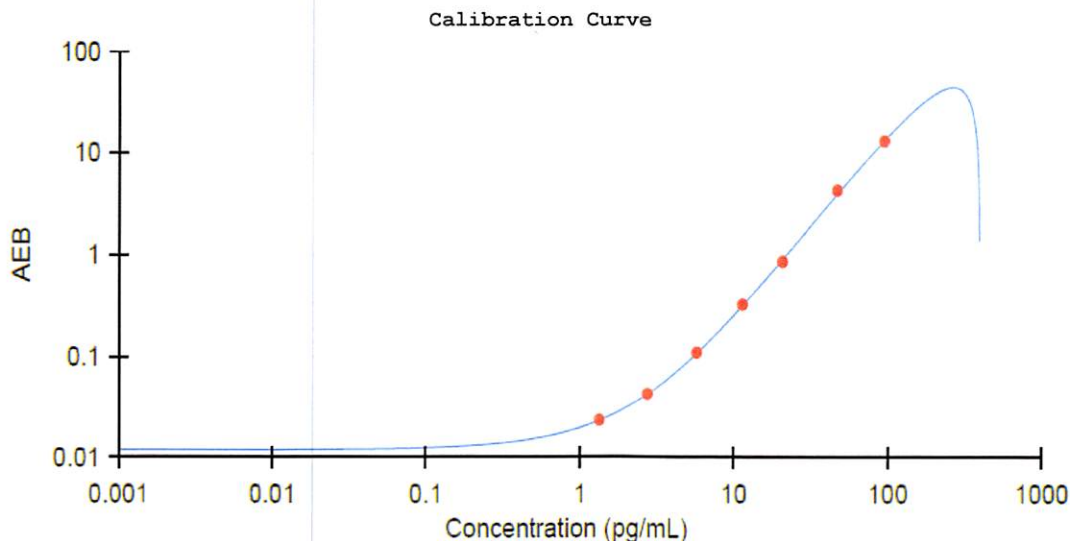
¹ Reagents are not Kit Lot Specific.

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



Data below represents results generated on the Quanterix SR-X™ Ultra-Sensitive Biomarker Detection System

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	909402	19.9	16.0-24.0	
Control 2	909402	175	133-200	
Panel 1	900701	36.8	29.8-44.7	
Panel 2	900702	43.9	35.2-52.8	
Panel 3	900703	22.9	20.7-31.0	



Review/Approval

Patrizia Stadler
Name

QC Assoc. Director
Title

p. Stodler

14 May 2019

Signature/Date