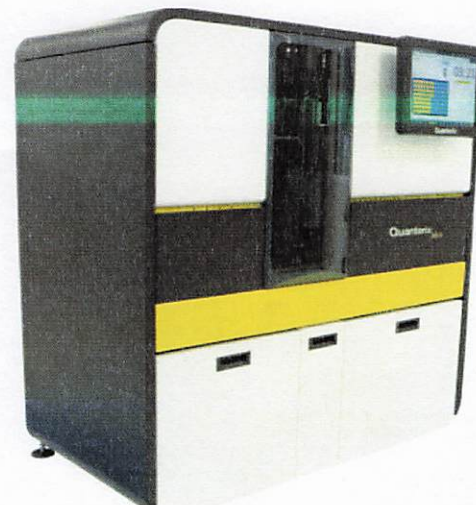


<b>Product Number:</b>	103293
<b>Lot Number:</b>	504000
<b>Expiration:</b>	22-Jan-2025
<b>Platform(s):</b>	SR-X and HD-1/HD-X



Component	Part Number	Lot Number
Bead Reagent	103281	401821
Detector Reagent	103283	402213
SBG Reagent	103285	402214
Sample Diluent	103286	401618
Calibrator Diluent	103288	402215
Calibrator Concentrate	103287	407801
SBG Diluent (SR-X)†	103158	N/A <sup>1</sup>
RGP Reagent	103159	N/A <sup>1</sup>

<sup>1</sup> Reagents are not Kit Lot Specific

† Specific to use on the SRX platform.

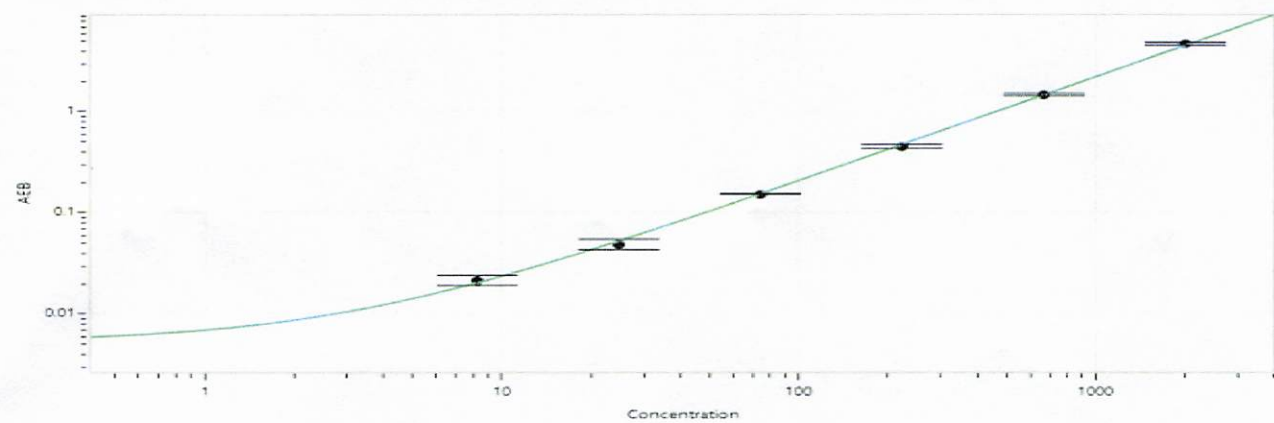
Refer to the Simoa TDP-43 Kit Instructions SR-X (KI-0132).

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

Release Materials*			
	Lot	Result (pg/mL)	Spec. Range (pg/mL)
Control 1	407801	199	150 - 250
Control 2	407801	1789	1500 - 2500
Panel 1	406803	76.2	59.4 - 103
Panel 2	406803	134	101 - 177
Panel 3	406803	256	207 - 311
Panel 4	406803	4218	3130 - 4695

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

<b>Calibrator Concentrate Value Assignment (pg/mL)</b>	26104
--	-------



Calibrator Levels (pg/mL)	
A	0.000
B	8.23
C	24.7
D	74.1
E	222
F	667
G	2000

Review/Approval

Cody Carlson  
Name

QC Manager  
Title

*Cody Carlson* 27 MAR 2024  
Signature/Date