

<b>Product Number:</b>	101736
<b>Lot Number:</b>	503784
<b>Expiration:</b>	14-Nov-2023



*Data below represents results generated on the Simoa® HD-1 / HD-X Analyzer.*

Component	Part Number	Lot Number
RGP Reagent	103159	23180

Means of Analysis	PASS/FAIL
HD-1/HD-X	Pass

Review/Approval

*Matt Kodouris*

*QC Supervisor*

*[Signature]* 10 Jan 2023

Name

Title

Signature/Date