

<b>Product Number:</b>	101736
<b>Lot Number:</b>	503466
<b>Expiration:</b>	28-Mar-2023



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

Component	Part Number	Lot Number
RGP Reagent	103159	20450
Means of Analysis	PASS/FAIL	
HD-1/HD-X	Pass	

Review/Approval

Richa Shrestha  
Name

QC Supervisor  
Title

*[Signature]* 28 Apr 2022  
Signature/Date