

Product Number:	101736
Lot Number:	503783
Expiration:	28-Nov-2023



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

Component	Part Number	Lot Number
RGP Reagent	103159	23320
Means of Analysis	PASS/FAIL	
HD-1/HD-X	Pass	

Review/Approval

<i>Matt Koulouris</i>	<i>QC Supervisor</i>	<i>[Signature]</i> 10Jan2023
Name	Title	Signature/Date