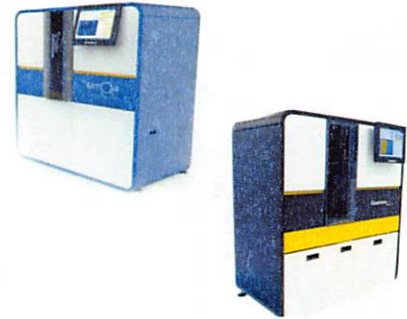


<b>Product Number:</b>	101736
<b>Lot Number:</b>	503425
<b>Expiration:</b>	27-Feb-2023



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

Component	Part Number	Lot Number
RGP Reagent	103159	20580

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

Review/Approval

Shivani Goel  
Name

QC Manager  
Title

Shivani Goel 04Apr 2022  
Signature/Date