

<b>Product Number:</b>	101736
<b>Lot Number:</b>	502776
<b>Expiration:</b>	11-Jan-2022
<b>Storage:</b>	2-8°C

Component	Part Number	Lot Number
RGP Reagent	103159	10110

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

Review/Approval

*Shivani Goel*

*QC Manager*

*Shivani Goel 10 Mar 2021*

Name

Title

Signature/Date