

<b>Product Number:</b>	101736
<b>Lot Number:</b>	503113
<b>Expiration:</b>	2-Aug-2022
<b>Storage:</b>	2-8°C

<b>Component</b>	<b>Part Number</b>	<b>Lot Number</b>
RGP Reagent	103159	12140

<b>Means of Analysis</b>	<b>PASS/FAIL</b>
HD-1/HD-X	Pass

Review/Approval

Shivani Goel  
Name

QC

Title

Shivani Goel 20 Aug 2021  
Signature/Date