

<b>Product Number:</b>	101736
<b>Lot Number:</b>	503052
<b>Expiration:</b>	30-Jun-2022
<b>Storage:</b>	2-8°C

<b>Component</b>	<b>Part Number</b>	<b>Lot Number</b>
RGP Reagent	103159	11810

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

Review/Approval

*Junpil Park*  
Name

QC Supervisor  
Title

*[Signature]* 16 Jul 2021  
Signature/Date