

<b>Product Number:</b>	101736
<b>Lot Number:</b>	502858
<b>Expiration:</b>	15-Dec-2021
<b>Storage:</b>	2-8 °C

Component	Part Number	Lot Number
RGP Reagent	103159	03500

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

Review/Approval

*Junpil Park*  
Name

QC Supervisor  
Title

*[Signature]* 08 Apr 2021  
Signature/Date