

<b>Product Number:</b>	101736
<b>Lot Number:</b>	502522
<b>Expiration:</b>	3-Aug-2021
<b>Storage:</b>	2-8°C

Component	Part Number	Lot Number
RGP Reagent	103159	02160

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

Review/Approval

Shivani Goel  
Name

QC Manager  
Title

Shivani Goel 02 Sep 2020  
Signature/Date