

Product Number:	101736
Lot Number:	502560
Expiration:	14-Sep-2021
Storage:	2-8°C

Component	Part Number	Lot Number
RGP Reagent	103159	02580

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

Review/Approval

Shivani Goel
Name

QC Manager
Title

Shivani Goel 05 Oct 2020
Signature/Date