

Product Number:	101736
Lot Number:	502880
Expiration:	6-Apr-2022
Storage:	2-8°C

Component	Part Number	Lot Number
RGP Reagent	103159	10960

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

Review/Approval

Shivani Goel
Name

QC Manager
Title

Shivani Goel 16Apr 2021
Signature/Date