

Product Number:	101736
Lot Number:	503192
Expiration:	16-Sep-2022
Storage:	2-8°C

Component	Part Number	Lot Number
RGP Reagent	103159	12590

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

Review/Approval

Shivani Goel
Name

QC

Title

Shivani Goel 22Oct 2021
Signature/Date