

Product Number:	101736
Lot Number:	503152
Expiration:	17-Jun-2022
Storage:	2-8 °C

Component	Part Number	Lot Number
RGP Reagent	103159	11680

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

Review/Approval Shivani Goel QC Shivani Goel 27 Sep 2021
Name Title Signature/Date