

Product Number:	101736
Lot Number:	502703
Expiration:	15-Dec-2021
Storage:	2-8°C

Component	Part Number	Lot Number
RGP Reagent	103159	03500

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

Review/Approval

Shivani Goel
Name

QC, Manager
Title

Shivani Goel 28 Jan 2021
Signature/Date