

Product Number:	101736
Lot Number:	503357
Expiration:	6-Dec-2022
Storage:	2-8°C

Component	Part Number	Lot Number
RGP Reagent	103159	13400

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

Review/Approval

Shivani Goel
Name

QC

Title

Shivani Goel 17 Feb 2022

Signature/Date