

Product Number:	101736
Lot Number:	503267
Expiration:	15-Nov-2022
Storage:	2-8°C

Component	Part Number	Lot Number
RGP Reagent	103159	13190

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

Review/Approval

Shivani Goel
Name

QC
Title

Shivani Goel 14 Dec 2021
Signature/Date