

Product Number:	101736
Lot Number:	504555
Expiration Date:	3-Sep-2026

Component	Part Number	Lot Number
RGP Reagent	103159	50620


Means of Analysis	PASS / FAIL
HD-X	PASS



Review/Approval

Dipesh Patel

Sr. Doc control

 17 MAR 2025

Name

Title

Signature/Date