

Product Number:	101736
Lot Number:	504620
Expiration Date:	9-Nov-2026

Component	Part Number	Lot Number
RGP Reagent	103159	51290

Means of Analysis	PASS / FAIL
HD-X	PASS



Review/Approval

Dipesh Patel Sr. Doc. Control - QA at 16 MAY 2025
 Name Title Signature/Date