

Product Number:	101736
Lot Number:	504469
Expiration Date:	3-Mar-2026



Component	Part Number	Lot Number
RGP Reagent	103159	42470

Means of Analysis	PASS / FAIL
HD-X	PASS

Review/Approval

Dhruvi Patel *Doc Control Specialist I, QA* *DP 04 Dec 2024*

Name

Title

Signature/Date