

Product Number:	101736
Lot Number:	504542
Expiration Date:	27-Jul-26



Component	Part Number	Lot Number
RGP Reagent	103159	50270

Means of Analysis	PASS / FAIL
HD-X	PASS

Review/Approval

Kyle Givoux

Senior Document Control Specialist, QA

Kyle Givoux 25 Feb 2025

Name

Title

Signature/Date