

Product Number:	101736
Lot Number:	504701
Expiration Date:	13-Dec-2026

Component	Part Number	Lot Number
RGP Reagent	103159	51640

Means of Analysis	PASS / FAIL
HD-X	PASS



Review/Approval

Dhruvi Patel

QA Specialist II, QA

EDP / 21 AUG 2025

Name

Title

Signature/Date