

Product Number:	101736
Lot Number:	504453
Expiration Date:	9-Jan-2026



Component	Part Number	Lot Number	Control Lot Number
RGP Reagent	103159	41910	41910A

Means of Analysis	PASS / FAIL
HD-X	PASS

Review/Approval

Dipesh Patel  
Name

Doc control QA  
Title

[Signature] 12 Nov 2024  
Signature/Date