

Product Number:	101736
Lot Number:	504380
Expiration Date:	2-Nov-2025

Component	Part Number	Lot Number
RGP Reagent	103159	41230

Means of Analysis	PASS / FAIL
HD-X	PASS



Review/Approval

Dipesh Patel
Name

Doc control-II QA
Title

cut / 10 sep 2024
Signature/Date