

Product Number:	101736
Lot Number:	536635
Expiration Date:	27-Feb-2027

Component	Part Number	Lot Number	Control Lot Number
RGP Reagent	103159	52390	52390A

Means of Analysis	PASS / FAIL
HD-X	PASS



Review/Approval	<u>Dilish Patel</u>	<u>Sr. QA Specialist</u>	<u>[Signature]</u> 06 Nov 2025
	Name	Title	Signature/Date