

Product Number:	103159
Lot Number:	50640
Control Lot Number:	50640A
Expiration Date:	5-Sep-2026
Storage:	2-8°C



Laboratory Analysis	
Means of Analysis	PASS / FAIL
HD-X	PASS

Review/Approval

Cody CarlsonQC ManagerSignature/Date22 Jul 2025