

Product Number:	103159
Lot Number:	40360
Expiration Date:	5-Aug-2025
Storage:	2-8°C



Laboratory Analysis

Means of Analysis	PASS / FAIL
HD-X	PASS

Review/Approval

ROBERT SONOGYI

QC MANAGER

[Handwritten Signature] 04/30/2024

Name

Title

Signature/Date