

Product Number:	103159
Lot Number:	41080
Expiration Date:	17-Oct-2025
Storage:	2-8°C



Laboratory Analysis

Means of Analysis	PASS / FAIL
HD-X	PASS

Review/Approval

ROBERT SOMOGYI QC MANAGER
 Name Title

TES 04 SEP 2024
 Signature/Date