

Product Number:	100457
Lot Number:	320109
Expiration:	20-Jul-2024
Storage:	2-8°C

Laboratory Analysis

Characteristic	Result	pH Target	PASS / FAIL
pH	7.36	7.35 - 7.45	PASS

Review/Approval

Sonia Khillan
Name

Senior QC Analyst
Title

Sonia Khillan 26 Oct 2023
Signature/Date