

Product Number:	100457
Lot Number:	206117
Expiration:	2-Mar-2023
Storage:	2-8°C

Laboratory Analysis

Characteristic	Result	pH Target
pH Measurement	PASS	7.35 - 7.45

Review/Approval Shivani Goel QC Manager Shivani Goel 08 Mar 2022
Name Title Signature/Date