

Product Number:	101631
Lot Number:	211101
Expiration:	21-Apr-2023
Storage:	2 - 8°C

Laboratory Analysis

Characteristic	Result	pH Target
pH	Pass	7.30-7.50

Review/Approval

Richa Shrestha
~~Shivani Goel~~ → *QC Supervisor*
Name *22Apr2022* Title *22Apr2022* Signature/Date *[Signature]* *22Apr2022*