

Product Number:	103847
Lot Number:	134211
Expiration:	8-Dec-2022
Storage:	2 - 8°C

Laboratory Analysis

Characteristic	Result	pH Target
pH	Pass	7.35-7.45

Review/Approval

Shivani Goel
Name

QC
Title

Shivani Goel 2Jan2022
Signature/Date ~~21Jan2022~~ SG
2Jan2022