

Product Number:	102002
Lot Number:	123412
Expiration:	22-Aug-2022
Storage:	2-8°C

Laboratory Analysis

Characteristic	Result	pH Target
pH Measurement	PASS	7.35-7.45

Review/Approval

Shivani Goel
Name

QC Manager
Title

Shivani Goel 28 Mar 2022
Signature/Date