

Product Number:	102002
Lot Number:	122311
Expiration:	11-Aug-2022
Storage:	2-8°C

Laboratory Analysis

Characteristic	Result	pH Target
pH Measurement	PASS	7.35-7.45

Review/Approval

Shivani Grel
Name

QC

Title

Shivani Grel 02 Feb 2022
Signature/Date