

Product Number:	102410
Lot Number:	120022
Expiration:	19-Jul-2022
Storage:	2-8°C

Laboratory Analysis

Characteristic	Result
pH Measurement	PASS

Review/Approval

Shivani Gael
Name

QC
Title

Shivani Gael 29 Jul 2021
Signature/Date