

Product Number:	101357
Lot Number:	127410
Expiration:	1-Oct-2022
Storage:	2-8°C

Laboratory Analysis

Characteristic	Result
pH Measurement	PASS

Review/Approval

Shivani Goel
Name

QC

Title

Shivani Goel 06Oct2021
Signature/Date