

Product Number:	101356
Lot Number:	127513
Expiration:	2-Oct-2022
Storage:	2-8°C

Laboratory Analysis

Characteristic	Result
pH Measurement	PASS

Review/Approval Shivani Goel QC Shivani Goel 29 NOV 2021
Name Title Signature/Date