

Product Number:	100376
Lot Number:	222007
Expiration:	8-Aug-2023
Storage:	2-8°C

**Laboratory Analysis**

Characteristic	Result	pH Target
pH Measurement	PASS	7.35 - 7.45

Review/Approval Matt Smart      QC      [Signature] 15 Aug 2022  
Name      Title      Signature/Date