

Product Number:	102252
Lot Number:	227220
Expiration:	29-Sep-2023
Storage:	2 - 8°C

Laboratory Analysis

Characteristic	Result	pH Target
pH	Pass	7.35 - 7.45

Review/Approval

Sonia Khillan QC manager

Name

Title

Soni Pill - 10/MAR/2023

Signature/Date