

Product Number:	102907
Lot Number:	503541
Expiration:	17-Mar-2023
Storage:	2 - 8°C

Component	Part Number	Lot Number	Means of Analysis	Pass / Fail
Lysate Diluent A Reagent	102908	215504	pH meter	Pass
Lysate Diluent B Reagent	102910	218102		Pass
Lysate Diluent C Reagent	102912	207611		Pass
Lysate Diluent D Reagent	102914	215603		Pass

Review/Approval

Sonia Khillan

QC

Sonia Khillan 07 JUL 2022

Name

Title

Signature/Date