

Laboratory Analysis			
Characteristic	Result	pH Target	PASS / FAIL
pH	6.19	6.18 - 6.22	PASS

Signature/Date