

Product Number:	101736 (103159)
Lot Number:	501455 (82690)
Expiration:	26-Sep-2019
Storage Condition:	2-8°C

	PASS/FAIL
Release Criteria	Pass

On behalf of:
Sivasankari Murugan

Sr. Director,
Manufacturing
Operations

Review/Approval

Bud Drummey

Name

Title

 12 Oct 2018

Signature/Date