

Product Number:	101736 (103159)
Lot Number:	502420 (01200)
Expiration:	29-Apr-2021
Storage:	2-8°C

	PASS/FAIL
Release Criteria	Pass

Review/Approval

Shivani Goel
Name

QC Manager
Title

Shivani Goel 09 Jun 2020
Signature/Date