

Product Number:	101736 (103159)
Lot Number:	502019 (92420)
Expiration:	30-Aug-2020
Storage:	2-8 °C

	PASS/FAIL
Release Criteria	Pass

Review/Approval

Patrizia Stadler

Assoc. Director, QC

P. Stadler

11 Sep 2019

Name

Title

Signature/Date