

Product Number:	101736 (103159)
Lot Number:	502117 (93010)
Expiration:	28-Oct-2020
Storage:	2-8°C

	PASS/FAIL
Release Criteria	Pass

Review/Approval

Patrizia Stadler Assoc. Director, QC
Name Title

P. Stadler

14 Nov 2019

Signature/Date