

Product Number:	101736 (103159)
Lot Number:	501924 (91690)
Expiration:	18-Jun-2020
Storage:	2-8°C

	PASS/FAIL
Release Criteria	Pass

Review/Approval

Patrizia Stadler
Name

Assoc. Director, QC
Title

P. Stadler

19 Jul 2019

Signature/Date