

Product Number:	103647
Lot Number:	107022
Expiration:	11-Jun-2021
Storage:	2-8°C

Laboratory Analysis

Characteristic	Means of Analysis	Result	Concentration
Ab Concentration	Nanodrop	PASS	1.01 mg/mL
Performance	HD-1 Analyzer	PASS	

Review/Approval

J. Park
Name

QC

Supervisor

[Signature] 29 Mar 2021
Signature/Date