

Product Number:	102010
Lot Number:	107002
Expiration:	11-Sep-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

Junpil Park

Name

QC Supervisor



29 Mar 2021

Signature/Date