

Product Number:	103647
Lot Number:	022709
Expiration:	14-Nov-2020
Storage:	2-8°C

Laboratory Analysis

Characteristic	Means of Analysis	Result	Concentration
Ab Concentration	Nanodrop	PASS	0.971 mg/mL
Performance	HD-1 Analyzer	PASS	

Review/Approval

Shivani Goel QC Manager
Name

Shivani Goel 16 Sep 2020.
Signature/Date