

<b>Product Number:</b>	103647
<b>Lot Number:</b>	012008
<b>Expiration:</b>	29-Jul-2020
<b>Storage:</b>	2-8°C

**Laboratory Analysis**

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

Review/Approval

Shivani Goel      QC Manager  
Name

*Shivani Goel*      06 May 2020  
Signature/Date