

<b>Product Number:</b>	100464
<b>Lot Number:</b>	122804
<b>Expiration:</b>	28-Jan-2024
<b>Storage:</b>	2 - 8°C

**Laboratory Analysis**

<b>Characteristic</b>	<b>Result</b>
Concentration	1.0 mg/mL

Review/Approval

Shivani Goel  
Name

QC  
Title

Shivani Goel 18 Aug 2021  
Signature/Date