

<b>Product Number:</b>	100464
<b>Lot Number:</b>	119304
<b>Expiration:</b>	28-Jul-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 03 Aug 2024

Signature/Date