

<b>Product Number:</b>	100463
<b>Lot Number:</b>	103515
<b>Expiration:</b>	13-Nov-2022
<b>Storage:</b>	2 - 8°C

**Laboratory Analysis**

<b>Characteristic</b>	<b>Result</b>
Concentration	5.2 mg/mL

Review/Approval

Junpil Park

Supervisor, QC

 04 Mar 2021

Name

Title

Signature/Date