

<b>Product Number:</b>	100463
<b>Lot Number:</b>	402922
<b>Expiration:</b>	31-Jul-2026
<b>Storage:</b>	2 - 8°C

**Laboratory Analysis**

<b>Characteristic</b>	<b>Result</b>
Concentration	5.1 mg/mL

<b>Characteristic</b>	<b>Result</b>
IgG Purity	> 95 %

Review/Approval

ROBERT SOMOGYI

Name

QC MANAGER

Title

 05 FEB 2024

Signature/Date