

Product Number:	100463
Lot Number:	414114
Expiration:	31-Jul-2026
Storage:	2 - 8°C

Laboratory Analysis


Characteristic	Result
Concentration	5.1 mg/mL

Characteristic	Result
IgG Purity	> 95 %

Review/Approval

Robert Somogyi

QC Manager

 28 MAY 2024

Name

Title

Signature/Date