

Product Number:	100464
Lot Number:	402605
Expiration:	22-Aug-2025
Storage:	2-8°C

Laboratory Analysis

Characteristic	Result
Concentration	1.1 mg/mL

Characteristic	Result
IgG Purity	> 95 %

Review/Approval

ROBERT SOMOGYI

Name

QC MANAGER

Title

*[Handwritten Signature]*

Signature/Date

30 JAN 2024