

Product Number:	100464
Lot Number:	406102
Expiration:	28-Jan-2027
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

[Signature] 05 MAR 2024

Signature/Date