

Product Number:	100464
Lot Number:	330605
Expiration:	22-Aug-2025
Storage:	2-8°C

Laboratory Analysis

Characteristic	Result
Concentration	1.1 mg/mL

Characteristic	Result
IgG Purity	>95 %

Review/Approval

Sonia Khilka Senior QC Analyst Sonia Khilka 14 Nov 2023
 Name Title Signature/Date