



CERTIFICATE OF ANALYSIS

**Product Description:** MSD GOLD SULFO-TAG Anti-Human IgG Antibody  
**Catalog Number:** D21APR-3  
**Lot Number:** D00V0044  
**Expiration Date:** 31 JUL 2028  
**Concentration:** 200X  
**Total Volume:** 200 µL  
**Storage:** 2-8°C

The above MSD GOLD™ SULFO-TAG Anti-Human IgG Antibody lot was tested according to MSD’s procedure and passed all specifications. Each lot of the MSD GOLD™ SULFO-TAG Anti-Human IgG Antibody was purified using size exclusion chromatography (SEC).

**Functional Testing Results:** This product was tested using Biotinylated Human IgG and MSD GOLD 96 well Streptavidin SECTOR Plate lot# Z0022134

**Precision, Accuracy, and Sensitivity Specifications:**

Metric	Specification	Result
Concentration Tested, Ratio of signal to reference lot signal	70% - 130%	Pass
Mean Intraplate CV	≤ 20%	Pass
Max Background Signal (counts)	≤ 250	Pass
Standard Deviation of Background Signal (counts)	≤ 20	Pass

This product is intended for Research Use Only. Not for use in Diagnostic Procedures.

	Name	Function	Signature	Date
<b>Review/Approval</b>	Karen Hamilla	Quality		27 MAR 2025