



CERTIFICATE OF ANALYSIS

Product Description: MSD GOLD SULFO-TAG Anti-Human IgG Antibody
Catalog Number: D21APR-3
Lot Number: D00V0050
Expiration Date: 30 SEP 2029
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# Z0022466

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
Review/Approval	Michelle Likiam	Quality		08 MAY 2026