



## CERTIFICATE OF ANALYSIS

Page 1 of 1

**Product Description:** MSD GOLD SULFO-TAG™ Anti-Human IgG Antibody

**Catalog Number:** D21APR-3

**Lot Number:** D00V0035

**Expiration Date:** 31 MAR 2027

**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# Z0022078.

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

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

	Name	Function	Signature	Date
Review/Approval	Karen Hamilla	Quality		08 NOV 2023