



CERTIFICATE OF ANALYSIS

Page 1 of 1

Product Description: MSD GOLD Read Buffer A
Catalog Number: R92TG-Series
Lot Number: Y0600046
Expiration Date: 31 OCT 2026

Statement: MSD GOLD Read Buffer A is verified to ensure that consistent signals are obtained across lots. Details on the quality control testing of MSD GOLD Read Buffer A are provided in the product insert available at www.mesoscale.com.

MSD GOLD Read Buffer A was tested using SULFO-TAG™ Conjugated Biotinylated IgG (BTI) lot # R0010830 and MSD GOLD™ 96-well Streptavidin SECTOR Plate lot # Z0022134

Concentration of BTI	Mean ECL Spec and Tolerance Range	Intraplate CV or SD spec	Pass/Fail
2.00 nM BTI (0.10 pmol)	17024 - 23032 counts	10 %	PASS
0 nM BTI (0 pmol)	48 - 68 counts	15 counts	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		05 JAN 2024