



CERTIFICATE OF ANALYSIS

Product Description: MSD GOLD Read Buffer A

CatalogNumber/Size: R92TG-Series

Lot Number: Y0600022

Expiration Date: 30 APR 2024

Storage Temperature: Room Temperature

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 # R0010698 and MSD GOLD™ 96-well Streptavidin SECTOR Plate lot # Z0021486

Concentration of BTI	Mean ECL spec and tolerance range	Intraplate CV or SD spec	Pass/Fail?
2.00 nM BTI (0.10 pmol)	21701 - 29360 counts	10 %	PASS
0 nM BTI (0 pmol)	48 - 68 counts	15 counts	PASS

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

	Name	Function	Signature	Date
Review/Approval	Karen Hamilla	Quality		28 MAY 2021