



CERTIFICATE OF ANALYSIS

Product Description: MSD GOLD Read Buffer
CatalogNumber/Size: R92TG-Series
Lot Number: Y0600013
Expiration Date: 31 JAN 2023
Storage Temperature: Room Temperature

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

MSD GOLD Read Buffer was tested using SULFO-TAG™ Conjugated Biotinylated IgG (BTI) lot # R0010605 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
0.05 nM BTI (0.0025 pmol)	474 - 642 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		16 JAN 2020