



CERTIFICATE OF ANALYSIS

Product Description: MSD GOLD™ Read Buffer

Catalog Number/Size: R92TG-2

Lot Number: Y0600008

Expiration Date: 31 JAN 2022

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 # Z0021294.

Concentration of BTI	Mean ECL spec and tolerance range	Intraplate CV or SD spec	Pass/Fail?
2.0 nM (0.10 pmol)	25,531 counts \pm 15%	10 % CV	PASS
0.05 nM (0.0025 pmol)	525 counts \pm 15%	10 % CV	PASS
0 nM (0 pmol)	58 \pm 10 counts	15 count SD	PASS

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

	Name	Function	Signature	Date
Review/Approval	Carissa Hubbard	Quality		08 OCT 2019