



CERTIFICATE OF ANALYSIS

Product Description: MSD GOLD™ Read Buffer

Catalog Number/Size: R92TG-Series

Lot Number: Y0600005

Expiration Date: 31 AUG 2021

Storage Temperature: Room Temp

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 # R0010578 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)	558 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
Preparer	Carissa Hubbard	Quality Control		29 AUG 2018
Review/Approval	Dawn Stollar	Quality Assurance		29 AUG 2018