



## CERTIFICATE OF ANALYSIS

**Product Description:** MSD GOLD™ Read Buffer

**Catalog Number/Size:** R92TG-Series

**Lot Number:** Y0600002

**Expiration Date:** 31 MAR 2020

**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](http://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 # Z0021179.

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 ± 15%	10 % CV	PASS
0.05 nM (0.0025 pmol)	525 counts ± 15%	10 % CV	PASS
0 nM (0 pmol)	58 ± 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
<b>Preparer</b>	Carissa Hubbard	Quality Control		13 APR 2017
<b>Review/Approval</b>	Dawn Stollar	Quality Assurance		13 APR 2017