



CERTIFICATE OF ANALYSIS

**Product Description:** MSD GOLD Read Buffer A  
**CatalogNumber/Size:** R92TG-Series  
**Lot Number:** Y0600016  
**Expiration Date:** 31 Jul 2023  
**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](http://www.mesoscale.com).

MSD GOLD Read Buffer A 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		06 AUG 2020