



## CERTIFICATE OF ANALYSIS

Page 1 of 1

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

**Catalog Number:** R60AM-Series

**Lot Number:** Y0800054


**Expiration Date:** 31 AUG 2027

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

MSD GOLD Read Buffer B was tested using SULFO-TAG™ conjugated biotinylated IgG (BTI) lot # R0010868 and MSD GOLD™ 96-well Streptavidin SECTOR Plate lot # Z0022134 .

Concentration of BTI	Specification Metric	Allowable Range	Pass/Fail?
2.00 nM BTI (0.10 pmol)	Test : Reference Signal Ratio	90 - 110 %	PASS
	Coefficient of Variation	≤ 10 %	PASS
0 nM BTI (0 pmol)	Test - Reference Signal Difference	± 10 counts	PASS
	Standard Deviation	≤ 15 counts	PASS

The above product is intended for Research Use Only. Not for use in Diagnostic Procedures.

	Name	Function	Signature	Date
Review/Approval	Karen Hamilla	Quality		10 APR 2025