


**CERTIFICATE OF ANALYSIS**

Product Description: MSD GOLD Read Buffer A
Catalog Number: R92TG-Series
Lot Number: Y0600064
Expiration Date 31 JUL 2028

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.

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	$\leq 10 \%$	Pass
0 nM BTI (0 pmol)	Test - Reference Signal Difference	± 10 counts	Pass
	Standard Deviation	≤ 15 counts	Pass

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

	Name	Function	Signature	Date
Review/Approval	Karen Hamilla	Quality		18 SEP 2025