


**CERTIFICATE OF ANALYSIS**

**Product Description:** MSD GOLD Read Buffer A  
**Catalog Number:** R92TG-Series  
**Lot Number:** Y0600069  
**Expiration Date** 30 NOV 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](http://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	$\pm 10$ counts	Pass
	Standard Deviation	$\leq 15$ 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		08 DEC 2025