


**CERTIFICATE OF ANALYSIS**

**Product Description:** MSD GOLD Read Buffer B  
**Catalog Number:** R60AM-Series  
**Lot Number:** Y0800061  
**Expiration Date** 30 JUN 2028

**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).

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		19 JAN 2026