



CERTIFICATE OF ANALYSIS

Product Description: MSD GOLD Read Buffer B
Catalog Number: R60AM-Series
Lot Number: Y0800062
Expiration Date 31 AUG 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.

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

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

	Name	Function	Signature	Date
Review/Approval	Karen Hamilla	Quality		02 APR 2026