



CERTIFICATE OF ANALYSIS

Page 1 of 1

Product Description: MSD GOLD™ Read Buffer B

Catalog Number: R60AM-Series

Lot Number: Y0800025

Expiration Date: 30 JUN 2024

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.

MSD GOLD™ Read Buffer B was tested using SULFO-TAG™ Conjugated Biotinylated IgG (BTI) lot # R0010698 and MSD GOLD™ 96-well Streptavidin SECTOR Plate Z0021725.

Concentration of BTI	Mean ECL Spec and Tolerance Range	Intraplate CV or SD spec	Pass/Fail
2.00 nM BTI (0.10 pmol)	21421 - 28982 counts	10 %	PASS
0 nM BTI (0 pmol)	24 - 44 counts	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		04 MAR 2022