



CERTIFICATE OF ANALYSIS

Page 1 of 1

Product Description: MSD® Read Buffer B

Catalog Number: R92BA-Series

Lot Number: Y0530003

Expiration Date: 31 DEC 2023

Statement: MSD Read Buffer B is verified to ensure that consistent signals are obtained across lots. Details on the quality control testing of Select Product Description are provided in the product insert available at www.mesoscale.com.

MSD Read Buffer B was tested using SULFO-TAG™ Conjugated Biotinylated IgG (BTI) lot # R0010698 and MSD GOLD™ 96-well Streptavidin SECTOR Plate Z0021486.

Concentration of BTI	Mean ECL Spec and Tolerance Range	Intraplate CV or SD spec	Pass/Fail
2.00 nM BTI	19354 – 26185	10%	Pass
Dil-100	16 - 56	15	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		28 JUL 2021

COA-14810-A

QA-FM-313-A