

# MSD GOLD™ Read Buffer B



## Ordering Information

MSD Customer Service  
Phone: 1-240-314-2795  
Fax: 1-301-990-2776  
Email: [CustomerService@mesoscale.com](mailto:CustomerService@mesoscale.com)  
[www.mesoscale.com/support](http://www.mesoscale.com/support)

## Scientific Support

Phone: 1-240-314-2798  
Email: [ScientificSupport@mesoscale.com](mailto:ScientificSupport@mesoscale.com)  
[www.mesoscale.com/support](http://www.mesoscale.com/support)

## Company Address

MESO SCALE DISCOVERY®  
A division of  
Meso Scale Diagnostics, LLC.  
1601 Research Boulevard  
Rockville, MD 20850-3173  
USA

[www.mesoscale.com](http://www.mesoscale.com)®

For Research Use Only.  
Not for use in  
diagnostic procedures.

<b>Contents</b>	MSD GOLD Read Buffer B (18 mL) MSD GOLD Read Buffer B (90 mL) MSD GOLD Read Buffer B (200 mL) MSD GOLD Read Buffer B (1000 mL)	Cat. No. R60AM-1 Cat. No. R60AM-2 Cat. No. R60AM-3 Cat. No. R60AM-4
<b>Storage</b>	Temperature range (15–30 °C)	

## Summary and Intended Use

Read buffer is an essential component of MSD® assays: it is required for the generation of electrochemiluminescence (ECL) signals. MSD GOLD Read Buffer B was optimized for use in all MSD assays and can be used as a replacement for MSD GOLD Read Buffer A and MSD Read Buffer T (4X). MSD GOLD Read Buffer B is also recommended for assays that measure whole-cells, vesicles, or other membrane-containing analytes.

**Note:** MSD GOLD Read Buffer B is not recommended for MSD assays developed using high-bind surface MSD plates.

MSD GOLD Read Buffer B has stringent quality control release criteria that ensure lot-to-lot reproducibility. The specifications are designed to meet the requirements of critical assays and longitudinal studies.

**Note:** MSD GOLD Read Buffer B is provided at the working concentration. We recommend using MSD GOLD Read Buffer B at the supplied concentration without any dilution.

## Reagent

MSD GOLD Read Buffer B is a Tris-based buffer containing a low volatile amine coreactant for light generation in ECL assays.

## Storage

To maximize consistency in ECL signals across assay runs, MSD GOLD Read Buffer B should be stored at the recommended temperature, away from direct sunlight, and handled per the instructions provided in this document. MSD GOLD Read Buffer B is stable for 30 months at the suggested temperature. Please refer to the certificate of analysis (COA) for more detail.

## Typical Protocol

After the final incubation step of the assay, wash the plate, add 150 µL/well of MSD GOLD Read Buffer B to each well, and then analyze the plate in an MSD® instrument. Unless necessary for optimal performance of your assay, incubation in MSD GOLD Read Buffer B is not required before reading the plate.

For the lot-specific specifications of MSD GOLD Read Buffer B, refer to the COA supplied with the product or available at [www.mesoscale.com](http://www.mesoscale.com).

## Safety

Use safe laboratory practices and wear gloves, safety glasses, and lab coats when handling MSD GOLD Read Buffer B. Handle and dispose of all hazardous samples properly in accordance with local, state, and federal guidelines. Additional product-specific safety information is available in the applicable safety data sheet(s) (SDS), which can be obtained from MSD Customer Service or at [www.mesoscale.com](http://www.mesoscale.com).

# MSD GOLD Read Buffer B

## MSD GOLD Read Buffer B Specifications

MSD GOLD Read Buffer B has stringent specifications for accuracy and precision. The quality control process is designed to ensure the reproducibility of ECL signals across experiments.

MSD GOLD Read Buffer B lots are tested on MSD GOLD 96-well Streptavidin plates at two levels of biotinylated IgG (BTI) labeled with SULFO-TAG™ BTI. The mean signal and %CV are calculated for each BTI concentration. The release specifications verify that the interlot signal performance at the high (2.0 nM BTI) and low signal (0 nM BTI) level is within the established tolerance range. Typical variability is below 10% at the high signal level and <7 counts at the low signal level. Quality control specifications for MSD GOLD Read Buffer B are shown below.

Concentration of BTI for 96-well Streptavidin SECTOR™ Plate	Signal Specification and Tolerance Range	Intra-plate CV or Standard Deviation (SD)
2.0 nM (0.1 pmole of IgG)	21421 - 28982 counts	CV ≤ 10%
0 nM ( 0 pmole of IgG)	24 - 44 counts	SD ≤ 15 counts

For the specifications of the actual lot, refer to the COA supplied with the read buffer.

## MSD GOLD Read Buffer B Reproducibility

The reproducibility of MSD GOLD Read Buffer B was tested on MSD GOLD Streptavidin plates. The data sets below demonstrate the signal reproducibility of six independent lots (color-coded data points) of MSD GOLD Read Buffer B when tested on a single plate lot across multiple test days. Each data point is the mean signal measured across one plate and the error bars represent the standard deviation across all the replicates (n = 48 replicates). The dotted lines represent the specification around the mean signal measured across the entire data set. The data demonstrate the uniform performance that is essential for critical assays and longitudinal studies.

