

MSD GOLD™ Read Buffer A



Ordering Information

MSD Customer Service
Phone: 1-240-314-2795
Fax: 1-301-990-2776
Email: CustomerService@mesoscale.com
www.mesoscale.com/support

Scientific Support

Phone: 1-240-314-2798
Email: ScientificSupport@mesoscale.com
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®

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

Contents	MSD GOLD Read Buffer A (200 mL)	Cat. No. R92TG-1
	MSD GOLD Read Buffer A (1,000 mL)	Cat. No. R92TG-2
	MSD GOLD Read Buffer A (18 mL)	Cat. No. R92TG-3
	MSD GOLD Read Buffer A (90 mL)	Cat. No. R92TG-4
Storage	Room temperature (15-30°C)	

Summary and Intended Use

MSD GOLD products provide a rapid and convenient method for the development of new immunoassays and the transfer of existing ELISAs utilizing the benefits of MSD's highly sensitive electrochemiluminescence (ECL) technology. Read buffer is an essential component of MSD® immunoassays and is required for the generation of ECL signals. MSD GOLD Read Buffer A has stringent quality control release criteria that ensure lot-to-lot reproducibility, consistent with other products in the MSD GOLD product line. The specifications are designed to meet the requirements of critical assays and longitudinal studies.

Note: MSD GOLD Read Buffer A is provided at the working concentration for most assays and we recommend using MSD GOLD Read Buffer A at the supplied concentration without any dilution. At this concentration, MSD GOLD Read Buffer A can be used as a replacement for 2X MSD Read Buffer T.

Reagent

MSD GOLD Read Buffer A is a tris-based buffer containing tripropylamine (TPA) as a co-reactant for light generation in ECL immunoassays.

Storage

In order to maximize consistency in ECL signals across assay runs, MSD GOLD Read Buffer A must be stored at the recommended temperature and handled per the instructions provided in this document.

Typical Protocol

After the final incubation step of the assay, wash the plate, add 150 µL/well of MSD GOLD Read Buffer A 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 A is not required before reading the plate.

For the lot-specific specifications of MSD GOLD Read Buffer A, refer to the certificate of analysis (COA) supplied with the product or available at www.mesoscale.com.

Safety

Use safe laboratory practices and wear gloves, safety glasses, and lab coats when handling MSD GOLD Read Buffer A. 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 safety data sheet (SDS), which can be obtained from MSD Customer Service or at www.mesoscale.com.

MSD GOLD Read Buffer A

MSD GOLD Read Buffer A Specifications

MSD GOLD designation certifies that the products meet stringent specifications for accuracy and precision. The quality control process is designed to ensure reproducibility of ECL signals across experiments.

MSD GOLD Read Buffer A lots are tested on MSD GOLD 96-well Streptavidin plates, with three levels of biotinylated IgG conjugated with SULFO-TAG™ label (BTI) at 0, 0.05, and 2.0 nM concentrations. The mean signal and %CV is calculated for each BTI concentration. The release specifications verify that the inter-lot signal performance at each level is within +/-15% of the established target, with typical variability below 10%. Quality control specifications for MSD GOLD Read Buffer A are shown below.

Concentration of BTI for 96-well Streptavidin SECTOR® Plate	Mean ECL Signal Specification and Tolerance Range	Intra-plate CV or Standard Deviation (SD)
2.0 nM	*25,500 +/- 15%	≤10%
0.05 nM	*520 +/- 15%	≤10%
0 nM	*58 +/- 10 counts	SD≤15 counts

*These signal specifications represent a specific lot of BTI reagent and may vary for different lots. For the actual specifications, refer to the COA supplied with the read buffer.

MSD GOLD Read Buffer A Reproducibility

The reproducibility of MSD GOLD Read Buffer A was tested on MSD GOLD Streptavidin plates. The data sets below demonstrate the signal reproducibility of three independent lots of MSD GOLD Read Buffer A when tested on a single plate lot across multiple test days. Each data point is the mean signal measured across three plates and the error bars represent the standard deviation across all the replicates (n=48 replicates). The dotted lines represent a +/-15% window around the mean signal measured across the entire data set. The data demonstrates the uniform performance that is essential for critical assays and longitudinal studies.

