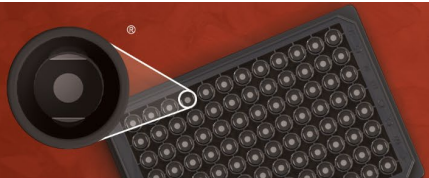


R-PLEX[®]

Human CD47



MSD Toxicology Assays

www.mesoscale.com®

Ordering Information

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

in rat serum and plasma

Scientific Support

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

Company Address

Meso Scale Discovery
A division of
Meso Scale Diagnostics, LLC.
1601 Research Boulevard
Rockville, MD 20850-3173 USA

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

MESO SCALE DISCOVERY, MESO SCALE DIAGNOSTICS, www.mesoscale.com, MSD, MSD (design), MSD GOLD, MESO, QuickPlex, QuickPlex Ultra, SECTOR, SULFO-TAG, R-PLEX, R-PLEX (design), U-PLEX, 96 WELL SMALL-SPOT (design), and Spot the Difference are trademarks and/or service marks of Meso Scale Diagnostics, LLC.
©2016-2025 Meso Scale Diagnostics, LLC. All rights reserved.

MK-DS-1191-v3-2025Jan

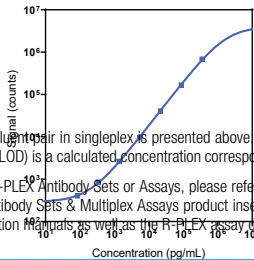


MSD Toxicology Assays

Product Options	Catalog Number	Description
Singleplex	K151ALZR-2	R-PLEX Human CD47 Assay with SECTOR™ plates
Antibody Set	F21ALZ-3	R-PLEX Human CD47 Antibody Set
Protocol	Assay Protocol 1 in the R-PLEX Antibody Sets Singleplex Assays Insert and the SECTOR and QuickPlex	

MESO SCALE DISCOVERY® R-PLEX Antibody Sets and Assays provide an expanded menu of electrochemiluminescence reagents for biomarker discovery and development. The Antibody Sets include a matched, biotinylated capture and SULFO-TAG™ conjugated detection antibody pair and Calibrator for the quick and easy development of highly sensitive immunoassays. R-PLEX® Assays add diluents, MSD GOLD™ Read Buffer B, and the appropriate plates, providing a complete set of components to develop and run an assay. R-PLEX Assays with QuickPlex Ultra plates can only be read on the MESO® QuickPlex Q 60MM instrument. R-PLEX Assays with SECTOR plates can be read on all MESO SECTOR® instruments and the MESO QuickPlex® SQ 120/120MM instrument.

Representative Calibration Curve and Sensitivity



Calibrator Stock (20X)	7,000,000 pg/mL
Calibrator Standard 1 (top of curve); 20-fold Calibrator stock dilution	350,000 pg/mL
LLOD	69 pg/mL

Representative data using the primary diluent pair in singleplex is presented above. The calibration curve was fitted with a 4-parameter logistic (or sigmoidal dose-response) model with a 1/Y² weighting. The lower limit of detection (LLOD) is a calculated concentration corresponding to the signal 2.5 standard deviations above the background (zero Calibrator).

For developing individual assays using R-PLEX Antibody Sets or Assays, please refer to the R-PLEX Antibody Sets & Singleplex Assays product insert. For multiplex assay development using R-PLEX Antibody Sets, use the R-PLEX Antibody Sets & Multiplex Assays product insert. For multiplexing a combination of R-PLEX and U-PLEX® Antibody Sets, refer to the U-PLEX Development Pack product insert. The detailed instruction Manuals as well as the R-PLEX assay datasheets are available at <https://www.mesoscale.com>.

Source

Calibrator	Recombinant Human CD47 expressed in a mouse cell line
Antibodies	Capture antibody - mouse monoclonal, detection antibody - sheep polyclonal
Assay Generation	A

Recommended Assay Components

The following items are recommended for developing assays using R-PLEX Antibody Sets.

Plates for Singleplex Assay Development

Plates are available in packs of 1, 5, 30, 120, and 510.

Description	Catalog No.	Included in:
MSD GOLD 96-well Small Spot Streptavidin SECTOR Plates	L45SA	K151ALZR-2
96-well Small Spot Streptavidin QuickPlex Ultra Plates	L4BLA	K151ALZR-22

Buffers

Description	Catalog No.	Included in:
-------------	-------------	--------------

MESO SCALE DISCOVERY, MESO SCALE DIAGNOSTICS, www.mesoscale.com, MSD, MSD (design), MSD GOLD, MESO, QuickPlex, QuickPlex Ultra, SECTOR, SULFO-TAG, R-PLEX, R-PLEX (design), U-PLEX, 96 WELL SMALL-SPOT (design), and Spot the Difference are trademarks and/or service marks of Meso Scale Diagnostics, LLC.
©2016-2025 Meso Scale Diagnostics, LLC. All rights reserved.

MK-DS-1191 -v3-2025Jan



Deleted: _____

MSD Toxicology Assays

Diluent 100 (50 mL)	R50AA-4	K151ALZR
MSD GOLD Read Buffer B (90 mL)	R60AM-2	K151ALZR
MSD Wash Buffer (20X, 100 mL)	R61AA-1	-
MSD Tris Wash Buffer (10X, 200 mL)	R61TX-2	-

in rat serum and plasma

Diluent Compatibility

The R-PLEX Human CD47 Antibody Set has been tested with the following diluent pair(s). The primary diluent pair was used for generating the data shown and is included in the R-PLEX Human CD47 Assay. Alternate diluent pairs may be considered for use with R-PLEX Antibody Sets, based on the sample type.

Tested Diluent Pair	Assay Diluent		Antibody Diluent	
	Description	Catalog No.	Description	Catalog No.
Primary	Diluent 57	R50BZ	Diluent 3	R50AP
Alternate 1	Diluent 12	R50JA	Diluent 11	R55BA

Diluents sufficient to run 40 samples per plate in duplicate with up to 10-fold sample dilution are provided with R-PLEX assays.

Additional assay diluent may be necessary for samples that are diluted greater than 10-fold and can be ordered separately. Diluent 100 may be used in place of assay diluent for samples that require high dilution. Testing of different diluents can help optimize assays for specific experimental conditions and sample types.

MSD offers a range of assay and antibody diluents for separate purchase. Please view our website at www.mesoscale.com/R-PLEX for more details on compatible diluents for R-PLEX assays.

