



www.mesoscale.com[®]

Ordering Information

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

Scientific Support

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

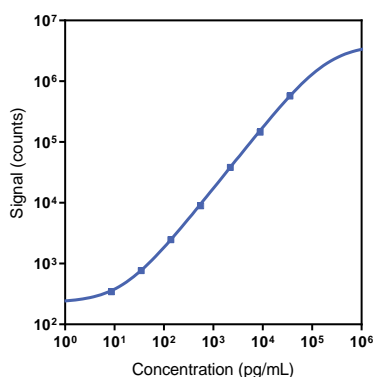
Company Address

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

Product Options	Catalog Number	Description
Singleplex	K151AMKR-2	R-PLEX Human TYRO-3 Assay with SECTOR™ plates
	K151AMKR-22	R-PLEX Human TYRO-3 Assay with QuickPlex [®] plates
Antibody Set	F21AMK-3	R-PLEX Human TYRO-3 Antibody Set
Protocol	R-PLEX Product Inserts are available at https://www.mesoscale.com	

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 either MSD GOLD small spot streptavidin QuickPlex or SECTOR plates, providing a complete set of components to develop and run an assay. R-PLEX Assays with QuickPlex plates can only be read on MESO[®] QuickPlex Q 60MM instruments. R-PLEX Assays with SECTOR plates can be read on all MESO SECTOR[®] and MESO QuickPlex[®] SQ 120/120MM instruments.

Representative Calibration Curve and Sensitivity



Calibrator Stock (20X)	700,000 pg/mL
Calibrator Standard 1 (top of curve); 20-fold Calibrator stock dilution	35,000 pg/mL
LLOD	6.1 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>.

Sample Dilution

The R-PLEX Human TYRO-3 Antibody Set has been tested with the following sample type(s) at the specified dilution in assay diluent.

Sample Type	Recommended Fold Dilution
Serum	4
EDTA Plasma	4

The fold dilution specified is a suggested dilution. We recommend that sample dilution and other assay parameters be optimized before running samples.

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

MSD® R-PLEX Assays

Source

Calibrator	Recombinant Human TYRO-3 expressed in a mouse cell line
Antibodies	Capture antibody - mouse monoclonal, detection antibody - goat 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	K151AMKR-2
MSD GOLD 96-well Small Spot Streptavidin QuickPlex Plates	L4BSA	K151AMKR-22

Buffers

Description	Catalog No.	Included in:
Diluent 100 (50 mL)	R50AA-4	K151AMKR
MSD GOLD Read Buffer B (90 mL)	R60AM-2	K151AMKR
MSD Wash Buffer (20X, 100 mL)	R61AA-1	-
MSD Tris Wash Buffer (10X, 200 mL)	R61TX-2	-

Diluent Compatibility

The R-PLEX Human TYRO-3 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 TYRO-3 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 12	R50JA	Diluent 11	R55BA
Alternate 1	Diluent 13	R56BB	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.

