



MMAF Antibody Drug Conjugate (ADC) ELISA Assay Kit

Catalog Number:

KTR-758 (1 x 96 wells)

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

v. 1.0

EAGLE BIOSCIENCES, INC.
20A Northwest Blvd., Suite 112, Nashua, NH 03063
Phone: 617-419-2019 Fax: 617-419-1110
WWW.EAGLEBIO.COM



INTENDED USE

MMAF Antibody Drug Conjugate (ADC) ELISA Assay kit is intended for use in the quantitative determination of antibody-MMAF-conjugate level in test sample. It is useful for pre-clinical and clinical pharmacology study of MMAF Antibody Drug Conjugate (ADC).

- Samples from tissue/cell culture and serum samples from human, rat, mouse, primate, etc. can be used directly with this MMAF Antibody Drug Conjugate (ADC) ELISA Assay kit.
- Both humanized monoclonal antibody based MMAF-ADC and mouse monoclonal antibody based MMAF-ADC can be measured with this MMAF Antibody Drug Conjugate (ADC) ELISA Assay kit.

PRINCIPLE OF THE ASSAY

This MMAF Antibody Drug Conjugate (ADC) ELISA Assay kit is designed, developed and produced for the quantitative measurement of antibody MMAF conjugate in serum. The assay utilizes the competitive immunoassay technique with an antibody that exclusively binds to MMAF.

Assay calibrators (antibody MMAF conjugate) and test serum samples are added directly to wells of a microtiter plate that is coated with specific anti-MMAF antibody. Subsequently, a horseradish peroxidase (HRP) conjugated MMAF is added to each well. During the incubation period, the antibody MMAF conjugate competes with the HRP conjugated MMAF for the limited binding sites of anti-MMAF antibody. An immune complex of well coated "anti-MMAF antibody – HRP conjugated MMAF" is formed. The unbound antibodies and buffer matrix are removed in the subsequent washing step. For the detection of this immunocomplex, the well is then incubated with a substrate solution in a timed reaction, which is terminated with an acidic reagent (i.e. ELISA stop solution). The absorbance is then measured in a spectrophotometric microplate reader. The enzymatic activity of the immunocomplex bound to the wall of each microtiter well is inversely proportional to the amount of antibody-MMAF conjugate in the test sample. A calibration curve is generated by plotting the absorbance versus the respective antibody-MMAF conjugate concentration for each calibrator on a 4-parameter or log-logit curve fitting. The concentration of antibody-MMAF conjugate in test samples is determined directly from this calibration curve.

REAGENTS: Preparation and Storage

The MMAF Antibody Drug Conjugate (ADC) ELISA Assay kit must be stored at 2 – 8°C upon receipt. For the expiration date of the kit refer to the label on the kit box. All components of the MMAF Antibody Drug Conjugate (ADC) ELISA Assay kit are stable until this expiration date.

Prior to use allow all reagents to come to room temperature. Regents from different kit lot numbers should not be combined or interchanged.

1. Anti-MMAF Antibody Coated Microplate (Cat. No. 30749)

One microplate with twelve by eight strips (96 wells total) coated with specific anti-MMAF antibody. The plate is framed and sealed in a foil zipper bag with a desiccant. This reagent should be stored at 2 – 8 °C and is stable until the expiration date on the kit box.



2. HRP Conjugated MMAF (Cat. No. 30719)

One vial containing 3 mL of ready to use HRP labeled MMAF in a stabilized protein matrix. This reagent should be stored at 2 – 8°C and is stable until the expiration date on the MMAF Antibody Drug Conjugate (ADC) ELISA Assay kit box.

3. ELISA Wash Concentrate (Cat. No. 10010)

One bottle containing 30 mL of 30-fold concentrate. Before use the contents must be diluted with 870 mL of demineralized water and mixed well. Upon dilution, this yields a working wash solution containing a surfactant in phosphate buffered saline with a non-azide, non-mercury preservative. The diluted wash solution may be stored at room temperature and is stable until the expiration date on the MMAF Antibody Drug Conjugate (ADC) ELISA Assay kit box.

4. ELISA HRP Substrate (Cat. No. 10020)

One bottle containing 15 mL of tetramethylbenzidine (TMB) with hydrogen peroxide. This reagent should be stored at 2 – 8°C and is stable until the expiration date on the MMAF Antibody Drug Conjugate (ADC) ELISA Assay kit box.

5. ELISA Stop Solution (Cat. No. 10030)

One bottle containing 15 mL of stop solution. This reagent may be stored at 2 – 8°C or room temperature and is stable until the expiration date on the MMAF Antibody Drug Conjugate (ADC) ELISA Assay kit box.

6. Antibody-MMAF Conjugated Zero Calibrator (Cat. No. 30721)

One vial containing 30mL zero calibrator (30721). This reagent is used for diluting the calibration stock to make assay calibrators, as well as for diluting test samples. This reagent should be stored at 2-8°C and is stable until the expiration date on the MMAF Antibody Drug Conjugate (ADC) ELISA Assay kit box.

7. Antibody-MMAF Conjugated Calibration Stock (Cat. No. 30722) –

Not provided in the kit (optional)

One vial (30722) containing the calibration stock of antibody-MMAF-conjugate in a lyophilized (0.5 mL) serum based matrix with a non-azide preservative. **Refer to the vial for exact concentration of the standard.** This reagent should be stored at 2 – 8°C and is stable until the expiration date on the kit box.

SAFETY PRECAUTIONS

The reagents must be used in professional laboratory. Source material for reagents containing bovine serum was derived in the contiguous 48 United States. It was obtained only from healthy donor animals maintained under veterinary supervision and found free of contagious diseases. Wear gloves while performing this assay and handle these reagents as if they are potentially infectious. Avoid contact with reagents containing TMB, hydrogen peroxide, or sulfuric acid. TMB may cause irritation to skin and mucous membranes and cause an allergic skin reaction. TMB is a suspected carcinogen. Sulfuric acid may cause severe irritation on contact with skin. Do not get in eyes, on skin, or on clothing. Do not ingest or inhale fumes. On contact, flush with copious amounts of water for at least 15 minutes. Use Good Laboratory Practices.



MATERIALS REQUIRED BUT NOT PROVIDED

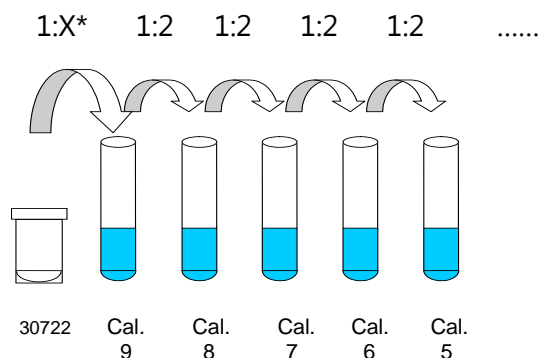
1. **Antibody-MMAF Conjugated Stock (Cat No.30722)**
2. Precision single channel pipettes capable of delivering 25 μ L, 50 μ L, 100 μ L, etc.
3. Disposable pipette tips suitable for above volume dispensing.
4. Aluminum foil.
5. Deionized or distilled water.
6. Plastic microtiter well cover or polyethylene film.
7. ELISA multichannel wash bottle or automatic (semi-automatic) washing system.
8. Spectrophotometric microplate reader capable of reading absorbance at 450/650 or 450/620 nm.

REAGENT PREPARATION

1. Reagent Preparation

- (1) Prior to use allow all reagents of the MMAF Antibody Drug Conjugate (ADC) ELISA Assay kit to come to room temperature. Reagents from different kit lot numbers should not be combined or interchanged.
- (2) ELISA Wash Concentrate must be diluted to working solution prior to use. Please see REAGENTS section for details.
- (3) Using EDI Calibrators:

Reconstitute calibration stock 30722 with **0.5 mL** DI-water. Dilute the reconstituted calibration stock (30722) 1:X* using the zero calibrator (30721) to obtain a level nine calibrator at 2 μ g/mL. Further create calibrator level eight to two by 1:2 serial dilutions to obtain these calibrators with concentrations of 1 μ g/mL, 0.5 μ g/mL, 0.25 μ g/mL, 0.125 μ g/mL, 0.063 μ g/mL, 0.032 μ g/mL and 0.016 μ g/mL. Assay calibrators should be used within 2 hours and should be stored below -20°C. Do not exceed 3 freeze-thaw cycles.



X^* = the concentration of 30722 / 2



(4) Test Configuration

ROW	STRIP 1	STRIP 2	STRIP 3	STRIP 4
A	STD 1	STD 5	STD 9	SAMPLE 4
B	STD 1	STD 5	STD 9	SAMPLE 4
C	STD 2	STD 6	SAMPLE 1	SAMPLE 5
D	STD 2	STD 6	SAMPLE 1	SAMPLE 5
E	STD 3	STD 7	SAMPLE 2	
F	STD 3	STD 7	SAMPLE 2	
G	STD 4	STD 8	SAMPLE 3	
H	STD 4	STD 8	SAMPLE 3	

NOTE:

The validation data of this test was generated by using EDI Antibody-MMAF Conjugated Stock (Cat. No. 30722). To order this calibrator stock, please order Ab-MMAF Conjugated Stock (Cat. No. 30722).

- (5) Place a sufficient number of Anti-MMAF antibody coated microwell strips in a holder to determine calibrators and unknown samples in duplicates.

ASSAY PROCEDURE

- (1) Add **100 µL** of calibrators and test samples into the designated microwells.
- (2) Seal the plate wells securely, cover with foil or other material to protect from light, and rotate on an ELISA plate shaker (small orbit radius) for **30 minutes** at 400 to 450 rpm.
- (3) Immediately add **25 µL** of HRP Conjugated MMAF (cat# 30719) to each well. (*Note: no wash step before add the HRP conjugated MMAF*)
- (4) Seal the plate wells securely, cover with foil or other material to protect from light, and rotate on an ELISA plate shaker (small orbit radius) for **2 hr. ± 10 minutes** at 400 to 450 rpm.
- (5) Wash each well 5 times by dispensing 350 µL of working wash solution into each well and then completely aspirating the contents. Alternatively, an automated microplate washer can be used.
- (6) Add **100 µL** of ELISA HRP Substrate into each of the wells.
- (7) Cover the plate with aluminum foil or other material to avoid exposure to light. Incubate plate static, at room temperature for **20 minutes**.
- (8) Immediately add **100 µL** of ELISA Stop Solution into each of the wells. Mix gently.
- (9) Read the absorbance at 450 nm.

PROCEDURAL NOTES

1. It is recommended that all calibrators and unknown samples be assayed in duplicate in the MMAF Antibody Drug Conjugate (ADC) ELISA Assay kit. The average absorbance reading of each duplicate should be used for data reduction and the calculation of results. It is recommended to add external controls to each assay.
2. Keep light sensitive reagents in the original amber bottles.
3. Store any unused antibody coated strips in the foil zipper bag with desiccant to protect from moisture.
4. Careful technique and use of properly calibrated pipetting devices are necessary to ensure reproducibility of the MMAF Antibody Drug Conjugate (ADC) ELISA Assay test.

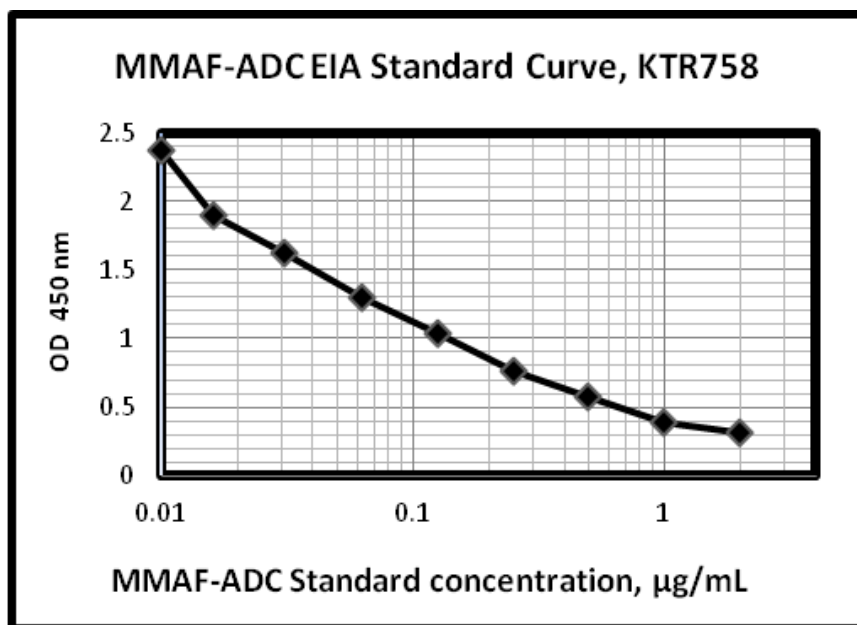
5. Incubation times or temperatures other than those stated in this insert may affect the results.
6. An orbital mixer with a larger orbit radius (e.g. > 1 cm) may be used at speeds of 150 to 200 rpm.
7. Avoid air bubbles in the microwell as this could result in lower binding efficiency and higher CV% of duplicate reading.
8. All reagents of the MMAF Antibody Drug Conjugate (ADC) ELISA Assay kit should be mixed gently and thoroughly prior to use. Avoid foaming.
9. If adapting this assay to automated ELISA system such as DS-2, DSX or Trituras, a procedural validation is necessary if there is any modification of the assay procedure.

INTERPRETATION OF RESULTS

It is recommended to use a 4-parameter or log-logit calibration curve fitting.

1. Calculate the average absorbance for each pair of duplicate test results.
2. The calibration curve is generated by the corrected absorbance of all calibrator levels on the ordinate against the calibrator concentration. Appropriate computer assisted data reduction programs should be used for the calculation of results.

The antibody-MMAF conjugate concentrations for the test samples are read directly from the calibration curve using their respective corrected absorbance.





EXAMPLE DATA AND STANDARD CURVE

A typical absorbance data and the resulting calibration curve from this MMAF Antibody Drug Conjugate (ADC) ELISA Assay kit are represented. This curve should not be used in lieu of calibration curve generated with each assay.

Well I.D.	OD 450 nm Absorbance		B/B ₀
	Readings	Average	
Cal-1: 0.000 µg/mL	2.362 2.373	2.368	100.0%
Cal-2: 0.016 µg/mL	1.899 1.899	1.899	80.2%
Cal-3: 0.032 µg/mL	1.623 1.626	1.624	68.6%
Cal-4: 0.063 µg/mL	1.320 1.265	1.293	54.6%
Cal-5: 0.125 µg/mL	1.030 1.034	1.032	43.5%
Cal-6: 0.250 µg/mL	0.721 0.799	0.760	32.1%
Cal-7: 0.500 µg/mL	0.598 0.561	0.580	24.5%
Cal-8: 1.000 µg/mL	0.413 0.373	0.393	16.6%
Cal-9: 2.000 µg/mL	0.291 0.331	0.311	13.1%

LIMITATION OF THE PROCEDURE

- (1) This assay requires serum or plasma sample for testing.
- (2) Serum or plasma samples from different species may show different matrix background. A modification of test procedure may be necessary for measuring rodent samples. Please contact Eagle Biosciences for technical support.
- (3) For sample values greater than 2 µg/mL, it is recommended to re-assay samples with dilution (i.e. 1:10 or 1:100 with zero calibrator matrix. This calibrator zero is available from kit manufacturer. Using a different buffer matrix for sample dilution may cause false high or low value because of matrix effect.). The best assay precision and most reliable test result is located between 15% B/B₀ to 85% B/B₀ of the standard curve.
- (4) Cell culture or tissue culture samples should be validated with total binding and other performance specifications before being used.
- (5) The kit calibrators are based on MMAF conjugated antibody or ADC concentration. It is not based on free MMAF concentration. The MMAF-ADC in different linker and DAR may give different curve shift.



QUALITY CONTROL

To assure the validity of the results each assay should include adequate controls.

PERFORMANCE CHARACTERISTICS

Sensitivity

The analytical sensitivity (LLOD) of this MMAF ADC EIA as determined by the 2 times standard deviation below the mean of B_0 on 12 duplicate determinations of zero standard (B_0) is approximately 0.019 $\mu\text{g/mL}$.

Specificity

This MMAF-ADC EIA does NOT show any cross reactivity to MMAE-ADC, DM1-ADC, DUO-3-ADC, and DUO-6-ADC.

High Dose "hook" effect

This assay has showed that it didn't have any high dose "hook" effect for MMAF ADC levels up to 1,000 $\mu\text{g/mL}$.

Precision

The intra-assay precision was validated by measuring three calibrators (L3, L4 and L7) in six replicate determinations. The CV% is 5.5%, 4.9% and 6.1%.

Linearity

Two samples were diluted with zero calibrator matrix and tested. The results of MMAF ADC dilution recovery value are as follows:

DILUTION	OBSERVED VALUE ($\mu\text{g/mL}$)	EXPECTED VALUE ($\mu\text{g/mL}$)	RECOVERY
Calibrator 5	0.125	0.125	-
20% + 80% buffer	0.024	0.025	96.0%
40% + 60% buffer	0.045	0.050	90.0%
60% + 40% buffer	0.049	0.075	105.3%
80% + 20% buffer	0.101	0.100	101.0%
Calibrator 7	0.500	0.500	-
20% + 80% buffer	0.109	0.100	109.0%
40% + 60% buffer	0.215	0.200	107.5%
60% + 40% buffer	0.347	0.300	111.3%
80% + 20% buffer	0.473	0.400	115.5%



Spike Recovery

Calibrator level 5 and 7 is equal volume mixed with calibrator level 4, 6, and 8 and tested. The results are as follows:

Spiked Sample	OBSERVED VALUE (µg/mL)	EXPECTED VALUE (µg/mL)	RECOVERY
Cal. 5	0.125	0.125	-
Cal. 5+Cal.-4 (0.063)	0.104	0.094	110.6%
Cal. 5+Cal.-6 (0.250)	0.208	0.188	110.9%
Cal. 5+Cal.-8 (1.000)	0.626	0.563	100.1%
Cal. 7	0.500	0.500	-
Cal. 7+Cal.-4 (0.063)	0.331	0.282	117.6%
Cal. 7+Cal.-6 (0.250)	0.407	0.375	108.5%
Cal. 7+Cal.-8 (1.000)	0.698	0.750	93.1%

REFERENCES

- (1) Sandhya Girish, et al. Clinical pharmacology of trastuzumab emtansine (T-DM1): an antibody–drug conjugate in development for the treatment of HER2-positive cancer. Cancer Chemother Pharmacol (2012) 69:1229–1240

Warranty Information

Eagle Biosciences, Inc. warrants its Product(s) to operate or perform substantially in conformance with its specifications, as set forth in the accompanying package insert. This warranty is expressly limited to the refund of the price of any defective Product or the replacement of any defective Product with new Product. This warranty applies only when the Buyer gives written notice to the Eagle Biosciences within the expiration period of the Product(s) by the Buyer. In addition, Eagle Biosciences has no obligation to replace Product(s) as result of a) Buyer negligence, fault, or misuse, b) improper use, c) improper storage and handling, d) intentional damage, or e) event of force majeure, acts of God, or accident.

Eagle Biosciences makes no warranties, either expressed or implied, except as provided herein, including without limitation thereof, warranties as to marketability, merchantability, fitness for a particular purpose or use, or non-infringement of any intellectual property rights. In no event shall the company be liable for any indirect, incidental, or consequential damages of any nature, or losses or expenses resulting from any defective product or the use of any product. Product(s) may not be resold, modified, or altered for resale without prior written approval from Eagle Biosciences, Inc.

For further information about this kit, its application or the procedures in this kit, please contact the Technical Service Team at Eagle Biosciences, Inc. at info@eaglebio.com or at 617-419-2019.