



EAGLE
BIOSCIENCES

DM1 Antibody Drug Conjugate (ADC) ELISA Assay Kit

Catalog Number:

KTR-756 (1 x 96 wells)

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

v. 1.2 (14 NOV 23)

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



INTENDED USE

The Eagle Biosciences DM1 Antibody Drug Conjugate (ADC) ELISA Assay Kit (enzyme-linked immunoassay kit) is intended for use in the quantitative determination of antibody-DM1-conjugate (Trastuzumab emtansine) level in Human/mouse/rat/primate serum.

*Both humanized monoclonal antibody based DM1-ADC and mouse monoclonal antibody based DM1-ADC can be measured with this ELISA Assay kit.

For further information about this kit, its application, or the procedures in this insert, please contact the Technical Service Team at Eagle Biosciences, Inc at www.EagleBio.com or at 866-411-8023.

PRINCIPLE OF THE ASSAY

The Eagle Biosciences DM1 Antibody Drug Conjugate (ADC) ELISA Assay Kit is designed, developed and produced for the quantitative measurement of antibody DM1 conjugate (Trastuzumab emtansine) in serum, tissue, and cell culture samples. The assay utilizes the competitive immunoassay technique with an antibody that exclusively binds to DM1.

Assay calibrators (antibody DM1 conjugate) and test serum samples are added directly to wells of a microtiter plate that is coated with specific anti-DM1 antibody. Subsequently, a horseradish peroxidase (HRP) conjugated DM1 is added to each well. During the incubation period, the antibody DM1 conjugate competes with the HRP conjugated DM1 for the limited binding sites of anti-DM1 antibody. An immune complex of well coated "anti-DM1 antibody – HRP conjugated DM1" 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-DM1 conjugate in the test sample. A calibration curve is generated by plotting the absorbance versus the respective antibody-DM1 conjugate concentration for each calibrator on a 4-parameter or log-logit curve fitting. The concentration of antibody-DM1 conjugate in test samples is determined directly from this calibration curve.

REAGENTS

Preparation and Storage

This DM1 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 are stable until the expiration date.

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

1. Anti-DM1 Antibody Coated Microplate

One microplate with twelve by eight strips (96 wells total) coated with specific anti-DM1 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 DM1

One vial containing **3 mL** of ready to use HRP labeled DM1 in a stabilized protein matrix. This reagent should be stored at 2 – 8°C and is stable until the expiration date on the kit box.

3. ELISA Wash Concentrate

One bottle containing **30 mL** of 30x 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 kit box.

4. ELISA HRP Substrate

One bottle containing **12 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 kit box.

5. ELISA Stop Solution

One bottle containing **12 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 kit box.

6. Antibody-DM1 Conjugated Zero Calibrator

One vial containing **30mL** calibrator zero. 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 kit box.

7. Antibody-DM1 Conjugated Calibration Stock

One vial containing the calibration stock of antibody-DM1-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 Human DM1 Antibody Drug Conjugate (ADC) Assay Kit reagents must be used in a professional laboratory environment and is for Research Use Only and is not to be used in diagnostic procedures. Only source material from which reagents of bovine serum was derived in the contiguous 48 United States. It was obtained only from donor health animals maintained under veterinary supervision and found free of contagious diseases. Wear gloves while performing this assay and handle these reagents as if they were 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.

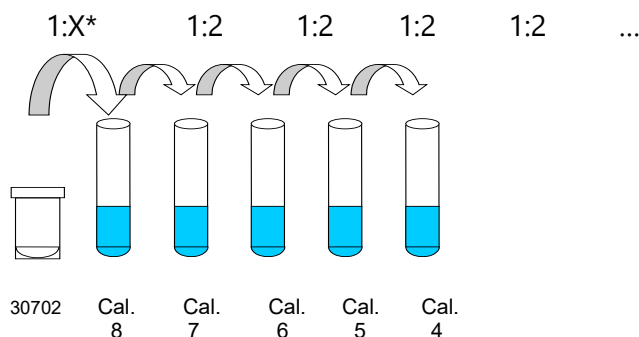


MATERIALS REQUIRED BUT NOT PROVIDED

- Antibody-DM1 Conjugated Stock
- Precision single channel pipettes capable of delivering 25 μ L, 50 μ L, and 100 μ L, etc.
- Disposable pipette tips suitable for above volume dispensing.
- Aluminum foil.
- Deionized or distilled water.
- Plastic microtiter well cover or polyethylene film.
- ELISA multichannel wash bottle or automatic (semi-automatic) washing system.
- Spectrophotometric microplate reader capable of reading absorbance at 450/650 nm or 450/620 nm.
- Calibrated Timer

REAGENT PREPARATION

1. Prior to use allow all reagents 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 with **0.5 mL** DI-water. Dilute the reconstituted calibration stock 1:X* using the zero calibrator to obtain a level eight calibrator at 1 μ g/mL. Further create calibrator level seven to two by 1:2 serial dilutions to obtain these calibrators with concentrations of 0.5 μ g/mL, 0.25 μ g/mL, 0.125 μ g/mL, 0.063 μ g/mL, 0.031 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 30702 / 1

4. Test Configuration

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



The validation data of this test was generated by using Antibody-DM1 Conjugated Stock

Assay Procedure

1. Place a sufficient number of Anti-DM1 antibody coated microwell strips/wells in a holder to run human Anti-DM1 standards, controls and unknown samples in duplicate.
2. Add 100 μL of calibrators and test samples into the designated microwells.
3. 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.
4. Immediately add **25 μL** of HRP Conjugated DM1 to each well. (*Note: no wash step before add the HRP Conjugated DM1*)
5. 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. \pm 10 minutes** at 400 to 450 rpm.
6. 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.
7. Add **100 μL** of ELISA HRP Substrate into each of the wells.
8. Cover the plate with aluminum foil or other material to avoid exposure to light. Incubate plate static, at room temperature for **20 minutes**.
9. Immediately add **100 μL** of ELISA Stop Solution into each of the wells. Mix gently.
10. Read the absorbance at 450 nm.

PROCEDURAL NOTES

1. It is recommended that all calibrators and unknown samples be assayed in duplicate. 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 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 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.

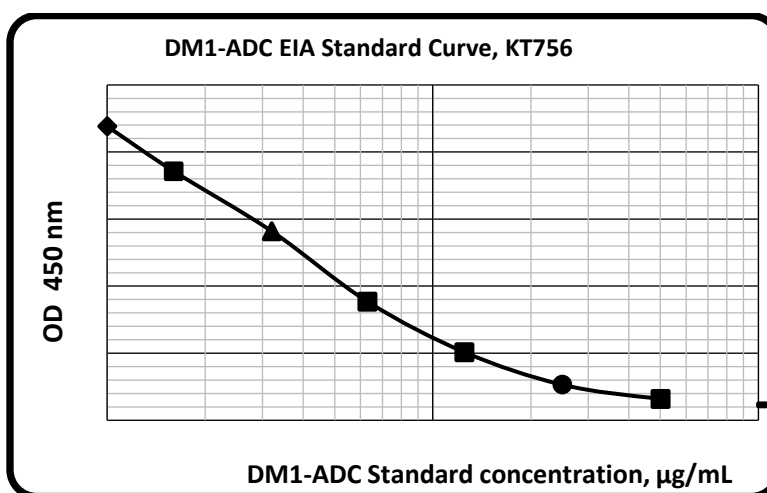


The antibody-DM1 conjugate concentrations for the test samples are read directly from the calibration curve using their respective corrected absorbance.

EXAMPLE DATA AND STANDARD CURVE

A typical absorbance data and the resulting standard curve from Human DM1 Antibody Drug Conjugate (ADC) ELISA Assay Kit are represented. This curve should not be used in lieu of standard curve run with each assay.

Well I.D.	OD 450 nm Absorbance		B/B ₀
	Readings	Average	
Cal-1: 0.000 µg/mL	2.234 2.150	2.192	100.0%
Cal-2: 0.016 µg/mL	1.753 1.960	1.857	84.7%
Cal-3: 0.032 µg/mL	1.446 1.375	1.410	64.3%
Cal-4: 0.063 µg/mL	0.902 0.868	0.885	40.4%
Cal-5: 0.125 µg/mL	0.495 0.520	0.508	23.2%
Cal-6: 0.250 µg/mL	0.260 0.270	0.265	12.1%
Cal-7: 0.500 µg/mL	0.157 0.165	0.161	7.3%
Cal-8: 1.000 µg/mL	0.111 0.119	0.115	5.2%



LIMITATION OF THE PROCEDURE

1. Serum or plasma samples from different species may show different matrix background. A modification of test procedure may be necessary for measuring rodent samples.
2. For sample values greater than 1 µg/mL, it is recommended to re-assay samples with dilution (i.e. 1:10 or 1:100 with calibrator zero). 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 10% B/B₀ to 85% B/B₀ of the standard curve.

3. Cell culture or tissue culture samples should be validated with total binding and other performance specifications before being used.
4. The kit calibrators are based on DM1 conjugated antibody or ADC concentration. It is not based on free DM1 concentration. The DM1-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 DM1 ADC EIA as determined by the 2 times calibrator deviation below the mean of B₀ on 8 duplicate determinations of zero calibrator (B₀) is approximately 0.024 µg/mL.

Specificity

This DM1-ADC EIA doesn't show any cross reactivity to MMAE-ADC, MMAF-ADC, DUO-3 ADC, or DUO-6 ADC.

High Dose "hook" effect

This assay has showed that it didn't have any high dose "hook" effect for DM1 ADC levels up to 1,000 µg/mL.

Precision

The intra-assay precision was validated by measuring three calibrators (L3, L6 and L8) in eight replicate determinations. The CV% is 6.8%, 6.2% and 8.9%.

Linearity

Two samples were diluted with calibrator zero and tested. The results of DM1 ADC dilution recovery value are as follows:

DILUTION	OBSERVED VALUE (µg/mL)	EXPECTED VALUE (µg/mL)	% RECOVERY
Calibrator 6	0.250	-	-
20% + 80% buffer	0.045	0.05	111.1%
40% + 60% buffer	0.104	0.1	96.2%
60% + 40% buffer	0.149	0.15	100.7%
80% + 20% buffer	0.219	0.2	91.3%
Calibrator 8	1	-	-
20% + 80% buffer	0.225	0.2	88.9%
40% + 60% buffer	0.488	0.4	82.0%
60% + 40% buffer	0.578	0.6	103.8%
80% + 20% buffer	0.735	0.8	108.8%

Spike Recovery

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

Spiked Sample	OBSERVED VALUE (µg/mL)	EXPECTED VALUE (µg/mL)	RECOVERY
Calibrator 5	0.125	0.125	-
Cal. 5+Cal.-4 (0.063)	0.085	0.094	90.4%
Cal. 5+Cal.-6 (0.250)	0.193	0.188	102.9%
Cal. 5+Cal.-8 (1.000)	0.637	0.563	113.2%
Calibrator 7	0.500	0.500	-
Cal. 5+Cal.-4 (0.063)	0.328	0.282	116.5%
Cal. 5+Cal.-6 (0.250)	0.378	0.375	100.8%
Cal. 5+Cal.-8 (1.000)	0.719	0.750	95.9%

REFERENCES

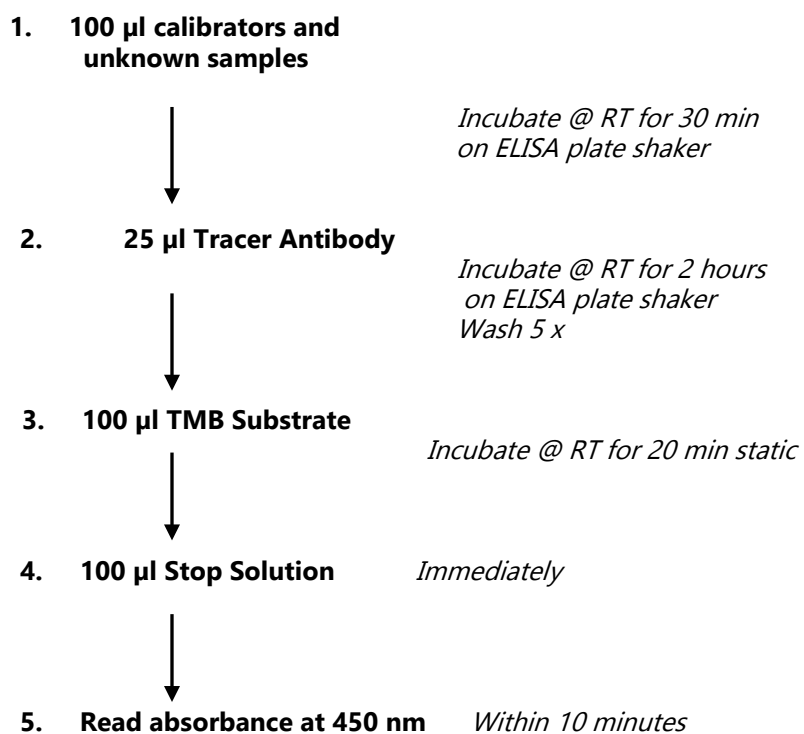
1. Sandhya Girish, et al. Clinical pharmacology of trastuzumab emtansine (T-DM1): an antibody–drug conjugate in development for the treatment of HER2-positive cancer. *Cancer Chemother Pharmacol* (2012) 69:1229–1240



PRECAUTIONS

- This kit is for research use only
- Compare contents and packing list, if there is breakage or shortage, notify Eagle Biosciences immediately
- Do not pipette reagents by mouth
- Do not smoke, eat or drink while performing assay
- Wear disposable gloves and proper lab protection and attire
- Treat all samples as potentially infectious
- Do not mix reagents from other lots
- Avoid contact with TMB and Stop solutions. If contact occurs, rinse thoroughly with water
- Eagle Biosciences is not responsible for outcomes as results of tampering with the reagents or using them unconventionally

DM1- ADC EIA: Condensed Assay Protocol





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 866-411-8023.