

# Product Information



## Anti-Endothelin EIA Strip Plate

Item No. 483152

### IMPORTANT

Water used to prepare all EIA reagents and buffers must be deionized and free of trace organic contaminants ('UltraPure'). Use activated carbon filter cartridges or other organic scavengers. Glass distilled water (even if double distilled), HPLC-grade water, and sterile water (for injections) are not adequate for EIA. UltraPure water may also be purchased (Item No. 400000).

### Laboratory Procedures

This plate is precoated with an Endothelin monoclonal antibody. Store at 4°C until ready to use. Immediately prior to use, open the plate packet and rinse the wells once with Wash Buffer. Be certain that all of the Wash Buffer is removed from each well before starting the assay. To accomplish this, the plate can be inverted and the last drops shaken out. *NOTE: If you do not need to use all the strips at once, place the unused strips back in the plate packet without rinsing and store at 2-4°C. Be sure the packet is sealed with the desiccant inside.*

### Buffer Preparation

#### 1. Sample Matrix Blank - Human plasma

The Sample Matrix Blank (SMB) provided (Item No. 400100) with this kit is IL-1 $\beta$ -free human plasma. If this matches your samples, reconstitute with 5 ml (100 dtn; 96-well kit) or 25 ml (500 dtn; 480-well kit) of UltraPure water. Store this solution at 4°C; it will be stable for approximately two weeks. If your SMB is not human plasma, you must obtain an IL-1 $\beta$ -free SMB that matches your samples.

*NOTE: The SMB provided is enough to run one standard curve once reconstituted. More SMB may be purchased by ordering Cayman's Human Plasma (Item No. 400100).*

#### 2. Phosphate Buffer

Prepare a 1.0 M phosphate buffer solution by combining 133 g K<sub>2</sub>HPO<sub>4</sub> and 32.15 g KH<sub>2</sub>PO<sub>4</sub> and diluting to a total volume of 1.0 liter with UltraPure water. The pH of this solution will be 7.4.

#### 3. EIA Buffer

Combine 100 ml of the phosphate buffer prepared above with 100 mg sodium azide, 23.4 g sodium chloride, 370 mg tetrasodium EDTA hydrate, and 1 g bovine serum albumin (Sigma A7030 or equivalent). Stir at room temperature until completely dissolved and dilute to a total volume of 1.0 liter with UltraPure water. Add 0.1 ml of Polysorbate 20 and mix thoroughly. This EIA Buffer may also be purchased from Cayman as a 10X concentrate buffer (Item No. 400060).

#### 4. Wash Buffer

Combine 10 ml of the phosphate buffer prepared above with 0.5 ml Polysorbate 20. Bring to a final volume of 1.0 liter with UltraPure water. This Wash Buffer may also be purchased from Cayman as a 400X concentrate buffer (Item No. 400062).

*NOTE: Polysorbate 20 is a viscous liquid and cannot be measured by a regular pipette. A positive displacement pipette or a syringe should be used to deliver small quantities accurately. It is normal for the concentrated buffer to contain crystalline salts after thawing. These will completely dissolve upon dilution with water.*

**WARNING: THIS PRODUCT IS FOR LABORATORY RESEARCH ONLY; NOT FOR ADMINISTRATION TO HUMANS. NOT FOR HUMAN OR VETERINARY DIAGNOSTIC OR THERAPEUTIC USE.**

#### MATERIAL SAFETY DATA

This material should be considered hazardous until information to the contrary becomes available. Do not ingest, swallow, or inhale. Do not get in eyes, on skin, or on clothing. Wash thoroughly after handling. This information contains some, but not all, of the information required for the safe and proper use of this material. Before use, the user must review the complete Material Safety Data Sheet, which has been sent *via* email to your institution.

#### WARRANTY AND LIMITATION OF REMEDY

Cayman Chemical Company makes **no warranty or guarantee** of any kind, whether written or oral, expressed or implied, including without limitation, any warranty of fitness for a particular purpose, suitability and merchantability, which extends beyond the description of the chemicals hereof. Cayman **warrants only** to the original customer that the material will **meet our specifications at the time of delivery.**

Cayman will carry out its delivery obligations with due care and skill. Thus, in no event will Cayman have **any obligation or liability**, whether in tort (including negligence) or in contract, for any direct, indirect, incidental or consequential damages, even if Cayman is informed about their possible existence.

This limitation of liability does not apply in the case of intentional acts or negligence of Cayman, its directors or its employees.

Buyer's **exclusive remedy** and Cayman's sole liability hereunder shall be limited to a **refund** of the purchase price, or at Cayman's option, the **replacement**, at no cost to Buyer, of all material that does not meet our specifications.

Said refund or replacement is conditioned on Buyer giving written notice to Cayman within thirty (30) days after arrival of the material at its destination. Failure of Buyer to give said notice within thirty (30) days shall constitute a waiver by Buyer of all claims hereunder with respect to said material.

For further details, please refer to our Warranty and Limitation of Remedy located on our website and in our catalog.

Copyright Cayman Chemical Company, 05/03/2011

## Cayman Chemical

### Mailing address

1180 E. Ellsworth Road  
Ann Arbor, MI  
48108 USA

### Phone

(800) 364-9897  
(734) 971-3335

### Fax

(734) 971-3640

### E-Mail

custserv@caymanchem.com

### Web

www.caymanchem.com

# Product Information



## Suggested Assay Protocol

This plate has been tested and formulated to work exclusively with ACE™ reagents. This plate may not perform as described if used with other assay reagents or protocols. The standard curve utilized in this assay should consist of a total of seven points made by 2-fold serial dilutions beginning at 1,000 pg/ml. The eighth point on the standard curve should only consist of sample matrix blank. *NOTE: This is an abbreviated protocol. If you are not familiar with this assay, please contact us for a complete protocol.*

1. Wash plate once with Wash Buffer.
2. Add 100 µl of Standard (Item No. 483154) or sample to the appropriate wells.
3. Add 100 µl Conjugate (Item No. 483150) to all wells except Blk.
4. Incubate for 90 minutes at 4°C.
5. Wash plate five times with Wash Buffer.
6. Add 200 µl Ellman's Reagent to each well.
7. Develop until S1 wells are visibly yellow.
8. Read absorbance between 405-420 nm.
9. Dilute samples may require longer development times.

Blk-Blank; S1-S8-Standards;  
1-7-Samples

	1	2	3	4
A	S1	S1	1	1
B	S2	S2	2	2
C	S3	S3	3	3
D	S4	S4	4	4
E	S5	S5	5	5
F	S6	S6	6	6
G	S7	S7	7	7
H	S8	S8	Blk	Blk

## Related Product

Endothelin EIA Kit - Item No. 583151

## Cayman Chemical

### Mailing address

1180 E. Ellsworth Road  
Ann Arbor, MI  
48108 USA

### Phone

(800) 364-9897  
(734) 971-3335

### Fax

(734) 971-3640

### E-Mail

[custserv@caymanchem.com](mailto:custserv@caymanchem.com)

### Web

[www.caymanchem.com](http://www.caymanchem.com)