

# Product Information



## STAT-8-Isoprostane EIA Standard

Item No. 400434

### 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

The STAT-8-Isoprostane (STAT-8-*iso* PGF<sub>2α</sub>) EIA Standard you have purchased contains approximately 500 μl of 8-isoprostane as a solution in ethanol. The concentration of this standard is 300 ng/ml. For long term storage, we suggest that the STAT-8-isoprostane EIA standard be stored as supplied at -20°C; it will be stable for at least one year. When ready to use, equilibrate a pipette tip in ethanol by repeatedly filling and expelling the tip with ethanol several times. Using the equilibrated pipette tip, transfer 100 μl of the 8-isoprostane EIA standard into a clean test tube, then dilute with 900 μl UltraPure water. The concentration of this solution (the bulk standard) will be 30 ng/ml. Store this solution at 4°C.

To prepare the standard for use in EIA: Obtain eight clean test tubes and number them #1 through #8. Aliquot 900 μl Tris Buffer to tube #1 and 500 μl Tris Buffer to tubes #2-8. Transfer 100 μl of the bulk standard (30 ng/ml) to tube #1 and mix thoroughly. Serially dilute the standard by removing 500 μl from tube #1 and placing in tube #2; mix thoroughly. Next, remove 500 μl from tube #2 and place it into tube #3; mix thoroughly. Repeat this process for tubes #4-8. The concentrations of these standards are 3,000, 1,500, 750, 375, 187.5, 93.8, 46.9, and 23.4 pg/ml. We recommend that you store these diluted standards for no more than 24 hours.

### Buffer Preparation

#### 1. Tris Buffer

Prepare a 0.1 M Tris-HCl buffer, pH 7.4, containing 0.15 M sodium chloride, 2 mM MgCl<sub>2</sub>, 0.1 mM ZnCl<sub>2</sub>, 0.05% azide, and 0.5% BSA solution by combining 12.11 g Tris base, 8.47 g sodium chloride, 0.406 g MgCl<sub>2</sub> • 6H<sub>2</sub>O, 0.014 g ZnCl<sub>2</sub>, 0.5 g sodium azide, and 5 g BSA in 800 ml UltraPure water and adjusting the pH to 7.4 with concentrated HCl. Stir at room temperature until completely dissolved and dilute to a total volume of 1.0 liter with UltraPure water. This buffer may also be purchased as a 10X concentrate buffer (Item No. 400080).

#### 2. Wash Buffer

Prepare a 0.01 M Tris-HCl buffer, pH 7.4, containing 0.05% Tween-20 by combining 1.3 g Tris-HCl and 0.194 g Tris base in 600 ml UltraPure water. Stir until dissolved and adjust the pH to 7.4 with HCl. Add 0.5 ml Tween 20 and dilute to a final volume of 1.0 liter with UltraPure water. This buffer may also be purchased as a 150X concentrated buffer (Item No. 411007).

#### 3. DEA Buffer

DEA Buffer may be purchased as a 10X concentrated buffer (Item No. 400082).

**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, 10/18/2010

### 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 standard has been tested and formulated to work exclusively with AP reagents. It is highly recommended that Cayman plates with Item No. 400004 or 400006 be used for this assay. This standard may not perform as described if used with other assay reagents or protocols. *NOTE: This is an abbreviated protocol. If you are not familiar with this assay, please contact us for a complete protocol.*

1. Add 100  $\mu$ l of Tris Buffer to NSB wells and 50  $\mu$ l to B<sub>0</sub> wells.
2. Add 50  $\mu$ l of Standard or sample to the appropriate wells.
3. Add 50  $\mu$ l Tracer (Item No. 400430) to all wells except Blk and TA.
4. Add 50  $\mu$ l Antiserum (Item No. 400432) to all wells except Blk, TA, and NSB.
5. Incubate one hour at room temperature on an orbital shaker.
6. Wash the plate five times with Wash Buffer.
7. Add 200  $\mu$ l pNPP to each well.
8. Add 5  $\mu$ l Tracer to the TA well.
9. Develop for approximately 60-90 minutes (B<sub>0</sub> = 0.3-1.0 AU) on an orbital shaker.
10. Read absorbance at a wavelength between 405 and 420 nm.

**Blk**-Blank; **NSB**-Non-specific Binding; **B<sub>0</sub>**-Maximum Binding; **TA**-Total Activity; **S1-S8**-Standards; **1-8**-Samples

	1	2	3	4
A	Blk	S1	S1	1
B	Blk	S2	S2	2
C	NSB	S3	S3	3
D	NSB	S4	S4	4
E	B <sub>0</sub>	S5	S5	5
F	B <sub>0</sub>	S6	S6	6
G	B <sub>0</sub>	S7	S7	7
H	TA	S8	S8	8

## Related Product

STAT-8-Isoprostane EIA Kit - Item No. 500431

## 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