



# Product Manual

## **LTB<sub>4</sub> ELISA kit**

Catalog #: ADI-900-068

96 Well Kit



# Product Manual

## **USE FOR RESEARCH PURPOSES ONLY**

Unless otherwise specified expressly on the packaging, all products sold hereunder are intended for and may be used for research purposes only and may not be used for food, drug, cosmetic or household use or for the diagnosis or treatment of human beings. Purchase does not include any right or license to use, develop or otherwise exploit these products commercially. Any commercial use, development or exploitation of these products or development using these products without the express written authorization of Enzo Life Sciences, Inc. is strictly prohibited. Buyer assumes all risk and liability for the use and/or results obtained by the use of the products covered by this invoice whether used singularly or in combination with other products.

## **LIMITED WARRANTY; DISCLAIMER OF WARRANTIES**

These products are offered under a limited warranty. The products are guaranteed to meet all appropriate specifications described in the package insert at the time of shipment. Enzo Life Sciences' sole obligation is to replace the product to the extent of the purchasing price. All claims must be made to Enzo Life Sciences, Inc., within five (5) days of receipt of order. THIS WARRANTY IS EXPRESSLY IN LIEU OF ANY OTHER WARRANTIES OR LIABILITIES, EXPRESS OR IMPLIED, INCLUDING WARRANTIES OF MERCHANTABILITY, FITNESS FOR A PARTICULAR PURPOSE, AND NONINFRINGEMENT OF THE PATENT OR OTHER INTELLECTUAL PROPERTY RIGHTS OF OTHERS, AND ALL SUCH WARRANTIES (AND ANY OTHER WARRANTIES IMPLIED BY LAW) ARE EXPRESSLY DISCLAIMED.

## **TRADEMARKS AND PATENTS**

Several Enzo Life Sciences products and product applications are covered by US and foreign patents and patents pending.

**FOR RESEARCH USE ONLY.  
NOT FOR USE IN DIAGNOSTIC PROCEDURES.**



Please read entire booklet before proceeding with the assay.



Carefully note the handling and storage conditions of each kit component.



Please contact Enzo Life Sciences Technical Support if necessary.

## Table of Contents

Description .....	2
Introduction .....	2
Precautions .....	2
Materials Supplied .....	3
Storage .....	3
Other Materials Needed .....	3
Sample Handling .....	4
Procedural Notes .....	5
Reagent Preparation .....	6
Assay Procedure .....	7
Calculation of Results .....	8
Typical Results .....	8
Typical Standard Curve .....	9
Typical Quality Control Parameters .....	9
Performance Characteristics .....	10
Sample Dilution Recommendations .....	12
References .....	13
Contact Information .....	14

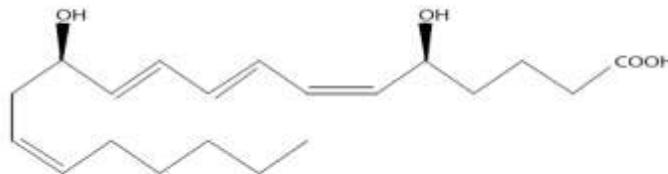
## DESCRIPTION

The LTB<sub>4</sub> ELISA kit is a competitive immunoassay for the quantitative determination of LTB<sub>4</sub> in biological fluids. Please read the complete kit insert before performing this assay. The kit uses a polyclonal antibody to LTB<sub>4</sub> to bind, in a competitive manner, the LTB<sub>4</sub> in the sample or an alkaline phosphatase molecule which has LTB<sub>4</sub> covalently attached to it. After incubations at room temperature the excess reagents are washed away and substrate is added. After a 2-hour incubation, the enzyme reaction is stopped and the yellow color generated is read on a microplate reader at 405nm. The intensity of the bound yellow color is inversely proportional to the concentration of LTB<sub>4</sub> in either standards or samples. The measured optical density is used to calculate the concentration of LTB<sub>4</sub>. For further explanation of the principles and practice of immunoassays please see the excellent books by Chard<sup>1</sup> or Tijssen<sup>2</sup>.

## INTRODUCTION

Leukotriene B<sub>4</sub> (LTB<sub>4</sub>) is a major product of arachidonic acid metabolism and is formed via the 5-lipoxygenase pathway<sup>3-5</sup>. LTB<sub>4</sub> stimulates leukocyte functions including lysosomal enzyme release<sup>6</sup>, adhesion<sup>7,8</sup>, and aggregation of polymorphonuclear leukocytes<sup>9</sup>. LTB<sub>4</sub> has been implicated as a potent mediator of inflammatory diseases<sup>10-13</sup> and immunoregulation<sup>14</sup>.

LTB<sub>4</sub>



## PRECAUTIONS

FOR RESEARCH USE ONLY. NOT FOR USE IN DIAGNOSTIC PROCEDURES.

1. Some kit components contain azide, which may react with lead or copper plumbing. When disposing of reagents always flush with large volumes of water to prevent azide build-up.
2. Stop Solution is a solution of trisodium phosphate. This solution is caustic; care should be taken in use.
3. The activity of the alkaline phosphatase conjugate is dependent on the presence of Mg<sup>2+</sup> and Zn<sup>2+</sup> ions. The activity of the conjugate is affected by concentrations of chelators (>10 mM) such as EDTA and EGTA.
4. We test this kit's performance with a variety of samples, however it is possible that high levels of interfering substances may cause variation in assay results.
5. The Leukotriene B<sub>4</sub> Standard provided, Catalog No. 80-0623, is supplied in ethanolic buffer at a pH optimized to maintain LTB<sub>4</sub> integrity. Care should be taken in handling this material because of the known and unknown effects of LTB<sub>4</sub>.



Handle  
with care

## MATERIALS SUPPLIED

1. **Goat anti-Rabbit IgG Microtiter Plate, One Plate of 96 Wells; Catalog No. 80-0060:**  
A plate using break-apart strips coated with goat antibody specific to rabbit IgG.
2. **LTB<sub>4</sub> ELISA Conjugate, 5 mL, Catalog No. 80-0066;**  
A blue solution of alkaline phosphatase conjugated with LTB<sub>4</sub>.
3. **LTB<sub>4</sub> ELISA Antibody, 5 mL, Catalog No. 80-0622:**  
A yellow solution of a rabbit polyclonal antibody to LTB<sub>4</sub>.
4. **Assay Buffer, 27 mL, Catalog No. 80-0010:**  
Tris buffered saline, containing proteins and detergents and sodium azide as preservative.
5. **Wash Buffer Concentrate, 27 mL, Catalog No. 80-1286:**  
Tris buffered saline containing detergents.
6. **Leukotriene B<sub>4</sub> Standard, 0.5 mL, Catalog No. 80-0623:**  
A solution of 120,000 pg/mL LTB<sub>4</sub>
7. **pNpp Substrate, 20 mL; Catalog No. 80-0075:**  
A solution of p-nitrophenyl phosphate in buffer. Ready to use.
8. **Stop Solution, 5 mL; Catalog No. 80-0247:**  
A solution of trisodium phosphate in water. Keep tightly capped. Caution: Caustic.
9. **LTB<sub>4</sub> Assay Layout Sheet, 1 each, Catalog No. 30-0115.**
10. **Plate Sealer, 1 each, Catalog No. 30-0012**



Reagents  
require  
separate  
storage  
conditions

## STORAGE

All components of this kit, except the conjugate, are stable at 4°C until the kit's expiration date. The conjugate must be stored frozen at -20°C.

## OTHER MATERIALS NEEDED

1. Deionized or distilled water.
2. Precision pipets for volumes between 5 µL and 1,000 µL.
3. Repeater pipets for dispensing 50 µL and 200 µL.
4. Disposable beakers for diluting buffer concentrates.
5. Graduated cylinders.
6. A microplate shaker.
7. Adsorbent paper for blotting.
8. A 37°C incubator.
9. Microplate reader capable of reading at 405nm, preferably with correction between 570 and 590nm.

## SAMPLE HANDLING

The LTB<sub>4</sub> enzyme-linked immunosorbent assay is compatible with LTB<sub>4</sub> samples in a wide range of matrices. Samples diluted sufficiently into Assay Buffer can be read directly from the standard curve. Please refer to the Sample Recovery recommendations for details of suggested dilutions. However, the end user must verify that the recommended dilutions are appropriate for their samples. Samples containing rabbit IgG may interfere with the assay.

Samples in the majority of tissue culture media, including those containing fetal bovine serum, can also be read in the assay, provided the standards have been diluted into the tissue culture media instead of Assay Buffer. There will be a small change in binding associated with running the standards and samples in media. Users should only use standard curves generated in media or buffer to calculate concentrations of LTB<sub>4</sub> in the appropriate matrix. For tissue, urine and plasma samples, prostaglandin synthetase inhibitors, such as, indomethacin or meclofenamic acid at concentrations up to 10 µg/mL should be added to either the tissue homogenate or urine and plasma samples.

Some samples normally have very low levels of LTB<sub>4</sub> present and extraction may be necessary for accurate measurement. A suitable extraction procedure is outlined below:

### Materials Needed

1. LTB<sub>4</sub> Standard to allow extraction efficiency to be accurately determined.
2. 2M hydrochloric acid, deionized water, ethanol, hexane and ethyl acetate.
3. 200 mg C18 Reverse Phase Extraction Columns.

## Procedure

1. Acidify the plasma, urine or tissue homogenate by addition of 2M HCl to pH of 3.5. Approximately 50  $\mu$ L of HCl will be needed per mL of plasma. Allow to sit at 4°C for 15 minutes. Centrifuge samples in a microcentrifuge for 2 minutes to remove any precipitate.
2. Prepare the C18 reverse phase column by washing with 10 mL of ethanol followed by 10 mL of deionized water.
3. Apply the sample under a slight positive pressure to obtain a flow rate of about 0.5 mL/minute. Wash the column with 10 mL of water, followed by 10 mL of 15% ethanol, and finally 10 mL hexane. Elute the sample from the column by addition of 10 mL ethyl acetate.
4. If analysis is to be carried out immediately, evaporate samples under a stream of nitrogen. Add at least 250  $\mu$ L of Assay Buffer to the dried samples. Vortex well, then allow to sit for five minutes at room temperature. Repeat twice more. If analysis is to be delayed, store samples as the eluted ethyl acetate solutions at -80°C until the immunoassay is to be run. Evaporate the organic solvent under a stream of nitrogen prior to running assay and reconstitute as above.

Please refer to references 15-18 for details of extraction protocols.

## PROCEDURAL NOTES

1. Do not mix components from different kit lots or use reagents beyond the kit expiration date.
2. Allow all reagents to warm to room temperature for at least 30 minutes before opening.
3. Standards can be made up in either glass or plastic tubes.
4. Pre-rinse the pipet tip with reagent, use fresh pipet tips for each sample, standard and reagent.
5. Pipet standards and samples to the bottom of the wells.
6. Add the reagents to the side of the well to avoid contamination.
7. This kit uses break-apart microtiter strips, which allow the user to measure as many samples as desired. Unused wells must be kept desiccated at 4°C in the sealed bag provided. The wells should be used in the frame provided.
8. Care must be taken to minimize contamination by endogenous alkaline phosphatase. Contaminating alkaline phosphatase activity, especially in the substrate solution, may lead to high blanks. Care should be taken not to touch pipet tips and other items that are used in the assay with bare hands.
9. Prior to addition of substrate, ensure that there is no residual wash buffer in the wells. Any remaining wash buffer may cause variation in assay results.

## REAGENT PREPARATION

### 1. LTB<sub>4</sub> Standard

Allow the 120,000 pg/mL LTB<sub>4</sub> standard solution to warm to room temperature. Label five 12x75 mm tubes #1 through #5. Pipet 975 µL of standard diluent (Assay Buffer or Tissue Culture Media) into tube #1. Pipet 750 µL of standard diluent into tubes #2 - #5. Add 25 µL of the 120,000 pg/mL standard to tube #1. Vortex thoroughly. Add 250 µL of tube #1 to tube #2 and vortex thoroughly. Add 250 µL of tube #2 to tube #3 and vortex thoroughly. Continue this for tubes #4 and #5.

The concentration of LTB<sub>4</sub> in tubes #1 through #5 will be 3,000, 750, 188, 46.9 and 11.7 pg/mL respectively. See LTB<sub>4</sub> Assay Layout Sheet for dilution details.

**Diluted standards should be used within 60 minutes of preparation.**

### 2. LTB<sub>4</sub> Conjugate

Allow the conjugate to warm to room temperature. Any unused conjugate should be dispensed into discrete aliquots and re-frozen at or below -20°C. Avoid repeated freeze-thaws of the aliquots.

### 3. Wash Buffer

Prepare Wash Buffer by diluting 5ml of the supplied concentrate with 95ml of deionized water. This can be stored at room temperature until the kit expiration, or for 3 months, whichever is earlier.

## ASSAY PROCEDURE

**Bring all reagents to room temperature for at least 30 minutes prior to opening. All standards and samples should be run in duplicate.**

1. Refer to the Assay Layout Sheet to determine the number of wells to be used and put any remaining wells with the desiccant back into the pouch and seal the ziploc. Store unused wells at 4°C.
2. Pipet 100 µL of standard diluent (Assay Buffer or Tissue Culture Media) into the NSB and the Bo (0 pg/mL Standard) wells.
3. Pipet 100 µL of Standards #1 through #5 into the appropriate wells.
4. Pipet 100 µL of the Samples into the appropriate wells.
5. Pipet 50 µL of Assay Buffer into the NSB wells.
6. Pipet 50 µL of blue Conjugate into each well, except the Total Activity (TA) and Blank wells.
7. Pipet 50 µL of yellow Antibody into each well, except the Blank, TA and NSB wells.

**NOTE:** Every well should be **Green** in color except the NSB wells which should be **Blue**. The Blank and TA wells are empty at this point and have no color.

8. Incubate the plate at room temperature on a plate shaker for 2 hours at ~500 rpm.\*
9. Empty the contents of the wells and wash by adding 400 µL of wash solution to every well. Repeat the wash 2 more times for a total of **3 Washes**.
10. After the final wash, empty or aspirate the wells, and firmly tap the plate on a lint free paper towel to remove any remaining wash buffer.
11. Add 5 µL of the blue Conjugate to the TA wells.
12. Add 200 µL of the pNpp Substrate solution to every well. Seal the plate and incubate at 37°C for 2 hours.
13. Add 50 µL of Stop Solution to every well. This stops the reaction and the plate should be read immediately.
14. Blank the plate reader against the Blank wells, read the optical density at 405 nm, preferably with correction between 570 and 590 nm. If the plate reader is not able to be blanked against the Blank wells, manually subtract the mean optical density of the Blank wells from all readings.

\* The plate shaker speed was based on a BellCo Mini Orbital Shaker (mod no. 7744-08096). The actual speed of the plate shaker should be such that the liquid in the plate wells mixes thoroughly, but does not splash out of the well.

## CALCULATION OF RESULTS

Several options are available for the calculation of the concentration of LTB<sub>4</sub> in the samples. We recommend that the data be handled by an immunoassay software package utilizing a 4 parameter logistic curve fitting program. If data reduction software (such as AssayBlaster™, catalog number ADI-28-0002) is not readily available, the concentration of LTB<sub>4</sub> can be calculated as follows:

1. Calculate the average Net Optical Density (OD) bound for each standard and sample by subtracting the average NSB OD from the average OD bound:

$$\text{Average Net OD} = \text{Average Bound OD} - \text{Average NSB OD}$$

2. Calculate the binding of each pair of standard wells as a percentage of the maximum binding wells (Bo), using the following formula:

$$\text{Percent Bound} = \frac{\text{Net OD}}{\text{Net Bo OD}} \times 100$$

3. Plot Percent Bound versus Concentration of LTB<sub>4</sub> for the standards. Approximate a straight line through the points. The concentration of LTB<sub>4</sub> in the unknowns can be determined by interpolation.

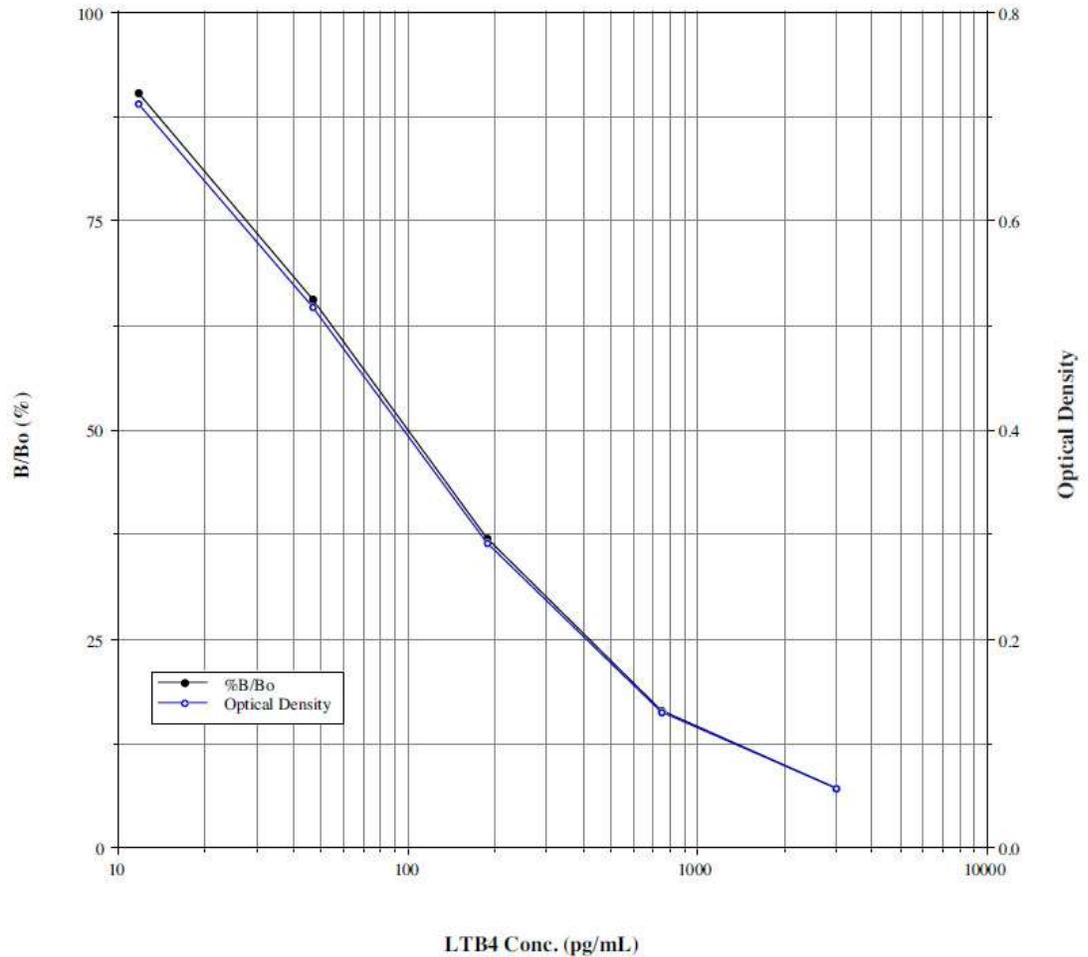
## TYPICAL RESULTS

The results shown below are for illustration only and should not be used to calculate results.

Sample	Mean OD (-Blank)	Average Net OD	Percent Bound	LTB <sub>4</sub> (pg/mL)
Blank	(0.145)			
TA	1.466			
NSB	-0.002	0.000		
Bo	0.786	0.788	100%	0
S1	0.054	0.056	7.0%	<b>3,000</b>
S2	0.127	0.129	16.4%	<b>750</b>
S3	0.289	0.291	36.9%	<b>188</b>
S4	0.515	0.517	65.6%	<b>46.9</b>
S5	0.710	0.712	90.4%	<b>11.7</b>
Unknown 1	0.465	0.467	59.3%	<b>63.5</b>
Unknown 2	0.155	0.157	19.9%	<b>556</b>

## TYPICAL STANDARD CURVES

A typical standard curve is shown below. This curve must not be used to calculate LTB<sub>4</sub> concentrations; each user must run a standard curve for each assay.



## Typical Quality Control Parameters

Total Activity Added =  $1.466 \times 10 = 14.66$   
 %NSB = 0.0%  
 %Bo/TA = 5.37%  
 Quality of Fit = 1.000

(Calculated from 4 parameter logistic curve fit)

20% Intercept = 552 pg/mL  
 50% Intercept = 98 pg/mL  
 80% Intercept = 23 pg/mL



## PERFORMANCE CHARACTERISTICS

The following parameters for this kit were determined using the guidelines listed in the National Committee for Clinical Laboratory Standards (NCCLS) Evaluation Protocols<sup>19</sup>.

### Sensitivity

Sensitivity was calculated by determining the average optical density bound for twenty (20) wells run as Bo, and comparing to the average optical density for twenty (20) wells run with Standard #5. The detection limit was determined as the concentration of LTB<sub>4</sub> measured at two (2) standard deviations from the zero along the standard curve.

Average Optical Density for the Bo = 0.802 ± 0.025 (3.2%)

Average Optical Density for Standard #5 = 0.696 ± 0.011 (1.6%)

Delta Optical Density (0-11.7 pg/mL) = 0.106

2 SD's of the Zero Standard = 2 x 0.025 = 0.051

Sensitivity =  $\frac{0.051}{0.106} \times 11.7 \text{ pg/mL}$  = 5.63 pg/mL

### Linearity

A sample containing 1,000 pg/mL LTB<sub>4</sub> was diluted 6 times 1:2 in the kit Assay Buffer and measured in the assay. The data was plotted graphically as actual LTB<sub>4</sub> concentration versus measured LTB<sub>4</sub> concentration.

The line obtained had a slope of 0.961 and a correlation coefficient of 0.999.

## Precision

Intra-assay precision was determined by taking samples containing low, medium and high concentrations of LTB<sub>4</sub> and running these samples multiple times (n=16) in the same assay. Inter-assay precision was determined by measuring three samples with low, medium and high concentrations of LTB<sub>4</sub> in multiple assays (n=8).

The precision numbers listed below represent the percent coefficient of variation for the concentrations of LTB<sub>4</sub> determined in these assays as calculated by a 4 parameter logistic curve fitting program.

	LTB <sub>4</sub> (pg/mL)	Intra-assay %CV	Inter-assay %CV
Low	305	6.0	
Medium	607	6.8	
High	1,078	5.9	
Low	99		15.7
Medium	308		16.5
High	507		5.0

## Cross Reactivities

The cross reactivities for a number of related compounds was determined by dissolving the cross reactant in Assay Buffer at concentrations from 40,000 to 0.4pg/mL. These samples were then measured in the LTB<sub>4</sub> assay and the measured LTB<sub>4</sub> concentration at 50% B/B<sub>0</sub> calculated. The % cross reactivity was calculated by comparison with the actual concentration of cross reactant in the sample and expressed as a percentage.

Compound	Cross Reactivity
LTB <sub>4</sub>	100%
6-trans-12-epi-LTB <sub>4</sub>	5.50%
6-trans-LTB <sub>4</sub>	4.90%
12-epi-LTB <sub>4</sub>	0.94%
PGE <sub>2</sub>	<0.2%
PGF <sub>2α</sub>	<0.2%
20-OH-LTB <sub>4</sub>	<0.2%
20-COOH-LTB <sub>4</sub>	<0.2%
LTC <sub>4</sub>	<0.2%
LTD <sub>4</sub>	<0.2%
LTE <sub>4</sub>	<0.2%
5(S)-HETE	<0.2%
12(S)-HETE	<0.2%
15(S)-HETE	<0.2%

## SAMPLE RECOVERIES

Please refer to pages 4 and 5 for Sample Handling recommendations and Standard preparation.

LTB<sub>4</sub> concentrations were measured in a variety of different samples including tissue culture media, human saliva and urine, and porcine plasma. For samples in tissue culture media, ensure that the standards have been diluted into the same media. LTB<sub>4</sub> was spiked into the undiluted samples of these media which were then diluted with the kit Assay Buffer and then assayed in the kit. The following results were obtained:

<b>Sample</b>	<b>% Recovery*</b>	<b>Recommended Dilution*</b>
Tissue Culture Media	97.3	None
Human Saliva	114.1	≥1:4
Human Urine	96.9	None
Porcine EDTA Plasma	109.6	1:2-1:4

\* See Sample Handling instructions on page 4 for details.

**REFERENCES**

1. Chard, T, "An Introduction. to Radioimmunoassay & Related Techniques 4th Ed.", (1990) Amsterdam:Elsevier
2. Tijssen, P, "Practice & Theory of Enzyme Immunoassays", (1985) Elsevier. Amsterdam:
3. B. Samuelsson & C.D. Funk, J. Biol. Chem., (1989) 264: 19469.
4. B. Samuelsson, et al., Science, (1987) 237: 1171.
5. P. Borgeat & P.H. Naccache, Clin. Biochem., (1990): 23: 459.
6. C.N. Serhan, et al., Biochem. Biophys. Res. Comm., (1982) 107: 1006.
7. S-E. Dahlen, et al., Proc. Natl. Acad. Sci., USA, (1981) 78: 3887.
8. P. Lindstrom, et al., Scand. J. Immunol., (1990) 31: 737.
9. F.M. Cunningham, et al., Agents Actions, (1981) 11: 583-4.
10. R.M. McMillan & S.J. Foster, Agents Actions, (1988) 24: 114.
11. S.D. Brain & T.J. Williams, Pharmacol. Ther., (1990) 46: 57.
12. W. König, et al., Eicosanoids, (1990) 3: 1.
13. A.W. Ford-Hutchinson, et al., Crit. Rev. Immunol., (1990) 10: 1.
14. M. Rola-Pleszczynski, J. Lipid Mediators, (1989) 1: 149.
15. K. Green, et al., Anal. Biochem, (1973) 54: 434.
16. J. Frolich, et al., J. Clin. Invest., (1975) 55: 763.
17. J.E. Shaw & P.W. Ramwell, Meth. Biochem. Anal., (1969) 17: 325.
18. K. Green, et al., Adv. Prostaglandin & Thromboxane Res., (1978) 5: 15.
19. National Committee for Clinical Laboratory Standards Evaluation Protocols, SC1, (1989) Villanova, PA:NCCLS.



# Product Manual

**Global Headquarters  
Enzo Life Sciences Inc.**

10 Executive Blvd  
Farmingdale, NY 11735  
(p) 1-800-942-0430  
(f) 1-631-694-7501  
(e) [info-usa@enzolifesciences.com](mailto:info-usa@enzolifesciences.com)

**Enzo Life Sciences (ELS) AG**

Industriestrasse 17, Postfach  
CH-4415 Lause / Switzerland  
(p) +41/0 61 926 89 89  
(f) +41/0 61 926 89 79  
(e) [info-ch@enzolifesciences.com](mailto:info-ch@enzolifesciences.com)

For local distributors and detailed product information visit us online:  
[www.enzolifesciences.com](http://www.enzolifesciences.com)