

Product Information



Interleukin-6 (human) EIA Standard

Item No. 483364

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 Interleukin-6 (IL-6) (human) EIA Standard you have purchased contains IL-6 (human) EIA Standard as a lyophilized powder. For long term storage, we suggest that the IL-6 (human) EIA Standard be stored as supplied at -20°C; it will be stable for at least one year. When ready to use, reconstitute the standard with 2 ml EIA Buffer (see buffer preparation instructions below). The concentration of this solution (the bulk standard) will be 5 ng/ml.

To prepare the standard for use in EIA: Obtain eight clean test tubes and number them #1 through #8. Aliquot 950 µl EIA Buffer to tube #1 and 500 µl EIA Buffer to tubes #2-8. Transfer 50 µl of the bulk standard (5 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 #4-7. Do not add any IL-6 to tube #8. This tube is the zero point vial, the lowest point on the standard curve. We recommend that you store these diluted standards for no more than 24 hours.

Buffer Preparation

1. Phosphate Buffer

Prepare a 1.0 M phosphate buffer solution by combining 133 g K_2HPO_4 and 32.15 g KH_2PO_4 and diluting to a total volume of 1.0 liter with UltraPure water. The pH of this solution will be 7.4.

2. 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, 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 Polysorbate 20 and mix thoroughly. This EIA Buffer may also be purchased from Cayman as a 10X concentrate buffer (Item No. 400060).

3. 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).

Suggested Assay Protocol

This standard has been tested and formulated to work exclusively with ACE™ reagents. 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. Wash plate once with Wash Buffer.
2. Add 100 µl of Standard or sample to the appropriate wells.
3. Add 100 µl Conjugate (Item No. 483360) 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

Interleukin-6 (human) EIA Kit - Item No. 583361

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, 03/21/2011

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