

PRODUCT INFORMATION



Resolvin D2 ELISA Antiserum

Item No. 401122

Overview

Contents:	This vial contains lyophilized Resolvin D2 Antiserum.
Storage:	-20°C
Stability:	≥18 months
Applications:	ELISA
Preparation:	Reconstitute the lyophilized antiserum with 6 ml of ELISA buffer (1X). Store reconstituted antiserum at 4°C and use within four weeks.

Suggested Competitive ELISA Protocol

Reagent Preparation

A source of ultrapure water, with a resistivity of 18.2 MΩ-cm and total organic carbon (TOC) levels of <10 ppb, is recommended. Pure water - glass-distilled or deionized - may not be acceptable. *NOTE: UltraPure Water is available for purchase from Cayman (Item No. 400000).*

1. **ELISA Buffer Concentrate (10X) (Item No. 400060)**

Dilute 10 ml of ELISA Buffer Concentrate (10X) with 90 ml of ultrapure water.

2. **Wash Buffer Concentrate (400X) (Item No. 400062)**

Dilute 5 ml of Wash Buffer Concentrate (400X) with ultrapure water to a total volume of 2 L and add 1 ml of Polysorbate 20 (Item No. 400035).

3. **Ellman's Reagent (Item No. 400050)**

Reconstitute 100 dtn vial with 20 ml of ultrapure water.

4. **Mouse Anti-Rabbit IgG Coated Plate (Item No. 400004)**

Ready to use as supplied.

5. **Resolvin D2 AChE Tracer (100 dtn vial; Item No. 401120)**

Reconstitute with 6 ml of ELISA Buffer (1X) before use. The reconstituted tracer will be stable for four weeks when stored at 4°C.

6. **Resolvin D2 ELISA Antiserum (100 dtn vial; Item No. 401122)**

Reconstitute with 6 ml of ELISA Buffer (1X) before use. The reconstituted antiserum will be stable for at least four weeks when stored at 4°C.

7. **Resolvin D2 ELISA Standard (Item No. 401076)**

Transfer 100 µl of the Resolvin D2 ELISA Standard into test tube and dilute with 900 µl of ultrapure water. The concentration of this bulk standard is 160 ng/ml. To prepare a standard curve (8,000 -1.24 pg/ml) for use in the assay obtain eight clean test tubes and number them #1-8. Aliquot 950 µl of ELISA Buffer (1X) to tube #1 and 500 µl to tubes #2-8. Transfer 50 µl of the bulk standard (160 ng/ml) to tube #1 and mix thoroughly. Serially dilute the standard by removing 200 µl from tube #1 and placing in tube #2; mix thoroughly. Next, remove 200 µl from tube #2 and place it into tube #3; mix thoroughly. Repeat this process for tubes #4-8. *NOTE: If assaying culture medium samples that have not been diluted with ELISA Buffer (1X), culture medium should be used in place of ELISA Buffer (1X) for dilution of the standard curve.*

WARNING

THIS PRODUCT IS FOR RESEARCH ONLY - NOT FOR HUMAN OR VETERINARY DIAGNOSTIC OR THERAPEUTIC USE.

SAFETY DATA

This material should be considered hazardous until further information becomes available. Do not ingest, inhale, get in eyes, on skin, or on clothing. Wash thoroughly after handling. Before use, the user must review the [complete](#) Safety Data Sheet, which has been sent via email to your institution.

WARRANTY AND LIMITATION OF REMEDY

Buyer agrees to purchase the material subject to Cayman's Terms and Conditions. Complete Terms and Conditions including Warranty and Limitation of Liability information can be found on our website.

Copyright Cayman Chemical Company, 12/18/2025

CAYMAN CHEMICAL

1180 EAST ELLSWORTH RD
ANN ARBOR, MI 48108 · USA

PHONE: [800] 364-9897

[734] 971-3335

FAX: [734] 971-3640

CUSTSERV@CAYMANCHEM.COM
WWW.CAYMANCHEM.COM

PRODUCT INFORMATION



Performing the Assay

1. Add 100 μ l of ELISA Buffer (1X) to the non-specific binding (NSB) wells and 50 μ l to B₀ wells. If culture media was used to dilute the standard curve, substitute 50 μ l of culture media for ELISA buffer (1X) in the NSB and B₀ wells (i.e., add 50 μ l culture medium to NSB and B₀ wells and 50 μ l ELISA Buffer (1X) to NSB wells).
2. Add 50 μ l of Standard or sample to the appropriate wells.
3. Add 50 μ l of Tracer (Item No. 401120) to all wells except blank (Blk) and total activity (TA).
4. Add 50 μ l Antiserum (Item No. 401122) to all wells except Blk, TA, and NSB.
5. Incubate for 2 hours at room temperature on an orbital shaker.
6. Wash the plate five times with Wash Buffer (1X).
7. Add 200 μ l of Ellman's Reagent to each well.
8. Add 5 μ l Tracer to the TA well.
9. To develop, incubate on an orbital shaker at room temperature for approximately 90 minutes.
10. Read absorbance at a wavelength between 405 and 420 nm.

Precaution

The suggested protocol above is supplied as a guideline. Users may need to optimize the assay conditions based on specific their applications.

Compound	Cross Reactivity
Resolvin D2	100%
12(S)-HHTrE	0.022%
5(S),6(R)-Lipoxin A ₄	0.02%
10(S),17(S)-DiHDoHE	0.012%
Docosahexaenoic Acid	<0.01%
17(R)-HDoHE	<0.01%
17(S)-HDoHE	<0.01%
17(S)-HpDoHE	<0.01%
Resolvin D1	<0.01%
Resolvin E1	<0.01%
17(R)-Resolvin D1	<0.01%
($\pm$)11(12)-DiHET	<0.01%
($\pm$)5(6)-DiHET	<0.01%
Prostaglandin A ₂	<0.01%
($\pm$)14(15)-DiHETE	<0.01%
($\pm$)8(9)-DiHET	<0.01%
($\pm$)8(9)-DiHETE	<0.01%
9(S),12(S),13(S)-TriHODE	<0.01%

	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 ₀	S5	S5	5
F	B ₀	S6	S6	6
G	B ₀	S7	S7	7
H	TA	S8	S8	8

Blk = Blank
 TA = Total Activity
 NSB = Non-Specific Binding
 B₀ = Maximum Binding
 S1-S8 = Standards
 1-8 - Samples

CAYMAN CHEMICAL

1180 EAST ELLSWORTH RD
ANN ARBOR, MI 48108 · USA

PHONE: [800] 364-9897
[734] 971-3335

FAX: [734] 971-3640

CUSTSERV@CAYMANCHEM.COM
WWW.CAYMANCHEM.COM