



PPAR α , δ , γ Complete Transcription Factor Assay Kit

Item No. 10008878

www.caymanchem.com

Customer Service 800.364.9897

Technical Support 888.526.5351

1180 E. Ellsworth Rd • Ann Arbor, MI • USA

TABLE OF CONTENTS

GENERAL INFORMATION	3	Materials Supplied
	4	Safety Data
	4	Precautions
	5	If You Have Problems
	5	Storage and Stability
	5	Materials Needed but Not Supplied
INTRODUCTION	6	Background
	7	About This Assay
PRE-ASSAY PREPARATION	9	Reagent Preparation
ASSAY PROTOCOL	11	Plate Set Up
	13	Performing the Assay
	17	Assay Procedure Summary
ANALYSIS	19	Performance Characteristics
RESOURCES	20	Interferences
	21	Troubleshooting
	23	References
	25	Plate Template
	26	Notes
	27	Warranty and Limitation of Remedy

GENERAL INFORMATION

Materials Supplied

Kit components may be stored at -20°C prior to use. For long term storage, the positive controls should be thawed on ice, aliquoted at 25 µl/vial, and stored at -80°C. After opening the kit, we recommend each kit component be stored according to the temperature listed below.

Item Number	Item	Quantity/Size	Storage
10006880	Transcription Factor Binding Assay Buffer (4X)	1 vial/3 ml	4°C
10007472	Transcription Factor Reagent A	1 vial/120 µl	-20°C
10008892	Transcription Factor Complete PPAR α Positive Control	1 vial/50 µl	-80°C
10008893	Transcription Factor Complete PPAR δ Positive Control	1 vial/50 µl	-80°C
10008894	Transcription Factor Complete PPAR γ Positive Control	1 vial/50 µl	-80°C
10006882	Transcription Factor Antibody Binding Buffer (10X)	1 vial/3 ml	4°C
10008895	Transcription Factor Complete PPAR α Primary Antibody	1 vial/50 µl	-20°C
10008896	Transcription Factor Complete PPAR δ Primary Antibody	1 vial/50 µl	-20°C
10008897	Transcription Factor Complete PPAR γ Primary Antibody	1 vial/50 µl	-20°C
400062	Wash Buffer Concentrate (400X)	1 vial/5 ml	RT
400035	Polysorbate 20	1 vial/3 ml	RT
10006885	Transcription Factor PPAR Competitor dsDNA	1 vial/120 µl	-20°C
10006884	Transcription Factor Goat Anti-Rabbit HRP Conjugate	1 vial/120 µl	-20°C
10006887	Transcription Factor PPAR 96-Well Strip Plate	1 plate	4°C
400012	96-Well Cover Sheet	1 cover	RT
10006888	Transcription Factor Developing Solution	1 vial/12 ml	4°C
10006889	Transcription Factor Stop Solution	1 vial/12 ml	RT

If any of the items listed above are damaged or missing, please contact our Customer Service department at (800) 364-9897 or (734) 975-3999. We cannot accept any returns without prior authorization.



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.

Precautions

Please read these instructions carefully before beginning this assay.

Kit components may be stored at -20°C prior to use. For long term storage, the Positive Controls should be thawed on ice, aliquoted at 25 µl/vial and stored at -80°C. If the assay will be used on multiple days, we recommend each kit component be stored according to the temperatures listed in the booklet.

If You Have Problems

Technical Service Contact Information

Phone: 888-526-5351 (USA and Canada only) or 734-975-3888
Fax: 734-971-3641
E-Mail: techserv@caymanchem.com

In order for our staff to assist you quickly and efficiently, please be ready to supply the lot number of the kit (found on the outside of the box).

Storage and Stability

This kit will perform as specified if stored as directed and used before the expiration date indicated on the outside of the box.

Materials Needed But Not Supplied

1. A plate reader capable of measuring absorbance at 450 nm
2. Adjustable pipettes and a repeating pipettor
3. An orbital microplate shaker
4. A source of UltraPure water; glass Milli-Q or HPLC-grade water are acceptable
5. 300 mM dithiothreitol (DTT)
6. Nuclear Extraction Kit available from Cayman (Item No. 10009277) or buffers for preparation of nuclear extracts

NOTE: The components in each kit lot have been quality assured and warranted in this specific combination only; please do not mix them with components from other lots.

Background

Peroxisome Proliferator-activated Receptors (PPARs) are ligand-activated transcription factors belonging to the large superfamily of nuclear receptors.^{1,2} They are activated by a variety of fatty acids and fatty acid derivatives such as prostaglandins and leukotrienes. PPARs play pivotal roles in the regulation of lipid metabolism and homeostasis and are important indirect as well as direct regulators of cellular insulin sensitivity.³ There are three major PPAR isotypes; PPAR α , PPAR γ , and PPAR δ/β which all bind to PPAR responsive elements (PPRE's) as heterodimers with RXR, another member of the nuclear receptor superfamily. PPAR α primarily activates genes encoding proteins involved in fatty acid oxidation, while PPAR γ primarily activates genes directly involved in lipogenic pathway and insulin signaling.^{1,4,5} Members of the PPAR family are important direct targets of many antidiabetic and hypolipidemic drugs.⁶

About This Assay

Cayman's PPAR α , δ , γ Complete Transcription Factor Assay is a non-radioactive, sensitive method for detecting specific transcription factor DNA binding activity in nuclear extracts and whole cell lysates. A 96-well enzyme-linked immunosorbent assay (ELISA) replaces the cumbersome radioactive electrophoretic mobility shift assay (EMSA). A specific double stranded DNA (dsDNA) sequence containing the PPAR response element is immobilized onto the bottom of wells of a 96-well plate (see Figure 1, on page 8). PPARs contained in a nuclear extract, bind specifically to the PPAR response element. PPAR α , δ , or γ are detected by addition of specific primary antibodies directed against the individual PPARs. A secondary antibody conjugated to HRP is added to provide a sensitive colorimetric readout at 450 nm. Cayman's PPAR α , δ , γ Complete Transcription Factor Assay comes with a single plate that measures all three isoforms of PPAR α , δ , and γ . There are enough reagents for one-third of a plate for each isoform.

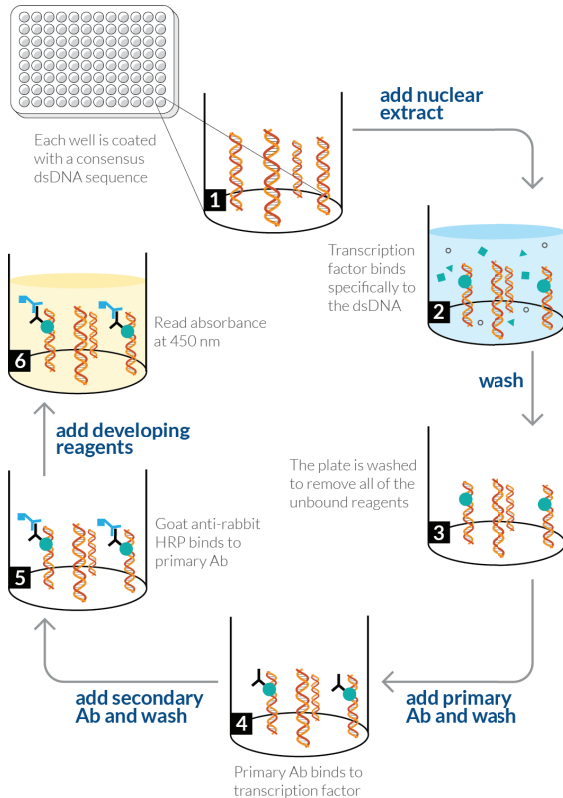


Figure 1. Schematic of the Transcription Factor Binding Assay

PRE-ASSAY PREPARATION

Reagent Preparation

1. Transcription Factor Antibody Binding Buffer (10X)

One vial (Item No. 10006882) contains 3 ml of a 10X stock of Transcription Factor Antibody Binding Buffer (ABB) to be used for diluting the primary and secondary antibodies. To prepare 1X ABB, dilute 1:10 by adding 27 ml of UltraPure water. Store at 4°C for up to six months.

2. Wash Buffer Concentrate (400X)

One vial (Item No. 400062) contains 5 ml of 400X Wash Buffer. Dilute the contents of the vial to a total volume of 2 liters with UltraPure water and add 1 ml of Polysorbate 20 (Item No. 400035). *NOTE: Polysorbate 20 is a viscous liquid and cannot be measured by a pipette. A positive displacement device such as a syringe should be used to deliver small quantities accurately. A smaller volume of Wash Buffer Concentrate can be prepared by diluting the Wash Buffer Concentrate 1:400 and adding Polysorbate 20 (0.5 ml/liter of Wash Buffer).* Store at 4°C for up to two months.

3. Complete Transcription Factor Binding Assay Buffer

Prepare 10 ml of Complete Transcription Factor Binding Assay Buffer (CTFB) by adding 2.5 ml of Transcription Factor Binding Buffer (4X) (Item No. 10006880), 0.1 ml of Transcription Factor Reagent A (Item No. 10007472), and 0.1 ml of 300 mM DTT to 7.3 ml of UltraPure water. Scale as necessary. *It is recommended that the CTFB be used the same day it is prepared.*

4. Transcription Factor PPAR α , δ , γ Positive Controls

Each vial (Item Nos. 10008892, 10008893, and 10008894) contains 50 μ l of clarified cell lysate. These lysates are provided as a positive controls for PPAR α , δ , and γ , respectively; they are not intended for plate to plate comparisons. The cell lysates provided are sufficient for 5 reactions and will provide a strong signal (>0.5 AU at 450 nm) when used at 10 μ l/well. When using this Positive Control, a decrease in signal may occur with repeated freeze/thaw cycles. It is recommended that the Positive Control be subaliquoted and stored at -80°C to avoid loss in signal from repeated freeze/thaw cycles.

ASSAY PROTOCOL

Plate Set Up

There is no specific pattern for using the wells on the plate. A typical layout of PPAR Positive Control (PC), Competitor dsDNA (C1), and samples of nuclear extracts (S1-S44) to be measured in duplicate is given below in Figure 2. We suggest you record the contents of each well on the template sheet provided (see page 25).

	1	2	3	4	5	6	7	8	9	10	11	12
A	S1	S1	S9	S9	S17	S17	S25	S25	S33	S33	S41	S41
B	S2	S2	S10	S10	S18	S18	S26	S26	S34	S34	S42	S42
C	S3	S3	S11	S11	S19	S19	S27	S27	S35	S35	S43	S43
D	S4	S4	S12	S12	S20	S20	S28	S28	S36	S36	S44	S44
E	S5	S5	S13	S13	S21	S21	S29	S29	S37	S37	NSB	NSB
F	S6	S6	S14	S14	S22	S22	S30	S30	S38	S38	PC	PC
G	S7	S7	S15	S15	S23	S23	S31	S31	S39	S39	Blk	Blk
H	S8	S8	S16	S16	S24	S24	S32	S32	S40	S40	C1	C1

S1-S44 - Sample Wells

NSB - Non-specific Binding Wells

PC - Positive Control Wells

Blk - Blank Wells

C1 - Competitor dsDNA Wells

Figure 2. Sample plate format

Pipetting Hints

- Use different tips to pipette each reagent.
- Before pipetting each reagent, equilibrate the pipette tip in that reagent (*i.e.*, slowly fill the tip and gently expel the contents, repeat several times).
- Do not expose the pipette tip to the reagent(s) already in the well.

General Information

- It is not necessary to use all the wells on the plate at one time; however a Positive Control should be run every time.
- For each plate or set of strips it is recommended that two Blk, two NSB, and two PC wells be included.

Performing the Assay

Binding of active PPAR α , δ , and/or γ to the consensus sequence

1. Equilibrate the plate and buffers to room temperature prior to opening. Remove the plate from the foil and select the number of strips needed. The 96-well plate supplied with this kit is ready to use.

NOTE: If you are not using all of the strips at once, place the unused strips back in the plate packet and store at 2-4°C. Be sure that the packet is sealed with the desiccant inside.

2. Prepare the CTFB as outlined on page 9
3. Add appropriate amount of reagent(s) listed below to the designated wells as follows:

Blk - add 100 μ l of CTFB to designated wells.

NSB - add 100 μ l of CTFB to designated wells. Do not add samples or Positive Control to these wells.

C1 - Add 80 μ l of CTFB prior to adding 10 μ l of Transcription Factor PPAR Competitor dsDNA (Item No. 10006885) to designated wells. Add 10 μ l of control cell lysate or sample.

NOTE: Competitor dsDNA must be added prior to adding the positive control or nuclear extracts.

S1-S44 - Add 90 μ l of CTFB followed by 10 μ l of Nuclear Extract to designated wells.

PC - Add 90 μ l of CTFB followed by 10 μ l of Positive Control to appropriate wells.

4. Use the 96-well cover sheet (Item No. 400012) provided to seal the plate. Incubate overnight at 4°C or one hour at room temperature on an orbital shaker.
5. Empty the wells and wash five times with 200 μ l of 1X Wash Buffer. After each wash empty the wells in the sink. After the final wash (wash #5), tap the plate on a paper towel to remove any residual Wash Buffer.

Addition of Transcription Factor PPAR Primary Antibody

1. Dilute the Transcription Factor PPAR Primary Antibody (Item Nos. 10008895, 10008896, and 10008897) 1:100 in 1X ABB as outlined in Table 1, below. Add 100 μ l of diluted PPAR Primary Antibodies to each well except the Blk wells.

Component	Volume/ Well	Volume/ Strip	Volume/ 96-well plate
1X ABB	99 μ l	792 μ l	9,504 μ l
PPAR Primary Antibody	1 μ l	8 μ l	96 μ l
Total required	100 μ l	800 μ l	9,600 μ l

Table 1. Dilution of Primary Antibody

2. Use the adhesive cover provided to seal the plate.
3. Incubate the plate for one hour at room temperature on an orbital shaker.
4. Empty the wells and wash each well five times with 200 μ l of 1X Wash Buffer. After each wash, empty the contents of the plate into the sink. After the final wash (wash #5), tap the plate three to five times on a paper towel to remove any residual Wash Buffer.

Addition of the Transcription Factor Goat Anti-Rabbit HRP Conjugate

1. Dilute the Transcription Factor Goat Anti-Rabbit HRP Conjugate (Item No. 10006884) 1:100 in 1X ABB as outlined in Table 2 below. Add 100 μ l of diluted secondary antibody to each well except the Blk wells.

Component	Volume/ Well	Volume/ Strip	Volume/ 96-well plate
1X ABB	99 μ l	792 μ l	9,504 μ l
Goat Anti-Rabbit HRP Conjugate	1 μ l	8 μ l	96 μ l
Total required	100 μ l	800 μ l	9,600 μ l

Table 2. Dilution of Secondary Antibody

2. Use the adhesive cover provided to seal the plate.
3. Incubate for one hour at room temperature on an orbital shaker.
4. Empty the wells and wash five times with 200 μ l of 1X Wash Buffer. After each wash, empty the contents of the plate into the sink. After the final wash (wash #5), tap the plate three to five times on a paper towel to remove any residual Wash Buffer.

Develop and Read the Plate:

1. To each well being used add 100 μ l of Transcription Factor Developing Solution (Item No. 10006888), which has been equilibrated to room temperature.
2. Incubate the plate for 30 minutes at room temperature on an orbital shaker protected from light. Allow the wells to turn medium to dark blue prior to adding Transcription Factor Stop Solution (Item No. 10006889).
3. Add 100 μ l of Stop Solution per well being used. The solution within the wells will change from blue to yellow after adding the Stop Solution.
4. Read absorbance at 450 nm within five minutes of adding the Stop Solution. Blank the plate reader according to the manufacturer's requirements using the blank wells.

Assay Procedure Summary

NOTE: This procedure is provided as a quick reference for experienced users. Follow the detailed procedure when initially performing the assay.

1. Prepare CTFB as described.
2. Add 90 μ l CTFB per sample well (80 μ l if adding Competitor dsDNA), 100 μ l to Blk and NSB wells).
3. Add 10 μ l of Competitor dsDNA (optional) to appropriate wells.
4. Add 10 μ l of Positive Control to appropriate wells.
5. Add 10 μ l of Sample containing PPAR α , δ , and/or γ to appropriate wells.
6. Incubate overnight at 4°C or one hour at room temperature on an orbital shaker.
7. Wash each well five times with 200 μ l of 1X Wash Buffer.
8. Add 100 μ l of diluted PPAR α , δ , and/or γ Primary Antibody per well (except Blk wells).
9. Incubate one hour at room temperature on an orbital shaker.
10. Wash each well five times with 200 μ l of 1X Wash Buffer.
11. Add 100 μ l of diluted Secondary Antibody (except Blk wells).
12. Incubate one hour at room temperature on an orbital shaker.
13. Wash each well five times with 200 μ l of 1X Wash Buffer.
14. Add 100 μ l of Developing Solution per well.
15. Incubate 30 minutes on an orbital shaker.
16. Add 100 μ l of Stop Solution per well.
17. Measure the absorbance at 450 nm.

Steps	Reagent	Blk	NSB	PC	C1	S1-S44
1. Add reagents	CTFB	100 μ l	100 μ l	90 μ l	80 μ l	90 μ l
	Competitor dsDNA				10 μ l	
	Positive Control			10 μ l	10 μ l	
	Samples					10 μ l
2. Incubate	Cover plate and incubate overnight at 4°C on an orbital shaker					
3. Wash	Wash all wells five times					
4. Add reagents	Primary Antibody		100 μ l	100 μ l	100 μ l	100 μ l
5. Incubate	Cover plate and incubate one hour at RT on an orbital shaker					
6. Wash	Wash all wells five times					
7. Add reagents	Secondary Antibody		100 μ l	100 μ l	100 μ l	100 μ l
8. Incubate	Cover plate and incubate one hour at RT on an orbital shaker					
9. Wash	Wash all wells five times					
10. Add reagents	Developer Solution	100 μ l	100 μ l	100 μ l	100 μ l	100 μ l
11. Incubate	30 minutes					
12. Add reagents	Stop Solution	100 μ l	100 μ l	100 μ l	100 μ l	100 μ l
13. Read	Read plate at wavelength of 450 nm					

Table 3. Quick Protocol Guide

ANALYSIS

Performance Characteristics

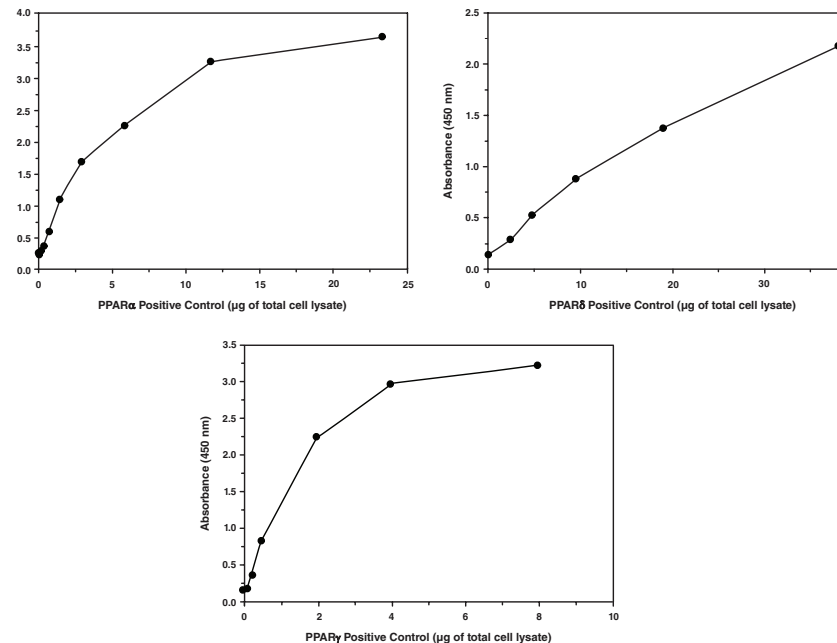


Figure 3. Schematic of the Transcription Factor Binding Assay
 Increasing amounts of positive control (total lysate) are assayed for PPAR α , δ , and γ DNA-binding activity using the Cayman's PPAR α , δ , γ Complete Transcription Factor Assay Kit.

Interferences

The following reagents were tested for interference in the assay.

Reagent	Will Interfere (Yes or No)
EGTA (≤ 1 mM)	No
EDTA (≤ 0.5 mM)	No
ZnCl (any concentration)	Yes
DTT (between 1 and 5 mM)	No
Dimethylsulfoxide ($\leq 1.5\%$)	No

Troubleshooting

Problem	Possible Causes	Recommended Solutions
No signal or weak signal in all wells	<ul style="list-style-type: none"> A. Omission of key reagent B. Plate reader settings not correct C. Reagent/reagents expired D. Salt concentrations affected binding between DNA and protein E. Developing reagent used cold F. Developing reagent not added to correct volume 	<ul style="list-style-type: none"> A. Check that all reagents have been added and in the correct order; perform the assay using the Positive Control B. Check wavelength setting on plate reader and change to 450 nm C. Check expiration date on reagents D. Reduce the amount of nuclear extract used in the assay, or reduce the amount of salt in the nuclear extracts (alternatively can perform buffer exchange) E. Prewarm the Developing Solution to room temperature prior to use F. Check pipettes to ensure correct amount of Developing Solution was added to wells
High signal in all wells	<ul style="list-style-type: none"> A. Incorrect dilution of antibody (too high) B. Improper/inadequate washing of wells C. Over-developing 	<ul style="list-style-type: none"> A. Check antibody dilutions and use amounts outlined in instructions B. Follow the protocol for washing wells using the correct number of times and volumes C. Decrease the incubation time when using the developing reagent
High background (NSB)	Incorrect dilution of antibody (too high)	Check antibody dilutions and use amounts outlined in the instructions

Problem cont.	Possible Causes cont.	Recommended Solutions cont.
Weak signal in sample wells	<ul style="list-style-type: none"> A. Sample concentration is too low B. Incorrect dilution of antibody C. Salt concentrations affecting binding between DNA and protein 	<ul style="list-style-type: none"> A. Increase the amount of nuclear extract used; loss of signal can occur with multiple freeze/thaw cycles of the sample; prepare fresh nuclear extracts and aliquot as outlined in product insert B. Check antibody dilutions and use amounts outlined in the instructions C. Reduce the amount of nuclear extract used in the assay or reduce the amount of salt in the nuclear extracts (alternatively can perform buffer exchange)

References

1. Desvergne, B. and Wahli, W. Peroxisome proliferator-activated receptors: Nuclear control of metabolism. *Endocr. Rev.* **20**, 649-688 (1999).
2. Clark, R.B. The role of PPARs in inflammation and immunity. *J. Leukoc. Biol.* **71**, 388-400 (2002).
3. Olefsky, J.M. Nuclear receptor minireview series. *J. Biol. Chem.* **276(40)**, 36863-36864 (2001).

4. Duplus, E., Glorian, M., and Forest, C. Fatty acid regulation of gene transcription. *J. Biol. Chem.* **275(40)**, 30749-30752 (2000).
5. Gervois, P., Torra, I.P., Fruchart, J.-C., *et al.* Regulation of lipid and lipoprotein metabolism by PPAR activators. *Clin. Chem. Lab. Med.* **38(1)**, 3-11 (2000).
6. Brown, P.J., Smith-Oliver, T.A., Charifson, P.S., *et al.* Identification of peroxisome proliferator-activated receptor ligands from a biased chemical library. *Chemistry & Biology* **4**, 909-918 (1997).

12								
11								
10								
9								
8								
7								
6								
5								
4								
3								
2								
1								
	A	B	C	D	E	F	G	H

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.

This document is copyrighted. All rights are reserved. This document may not, in whole or part, be copied, photocopied, reproduced, translated, or reduced to any electronic medium or machine-readable form without prior consent, in writing, from Cayman Chemical Company.

©12/14/2020, Cayman Chemical Company, Ann Arbor, MI, All rights reserved. Printed in U.S.A.

