



6042 Cornerstone Court W, Ste B
San Diego, CA 92121
Tel: 1.858.829.3082
Fax: 1.858.481.8694
Email: info@bpsbioscience.com

Data Sheet

Acetyl-Coenzyme A Carboxylase 2 (ACC2) Assay Kit

Cat # 79282

DESCRIPTION: Acetyl-Coenzyme A Carboxylase (ACC) plays an important role in fatty acid metabolism so it has been proposed to be a drug target for the fatty acid-related metabolic diseases including obesity and diabetes. The *Acetyl-Coenzyme A Carboxylase 2 (ACC2) assay kit* is designed to measure ACC2 activity for screening and profiling applications using ADP-Glo™ reagents as a detection reagent. The *Acetyl-coenzyme A carboxylase 2 (ACC2) assay kit* comes in a convenient 96-well format, with enough purified recombinant ACC2 enzyme, ACC2 substrate, ATP and ACC assay buffer for 100 enzyme reactions.

COMPONENTS:

Catalog #	Reagent	Amount	Storage	
50101	ACC2	5 µg	-80°C	Avoid multiple freeze/thaw cycles!
79283	5x ACC assay buffer	1 ml	-20°C	
	ATP (100 µM)	500 µl	-20°C	
	Acetyl-CoA (2 mM)	25 µl	-20°C	
	Sodium Bicarbonate (400 mM)	75 µl	-20°C.	
	96-well plate, white	1	Room Temp.	

MATERIALS OR INSTRUMENTS REQUIRED BUT NOT SUPPLIED:

ADP-Glo™ Kinase Assay (Promega #V6930)
Microplate reader capable of reading luminescence
Adjustable micropipettor and sterile tips

APPLICATIONS: Useful for studying enzyme kinetics and screening small molecular inhibitors for drug discovery and HTS applications.

STABILITY: Up to 6 months when stored as recommended.

REFERENCE:

Ting, L. *J. Cell. Biochem.* **99(6)**:1476-1488 (2006)

OUR PRODUCTS ARE FOR RESEARCH USE ONLY. NOT FOR DIAGNOSTIC OR THERAPEUTIC USE.

To place your order, please contact us by Phone **1.858.829.3082** Fax **1.858.481.8694**

Or you can Email us at: info@bpsbioscience.com

Please visit our website at: www.bpsbioscience.com



6042 Cornerstone Court W, Ste B
San Diego, CA 92121
Tel: 1.858.829.3082
Fax: 1.858.481.8694
Email: info@bpsbioscience.com

ASSAY PROTOCOL:

All samples and controls should be tested in duplicate.

- 1) Thaw 5x ACC assay buffer, ATP and Acetyl-CoA and Sodium Bicarbonate.
- 2) Prepare the master mixture (15 μ l per well): N wells x (3.5 μ l 5x ACC assay buffer + 5 μ l ATP (100 μ M) + 0.25 μ l Acetyl-CoA + 0.75 μ l Sodium Bicarbonate + 5.5 μ l water). Add 15 μ l to every well except the Blank well.
- 3) For the Blank well, prepare 'no substrate master' (15 μ l per well): N wells x (3.5 μ l 5x ACC assay buffer + 5 μ l ATP (100 μ M) + 6.5 μ l water). Add 15 μ l to the Blank well. (Alternatively, 'master mix + no ACC2' can be used as a Blank well)

	Positive Control	Test Inhibitor	Blank (no substrate)	Blank (no ACC2)
5x ACC assay buffer	3.5 μ l	3.5 μ l	3.5 μ l	3.5 μ l
ATP (100 μ M)	5 μ l	5 μ l	5 μ l	5 μ l
Acetyl-CoA (2 mM)	0.25 μ l	0.25 μ l	–	0.25 μ l
Sodium Bicarbonate (400 mM)	0.75 μ l	0.75 μ l	–	0.75 μ l
Water	5.5 μ l	5.5 μ l	6.5 μ l	5.5 μ l
Test Inhibitor	–	2.5 μ l	–	–
Inhibitor Diluent (no inhibitor)	2.5 μ l	–	2.5 μ l	2.5 μ l
1x ACC buffer	–	–	–	7.5 μ l
ACC2 (5.4 ng/ μ l)	7.5 μ l	7.5 μ l	7.5 μ l	–
Total	25 μ l	25 μ l	25 μ l	25 μ l

- 4) Add 2.5 μ l of Inhibitor solution of each well labeled as "Test Inhibitor". For the "Positive Control" and "Blank", add 2.5 μ l of the same solution without inhibitor (Inhibitor Diluent; e.g. 10% DMSO (aqueous) is recommended for inhibitor diluent, resulting in 1% DMSO at final concentration)
 - 5) Prepare 2 ml of 1x ACC assay buffer by mixing 400 μ l of 5x ACC assay buffer with 1600 μ l water. 2 ml of 1x Kinase assay buffer is sufficient for 100 reactions.
 - 6) To the wells designated as "Blank (no ACC2)", add 7.5 μ l of 1x ACC assay buffer.
 - 7) Thaw ACC2 enzyme on ice. Upon first thaw, briefly spin tube containing enzyme to recover full content of the tube. Calculate the amount of ACC2 required for the
- OUR PRODUCTS ARE FOR RESEARCH USE ONLY. NOT FOR DIAGNOSTIC OR THERAPEUTIC USE.

To place your order, please contact us by Phone **1.858.829.3082** Fax **1.858.481.8694**

Or you can Email us at: info@bpsbioscience.com

Please visit our website at: www.bpsbioscience.com

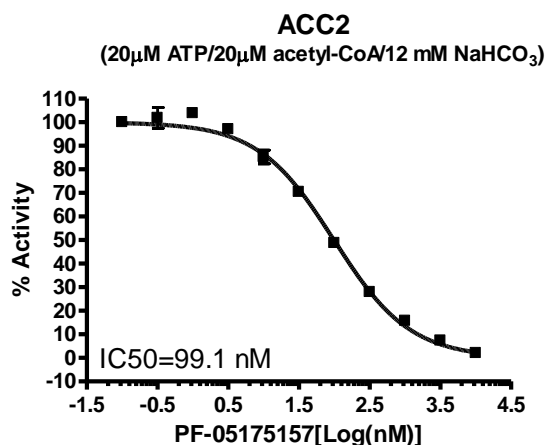


6042 Cornerstone Court W, Ste B
San Diego, CA 92121
Tel: 1.858.829.3082
Fax: 1.858.481.8694
Email: info@bpsbioscience.com

assay and dilute enzyme to ~5.4 ng/μl with 1x ACC assay buffer. Store remaining undiluted enzyme in aliquots at -80°C. *Note: ACC2 enzyme is sensitive to freeze/thaw cycles. Avoid multiple freeze/thaw cycles. Do not re-use thawed aliquots or diluted enzyme.*

- 8) Initiate reaction by adding 7.5 μl of diluted ACC2 enzyme to the wells designated "Positive Control", "Test Inhibitor Control" and "Blank (no substrate)". Incubate at room temperature for 40 minutes.
- 9) Thaw ADP-Glo reagent (Promega).
- 10) After the 40 minutes reaction, add 20 μl of ADP-Glo reagent to each well. Cover plate and incubate at room temperature for 45 minutes.
- 11) Thaw Kinase Detection reagent (Promega).
- 12) After 45 min incubation with ADP-Glo reagent, add 50 μl of Kinase-Detection reagent to each well. Cover plate with aluminum foil and incubate at room temperature for 45 ~ 90 minutes
- 13) Measure luminescence using the microplate reader.

Example of Assay Results:



Inhibition of ACC2 enzyme by PF-05175157, measured using the *Acetyl-Coenzyme A Carboxylase 2 (ACC2) assay kit* (Cat. #xxxxxx). *Data shown is lot-specific. For lot-specific information, please contact BPS Bioscience, Inc. at info@bpsbioscience.com*

OUR PRODUCTS ARE FOR RESEARCH USE ONLY. NOT FOR DIAGNOSTIC OR THERAPEUTIC USE.

To place your order, please contact us by Phone **1.858.829.3082** Fax **1.858.481.8694**

Or you can Email us at: info@bpsbioscience.com

Please visit our website at: www.bpsbioscience.com



6042 Cornerstone Court W, Ste B
San Diego, CA 92121
Tel: 1.858.829.3082
Fax: 1.858.481.8694
Email: info@bpsbioscience.com

RELATED PRODUCTS:

<u>Product Name</u>	<u>Catalog #</u>	<u>Size</u>
ACC2, His-tag	50201	10 µg
ACC2, His-tag, Strep-tag	50203	10 µg
ACC1, His-tag	50200	10 µg
ACC1, FLAG-His-tags	50202	10 µg

OUR PRODUCTS ARE FOR RESEARCH USE ONLY. NOT FOR DIAGNOSTIC OR THERAPEUTIC USE.

To place your order, please contact us by Phone **1.858.829.3082** Fax **1.858.481.8694**

Or you can Email us at: info@bpsbioscience.com

Please visit our website at: www.bpsbioscience.com