

# Acetyl CoA Assay

A plate based fluorometric assay to quantify Acetyl CoA levels in biological samples.

Acetyl CoA is a key molecule in cellular metabolism and energy production. A central hub, connecting different metabolic pathways within the cell. Acetyl CoA plays a pivotal role in the citric acid cycle resulting in the production of energy-rich molecules such as ATP and is a precursor for the synthesis of fatty acids, cholesterol, and other important cellular components. AkrivisBio's Acetyl CoA Assay is a simple, sensitive method of measuring Acetyl CoA levels in a variety of samples with a detection range of 0 - 400 picomoles. In the assay, Acetyl CoA in samples is hydrolyzed to CoA and acetate. CoA is then converted to succinyl CoA forming NADH, which is utilized by diaphorase to reduce resazurin to resorufin with an intense fluorescence measured at Ex/Em = 535/587 nm.

## Ordering Information

[Order Online »](#)

AKR-MA-0127	100 wells
-------------	-----------

## Manuals, SDS & CofA

[View Online »](#)

## Handling & Storage

Long Term Storage      -20°C

Shipping                      Blue Ice

**Regulatory Status**      RUO - Research Use Only

## Product Details

Application                      Fluorescent detection

Species Reactivity              Species independent

**Technical Info / Product Notes**      For the Original Manufacturer's data sheet please [click here](#).



ENZO LIFE SCIENCES,  
INC.  
Phone: 800.942.0430  
[info-usa@enzolifesciences.com](mailto:info-usa@enzolifesciences.com)

European Sales Office  
ENZO LIFE SCIENCES  
(ELS) AG  
Phone: +41 61 926 8989  
[info-eu@enzolifesciences.com](mailto:info-eu@enzolifesciences.com)

Belgium, The Netherlands  
& Luxembourg  
Phone: +32 3 466 0420  
[info-be@enzolifesciences.com](mailto:info-be@enzolifesciences.com)

France  
Phone: +33 472 440 655  
[info-fr@enzolifesciences.com](mailto:info-fr@enzolifesciences.com)

Germany  
Phone: +49 7621 5500 526  
[info-de@enzolifesciences.com](mailto:info-de@enzolifesciences.com)

UK & Ireland  
Phone (UK customers):  
0845 601 1488  
Phone: +44 1392 825900  
[info-uk@enzolifesciences.com](mailto:info-uk@enzolifesciences.com)