## Acetyltransferase activity kit

The Acetyltransferase activity kit is a homogeneous mix-and-read fluorescent assay for the determination of any acetyl-CoA dependent acetyltransferase activity. It is suitable for end-point or kinetic read options, which is ideal for determining mechanism of action, kinetics, and screening candidate compounds. The assay is amendable to HTS and miniaturization.

Acetylation of proteins and drugs is a regulatory event that can result in the activation, inactivation, or detoxification of molecules including histones, kinases, transcription factors, anti-microbial agents, and neurotransmitters. While acetylation can be carried out by a variety of acetyltransferase enzymes using diverse substrates, their common cofactor or acetyl donor is acetyl-CoA.

Citations: 3

**View Online »** 

**Ordering Information** 

Order Online »

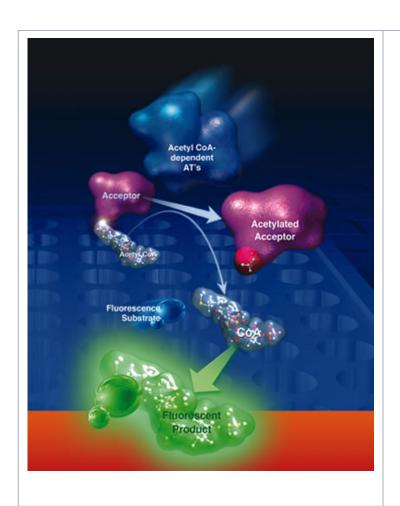
ADI-907-026

96 wells

Manuals, SDS & CofA

View Online »

- Universal suitable for any AcetylCoA-dependent acetyltransferase
- Robust Z' factor = 0.87
- Sensitive fluorescent detection allows for both DNA and protein assays
- Higher throughput homogeneous mix-and-read assay



## **Handling & Storage**

**Use/Stability** Store all components at -20°, except positive control at -80°C.

Shipping Dry Ice

Regulatory Status RUO - Research Use Only

**Product Details** 

**Application** Activity assay, Fluorescent detection, HTS

**Application Notes** For the measurement of acetyl-CoA dependent

acetyltransferase activity in purified in vitro samples that

do not contain reductants from any species.

**Contents** Black Microtiter plate, Reaction buffer, Assay buffer,

Detection solution, Positive control

Species Reactivity Species independent

Last modified: May 29, 2024



eu@enzolifesciences.com