

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 »](#)

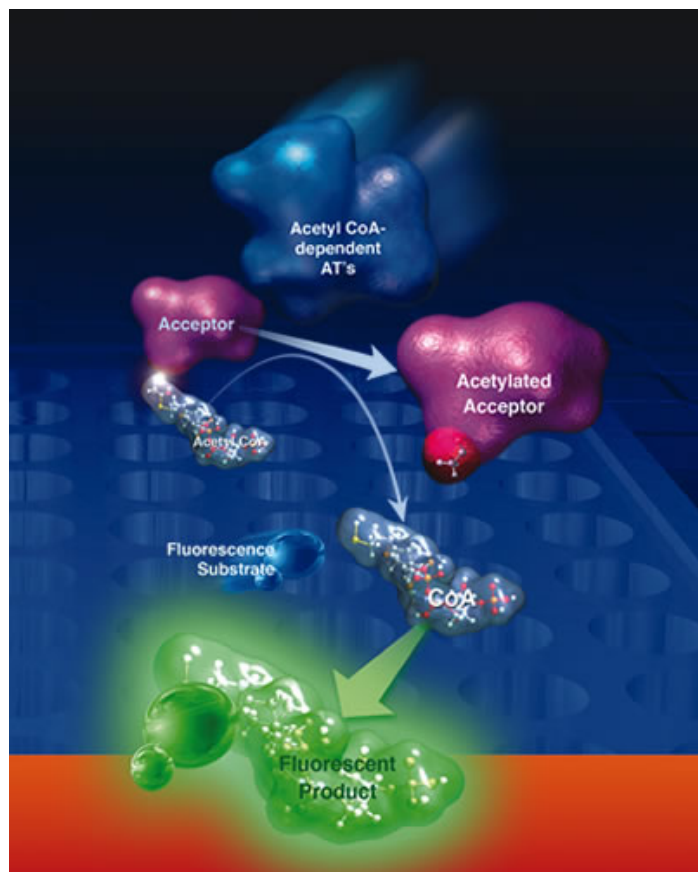
ADI-907-026

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



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)