

# p300 Inhibitor Screening Assay

A plate based fluorometric assay to screen potential p300 inhibitors in a rapid and high-throughput adaptable way.

p300, also known as KAT3B, is a histone acetyltransferase. One of the primary functions of p300 is its role in epigenetic regulation. By acetylating histones, p300 alters the structure of chromatin, promoting transcriptional activation and accessibility of genes. This process allows the cells to respond dynamically to external cues and ensure appropriate gene expression during development, tissue homeostasis, and cellular response to stress and environmental changes. Dysregulation of p300 activity has been implicated in various diseases, including cancer, neurodegenerative disorders, and cardiovascular diseases, making it an attractive target for therapeutic interventions. AkrisBio's p300 Inhibitor Screening Assay is a simple method of assessing potential p300 inhibitors in a rapid and high-throughput adaptable assay.

## Ordering Information

[Order Online »](#)

AKR-MA-0146	100 wells
-------------	-----------

## Manuals, SDS & CofA

[View Online »](#)

## Handling & Storage

Long Term Storage -80°C

Shipping Dry 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)