

Pyruvate Assay

A plate based colorimetric/fluorometric assay to quantitate pyruvate in the range of 1 μ M - 10 mM in a variety of biological samples.

Pyruvate is a crucial molecule in cellular metabolism. It also serves as a key precursor for various biosynthetic pathways, including the synthesis of amino acids and fatty acids. Pyruvate's central role in energy production and biosynthesis makes it a fundamental molecule in cellular processes. Measuring pyruvate levels can provide insights into various pathological conditions, including metabolic disorders, liver diseases, cancer etc.

AkrivisBio's Pyruvate Assay is a simple, sensitive method of quantitating pyruvate in a variety of biological sample types. In the assay, Pyruvate in the sample is oxidized by pyruvate oxidase with the formation of hydrogen peroxide. Hydrogen peroxide is then utilized by the peroxidase to oxidize ADHP forming an intense color or fluorescence measured at OD = 570 nm or at Ex/Em = 538/587 nm respectively.

Ordering Information

[Order Online »](#)

| | |
|-------------|-----------|
| AKR-MA-0129 | 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 Colorimetric 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

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

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

France
Phone: +33 472 440 655
info-fr@enzolifesciences.com

Germany
Phone: +49 7621 5500 526
info-de@enzolifesciences.com

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