

α -Ketoglutarate Assay

A plate based colorimetric/fluorometric assay to measure the concentration of α -Ketoglutarate in a wide range of sample types.

α -Ketoglutarate is a key intermediate in multiple pathways, playing essential roles in energy production, amino acid metabolism, and cellular homeostasis. In the TCA cycle, α -ketoglutarate is converted into succinyl-CoA, producing ATP. α -Ketoglutarate is a precursor for the synthesis of several amino acids, including glutamate and glutamine which are critical for neurotransmission, immune function, and cellular signaling. α -Ketoglutarate also acts as a signaling molecule, influencing gene expression and processes such as oxidative stress response and protein synthesis. AkrisBio's α -Ketoglutarate Assay is a simple, sensitive way to measure this important molecule in a wide range of sample types, with a useful measurement range from 10 nmoles down to below 10 pmoles per well. In the assay, α -ketoglutarate is transaminated with the generation of pyruvate, which is oxidized by pyruvate oxidase forming hydrogen peroxide. Hydrogen peroxide is then utilized by the peroxidase to oxidize ADHP resulting in the formation of intense color or fluorescence measured at OD 570 nm or at Ex/Em = 535/587 nm respectively.

Ordering Information

[Order Online »](#)

AKR-MA-0141	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