

Total Carbohydrate Assay

A plate based colorimetric assay to determine total carbohydrate content in various samples.

Carbohydrates are one of the fundamental building blocks of biochemistry. The total carbohydrate assay is a crucial analytical technique used to determine overall carbohydrate content in each sample. This assay provides valuable information for a wide range of applications in the food and beverage industries, pharmaceutical research, and basic biochemistry. The assay involves the oxidation of sugars to aldehydes which react with phenol to give color around 490 nm. Scientists utilize the assay to assess the nutritional value of food products, verify the efficiency of carbohydrate-based pharmaceutical formulations, and monitor carbohydrate metabolism in biological samples. AkrisBio's Total Carbohydrate Assay's accuracy and reliability make it an invaluable tool in scientific research and commercial industries, contributing to improved product development and a deeper understanding of carbohydrate-related phenomena. Moreover, it serves as an essential tool in quality control processes, enabling manufacturers to ensure consistency and adherence to standards.

Ordering Information

[Order Online »](#)

AKR-MA-0144	100 wells
-------------	-----------

Manuals, SDS & CofA

[View Online »](#)

Handling & Storage

Long Term Storage Ambient

Shipping Ambient Temperature

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](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)