

# Choline/Acetylcholine Assay

**A plate based colorimetric/fluorometric assay to measure choline in a variety of biological samples with a sensitivity from the low picomole to the multi-nanomole range.**

Choline plays a vital role in various metabolic processes. As a precursor to acetylcholine, it's crucial for maintaining healthy neurological function. Choline is involved in lipid metabolism both aiding in the transport and metabolism of fats in the liver and as a precursor to phosphatidylcholine, a major component of membranes. Choline is also highly important as a methyl donor, contributing to the formation of S-adenosylmethionine, a molecule with diverse functions. Small amounts of choline can be synthesized but dietary intake is important to maintain the amounts required for optimal health. AkrisBio's Choline Assay is a simple sensitive method of measuring choline in a variety of biological samples with a sensitivity from the low picomole to the multi-nanomole range.

## Ordering Information

[Order Online »](#)

|             |           |
|-------------|-----------|
| AKR-MA-0152 | 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](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)