

# Cholesterol- Cholesterol Ester Assay

## A plate based colorimetric/fluorometric assay to quantify free, esterified or total

**cholesterol in biological samples.**  
Cholesterol is a lipid molecule essential to all animal life. It is the precursor of steroid hormones and vitamin D. It makes up about 30% of cell membranes. It is transported via lipoproteins including HDL, LDL, VLDL in the blood, as free cholesterol and as cholesteryl esters. AkrivisBio's Cholesterol/Cholesteryl Ester Assay provides a simple, sensitive method of quantifying free, esterified or total cholesterol (cholesterol and cholesteryl esters). In this assay, cholesteryl esters are hydrolyzed to cholesterol by cholesterol esterase, which is then oxidized by cholesterol oxidase to yield hydrogen peroxide. Hydrogen peroxide is utilized by the peroxidase to oxidize ADHP thereby 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-0119	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)