

HDL-LDL Cholesterol Assay

A plate based colorimetric/fluorometric assay to quantify HDL and LDL/VLDL cholesterol in samples separately.

HDL and LDL play a role in various pathological conditions. Low levels of HDL combined with high level of LDL are associated with an increased risk of cardiovascular events including plaque rupture, stroke, hypertension, etc. AkrisBio's HDL/LDL Assay provides a simple, sensitive method for quantifying HDL and LDL/VLDL cholesterol separately. In this assay, cholesterol oxidase oxidizes free cholesterol with the formation of hydrogen peroxide. Hydrogen peroxide is then 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 = 538/587 nm respectively. In an optional step, cholesterol esterase cleaves cholesterol esters to free cholesterol, permitting the assessment of free and total cholesterol separately. Thus, if cholesterol esterase is included in the assay, total cholesterol is measured. If it is not included, free cholesterol is measured. The amount of cholesterol ester can be calculated by subtracting free cholesterol from total cholesterol.

Ordering Information

[Order Online »](#)

AKR-MA-0124	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