

Total Cholesterol Assay Kit (Colorimetric)

Cell Biolabs' Total Cholesterol Assay Kits measure the total cholesterol within serum, plasma, cell lysate, or tissue samples. The assays will detect total cholesterol (cholesteryl esters plus free cholesterol) in the presence of cholesterol esterase or only free cholesterol in the absence of the esterase enzyme.

Ordering Information

[Order Online »](#)

CBI-STA-384	192 assays
-------------	------------

Manuals, SDS & CofA

[View Online »](#)

- Suitable for use with serum, plasma, lysate, or tissue samples
- Cholesterol standard included

Handling & Storage

Use/Stability Upon receipt, store the Cholesterol Standard, 50X Colorimetric Probe, HRP, Cholesterol Oxidase, and Cholesterol Esterase at -20°C. The 50X Colorimetric Probe is light sensitive and must be stored accordingly. Avoid multiple freeze/thaw cycles. Store the remaining kit components at 4°C.

Shipping Ambient Temperature

Regulatory Status RUO - Research Use Only

Product Details

Application Colorimetric detection

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