

EN

Cholesterol Food Std

REF 984391

3 x 3 ml

INTENDED USE

Standard for determination of Cholesterol in homogenous liquid samples using automated Thermo Scientific™ Arena™ or Gallery™ analyzer. The Cholesterol standard is a stable aqueous solution.

STORAGE AND STABILITY

The standard is stable up to the end of the indicated month of expiry, if stored unopened at 2...8 °C and contamination is avoided.

WARNINGS AND PRECAUTIONS

Standard is hazardous.

See separate sheet inside the kit for Hazardous- and Precautions-phrases: H315, H317, H319, P280, P302 + P352, P305 + P351 + P338.

Standard contains sodium azide as preservative.

Exercise the normal precautions required for handling all laboratory reagents.

The product has to be disposed of as laboratory chemical in accordance with local regulations.

STANDARD PREPARATION

Standard concentration is high for the application and therefore should be diluted manually before use with isopropanol. Use 1 part of Standard and 4 parts of isopropanol.

PROCEDURE

The assay procedure is indicated in the package insert of the respective test-kit.

Notes for calibration are indicated in the applications for automated systems.

COMPONENTS AND CONCENTRATIONS

Lot.	See the CoA/Value sheet	
Parameter	Concentration	Unit
Cholesterol	See the CoA/Value sheet	mg/l

Concentrations are given in the CoA/Value sheet of the product. The concentrations were measured with Thermo Scientific reagents.

Differences to the nominal values could be caused by using other reagents and methods.

WASTE MANAGEMENT

Please refer to local legal requirements. It is recommended to empty the analyzer cuvette waste bin and waste water daily. Emptying should be done immediately after the analysis when using hazardous reagents/solutions.

Note: If using reagents/solutions that react with each other, cuvette waste bin and waste water should be emptied and washed between use of these reagents.

ADDITIONAL MATERIAL

Certificate of analysis and SDS are available at

www.e-labeling.eu/TSF

Applications for Gallery and Arena automated analyzers are available upon request from the local sales representative. Information in the Application note can change without prior notice.

MANUFACTURER

Thermo Fisher Scientific Oy
Ratastie 2, P.O. Box 100, FI-01621 Vantaa, Finland
Tel. +358 10 329200

CONTACT INFORMATION

www.thermoscientific.com

e-mail: system.support.fi@thermofisher.com

Date of revision (yyyy-mm-dd)

2015-05-22

Changes from previous version

Warnings and precautions updated.