

EN

Lactose standard

REF 984390

3 x 3 ml

INTENDED USE

Standard for determination of Lactose in in homogenous liquid samples using automated Thermo Scientific™ Arena™ or Gallery™ analyzer. The Lactose 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

The standard contains sodium azide (< 0.1 %) as preservative. Do not swallow. Avoid contact with skin and mucous membranes. Take the necessary precautions for the use of laboratory reagents.

STANDARD PREPARATION

The standard is ready-to-use.

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
Lactose	See the CoA/Value sheet	g/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)
2015-02-03

Changes from previous version

Components and concentrations updated.
General update.