



# Galactosemia

## Quantase™ Neonatal Total Galactose

### Newborn Screeners Are Talking

*“In my opinion, quality is the most important feature of any product.”*

### Bio-Rad Is Listening

Quantase™ legacy in newborn screening assays demonstrate our commitment to quality. We manufacture in compliance to international standards to provide the performance and reproducibility that you demand. Bio-Rad's commitment to quality gives confidence in your results.

### Galactosemia Assay

Bio-Rad Quantase™ offers a robust total Galactose newborn screening test providing quality results you can have confidence in.

### The Bio-Rad Assay

- Not subject to the interferences commonly observed with fluorescent and bacterial inhibition methods
- Simplified protocol can be used on most plate readers
- Rapid results in under 2 hours

**BIO-RAD**

## Galactosemia Assays

### Clinical Background

Galactosemia is an inherited disorder of galactose metabolism, caused by a deficiency in one of three enzymes, the most important being galactose-1-phosphate uridyl transferase. Galactosemia may also be caused by a lack of galactokinase or UDP-galactose-4-epimerase activity.

Long-term consequences of galactosemia include mental retardation, developmental abnormalities, and even death.

The worldwide prevalence of galactosemia is estimated to be 1:60,000-1:80,000. The Quantase™ Neonatal Total Galactose Screening Assay is an enzymatic colorimetric end-point method for the determination of total D(+) galactose in dried-blood spot specimens.

### Ordering Information

Catalog No.	Description	
532-6054	Quantase™ Neonatal Total Galactose Assay .....	500 test kit
532-6050	Quantase™ Neonatal Total Galactose Assay .....	2000 test kit
532-5038	Quantase™ Accessory Kit 3, filter plates .....	2000 test kit
532-5039	Quantase™ Accessory Kit 4, assay plates.....	2000 test kit
532-5040	Quantase™ Accessory Kit 5, elution buffer.....	2000 test kit
532-5025	Vacuum Manifold System.....	1 unit



**Bio-Rad  
Laboratories**

*For further information, please contact the Bio-Rad office nearest you or visit our website at [www.bio-rad.com/diagnostics](http://www.bio-rad.com/diagnostics)*

*Clinical  
Diagnostics Group*

**Website** [www.bio-rad.com/diagnostics](http://www.bio-rad.com/diagnostics) **U.S.** 1-800-224-6723 **Australia** 61-2-9914-2800 **Austria** 43-1-877-8901 **Belgium** +32 (0)710-53-00 **Brazil** +55 (31)3689-6600  
**Canada** 1-514-334-4372 **China** 86-21-61698500 **Czech Republic** 420-241-430-532 **Denmark** +45-4452-1000 **Finland** 358-9-804-22-00 **France** 33-1-47-95-60-00  
**Germany** +49 (0)89-318-840 **Greece** 30-210-7774396 **Hong Kong** 852-2789-3300 **Hungary** +36-1-459-6100 **India** 1800-180-1224 **Israel** 972-3-9636050 **Italy** +39-02-216091  
**Japan** 81-3-6361-7070 **Korea** 82-2-3473-4460 **Mexico** +52 (55)5488-7670 **The Netherlands** +31-318-540686 **New Zealand** 64-9-415-2280 **Norway** +47-23-38-41-30  
**Poland** 48-22-3319999 **Portugal** 351-21-472-7700 **Russia** +7-495-721-1404 **Singapore** 65-6415-3170 **South Africa** 27-11-442-85-08 **Spain** 34-91-590-5200  
**Sweden** +46-8-555-127-00 **Switzerland** +41 (0)26-674-55-05/06 **Taiwan** 886-2-2578-7189 **Thailand** 662-651-8311 **United Kingdom** +44 (0)20-8328-2000