



Galactosemia

Quantase™ Neonatal Total Galactose and Neonatal GALT Assays

Newborn Screeners Are Talking

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

Bio-Rad Is Listening

Quantase™ newborn screening assays demonstrate our commitment to quality. We manufacture under the most stringent standards to provide the performance and reproducibility that you demand. Bio-Rad’s quality gives you confidence in your results.

Galactosemia Assay

Bio-Rad Quantase™ offers a complete package for galactosemia testing. Whether screening with Total Galactose or with GALT (transferase), Bio-Rad provides results you can have confidence in.

The Bio-Rad Assay

- Not subject to the interferences commonly observed with fluorescent and bacterial inhibition methods
- Simplified protocols can be used on most plate readers
- Rapid results in under 2 hours (Total Galactose) and 3.5 hours (GALT)

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.

The Quantase™ Neonatal GALT Assay measures galactose-1-phosphate uridyl transferase activity in neonatal dried-blood spots to screen for galactosemia.

Ordering Information

Catalog No.	Description	
532-6000	Quantase™ Neonatal GALT Assay.....	192 test kit
532-6001	Quantase™ Neonatal GALT Assay.....	960 test kit
532-6002	Quantase™ Neonatal GALT Assay Controls, bi-level.....	1 unit
532-6054	Quantase™ Neonatal Total Galactose Assay	500 test kit
532-6050	Quantase™ Neonatal Total Galactose Assay	2000 test kit
532-5036	Quantase™ Accessory Kit 1, elution buffer, assay plates, filter plates	2000 test kit
532-5037	Quantase™ Accessory Kit 2, elution buffer, assay plates	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

*Clinical
Diagnostics Group*

Website www.bio-rad.com/diagnostics **U.S.** 1-800-2BIO-RAD **Australia** 61-2-9914-2800 **Austria** 43-1-877-8901 **Belgium** 32-9-385-5511 **Brazil** 5521-3237-9400 **Canada** 1-514-334-4372 **China** 86-21-64260808 **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** 1-800-180-1224 **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** 7-495-721-14-04 **Singapore** 65-6415-3170 **South Africa** 27-11-442-85-08 **Spain** 34-91-590-5200 **Sweden** 46-8-555-127-00 **Switzerland** 41-61-717-95-55 **Thailand** 662-651-8311 **United Kingdom** +44-(0)20-8328-2000