



G6PD Deficiency

Quantase™ Neonatal G6PD* Assay

Newborn Screeners Are Talking

“Simpler is better.”

Bio-Rad Is Listening

Quantase™ newborn screening assays feature streamlined protocols that are readily adapted to any laboratory environment. Whether you are looking for manual or automated solutions, Quantase™ assays provide the efficiency you demand. When you are ready to simplify your workflow, choose Bio-Rad.

Quantase™ Neonatal G6PD Assay

Bio-Rad introduces a convenient and precise assay for the screening of G6PD deficiency. The Quantase™ G6PD Assay utilizes a rapid colorimetric procedure designed specifically for dried-blood spot screening, free of a red blood cell lysis step.

* Not available in the US

The Bio-Rad Assay

- Simplified protocol for use on most microplate systems
- Designed for use with dried-blood spots
- Established ranges for newborn screening
- Room temperature procedure incubator not required
- Numerical results eliminate subjective interpretations
- Rapid results in about 45 minutes

BIO-RAD

G6PD Deficiency

Clinical Background

Glucose-6-phosphate dehydrogenase, an enzyme found in most body cells, is part of the pentose phosphate pathway that metabolizes glucose. G6PD deficiency is a hereditary, sex-linked condition carried on the X chromosome, with clinical disease found mostly in males. G6PD deficiency impairs the stability of red cell membranes and makes red cells susceptible to destruction by strong oxidizing agents. Affected patients are at risk for hemolytic anemia, which can be induced by exposure to certain drugs and foods.

G6PD is one of the most common hereditary enzyme disorders, affecting over 200 million people worldwide. About 10% of all African American males inherit mild forms of G6PD deficiency. Some people of Mediterranean origin inherit severe forms of the disease. The ingestion of fava beans may also elicit hemolytic episodes in some affected Caucasians.

Ordering Information

Catalog No.	Description
532-6100	Quantase™ Neonatal G6PD Assay 200 test kit
532-6101	Quantase™ Neonatal G6PD Assay 500 test kit
532-6102	Quantase™ Neonatal G6PD Assay 2000 test kit



**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 91-124-4029300 Israel 972-3-9636050 Italy +39-02-216091 Japan 81-3-6361-7070 Korea 82-2-3473-4460 Mexico 52(55)5200-0520 The Netherlands +31-318-540666 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 65-6415-3188 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