

HUMAN HEALTH

ENVIRONMENTAL HEALTH

RELIABLE
RESULTS
– EFFICIENTLY



GSP™ NEONATAL ASSAY
for congenital hypothyroidism screening


PerkinElmer®
For the Better



GSP™ NEONATAL KITS



Results you can rely on

GSP™ is the new, automated neonatal screening system from PerkinElmer. It has multi-technology capability, which means that both DELFIA® and prompt fluorescence assays can now be run on a common platform. Its versatility and speed help it to accommodate present and future screening needs.

The first kit, already available for the GSP instrument is GSP Neonatal TSH, for congenital hypothyroidism (CH) screening. The new kit employs the DELFIA chemistry, acclaimed for the analytical sensitivity and robustness of assays based on it.

In addition, several new kits are currently in development as we work to fill the GSP assay portfolio.

Reduced risk of error

All contents of GSP Neonatal kits, including reagents, plates, QC material and lot specific QC certificates are barcoded to reduce the risk of errors.

Less work – faster results

To minimize the pre-analytical work all reagents are ready to use. On the instrument itself, there are direct access water and waste lines, with automatic dilution of wash concentrate. The reagent storage compartment is temperature controlled to provide extended on board stability, and relieve the need to unload reagents once loaded.

For further time efficiency, a single calibration curve can be used for up to 24 hours. This enables more samples to be run on plates and easier reruns. Additionally, all calibrators and controls come in cassette format to support automatic punching.



SCREENING FOR CONGENITAL HYPOTHYROIDISM

using the GSP Neonatal TSH assay

Severe congenital hypothyroidism (CH) can be found in 1 in 4000 newborns

while milder forms of the disease are more frequent. CH results from a failure of the thyroid glands to produce thyroid hormones in adequate amounts. The condition can easily be treated with daily doses of thyroid hormones but clinical diagnosis is difficult to establish and the disease may continue unrecognized for a long time causing irreversible brain damage. However, increased thyroid stimulating hormone (TSH) and decreased thyroxine are clear signs of CH, which have led to the establishment of large scale screening programs.

The PerkinElmer GSP Neonatal TSH assay is used with dried blood spot specimens as an aid in screening newborns for congenital hypothyroidism.

- The incubation time is only 3.5 h
- Sensitive and robust DELFIA assay
- The kit contains reagents and plates for 1152 tests (12 plates)



Work with the leader in Newborn Screening

PerkinElmer is the global market leader in neonatal screening, currently serving customers in some 74 countries.

The company is a total solution provider offering complete systems based on a broad range of high quality, validated products, including newborn screening kits, consumables, instruments and software. Our global presence and comprehensive support philosophy mean that our expertise is available to you at all times.

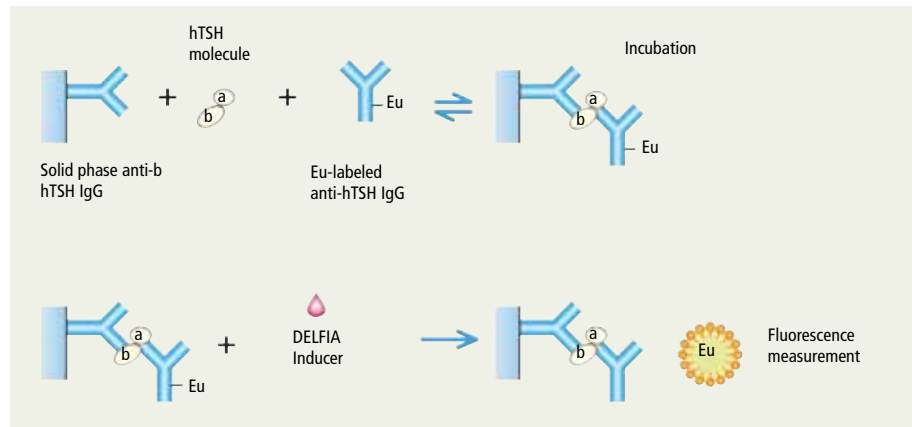
125,000 babies saved so far

The first DELFIA TSH kit was developed in 1985, to allow dried blood spot measurement of TSH from neonatal samples. In 2008, for this analyte alone PerkinElmer sold kits in 64 countries.

By 2008, almost 300 million babies had been screened with PerkinElmer products. More than 125,000 of them were identified as having a disorder.

GSP Neonatal TSH assay

The GSP Neonatal TSH assay is based on a direct sandwich technique where two monoclonal antibodies recognize separate antigenic determinants on the TSH molecule. The fluorescence signal is proportional to the analyte concentration in the sample.



ORDERING INFORMATION

3301-001U	GSP Neonatal hTSH kit
3304-0010	DELFLIA Inducer
4080-0010	GSP Wash concentrate

All PerkinElmer neonatal products may not be available in all countries.

For the better

At PerkinElmer, we're working to improve the health and safety of people and their environment. From safer water to cleaner air to healthier babies, our solutions touch every part of your life.

For information on availability of PerkinElmer diagnostic products please contact your local representative.

PerkinElmer, Inc.
940 Winter Street
Waltham, MA 02451 USA
Phone: (800) 762-4000 or
(+1) 203-925-4602
www.perkinelmer.com

PerkinElmer Genetic Screening
Center of Excellence
Wallac Oy, PO Box 10
20101 Turku, Finland
Phone: + 358 2 2678 111
Fax: + 358 2 2678 357



For a complete listing of our global offices, visit www.perkinelmer.com/ContactUs

©2009 PerkinElmer, Inc. All rights reserved. PerkinElmer is a registered trademark of PerkinElmer, Inc. All other trademarks depicted are the property of their respective holders or owners. PerkinElmer reserves the right to change this document at any time and disclaims liability for editorial, pictorial or typographical errors. All PerkinElmer diagnostic products may not be available in all countries. For information on availability please contact your local representative.