

Cystic fibrosis screening



PerkinElmer's Neonatal IRT assay for its DELFIA® and AutoDELFLIA® platforms allows easy-to-perform and accurate cystic fibrosis (CF) screening to support every customer's needs. It is globally the most widely used assay for 1st tier CF screening, and in 2008 either the manual or automatic version of the assay was used in 34 countries.

Top performance based on robust DELFIA® chemistry

The Neonatal IRT assay provides high analytical sensitivity with the limit of blank of 0.43 ng/mL for DELFIA and 0.53 ng/mL for AutoDELFLIA platforms. Intra-assay imprecision for the clinically relevant area is close to 9% for both assays. All calibrators and controls are lot-specific to ensure that every lot is of the same level.

Fully automated AutoDELFLIA® system leads the way in terms of results quality

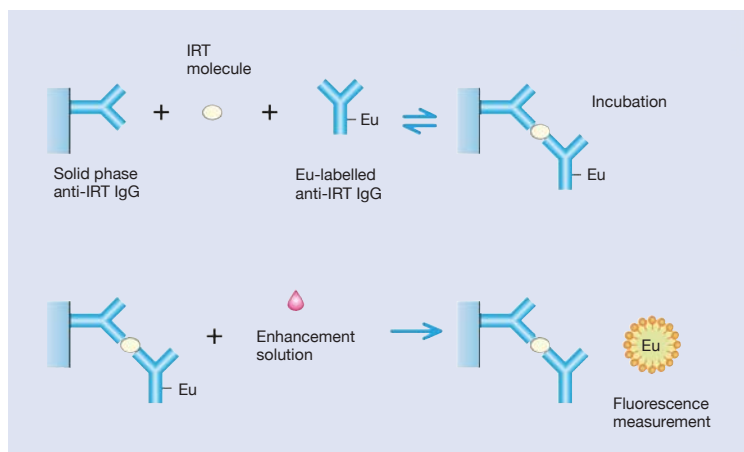
The automatic system allows for the best assay performance in terms of precision, and brings robustness, reliability and cost-effectiveness to your CF screening program.

Early diagnosis for better quality of life

Cystic fibrosis is a common genetic disorder affecting 1 in 2,500 Caucasians. The diagnosis is often based on the symptoms which may cause considerable delays in the disease intervention, and evidence indicates that early attention may be important in determining the clinical outcome. The amount of the pancreatic enzyme, immunoreactive trypsin (IRT) has been shown to be increased in blood of CF patients especially during the first weeks after birth.

PerkinElmer's Neonatal IRT assay allows the quantitative determination of IRT from dried blood spot specimens, and is intended as an aid in screening for CF.

Robust DELFIA® “sandwich-type” assay



The assay is based on the direct sandwich technique in which two monoclonal antibodies bind to different epitopes on the target molecule. The enhanced signal is measured with high analytical sensitivity using time-resolved fluorometry.

Ordering information

- A005-204 DELFIA® Neonatal IRT kit (4 plates)
- A005-210 DELFIA® Neonatal IRT kit (10 plates)
- B005-204 AutoDELFLIA® Neonatal IRT kit (4 plates)
- B005-212 AutoDELFLIA® Neonatal IRT kit (12 plates)

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

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, DELFIA and AutoDELFLIA are registered trademarks of PerkinElmer, Inc. All 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.