



HUMAN HEALTH

ENVIRONMENTAL HEALTH

CONTROLS OF DEMONSTRATED STABILITY FOR YOUR HB ASSAYS

HB610 extendSURE™ HbFASC Control

In the detection of hemoglobin variants using gel electrophoresis, immunoassay or other technologies, HB610 extendSURE™ HbFASC controls provide the stability needed for reliable assay performance monitoring.

Covering four key Hb variants

HB610 extendSURE™ HbFASC controls represent an optimal control solution for laboratories performing hemoglobinopathy screening or diagnostic testing. The product comprises controls for the key hemoglobin variants, HbF, HbA, HbS and HbC.

Compatible with most detection technologies

HB610 extendSURE HbFASC controls are ideal for laboratories using PerkinElmer's RESOLVE® and AutoDELFIA® platforms. They are also well-suited for use with other platforms, and with other technologies, such as HPLC.

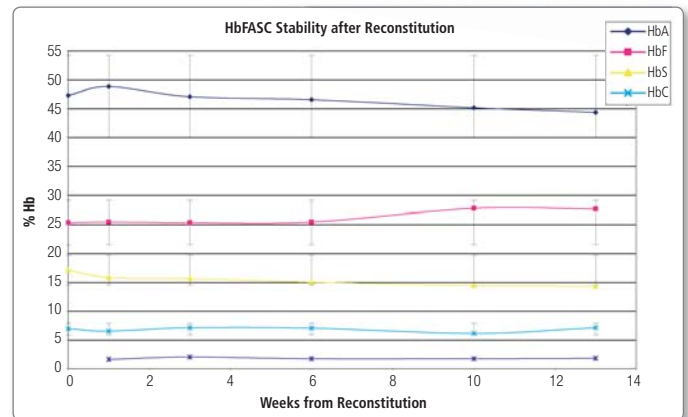
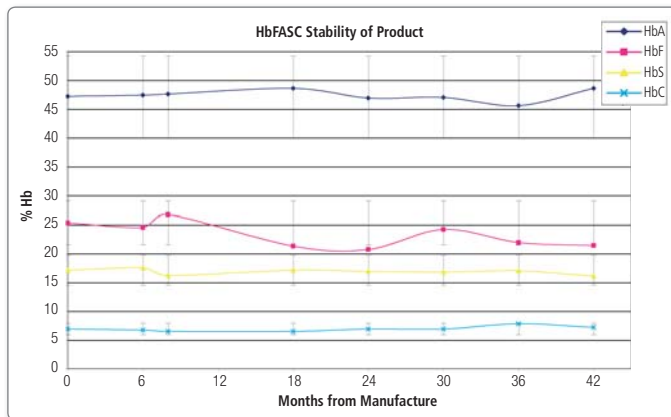
- Optimal controls for use with RESOLVE® or AutoDELFIA®
- Ideal for HPLC, capillary IEF, cellulose acetate and agar/agarose gels, and immunoassays
- CE-marked and FDA cleared
- Very good stability after reconstitution - six weeks shelf life

Convenient and cost-effective

Four individual vials of lyophilized preparation each reconstitute to 1.0mL of control material in which all four variants are present. After reconstitution, the controls have a stated shelf life of 6 weeks when stored at 2-8°C. This makes them particularly cost effective and easy to use. No aliquoting and freezing is required and wastage is minimized. Unopened, the controls have a lifetime of 30 months from the date of manufacture.

Kit contains:

- 4 vials of control
- 2 vials of reconstitution buffer
- Information and instruction sheet in eight languages (Brazilian Portuguese, Danish, English, French, German, Italian, Spanish, Swedish)



The graph on the left shows results for the four Hb variants assayed after reconstitution at intervals over three and a half years. The graph on the right shows results over a period of 13 weeks following reconstitution. In both graphs the individual points' error bars represent the stability specification of $\pm 15\%$ of the initial concentration. The graphs provide strong evidence of the product's excellent stability.

ORDERING INFORMATION

Part Number	Product Description
HB610	extendSURE™ HbFASC Control
B019-104	B019-104 AutoDELFLIA® Hb Immunoassay
FR-9120	RESOLVE® Hemoglobin kit (135 tests)
FR-9400	RESOLVE Hemoglobin kit (360-440 tests) – recommended for neonatal use
FR-9360	RESOLVE Hemoglobin kit (3600-4400 tests) – recommended for neonatal use

Manufacturer: Canterbury Scientific Ltd.

PerkinElmer is a global distributor of the product for all applications, and the exclusive supplier for IEF use.

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, AutoDELFLIA and RESOLVE are registered trademarks of PerkinElmer, Inc. ExtendSURE is a trademark of Canterbury Scientific Ltd. 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.