

HUMAN HEALTH

ENVIRONMENTAL HEALTH

24 CHROMOSOMES IN 24 HOURS

KARYOLITE™ BoBs™*

Part of the BACs-on-Beads™ product family

* For Research Use Only. Not for Use in Diagnostic Procedures.


PerkinElmer®
For the Better

WHAT IS KARYOLITE™ BoBs™?

KaryoLite™ BoBs™ is a new member of the BACs-on-Beads™ family of products.

KaryoLite BoBs has been developed to detect aneuploidies, gains and losses in all 24 chromosomes in a single assay. This assay is being offered for research use only. The product covers both p and q arms of all chromosomes 1-22, X and Y. It is a bead-based molecular karyotyping assay utilizing

probes derived from BAC DNA and coupled to encoded microspheres. It consists of 90 DNA probes immobilized onto polystyrene microspheres distinguishable by the Luminex® instrument system.



The kit is easy to use with objective and fast result interpretation. The protocol is very robust, assuring results that can be relied on. The kit contains sample labeling reagents, the BACs-on-Beads mix, hybridization reagents, streptavidin-phycoerythrin reporter and wash reagents.

Research may be performed with KaryoLite BoBs using minute amounts of genomic DNA extracted directly from amniotic fluid, chorionic villae or fetal tissue. In fact only 50 ng of DNA is needed, and no cell culturing is required.

For True Rapid Aneuploidy Detection

- **Results in less than 24 hours**

Complete procedure from sample to result takes less than 24 hours, allowing your laboratory to obtain results the following day.

- **Quick and easy implementation**

The KaryoLite BoBs reagents are provided in a single kit.

- **Cost-efficiency**

Tens of samples can be run simultaneously reducing the hands-on time required.

- **Easy interpretation**

Intuitive BoBsoft™ analysis software presents results that are clear and easy to interpret.

- **More information**

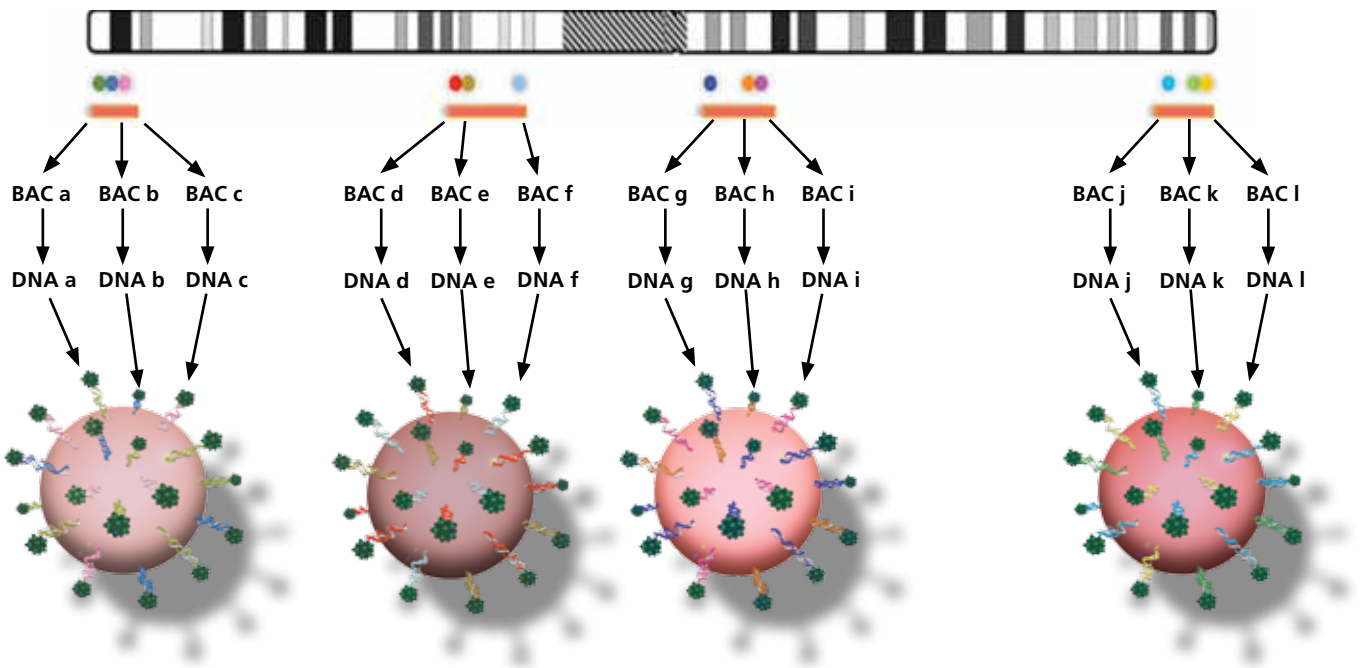
Aneuploidies, gains and losses in all 24 chromosomes in one assay.



Results are obtained in less than 24 hours and the hands-on-time is only about 3.5 hours.

The probes used have been carefully selected to provide information about the whole chromosome including the subtelomere region of the chromosome. Separate bead types provide coverage of at least two separate regions of each arm of the chromosome. In acrocentric chromosomes (13, 14, 15, 21 and 22) three beads cover the q arm of the chromosome.

KARYOLITE™ BoBs™



KaryoLite BoBs utilizes a new concept of composite beads having three different BAC clones on each bead type. The composite clone format expands the region of chromosomal DNA interrogated by each bead. This will result in an averaging of the signal thus minimizing the effect of allelic dropouts or bias in sample labeling.

ORDERING INFORMATION

Product	Product Name	Comment
4501-0010	KaryoLite™ BoBs™	Research Use Only, Available in May 2011
4500-0010	Constitutional BoBs™	Research Use Only
3100-0010	Prenatal BoBs™	Not available in the USA
1014-0020	Luminex® 200 w/ xPONENT®	
5012-0010	BoBsoft™ Analysis Software	Research Use Only
5012-0020	BoBsoft™ Analysis Software	Not available in the USA
4502-0010	EasyAmp™ WGA kit	Research Use Only

All PerkinElmer products may not be available in all countries.
For information on availability 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, Inc.
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

©2009-2011 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.