



With all the changes
in nylon fashion,
our style has stayed
the same.

**Genomic Tested Zeta-Probe® GT nylon membranes.
No formula change, no compromise.**

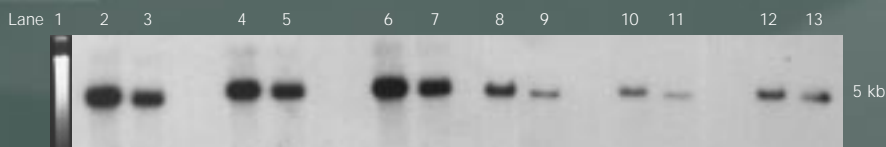
Low background and high signal-to-noise ratios are critical for nucleic acid blotting. That's why Zeta-Probe GT membranes have the same formula they've always had, with the same quality you expect from Bio-Rad. We don't follow the changing trends; we set them and keep our standards high. Zeta-Probe GT membranes are positively charged nylon with quaternary amine groups for maximum binding and retention. A high tensile strength makes them resistant to tearing and shrinking. Most importantly, each lot is tested for picogram level detection of a single copy gene on a genomic Southern and alkaline transfer blot. For assurance, a lot-specific result certificate is always enclosed.

If your old nylon membranes just don't seem the same, they probably aren't. Call us when you're ready to change into something a little more comfortable — Zeta-Probe GT. For more information, contact your Bio-Rad representative.

BIO-RAD

Zeta-Probe GT Recommended Uses

Capillary, vacuum, or electrophoretic transfer of single and double-stranded DNA and RNA of all sizes for subsequent hybridization. Alkaline blotting for single step binding of nucleic acids.



Superior Lot Results: *Bgl* I digested human genomic DNA was transferred to Zeta-Probe GT by Southern or alkaline method. Membranes were hybridized with ³²P labeled plasmid containing a 1.8 kb fragment from a human factor VIII cDNA. Probe hybridizes to a 5 kb *Bgl* I restriction fragment. Blots were exposed to x-ray film overnight with intensifying screens at -70 °C.

Lane 1: ethidium bromide stain of genomic DNA digest; alternating 5 µg and 2.5 µg human genomic DNA digested with *Bgl* I; lanes 2 and 3 – Southern control; lanes 4 and 5 – Southern test of side 1; lanes 6 and 7 – Southern test of side 2; lanes 8 and 9 – alkaline control; lanes 10 and 11 – alkaline test of side 1; lanes 12 and 13 – alkaline test of side 2.

Ordering Information

Catalog #	Product Description
162-0190	9 x 12 cm, 15 sheets per box
162-0191	10 x 15 cm, 15 sheets per box
162-0192	15 x 15 cm, 15 sheets per box
162-0193	15 x 20 cm, 15 sheets per box
162-0194	20 x 20 cm, 15 sheets per box
162-0195	20 x 25 cm, 15 sheets per box
162-0196	30 cm x 3.3 m, 1 roll
162-0197	20 cm x 3.3 m, 1 roll

BIO-RAD

Molecular
Bioscience Group

U.S. (800) 4BIORAD • California (510) 741-1000 • Australia 02-9914-2800 • Austria (1)-877 89 01 • Belgium 09-385 55 11 • Canada (905) 712-2771 • China (01) 2046622 • Denmark 39 17 9947 • Finland 90 804 2200 • France (1) 49 60 68 34 • Germany 089 318 84-0 • India 91-11-461-0103 • Italy 02-21609 1 • Japan 03-5811-6270 • Hong Kong 7893300 • The Netherlands 0318-540666 • New Zealand 09-443 3099 • Singapore (65) 272-9877 • Spain (91) 661 70 85 • Sweden 46 (0) 8-735 83 00 • Switzerland 01-809 55 55 • United Kingdom 0800 181134

SIG 020996 Printed in USA

With all the changes
in nylon fashion,
our style has stayed the same.

**Genomic Tested Zeta-Probe® GT nylon membranes.
No formula change, no compromise.**

Low background and high signal-to-noise ratios are critical for nucleic acid blotting. That's why Zeta-Probe GT membranes have the same formula they've always had, with the same quality you expect from Bio-Rad. We don't follow the changing trends; we set them and keep our standards high. Zeta-Probe GT membranes are positively charged nylon with quaternary amine groups for maximum binding and retention. A high tensile strength makes them resistant to tearing and shrinking. Most importantly, each lot is tested for picogram level detection of a single copy gene on a genomic Southern and alkaline transfer blot. For assurance, a lot-specific result certificate is always enclosed.

If your old nylon membranes just don't seem the same, they probably aren't. Call us when you're ready to change into something a little more comfortable — Zeta-Probe GT. For more information, contact your Bio-Rad representative.

For a Free Sample, Contact:

BIO-RAD