



226PR

G-Biosciences ♦ 1-800-628-7730 ♦ 1-314-991-6034 ♦ technical@GBiosciences.com

A Geno Technology, Inc. (USA) brand name

Hyb-LINK™

Hybridization Buffer

(Cat. # 786-405, 786-406)



think proteins! think G-Biosciences www.GBiosciences.com

INTRODUCTION

Hyb-LINK™ is a unique high performance proprietary hybridization buffer for Southern, Northern blots or Southwestern blots. It provides researchers with unprecedented simplicity, reliability, accuracy for hybridization experiments. Hyb-LINK™ is suitable for all types of probes and protocol, which involves incubation of blots with denatured probes in Hyb-LINK™ followed by washing of the blot to remove excess unhybridized probes. It also provides high resolution and sensitivity and is suitable for Southern/Northern or Southwestern hybridization analysis, hybridization with oligonucleotides or other types of probes, and for hybridization screening of gene libraries.

ITEM(S) SUPPLIED

Description	Cat. # 786-405	Cat. # 786-406
Hyb-LINK™	125ml	4 x 125ml

STORAGE CONDITIONS

Shipped at ambient temperature. Upon arrival, store Hyb-LINK™ at room temperature. Stable for 1 year, when stored and handled properly. To avoid possible microbial contamination, use aseptic techniques when handling.

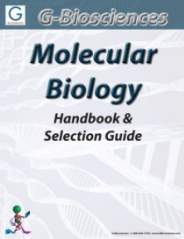
NOTE: Reagents must not be stored in the cold (4°C).

PROTOCOL

1. Incubate the blot in a plastic bag with appropriate volume (0.1ml/cm²) of Hyb-LINK™ for 1 hour at 65°C.
2. Add fresh (0.1ml/cm²) volume of Hyb-LINK™ to the bag with the denatured probe and incubate overnight at 65°C.
3. After incubation, wash the blot with SSC or sodium phosphate buffer to remove excess probe.
4. Proceed further with your experiment.

RELATED PRODUCTS

Download our Molecular Biology Handbook.



<http://info.gbiosciences.com/complete-molecular-biology-handbook>

For other related products, visit our website at www.GBiosciences.com or contact us.

Last saved: 8/21/2012 CMH



www.GBiosciences.com