

Amersham ECL gold hybridization buffer

Product Specification Sheet

Introduction

Product code

RPN3006

About

For use with Horseradish Peroxidase-labelled probes, using the ECL™ direct labelling and detection systems. Sufficient for hybridization of 4000 cm² membrane.

Important

Read these instructions carefully before using the products.

Intended use

The products are intended for research use only, and shall not be used in any clinical or *in vitro* procedures for diagnostic purposes.

Safety

For use and handling of the products in a safe way, refer to the Safety Data Sheets.

Storage

Store at 15°C to 25°C. The buffer and blocking agent are stable at 15°C to 25°C for at least twelve months when stored under the recommended conditions.

Expiry

See outer packaging.

Components

Hybridization buffer: 2 x 500 ml

Blocking reagent: 2 x 25 g

Quality control

The ECL gold hybridization buffer is tested by our quality control group using the ECL direct system (RPN3000/3001), to ensure detection of a single copy gene in human genomic DNA blots, representing a minimum sensitivity of 0.5 pg target in 1 µg genomic DNA.

Preparation of buffer for hybridization

Step	Action
1	Take the required amount of hybridization buffer and add solid Sodium Chloride (analytical grade) to a concentration that is most suitable for effective probe hybridization.

Note:

The optimum concentration may be different for different probes, 0.5 M NaCl generally gives acceptable results.

Step	Action
2	Add the blocking agent to a final concentration of 5% (w/v). Immediately mix thoroughly to get the blocking agent into a fine suspension.
3	Continue mixing at room temperature for 1 hour on a magnetic stirrer or roller mixer, then preheat to 42°C for 0.5–1 hour.

The buffer is then ready for use. (A few small particles of undissolved blocking agent will not affect the hybridization).

If excess buffer is prepared this should be dispensed into suitable aliquots into sterile plastic containers and stored at -15°C to -30°C. Prepared buffer may be stored for at least three months. Do not store in a frost-free freezer.

Hybridization

Hybridization should be carried out as described in the protocol booklet supplied with the ECL direct systems.

Note: *Do not exceed a temperature of 42°C during hybridizations.*

Related products

ECL direct nucleic acid labelling and detection systems	
Labelling reagent for 5 µg DNA Hybridization buffer and detection reagents sufficient for 2000 cm ² membrane	RPN3000
Labelling reagents for 10 µg DNA Hybridization buffer and detection reagents sufficient for 4000 cm ² membrane	RPN3001
ECL direct nucleic acid labelling system	
Labelling reagents for 5 µg DNA Hybridization buffer sufficient for 2000 cm ² membrane	RPN3005
ECL detection reagents	
Sufficient for 2000 cm ² membrane	RPN3004
Sufficient for 4000 cm ² membrane	RPN2105
Hybond™-N+ and Hybond-ECL	

cytiva.com

Cytiva and the Drop logo are trademarks of Global Life Sciences IP Holdco LLC or an affiliate.

Amersham™, ECL and Hybond are trademarks of Global Life Sciences Solutions USA LLC or an affiliate doing business as Cytiva.

All other third-party trademarks are the property of their respective owners.

© 2020–2021 Cytiva

All goods and services are sold subject to the terms and conditions of sale of the supplying company operating within the Cytiva business. A copy of those terms and conditions is available on request. Contact your local Cytiva representative for the most current information.

For local office contact information, visit [cytiva.com/contact](https://www.cytiva.com/contact)

29460175 AA V:5 01/2021

