

GelBond PAG film

Introduction

Product codes

80112936

80112937

Important

Read these instructions carefully before using the products.

Intended use

GelBond PAG film is used as a support for moulding polyacrylamide gels.

Storage

GelBond PAG film is sensitive to light and elevated temperatures. Store in resealable light-proof package at room temperature, and at a dry dark location.

Contents

80112936 Pack of 50 films, 124 x 258 mm.

80112937 Pack of 50 films, 203 x 260 mm.

Instructions

Each film has one hydrophilic side and one hydrophobic side. The gel must be cast on the hydrophilic side, which is the side protected by a paper sheet. The two sides can also be distinguished by noting that a small drop of water will "spread" on the hydrophilic surface whereas it will "bead up" on the hydrophobic one. GelBond PAG film can be used for gels containing sodium dodecyl sulfate or urea. Nonionic detergents such as Nonidet P-40 have been found to interfere with the adhesion of polyacrylamide gels to GelBond PAG film.

Certain basic ampholytes (pH \geq 8.0) have been found to interfere with the adhesion of polyacrylamide gel to GelBond PAG film. To overcome this problem, increase the concentration of ammonium persulfate to 0.07% and TEMED to 0.1% in the final acrylamide gel solution. For isoelectric focusing the recommended carrier ampholyte concentrations are 2.5% w/v (Ampholine) or dilution 1:16 (Pharmalyte™).

Release the polyacrylamide gel from the opposing glass plate can be facilitated by rinsing the glass with Repel-Silane ES, Product code 17133201. Note that all glass plates must be meticulously cleaned before use.

Prefocus the gel to remove the ammonium persulfate before applying the samples.

Page intentionally left blank



cytiva.com

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

Pharmalyte is a trademark 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)

29645873 AA V:5 02/2021