

## **Nitrocellulose Membrane**

Catalog nos. LC2000, LC2001

## Instructions for Use

Description	Nitrocellulose membrane is a high quality membrane ideal for blotting of proteins and nucleic acids. The Nitrocellulose membrane is available in two pore sizes:				
	Catalog no. P	ore Size	Application		
	LC2000 0.	.2 μm	Transfer of low molecular weight proteins (<20 kDa) and nucleic acids (<300 bp)		
	LC2001 0.	.45 µm	Transfer of most proteins (>20 kDa) and nucleic acids (>300 bp)		
	The major features of the Nitrocellulose membrane available from Invitrogen are:				
	Composed of 100% pure nitrocellulose to provide high-quality transfer				
	Contains no support fabric or detergents				
	• Compatible with commonly used transfer conditions and detection methods such as staining, immunodetection, fluorescence, or radiolabeling				
	Provides high-sensitivity with low background				
	• Supplied in a pre-cut, pre-assembled membrane/filter paper sandwich for convenience				
Specifications	The specifications of Nitrocellulose membrane are listed below.				
	Quantity:		0 membrane/filter paper sandwiches (sufficient for 20 transfers)		
	Pore Size:	0	.2 μm (LC2000); 0.45 μm (LC2001)		
	Dimensions:	8	.3 cm x 7.3 cm		
	Filter Media:		100% pure nitrocellulose		
			209 μg/cm² of proteins (0.2 μm) 30 μg/cm² of proteins (0.45 μm)		
	Binding Interaction	n: F	Hydrophobic and electrostatic		
	Re-probe Characte	ristics: N	Jo		
Applications	Nitrocellulose mer	nbrane is s	uitable for the following applications:		
	Western, Southern, and Northern Transfers				
	Amino Acid Analysis				
	Dot/Slot Blotting				

Continued on next page

Part No: IM1600

Rev. Date: 032403

This product is distributed for laboratory research use only. CAUTION: Not for diagnostic use. The safety and efficacy of this product in diagnostic or other clinical uses has not been established.

For technical questions about this product, call the Invitrogen TECH-LINE<sup>SM</sup> 800 955 6288

## Instructions for Use, Continued

	Product	Quantity	Catalog no.
	Novex <sup>®</sup> Tris-Glycine Transfer Buffer (25X)	500 ml	LC3675
	NuPAGE <sup>®</sup> Transfer Buffer (20X)	1 L	NP0006-1
	XCell II <sup>™</sup> Blot Module	1 unit	EI9051
	WesternBreeze® Chromogenic Kit-Anti-Mouse	1 kit	WB7103
	WesternBreeze® Chromogenic Kit-Anti-Rabbit	1 kit	WB7105
	WesternBreeze® Chromogenic Kit-Anti-Goat	1 kit	WB7107
	WesternBreeze® Chemiluminescent Kit-Anti-Mouse	1 kit	WB7104
	WesternBreeze <sup>®</sup> Chemiluminescent Kit-Anti-Rabbit	1 kit	WB7106
	WesternBreeze® Chemiluminescent Kit-Anti-Goat	1 kit	WB7108
	MagicMark <sup>™</sup> Western Protein Standard	250 µl	LC5600
	PVDF Membrane (0.2 μm pore size)	20 membrane/filter papers	LC2002
	Invitrolon <sup>™</sup> PVDF Membrane (0.45 µm pore size)	20 membrane / filter papers	LC2005

goal is to ensure that every customer is 100% satisfied with our products and our services. Our goal is to ensure that every customer is 100% satisfied with our products and our service. If you should have any questions or concerns about an Invitrogen product or service, contact our Technical Service Representatives. Invitrogen warrants that all of its products will perform according to specifications stated on the certificate of analysis. The company will replace, free of charge, any product that does not meet those specifications. This warranty limits Invitrogen Corporation's liability only to the cost of the product. No warranty is granted for products beyond their listed expiration date. No warranty is applicable unless all product components are stored in accordance with instructions. Invitrogen reserves the right to select the method(s) used to analyze a product unless Invitrogen agrees to a specified method in writing prior to acceptance of the order. Invitrogen makes every effort to ensure the accuracy of its publications, but realizes that the occasional typographical or other error is inevitable. Therefore Invitrogen makes no warranty of any kind regarding the contents of any publications, please report it to our Technical Service Representatives.

Invitrogen assumes no responsibility or liability for any special, incidental, indirect or consequential loss or damage whatsoever. The above limited warranty is sole and exclusive. No other warranty is made, whether expressed or implied, including any warranty of merchantability or fitness for a particular purpose.

©1997-2003 Invitrogen Corporation. All rights reserved.