

Glogos® II Autorad Markers

Catalog #420201, #420202, #420203



For Research Use Only. Not for use in diagnostic procedures.

Storage Store Glogos® II Autorad Markers at room temperature away from direct sunlight.

INTRODUCTION

Stratagene's Glogos® II autorad markers are luminescent peel-off stickers used for accurately marking and aligning autoradiographs. Easily affixed to plastic wrap or paper, these convenient markers emit low-level luminescence rather than radioactivity, making the Glogos II autorad markers a safe, nonhazardous alternative to radioactive ink. The asymmetric design of the Glogos II autorad markers ensures accurate orientation of an autoradiograph to its membrane. To assist in the alignment of a blot or membrane to its autoradiograph, a ruler is included with the markers. The Glogos II autorad markers are adaptable for long or short exposures, and their usage range is from -70°C to room temperature.

MATERIALS PROVIDED

Material Provided	Catalog #420201	Catalog #420202	Catalog #420203
Glogos® II Autorad Markers	100 markers	300 markers	1000 markers

INSTRUCTIONS

After probing, washing, and wrapping the membrane, peel off a Glogos II autorad marker and affix the marker to the wrapping in a place where the marker will not cover the data. Expose and develop the film as usual and then align the image on the film with the Glogos II autorad marker to achieve correct orientation.

Extent of Light Activation

The Glogos II autorad markers are calibrated, so exposure to ambient levels of room light for at least 1 minute will provide sufficient activation for autoradiograph exposures lasting longer than ~45 minutes. Because most light decay occurs within the first hour, a Glogos II autorad marker exposed for 1 hour or for several days will appear similar.

If exposing data to film for a short period of time (e.g., 15 minutes) or for quick exposures, hold the Glogos II autorad marker close to a bright light for just a few seconds to activate the marker.

Decay

Because luminescence decays rapidly and to avoid light-colored images, put the Glogos II autorad marker on film within 3 minutes after placing the marker in the dark following light activation.

Exposure to Film

Place the Glogos II autorad marker directly facing the film. If superactivated with *bright* light, minimize the movement of the Glogos II autorad marker once the marker is adjacent to the film to avoid double exposures. Sharper images are obtained when the marker is affixed to clear areas of the wrapping, rather than against a background of white, reflective material (e.g., Whatman® 3MM paper). If *extreme* accuracy in autoradiograph alignment is required and/or if data is large in size, use two Glogos II autorad markers, placing one on each side of the longest diagonal.

Temperature

The Glogos II autorad markers are effective for exposures ranging from room temperature to -70°C .

Intensifying Screen

The use of an intensifying screen will increase the intensity of the image by two- to threefold, but will *slightly* decrease the resolution of the image.

LIMITED PRODUCT WARRANTY

Stratagene warrants to the original purchaser that this product will be free from defects in workmanship or materials for a period of one (1) year from the date of purchase, as indicated on the original invoice, under normal use.

The sole and exclusive remedy under this limited warranty is repair or replacement of defective products or parts thereof. Repair or replacement products or parts thereof will be furnished solely on an exchange basis and are obtainable only by the original purchaser. The original purchaser shall return the defective product, or part thereof, properly packaged, postage or shipping costs prepaid to Stratagene. Loss or damage during shipment shall be at the risk of the purchaser.

This limited warranty does not include repair or replacement necessitated by accident, neglect, misuse, relocation, unauthorized repair, or modification of the product.

THE ABOVE LIMITED WARRANTY IS THE SOLE WARRANTY PROVIDED BY STRATAGENE. NO OTHER WARRANTIES, EXPRESS OR IMPLIED, INCLUDING WARRANTIES OF MERCHANTABILITY OR FITNESS FOR A PARTICULAR PURPOSE, ARE PROVIDED BY STRATAGENE.

In no event will Stratagene be liable for any direct, indirect, consequential, or incidental damages, including lost profits, or for any claim by any third party, arising out of the use, the results of use, or the inability to use this product.

ORDERING INFORMATION AND TECHNICAL SERVICES

United States and Canada

Stratagene

11011 North Torrey Pines Road
La Jolla, CA 92037

Telephone (858) 373-6300

Order Toll Free (800) 424-5444

Technical Services (800) 894-1304

Internet tech_services@stratagene.com

World Wide Web www.stratagene.com

Stratagene European Contacts

Location	Telephone	Fax	Technical Services
Austria	0800 292 499	0800 292 496	0800 292 498
Belgium	00800 7000 7000	00800 7001 7001	00800 7400 7400
	0800 15775	0800 15740	0800 15720
France	00800 7000 7000	00800 7001 7001	00800 7400 7400
	0800 919 288	0800 919 287	0800 919 289
Germany	00800 7000 7000	00800 7001 7001	00800 7400 7400
	0800 182 8232	0800 182 8231	0800 182 8234
Netherlands	00800 7000 7000	00800 7001 7001	00800 7400 7400
	0800 023 0446	+31 (0)20 312 5700	0800 023 0448
Switzerland	00800 7000 7000	00800 7001 7001	00800 7400 7400
	0800 563 080	0800 563 082	0800 563 081
United Kingdom	00800 7000 7000	00800 7001 7001	00800 7400 7400
	0800 917 3282	0800 917 3283	0800 917 3281

All Other Countries

Please contact your local distributor. A complete list of distributors is available at www.stratagene.com.

ENDNOTES

Glogos® is a registered trademark of Stratagene in the United States.

Whatman® is a registered trademark of Whatman Ltd.