



Molecular Analyst 2-D PAGE Software

Software Instruction Manual and Tutorial

Catalog Number
170-7560

For Technical Service
Call Your Local Bio-Rad Office or
in the U.S. Call 1-800-4BIORAD
(1-800-424-6723)

Information contained in this document is subject to change without notice and does not represent a commitment on the part of Bio-Rad Laboratories. The software described in this document is furnished under a licence agreement. The software may be used and copied only in accordance with the terms of the agreement. It is against the law to copy the software on any medium except as specifically allowed in the licence. No part of this manual may be reproduced or transmitted in any form or by any means, electronic or mechanical, including photocopying and recording, for any purpose without the express written permission of Bio-Rad Laboratories.

Microsoft, MS, and MS-DOS are registered trademarks.
Windows is a trade mark of the Microsoft Corporation.

All other registered trademarks acknowledged.

Licence Agreement

This Licence Agreement sets out the terms and conditions relating to the acquisition and use by you of the Molecular Analyst 2D package ("the Product"). The Licensee acknowledges that these terms form part of the contract of sale for the product and the purchase by the Licensee is evidence of his acceptance and agreement of these terms.

When the Product is purchased by you the disks contained in the Product become your property. All software on the disks (including the Molecular Analyst 2D application and supporting files), "the Software", shall remain the sole property of Bio-Rad Laboratories, the Licensor, and you, the Licensee, are permitted to use the Software only in accordance with the terms of this Licence Agreement.

These terms are as follows:

1. The Licensee may make one backup copy on floppy disk of the Software on each of the original disks for use by the Licensee only. No other backup copies may be made.
2. The Software must not be installed or used on more than one computer workstation at any one time.
3. The Software and any documentation or other material supplied in the Product shall not be adapted, modified, decompiled or disassembled in any way, or translated into any other computer or foreign language either in whole or in part, or be reproduced in any way whatsoever.
4. The Software or any part of the Product shall not be lent out on loan, rent or lease or supplied to any third party in any manner.
5. The Licensee is responsible for and must use his best efforts to ensure that no other person breaks the terms and conditions of this agreement.
6. If the Licensee wishes to transfer the right to use the product to any other party he must transfer the whole Product and its original contents together to the same party and may not transfer part only. Upon transfer the Licensee shall destroy all backup copies made by him and shall delete from his computer all files installed from the Product. The Licensee shall ensure that the terms of this agreement are drawn to the attention of any transferee and any transferee shall, upon acceptance of the product, become subject to this Agreement.
7. Warranty - In the event that there exists any material defect in the Product as supplied to the Licensee, which affects the performance of the Product as set out in the manual, Bio-Rad Laboratories undertake to replace the Product supplied to the Licensee free of charge to the Licensee.
8. Disclaimer - The sole liability of Bio-Rad Laboratories in respect of the Product is as set out in paragraph 7 and all warranties and conditions (whether express or implied by statute or otherwise) relating to the merchantability or fitness for purpose of the Product are expressly excluded. Bio-Rad Laboratories shall not be liable (whether in contract or tort) in any manner whatsoever for any indirect or consequential loss or damage caused by or suffered by reason of any loss or corruption of data arising in the use of the Product.

PLEASE REMEMBER

Illicit copying of software increases the cost of products for bona fide software users and also prevents us from creating more and better products for you.