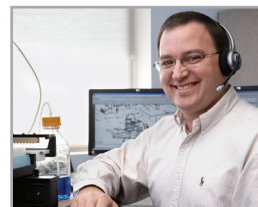
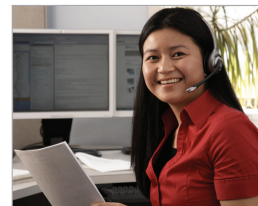


Our teams are committed to providing the service and support you need to sustain optimal performance of your BioTek products. Contact your local BioTek Service Center to discuss available warranty and contract coverage options for your new instrument.

In order to ensure trouble free operation of your instrument, BioTek recommends:

- ▶ Routine user maintenance per the operator's manual
- ▶ Regular Preventive Maintenance service
- ▶ Utilizing Absorbance / Fluorescence Test Plates to verify the continued accuracy of your microplate reader

Remember, Test Plates should be certified every year. The BioTek Global Technical Support Team is available in the United States to answer your instrument, software, operation and applications questions.



Guter Service und Support ist unser Selbstverständnis für die optimale Nutzung der BioTek Produkte. Regionale Service Center sind gerne für Sie da, wenn es um Dienstleistungsprogramme oder individuelle Lösungen geht.

Für eine lange Lebensdauer Ihres BioTek Gerätes empfehlen wir:

- ▶ Routinemäßige Durchführung der Wartungsintervalle (wie im Handbuch angegeben)
- ▶ Eine regelmäßige vorbeugende Wartung durch BioTek Service-Mitarbeiter
- ▶ Verwendung der entsprechenden Absorptions-oder Fluoreszenz-Testplatte für die kontinuierliche Überprüfung der Messgenauigkeit Ihres Mikroplatten-Readers

(Bitte berücksichtigen Sie, dass Testplatten einmal jährlich rezertifiziert werden müssen).

我们的团队承诺为您提供优异的售后服务和技术支持。以保证您所拥有的BioTek产品保持良好的工作状态。请联系当地的BioTek服务中心了解新产品的质保和相关服务条款。为了保证您的产品无障碍操作，BioTek推荐：

- ▶ 请按照操作手册要求来进行日常维护
- ▶ 定时进行预防性维护保养
- ▶ 使用吸收光/荧光测试板，确保您的微孔板检测仪保持检测的精准性

注意，测试板每年都需要进行校准。BioTek在美国的全球技术支持团队随时准备回答您有关产品、软件、操作和应用的相关问题。

Nuestros equipos están comprometidos en proveer el servicio y soporte que usted necesita para un desempeño óptimo de su producto BioTek. Contacte a su Centro Local de Servicio BioTek, para discutir las opciones de garantía y cobertura por contrato, disponibles para su nuevo instrumento.

Para asegurar una operación libre de problemas, BioTek recomienda para su equipo:

- ▶ Realizar los mantenimientos de rutina descritos en el manual del usuario
- ▶ Realizar servicios regulares de mantenimiento preventivo
- ▶ Utilizar Placas de Pruebas de Absorbancia/ Fluorescencia, para verificar se mantiene la exactitud de su lector de microplacas

Recuerde, las Placas de Prueba deben ser certificadas cada año. El equipo del Centro de Soporte Técnico Global de BioTek está disponible en los Estados Unidos, para responder a sus consultas sobre su equipo, software, operación y aplicaciones.

Available Service	Field Service	Service Center
Installation	•	
Installation Qualification	•	
Operational Qualification	•	
Training	•	•
Preventive Maintenance	•	•
Test Plate Certification		•
Dispense Cassette Calibration		•
Hardware Upgrades	•	•
Software Upgrades	•	•
Repair	•	•

BioTek Instrument Service and Repair is available worldwide at several of BioTek's International Service Centers and in the field at your location. Visit www.biotek.com/sales for a list of Service Centers by country.

BIOTEK DIRECT OFFICES

BioTek World Headquarters US Service Center

Service Shipping Address:
15 Tigan Street
Winooski, VT 05404, United States
Phone: (800) 242-4685
Email: TAC@biotek.com
Website: www.biotek.com

BioTek China - Beijing Service Center

Room 304, Tower D
Ocean International Center
62 Middle 4th East Ring Road
Chaoyang District
Beijing 100025, P.R. China
Phone: +86 (10) 85865569
Email: infochina@biotek.com
Website: www.biotekchina.com.cn

BioTek China - Shanghai Service Center

Room 407, Tower 2,
Kaixin International Plaza
1299 Zhangheng Road
Pudong New District
Shanghai 201203, P.R. China
Phone: +86 (10) 85865569
Email: infochina@biotek.com
Website: www.biotekchina.com.cn

BioTek Instruments SAS

Bureau de Liaison France
50 avenue d'Alsace
68025 Colmar Cedex, France
Phone: +33 (0) 389206329
Email: info@biotek.fr
Website: www.biotek.fr

BioTek Germany Service Center & European Coordination Center

Kocherwaldstr. 34
D-74177 Bad Friedrichshall, Germany
Phone: +49 (0) 71369680
Email: info@biotek.de
Website: www.biotek.de

BioTek India Service Center

A Wing, 601 Lotus Corporate Park
Graham Firth Compound,
Off W.E. Highway, Goregaon East
Mumbai 400063, India
Phone: +91 (77) 18810038
Email: biotek@biotek.in
Website: www.biotek.in

BioTek Japan Service Center

8F Andoh Building
2-7-1 Taito, Taito-ku
Tokyo 110-0016, Japan
Phone: +81 (0)3 5812 8109
Email: infojapan@biotek.com
Website: www.biotek.jp

BioTek South Korea Service Center

3F, Gyungnam building, 830-48
Yeoksam-dong, Gangnam-gu
Seoul, South Korea (135-936)
Phone: +82 (0) 2-562-4740
Email: korea@biotek.com
Website: www.biotekinstruments.co.kr

BioTek Switzerland Service Center

Zentrum Fanghöfli 8
6014 Luzern, Switzerland
Phone: +41 (41) 2504060
Email: info@biotek.ch
Website: www.biotek.ch

BioTek Taiwan Service Center

5F-4, No. 15, Lane 360, Sec. 1
Neihu Road, Neihu Dist.
Taipei, Taiwan R.O.C.
Phone: +886 (2) 26277725
Email: infotaiwan@biotek.com
Website: www.biotek.com

BioTek UK Service Center

Papermakers House
Rivenhall Road
Swindon SN5 7BD, UK
Phone: +44 (1793) 575020
Email: infouk@biotek.com
Website: www.biotek.co.uk