










Quick start guide | Kurzanleitung | Guide de démarrage rapide |
Guía de inicio rápido | Guida rapida

Biosealer® TC



SARTORIUS

English	page	4
Deutsch	page	6
Français	Seite	8
Español	página	10
Italiano	pagina	12

Symbol	Meaning
	Read "Operating Manual" before using the device!
	Hand injury warning Never reach into openings with your hands during operation. Noncompliance can lead to serious injury.
	Remove power cord before cleaning. Never open up the device!
	Warning of hot surface Inside the device, the surface can be hot and cause burns injury. These surfaces must not be touched. The device must first be cooled before maintenance cleaning.
	Ratings on marking plate must be followed!
	Only original equipment may be used!
	The device can be operated between +5°C and +40°C. The sealing parameters are qualified for room temperature operation only.
	The device was developed for laboratory use and must be protected from every type of fluid or vapour.
	The device is intended to use for technical qualified personnel only!

Operating Manual

Please download the latest operating manual and read it carefully!
www.sartorius.com/biosealer-tc

Installation

- Place the Biosealer® TC on a flat, stable surface
- Connect the power cord into the device socket from the power supply and to the according mains supply
- Connect the power supply with the Biosealer® TC
- Switch Biosealer® TC on

Initialization

- Remove everything from the sealing area
- Close cover to start the initialization process

Sealing

- Choose the right tubing type and diameter
- Insert the correct tubing
- Close cover
- Make sure that there is no pressure and tension on the tubing
- Press start button to start the run
- Follow carefully the instructions on the display to avoid any issues due to overpressure

After the sealing is complete:

- Remove tubing
- Check sealing visually according to operating manual before proceeding

Symbol	Bedeutung
	Vor Gebrauch des Gerätes "Bedienungsanleitung" lesen!
	Warnung vor Handverletzungen Greifen Sie während des Betriebs niemals mit den Händen in Öffnungen. Nichtbeachtung kann schwere Verletzungen verursachen.
	Netzkabel vor dem Reinigen entfernen. Niemals das Gerät öffnen!
	Die Oberflächen in diesem Bereich können sehr hohe Temperaturen aufweisen. Hautkontakt kann zu Verbrennungen Verletzungen führen. Vermeiden Sie direkten Hautkontakt in diesem Bereich und tragen Sie entsprechende Schutzkleidung -ausrüstung.
	Werte auf dem Typenschild müssen eingehalten werden!
	Es dürfen nur Originalteile verwendet werden!
	Stellen Sie sicher, dass die Umgebungstemperatur innerhalb +5 °C und +40 °C ist. Die Schweißparameter sind für Raumtemperatur qualifiziert.
	Das Gerät wurde für den Einsatz im Labor entwickelt und muss vor jeder Art von Flüssigkeiten oder Dämpfen geschützt werden.
	Das Gerät darf nur von qualifiziertem Fachpersonal benutzen werden!

Bedienungsanleitung

Bitte laden Sie die aktuelle Betriebsanleitung herunter und lesen Sie sie sorgfältig durch! www.sartorius.com/biosealer-tc

Installation

- Stellen Sie den Biosealer® TC auf eine flache, stabile Oberfläche
- Schliessen Sie das Netzkabel an der Gerätebuchse des Netzteils und der Steckdose an
- Verbinden sie das Netzteil mit dem Biosealer® TC
- Schalten Sie den Biosealer® TC ein

Initialisierung










- Entfernen Sie alles aus dem Versiegelungsbereich
- Schliessen Sie die Haube um die Initialisierung zu starten

Versiegelung

- Wählen Sie den richtigen Schlauchtyp und Durchmesser
- Legen Sie den richtigen Schlauch ein
- Schliessen Sie die Haube
- Achten Sie darauf, dass während des gesamten Versiegelungsvorgangs keine Zugkraft und kein Druck auf den Schlauch einwirkt
- Drücken Sie den Startknopf um die Schweissung zu starten
- Befolgen Sie die Anweisungen auf dem Display sorgfältig, um Probleme durch Überdruck zu vermeiden

Nach der Versiegelung:

- Entfernen Sie den Schlauch
- Überprüfen Sie die Versiegelung visuell gemäss Betriebsanleitung, bevor Sie fortfahren

Symbole	Explication
	Lisez "Manuel d'utilisation" avant d'utiliser la machine!
	Avertissement de blessure à la main Ne jamais introduire les mains dans les ouvertures pendant le fonctionnement. La non-conformité peut entraîner des blessures graves.
	Retirez le cordon d'alimentation avant de le nettoyer. N'ouvrez jamais l'appareil!
	La surface dans cette zone peut avoir des températures très élevées. Le contact avec la peau peut entraîner des brûlures blessures. Eviter le contact direct avec la peau dans cette zone et porter un vêtement équipement de protection correspondant. Utilisez toujours l'outil de décolleté (accessoire) pour enlever la lame.
	Les notes sur la plaque de marquage doivent être suivies !
	Seul l'équipement d'origine peut être utilisé !
	L'appareil peut fonctionner entre +5 °C et +40 °C. Les paramètres de soudage ne sont adaptés qu'à la température ambiante.
	L'appareil a été développé pour une utilisation en laboratoire et doit être protégé contre tout type de fluide ou de vapeur.
	L'appareil est destiné à être utilisé uniquement pour le personnel qualifié technique.

Manuel d'utilisation

Veuillez télécharger la dernière notice d'utilisation et la lire attentivement !
www.sartorius.com/biosealer-tc

Installation

- Placer le Biosealer® TC sur une surface plane et stable
- Brancher le cordon d'alimentation dans la prise de l'appareil à partir du bloc d'alimentation et sur le secteur correspondant.
- Raccorder l'alimentation électrique avec le Biosealer® TC
- Allumer le Biosealer® TC

Initialisation










- Enlever tout ce qui se trouve dans la zone de soudure
- Fermer le capot pour lancer le processus d'initialisation

Soudure

- Choisir le bon type et le bon diamètre de tube
- Insérez le bon tube
- Fermer le capot
- S'assurer qu'il n'y a pas de pression ou de tension sur le tube
- Appuyez sur le bouton de démarrage pour démarrer la course.
- Suivez attentivement les instructions sur l'écran pour éviter tout problème dû à une surpression

Une fois le soudage terminé:

- Retirer le tube
- Vérifier visuellement la soudure selon le manuel d'utilisation avant de procéder

Símbolo	Explicación
	Lea el “Manual de instrucciones” antes de utilizar la máquina!
	Advertencia de lesiones Nunca introduzca las manos en las aberturas durante el funcionamiento. El incumplimiento puede causar lesiones graves.
	Retire el cable de alimentación antes de limpiarlo. ¡Nunca abra el dispositivo!
	La superficie en esta área puede tener temperaturas muy altas. El contacto con la piel puede causar quemaduras lesiones. Evite el contacto directo con la piel en esta área y use la ropa equipo de protección correspondiente. Utilice siempre la herramienta removedora de cuchillas (accesorio) para quitar la cuchilla.
	¡Las calificaciones en la placa de marcado deben ser seguidas!
	¡Sólo se puede utilizar el equipo original!
	El dispositivo puede funcionar entre +5 °C y +40 °C. Los parámetros de soldadura están calificados para operación a temperatura ambiente solamente.
	El dispositivo fue desarrollado para uso en laboratorio y debe estar protegido de cualquier tipo de líquido o vapor.
	El dispositivo está diseñado para ser utilizado únicamente por personal técnico cualificado!

Manual de instrucciones

Por favor, descargue el manual de instrucciones más reciente y léalo detenidamente! www.sartorius.com/biosealer-tc

Instalación

- Coloque el Biosealer® TC sobre una superficie plana y estable.
- Conecte el cable de alimentación a la toma de corriente del aparato desde la fuente de alimentación y a la fuente de alimentación correspondiente.
- Conecte la fuente de alimentación con el Biosealer® TC
- Conectar Biosealer® TC

Inicialización










- Retire todo del área de sellado
- Cierre la cubierta para iniciar el proceso de inicialización.

Sellado

- Elija el tipo y diámetro de tubo adecuado
- Inserte el tubo correcto
- Cerrar la tapa
- Asegúrese de que no haya presión ni tensión en el tubo.
- Presione el botón de inicio para iniciar la ejecución
- Siga cuidadosamente las instrucciones de la pantalla para evitar cualquier problema debido a la sobrepresión.

Después de que se complete el sellado:

- Remueva la tubería
- Compruebe visualmente el sellado según el manual de instrucciones antes de proceder

Simbolo	Spiegazione
	Leggere "Manuale d'uso" prima di utilizzare la macchina!
	Avvertimento per lesioni alle mani Non toccare mai le aperture con le mani durante il funzionamento. La non conformità può portare a lesioni gravi.
	Rimuovere il cavo di alimentazione prima della pulizia. Non aprire mai il dispositivo!
	La superficie in questa zona può avere temperature elevate. Il contatto con la pelle può provocare ustioni lesioni. Evitare il contatto diretto con la pelle in questa zona e indossare indumenti protettivi attrezzature corrispondenti. Utilizzare sempre lo strumento di rimozione della lama (accessorio) per rimuovere la lama.
	Devono essere rispettati i punteggi sulla piastra di marcatura!
	Possono essere utilizzati solo apparecchi originali!
	Il dispositivo può essere azionato tra +5 °C e +40 °C. I parametri di saldatura sono qualificati solo per il funzionamento della temperatura ambiente.
	Il dispositivo è stato sviluppato per uso di laboratorio e deve essere protetto da ogni tipo di fluido o vapore.
	Il dispositivo è destinato esclusivamente a personale tecnico qualificato!

Manuale operativo

Si prega di scaricare le ultime istruzioni per l'uso e di leggerle attentamente!
www.sartorius.com/biosealer-tc

Installazione

- Posizionare il Biosealer® TC su una superficie piana e stabile.
- Collegare il cavo di alimentazione alla presa dell'apparecchio dall'alimentatore e alla relativa alimentazione di rete.
- Collegare l'alimentazione elettrica con il Biosealer® TC
- Accendere il Biosealer® TC

Inizializzazione

- Rimuovere tutto dall'area di sigillatura
- Chiudere la copertura per avviare il processo di inizializzazione

Guarnizione

- Scegliere il tipo e il diametro del tubo giusto
- Inserire il tubo corretto
- Chiudere il coperchio
- Assicurarsi che non ci sia pressione e tensione sul tubo.
- Premere il pulsante di avvio per avviare la corsa
- Seguire attentamente le istruzioni sul display per evitare problemi di sovrappressione.

Dopo che la sigillatura è stata completata:

- Rimuovere il tubo
- Controllare visivamente la sigillatura secondo le istruzioni per l'uso prima di procedere

Sartorius Stedim Biotech GmbH
August-Spindler-Strasse 11
37079 Goettingen, Germany

Phone: +49 551 308 0
www.sartorius.com

The information and figures contained in these instructions correspond to the version date specified below.

Sartorius reserves the right to make changes to the technology, features, specifications and design of the equipment without notice.

Masculine or feminine forms are used to facilitate legibility in these instructions and always simultaneously denote the other gender as well.

Copyright notice:

This instruction manual, including all of its components, is protected by copyright.

Any use beyond the limits of the copyright law is not permitted without our approval.

This applies in particular to reprinting, translation and editing irrespective of the type of media used.

© Sartorius Germany

Last updated:
10 | 2020