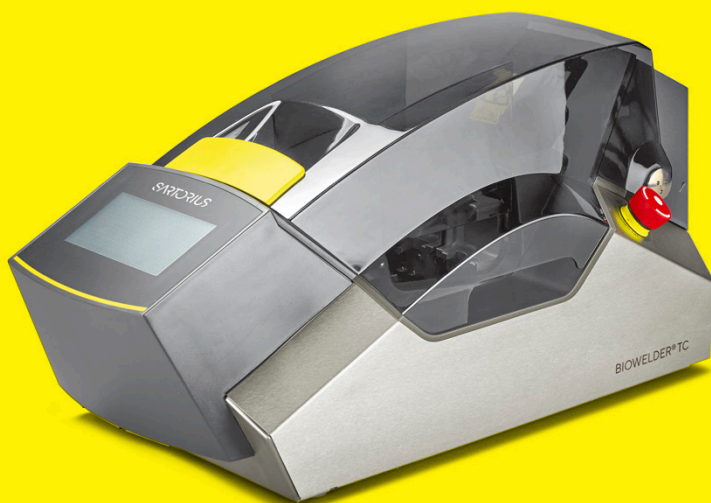





Quick start guide | Kurzanleitung | Guide de démarrage rapide |
Guía de inicio rápido | Guida rapida

Biowelder® 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 grab hold of anything inside openings during operation. Noncompliance can lead to serious injury
	In case of emergency, press emergency off button!
	Remove power cord before cleaning. Never open up the device!
	The surface in this area can have very high temperatures. Skin contact can result in burns injury. Avoid direct contact with skin in this area and wear corresponding protective clothing equipment. Always use the blade remover tool (accessory) to remove the blade.
	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 welding 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/biowelder-tc

Installation

- Place the Biowelder® TC on a flat, stable surface
- Connect the power cord into the device socket and to the according mains supply
- Release emergency off and switch Biowelder® TC on

Initialization

- Remove everything from the welding area
- Close cover to start the initialisation process

Weld

- Choose the right tube type (use minimum 45cm each tube)
- Insert tube holders
- Insert both tubes
- Insert a new blade
- Close cover
- Make sure that during the entire welding process there is no tension and no pressure on the tube, as well as free rotation on the left side
- Press start button to start the run

After the weld is complete:

- Remove the blade with the blade removing tool
- Remove both tubes
- Check welding visually before proceeding

Symbol	Bedeutung
	Vor Gebrauch des Gerätes "Bedienungsanleitung" lesen!
	Warnung vor Handverletzungen Während des Betriebes darf niemals mit den Händen in Öffnungen gegriffen werden. Nichtbeachtung kann schwere Verletzungen verursachen.
	Im Notfall die Not-Aus-Taste drücken!
	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/biowelder-tc

Installation

- Stellen Sie den Biowelder® TC auf eine flache, stabile Oberfläche
- Schliessen Sie das Netzkabel an der Gerätebuchse und der Steckdose an
- Entriegeln Sie den Not-Aus und schalten Sie den Biowelder® TC ein

Initialisierung

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

Schweissung

- Wählen Sie den richtigen Schlauchtyp (verwenden Sie mindestens 45 cm pro Schlauch)
- Setzen Sie die Schlauchhalter ein
- Legen Sie beide Schläuche ein
- Legen Sie eine neue Klinge ein
- Schliessen Sie die Haube
- Achten Sie darauf, dass während des gesamten Schweißvorganges keine Zugkraft und kein Druck auf den Schlauch einwirkt sowohl freie Rotation auf der linken Seite vorhanden ist
- Drücken Sie den Startknopf um die Schweissung zu starten

Nach der Schweissung:

- Entfernen Sie die Klinge mit dem Klingenentfernungstool
- Entfernen Sie beide Schläuche
- Überprüfen Sie die Schweissung visuell, bevor Sie fortfahren

Symbole	Explication
	Lisez "Manuel d'utilisation" avant d'utiliser la machine!
	Avertissement de blessure à la main Ne jamais saisir rien dans les ouvertures pendant l'opération. La non-conformité peut entraîner des blessures graves.
	En cas d'urgence, appuyez sur le bouton d'urgence!
	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. Éviter 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

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

Installation

- Placez le Biowelder® TC sur une surface plane et stable
- Connectez le cordon d'alimentation dans la prise du périphérique et à l'alimentation secteur
- Relâchez le bouton d'arrêt d'urgence et allumez le Biowelder® TC

Initialisation











- Enlever tout ce qui se trouve dans la zone de soudage
- Fermer la couverture pour commencer le processus d'initialisation

Souder

- Choisissez le type de tube correct (utiliser au minimum 45 cm chaque tube)
- Insérer des supports de tubes
- Insérer les deux tubes
- Insérer une nouvelle lame
- Fermer la couverture
- Veillez à ce que pendant tout le processus de soudage, il n'y ait pas de force de traction et de pression sur le flexible, ainsi qu'une rotation libre du côté gauche
- Appuyez sur le bouton de démarrage pour lancer la course

Une fois la soudure terminée:

- Retirer la lame avec l'outil d'enlèvement de lame
- Retirer les deux tubes
- Vérifier visuellement la soudure avant de procéder

Símbolo	Explicación
	Lea el “Manual de instrucciones” antes de utilizar la máquina!
	Advertencia de lesiones Nunca agarre nada dentro de las aberturas durante el funcionamiento. El incumplimiento puede causar lesiones graves.
	En caso de emergencia, presione el botón de apagado de emergencia!
	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/biowelder-tc

Instalación

- Coloque el Biowelder® TC sobre una superficie plana y estable
- Conecte el cable de alimentación en la toma del dispositivo y en la fuente de alimentación correspondiente
- Suelte el botón de parada de emergencia y encienda el Biowelder® TC

Inicialización

- Retirar todo del área de soldadura
- Cierre la cubierta para iniciar el proceso de inicialización

Soldar

- Elija el tipo de tubo correcto (usar un mínimo de 45cm cada tubo)
- Soportes de tubo de inserción
- Inserte ambos tubos
- Inserte una nueva pala
- Cierre la cubierta
- Asegúrese de que durante todo el proceso de soldadura no haya fuerza de tracción ni presión en la manguera, así como rotación libre en el lado izquierdo
- Pulse el botón de inicio para iniciar la ejecución

Una vez terminada la soldadura:

- Retire la cuchilla con la herramienta de extracción de la cuchilla
- Retire ambos tubos
- Compruebe visualmente la soldadura antes de proceder

Simbolo	Spiegazione
	Leggere "Manuale d'uso" prima di utilizzare la macchina!
	Avvertimento per lesioni alle mani Non tirate mai dentro le aperture durante il funzionamento. La non conformità può portare a lesioni gravi.
	In caso di emergenza, premere il pulsante di emergenza!
	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/biowelder-tc

Installazione

- Posizionare il Biowelder® TC su una superficie piana e stabile
- Collegare il cavo di alimentazione alla presa del dispositivo e alla rete di alimentazione corrispondente
- Rilasciare l'emergenza e accendere il Biowelder® TC

Inizializzazione

- Rimuovere tutto dall'area di saldatura
- Chiudere il coperchio per avviare il processo di inizializzazione

Saldare

- Scegliere il tipo di tubo corretto (utilizzare un minimo di 45 cm per ogni tubo)
- Inserire i porta tubi
- Inserire entrambi i tubi
- Inserire una nuova lama
- Chiudere il coperchio
- Accertarsi che durante l'intero processo di saldatura non vi siano forze di trazione e pressione sul tubo flessibile, nonché rotazione libera sul lato sinistro
- Premere il pulsante di avvio per avviare l'esecuzione

Dopo che la saldatura è completa:

- Rimuovere la lama con l'utensile di rimozione della lama
- Rimuovere entrambi i tubi
- Controllare visivamente la saldatura 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