

# Isotemp Advanced Hot Plates, Stirrers and Stirring Hotplates

OPERATING MANUAL  
AND PARTS LIST

CIC0000812 V11 11/29/17

---

**This manual covers the models shown below**

<b>NA Model</b>	<b>EU Model</b>	<b>Voltage</b>	<b>Description</b>
HP88857290	N/A	100-120V	7x7 Ceramic Hot Plate
HP88857295	15353528	220-240V-EU	7x7 Ceramic Hot Plate
HP88857296	15363528	220-240V-AUS/CHN/UK	7x7 Ceramic Hot Plate
S88857290	N/A	100-120V	7x7 Ceramic Stirrer
S88857295	15373528	220-240V-EU	7x7 Ceramic Stirrer
S88857296	15383528	220-240V-AUS/CHN/UK	7x7 Ceramic Stirrer
SP88857290	N/A	100-120V	7x7 Ceramic Stirring Hotplate
SP88857295	15393528	220-240V-EU	7x7 Ceramic Stirring Hotplate
SP88857296	15303538	220-240V-AUS/CHN/UK	7x7 Ceramic Stirring Hotplate
HP88857294	N/A	100-120V	7x7 Aluminum Hot Plate
SP88857294	N/A	100-120V	7x7 Aluminum Stirring Hotplate



Important Before using this product, read this entire operation manual carefully. Users should follow all of the operational guidelines contained in this manual and take all necessary safety precautions while using this product. Failure to follow these guidelines could result in potentially irreparable bodily harm and/or property damage.

Caution all internal adjustments and maintenance must be performed by qualified service personnel.

Material in this manual is for information purposes only. Fisher Scientific is committed to a continuing program of product development and improvement, and reserves the right to change information, such as specifications, appearance, and dimensions, described in this document without notice. Fisher Scientific makes no representations or warranties with respect to this manual. In no event shall Fisher Scientific be held liable for any damages, direct or incidental, arising out of or related to the use of this manual.

No part of this manual may be reproduced or transmitted in any form or by any means, including photocopying, recording, or using information storage and retrieval systems, for any purpose other than the purchaser's own use, without the express written permission of the manufacturer.

Any other product names and services identified in this manual are trademarks or registered trademarks of their respective owners. No such use, or the use of any trade name, is intended to convey endorsement or other affiliation with Fisher Scientific.

©2017 Fisher Scientific. All rights reserved.

---

# Table of Contents

Safety Information .....	6
Alert Signals .....	6
Warnings .....	6
Información de seguridad .....	8
Señales de alerta .....	8
Advertencias .....	9
Sicherheitshinweise .....	10
Gefahrensymbole .....	10
Warnungen .....	10
Informations de sécurité .....	12
Signaux d'alerte .....	12
Avertissements .....	13
Informação de Segurança .....	14
Sinais de Aviso .....	14
Avisos .....	14
Информация по безопасности .....	16
Сигналы тревоги .....	16
Предупреждения .....	17
General Specification .....	18
Ceramic Top Hot Plate Specifications .....	18
Ceramic Top Stirrer Specifications .....	18
Ceramic Top Stirring Hotplate Specifications .....	19
Aluminum Top Hot Plate Specifications .....	19
Heating Specifications* .....	20
Stirring Speed Specifications .....	20
Declaration of Conformity .....	21
Environmental Conditions .....	21
Introduction .....	22
General Cleaning Instructions .....	22
Unpacking and Installation .....	23
Unpacking .....	23
Installation .....	23
Isotemp Advanced Hot Plate Control Panel .....	24
Isotemp Advanced Stirrer Control Panel .....	24
Isotemp Advanced Stirring Hot Plate Control Panel .....	24
Control Panel .....	25
Principles of Operation .....	26

---

Operation .....	27
Power Switch .....	27
Setting the Stirring Speed .....	27
Setting the Temperature .....	28
Controlling Solution Temperature Using External Probe .....	29
Using Preset Programs .....	29
Heating Metal Vessels and Sand Baths .....	30
Setting the Time .....	30
Hardware Over-Temperature-Protection .....	30
System Configuration .....	31
Temp Unit .....	32
Over-temperature Set Point –OTP .....	32
Preset Continue .....	33
Temperature Probe Calibration .....	33
Troubleshooting .....	34
Error Codes .....	34
Accessories .....	35
Warranty .....	36
WEEE Compliance .....	37

---

# Safety Information

---

## Alert Signals



### Warning

Warnings alert you to a possibility of personal injury.



### Caution

Cautions alert you to a possibility of damage to the equipment.



### Note

Notes alert you to pertinent facts and conditions.



### Hot Surface

Hot surfaces alert you to a possibility of personal injury if you come in contact with a surface during use or for a period of time after use.



### Note

The Isotemp Advanced hot plates are not explosion proof. If explosion proof models are required, contact Customer Service at 1-800-553-0039 for more information.



### Warning

Refer servicing to qualified personnel.

Your Fisher Scientific Isotemp Advanced hot plate, stirrer or stirring hot plate has been designed with function, reliability, and safety in mind. It is your responsibility to install it in conformance with local electrical codes. For safe operation, please pay attention to the alert signals throughout the manual.

Warning: These products should be used only under the operating conditions specified in the Operating Manual.

- If the equipment is used in a manner not specified by the manufacturer, protection provided by the equipment may be impaired.
- Always use safe laboratory practices and do not leave the hotplate in operation while unattended as product functionality or laboratory practice failures could occur that might lead to uncontrolled or excessive heating of the top surface.
- Safety procedures (including, but not limited to, unplugging when not in use) and emergency response plans should be put in place to address the worst case possibility.
- If an over-temperature failure occurs, the top surface temperature could rise to the maximum temperature (300-450°C depending on your model's specification) and remain at that temperature indefinitely. Under these conditions, the material being heated on the surface of the hotplate could reach levels in excess of the maximum temperature.

This manual contains important operating and safety information. The user must carefully read and understand the contents of this manual prior to the use of this equipment.

---

## Warnings

### To avoid electrical shock, always:

1. Use a properly grounded electrical outlet of correct voltage and current handling capacity.

2. Disconnect from the power supply prior to maintenance and servicing.

### To avoid personal injury:

1. Do not use in the presence of flammable vapors outside of a fume hood or near combustible materials where heating could result in a fire or explosion. This device contains components which may ignite such materials. This device is not rated for use in designated hazardous locations where flammable vapors may reach the lower flammability limit.
2. Use caution when heating volatile materials; top surface and element can reach the "Flash Point Temperature" of many chemicals. These hot plates are not explosion proof. Fire or explosion may result. Unit contains components which may ignite such materials.
3. Keep top surface clean. Use a non-abrasive cleaner. Alkali spills, hydrofluoric acid spills or phosphoric acid spills may damage top and lead to thermal failure. Unplug unit and remove spills promptly. Do not immerse unit for cleaning.
4. Replace the top immediately if damaged by etching, scratching or chipping. A damaged top can break in use.
5. Do not use metal foil on hot plate which may block air flow. Overheating will result.
6. Check and tighten the removable cord periodically making sure it is secure. If loosened, the cord could become hot and/or spark and be a potential fire hazard. If cord appears damaged, replace immediately. If cord is repeatedly loosened it is recommended to purchase the power cord.
7. Do not remove or modify grounded power plug. Use only properly grounded outlets to avoid shock hazard.
8. Use appropriate hand and eye protection when handling hazardous chemicals.
9. Gross weight of items placed on top of hot plates should not exceed 35 lbs. (15.9 kg.) on the 10" x 10" models and 25 lbs. (11.3 kg) on the 7" x 7" models.
10. "Caution: Hot Top. Avoid Contact." The top plate of the unit can remain hot for some time after use. A "CAUTION-HOT TOP" light will remain on until top plate temperature cools to below 50°C.
11. Do not leave an active probe out of the fluid. This may cause uncontrolled heating of the fluid on the hot plate and unintentional boiling or an explosion could occur.
12. Localized heater element temperature can be significantly higher than the temperature indicated on the display. If flammable concentrations reach internal element, a fire could result.
13. Note that the exterior housing will be hot during and for a period of time after use.
14. Refer servicing to qualified personnel.

# Información de seguridad

## Señales de alerta



### Advertencia

Las advertencias le alertan de la posibilidad de lesiones personales.



### Precaución

Precauciones le alerta de posibilidades de daños en el equipo.



### Nota

Las notas le alertan sobre información y condiciones pertinentes.



### Superficie caliente

Las superficies calientes le alertan de la posibilidad de lesiones personales si tiene contacto con la superficie durante su uso o después de un determinado periodo de tiempo.

Su Plancha Caliente de Agitación Fisher Scientific Isotemp Advanced fue diseñada con la función, la fiabilidad y la seguridad en mente. Para su instalación es su responsabilidad hacerlo en conformidad con los códigos eléctricos locales. Para su seguro funcionamiento, por favor, preste atención a las señales de alerta especificados en el manual.

Advertencia: Estos productos deben ser usados siguiendo las condiciones de operación especificadas en el Manual de Operación.

- Si el equipo se utiliza en una manera distinta a la indicada por el fabricante, la protección del equipo puede verse afectada.
- Siempre use prácticas seguras de laboratorio y no deje la plancha de calentamiento en funcionamiento sin vigilancia ya que puede haber cualquier falla en la funcionalidad del producto o en la práctica de laboratorio que podría conducir a un calentamiento incontrolado o excesivo de la superficie superior.
- Se deben poner en marcha los procedimientos de seguridad (incluyendo, pero no limitado a, desconectar cuando no está en uso) y los planes de respuesta para hacer frente en el peor de los escenarios.
- Si se produce un fallo de exceso de temperatura, la temperatura de la superficie superior podría elevarse a una temperatura máxima de (300-540°C dependiendo de las especificaciones de su modelo) y permanecer a esa temperatura indefinidamente. En estas condiciones, el material se calienta en la superficie de la plancha calefactora que podría alcanzar niveles que superen la temperatura máxima permitida.

Este manual contiene información importante acerca del funcionamiento y seguridad. Antes de utilizar el equipo el usuario debe leer cuidadosamente y entender el contenido de este manual.



### Advertencias

#### Para evitar descarga eléctrica, siempre:

1. Utilice una toma eléctrica con polo a tierra con la correcta tensión y la capacidad de manejo de corriente.
2. Desconecte de la red eléctrica antes de hacer cualquier mantenimiento y/o reparación.

#### Para evitar lesiones personales:

1. No utilizar cerca de materiales inflamables o combustibles - podría conllevar a un incendio o una explosión. Este dispositivo contiene componentes que pueden encender estos materiales. No está clasificado para ser usado en entornos peligrosos.
2. Tenga cuidado al calentar materiales volátiles; la superficie superior y el elemento puede alcanzar a una "Temperatura Punto de inflamación" de muchos productos químicos. Estas planchas no son a prueba de explosiones. Podría producirse un incendio o una explosión. La unidad contiene componentes que pueden encender estos materiales.
3. Mantenga limpia la parte superior superficie. Use un limpiador no abrasivo. Los vertidos alcalinos, los derrames de ácido fluorhídrico o derrames de ácido fosfórico pueden dañar la parte superior y conllevar a una insuficiencia térmica. Desconecte la unidad y elimine los derrames de inmediato. No sumerja la unidad en líquidos para su limpieza.
4. Reemplace la tapa inmediatamente si está dañada por el efecto de los químicos, rayado o astillado. Una tapa dañada puede romperse durante su uso.
5. No utilice papel de aluminio en la plancha caliente este puede bloquear el flujo de aire dando lugar a un sobrecalentamiento.
6. Se debe revisar y apretar el cable desmontable periódicamente para asegurarse de que es seguro.

Si esta aflojado, el cable se podría calentar y/ o provocar chispas siendo un peligro potencial de incendio. Sustituya el cable inmediatamente si el cable está dañado. Si el cable se afloja continuamente se recomienda comprar el cable de energía eléctrica.

7. No retire o modifique el enchufe con polo a tierra. Utilice sólo enchufes con polo a tierra para evitar el riesgo de descarga eléctrica.
8. Utilice la protección adecuada para manos y ojos al manipular productos químicos peligrosos.
9. El peso bruto de los objetos que se pongan en la parte superior de las planchas calientes no debe exceder las 35 libras. (15,9 kg.) En los modelos 10"x 10", 25 libras. (11.3 kg) en los modelos 7"x 7" .
10. "Precaución: Hot Top. Evite el contacto. "La placa superior de la unidad puede permanecer caliente durante algún tiempo después de su uso. Una luz de "PRECAUCIÓN - HOT TOP" permanecerá encendida hasta que la temperatura plancha superior se enfríe por debajo de los 50°C.
11. No deje una sonda activa fuera del fluido. Esto puede causar un calentamiento incontrolado del fluido en la plancha caliente y podría producirse una ebullición o una explosión.
12. El elemento de temperatura del calentador localizado puede ser significativamente mayor que la temperatura indicada en la pantalla. Si las concentraciones inflamables alcanzan algún componente interno, podrían provocar un incendio.
13. Debe tener en cuenta que la carcasa exterior permanecerá caliente durante su uso y por un periodo de tiempo después de su uso.
14. Siempre solicite los servicios del personal cualificado.

# Sicherheitshinweise

## Gefahrensymbole



### Warnung

Eine Warnung weist Sie auf die Möglichkeit einer -Verletzung hin.



### Vorsicht

Das Symbol Vorsicht weist Sie auf die mögliche Beschädigung der Ausrüstung hin.



### Hinweis

Hinweise weisen Sie auf einschlägige Fakten und Bedingungen hin.



### Heiße Oberfläche

Das Symbol heiße Oberfläche weist Sie auf das Risiko von Verletzungen hin, sollten Sie während des Betriebs oder für einen bestimmten Zeitraum danach in Kontakt mit einer Oberfläche kommen.

Ihre Fisher Scientific Isotemp Advanced Rühr- Heizplatte wurde unter den Gesichtspunkten der Funktionalität, Zuverlässigkeit und Sicherheit entworfen. Es obliegt Ihrer Verantwortung diese in Übereinstimmung mit den örtlichen elektrischen Vorschriften zu installieren. Für einen sicheren Betrieb beachten Sie bitte die Gefahrensymbole in der gesamten Bedienungsanleitung.

Warnung: Diese Produkte sollten nur unter den in der Betriebsanleitung angegebenen Betriebsbestimmungen eingesetzt werden.

- Falls die Ausrüstung in einer vom Hersteller nicht angegebenen Weise benutzt wird, kann dadurch der Schutz, den das Gerät bietet, beeinträchtigt werden.
- Wenden Sie immer sichere Laborpraktiken an und lassen Sie die Heizplatte nicht eingeschaltet, während diese unbeaufsichtigt ist, da das Fehler der Produktfunktionalität oder der Laborpraktiken zu unkontrollierter oder übermäßiger Erwärmung der Oberfläche führen kann.
- Sicherheitsverfahren (einschliesslich, aber nicht beschränkt auf das Ziehen des Netzsteckers bei Nichtgebrauch) und Reaktionspläne sollten erstellt werden, um auf den Ernstfall vorbereitet zu sein.
- Wenn eine Störung zu einer Überhitzung führt, könnte die Oberflächentemperatur auf die Maximaltemperatur ansteigen (300-540°C, je nach Modellspezifikation) und für unbestimmte Zeit auf dieser Temperatur verbleiben. Unter diesen Bedingungen kann das zu erhitzende Material an der Oberfläche der Heizplatte ein höhere Temperatur als die Höchsttemperatur erreichen.

Dieses Handbuch enthält wichtige Bedienungs- und Sicherheitshinweise. Der Benutzer muss dieses Handbuch sorgfältig durchlesen und die Inhalte vor der Verwendung dieses Geräts verstehen.

## Warnungen

### Zur Vermeidung von Stromschlag, immer:

1. Eine ordnungsgemäß geerdete Steckdose verwenden, welche auf die korrekte Spannung und Strombelastbarkeit ausgelegt ist.

2. Das Gerät vor der Wartung und Reparatur von der Stromversorgung abtrennen.

## Um Verletzungen zu vermeiden:

1. Nicht in der Nähe von entzündlichen oder brennbaren Materialien verwenden - es besteht Brand- oder Explosionsgefahr. Dieses Gerät enthält Komponenten, die solche Materialien entzünden können. Nicht für den Einsatz in Bereichen freigegeben, in denen Explosionsgefahr besteht.
2. Vorsicht beim Erhitzen flüchtiger Materialien. Die Oberfläche und das Element können die Temperatur des "Entflammungspunkts" vieler Chemikalien erreichen.  
Die Heizplatten sind nicht explosionsgeschützt. Dies kann zu einem Brand oder einer Explosion führen. Dieses Gerät enthält Komponenten, die solche Materialien entzünden können.
3. Halten Sie die Oberfläche sauber. Verwenden Sie ein nicht scheuerndes Reinigungsmittel. Leckagen von Alkali, Flußsäure oder Phosphorsäure können die Oberfläche beschädigen und zum Ausfall der Heizkomponenten führen. Ziehen Sie den Netzstecker der Einheit und entfernen Sie Leckagen unverzüglich. Nicht zur Reinigung in Flüssigkeiten eintauchen.
4. Ersetzen Sie die Oberfläche sofort, falls es durch Verätzungen, Kratzer oder Absplitterungen beschädigt ist. Eine beschädigte Oberfläche kann während dem Gebrauch brechen.
5. Verwenden Sie keine Metallfolie auf der Heizplatte, welches den Luftstrom blockieren kann. Das kann Überhitzung zur Folge haben.
6. Überprüfen und ziehen Sie das abnehmbare Kabel regelmäßig an und gewährleisten, dass es sicher ist. Wenn gelockert, kann das Kabel heiß werden und/oder Funken schlagen und eine potentielle Brandgefahr darstellen. Wenn das Kabel beschädigt erscheint, sofort ersetzen. Wenn das Kabel sich wiederholt lockert, wird es empfohlen, das Netzkabel zu kaufen.
7. Den geerdeten Netzstecker nicht entfernen oder modifizieren. Verwenden Sie nur geerdete Steckdosen, um einen Stromschlag zu vermeiden.
8. Benutzen Sie geeigneten Hand und Augenschutz beim Umgang mit gefährlichen Chemikalien.
9. Das Bruttogewicht der Gegenstände, die auf die Heizplatten gegeben werden sollte 15,9 kg nicht überschreiten (35 lbs). Auf den 10 Zoll x 10 Zoll Modellen, 11,3 kg (25 lbs) auf den 7 Zoll x 7 Zoll Modellen.
10. "Vorsicht: Heiße Oberfläche. Kontakt vermeiden."  
Die obere Heizplatte des Geräts kann für einige Zeit nach der Benutzung heiß bleiben. Eine Warnleuchte "ACHTUNG - HEISSE OBERFLÄCHE" wird aufleuchten bis die Temperatur der oberen Heizplatte auf unter 50°C abkühlt.
11. Lassen Sie keine Flüssigkeit aus der aktiven Probe entweichen. Dies kann eine unkontrollierte Erwärmung der Flüssigkeit auf der Heizplatte zur Folge haben und ein unbeabsichtigtes Sieden oder eine Explosion verursachen.
12. Die lokalisierte Heizelementtemperatur kann wesentlich höher sein als die Temperatur die auf dem Display angegeben ist. Wenn brennbare Mischungen innenliegende Elemente erreichen, könnte dies zu einem Brand führen.
13. Beachten Sie, dass das Außengehäuse während dem Gebrauch und eine gewisse Zeit danach heiß ist.
14. Kontaktieren Sie für alle Wartungsarbeiten qualifiziertes Servicepersonal.

# Informations de sécurité

## Signaux d'alerte



### Avertissement

Les avertissements vous alertent de la possibilité de blessures corporelles.



### Mise en garde

Les mises en garde vous avertissent de la possibilité d'endommagement de l'équipement.



### Remarque

Les remarques vous alertent de conditions et de faits pertinents.



### Surface chaude

Les surfaces chaudes vous avertissent de la possibilité de blessures corporelles, si vous entrez en contact avec une surface en cours d'utilisation ou pour une période de temps après l'utilisation.

Votre Plaque Chauffante d'Agitatrice Isotemp de Fisher Scientific a été conçue, tout en ayant à l'esprit: la fonction, la fiabilité et la sécurité. Il est de votre responsabilité de l'installer en conformité aux codes électriques locaux. Pour un fonctionnement en toute sécurité, veuillez prêter attention aux signaux d'alerte mentionnés à travers le manuel.

Avertissement: Ces produits ne devraient être utilisés que dans des conditions d'utilisation spécifiées dans le manuel d'utilisation.

- Si l'équipement est utilisé d'une manière non-spécifiée par le fabricant, la protection fournie par l'équipement peut en être réduite.
- Employer toujours des pratiques de laboratoire sûres et ne laissez pas la plaque chauffante fonctionner sans surveillance puisque des défaillances de fonctionnement du produit ou de pratique de laboratoire pourraient survenir et conduire à un chauffage incontrôlé ou excessif de la surface supérieure.
- Des procédures de sécurité (y compris, mais non limité au débranchement lorsqu'il n'est pas en cours d'utilisation) et des plans d'intervention devraient être mis en place pour résoudre le pire des scénarios.
- Si une panne de surchauffe se produit, la température de la surface supérieure pourrait augmenter jusqu'à la température maximale (300-540°C, en fonction des spécifications de votre modèle) et demeurer indéfiniment à cette température. Dans ces conditions, le matériel chauffé sur la surface de la plaque chauffante pourrait atteindre des niveaux au-delà de la température maximale.

Ce manuel contient d'importantes informations relatives au fonctionnement et à la sécurité. L'utilisateur doit attentivement lire et comprendre le contenu de ce manuel avant d'utiliser cet équipement.

### **Avertissements**

#### **Pour éviter tout risque de choc électrique, toujours:**

1. Utiliser une prise électrique correctement mise à la terre avec une tension et capacité de gestion du courant correctes.
2. Débrancher de la source d'alimentation, avant de procéder à l'entretien et la maintenance.

#### **Pour éviter les blessures:**

1. Ne pas utiliser en présence de matériaux inflammables ou combustibles - cela peut entraîner un incendie ou une explosion. Cet appareil contient des composants qui peuvent enflammer ces matériaux. Ne pas utiliser dans des lieux dangereux.
2. Faire preuve de prudence lors du chauffage de matériaux volatils; la surface supérieure et l'élément.
3. Garder propre la surface supérieure. Utiliser un produit de nettoyage non abrasif. Les déversements alcalis, les déversements d'acide fluorhydrique ou les déversements d'acide phosphorique peuvent endommager la partie supérieure et conduire à une défaillance thermique. Débrancher l'appareil et enlever rapidement les déversements. Ne pas plonger l'unité dans de l'eau pour le nettoyer.
4. Remplacer immédiatement la partie supérieure, si elle est endommagée par une attaque chimique, une rayure ou un écaillage. Une partie supérieure endommagée peut tomber en panne en cours d'utilisation.
5. Ne pas utiliser la feuille de métal sur la plaque chaude, cela peut bloquer la circulation d'air. Une surchauffe pourra se produire.
6. Vérifiez et serrez périodiquement la corde amovible afin d'assurer sa sécurisée. Si la corde se desserre, elle va devenir chaude et/ou produire de l'étincelle et aussi devenir un risque d'incendie potentiel. S'il y a des endommagements, on doit remplacer la corde immédiatement. Si la corde se desserre plusieurs fois, on vous recommande d'acheter la corde d'alimentation.
7. Ne pas retirer ou modifier la fiche d'alimentation mise à la terre. Utiliser uniquement des prises mises à la terre correctement, afin d'éviter tout risque de choc électrique.
8. Utiliser des lunettes et des gants de protection, lors de la manipulation de produits chimiques dangereux.
9. Le poids brut des éléments placés sur la partie supérieure des plaques chauffantes ne doit pas dépasser 35 lbs. (15,9 Kg.) sur les modèles 10" x 10", 25 lbs. (11,3 Kg) sur les modèles 7" x 7" .
10. "Mise en garde: Partie supérieure Chaude. Éviter tout contact." La partie supérieure de la plaque de l'unité peut rester chaude pendant un certain temps après utilisation. Un témoin "ATTENTION - PARTIE SUPÉRIEURE CHAUDE" restera allumé jusqu'à ce que la température de la partie supérieure de la plaque descende en-dessous de 50°C.
11. Ne pas laisser de sonde active hors du liquide. Cela peut provoquer un réchauffement incontrôlé du liquide sur la plaque chaude et une ébullition involontaire ou une explosion pourraient se produire.
12. La température de l'élément chauffant localisé peut être sensiblement plus élevée que la température indiquée sur l'affichage. Si des concentrations inflammables atteignent l'élément interne, un incendie pourrait se produire.
13. Notez que le boîtier extérieur sera chaud pendant et pour une période de temps après utilisation.
14. S'adresser à un personnel qualifié pour la maintenance.

# Informação de Segurança

## Sinais de Aviso



### Aviso

Os avisos alertam para a possibilidade de ferimentos pessoais.



### Cuidado

Os sinais de cuidados alertam para a possibilidade de danos no equipamento.



### Nota

As notas alertam para factos e condições pertinentes.



### Superfície Quente

Os sinais de superfície quente alertam para a possibilidade de ferimentos pessoais se entrar em contacto com uma superfície durante a utilização ou por um período de tempo após a utilização.

A Chapa Eléctrica Isotemp Térmica de Fisher Scientific foi desenhada com funcionalidade, confiança e segurança em mente. É sua responsabilidade instalar de acordo com os regulamentos eléctricos locais. Para operação segura, preste atenção aos sinais de alerta presentes no manual.

Aviso: Estes produtos devem ser unicamente utilizados sob as condições de operação especificadas no Manual de Utilização.

- Se o equipamento for utilizado de forma não especificada pelo fabricante, a protecção fornecida pelo equipamento poderá ser anulada.
- Siga sempre todas as práticas de laboratório seguras e não deixe a chapa eléctrica em funcionamento sem supervisão, uma vez que a funcionalidade do produto ou falhas de práticas de laboratório podem ocorrer e conduzir ao aquecimento não controlado ou excessivo na superfície superior.
- Os procedimentos de segurança (incluindo, mas não limitado a, desligar da ficha quando não estiver em utilização) e planos de resposta devem ser postos em prática no pior caso possível.
- Em caso de falha de sobreaquecimento, a temperatura da superfície superior pode aumentar para a temperatura máxima (300 - 540°C dependendo das especificações do seu modelo) e manter-se na mesma durante um tempo indefinido. Sob estas condições, o material a ser aquecido na superfície da chapa pode alcançar níveis excessivos de temperatura.

Este manual contém informação de operação e segurança importante. O utilizador deve ler atentamente e compreender os conteúdos deste manual antes de utilizar o equipamento.

## Avisos

### Para evitar choque eléctrico:

1. Utilize sempre uma tomada eléctrica aterrada com a voltagem e capacidade de manuseamento de corrente correctas.

2. Desligue sempre a alimentação antes da manutenção e assistência.

### **Para evitar ferimentos pessoais:**

1. Não utilize na presença de materiais combustíveis ou inflamáveis - poderá resultar em incêndio e explosão. Este dispositivo contém componentes que podem incendiar esses materiais. Não aprovado para utilização em atmosferas perigosas.
2. Tenha cuidado quando aquecer materiais voláteis; a superfície e elemento superiores podem alcançar a "Temperatura de Ponto de Incêndio" de muitos químicos. Estas chapas eléctricas não são à prova de explosão. Poderá resultar em incêndio ou explosão. A unidade contém componentes que.
3. Mantenha a superfície superior limpa. Utilize um detergente não abrasivo. Derramamentos de alcalinos, ácidos hidrofluóricos ou ácidos fósforos podem danificar a superfície e conduzir a falha termal. Desligue a unidade da tomada e limpe os derramamentos de imediato. Não mergulhe a unidade para limpeza.
4. Substitua a superfície de imediato se danificada por cauterização, arranhões ou lascas. A superfície pode partir durante a utilização.
5. Não utilize folha de metal na chapa, pois poderá bloquear o fluxo de ar. Poderá resultar em sobreaquecimento.
6. Verifique e aperte o cabo removível periodicamente para ter certeza que está seguro. Caso esteja frouxo, o cabo poderá se aquecer e/ou soltar faíscas, gerando perigo de incêndio. Se o cabo parecer danificado, substitua-o imediatamente. Se o cabo se afrouxar com frequência, é recomendada a aquisição do cabo de alimentação.
7. Não retire nem modifique a ficha de alimentação aterrada. Utilize unicamente tomadas correctamente aterradas para evitar perigo de choque eléctrico.
8. Utilize protecção de mãos e olhos quando manusear químicos perigosos.
9. O peso bruto dos produtos colocados em cima da chapa não deve exceder os 35 lbs. (15.9 kg.) em modelos 10" x 10", 25 lbs. (11.3 kg.) em modelos 7" x 7".
10. "Cuidado: Topo Quente. Evite Contacto." A chapa superior da unidade pode permanecer quente durante algum tempo após a utilização. A luz "CUIDADO - TOPO QUENTE" irá permanecer ligada até a temperatura descer abaixo dos 50°C.
11. Não deixe um estilete activo fora do fluído. Isto poderá causar aquecimento descontrolado do fluído na chapa e poderá resultar em fervura ou explosão não intencional.
12. A temperatura de elemento de aquecimento localizada pode ser significativamente superior à temperatura indicada no ecrã. Se as concentrações inflamáveis alcançarem o elemento interno, poderá resultar em incêndio.
13. Note que o revestimento exterior estará quente durante a utilização e após a mesma.
14. Refira a assistência ao pessoal qualificado.

# Информация по безопасности

## Сигналы тревоги



### Предупреждение

Символы предупреждения указывают на возможность получения травмы.



### Осторожно

Символы осторожно указывают на возможность повреждения оборудования.



### Примечание

Примечания указывают на соответствующие факты и условия.



### Горячая поверхность

Символ горячей поверхности предупреждает о возможности получения травмы, если вы вступите в контакт с поверхностью во время использования или в течение периода времени после использования.

Термо нагревательная плита Термо Фишер Сайентифик была разработана с функцией надежностью и безопасности. Вам необходимо установить ее в соответствии с местными нормативами. Для безопасной эксплуатации, пожалуйста, обратите внимание на сигналы тревоги по всему руководству.

Предупреждение: Данные продукты следует использовать только в рабочих условиях, указанных в руководстве по эксплуатации.

- Если устройство используется отличным от указанного производителем образом, предусмотренная защита оборудования может быть повреждена.
- Всегда следуйте технике безопасности и не оставляйте работающую плиту без присмотра, так как функциональность продукта или несоблюдение техники безопасности может привести к неконтролируемому или чрезмерному нагреву верхней поверхности.
- Техника безопасности (включая, но не ограничиваясь, отсоединение от розетки, во время, когда плита не используется) и планы реагирования должны быть приняты для решения крайней необходимости.
- Если происходит сбой из-за перегрева, температура верхней поверхности может подняться до максимальной температуры (300-540°C в зависимости от спецификации вашей модели) и оставаться при этой температуре в течение длительного времени. В этих условиях, продукты, нагреваемые на поверхности плиты, могут достичь уровня выше максимальной температуры.

Это руководство содержит важную операционную информацию и информацию по безопасности. Пользователь должен внимательно прочитать и понять содержание данного руководства перед началом использования данного оборудования.



### Предупреждения

#### Во избежание поражения электрическим шоком, всегда:

1. Используйте правильно заземленную розетку соответствующего напряжения и пропускной способностью тока.
2. Отсоединяйте от источника питания перед обслуживанием и сервисом.

#### Во избежание получения травмы:

1. Не используйте возле легковоспламеняющихся или горючих материалов – это может привести к пожару или взрыву. Это устройство содержит компоненты, которые могут воспламенить эти материалы. Не предназначено для использования в опасной атмосфере.
2. Будьте внимательны при нагревании летучих веществ; верхняя поверхность и элемент могут достигнуть "температуры вспышки" многих химических веществ. Данные нагревательные плиты не взрывозащищенные. Это может привести к пожару или взрыву. Блок содержит компоненты, которые могут воспламенить эти материалы.
3. Держите верхнюю поверхность в чистоте. Используйте неабразивный очиститель. Щелочные растворители, растворители с содержанием фтористоводородной кислоты или фосфорной кислоты могут повредить верхний слой и привести к термическим повреждениям. Отключайте устройство и тщательно удаляйте растворители. Не опускайте прибор в воду для чистки.
4. Немедленно замените верхнюю поверхность, если она повреждена в результате травления, царапин или обкалывания. Поврежденная верхняя поверхность может сломаться во время использования.
5. Не используйте металлическую фольгу на нагревающей плите, это может блокировать поток воздуха и привести к перегреву.

6. Проверить и заменить съемный кабель, в процессе проверяя безопасность. Если кабель слабо затянут, то он может нагреваться или искрить, что может быть причиной возгорания. Если кабель поврежден, его необходимо немедленно заменить. Если кабель постоянно слабозатянут, то рекомендуется приобрести кабель питания.
7. Не удаляйте или не изменяйте заземленную вилку. Используйте только правильно заземленные розетки, чтобы избежать поражения электрическим током.
8. Используйте соответствующие защитные средства для рук и глаз при работе с опасными химическими веществами.
9. Общий вес предметов, помещаемых на плиту, не должен превышать 35 фунтов (15,9 кг) на 10" x 10" модели, 25 фунтов (11,3 кг) на 7" x 7" модели.
10. "Осторожно: Горячая поверхность. Избегайте контакта." Верхняя плита блока может оставаться горячей некоторое время после использования. Надпись "ОСТОРОЖНО – ГОРЯЧАЯ ПОВЕРХНОСТЬ" будет гореть до тех пор, пока температура верхней плиты не остынет до температуры ниже 50°C.
11. Не оставляйте активную емкость с чрезмерным количеством жидкости. Это может привести к неконтролируемому нагреву жидкости на горячей плите и непреднамеренному кипению или взрыву.
12. Локализованная температура нагревательного элемента может быть значительно выше, чем температура, указанная на дисплее. Если горючие концентрации достигают внутреннего элемента, это может привести к пожару.
13. Необходимо принять во внимание, что внешний корпус будет горячим во время и в течение периода времени после использования.
14. За сервисным обслуживанием обращайтесь к квалифицированному персоналу.

# General Specification

## Ceramic Top Hot Plate Specifications

Overall Dimensions	HP88857290 (NA)	HP88857295 (NA)	HP88857296 (NA)
		15353528 (EU)	15363528 (EU)
Width - in.(cm)	8.2 (20.8)	8.2 (20.8)	8.2 (20.8)
Height - in.(cm)	4.1 (10.5)	4.1 (10.5)	4.1 (10.5)
Depth - in.(cm)	14.2 (36)	14.2 (36)	14.2 (36)
Weight - lbs.(kg)	7.5 (3.4)	7.5 (3.4)	7.5 (3.4)
<b>Top Plate Dimensions</b>			
Width - in.(cm)	7.25 (18.4)	7.25 (18.4)	7.25 (18.4)
Height - in.(cm)	1.0 (2.5)	1.0 (2.5)	1.0 (2.5)
Depth - in.(cm)	7.25 (18.4)	7.25 (18.4)	7.25 (18.4)
Volts	100-120	220-240	220-240
Amps	12.3	4.4	4.4
Watts	1480	1056	1056
Freq.	50/60	50/60	50/60
Phase	1	1	1
Max. Temp °C (°F)	450°C(842°F)	450°C(842°F)	450°C(842°F)

## Ceramic Top Stirrer Specifications

Overall Dimensions	S88857290 (NA)	S88857295 (NA)	S88857296 (NA)
		15373528 (EU)	15383528 (EU)
Width - in. (cm)	8.2 (20.8)	8.2 (20.8)	8.2 (20.8)
Height - in. (cm)	4.1 (10.5)	4.1 (10.5)	4.1 (10.5)
Depth - in. (cm)	14.2 (36)	14.2 (36)	14.2 (36)
Weight - lbs. (kg)	9.0 (4.1)	9.0 (4.1)	9.0 (4.1)
<b>Top Plate Dimensions</b>			
Width - in. (cm)	7.25 (18.4)	7.25 (18.4)	7.25 (18.4)
Height - in. (cm)	1.0 (2.5)	1.0 (2.5)	1.0 (2.5)
Depth - in. (cm)	7.25 (18.4)	7.25 (18.4)	7.25 (18.4)
Volts	100-120	220-240	220-240
Amps	0.2	0.1	0.1
Watts	24	24	24
Freq.	50/60	50/60	50/60
Phase	1	1	1

## Ceramic Top Stirring Hotplate Specifications

Overall Dimensions	SP88857290 (NA)	SP88857295 (NA)	SP88857296 (NA)
		15393528 (EU)	15303538 (EU)
Width - in. (cm)	8.2 (20.8)	8.2 (20.8)	8.2 (20.8)
Height - in. (cm)	4.1 (10.5)	4.1 (10.5)	4.1 (10.5)
Depth - in. (cm)	14.2 (36)	14.2 (36)	14.2 (36)
Weight - lbs. (kg)	9.0 (4.1)	9.0 (4.1)	9.0 (4.1)
<b>Top Plate Dimensions</b>			
Width - in. (cm)	7.25 (18.4)	7.25 (18.4)	7.25 (18.4)
Height - in. (cm)	1.0 (2.5)	1.0 (2.5)	1.0 (2.5)
Depth - in. (cm)	7.25 (18.4)	7.25 (18.4)	7.25 (18.4)
Volts	100-120	220-240	220-240
Amps	12.3	4.4	4.4
Watts	1480	1056	1056
Freq.	50/60	50/60	50/60
Phase	1	1	1
Max. Temp °C (°F)	450°C(842°F)	450°C(842°F)	450°C(842°F)

## Aluminum Top Hot Plate Specifications

Overall Dimensions	HP88857294 (NA)	SP88857294 (NA)
Width - in. (cm)	8.2 (20.8)	8.2 (20.8)
Height - in. (cm)	4.1 (10.5)	4.1 (10.5)
Depth - in. (cm)	14.2 (36)	14.2 (36)
Weight - lbs. (kg)	7.5 (3.4)	9.0(4.1)
<b>Top Plate Dimensions</b>		
Width - in. (cm)	7.25 (18.4)	7.25 (18.4)
Height - in. (cm)	1.0 (2.5)	1.0 (2.5)
Depth - in. (cm)	7.25 (18.4)	7.25 (18.4)
Volts	100-120	100-120
Amps	6.0	6.0
Watts	720	720
Freq.	50/60	50/60
Phase	1	1
Max. Temp °C (°F)	300°C(572°F)	300°C(572°F)

## GENERAL SPECIFICATIONS

### Heating Specifications\*

#### Top Plate Surface - Solid Ceramic

Models	7" x 7"	10" x 10"
Temperature range	30°C - 450°C (86°F - 842°F)**	30°C - 400°C (86°F - 752°F)**
Heat-up time to maximum temperature (unloaded top plate)	5 minutes	7 minutes
Accuracy of the temperature display vs. the actual average temperature of a 2" diameter of setting area at the center of the top plate (setpoint 100°C unloaded)	± 5.0°C	
Temperature stability at the center of the top plate surface (@ 100°C unloaded)	± 1.0°C	
Typical accuracy of remote probe at user selected calibration temperature after calibration procedure	± 0.5°C	
Temperature stability using remote probe (500 ml of water in a 1000 ml flask at 70°C)	± 0.5°C	

#### Top Plate Surface - Solid Aluminum

Models	7" x 7"
Temperature range	30°C - 300°C (86°F - 572°F)**
Heat-up time to maximum temperature (unloaded top plate)	5 minutes
Accuracy of the temperature display vs. the actual average temperature of a 2" diameter of setting area at the center of the top plate (setpoint 100°C unloaded)	± 5.0°C
Temperature stability at the center of the top plate surface (@ 100°C unloaded)	± 1.0°C
Typical accuracy of remote probe at user selected calibration temperature after calibration procedure	± 0.5°C
Temperature stability using remote probe (500 ml of water in a 1000 ml flask at 70°C)	± 0.5°C

### Stirring Speed Specifications

Speed Range 50 to 1500 RPM (Maximum speed is dependent on the viscosity of the solution).

Stability of the stirring speed setpoint (600 ml of water in a 1000 ml glass flask) ± 2.0% at 1000 RPM

Top Plate Size	Max Recommended Flask Size	Max Weight on Top Plate
7" x 7"	4 liters	25 lbs
10" x 10"	6 liters	35 lbs

\* The specification defined in this section is derived from testing data statistics of Thermo Fisher Scientific lab. We reserve the right to make changes in design and specification without prior notice.

\*\* As the top plate becomes dirty, the maximum temperature will decrease. To return the unit to its maximum temperature performance, use a mild abrasive to remove stained areas.

---

## Environmental Conditions

Operating: 5°C to 35°C; 20% to 80% relative humidity, non-condensing. Installation category II (overvoltage) in accordance with IEC 664. Pollution degree 2 in accordance with IEC 664. Altitude Limit: 2,000 meters.

Storage: -25°C to 65°C  
10% to 85% relative humidity

---

## Declaration of Conformity

We hereby declare under our sole responsibility that this product conforms with the technical requirements of the following standards:

EMC:	EN 61000-3-2	Limits for harmonic current emissions
	EN 61000-3-3	Limits for voltage fluctuations and flicker
	EN 61326-1	Electrical equipment for measurement, control, and laboratory use; Part I: General Requirements
Safety:	EN 61010-1	Safety requirements for electrical equipment for measurement, control, and laboratory use; Part I: General Requirements
	EN 61010-2-010	Part II: Particular requirements for laboratory equipment for the heating of materials
	EN 61010-2-051	Part II: Particular requirements for laboratory equipment for mixing and stirring

*per the provisions of the Electromagnetic Compatibility Directive 2014/30/EU, and per the provisions of the Low Voltage Directive 2014/35/EU.*

Copies of the Declaration of Conformity are available upon request.

---

# Introduction



Isotemp Advanced Stirring Hotplate

Please read all the information in this manual before operating the unit.

Your Isotemp Advanced hot plate, stirrer or stirring hot plate is a heating and/or stirring plate designed for laboratory procedures requiring precise control of temperature and/or stirring speed. Each Isotemp Advanced model includes a liquid crystal display for monitoring actual and setting temperature and/or stirring speed. The hot plate is capable of producing accurately controlled top plate temperatures from 30°C through 300°C, 400°C or 450°C depending on models. The temperature is controlled at the plate surface by an internal Type K thermocouple sensor, or the solution temperature may be controlled by utilizing the included PT100 probe. A PT1000 probe may be ordered separately. The 7 x 7 or 10 x 10 top plate on the Isotemp Advanced units are solid ceramic or solid aluminum, and are suitable for use with glass or metal vessels and sand baths.

Other features of the Isotemp Advanced include a time feature that can be used to shut off heating. Specific temperature setpoints may be set using the PROGRAM buttons on the Isotemp Advanced unit. An over temperature protection (OTP) can be displayed in the system configuration menus. See details in “System Configuration”.

Your Isotemp Advanced model may be used for general purpose heating applications including sample preparation, heating reagents, melting paraffin, warming resinous chemicals, content analysis, solvent evaporations, digestions, media preparation and sterilization, titrations, sand baths, and micro-scale chemistry applications.

---

## General Cleaning Instructions

Keep top surface clean. Use a non-abrasive cleaner. Alkali spills, hydrofluoric acid spills or phosphoric acid spills may damage top and lead to thermal failure. Unplug unit and remove spills promptly. Do not immerse unit for cleaning. Wipe exterior housing with lightly dampened cloth containing mild soap solution.

---

# Unpacking and Installation



## Warning

Use a properly grounded electrical outlet of correct voltage and current handling capacity.

Do not remove or modify grounded power plug. Use only properly grounded outlets to avoid shock hazard. Not rated for use in hazardous atmospheres.

Do not use in the presence of flammable or combustible materials; fire or explosion may result. This device contains components which may ignite such materials. If explosion proof models are required, contact Customer Service at 1-800-553-0039 for more information.

Do not use in highly corrosive atmospheres; corrosive fumes and spills may damage top and internal components, creating shock hazard.



## Note

Some misalignment of the motor bearings in this product may have occurred during shipping. Prior to using this product, run the stirrer at maximum speed for ten minutes to realign the bearings.

---

## Unpacking

Remove your Isotemp Advanced unit from the carton. Inspect to ensure that the unit has not been damaged during shipment. If the unit appears to have sustained shipping damage contact the distributor from whom you purchased this product or Customer Service at 800-553-0039. Check for remote probe and thumbscrew prior to discarding packaging.

### ***The following items are included in the shipment:***

Isotemp Advanced unit

Cord

Remote Probe – PT100

Stir Bar (Stirring models only)

Knob

Operating Manual

If any of these items are missing from the carton, contact Customer Service.

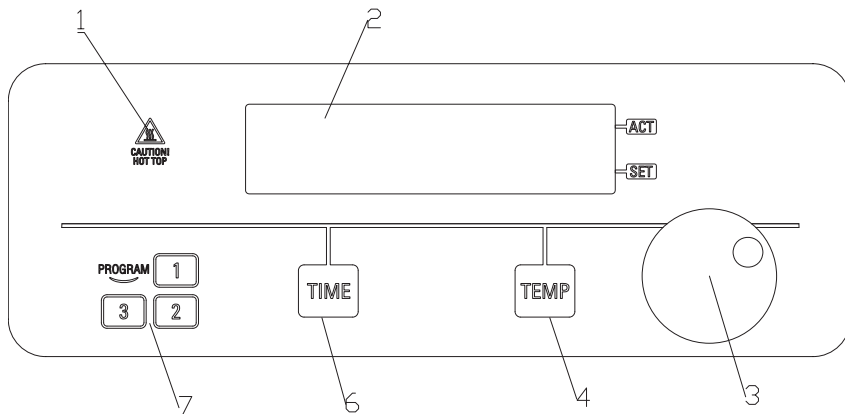
---

## Installation

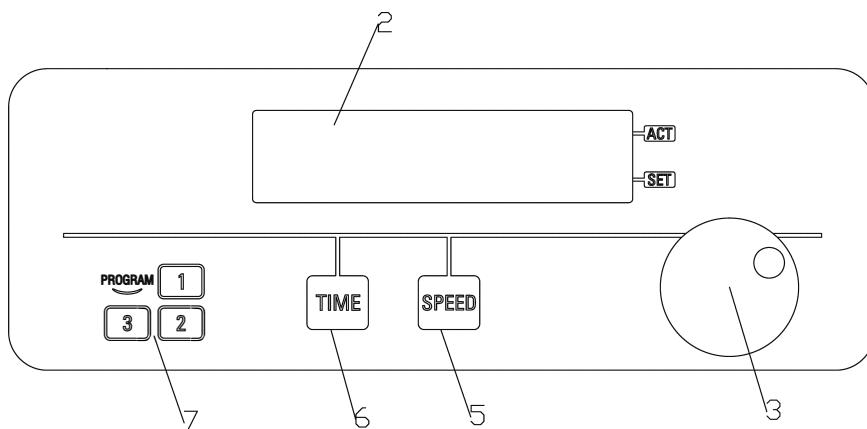
Set the unit on a flat stable surface at least 12" away from combustible materials, and plug the cord set into a properly grounded electrical outlet of correct voltage and current handling capacity.

# UNPACKING AND INSTALLATION

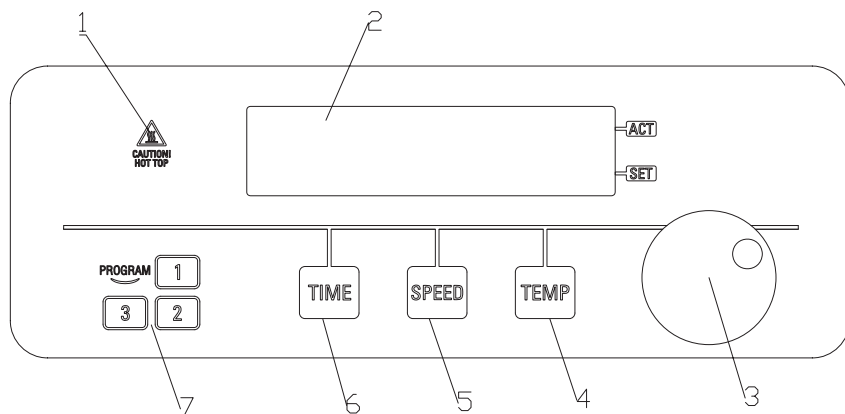
## Isotemp Advanced Hot Plate Control Panel



## Isotemp Advanced Stirrer Control Panel



## Isotemp Advanced Stirring Hot Plate Control Panel





---

## Control Panel

<b>Key</b>	<b>Description</b>
1	CAUTION! HOT TOP Indicator
2	Display
3	Knob
4	Temperature Set Button
5	Speed Set Button
6	Time Set Button
7	Program Buttons

---

# Principles of Operation

Each Isotemp Advanced unit utilizes the latest microprocessor technology to deliver the most reliable, accurately controlled, ceramic top or aluminum top hot plate on the market.

Your Isotemp Advanced unit with heating functionality has an electronic, closed-loop feedback control which will accurately maintain temperature setpoints from 30°C through 300°C, 400°C or 450°C depending on models.

The Isotemp Advanced units use a Type K thermocouple for measuring the top plate temperature, OTP and the external probe. Both top plate surface and OTP temperature are measured with independent analog circuits. This provides a significant increase in safety due to the redundancy of the circuits. It is possible to heat certain types of metal vessels and sand baths on the Isotemp Advanced without damaging the ceramic top or aluminum top.

The electronic stirring speed control will maintain the speed setpoint when the unit is loaded. The motor is combined with a powerful magnet to provide exceptional magnetic coupling with a stir bar.

The time feature on the Isotemp Advanced can be programmed to turn off the stirring, heating or both. See “Setting the Time” section of this manual.

# Operation



## Warning

Use caution when heating volatile materials; top surface and element can reach the "Flash Point Temperature" of many chemicals. These hot plates are not explosion proof. Fire or explosion may result. Unit contains components which may ignite such materials.

Use appropriate hand and eye protection when handling hazardous chemicals.

"Caution: Hot Top. Avoid Contact." The top plate of the unit can remain hot for some time after use. A 'CAUTION - HOT TOP' light will remain on until top plate temperature cools to below 50°C.



## Note

Boiling times are dependent on solution volume and the surface area of the flask that is exposed to the hot plate. For example, when heating the same amount of solution in a 2L flask vs. a 1L flask, the solution will heat about 20% faster.



## Note

The top plate temperature will be higher than the solution temperature.

## Power Switch

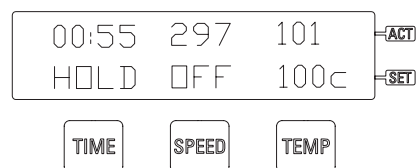
The Isotemp Advanced unit has Power Switch on the left side of the unit. The unit is powered on by pressing the Power Switch. When the unit is turned on, the unit will initialize and then model name will be displayed on the first row of display screen while software version in second row. To turn off power to the unit, press the Power Switch off. It is recommended to unplug the unit when not in use for 24 hours or longer.



## Setting the Stirring Speed

The Isotemp Advanced stirring unit has an electronic feedback speed control which will maintain a precise speed setpoint from 50 rpm through 1500 rpm. (Maximum speed is dependent on viscosity of the solution.) The Isotemp Advanced stirring hot plates are equipped with a strong magnet and high torque motor.

To set the stirring speed, press "SPEED" key located under display. Then the speed setpoint will flash. Use the knob on the front panel to set a speed. Turn the knob clockwise to increase the speed or counterclockwise to decrease the speed. When desired setpoint has been reached, press the "SPEED" key or wait 10 seconds, the speed setpoint will no longer flash indicating that it has been set.



Push the knob to turn on stirring, once the unit begins stirring, the display will indicate both actual speed and setpoint. To stop the stirring action without changing your setpoint press the knob once. To resume stirring at the set speed, press the knob again.

To turn stirring off, press "SPEED" key under the display and turn the knob counterclockwise until the display reads "OFF".



## Note

Boiling times are dependent on solution volume and the surface area of the flask that is exposed to the hot plate. For example, when heating the same amount of solution in a 2L flask vs. a 1L flask, the solution will heat about 20% faster.

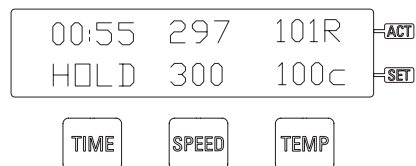


## Note

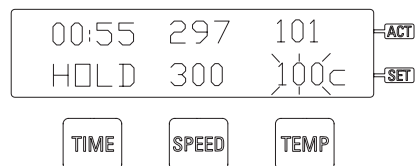
The top plate temperature will be higher than your solution temperature.

## Setting the Temperature

The Isotemp Advanced heating unit has an electronic closed-loop feedback control which will accurately maintain temperature setpoints from 30°C through 300°C, 400°C or 450°C (depending on the model). An unloaded hot plate will heat to maximum temperature in just 5 minutes. A "CAUTION HOT TOP" light on the front panel will illuminate whenever the top surface temperature exceeds 50°C. The temperature may be controlled either at the top plate by the internal Type K thermocouple sensor, or in the solution using an accessory ungrounded PT100 probe. If a probe is used, the display will indicate "R" after actual temperature. The Isotemp Advanced heating units will accept any ungrounded PT100 (PT1000) probe, however the accuracy of probes other than those tested and specified may not reach the probe accuracy rating designated.



To set the temperature, press "TEMP" key located under display. Then the temp setpoint will flash. Use the knob on the front panel to select a temp. Turn the knob clockwise to increase the temp or counterclockwise to decrease the temp. When desired setpoint has been reached, press the "TEMP" key or wait 10 seconds, the temp setpoint will no longer flash indicating that it has been set, then push the knob to start heating.



## Controlling Solution Temperature Using External Probe

To control the solution temperature plug a PT100 probe into the probe receptacle located on the back side of the unit. When the probe is connected, the adjustable temperature range is 30-250°C. Place the probe into the solution. The display will indicate the actual temperature of the solution as measured by the probe.

The external probe offers more exact temperature control than regulating the top plate by the internal sensor. If you need to maintain a precise setpoint it is recommended to use a probe to control the solution temperature instead of controlling by the top plate temperature.

When using a probe with the Isotemp Advanced unit it is recommended that a clamp on a support rod be used to hold the probe in the solution.

To ensure accurate probe readings, the terminal of probe should not touch the container side or bottom, and as much of the probe sheath as possible should be immersed in the solution. Make sure the probe is immersed in the liquid and is not located in air or outside of the solution. If the probe is plugged into the hot plate, but is not in solution while the heat control is operating, the temperature display will continue to indicate an ambient temperature, and a probe out of solution error will occur. Because the set point cannot be reached the heating element will continue to supply heat to the top plate, and the maximum top plate temperature(300°C, 400°C or 450°C) may be reached. If the remote probe does not sense a temperature change in 15 minutes the unit will display ExProbe Err and shut down.

## Using Preset Programs

The Isotemp Advanced unit features the option of presetting 3 setpoints into its memory. To set a preset, press any "preset" button, and the indicator LED on the number button will be illuminated. Then adjust the time, speed and heat features to your desired setpoints. Press the number button again to exit the preset program. The setpoints are stored in memory automatically. The presets will be held in memory, even if the unit is unplugged, until you choose to override it with another value.

To activate one of the presets, press the corresponding number then press the knob, the time, speed and heat features will automatically adjust itself.

The Isotemp Advanced unit has a preset continuous mode. In continuous mode, the preset setting value can run continuously.

By pressing 1 and then pressing the knob the unit will run by continuous presets by order 1, 2, 3. By selecting 2 and then pressing the knob the unit will run by continuous presets by order 2, 3.

The unit will stop running when the next preset value of temperature and speed are both OFF, or preset 3 is completed.

The default setting is OFF. See "System Configuration" section to active or disable the preset continuous mode.

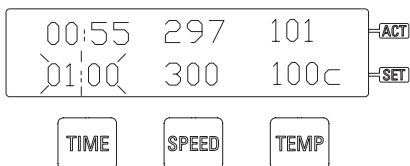
# OPERATION

## Heating Metal Vessels and Sand Baths

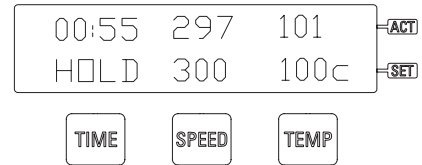
Metal vessels and sand baths cannot be heated on most solid ceramic tops because of the tendency for the metal and the sand to reflect heat back into the top, eventually exceeding the maximum temperature rating for a ceramic top, which causes it to break. Because of its advanced electronic control, the Isotemp Advanced is capable of precisely regulating the top plate temperature. Metal vessels and sand baths may be heated safely without the danger of the ceramic top breaking.

## Setting the Time

Isotemp advanced provides two timer modes. One is countdown mode. the timer function on the Isotemp advanced can shut off heating when time is reached. To set the time, press "TIME" key located under display. Hold will begin to flash. Turn the knob until time shows then press "TIME" key again. Use the knob on the front panel to select a time. Turn the knob clockwise to increase the time or counterclockwise to decrease the time. When desired setpoint has been reached, press the "TIME" key or wait 10 seconds, the time setpoint will no longer flash indicating that it has been set. In the temperature rising mode, when the temperature is more than 2 degree below the temperature setpoint, the timer will start to count. In the temperature declining mode, when the temperature is less than 2 degree above the temperature setpoint, the timer will start to count. When the timer counts down to 0, the unit will turn heating off.



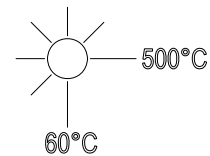
Another mode is timing mode. when the display shows "HOLD", the timer records the total running duration.



## Hardware Over-Temperature-Protection

A hardware Over-Temperature-Protection(OTP) analog circuit is designed separately to protect unexpected heating.

A small straight blade screwdriver can be used to adjust the scale at right side of the unit. The protection value can be set from 60°C to 500°C(450°C depending on models). When the actual temperature is higher than the Hardware OTP setting, an error message occurs and the heating is cut off, unplug the power cord can clear the error message.



Please notice that this function is only purposed to avoid abnormal heating if unit is out of control. Users should not use this function to limit the temperature in normal applications. Refer to the section "Over-temperature Set Point – OTP" for accurate temperature protection.

Default set is at maximum scale.

# System Configuration

The System Configuration has four features that will allow a user to customize their unit.

A table of the features available in the System Configuration is given below. The features are given in order of their appearance in the System Configuration when rotating the knob in a clockwise rotation. The

Feature column is the name of the feature, the Display column is the designation of that feature as shown on the display of the unit, the Availability column gives which type of unit the feature is applicable (HP = hot plate SP = stirring hot plate S=stir plate), and the last column gives a brief description of the purpose of the feature.

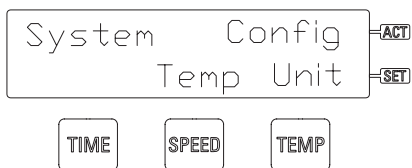
Feature	Display	Availability	Brief Description
Temperature Unit Conversion Over-temperature Set Point	Temp Unit	HP , SP	Allows the user to converse temperature unit between °C and °F.
Over-temperature Set Point	User OTP	HP , SP	Allows the user to see the setting of the Over Temperature Protection control.
Preset Continue Setting	Preset Continue	HP , S, SP	Allows user to activate or enable the preset continuous running
Temperature Probe Calibration	Calibrate Temp	HP , SP	Allows calibration of the temperature probe and associated circuitry only when the probe is connected.

# System Configuration

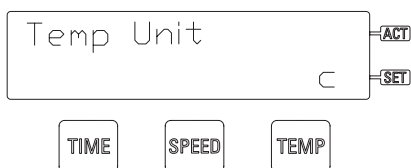
## Temp Unit

This option is only available for units that have heating functionality.

Plug the unit into the appropriate power, enter the System Configuration by pressing and holding the knob. After about 5 seconds you may remove your finger from the knob. The display indicate "System Config" on the first row, the "Temp Unit" will show on the second row of the display.



Press the knob to enter the Temp Unit setting interface. (Long press will return to operation menu.) The "Temp Unit" will display on the first row, the Actual temperature and unit "c" or "F" flash on the second row. Turn the knob clockwise or counterclockwise to change the temperature unit. When the desired temperature unit has reached, wait 10 seconds, the display will return to operation menu indicating that it has been set.



## Over-temperature Set Point – OTP

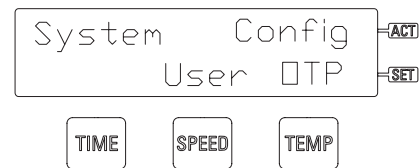
This option is only available for units that have heating functionality. The Over-temperature Set Point is the set point of the independent temperature control. The purpose of the control is to limit the top temperature of the element independently of the standard heating control.

This is helpful if there is a failure of the main control that would cause the top to heat excessively. Typically the set point will be adjustable from 60 to 350°C, 450°C or 500°C(depending on the model).

The Over-temperature can be set or viewed with this menu option. To do so, follow the instructions given below:

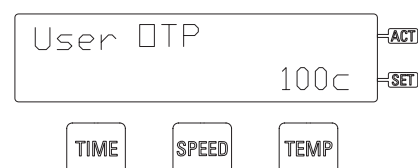
Enter the System Configuration by pressing and holding the knob. After about 5 seconds a single you may remove your finger from the knob. The display indicate "system config" on the first row, the "Temp Unit" will show on the second row of the display.

Turn the knob clockwise until the "User OTP" show on the second row of the display. If you do no action, the display will return to the operation status.



Press the knob to enter the Over-temperature setting interface.(Long press will return to operation menu.) The "User OTP" will display on the first row, the Over-temperature setpoint flash on the second row. Turn the knob clockwise to increase the temp or counterclockwise to decrease the temp. When the desired Over-temperature setpoint has been reached, wait 10 seconds, the display will return to operation menu indicating that it has been set.

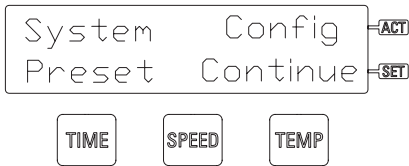
When the actual top plate temperature is higher than the OTP setting, the heating will be shut OFF, and error message "OVER TEMP" occurs, unplug the power cord can clear the error message.





## Preset Continue

Enter the System Configuration by pressing and holding the knob. After about 5 seconds a single you may remove your finger from the knob. The display indicate "System Config" on the first row, the "Temp Unit" will show on the second row of the display. Turn the knob clockwise until the "Preset Continue" show on the second row of the display. If you do no action, the display will return to the operation status.



Press the knob to enter the Preset Continue setting interface.(Long press will return to operation menu.)

The "User OTP" show on the first row, "OFF" flash on the second row. Turn the knob clockwise to change the status. When "On" has reached, wait 10 seconds, the display will return to operation menu indicating that it has been set.



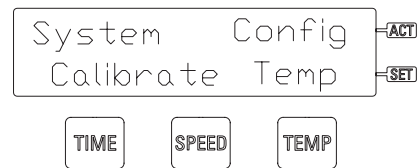
## Temperature Probe Calibration

This section applies only to the calibration of the temperature probe system. The hot plate surface and the motor speed cannot be calibrated. This method is only applicable to units that have heating functionality and with

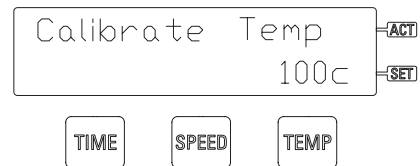
the probe connected.

Before calibrate the probe, connect the probe on the unit and put the probe sensor into the temperature reference.

Enter the System Configuration by pressing and holding the knob. After about 5 seconds a single you may remove your finger from the knob. The display indicate "System Config" on the first row, the "Temp Unit" will show on the second row of the display. Turn the knob clockwise until the "Calibrate Temp" show on the second row of the display. If you do no action, the display will return to the operation status. The option "Calibrate Temp" will only show when there is a remote probe PT100 or PT1000 adopted.



Press the knob to enter the Calibrate Temp setting interface. The "Calibrate Temp" show on the first row, the measured temperature flash on the second row. If the data is not matched the temperature you measure by external thermo detector, you can turn the knob to calibrate. When it's matched, press the knob to complete the calibration.



When setting and/or viewing is complete, the knob can be pressed to return to the previous operation menu.

# Troubleshooting

## Error Codes

Error Handler will lock out heating functions if heating error is detected. Stirring functionality is unaffected. Error

Handler will lock out stirring functions if stirring error is detected. Heating functionality is unaffected. If the condition that caused the error is no longer present, the unit will clear Errors automatically.

Displayed Message	Intended to Detect	Cause	Solution
InProbe Err	Internal thermocouple out of range.	Internal thermocouple not connected. Thermocouple open. Thermocouple connected backwards (reversed polarity).	Call technical service
Check Heater	Excessive top heat-up time	Internal thermocouple short circuit. Failure in Internal thermocouple. Failure in Element. Failure in optocoupler/triac circuit.	Call technical service
ExProbe Err	External Probe left out of solution.	External Probe left out of solution. External probe connected backwards.	Place external probe into solution.
Over Temp	OTP thermocouple out of range.	OTP Circuit failure. OTP potentiometer set too low.	Increase OTP setting.
Check Motor	Motor system failure.	Locked rotor condition. Failure of motor. Failure of motor circuit. Failure of motor sensor.	Free locked rotor. Call technical service

# Accessories

NA Order No.	EU Order No.	Description
CIC0001462	15373538	Splash guard protection shield 4", for 4 x 4 Stirrers and Hotplate Stirrers
CIC0001463	15383538	Splash guard protection shield 7", for 7 x 7 Stirrers and Hotplate Stirrers
CIC0001464	15393538	Splash guard protection shield 10", for 10 x 10 Stirrers and Hotplate Stirrers
CIC0001465	15303548	PT100 External probe, Stainless steel, for all sizes of Hotplates and Hotplate Stirrers
CIC0001466	15313548	PT1000 External probe, Stainless steel, for all sizes of Hotplates and Hotplate Stirrers
CIC0001467	15323548	Hastelloy PT1000 External probe, Hastelloy, for all sizes of Hotplates and Hotplate Stirrers
CIC0001472	N/A	Power cord assembly China plug 10A 250V
CIC0001474	N/A	Power cord assembly British BS 1363 plug 10A 250V
CIC0001473	N/A	Power cord assembly EURO plug 10A 250V
CIC0001475	N/A	Power cord assembly US&JP plug 16A 125V
CIC0001832	N/A	Fisher Scientific supporting Rod to be used together with clamps and clamps holder to hold temperature probe and flask, suitable for all sizes of hotplates, stirrers, stirring hotplates
CIC0001833	N/A	Fisher Scientific clamps and clamp holder to attach on supporting rod, suitable for all sizes of hotplates, stirrers, stirring hotplates

---

# Warranty

When used in laboratory conditions and according to these operation instructions and maintenance, this product is warranted for 36 months against defective materials or workmanship. The 36 month warranty period begins from the delivery date of this product.

For product quality or performance issues, contact Fisher Scientific Customer Service.

# WEEE Compliance

WEEE Compliance. This product is required to comply with the European Union's Waste Electrical & Electronic Equipment (WEEE) Directive 2002/96EC. It is marked with the following symbol. Fisher Scientific has contracted with one or more recycling/disposal companies in each EU Member State, and this product should be disposed of or recycled through them. Further information on our compliance with these Directives, the recyclers in your country, and information on Fisher Scientific products which may assist the detection of substances subject to the RoHS Directive are available at [www.fishersci.com/](http://www.fishersci.com/)

## Great Britain



WEEE Konformität. Dieses Produkt muss die EU Waste Electrical & Electronic Equipment (WEEE) Richtlinie 2002/96EC erfüllen. Das Produkt ist durch folgendes Symbol gekennzeichnet. Fisher Scientific hat Vereinbarungen getroffen mit Verwertungs-/Entsorgungsanlagen in allen EU-Mitgliedstaaten und dieses Produkt muss durch diese Firmen wiederverwertet oder entsorgt werden. Mehr Informationen über die Einhaltung dieser Anweisungen durch Fisher Scientific, die Verwerter und Hinweise die Ihnen nützlich sein können, die Fisher Scientific Produkte zu identifizieren, die unter diese RoHS-Anweisung fallen, finden Sie unter [www.fishersci.com/](http://www.fishersci.com/)

## Deutschland



Conformità WEEE. Questo prodotto deve rispondere alla direttiva dell'Unione Europea 2002/96EC in merito ai Rifiuti degli Apparecchi Elettrici ed Elettronici (WEEE). È marcato col seguente simbolo. Fisher Scientific ha stipulato contratti con una o diverse società di riciclaggio/smaltimento in ognuno degli Stati Membri Europei. Questo prodotto verrà smaltito o riciclato tramite queste medesime. Ulteriori informazioni sulla conformità di Fisher Scientific con queste Direttive, l'elenco delle ditte di riciclaggio nel Vostro paese e informazioni sui prodotti Fisher Scientific che possono essere utili alla rilevazione di sostanze soggette alla Direttiva RoHS sono disponibili sul sito [www.fishersci.com/](http://www.fishersci.com/)

## Italia



Conformité WEEE. Ce produit doit être conforme à la directive euro-péenne (2002/96EC) des Déchets d'Equipements Electriques et Electroniques (DEEE). Il est marqué par le symbole suivant. Fisher Scientific s'est associé avec une ou plusieurs compagnies de recyclage dans chaque état membre de l'union européenne et ce produit devrait être collecté ou recyclé par celles-ci. Davantage d'informations sur la conformité de Fisher Scientific à ces directives, les recycleurs dans votre pays et les informations sur les produits Fisher Scientific qui peuvent aider la détection des substances sujettes à la directive RoHS sont disponibles sur [www.fishersci.com/](http://www.fishersci.com/)

## France



## North America

United States  
1-800-766-7000  
fishersci.com

Canada  
1-800-234-7437  
fishersci.ca

## Europe

Austria:  
+43(0)800-20 88 40  
at.fishersci.com

Netherlands:  
+31(0)20 4887 70 00  
nl.fishersci.com

Belgium:  
+32(0)56 260 260  
be.fishersci.com

Norway:  
+47 22 95 59 59  
fishersci.no

Denmark:  
+45 70 27 99 20  
fishersci.dk

Portugal:  
+351 21 425 33 50  
pt.fishersci.com

Germany:  
+49(0)180 5258221  
de.fishersci.com

Spain:  
+34 002 239 303  
es.fishersci.com

Ireland:  
+959(0)1 885 5854  
ie.fishersci.com

Sweden:  
+46 31- 68 94 30  
fishersci.se

Italy:  
+39 02 950 59 478  
it.fishersci.com

Switzerland:  
+41(0)56 618 41 11  
ch.fishersci.com

Finland:  
+358(0)9 8027 6280  
fishersci.fi

UK:  
+44(1)1509 555 500  
fisher.co.uk

France:  
+33(0)388 67 14 14  
fishersci.fr

Find out more at **fishersci.com**

© 2017 Thermo Fisher Scientific Inc. All rights reserved.  
Trademarks used are owned as indicated at fishersci.com/trademarks.