**Thermo Scientific** 

#### Heratherm Hornos Advanced Protocol y Advanced Protocol Security OGH 60/100/180

OGH 60/100/180 OGH 60-S/100-S/180-S OMH 60/100/180/400/750 OMH 60-S/100-S/180-S

50129640 A

Manual de instrucciones

14.06.12





© 2012 Thermo Fisher Scientific Inc. Todos los derechos reservados.

Estas instrucciones de operación están protegidas por derechos de autor. Los derechos resultantes del mismo, en particular, de reimpresión, de posprocesamiento o reproducción por métodos fotomecánicos o digitales, incluso de manera parcial, están permitidos sólo con el consentimiento escrito de Thermo Fisher Scientific. Esta disposición no rige para las reproducciones destinadas al uso en la planta.

El contenido de este manual de operación puede ser modificado en cualquier momento y sin aviso previo. En el caso de versiones confusas en lenguas extranjeras, la versión en idioma alemán de estas instrucciones de operación será vinculante.

#### Marcas registradas

Heratherm<sup>®</sup> es una marca registrada de Thermo Scientific.

Thermo Scientific es una marca propiedad de Thermo Fisher Scientific, Inc.

Todas las demás marcas registradas que se mencionan en las instrucciones de operación son de propiedad exclusiva de los fabricantes respectivos.

Thermo Electron LED GmbH Robert-Bosch-Straße 1 D - 63505 Langenselbold Alemania

Thermo Electron LED GmbH es filial de: Thermo Fisher Scientific Inc. 81 Wyman Street Waltham, MA 02454 EEUU

Thermo Fisher Scientific Inc. suministra este documento a sus clientes con la compra de un producto, para ser usado en la operación de este último. Este documento está protegido por derechos de copia y cualquier reproducción total o parcial del mismo queda estrictamente prohibida, salvo autorización escrita de Thermo Fisher Scientific Inc. El contenido de este documento es pasible de modificaciones sin previo aviso.

Toda la información técnica en este documento sirve solamente para propósitos de referencia. Las configuraciones y especificaciones de sistema en el presente documento reemplazan toda información previa recibida por el comprador. Thermo Fisher Scientific Inc. no declara que este documento sea completo, preciso o sin errores no asumiendo responsabilidad, ni responderá por ningún error, omisión, daño o pérdida que podría resultar del uso de este documento, aún en el caso de que se siga, de manera adecuada, la información incluida en el mismo. Este documento no constituye ningún contrato de ventas entre Thermo Fisher Scientific Inc. y un comprador. Este documento no condicionará ni modificará Términos y Condiciones de Venta, dichos Términos y Condiciones de Venta prevalecerán sobre toda información conflictiva entre ambos documentos.

## Tabla de contenido

Capítulo 1	Notas de seguridad	1-1
-	Precauciones básicas de operación	1-1
	Reglas de seguridad operativa	1-2
	Garantía	1-2
	Aclaración de la información y símbolos de seguridad	1-3
	Notas y símbolos de seguridad utilizados a lo largo de estas instrucciones de operación	1-3
	Símbolos adicionales para la información de seguridad	1-4
	Símbolos en el horno	1-5
	Utilización prevista para el horno	1-5
	Uso correcto	1-5
	Uso incorrecto	1-6
	Normas y directivas	1-6
Capítulo 2	Envío del horno	2-i
	Embalaje	2-i
	Inspección de aceptación	2-i
	Alcance del suministro	2-ii
Capítulo 3	Instalación	3-1
	Condiciones ambientales	
	Requisitos de ubicación	
	Uso con sistemas de escape de aire	
	Almacenamiento intermedio	
	Ventilación del recinto	
	Uso con sistemas de escape de aire	
	Requisitos de espacio	
	Instalación, unidades empotradas	
	Unidades de mesa	
	Unidades para piso	
	Transporte	
	Unidades de mesa	
	Juego de apilado	
	Instalación de los pies de apilado	
	Instalación del adaptador de apilado	
	Instalación del anclaje antivuelco	
	Unidades de mesa	
	Unidades para piso	
	Espaciadores para unidades de montaje sobre el piso	
Capítulo 4	Descripción de producto	4-1
	Vista general de la serie OGH de hornos Heratherm	
	Vista general de los hornos Heratherm OGH-S	
	Vista general del horno Heratherm OMH	

T

	Vista general de los hornos Heratherm OMH-S	
	Dispositivos de seguridad	
	Entorno operativo	
	Interruptor de puerta	
	Sistema de sensórica y control	
	Comunicación de datos e interfaz de alarma	
	Interfaz RS 232	
	Contacto de alarma	
	Enchufe eléctrico	
	Fusibles	
	Componentes de la cámara	
	Cámara interna	
	Opciones de producto en las unidades de mesa	
	Puerta provista de ventanas para observación, e iluminación de la cámara	
	Puertos para acceso de ductos en unidades de mesa	
Capítulo 5	Puesta en marcha	5-1
	Unidades de mesa	
	Instalación del sistema de bandejas	
	Instalación inicial	
	Instalación del sistema de bandejas	
	Preparación de la cámara	
	Ûnidades de mesa	
	Instalación o extracción de los rieles de soporte (sólo series OMH y OMH-S)	
	Instalando y desinstalando el panel de aire posterior (sólo series OMH y OMH-S)	
	Instalando y desinstalando los paneles de aire laterales (sólo series OGH y OGH-S)	
	Nivelación del horno	
	Puesta en servicio de unidades para piso	
	Colocación y remoción de compartimientos de aire (Serie OMH)	
	Puesta en servicio, general	5-9
	Instalación de las abrazaderas de soporte de bandeja	5-9
	Instalación de las bandejas de malla metálica	
	Conexión de energía eléctrica	
	Conexión a la fuente suministro eléctrico	
	Conectar la Interfaz RS 232	
	Interconectando el horno con una computadora	
	Cableado del contacto de alarma	5-14
	Descripción funcional	5-14
	Especificaciones del relé de alarma	
	Ejemplo de conexión	
Capítulo 6	Operación	6-1
	Preparación del horno	6-1
	Inicio de la operación	6-1
Capítulo 7	Manipulación y control	7-1
	Encendido	7-6
	Conmutar el horno en Off / Apagado	7-6
	Valor teórico de temperatura	7-7
	Temporizador	7-9
	Configurar un temporizador de encendido o apagado de hora fija	7-11
	Configurar un temporizador semanal	7-12

	Detener un temporizador	
	Luz	
	Boost	
	Ventilador	
	Unidades de mesa	
	Unidades para piso	
	Regulador	
	Ajustes	
	Registro de errores (log)	
	Calibración	
	Fecha y Hora	
	Unidad de visualización de temperatura	
	Protección de proceso	
	Preferencias de programa	
	Configuración	
	Programación	
	End of Program	
	Programa de conveniencia ECO (sólo para unidades de mesa)	
	Instrucciones	
Canítulo 8	Anagado	8-1
Capitalo C	Apagado del horne	0 - 1 0 1
	Apagado del normo	
Capítulo 9	Limpieza y Desinfección	
	Limpieza	
	Limpieza de las superficies externas	
	Desinfección por limpieza y rociado	
	Preparación de la desinfección manual por limpieza y rociado	
	Predesinfección	
Capítulo 10	) Mantenimiento	
	Inspecciones y controles	10-1
	Intervalos de servicio	10-2
	Preparación para la calibración de la temperatura	10-2
	Procedimiento de medición por comparación	
	Procedimiento de calibración de la temperatura	10-3
	Reemplazo de la junta de sellado de la puerta	
	Reemplazo del cable de alimentación eléctrica	
	Devoluciones para reparación	
	1 1	-
Capítulo 11	Eliminación	11-1
	Vista general de los materiales usados	
Capítulo 12	2 Códigos de error	
Capítulo 13	B Datos técnicos	13-1
Capítulo 14	Piezas de repuesto y accesorios	
Capítulo 15	5 Dispositivo de registro	15-1
Capítulo 16	o Contacto	

Tabla de contenido

## **Figuras**

Figura 3-1 Unidades de horno para mesa, dimensiones y espacios libres necesarios	3-5
Figura 3-2 Unidades de horno para piso, dimensiones y espacios libres necesarios	3-6
Figura 3-3 Unidades de horno para piso, dimensiones y espacios libres necesarios	3-7
Figura 3-4 Puntos de izamiento	3-8
Figura 4-1 Serie Heratherm OGH 60 / OGH 100 / OGH 180, vista frontal 4	í- 2
Figura 4-2 Serie Heratherm OGH 60 / OGH 100 / OGH 180, vista posterior 4	í- 3
Figura 4-3 Vista frontal del horno Heratherm OGH-S 4	í- 5
Figura 4-4 Vista posterior del horno Heratherm OGH-S 4	í- 6
Figura 4-5 Serie Heratherm OMH 60/ OMH 100/ OMH 180: vista frontal 4	<b>í-</b> 7
Figura 4-6 Serie Heratherm OMH 60/ OMH 100/ OMH 180: vista posterior	í- 8
Figura 4-7 Serie Heratherm OMH 400 vista frontal	í- 9
Figura 4-8 Serie Heratherm OMH 400: vista posterior	<del>í</del> - 10
Figura 4-9 Serie Heratherm OMH 750 vista frontal	<b>í</b> - 11
Figura 4-10 Serie Heratherm OMH 750: vista posterior	<del>í</del> - 12
Figura 4-11 Vista frontal del horno Heratherm OMH-S 4	<b>í</b> - 14
Figura 4-12 Vista posterior del horno Heratherm OMH-S	<del>í</del> - 15
Figura 4-13 Ubicación del montaje del sensor - Serie OGH - Unidades de mesa	<del>í</del> - 16
Figura 4-14 Ubicación del montaje del sensor - Serie OMH - Unidades de mesa 4	<b>í</b> - 17
Figura 4-15 Ubicación del montaje del sensor - Serie OMH - Unidades de mesa 4	<b>í-</b> 17
Figura 4-16 Interfaces de señal y enchufe eléctrico 4	í- 18
Figura 4-17 Sistema de bandejas - Hornos OGH y OGH-S 4	í- 19
Figura 4-18 Sistema de bandejas - Hornos OMH y OMH-S 4	í- 20
Figura 4-19 Vista frontal del Paquete de visualización 4	<b>í-</b> 21
Figura 4-20 Puertos de acceso para conductos 4	<b>í</b> - 22
Figura 5-1 Deslizando el resorte de retención en el riel de soporte	5-2
Figura 5-2 Series OGH y OGH-S - Instalación del sistema de bandejas5	5-2
Figura 5-3 Series OMH y OMH-S - Instalación del sistema de bandejas	5-3
Figura 5-4 Instalación de riel de soporte	5-4
Figura 5-5 Instalando el panel de aire posterior	5-5
Figura 5-6 Desmontaje del panel inferior	5-6
Figura 5-7 Panel inferior retirado	5-6
Figura 5-8 Desmontaje el panel de aire lateral	5-7
Figura 5-9 Remoción de los perfiles de soporte izquierdo y derecho	5-8
Figura 5-10 Remoción del compartimiento de aire posterior, OMH 400/750 5	5-8
Figura 5-11 Instalación de los soportes de las bandejas5	5-9
Figura 5-12 Bandeja de malla metálica	5-10
Figura 5-13 Receptáculo para enchufe de red de CA5	5-12
Figura 7-1 Panel de control para los hornos Heratherm OGH, OMH, OGH-S y OMH-S	7-1
Figura 10-1 Reemplazo de la junta de la puerta (el ejemplo muestras un horno de la serie OMH) 1	10-4

F

#### Figuras

## Notas de seguridad

## Precauciones básicas de operación

Estas instrucciones de operación describen a los hornos Heratherm.

Los hornos Heratherm han sido fabricados con las técnicas más avanzadas habiéndoselas controlado minuciosamente antes de su envío para un funcionamiento perfecto. No obstante, el horno puede implicar peligros potenciales, particularmente al ser operada por personal entrenado inadecuadamente o al utilizarla con un propósito diferente al previsto. Por lo tanto, se debe observar lo siguiente con el objetivo de evitar accidentes:

- Los hornos Heratherm deben ser operados por personal profesional autorizado y adecuadamente entrenado.
- Los hornos Heratherm no deben ser operados hasta haberse leído y entendido completamente estas instrucciones de operación.
- Las siguientes instrucciones de uso, hojas de datos de seguridad aplicables, lineamientos de higiene en planta y las correspondientes normativas técnicas emitidas por el operador deben usarse para crear procedimientos escritos orientados al personal que trabaja con el dispositivo en cuestión, detallando:
  - las medidas de seguridad a implementarse al procesar agentes específicos,
  - las medidas de seguridad a implementarse en caso de accidente.
- Las tareas de reparación en el horno deben encargarse solamente a personal experto autorizado y entrenado.
- El contenido de estas instrucciones de operación está sujeto a cambios en cualquier momento y sin previo aviso.
- En cuanto a traducciones en lenguas extranjeras, la versión alemana de estas instrucciones de operación se considera vinculante.
- Mantenga estas instrucciones de uso cerca del horno de manera tal que, tanto las instrucciones de seguridad como la información importante estén siempre accesibles.
- Si usted encuentra problemas que no se hubiesen detallado adecuadamente en estas instrucciones de uso, póngase en contacto inmediatamente con Thermo Electron LED GmbH, por su propia seguridad.

## Reglas de seguridad operativa

Se deben considerar las siguientes normas al trabajar con hornos Heratherm:

- Respete los límites de peso de la muestra especificados para su horno Heratherm en su totalidad y especialmente en sus bandejas; ver "Datos técnicos" en página 13-1).
- No cargue la parte inferior del espacio de trabajo interior para evitar el riesgo de sobrecalentamiento de las muestras ubicadas allí.
- Ubique las muestras de manera uniforme en el espacio de trabajo, controlando de no ubicarlas muy cerca de las paredes internas para garantizar una distribución térmica uniforme.
- No cargue su horno Heratherm con sustancias que excedan la capacidad del dispositivo de laboratorio y del equipo de protección personal disponibles para brindar el grado de protección suficiente para usuarios y terceros.
- Controle la junta de goma de la puerta cada mes semana para una adecuada eficiencia de sellado y ante posibles averías.
- No procese muestras que contengan sustancias químicas peligrosas, que podrían liberarse en el aire ambiental a través de sellados defectuosos o que podrían causar corrosión u otras fallas en componentes de su horno Heratherm.

## Garantía

Thermo Electron LED GmbH garantiza la seguridad operativa y las funciones de los hornos únicamente bajo la siguiente condición:

- que el horno sea operado y mantenido exclusivamente acorde a su propósito previsto y tal como se describe en estas instrucciones operativas,
- que el horno no sea modificado,
- que solamente se usen repuestos y accesorios originales que hayan sido aprobados por Thermo Electron LED GmbH (los repuestos de terceros sin la aprobación de Thermo Scientific conllevan la anulación de la garantía limitada),
- que se realicen las inspecciones y el mantenimiento en los intervalos especificados,
- se realiza una prueba de verificación de instalación al poner en servicio el horno por primera vez, y luego reiteradamente después de cada tarea de inspección y reparación.

La garantía tiene validez a partir de la fecha de entrega del horno al operador.

## Aclaración de la información y símbolos de seguridad

# Notas y símbolos de seguridad utilizados a lo largo de estas instrucciones de operación

	A PELIGRO
	Indica una situación peligrosa la cual, de no evitársela, producirá la muerte o heridas de gravedad.
	ADVERTENCIA Indica una situación peligrosa la cual, de no evitársela, podría producir la muerte
	Indica una situación la cual, de no evitársela, podría dañar el equipo o la propiedad.
	<b>NOTA</b> Se usa para sugerencias y datos útiles referidos al uso.

### Símbolos adicionales para la información de seguridad

	¡Utilice guantes de seguridad!
	¡Utilice gafas protectoras de seguridad!
2	¡Líquidos perjudiciales!
Z	¡Choque eléctrico!
	¡Superficies calientes!
*	¡Peligro de incendio!
	¡Peligro de explosión!
	¡Peligro de sofocación!
	¡Riesgo de vuelco!

#### Símbolos en el horno



Respete las instrucciones operativas



Sello de la prueba VDE



Sello de conformidad de la CE: confirma la conformidad según los lineamientos de la UE

NC PE NO COM

Contacto de alarma

## Utilización prevista para el horno

#### Uso correcto

Las estufas y hornos de secado empotrados Heratherm OGH, OGH-S, OMH y OMH-S deben hacerse funcionar hasta la temperatura de trabajo máx. de 250 °C (482 °F), únicamente.

Los hornos son equipos de laboratorio para aplicaciones de calentamiento, equipados con control de temperatura de precisión.

Están diseñados para el tratamiento térmico de muestras o materiales a temperaturas de trabajo entre 50 °C (122 °F) y 330 °C (626 °F) (unidades de montaje sobre el piso: hasta 250 °C/482 °F) como por ejemplo: secado, envejecimiento, análisis, descomposición, incineración, oxidación, reducción y precalentamiento.

Los hornos se han diseñado para su instalación y operación en los siguientes entornos:

- tratamiento térmico;
- secado de materiales.

#### Uso incorrecto

Para evitar el riesgo de explosión, no cargue el horno con tejido, material o líquidos los cuales:

- sean fácilmente inflamables o explosivos;
- liberen vapores o polvo que formen mezclas combustibles o explosivas al contacto con el aire;
- desprendan venenos;
- crean una atmósfera húmeda;
- liberen polvillo;
- muestren reacciones exotérmicas;
- sean substancias pirotécnicas;
- superan la carga total especificada.

## Normas y directivas

El horno cumple con las siguientes normas y lineamientos:

- IEC EN 61010 1, IEC EN 61010 2 010
- Directiva para baja tensión 2006/95/CE
- Directiva CEM 2004/108/CE

Además, el horno cumple con muchas otras normas, reglamentos y directivas internacionales no incluidas aquí. En caso de dudas relacionadas con el cumplimiento de las normas, reglamentos y directivas nacionales vigentes en su país, contáctese con su representación comercial de Thermo Fisher Scientific.

## Envío del horno

## Embalaje

Los hornos Heratherm se entregan en una caja de embalaje resistente. Todos los materiales de embalaje pueden ser separados y son reutilizables:

Materiales de embalaje

Cartón de embalaje:	papel reciclable
Elementos de espuma:	espuma de poliestireno (sin CFC)
Pallet:	madera sin tratamiento químico
Lámina de embalaje:	polietileno
Cintas de embalar:	polipropileno

## Inspección de aceptación

Tras recibir el horno, controle inmediatamente el envío en cuanto a:

- integridad del suministro,
- posibles averías.

Si faltan componentes o se registran averías en el horno o en el embalaje, en particular daño causado por humedad y/o agua, informe inmediatamente al transportista así como al Soporte Técnico.

	Riesgo de lesiones		
En caso de haberse generado bordes cortantes en las zonas dañadas o en cualquier otro lugar del dispositivo, implemente todas las medidas precautorias necesarias para proteger la personal a cargo de manipular el horno. Por ejemplo, oblíguelos a que usen guantes protectores y todo otro equipo de protección personal.			

2

## Alcance del suministro

#### Hornos

Cantidad de componentes suministrados (piezas)	Serie OGH Serie OGH-S	Serie OMH-S Serie OMH
Bandeja de malla metálica	2	2
Riel de soporte para bandeja (sólo para unidades de mesa)	0	2
Soportes de las bandejas	4	4
Cordón de alimentación	1	1
Conector, contacto con separación galvánica	1	1
Resortes de sujeción (sólo para unidades de mesa)	0	2
Manual de operación	1	1
Guía de referencia resumida	1	1

## Instalación

## **Condiciones ambientales**

### Requisitos de ubicación

Las unidades de incubadoras, así como estufas y hornos de secado empotrados, deben hacerse funcionar con un sistema de escape de aire y una manguera de escape (únicamente utilice accesorios de Thermo originales).

Por razones de seguridad, el lugar de emplazamiento debe construirse de materiales no combustibles, según la norma DIN 4102.



Al instalar unidades empotradas, asegúrese de que el aire saliente se descargue correctamente fuera del lugar de emplazamiento.

#### Uso con sistemas de escape de aire

Para unidades empotradas, debe utilizarse una manguera de escape para alta temperatura y resistente a la corrosión (únicamente utilice accesorios de Thermo originales) la cual deberá conectarse al puerto de salida de aire con una llave de paso.

Al conectar varias unidades en fila empotradas a un sistema central de escape de aire (ver figura), debe preverse una llave de paso.



3

El horno debe operarse solamente en un lugar que reúna todos los requisitos de las condiciones ambientales abajo listados:

- Lugar interiorsin corrientes de aire y seco.
- La carga de polvo no debe exceder la categoría de contaminación 2 definida en la norma EN 61010-1. Se prohíbe usar el horno en una atmósfera con polvo conductor eléctrico.
- Debe respetarse la distancia mínima a superficies adyacentes en todos los lados (ver sección "Requisitos de espacio" en página 3-4.
- El cuarto de operaciones debe estar equipado con la ventilación adecuada.
- Materiales sólidos, nivelados, de superficie antiflama e ignífugos detrás del panel posterior del horno.
- La subestructura debe ser antivibratoria (apoyo sobre del piso, mesa de laboratorio) y capaz de soportar el peso muerto del horno y sus accesorios (en particular si se apilan dos equipos).
- Los hornos se han diseñados para operar a una altitud de hasta 2000 m sobre el nivel del mar.
- El rango de temperaturas es de 18 °C a 32 °C / 64,4° F a 89,6° F.
- Humedad relativa de hasta el 80% (máximo; preferentemente entre 60 y 70%), sin condensación.
- Debe evitarse condensación, por ejemplo, después de mover o transportar el equipo. En caso de condensación, espere hasta que la humedad se haya evaporado completamente antes de conectar el horno al suministro eléctrico y encenderlo.
- Evite la exposición solar directa.
- No deben ubicarse equipos que generen calor excesivo cerca del horno.
- Para evitar la operación de secado sin la provisión adecuada de aire puro controle que la entrada de aire (la cual puede estar equipada con un filtro de aire puro opcional) no se vea obstruida o bloqueada por ningún objeto cercano.
- Las fluctuaciones de la tensión de línea no deben exceder ±10 % del voltaje nominal.
- Los picos transitorios de tensión no deben exceder los valores usuales de la red de suministro eléctrico. El nivel de sobretensión transitoria nominal debe ser la tensión impulsiva resistiva de categoría de sobretensión II de IEC 60364-4-443.
- Considere instalar un interruptor dedicado por horno, aguas arriba de la instalación, para evitar múltiples caídas de equipos ante una falla eléctrica.

## Almacenamiento intermedio

Si el horno se ubica en almacenamiento intermedio, lo cual puede extenderse por un tiempo máximo de cuatro semanas, asegúrese de que la temperatura ambiente se mantenga entre los 20 °C y los 60 °C (68 °F a 140 °F) y que la humedad relativa máxima no exceda el 90%, sin condensación.

## Ventilación del recinto

El calor disipado por el horno en operación continua puede causar un cambio en el clima del cuarto.

- Por lo tanto sólo debe instalarse el horno en recintos con suficiente ventilación.
- No instale el horno en recesos del recinto carentes de ventilación.
- Cuando varios dispositivos deban colocarse en el mismo cuarto, se deberá proveer ventilación adicional en caso de ser necesario.
- Para evitar todo impacto del calor disipado por el horno en el clima ambiental, se debe ventilar el recinto por medio de un sistema de ventilación de grado de laboratorio, conforme a las normas de seguridad y sanitarias locales y nacionales vigentes, y de suficiente capacidad.
- Si hay una tendencia a las temperaturas excesivas en el cuarto de operaciones, asegúrese de prever una protección térmica que seccione el suministro eléctrico para mitigar el impacto de las condiciones de la temperatura sobreelevada.

## Uso con sistemas de escape de aire

Si se conecta el horno a un sistema de escape de aire, asegúrese de ajustar su flujo de aire como para que el patrón de distribución de temperatura en todos los puntos del horno no se modifique y se conserve el control de temperatura de precisión.

Antes de utilizar los ductos existentes en el edificio para extraer el aire caliente del horno, cerciórese de que dicha tubería de escape esté construida de materiales resistentes al calor, de tal modo que no se recaliente, funda y/o inflame.

Las superficies externas del horno y su tubería de escape de aire pueden calentarse, por lo que deberá conservarse el espaciado adecuado con respecto a paredes y techos, conforme lo especifique la norma constructiva y/o el reglamento de protección contra incendios.

• Marque el ducto de salida con señales de advertencia de superficie caliente adecuadas restrinja el acceso a tales conductos para evitar riesgos de lesiones por contacto con superficies calientes.

## Requisitos de espacio

En las unidades empotradas debe dejarse el siguiente espacio libre:



A, B, C y D, ver Instrucciones de Operación de la unidad.

E (mm/inch)	F (mm/inch)	G (mm/inch)	H (mm/inch)
100 / 4	50 / 2	80 / 3,2	30 / 1,2

### Instalación, unidades empotradas

IGS	Espacios libres, según Fig. 3, sin manguera de escape.
IMH/IMH-S	Espacios libres, según Fig. 3, cierre el puerto de aire con el tapón suministrado con la unidad.
OGS/OMS	Espacios libres, según Fig. 3, con espacio adicional para el funcionamiento con esclusa de aire, con sistema de escape de aire y manguera de escape de $Ø$ 40 mm (1,58")/1,5 m (59") (accesorio Thermo original), acortada a la longitud necesaria, instalada según Fig. 1 y 2.
OGH/OGH-S OMH/OMH-S	Espacios libres, según Fig. 3, con sistema de escape de aire y manguera de escape de Ø 40 mm (1,58")/1,5 m (59") (accesorio Thermo original), acortada a la longitud necesaria, instalada según Fig. 1 y 2, temperatura de trabajo máx. 250 °C (482 °F).



Una vez conectada la unidad empotrada a la alimentación eléctrica, evite dañar el cable eléctrico, deslizándolo a la posición de instalación.

Al instalar el horno, asegúrese de que la instalación y las conexiones de suministro queden fácilmente accesibles.

Los espacios laterales especificados representan distancias mínimas.

### Unidades de mesa



## Figura 3-1 Unidades de horno para mesa, dimensiones y espacios libres necesarios

#### Tabla 3-1 Dimensiones del horno Heratherm

Modelo	A (mm/inch <sup>*</sup> )	B (mm/inch)	C (mm/inch	D (mm/inch)
OMH 60	530 / 20,8	565 / 25,2	720 / 28,3	540 / 21,3
OMH 100	640 / 25,2	565 / 25,2	820 / 32,3	650 / 25,6
OMH 180	640 / 25,2	738 / 29,1	920 / 36,2	650 / 25,6
OMH 60-S	530 / 20,8	565 / 25,2	720 / 28,3	540 / 21,3
OMH 100-S	640 / 25,2	565 / 25,2	820 / 32,3	650 / 25,6
OMH 180-S	640 / 25,2	738 / 29,1	920 / 36,2	650 / 25,6

\*Las dimensiones en pulgadas son equivalentes redondeados, especificados solamente para información. Profundidad de manija /display (66 mm/2.6") no se incluye en la profundidad total especificada; altura de los pies ajustables (36 mm/1.4") no se incluye en la altura total especificada.

E (mm/inch)	F (mm/inch)	G (mm/inch)	H (mm/inch)
80 / 3,2	50 / 2	300 / 12	300 / 12

#### Tabla 3-2 Espacios libres requeridos

### Unidades para piso

#### Unidades de 400 litros



## Figura 3-2 Unidades de horno para piso, dimensiones y espacios libres necesarios

Tabla 3-3 Dimensiones del horno

Modelo	A (mm/inch)	B (mm/inch)	C (mm/inch)	D (mm/inch)
OMH 400	755 / 29,7	770 / 30,3	1655 / 65,2	810 / 31,9

\* Profundidad de manija /display (66 mm/2.6") no se incluye en la profundidad total especificada. Ancho de bisagra (23 mm) no se incluye en el ancho total especificado.

Tabla 3-4 Espacios libres requeridos

E (mm/inch)	F (mm/inch)	H (mm/inch)	l (mm/inch)
120 / 4,7	50 / 2,0	200 / 7,9	200 / 7,9

#### Unidades de 750 litros



Figura 3-3 Unidades de horno para piso, dimensiones y espacios libres necesarios

#### Tabla 3-5 Dimensiones del horno

Modelo	A (mm/inch)	B (mm/inch)	C (mm/inch)	D (mm/inch)
OMH 750	1215 / 47,8	770 / 30,3	1655 / 65,2	670 / 26,4

\* Profundidad de manija /display (66 mm/2.6") no se incluye en la profundidad total especificada. Ancho de bisagra (23 mm) no se incluye en el ancho total especificado.

Tabla 3-6	Espacios	libres	requeridos
-----------	----------	--------	------------

E (mm/inch)	F (mm/inch)	H (mm/inch)	l (mm/inch)
120 / 4,7	50 / 2,0	200 / 7,9	350 / 13,8

## Transporte

### Unidades de mesa

Al transportarla, no levante el horno de las puertas o de los componentes adosados a el horno como si se tratasen de puntos de izamiento.



#### Figura 3-4 Puntos de izamiento



#### Unidades para piso



Las unidades para piso se proveen equipadas con cuatro (4) ruedas. La palanca de liberación de rueda se sitúa encima de la palanca de bloqueo. Luego de posicionar la unidad en su lugar de ubicación, asegúrese de que las palancas de bloqueo estén aprisionando las ruedas.

Para garantizar el grado de estabilidad especificado por los requisitos de seguridad, las ruedas frontales deben girarse de tal modo que se orienten hacia adelante después de haber ubicado la unidad en su posición de instalación, con las palancas de bloqueo presionadas sobre dichas ruedas.



## Juego de apilado

El adaptador de apilado está disponible sólo para unidades de mesa.

Nº de Material	Descripción
50126665	Adaptador de apilado Heratherm 60L
50126666	Adaptador de apilado Heratherm 100L
50126667	Adaptador de apilado Heratherm 180 L

#### Alcance de suministro:

1 adaptador de apilado

1 anclaje antivuelco

1 Bolsa plástica con 2 pies de apilado y 2 tornillos M4x16 Torx.

#### Herramientas necesarias:

Destornillador para ranura 5,5x100 o destornillador Torx 20x100.

## Instalación de los pies de apilado

Quite los tapones ciegos izquierdo y derecho en la parte superior.

Instale los pies de apilado con los tornillos suministrados, usando un destornillador de ranura o Torx.

## Instalación del adaptador de apilado



Apile los equipos como sigue si utiliza un adaptador de apilado (las cifras indican el volumen del horno en litros):

- para 60/100/180 sobre 180 use el adaptador de apilado Heratherm 180L,
- para 60/100/100 sobre 180 use el adaptador de apilado Heratherm 100L,
- para 60/60/100 sobre 180 use el adaptador de apilado Heratherm 60L.

Para evitar que el horno superior se deslice o se caiga, deben cumplimentarse los siguientes requisitos antes de proceder a apilar los equipos:

- Solamente se pueden superponer dos unidades. Al apilar equipos con el mismo tipo de gabinete pero con diferente plantilla, el equipo con la plantilla más pequeña debe ubicarse arriba del otro que posee plantilla más grande.
- El incubador inferior debe estar correctamente nivelado.
- Asegúrese de usar el adaptador de apilado apropiado.
- Los pies niveladores en el incubador superior deben estar atornillados por completo.
- Los pies niveladores del dispositivo superior deben estar alineados con y ubicados exactamente sobre los receptáculos del adaptador de apilado.
- El anclaje antivuelco debe instalarse en la parte superior del equipo.

## Instalación del anclaje antivuelco

### Unidades de mesa

El anclaje antivuelco asegura el equipo superior de la pila a una parte firme del edificio. El anclaje antivuelco debe montarse del lado opuesto a las bisagras de la puerta.

Doble ambas pestañas de fijación del anclaje antivuelco hacia arriba y hacia abajo, respectivamente, en un ángulo aprox. 90°.



- 1. No use esta posición si la puerta pivota de ese lado. Las bisagras a mano derecha son la configuración estándar.
- 2. Posición preferida.
- 3. Posición alternativa. No use esta posición si la puerta pivota de ese lado.

Quite los tornillos del soporte. Use la posición preferida, dentro de lo posible.

Fije el anclaje antivuelco con la cara del soporte mirando hacia la unidad, hacia abajo (ver figura).

Posicione la unidad con un ángulo de aprox. 90° +/- 20% con respecto al anclaje antivuelco.

Cuide que los pies de apilado de la unidad ya estén ubicados en su lugar correcto sobre la unidad inferior, sobre el adaptador de apilado.

Fije el anclaje antivuelco a una parte firme del edificio.

### Unidades para piso



Las unidades Heratherm de montaje sobre el piso siempre deben estar fijadas a la pared mediante dos (2) soportes de retención por el lado externo izquierdo y derecho de la parte posterior de la misma.



Quite los tornillos.

Adose el extremo del soporte de retención de tal modo que mire hacia abajo de la unidad.

Alinee el dispositivo a aprox. 90°, +/-20° con respecto al soporte de retención.

Fije el soporte de retención a la pared.



¡Parte insegura del edificio!

Instale el anclaje antivuelco contra una parte firme del edificio, capaz de soportar cargas.

La instalación debe ser realizada únicamente por personal cualificado.

La vinculación con el edificio debe prever tornillos y tarugos apropiados para la consistencia del componente edilicio comprometido.

Adicionalmente, hay que tener en cuenta las siguientes indicaciones de precaución:

		Riesgo de sobrecalentamiento con equipos apilados
	¡No exceda la altura especificada de estibación para evitar así riesgo de sobrecalentamiento del gabinete externo o la pérdid control de temperatura a causa de la ventilación insuficiente!	

		Riesgo de inclinación o de caída de equipos apilados
	Debe tener cuidado cada vez conformen una unidad estab de apilado y los pies estén co superior puede inclinarse y o evitar lesiones personales y dispositivos apilados como u uno por uno, y luego superp	e que los dispositivos apilados no le, aún en el caso de que los receptáculos orrectamente acoplados. El dispositivo caerse si se lo transporta apilado. ¡Para del equipo, no intente trasladar los una unidad! Separe y traslade cada equipo, óngalos.
	Thermo Scientific no asume de otra procedencia que se h riesgo del usuario.	responsabilidad por cualquier dispositivo nubiese apilado; esto se realiza a cuenta y

## Espaciadores para unidades de montaje sobre el piso

El espaciador sobre el módulo eléctrico debe extraerse y fijarse en su lugar antes de colocar la unidad para piso en su lugar de emplazamiento.



- 1 Afloje los 2 tornillos
- 2 Extraiga el espaciador y deslice los tornillos dentro de los recesos
- 3 Ajuste ambos tornillos

## Descripción de producto

Esta sección describe los hornos Heratherm Advanced Protocol para aplicaciones de laboratorio sofisticadas, las que están disponibles en dos versiones diferentes según el nivel de seguridad:

- Hornos Heratherm Advanced Protocol serie OGH de convección por efecto de la gravedad (ver "Vista general de la serie OGH de hornos Heratherm" en página 4-1)
- Los hornos Heratherm de seguridad de convección natural y de protocolo avanzado de la serie OGH-S se proveen con características coincidentes con los equipos de la serie OGH, incluyendo funciones de seguridad adicionales para el cliente (ver "Vista general de los hornos Heratherm OGH-S" en página 4-3):
- Serie Heratherm OMH (OMH es la abreviatura de Oven with Mechanical convection); el horno Advanced Protocol de convección a ventilador viene con las siguientes funciones (ver "Vista general del horno Heratherm OMH" en página 4-6):
- Los hornos Heratherm de convección mecánica y de protocolo avanzado de la serie OMH-S se proveen con las siguientes características, incluyendo funciones de seguridad adicionales para el cliente (ver "Vista general de los hornos Heratherm OMH-S" en página 4-13):

## Vista general de la serie OGH de hornos Heratherm

La serie OGH (OGH es la abreviatura de <u>O</u>ven with <u>G</u>ravity convection) de hornos de convección natural y Advanced Protocol Heratherm viene con las siguientes funciones:

- control térmico de cámara de alta precisión, ajustable en pasos de un décimo de grado hasta los 330 °C / 626 °F);
- función aceleradora de calentamiento para un horno frío y vacío;
- un regulador eléctrico para ventilar la cámara, ajustable desde el panel de control;
- temporizador semanal, de hora fija y de cuenta regresiva para control de procesos cronometrados;
- dos bandejas de malla metálica;
- soporte de programa de usuario para control de proceso automático;
- un puerto de acceso para tubería, cables de sensores, etc.

Las características individuales de los hornos OGH se incluyen a continuación en las figuras.



#### Figura 4-1 Serie Heratherm OGH 60 / OGH 100 / OGH 180, vista frontal

- [1] Puerta externa
- [2] Ranura para el pestillo
- [3] Pestillo y manija de puerta
- [4] Bisagra de puerta, inferior
- [5] Pie nivelador
- [6] Placa de características
- [7] Deflector de aire, abajo
- [8] Deflector de aire, lateral
- [9]
- [10] Gancho de cierre de la puerta
- [11] Bandeja de malla metálica
- [12] Riel de soporte para bandeja de malla metálica
- [13] Junta estanca de la puerta
- [14] Receptáculo de apilado
- [15] Sensor de temperatura
- [16] Conducto de escape de aire
- [17] Puerto de acceso



#### Figura 4-2 Serie Heratherm OGH 60 / OGH 100 / OGH 180, vista posterior

- [1] Pestillo y manija de puerta
- [2] Panel de control
- [3] Receptáculo de apilado
- [4] Compuerta del puerto de acceso
- [5] Puerto de acceso
- [6] Espaciador
- [7] Escape de aire
- [8]
- [9] Deflector de escape de aire, de accionamiento eléctrico
- [10]
- [11] Compartimiento de electrónica
- [12] Pie nivelador

## Vista general de los hornos Heratherm OGH-S

Serie Heratherm OGH-S (OGH-S es la abreviatura de Oven with Gravity convection with exacting Safety requirements); el horno Advanced Protocol Security de convección natural viene equipado con las funciones que también se encuentran en los equipos de la serie OGH, incluyendo:

- control térmico de cámara de alta precisión, ajustable en pasos de un décimo de grado hasta los 330 °C / 626 °F);
- función aceleradora de calentamiento para un horno frío y vacío;
- un regulador eléctrico para ventilar la cámara, ajustable desde el panel de control;
- temporizador semanal, de hora fija y de cuenta regresiva para control de procesos cronometrados;
- dos bandejas de malla metálica;
- un puerto de acceso para tubería, cables de sensores, etc.;
- soporte de programa de usuario para control de proceso automático;

Adicionalmente, los hornos de la serie OGH-S porporcionan la siguiente funcionalidad extra:

- una puerta bloqueable para asegurar un proceso en curso contra el acceso no autorizado;
- un interruptor e indicador de puerta en el panel de control que indica si la puerta está abierta;

- una funcionalidad de protección de muestras que reduce la temperatura de la cámara a un nivel seguro cuando el controlador interno está sujeto a una condición de error;
- monitorización de temperatura inferior;
- está previsto el accesorio sensor de muestras;
- soporte de secado ECO (requiere del sensor de muestras opcional).



Las características individuales de los hornos OGH-S se incluyen a continuación en figura 4-5 y figura 4-6.

#### Figura 4-3 Vista frontal del horno Heratherm OGH-S

- [1] Puerta externa
- [2] Interruptor de puerta
- [3] Pestillo y manija de puerta
- [4] Bisagra de puerta, inferior
- [5] Pie nivelador
- [6] Placa de características
- [7] Deflector de aire, abajo
- [8] Deflector de aire, lateral

[9]

- [10] Gancho de cierre de la puerta
- [11] Bandeja de malla metálica
- [12] Riel de soporte para bandeja de malla metálica
- [13] Junta estanca de la puerta
- [14] Receptáculo de apilado
- [15] Sensor de temperatura
- [16] Conducto de escape de aire
- [17] Puerto de acceso



#### Figura 4-4 Vista posterior del horno Heratherm OGH-S

- [1] Pestillo y manija de puerta
- [2] Panel de control
- [3] Receptáculo de apilado
- [4] Compuerta del puerto de acceso
- [5] Puerto de acceso
- [6] Espaciador
- [7] Escape de aire
- [8] Conexión del sensor de muestras
- [9] Deflector de escape de aire, de accionamiento eléctrico

[10]

- [11] Compartimiento de electrónica
- [12] Pie nivelador

## Vista general del horno Heratherm OMH

La serie OMH (OMH es la abreviatura de Oven with Mechanical convection) de hornos Heratherm Advanced Protocol y convección mecánica viene con las siguientes funciones:

- control térmico de cámara de alta precisión, ajustable en pasos de un grado hasta 330 °C/626 °F (unidad de montaje sobre piso: hasta 250 °C (482 °F);
- función aceleradora de calentamiento para un horno frío y vacío (sólo para unidades de mesa);
- un ventilador de velocidad variable para optimizar la uniformidad térmica en todos los puntos de la cámara;
- un regulador eléctrico para ventilar la cámara, ajustable desde el panel de control;
- temporizador semanal, de hora fija y de cuenta regresiva para control de procesos cronometrados;
- dos bandejas de malla metálica;
- un puerto de acceso para tubería, cables de sensores, etc.
- soporte de programa de usuario para control de proceso automático;
- interruptor de puerta para unidades para piso
Las características individuales de los hornos OMH se incluyen a continuación en la figuras.



### Figura 4-5 Serie Heratherm OMH 60/ OMH 100/ OMH 180: vista frontal

- [1] Puerta externa
- [2] Ranura para el pestillo
- [3] Pestillo y manija de puerta
- [4] Bisagra de puerta, inferior
- [5] Pie nivelador
- [6] Placa de características
- [7] Deflector de aire, pieza superior
- [8] Riel de soporte para bandeja de malla metálica
- [9]
- [10] Gancho de cierre de la puerta
- [11] Bandeja de malla metálica
- [12] Riel de soporte para bandeja de malla metálica
- [13] Junta estanca de la puerta
- [14] Receptáculo de apilado
- [15] Resorte para riel de soporte
- [16] Sensor de temperatura
- [17] Conducto de escape de aire
- [18] Puerto de acceso



### Figura 4-6 Serie Heratherm OMH 60/ OMH 100/ OMH 180: vista posterior

- [1] Pestillo y manija de puerta
- [2] Panel de control
- [3] Receptáculo de apilado
- [4] Compuerta del puerto de acceso
- [5] Puerto de acceso
- [6] Conducto de escape de aire
- [7] Escape de aire
- [8] Ventilador
- [9] Deflector de escape de aire, de accionamiento eléctrico
- [10] -
- [11] Compartimiento de electrónica
- [12] Pie nivelador



#### Figura 4-7 Serie Heratherm OMH 400 vista frontal

- [1] Puerta externa
- [2] Pestillo de la puerta
- [3] Rueda de la unidad
- [4]

\_

- [5] Bandeja de malla metálica
- [6] Gancho de cierre de la puerta
- [7]
- [8] Sensor de temperatura
- [9] Puerto de acceso
- [10] -
- [11] -
- [12] -
- [13] \_
- [14] \_
- [15] \_
- [16] --
- [17] [18] -
- [19] -
- [20] -[21] Apertura del ventilador, deflector de aire



### Figura 4-8 Serie Heratherm OMH 400: vista posterior

- [1] Puerta externa
- [2]
- [3] Rueda de la unidad
- [4]

\_

-

-

\_

\_

- [5]
- [6]
- [7]
- [8]
- [9]
- [10] Conducto de escape de aire
- [11] Puerto de acceso
- [12] Anclaje antivuelco
- [13] Compartimiento de electrónica
- [14] Deflector de escape de aire, de accionamiento eléctrico
- [15] Bisagra, derecha
- [16] Manija
- [17] Display
- [18] Placa de características
- [19] -
- [20] -
- [21] -
- [22] Ventilador



### Figura 4-9 Serie Heratherm OMH 750 vista frontal

- [1] Puerta externa
- [2] Gancho de cierre de la puerta
- [3] Rueda de la unidad
- [4] Deflector de aire
- [5] Bandeja de malla metálica
- [6] Pestillo de la puerta
- [7]
- [8] Sensor de temperatura
- [9] Puerto de acceso
- [10] -
- [11] -
- [12] -
- [13] -
- [14] -
- [15] -
- [16] -
- [17] -
- [18] -
- [19] -
- [20] -
- [21] Apertura del ventilador, deflector de aire



### Figura 4-10 Serie Heratherm OMH 750: vista posterior

- [1] Puerta externa
- [2]
- [3] Rueda de la unidad
- [4]

-

-

-

\_

\_

- [5]
- [6]
- [7]
- [8]
- [9] -
- [10] -
- [11] Puerto de acceso
- [12] Anclaje antivuelco
- [13] Compartimiento de electrónica
- [14]
- [15] Bisagra, derecha
- [16] Manija

\_

- [17] Display
- [18] Placa de características
- [19] Bisagra, izquierda
- [20] Deflector de escape de aire, de accionamiento eléctrico
- [21] -
- [22] Ventilador

## Vista general de los hornos Heratherm OMH-S

La serie OMH-S (OMH-S es la abreviatura de Oven with Mechanical convection for High-end laboratory applications with exacting Safety requirements) de los hornos Heratherm Advanced Protocol Security de convección mecánica viene con las funciones que también se encuentran en los equipos OMH, incluyendo:

- control térmico de cámara de alta precisión, ajustable en pasos de un décimo de grado hasta los 330 °C / 626 °F);
- función aceleradora de calentamiento para un horno frío y vacío;
- un ventilador de velocidad variable para optimizar la uniformidad térmica en todos los puntos de la cámara;
- un regulador eléctrico para ventilar la cámara, ajustable desde el panel de control;
- temporizador semanal, de hora fija y de cuenta regresiva para control de procesos cronometrados;
- dos bandejas de malla metálica;
- un puerto de acceso para tubería, cables de sensores, etc.;
- soporte de programa de usuario para control de proceso automático;

Adicionalmente, los hornos de la serie OMH-S porporcionan la siguiente funcionalidad extra:

- una puerta bloqueable para asegurar un proceso en curso contra el acceso no autorizado;
- un interruptor e indicador de puerta en el panel de control que indica si la puerta está abierta;
- una funcionalidad de protección de muestras que reduce la temperatura de la cámara al punto de ajuste cuando el controlador interno está sujeto a una condición de error.
- Monitorización de temperatura inferior;
- está previsto el accesorio sensor de muestras;
- Soporte de secado ECO (requiere del sensor de muestras opcional).

Las características individuales de los hornos OMH-S se incluyen a continuación en figura 4-16 y figura 4-17.



### Figura 4-11 Vista frontal del horno Heratherm OMH-S

- [1] Puerta externa
- [2] Interruptor de puerta
- [3] Pestillo y manija de puerta
- [4] Bisagra de puerta, inferior
- [5] Pie nivelador
- [6] Placa de características
- [7] Deflector de aire, pieza superior
- [8] Soportes de las bandejas

[9]

\_

- [10] Gancho de cierre de la puerta
- [11] Bandeja de malla metálica
- [12] Riel de soporte para bandeja de malla metálica
- [13] Junta estanca de la puerta
- [14] Receptáculo de apilado
- [15] Resorte para riel de soporte
- [16] Sensor de temperatura
- [17] Conducto de escape de aire
- [18] Puerto de acceso, puerto de salida de aire



### Figura 4-12 Vista posterior del horno Heratherm OMH-S

- [1] Pestillo y manija de puerta
- [2] Panel de control
- [3] Receptáculo de apilado
- [4] Compuerta del puerto de acceso
- [5] Puerto de acceso
- [6] Espaciador
- [7] Escape de aire
- [8] Conexión del sensor de muestras
- [9] Deflector de escape de aire, de accionamiento eléctrico
- [10] Ventilador
- [11] Compartimiento de electrónica
- [12] Pie nivelador

## Dispositivos de seguridad

Los hornos están equipados con las siguientes características de seguridad:

- una funcionalidad de protección de muestra, que salvaguarda las muestras de su destrucción por sobrecalentamiento en caso de falla del controlador;
- una funcionalidad de corte para protección contra sobrecalentamiento, que detiene por completo el funcionamiento del horno cuando hay temperaturas excesivas en la cámara;
- fusibles duales calibrados a 16 amperes.

## Entorno operativo

Para asegurarse una operación sin problemas, la temperatura ambiente en el cuarto de operaciones debe ser de por lo menos 18 °C (64,4 °F). El sistema de calentamiento controla la temperatura en la cámara del horno de 50 °C/122 °F hasta el máximo de 330 °C/626 °F. (unidades para piso: 250 °C/482 °F)

## Interruptor de puerta

Los hornos Heratherm OGH 60/100/180-S y Heratherm OMH 60/100/180/400/750 vienen con un interruptor de puerta integrado dentro del mecanismo de cerrojo. Si se activa el interruptor de la puerta al abrirla, se suspenden las operaciones de calentamiento en la cámara y se ilumina un ícono en la ventana del display (ver D4 en figura 7-1 en página 7-1).

Si la puerta permanece abierta durante más de 30 segundos, se enciende una alarma acústica breve además del ícono en la ventana del visualizador. Si se deja la puerta abierta por más de 10 minutos, se activa una alarma audible, apareciendo en el display un mensaje de alarma diciendo "door open" (E001) y que se emite a través de la interfaz RS-232 con activación del relé de la alarma.

## Sistema de sensórica y control

El sensor tipo PT 100 para el control de la temperatura de la cámara y para la protección térmica [1] está montado en el panel superior del panel de cámara.



Figura 4-13 Ubicación del montaje del sensor - Serie OGH - Unidades de mesa



Figura 4-14 Ubicación del montaje del sensor - Serie OMH - Unidades de mesa



Figura 4-15 Ubicación del montaje del sensor - Serie OMH - Unidades de mesa

El sensor de temperatura de cámara provee de información para el controlador integrado del horno, el cual compara continuamente los valores medidos con el valor teórico especificado por el usuario y regula los calefactores según el resultado.



La unidad incluye una función de protección térmica que está preprogramada en fábrica y no es ajustable. La misma protege los cultivos dentro de la cámara contra el sobrecalentamiento: La protección térmica interviene ante la mínima violación del límite superior de temperatura, basándose en la temperatura del punto teórico definido, reduciendo automáticamente la temperatura en la cámara al valor teórico establecido por el usuario y permitiendo que el proceso de incubación continúe aún en caso de una falla en el controlador. Si se activa la protección térmica, aparece el mensaje (E111) "Temperatura excesiva" en la ventana del display y suena una alarma acústica.

Cuando el usuario recibe el mensaje de error, el ícono de alarma rojo (D4 en figura 7-1 en página 7-1) se ilumina y el ícono del valor teórico de temperatura (ver tabla 7-3 en página 7-4) es resaltado con un borde rojo que indica que ha intervenido la protección térmica.

## Comunicación de datos e interfaz de alarma

Todas las conexiones de señal se instalan en el panel de interfaz eléctrica en la parte posterior del horno.

### Interfaz RS 232

La interfaz RS 232 (a la izquierda en figura 4-16 abajo) puede ser usada para conectar las hornos al puerto de interfaz serie de una computadora para permitir la adquisición y documentación computarizada de los principales parámetros de funcionamiento (temperatura, códigos de error, etc.).



Figura 4-16 Interfaces de señal y enchufe eléctrico

## Contacto de alarma

Se puede conectar el horno a un sistema de alarma externa in situ, (como puede ser una central telefónica privada, un sistema de monitoreo de instalación, indicadores de alarma acústicos o visuales; ver "Cableado del contacto de alarma" en página 5-14). A tal fin, los equipos proporcionan un contacto de alarma con separación galvánica precableado (ver figura 4-16). El contacto de alarma se activa toda vez que ocurra un error en un lazo de control interno, en los circuitos eléctricos del horno o en el hardware.

## Enchufe eléctrico

El equipo se conecta a las líneas de suministro de CA a través del receptáculo IEC el cual acepta un cable con un enchufe estándar IEC (ver figura 4-16).

### **Fusibles**

Dos fusibles de 16 A de fusión lenta montados en el tablero electrónico principal del horno protegen al circuito interno del impacto de un consumo energético excesivo.

En la opción de producto "Ventana visora con iluminación de cámara de trabajo", las lámparas está protegidas con un fusible de 5 A en el compartimiento electrónico.



## NOTA

### Reemplazo del fusible

Los fusibles del dispositivo no admiten intervención del usuario para su reemplazo. Cuando el horno muestra los signos típicos de un fusible quemado (sin respuesta al presionar el botón On/Off, el panel de control no se enciende, no hay operación de calentamiento), contacte al servicio de atención de cliente para que le reemplacen los fusibles.

## Componentes de la cámara Cámara interna

Todos los componentes de la cámara se fabrican de acero inoxidable resistente a la corrosión y tienen una superficie lisa y fácil de limpiar. Todos los relieves tienen un radio amplio.

## Sistema de bandejas

Se provee el horno con dos bandejas de malla metálica. Los rieles de soporte de bandejas tienen una patrón alternativo de perforaciones ovales y circulares espaciadas regularmente cada 30 mm, lo que permite que los soportes de bandejas se inserten sin margen de error y con gran flexibilidad para alojar cualquier altura de recipiente de muestra que se requiera. Las bandejas tienen una protección integrada antivuelco y tope de extracción. Por detalles en cuanto al uso del sistema de bandejas, vea la sección"Puesta en marcha" en página 5-1.



- [3] Riel de soporte para bandeja de r[4] Bandeja de malla metálica
- [5] Deflector de aire, abajo

Figura 4-17 Sistema de bandejas - Hornos OGH y OGH-S



## Opciones de producto en las unidades de mesa

Esta sección describe las opciones disponibles para las estufas y hornos de secado Heratherm para aplicaciones avanzadas de laboratorio.

# Puerta provista de ventanas para observación, e iluminación de la cámara

Las estufas y hornos de secado Heratherm pueden equiparse con un paquete de visualización. En equipos de 60 y 100 litros de capacidad de cámara, este paquete consiste en una ventana visora junto a la iluminación interior para la cámara, mientras que los equipos de 180 litros se completan con dos ventanas visoras e iluminación de la cámara.

La iluminación de cámara no está diseñada para usarse como luminaria permanente. Úsela únicamente para realizar controles visuales breves del interior del equipo. El Paquete de visualización limita el rango operativo de temperaturas de la estufa y horno de secado a 250 °C (482 °F).

La sección "Luz" en página 7-16da instrucciones para encender y apagar la iluminación interior de la cámara.



### Figura 4-19 Vista frontal del Paquete de visualización

- [1] Panel de control
- [2] Pestillo y manija de puerta
- [3] Pie nivelador
- [4] Placa de características
- [5] Bisagra de puerta, inferior
- [6] Ventana visora

## Puertos para acceso de ductos en unidades de mesa

Las estufas y hornos de secado Heratherm pueden equiparse con puertos adicionales de acceso para conductos, tanto en paneles laterales como superiores.

Las opciones disponibles en cuanto a puertos de acceso para conductos se listan a continuación en tabla 4-1.

### Tabla 4-1 Puertos de acceso para conductos para estufas y hornos de secado Heratherm (opcional)

Modelo	Puerto montado en panel lateral, diám. en mm	Puerto montado en panel lateral, diám. en mm
OGH y OGH-S	19 o 53	24 o 58
OMH y OMH-S	24 o 58	24 o 58

Los puertos de acceso de conductos se instalan en lugares fijos sobre los paneles laterales y superior (ver figura 4-20).



### Figura 4-20 Puertos de acceso para conductos

- [1] Puerto de acceso de conductos en panel superior
- [2] Puerto de acceso de conductos en panel lateral
- [3] Tapón de sellado para puerto de acceso de conductos del panel lateral

Una vez insertados los cables, mangueras u otros conductos, estos puertos de acceso deben envolverse con las almohadillas de fibra resistentes al calor suministradas con el equipo, colocándole luego la tapa, para así sellar el puerto dentro de lo posible.



## Puesta en marcha

## Unidades de mesa

## Instalación del sistema de bandejas

No se necesitan herramientas para la instalación del sistema de bandejas. Los rieles de soporte se mantienen en posición por la acción de un resorte. Una vez insertas las abrazaderas de soporte de bandeja en los rieles, las bandejas perforadas simplemente se pueden empujar sobre los ganchos de soporte para completar la instalación.



## Instalación inicial

Los hornos de las series Heratherm OGH y OGH-S poseen los rieles de soporte de las bandejas integrados en los paneles de aire, los cuales se suministran preinstalados de fábrica.

Los hornos de las series Heratherm OMH y OMH-S están provistos de rieles de soporte separados, los cuales deberán instalarse como sigue:

- Quite la lámina protectora de los rieles de soporte.
- Introduzca el resorte de retención [1] dentro de la guía sobre el riel de soporte [2], asegurando que la protuberancia de bloqueo [3] del resorte de retención engrane firmemente con el orificio coincidente del riel de soporte.



Figura 5-1 Deslizando el resorte de retención en el riel de soporte

## Instalación del sistema de bandejas

Las siguientes ilustraciones muestran la ubicación de los elementos del sistema de bandejas.



[4] Bandeja de malla metálica[5] Deflector de aire, abajo

Figura 5-2 Series OGH y OGH-S - Instalación del sistema de bandejas



Figura 5-3 Series OMH y OMH-S - Instalación del sistema de bandejas

## Preparación de la cámara

Debe controlarse la limpieza de los siguientes componentes de la cámara y se los debe limpiar antes de su uso:

- rieles de soporte,
- soporte de las bandejas,
- bandejas de malla metálica,
- superficies de la cámara,
- sellos y juntas estancas de la cámara.



## Unidades de mesa

# Instalación o extracción de los rieles de soporte (sólo series OMH y OMH-S)



Figura 5-4 Instalación de riel de soporte

Los relieves en [2] y [7] sirven de guías laterales para los rieles de soporte, mientras que los relieves en [1] y [6] fijan los rieles de soporte en su lugar. Para una correcta instalación del riel de soporte, el resorte de retención [3] debe estar con su cara hacia arriba.

- 1. Posicione el riel de soporte [4] en el relieve inferior [6] e inclínelo, elevándolo hacia la pared lateral de la cámara de manera tal que el riel se posicione por encima de los dos relieves en [5] y [2].
- 2. Trabe el resorte de retención [3] detrás del relieve superior [1].
- 3. Para extraer los rieles de soporte, tire hacia abajo el botón del resorte de bloqueo retirándolo del relieve y extraiga el riel.

# Instalando y desinstalando el panel de aire posterior (sólo series OMH y OMH-S)

Los hornos de las series Heratherm OMH y OMH-S se entregan de fábrica con un panel de aire preinstalado en la pared posterior de la cámara (mientras que los hornos de las series OGH y OGH-S no poseen dicha prestación). Antes de poder retirar el panel de aire de la pared posterior, es necesario desmontar los rieles de soporte como se explicase más arriba.



Figura 5-5 Instalando el panel de aire posterior

1. Afloje los dos tornillos [1] que sostienen la chapa del panel de aire contra la pared posterior de la cámara.



- 2. Apriete los dos resortes de retención [2] por sus pestañas y tire de ellos hacia abajo, extrayéndolos del relieve, luego retire el panel de aire posterior.
- 3. Para una correcta instalación del panel de aire lateral, ambos resortes de retención [2] deben estar con su cara hacia arriba. Coloque el panel de aire sobre los relieves inferiores e inclínelo contra la pared posterior de la cámara.
- 4. Asegure ambos resortes de retención [2] en los relieves superiores.
- 5. Asegure el panel de aire contra la pared posterior de la cámara ajustando sendos tornillos en [1].

# Instalando y desinstalando los paneles de aire laterales (sólo series OGH y OGH-S)



La siguiente sección describe la instalación y desinstalación de los paneles de aire laterales.

### Figura 5-6 Desmontaje del panel inferior

1. Afloje sendos tornillos [1] en el panel inferior, luego retire todo el panel inferior levantándolo de sus dos relieves.



### Figura 5-7 Panel inferior retirado

2. Apriete los dos resortes de retención [1] por sus pestañas y tire de ellos hacia abajo, fuera del relieve, luego retire el panel de aire lateral.





Figura 5-8 Desmontaje el panel de aire lateral

- 3. Para una correcta instalación de los paneles de aire laterales, ambos resortes de retención [1] deben estar con su cara hacia arriba. Coloque el panel de aire lateral sobre los relieves inferiores e inclínelo contra la pared lateral de la cámara.
- 4. Asegure ambos resortes de retención [1] en los relieves superiores.
- 5. Vuelva a colocar el panel inferior en los relieves y asegúrelo ajustando sendos tornillos en [1].

### Nivelación del horno

- 1. Ubique un nivel de burbuja en el centro de la bandeja.
- 2. Ajuste manualmente los pies niveladores hasta que la bandeja de malla metálica quede alineada horizontalmente en todas las direcciones. Realice el ajuste de los pies niveladores desde la izquierda hacia la derecha y desde atrás hacia adelante.

## Puesta en servicio de unidades para piso

Colocación y remoción de compartimientos de aire (Serie OMH)



Figura 5-9 Remoción de los perfiles de soporte izquierdo y derecho

Afloje y extraiga los ocho (8) tornillos de los perfiles de soporte izquierdo y derecho y luego quite los compartimientos de aire laterales.





En el modelo OMH 400 afloje y quite los seis (6) tornillos del compartimiento de aire posterior, mientras que en el modelo OMH 750 afloje y quite los seis (6) tornillos de los compartimientos de aire superior e inferior posteriores, y a continuación retírelo/s.

## Puesta en servicio, general

## Instalación de las abrazaderas de soporte de bandeja

1. Inserte las abrazaderas de soporte de bandeja [3] en las perforaciones [1] del riel de soporte y del panel de aire, e inclínelo hacia abajo.



2. Asegúrese de que ambos elementos verticales [2] del soporte de estantería estén al ras de la superficie del riel de soporte y el panel de compartimiento de aire.



Figura 5-11 Instalación de los soportes de las bandejas

## Instalación de las bandejas de malla metálica



Figura 5-12 Bandeja de malla metálica

- 1. Empuje la bandeja de malla metálica sobre los soportes de la misma con el elemento de protección antivuelco [1] orientado hacia el panel trasero del horno. Los dispositivos de protección antivuelco [1] también sirven de guía para las bandejas de malla metálica.
- 2. Levante ligeramente la bandeja de malla metálica para que los topes de extracción [2] puedan deslizarse sobre los soportes de la bandeja.
- 3. Asegúrese de que las bandejas y sendos dispositivos antivuelco tengan libertad de movimiento sobre los soportes de bandeja.

## Conexión de energía eléctrica



El horno posee una carcasa con protección a tierra de clase I. Para reducir al mínimo el riesgo de choque eléctrico, utilice el cable de CA incluido para conectar el horno a una toma de energía correctamente instalada y con protección de tierra, con las siguientes funciones colocadas para cada horno:

- fusibles de fusión lenta T 16 A
- interruptor B 16



Ventajas de utilizar líneas de alimentación eléctrica independientes

Si bien se pueden hacer funcionar varios equipos desde una alimentación eléctrica común si no se excede la corriente nominal, recomendamos especialmente prever una línea de alimentación con su interruptor aguas arriba por horno para evitar la falla de múltiples equipos en caso de una avería eléctrica.

## Conexión a la fuente suministro eléctrico

- Antes de conectar el horno al suministro eléctrico, verifique si el voltaje de la red corresponde a las especificaciones de la placa de denominación en el frente del horno. ¡Si los niveles de tensión (V) y corriente (A) no son los requeridos, no conecte el horno a la fuente de energía!
- 2. Asegúrese de que el contacto de la alarma permanezca desconectado en ese momento. Si está conectado, desconéctelo ahora para evitar una falsa alarma en el terminal receptor. Volverá al contacto de la alarma más tarde una vez que haya avanzado en el procedimiento de inicio.
- 3. Conecte el conector IEC en el receptáculo de la parte trasera del horno.
- 4. Lleve el cable por un espacio que no atraviese tuberías de escape de aire, o conductos y pasillos. Al apilar equipos, mantenga el cable lejos de los puntos calientes del otro horno en la pila.
- 5. Conecte el enchufe con terminal de tierra del cable a un enchufe de pared correctamente puesto a tierra y protegido mediante disyuntor diferencial.
- 6. Asegúrese de que el cable no esté sujeto a fuerzas de tensión o compresión.



¡Asegúrese de que los enchufes de alimentación eléctrica permanezcan accesibles en todo momento!

Debe desconectarse rápidamente el cordón de alimentación ante una emergencia; ¡asegúrese de que sus tomas de corriente siempre estén accesibles y despejadas!



- [4] Enchufe eléctrico IEC
- [5] Fusible de 5 A (con la opción "ventanas para observación e iluminación de la cámara", solamente)

### Figura 5-13 Receptáculo para enchufe de red de CA



## **Conectar la Interfaz RS 232**



La interfaz de comunicación de datos RS 232 permite consultar la información de estado y los datos de temperatura del horno introduciendo comandos básicos en una ventana de terminal estándar del sistema operativo de su computadora. La interconexión requiere un cable RS 232 estándar con conectores de 9 pines y un pinout directo "1:1", sin cruce de líneas, el cual no se suministra con el horno.

## Interconectando el horno con una computadora

- 1. Apague la computadora.
- 2. Tienda el cable de interfaz serie a lo largo de un espacio que no cruce tuberías de escape de aire caliente, mesas, pasillos o conductos. Con equipos apilados, mantenga el cable de interfaz serie lejos de los puntos calientes del otro horno en la pila.
- Conecte un conector del cable de interfaz serie (longitud del cable, desde 5 hasta un máximo de 10 metros, no suministrado como componente estándar) hasta el enchufe etiquetado RS 232 en la sección de interfaz en la parte posterior del horno (ver "Interfaz RS 232" en página 4-18).
- 4. Conecte el segundo conector a un COM 1 /COM 2 no utilizado u otro puerto serie en la computadora.
- 5. Inicie el sistema de la computadora.
- 6. Inicie su programa de terminal estándar y parametrice la conexión como se indica a continuación:
  - 57600 bits por segundo

- 8 bits de datos
- 1 bit de stop
- Sin paridad
- 7. Una vez que su terminal indique que la comunicación serie se ha establecido exitosamente, introduzca cualquiera de los comandos en la lista tabla 5-1 de abajo, según el tipo de información que quiera averiguar.



8. Utilice la siguiente sintaxis genérica para los comandos:

#### ?:aaaa:bb::cc<CR>, donde:

- --- ?: identifica la línea de comando como consulta;
- **aaaa:** es la dirección del parámetro;
- **bb::** es un campo de respuesta que debe ser dejado en"00 por razones técnicas;
- cc es una suma de verificación para los comandos indicados en la tabla siguiente
- **<CR>** simboliza retorno de carro (Enter).

Recibirá una repuesta con el siguiente formato general:

#### !:aaaa:bb:XXXXX:cc<CR>, donde:

- l: identifica la línea de comando como una respuesta a una consulta;
- aaaa: es la dirección del parámetro introducida con la consulta;
- bb: es el número de bytes de dato en código hexadecimal, por ejemplo, 1F para el valor decimal 31;
- XXXXXX: es la información significativa de estado consultada;
- cc: es un check-sum (técnicamente, una XOR negada de todos los bytes retornados, excluyendo los bytes del check-sum y el carácter <CR>);
- **<CR>** simboliza retorno de carro (Enter).

Sintaxis del comando	Ejemplo de respuesta
Fecha y hora combinada	S
?:0010:00::c1	!:0010:11: <b>31.07.10;01:02:23</b> :e2 <b>Fecha Hora</b>
Sólo fecha	
?:0011:00::c0	!:0011:08: <b>31.07.10</b> :d2 <b>Fecha</b>
Sólo Hora	
?:0012:00::c3	l:0012:08: <b>01:02:23</b> :dc <b>Hora</b>
Valor Fijo de temperatura Temperatura de referenc	a (T1); Temperatura de cámara habitual (T2); ia (T3); Temperatura del sensor de muestra (T4)
?:3010:00::c2	!:3010:1f:+125.00;+124.96;+000.000;+000.00:b0 T1 T2 T3 T4

### Tabla 5-1 Comandos de terminal para los datos de consulta

## Cableado del contacto de alarma

NOTA	Trabajo especializado
Thermo Electron LED GmbH operatividad del equipo única son realizadas por personal	garantiza la seguridad operativa y la amente si la instalación y las reparaciones especializado.
¡La conexión del horno a u encomendarse únicamente telecomunicaciones, adecu	n sistema de alarma externo debe a personal experto de ingeniería o ladamente entrenado y autorizado!

## **Descripción funcional**

Cuando los errores y fallas del sistema tiene lugar en los circuitos de control térmico, se emite un mensaje de alarma al sistema de monitoreo de alarma conectado. El contacto con separación galvánica (contacto tipo changeover) ha sido diseñado para la configuración circuital especificada a continuación.



### Especificaciones del relé de alarma

Circuit	Voltage	External fusing
Circuits with system voltage	max. 250 V ~	max. 6 A
SELV circuits (cf.	25 V ~	max. 2 A
VDE 0100, Part 410)	60 V =	max.1A
SELV-E circuits (cf. VDE 0100,	50 V ~	max. 1 A
Part 410)	120 V =	max. 0.5 A

PRECAUCION

Consideraciones de compatibilidad eléctrica del contacto de alarma

Para evitar la sobrecarga y avería del contacto de alarma, verifique los parámetros de interfaz eléctrica del sistema de recepción de alarma en cuanto a su compatibilidad con las especificaciones del relé de alarma expuestas arriba.

## Ejemplo de conexión

Se provee el conector [5] para el cable de interfaz con el horno, como componente estándar. Las especificaciones para la tensión de trabajo y la protección fusible del circuito de alarma externo se incluyen en la tabla de la página precedente.

- 1. Cablee los conductores individuales [1] del cable de interfaz [4] como se muestra en el diagrama de cableado.
- 2. Lleve el cable de alarma por un espacio que no cruce tuberías de escape de aire caliente, mesas, pasillos o conductos. Con equipos apilados, mantenga el cable de interfaz serie lejos de los puntos calientes del otro horno en la pila.
- 3. Enchufe el conector del cable de interfaz del sistema de alarma en el puerto de interfaz [5] en el panel posterior del horno.



Ejemplo de conexión del relé de alarma

El diagrama circuital mostrado arriba representa la condición operativa sin fallas. En caso de una condición de error - incluyendo un corte energético - se cierran los contactos en el tramo entre contactos 1-4.

## Operación

## Preparación del horno

El horno no debe ser habilitado para su operación antes de haberse completado todas las tareas de arranque principales (ver "Puesta en marcha" en página 5-1).

### **Control del equipo**

Antes de la puesta en marcha inicial, controle el correcto funcionamiento de los siguientes componentes del horno:

- El sello de la puerta en el marco frontal no debe estar dañado.
- Los componentes de bandejas se deben instalar de manera segura.

## Inicio de la operación

1. Encienda el horno desde el panel de control.

- 2. Ajuste el valor de temperatura en el panel de control.
- 3. El controlador de temperatura comienza ahora a regular la cámara al valor teórico de temperatura especificado por el usuario.

Pera evitar cualquier riesgo de explosión o incendio		
<ul> <li>absténgase de cargar el horno con cualquiera de las sustancias incluidas en la lista de la sección "Uso incorrecto" en página 1-6</li> </ul>		
<ul> <li>compruebe que el aire ambiental esté libre de todo solvente</li> </ul>		
<ul> <li>no opere el horno en áreas con peligro de explosión</li> </ul>		

4. Cargue la cámara con muestras.

6

			_
Γ			
	7	1	
Ŀ			

#### 

### **Superficies calientes**

Al abrir la puerta, ésta puede volverse automáticamente.

El panel interno de la puerta externa así como las superficies de la cubierta externa, la estantería y la cámara se calientan extremadamente durante los ciclos de calentamiento del horno.

¡Al extraer muestras de un ciclo de calentamiento en proceso o de uno recientemente concluido, use siempre guantes de seguridad y cualquier otro equipo de protección de personal apropiado para evitar quemaduras con las superficies calientes!

	Riesgo de sobrecarga
La sobrecarga puede daña y/o el horno se inclinen cua destruyendo las muestras. de su bandeja, respete los especificados en "Datos té	r las bandejas o hacer que las mismas ando se extraen las bandejas, incluso Para evitar la sobrecarga del horno o límites de peso de muestras cnicos" en página 13-1.

NOTA	Carga correcta
Para asegurar suficiente circo de las muestras, no use más cámara. Los objetos volumino impedir la distribución del mis	ulación de aire y el calentamiento uniforme del 70% del área superficial máxima de la osos en la cámara pueden disipar el calor e smo.

## Manipulación y control

Los hornos de las series Heratherm OGH, OMH, OGH-S y OMH-S tienen una unidad de control montada en el panel frontal que consiste en un display multifuncional, cuatro botones de control y un botón on/off. Los cuatro botones de control interactúan con la ventana del display para permitir que los usuarios tengan acceso a todas las funciones de control y de ajuste de usuario del horno, incluyendo, por ejemplo, el valor de temperatura teórico, temporizador, así como una variedad de otras funciones.

En condiciones operativas normales el display muestra al usuario la temperatura de la cámara. La pantalla retorna a su modo predeterminado al completarse los ajustes o cuando no se introduzcan datos por un lapso de 30 segundos.

El gráfico siguiente muestra el panel de control del Heratherm OGH 60/100/180, Heratherm OMH 60/100/180/400/750, Heratherm OGH 60/100/180-S, y Heratherm OMH 60/100/180-S con todos sus elementos de visualización y control.





La tabla siguiente contiene descripciones breves de los botones en el panel de control (ítems K2 a K5 en figura 7-1).

Ícono	Ítem	Función
MENU	K1	<ul> <li>Botón de menú/Enter</li> <li>Presionando una vez la tecla: Activa el menú, resaltando el primer componente del menú con un contorno rojo.</li> <li>Presionando dos veces la llave: Seleccione el ítem de menú actualmente activo (resaltado con un borde rojo) y, según la función actualmente seleccionada, este botón permitirá cargar datos con ítem D2, D5 o D6.</li> <li>Presionando tres veces la tecla (habiendo cambiado un ajuste): Confirma una entrada o selección previa.</li> </ul>
<ul> <li></li> </ul>	К2	<ul> <li>Botón de flecha izquierda</li> <li>Después de la primera presión del botón Menú/Enter: <ul> <li>Desplaza la selección en el menú (ver ítem D3) al próximo ícono de la izquierda.</li> </ul> </li> <li>Habiendo seleccionado un ítem de menú: <ul> <li>Disminuye el valor de un parámetro ajustable, por ejemplo, el valor teórico de la temperatura en D5 o el nivel de velocidad del ventilador a D6. Manteniendo este botón presionado durante algunos segundos, cambia el valor seleccionado al modo rápido.</li> <li>Traslada la selección en el panel de display multifunción en D2 en el componente del menú actualmente activado a la opción siguiente a la izquierda, por ejemplo, del estado Off del temporizador al estado On.</li> </ul> </li> </ul>
	КЗ	<ul> <li>Botón On/Off</li> <li>Manteniendo presionado este botón durante 2 segundos conmuta el horno a off. La ventana del display desaparece, excepto el ícono del indicador de disposición en el sector de visualización de estado del ítem D4.</li> <li>El campo de visualización de temperatura D1 ofrece una lectura atenuada de la temperatura de la cámara, siempre que la temperatura exceda los 50 °C (122 °F).</li> </ul>
>	К4	<ul> <li>Botón de flecha derecha</li> <li>Después de la primera presión del botón Menú/Enter: <ul> <li>Traslada la selección en el menú (ver ítem D3) al ícono siguiente a la derecha.</li> </ul> </li> <li>Habiendo seleccionado un ítem de menú: <ul> <li>Incrementa el valor de un parámetro ajustable, por ejemplo, el valor teórico de la temperatura en D5 o el nivel de velocidad del ventilador a D6. Manteniendo este botón presionado durante algunos segundos, cambia el valor seleccionado al modo rápido.</li> <li>Mueva la selección del panel de diplay multifunción en D2 a la siguiente opción a la derecha - por ejemplo, del estado On del temporizador a Off.</li> </ul> </li> </ul>
ESC	K5	<b>Botón de escape</b> Retorna al nivel previo del menú o a la visualización estándar. Al abandonar el ítem de menú actual, se le puede sugerir al usua- rio guardar alguna configuración realizada previamente.

Tabla 7-1 Botones de control
La tabla siguiente contiene descripciones breves de las funciones de visualización del panel de control (ítems D1 a D6 en figura 7-1; los identificadores K1 a K4 se refieren a los botones mostrados en esta figura).

Función	Ítem	Función
24.0° 12:30°	D1	Campo de visualización que muestra la lectura permanente de la temperatura actual en la cámara ya sea en °C o °F (depen- diendo de las preferencias del usuario, ver "Unidad de visualiza- ción de temperatura" en página 7-34). Alternativamente, cuando el usuario configura el reloj integrado en el horno aparece en este lugar un mensaje de entrada tem- poral intermitente en el formato general <b>hh:mm</b> (horas:minutos, ambos con dos dígitos).
2010-03-29 12:59rm (On Off) 29.03.2010 12:59 heating relay error (E109)	D2	Panel de display multifuncional de cuatro líneas con campos para fecha y hora, un área de visualización de las opciones específicas del ítem del menú seleccionado, mensajes de alarma detallados con códigos de alarma, indicadores de avance para procesos continuos (por ejemplo, rampa de tempe- ratura controlada por programa), etc.
	D3	Barra de menú con representaciones iconizadas de parámetros ajustables. Se utiliza un contorno rojo para resaltar el ítem actual del <b>menú</b> , seleccionado con los botones Menú (K1) y de flechas <b>izquierda</b> (K2) y <b>derecha</b> (K4). Más abajo, en tabla 7-3 se incluyen descripciones breves de los ítems de menú individuales. <b>Nota</b> Si no se puede seleccionar un ítem, significa que la función no forma parte de la configuración equipada en su unidad.
	D4	<ul> <li>Área de visualización de estado con tres íconos representando estados específicos del horno (de izquierda a derecha):</li> <li>El ícono de puerta abierta aparece cuando la puerta frontal del horno está abierta o no ha sido cerrada correctamente (ver "Interruptor de puerta" en página 4-16).</li> <li>Nota El ícono de puerta abierta solamente es funcional con los equipos de la serie OGH-S y de la serie OMH-S y unidades para piso de la serie OMH.</li> <li>Al presentarse una condición de error, se encenderá el ícono de alarma rojo. Al mismo tiempo, el código de error actual parpadeará en el campo de visualización D2. Puede resetearse la alarma presionando el botón </li> </ul>
	<b>D</b> 5	puesto en off mediante el botón <b>On/Off</b> (ítem K3 en figura 7-1).
Set 899.3 °C	D5	Panel de ajustes etiquetado Set para establecer el valor de tem- peratura tanto en ℃ como en ℉ (según preferencia del usuario; ver "Unidad de visualización de temperatura" en página 7-34).

Tabla 7-2 Funciones del visualizador

Función	Ítem	Función
	D6	El gráfico izquierdo de los dos gráficos de barra verticales perte- nece al ícono de <b>ventilador</b> inmediato inferior y muestra el nivel actual de velocidad del ventilador. Gráfico de barra para ajuste de la velocidad del ventilador (en 5 pasos: 1 – 5) - 20% (1 signo encendido) - 40% (signos 1 y 2 encendidos) - 60% (signos 1 a 3 encendidos) - 80% (signos 1 a 4 encendidos) - 100% (signos 1 a 5 encendidos)
	D6	El gráfico derecho de los dos gráficos de barra verticales perte- nece al ícono del regulador inmediato inferior y muestra la posi- ción actual del regulador. Gráfico de barra de la posición del regulador (en cuatro pasos: 0 - 3) - posición 0 (regulador cerrado, todos los chevrones están apagados) - posición 1 (signos 1 y 2 encendidos) - posición 2 (signos 1 a 4 encendidos) - posición 3 (regulador totalmente abierto; signos 1 a 5 encendidos

Tabla 7-2 Funciones del visualizador

La tabla siguiente contiene descripciones breves de los íconos de barra de menú (ítem D3 en figura 7-1)

Tabla 7-3 Íconos de la barra de menú

Ícono	Función
	Valor teórico de temperatura Permite el cambio del valor teórico de la temperatura dentro del rango de temperaturas aceptable. El valor teórico puede cambiarse al presionar izquierda y derecha (ítem K2 o K4) y se puede, después de confirmar los cambios con el botón Menu/Enter (ítem K1), observar su efecto sobre la temperatura actual en el campo de visualización multifuncional de D1. Instrucciones: "Valor teórico de temperatura" en página 7-7,
	<b>Temporizador</b> Permite tener el horno encendido y/o apagado al finalizar un período de cuenta regresiva especificado por el usuario o por un tiempo fijo de encendido o apagado, o hacerla operar según un cronograma semanal completo de horarios de encendido y apagado. Cuando el usuario habilita un "on timer" el horno se apaga. Una mano giratoria en el ícono de <b>Temporizador</b> y el ícono encendido del indicador de disposición en el área de visualización de estado indica que el temporizador está funcionando. <b>Instrucciones:</b> "Temporizador" en página 7-9,
	Booster (refuerzo) (sólo para unidades de mesa) (Reforzador) una funcionalidad muy útil, pensada para lograr espontá- neamente un calentamiento rápido de una unidad fría y vacía (requiere una temperatura programada de al menos 150 ℃/302 ℉). Instrucciones: "Boost" en página 7-18,

Tabla 7-3	Íconos	de la	barra	de	menú

lcono	Función
	Ventilador Enciende el ventilador y habilita la elección de los niveles de velocidad descritos en D6. La configuración actual se muestra en el gráfico de barra situado directamente sobre el ícono y expresado como valor numérico en el panel de display en D2. Instrucciones: "Ventilador" en página 7-20.
	<b>Regulador</b> Abre y cierra el regulador en pasos graduados. La posición actual del regulador se muestra en el gráfico de barra situado directamente sobre el ícono y expresado como porcentaje en el panel de display en D2. <b>Instrucciones:</b> "Regulador" en página 7-24,
	<ul> <li>Ajustes</li> <li>Llama a un submenú con las siguientes funciones: <ul> <li>Acceso a la lectura del log de errores</li> <li>Calibración del horno</li> <li>Configurar la fecha y hora</li> <li>Conmuta el unidad del display de temperatura entre °C y °F</li> <li>Preferencias de ajustes para programas del usuario (program cycles / mode after end)</li> <li>Selecciona el modo de operación del temporizador (cuenta regresiva / hora fija en el día / temporizador semanal)</li> <li>Introducción de un código de control de configuración</li> </ul> </li> </ul>
/ <b>-</b> \_	<b>Programa (Instrucciones:</b> "Programación" en página 7-40) Permite ejecutar, crear, borrar, copiar y editar programas de usuario.

NOTA	Monitorización de inactividad del botón
Si no se presiona un botór ítem seleccionado del men seleccionada y retorna al n	) por más de 30 segunos en cualquier ú, el display abandona la pantalla nodo de visualización predeterminado.

# Encendido

1. Enchufe el cable eléctrico del horno en una toma de CA con adecuada puesta a tierra.

En la ventana del display, en el panel frontal, el ícono del indicador de disposición (ícono más a la derecha en el área de visualización de estado en D4 en figura 7-1 en página 7-1) se enciende.

(1)

2. Mantenga el botón **On/Off** presionado durante dos segundos.

Una rutina de inicialización se ejecutará después de que el horno haya sido encendido. Al completarse el test del sistema de inicialización, se encenderá el display y la temperatura actual de la cámara aparecerá en el campo de visualización de temperatura (ítem D1 en figura 7-1 en página 7-1). El horno ahora está listo para su uso.

# Conmutar el horno en Off / Apagado

1. Mantenga el botón **On/Off** presionado durante dos segundos.

La ventana del display desaparece, excepto el ícono del indicador de disposición (ícono más a la derecha en el área de visualización de estado en D4 en figura 7-1 en página 7-1) y la lectura de la temperatura de calor residual en caso de que la temperatura de la cámara sea todavía  $\geq$  50 °C/122 °F. El horno ahora se apaga.

2. Si es necesario, desenchufe el cable de alimentación de CA para apagar completamente el horno.

# Valor teórico de temperatura

Los hornos Heratherm permiten configurar la temperatura deseada de la cámara directamente con unas pocas maniobras con los botones. Después de confirmar el nuevo valor de temperatura en el panel de ajustes **Set** (ítem D5 en figura 7-1 en página 7-1), el usuario puede consultar el cambio de temperatura resultante en el campo de visualización de temperatura (ítem D1 en figura 7-1 en página 7-1).

Tras finalizar un proceso de secado o de calefacción Ud. también puede utilizar el valor teórico de temperatura para acelerar el proceso de enfriamiento. Al seleccionar temperaturas de una temperatura ambiente de +10 °C, la válvula de alimentación de aire debe estar abierta y no debe entrar calor adicional al aparato. No obstante, usted puede ajustar el horno para una temperatura deseada de 0 °C (32 °F), para acelerar el enfriamiento. Esto se logra en un solo paso, como se describe a continuación.

	Presione para activar la barra del menú, luego emplee para seleccionar el ícono <b>Temperatura</b> y presione para confirmar.
Set Set Set Set Set Set Set Set Set Set	En el panel de ajuste intermitente Set, presione o <>>>>>>>>>>>>>>>>>>>>>>>>>>>>>>>>>>>
	La pantalla regresa a su modo predeterminado. La temperatura real medida en la cámara y mostrada en el área de visualización de temperatura comienza a cambiar hasta que alcanza el valor teórico recientemente programado.

Tabla 7-4	Programación	del valor teórico	de la te	mperatura



Las estufas y hornos de secado empotrados Heratherm OGH, OGH-S, OMH y OMH-S deben hacerse funcionar hasta una temperatura de trabajo máx. de 250 °C (482 °F), únicamente.

	Al terminar un proceso de secado o calefacción,
	presione 🗮 para activar la barra del menú, luego
	utilice ≥ para seleccionar el ícono Temperatura y
	presione ਦ para confirmar.
Set	En el panel de ajustes intermitente Set, presione < para reducir el valor teórico de temperatura a 50 ºC
000.00	(122 °F), luego en < a 0 °C (32 °F) en un paso adicional. Cuando el display muestre 0 °C (32 °F),
	presione 🗮 para confirmar sus ajsutes.
	La pantalla regresa a su modo predeterminado.
	La temperatura real medida en la cámara y mostrada
	en el área de visualización de temperatura comienza a
	descender, mostrando el avance del proceso de
	enfriamiento.

#### Tabla 7-5 Uso del valor teórico de temperatura para acelerar el enfriamiento



# Temporizador

La función **Temporizador** de la barra de menú permite al usuario encender y apagar el horno según un cronograma. El temporizador acepta tres modos diferentes de operación, según las preferencias del usuario:

- Temporizador del tipo cuenta regresiva "on" u "off": Enciende (on) o apaga (off) el horno después de un tiempo especificado por el usuario. Las instrucciones para la configuración de las preferencias para esta opción se incluyen en tabla 7-6 más abajo, mientras que su uso como temporizador de encendido y de apagado se describe en tabla 7-7 en página 7-10 y tabla 7-8 en página 7-10, respectivamente.
- Temporizador de encendido y apagado de hora fija: Enciende o apaga el horno en una hora programada. Las instrucciones para la configuración de las preferencias para esta opción se incluyen en tabla 7-9, mientras que su uso como temporizador de encendido y de apagado se describe en tabla 7-10 en página 7-11 y tabla 7-11 en página 7-12, respectivamente.
- Temporizador semanal: Enciende o apaga el horno en horas programadas en días específicos de la semana. El proceso para la configuración de las preferencias para esta opción se describe en tabla 7-12, mientras que las instrucciones para la programación de las horas diarias de encendido y apagado se incluyen en tabla 7-13 en página 7-13.

La programación de un tiempo de encendido hace que el horno permanezca en modo standby hasta el horario de reinicio, mientras que el tiempo de apagado mantiene al dispositivo en funcionamiento antes de apagarse a la hora especificada por el usuario. El temporizador comienza a funcionar apenas el usuario confirma sus datos.

Las rutinas preprogramadas (programa de usuario, descontaminación) omiten el cronograma temporizado. Un temporizador no arrancará antes de que un programa concurrente termine de ejecutarse.

# Tabla 7-6 Programación del modo de operación del temporizador en modo cuenta regresiva

<b></b>	Presione 🗮 para activar la barra del menú, luego use
	para seleccionar el ícono de Ajustes y presione
	para confirmar.
2010-04-12 10:14рм	Presione ⋗ para conmutar al ítem de menú
Settings ∢ Timer ▶	Temporizador y confirme la selección con 💭.

# Tabla 7-6 Programación del modo de operación del temporizador en modo cuenta regresiva

2010-04-12 10:14 <sub>PM</sub>	Presione >>>>>>>>>>>>>>>>>>>>>>>>>>>>>>>>>>>>	
Tabla 7-7 Configurar u	un temporizador de apagado del tipo cuenta regresiva	
	Presione $\stackrel{\blacksquare}{\longleftarrow}$ para activar la barra del menú, luego use para seleccionar el ícono <b>Temporizador</b> y presione para confirmar.	
2010-03-29 12:59⊧м ∢On Off ▶	Presione — para seleccionar la opción de temporizador de apagado Off.	
2010-03-29 1:05 <sub>РМ</sub> Off-Timer 00:00	Establezca las horas y minutos que transcurrirán hasta que el horno se apague presionando 🔊 o < y luego confirme con 🗮.	
	La pantalla regresa a su modo predeterminado. En la barra de menú, el ícono de temporizador se ilumina y una mano gira sobre la cara del ícono.	
Tabla 7-8 Configurar un temporizador de encendido del tipo cuenta regresiva		

	Presione 💭 para activar la barra del menú, luego use para seleccionar el ícono Temporizador y presione para confirmar.
2010-03-29 12:59⊧м	Presione $\checkmark$ para seleccionar la opción de temporizador
∢On Off ▶	On y confírmela con $\overset{\text{MENU}}{\longleftarrow}$ .

2010-03-29 1:05թм	Establezca las horas y minutos que el horno esperará
On-Timer	para encenderse presionando Do I (luego
00:00	presione para confirmar.
	El horno se apaga. La pantalla se apaga, el ícono del temporizador se ilumina en la barra de menú, con una mano giratoria en la cara del ícono. Adicionalmente, el ícono del indicador de disposición se enciende.

#### Tabla 7-8 Configurar un temporizador de encendido del tipo cuenta regresiva

### Configurar un temporizador de encendido o apagado de hora fija

Tabla 7-9 Preconfigurar el modo de operación de temporizador de "Hora fija"

F	Presione 🗮 para activar la barra del menú, luego use
	para seleccionar el ícono de Ajustes y presione
	para confirmar.
	Presione ⋗ para conmutar al ítem de menú
	Temporizador y confirme la selección con 🛁.
2010-04-12 10:14⊧м Timer ∢Absolute ▶	Presione >>>>>>>>>>>>>>>>>>>>>>>>>>>>>>>>>>>>
	La pantalla regresa a su modo predeterminado.

#### Tabla 7-10 Configurar un temporizador de apagado de hora fija

	Presione ara activar la barra del menú, luego use para seleccionar el ícono <b>Temporizador</b> y presione
2010-03-29 12:59⊧м ∢On Off ▶	Presione para seleccionar la opción de temporizador de apagado Off.

2010-03-29 1:05рм	Configure el año, mes, día, horas y minutos mediante	
Off-Timer	🔊 o < , seguido de 🗮 para confirmar.	
2010-03-29 1:05PM		
	En la barra de menú, el ícono de temporizador se	
	ilumina y una mano gira sobre la cara del ícono.	
Tabla 7-11 Configurar (	Tabla 7-11 Configurar un temporizador de encendido de hora fija	
<u>(</u> )	Presione ਦ para activar la barra del menú, luego use	
	para seleccionar el ícono Temporizador y presione	
	para confirmar.	
2010-03-29 12:59рм	Presione < para seleccionar la opción de	
∢On Off ▶	temporizador $On$ y confírmela con $$ .	
2010-03-29 1:05рм	Configure el año, mes, día, horas y minutos mediante	
On-Timer 2010-03-29 1:05⊳м	🔊 o < , seguido de 🗮 para confirmar.	
	El horno se apaga.	
$\bigcirc$	La pantalla se apaga, el ícono del temporizador se ilumina	

#### Tabla 7-10 Configurar un temporizador de apagado de hora fija

### Configurar un temporizador semanal

#### Tabla 7-12 Programación del modo temporizador semanal

disposición se enciende.

F	Presione $\stackrel{\blacksquare}{\longleftarrow}$ para activar la barra del menú, luego use para seleccionar el ícono de Ajustes y presione para confirmar.
2010-04-12 10:14թм Settings ∢ Timer ▶	Presione <i>para conmutar al ítem de menú</i> Temporizador y confirme la selección con <i>.</i>

en la barra de menú, con una mano giratoria en la cara

del ícono. Adicionalmente, el ícono del indicador de

2010-04-12 10:14 <sub>РМ</sub> Timer	Presione 🔊 o < para preseleccionar el modo de operación de temporizador Semanal como su modo
♦ Weekly ►	elegido, luego confirme la selección con 🗮.
	La pantalla regresa a su modo predeterminado.

#### Tabla 7-12 Programación del modo temporizador semanal

#### Tabla 7-13 Configurar un temporizador semanal

	Presione 🗮 para activar la barra del menú, luego use
	para seleccionar el ícono Temporizador y presione
	para confirmar.
2010-03-29 12:59рм	En la pantalla de selección mostrada a la izquierda,
Timer ∢ On ∢	presione ⋗ para conmutar de On a la opción Edit.
2010-03-29 12:59⊧м Timer ∢ Edit ▶	Seleccione la opción Edición presionando
2010-03-29 12:59рм	Presione 🗮 para seleccionar la hora de encendido
♦ Monday ► On: 08:00AM	del lunes, la cual debe empezar a parpadear al
Оff: 06:00рм	seleccionaria (o continue nasta el día de la semana
	deseado presionando N, lo cual hará que la hora de
	Presione 🕐 o 🔇 para fijar las noras, luego continue
	con los minutos presionando 🛁.
	Use 🔊 o < para fijar los minutos 든 Luego, siga
	con el martes o cualquier otro día de la semana
	deseado presionando 🗮.
	Para evitar que el horno se encienda y vuelva a
	apagarse en un día específico, configure tanto las
	horas como los minutos en:

	Fije las horas programadas de encendido y apagado
2010-03-29 12:59рм	Tije las horas programadas de encendido y apagado
♦ Sunday ►	para cada dia de la semana por separado hasta el
Оп::ам Оff::РМ	domingo y presione . (Las horas de encendido y apagado para sábado y domingo están deshabilitadas por defecto.) Si no hay necesidad de fijar o cambiar las horas de encendido y/o apagado para un día específico,
	presione 💷 y ⋗ o < para regresar y avanzar al
	día previo y al siguiente respectivamente.
2010-03-29 12:59рм Save ?	Cuando se le sugiera guardar los cambios, presione
∢ No Yes ▶	<b>Nota</b> Este mensaje para guardar también aparece cuando presiona el botón Esc mientras trabaja en las configuraciones del temporizador semanal.
2010-03-29 12:59рм	Para habilitar el temporizador semanal presione ≕.
ilmer ∢ On ∢	Alternativamente, presione 🔊 y 🗮 para seleccionar
	la opción Apagado (Off) si el temporizador semanal recién
	guardado debiera activarse en una fecha posterior.
	En la barra de menú, el ícono de temporizador se
(-)	ilumina y una mano gira sobre la cara del ícono.

### Detener un temporizador

#### Tabla 7-14 Detener un temporizador de apagado antes de que expire

	Presione $\stackrel{\blacksquare}{\longleftarrow}$ para activar la barra del menú, luego use para seleccionar el ícono <b>Temporizador</b> y presione para confirmar.
2010-03-29 1:12PM Stop Timer ? ◀ No Yes ▶	Presione ara confirmar la selección Yes (Sí) por defecto.
	En la barra de menú, el ícono del temporizador desaparece. La pantalla regresa a su modo predeterminado.

#### Tabla 7-15 Detener un temporizador de encendido antes de que expire

	Para cancelar un temporizador de encendido preprogramado mientras el horno está apagado, mantenga el botón On/Off presionado por algunos segundos.
2010-03-29 1:12PM Stop Timer ? ◀ No Yes ▶	Al aparecer la pregunta Stop Timer ? para detener el temporizador, confirme la selección por defecto Yes presionando
<b>-</b>	En la barra de menú, el ícono del temporizador desaparece. La pantalla regresa a su modo predeterminado.

#### Tabla 7-16 Detener un temporizador semanal antes de que expire

()	Presione 💭 para activar la barra del menú, luego use
	para seleccionar el ícono Temporizador y presione
	para confirmar.
2010-03-29 12:59рм	La palabra encendida On aparece en el panel de
ilmer ∢ On ∢	Presione Presione
	presione 🛁.
2010-03-29 12:59թм Timer ∢ Off ▶	
2010-03-29 4:05 <sub>РМ</sub> Timer stopped!	El mensaje Timer stopped aparece como confirmación.
	En la barra de menú, el ícono del temporizador desaparece. La pantalla regresa a su modo predeterminado.



### Luz

Este ítem de menú apaga o enciende la luz interior en la cámara del equipo (requiere paquete de inspección opcional; sólo para unidades de mesa). Mientras la luz está encendida, su ícono **Luz** se ilumina en la barra del menú.

#### Tabla 7-17 Encender la luz

-œ́-	Presione ਦ para activar la barra del menú, luego use
Į	🔊 para seleccionar el ícono de Luz y presione 🗮
	para confirmar.
2010-03-31 1:34рм	En la pantalla de selección emergente, confirme la
∢ On Off ▶	opción seleccionada presionando 🗮.
2010-03-31 1:34рм	La opción seleccionada $\mathbf{On}$ permanece en el display por
On	algunos segundos para permitirle hacer el seguimiento de
	la acción.
` <b>`</b> <u><u><u></u><u></u><u></u><u></u><u></u><u></u><u></u><u></u><u></u><u></u><u></u><u></u><u></u><u></u><u></u><u></u><u></u><u></u></u></u>	La pantalla regresa a su modo predeterminado.
	El ícono Luz en la barra de menú se ilumina para
	indicar que la luz está encendida dentro del horno.

#### Tabla 7-18 Apagar la luz

-``@`-	Presione 🗮 para activar la barra del menú, luego use
	para seleccionar el ícono de Luz y presione
	para confirmar.
2010-03-31 1:34рм	En la pantalla de selección emergente, confirme la
∢On Off ▶	opción seleccionada presionando 🚛.
	La opción seleccionada Off permanece en el display por
2010-03-31 1:35рм	algunas sogundas para parmitirla basar al soguimiento do
Off	la acción.
	La pantalla regresa a su modo predeterminado.
[- <u>_</u> ]	El ícono $\mathbf{Luz}$ en la barra de menú se extingue para
	indicar que la luz está apagada dentro del horno.



# Restricciones en cuanto a la iluminación

La iluminación de cámara no está diseñada para usarse como luminaria permanente. Úsela únicamente para realizar controles visuales breves del interior del equipo.



#### Temperatura de trabajo máxima

El Paquete de visualización con ventanas visoras e iluminación de cámara limita el rango operativo de temperaturas de la estufa y horno de secado a 250°C (482 °F).



## Boost

La funcionalidad "boost" (sólo para unidades de mesa) permite acelerar el precalentamiento de un horno frío y vacío a un valor teórico de temperatura previamente ajustado (ver "Valor teórico de temperatura" en página 7-7) no inferior a 150 °C (302 °F). En modo boost, los calefactores de la unidad trabajan con su capacidad de calefacción máxima para aproximarse a la temperatura deseada con la mayor rapidez posible. Esta funcionalidad puede desactivar la protección por sobretemperatura por hasta 30 minutos y hacer que el calefactor trabaje con temperaturas pico transitorias. Una vez alcanzada la temperatura de preajuste, la unidad retorna al control normal de temperatura y el usuario puede proceder a cargar el horno con las muestras.

PRECAUCION	<b>i</b> ]

Riesgo de dañar las muestras!

Asegúrese de quitar de la cámara del horno el material a secar y todo otro ítem sensible a la temperatura antes de acelerar el calentamiento del horno. Temporariamente, tras haber calentado el horno rápidamente con la función boost, podrían permanecer elevadas temperaturas en la cámara, capaces de dañar las muestras.

#### Tabla 7-19 Activación de la función Boost

	Establezca el valor teórico de temperatura en al menos 150 °C/302 °F (ver "Valor teórico de temperatura" en pág. 7-9).
	Presione 🚅 para activar la barra del menú, luego use para seleccionar el ícono de Boost y presione 🚝 para confirmar.
2010-03-31 1:34 <sub>РМ</sub>	En la pantalla de selección emergente, confirme la opción seleccionada <b>On</b> presionando 🗮.
2010-03-31 1:34 <sub>РМ</sub> Оп	La opción seleccionada <b>On</b> permanece en el display por algunos segundos para permitirle hacer el seguimiento de la acción.

#### Tabla 7-19 Activación de la función Boost

	La pantalla regresa a su modo predeterminado.
	El ícono <b>Boost</b> en la barra de menú se ilumina para
	indicar que el modo de refuerzo está activado.
	Una vez alcanzado el valor teórico de temperatura
	preestablecido, el calefactor de refuerzo se apaga
	automáticamente, y el ícono Boost se va.
	¡Riesgo de quemaduras! Dependiendo del valor teórico
PRECAUCION	de temperatura elegido, pueden imperar elevadas
	temperaturas en la cámara después de que el horno
	estuvo operando con la funcionalidad "boost".
	¡Para evitar riesgos de quemadura los usuarios deben
	utilizar guantes protectores u otro equipamiento de
	protección personal al cargar la cámara del horno!

#### Tabla 7-20 Apagado prematuro del "booster"

	Presione 💭 para activar la barra del menú, luego use
	para seleccionar el ícono de Boost y presione para confirmar.
2010-03-31 1:34рм	En la pantalla de selección emergente, confirme la
<b>∢</b> On Off <b>}</b>	opción seleccionada Off presionando 🗮.
2010-03-31 1:35рм	La opción seleccionada Off permanece en el display
Off	por algunos segundos para permitirle hacer el
	seguimiento de la acción.
	La pantalla regresa a su modo predeterminado.
	El ícono <b>Boost</b> en la barra de menú se extingue para
	indicar que el modo de refuerzo fue apagado.



# Ventilador

Este componente del menú permite encender el ventilador que ventila la cámara, al mismo tiempo que se ajusta gradualmente su velocidad. El estado actual del ventilador se indicará a través del ícono iluminado **Ventilador** en la barra del menú y el gráfico de barra de cinco niveles (ver ítem D6 en figura 7-1 en página 7-1) para la regulación de la velocidad del ventilador, situado directamente sobre el ícono.

# Unidades de mesa

La unidad contiene un ventilador regulable; puede reducirse su velocidad (ajustarse) en 5 niveles:

- 20% (1 signo encendido)
- 40% (signos 1 y 2 encendidos)
- 60% (signos 1 a 3 encendidos)
- 80% (signos 1 a 4 encendidos)
- 100% (signos 1 a 5 encendidos)

Para proteger el horno del sobrecalentamiento, no obstante, el controlador integrado puede omitir el ajuste del usuario e incrementar la velocidad del ventilador automáticamente una vez que la temperatura de la cámara alcance el valor teórico. Una pérdida de velocidad puede disparar un error de ventilador (fan error).

#### Tabla 7-21 Encender el ventilador

	Presione $\stackrel{\texttt{MENU}}{\longleftarrow}$ para activar la barra del menú, luego use o o para seleccionar el ícono Ventilador y presione $\stackrel{\texttt{MENU}}{\longleftarrow}$ para confirmar.
31.03.2010 14:15	El diálogo de configuración que se muestra a la izquierda aparece en el panel de la pantalla multifunción, con el valor 20% encendido.

#### Tabla 7-21 Encender el ventilador

31.03.2010       14:15         ▲ 80%       ▲         ●       ●       ●         ●       ●       ●       ●         ●       ●       ●       ●         31.03.2010       14:15       ▲         ↓       100%       ▶       ▲         ●       ●       ●       ●       ●         ●       ●       ●       ●       ●	<ul> <li>Mantenga este ajuste o presione  las veces que sean necesarias para alcanzar el nivel deseado de velocidad del ventilador, confirmando con .</li> <li>El panel de pantalla multifuncional mostrará el nivel de velocidad del ventilador actual como porcentaje (20%, 40%, 60%, 80% o 100%). Adicionalmente, el número coincidente con el patrón en forma de V se iluminará en el gráfico de barras da la derecha.</li> </ul>
	La pantalla regresa a su modo predeterminado. El ícono de ventilador en la barra de menú se iluminará para indicar que el ventilador está funcionando.

#### Tabla 7-22 Ajuste de la velocidad del ventilador

	Presione para activar la barra del menú, luego use o para seleccionar el ícono Ventilador y presione para confirmar.
31.03.2010 14:15	El diálogo de ajustes aparece a la izquierda del panel multifunción, con la configuración de la velocidad del ventilador actual, intermitente.
31.03.2010 14:15	Cambie la velocidad del ventilador con 🔊 o < , luego presione 🗮 para confirmar.

#### Tabla 7-22 Ajuste de la velocidad del ventilador

ΝΟΤΑ	Si aparece un mensaje Heater prot. (protecc. de calefactor) al tratar de apagar el ventilador o bajar su velocidad:
	Los hornos de las series OMH y OMH-S emplean ventilación forzada, esto es, el ventilador no puede apagarse por completo y funcionará a una velocidad minima variable cada vez que se sobrepase cierto límite de temperatura. Todo intento de retornar la velocidad del ventilador al nivel de 0% será denegado como se indica en el mensaje <b>heater prot</b> en el panel visualizador multifunción.
	Idénticamente, el mensaje de protecc. de calefactor puede aparecer al presionar una vez el botón <a>o</a> o en cualquier paso sucesivo, al intentar reducir la velocidad del ventilador. Esto indica que la función de protección por temperatura ha tomado el control del ventilador, regulando su velocidad automáticamente hasta un nivel seguro que impida el daño térmico de los circuitos calefactores del horno.
	De otro modo, el control de velocidad de ventilador se ve limitado al nivel de velocidad admitido por el controlador o permanece inhabilitado por completo hasta que la temperatura de la cámara haya retornado a un nivel seguro.
$\leq$	La pantalla regresa a su modo predeterminado. El ícono de <b>Ventilador</b> permanece iluminado en la barra de menú.

## Unidades para piso

La unidad contiene un ventilador regulable; puede reducirse su velocidad (ajustarse) en 2 niveles:

Velocidad mínima del ventilador 60% (chevrones 1 a 3 encendidos)

Velocidad máxima del ventilador (chevrones 1 a 5 encendidos)

#### Tabla 7-23 Encender el ventilador

$\leq$	Presione 💭 para activar la barra del menú, luego use
	🔊 o < para seleccionar el ícono Ventilador y
	presione 🗮 para confirmar.
03 04 2012 12:45	El diálogo de ajustes aparece a la izquierda del panel
60%	multifunción, con la configuración de la velocidad del
	ventilador actual, intermitente.

#### Tabla 7-23 Encender el ventilador

03.04.2012 12:46 100%	<ul> <li>Presione Nas veces que sean necesarias para alcanzar el nivel 2 de velocidad del ventilador, confirmando con National.</li> <li>El panel de pantalla multifuncional mostrará el nivel de velocidad del ventilador actual como porcentaje (60% o 100%). Adicionalmente, el número coincidente con el patrón en forma de V se iluminará en el gráfico de barras da la derecha.</li> </ul>
	La pantalla regresa a su modo predeterminado. El ícono de ventilador en la barra de menú se iluminará para indicar que el ventilador está funcionando.

#### Tabla 7-24 Ajuste del ventilador

	Presione vara activar la barra del menú, luego use o vara seleccionar el ícono Ventilador y presione vara confirmar.
03.04.2012 12:45 100%	El diálogo de ajustes aparece a la izquierda del panel multifunción, con la configuración de la velocidad del ventilador actual, intermitente.
03.04.2012 12:45	Cambie la velocidad del ventilador con 🔊 o < , luego presione 🗮 para confirmar.
$\mathbf{x}$ / $\mathbf{x}$	La pantalla regresa a su modo predeterminado. El ventilador sigue funcionando hasta que la unidad se apaga.



# Regulador

Este ítem de menú permite abrir y cerrar gradualmente el regulador que ventila la cámara del horno. El estado actual del regulador se indicará a través del ícono iluminado **Regulador** en la barra del menú y el gráfico de barra de cuatro niveles (ver ítem D6 en figura 7-1 en página 7-1) para la posición angular del regulador, situada directamente por encima del ícono. Se puede ajustar el regulador en 4 niveles:

- Posición **0**: regulador cerrado todos los chevrones están apagados
- Posición 1: signos 1 y 2 encendidos
- Posición 2: signos 1 a 4 encendidos
- Posición 3: regulador totalmente abierto; signos 1 a 5 encendidos

#### Tabla 7-25 Apertura del regulador

	Presione 🗮 para activar la barra del menú, luego use
	🔊 o < para seleccionar el ícono Regulador y
	presione 🗮 para confirmar.
	El diálogo de configuración que se muestra a la izquierda aparece en el panel de la pantalla multifunción, con el valor inicial <b>0</b> encendido.
2010-04-10 2:37PM	<ul> <li>Presione Nas veces que sean necesarias para alcanzar el ángulo del regulador deseado, luego confirme con New.</li> <li>El panel de pantalla multifunción mostrará el ajuste actual del regulador como un valor entre 0 y 3. Adicionalmente, el número coincidente con el patrón en forma de V se iluminará en el gráfico de barras da la derecha.</li> </ul>
	La pantalla regresa a su modo predeterminado. El ícono <b>Regulador</b> en la barra de menú se ilumina para indicar que el regulador está abierto y cuánto.

×	Presione 💭 para activar la barra del menú, luego use
	▶ o 🔇 para seleccionar el ícono Regulador y
	presione 🛁 para confirmar.
10.04.2010 14:37	El diálogo de ajustes aparece a la izquierda del panel multifunción, con la configuración de la posición del regulador actual, intermitente.
2010-04-10 2:37рм	Use ⋗ o < para cambiar la posición del regulador,
	luego confírmela con 🛁.
	Para cerrarlo, fije la posición del mismo en 0 mediante
	< , luego confirme con 🗮.
	La pantalla regresa a su modo predeterminado.
	Si acaba de reajustar la posición del regulador, el ícono
	Regulador quedará iluminado en la barra del menú.
	Si decide cerrar por completo el regulador, el ícono de
	Regulador en la barra del menú se apagará.

#### Tabla 7-26 Ajuste de la posición del regulador o cierre del regulador



## Ajustes

El componente del menú **Ajustes** abre un submenú con varios comandos para visualizar la información de estado general en la unidad y para realizar preferencias de ajustes de funcionamiento del horno o su ventana visualizadora:

- Acceso a la lectura del log de errores
- Calibración del horno
- Configurar la fecha y hora
- Conmuta el unidad del display de temperatura entre °C y °F
- Encendiendo y apagando la protección de proceso
- Preferencias de ajustes para programas del usuario (program cycles / mode after end)
- Selecciona el modo de operación del temporizador (cuenta regresiva / hora fija en el día / temporizador semanal)
- Introducción de un código de control de configuración

Las instrucciones para el uso de estas funcionalidades se brindan en lo sucesivo.

### Registro de errores (log)

Cuando los usuarios llaman a Atención al Cliente para hacer consultas, el agente de les puede solicitar información del registro de errores del horno. Se accede a esta memoria interna con el ítem de menú **Settings -> Error**. Esto permite al usuario recorrer los mensajes de alarma más recientes causados por errores de hardware o de los lazos de control. Se muestra cada error con la fecha y la hora de su ocurrencia, una breve descripción en texto legible y un código de error interno.

Los códigos de error e instrucciones para despejar las condiciones de alarma aparecen en la sección "Códigos de error" en página 12-1.

F	Presione 逆 para activar la barra del menú, luego use o 🔇 para seleccionar el ícono Ajustes y presione para confirmar.
2010-04-06 1:33рм ∢ Error ▶	Presione ਦ para seleccionar el ítem de Error con el submenú Ajustes.
2010-04-06 1:36рм Error 0 2010-04-06 1:31рм Fan Error (E009)	La primera entrada del registro de error se muestra, numerada "0". Esta entrada representa el error registrado más recientemente.

#### Tabla 7-27 Leer el Registro de errores

#### Tabla 7-27 Leer el Registro de errores

2010-04-06 1:37рм Error 1 2010-04-06 1:34рм Heat Relay (E109)	Presione >>>>>>>>>>>>>>>>>>>>>>>>>>>>>>>>>>>>
F	Para salir del registro de errores y retornar al modo de pantalla normal presione Esc. El ícono Ajustes en la barra del menú desaparecerá.

### Calibración

El componente del menú **Ajustes -> Calibración** permite al usuario iniciar un proceso de calibración de temperatura (ver "Procedimiento de calibración de la temperatura" en página 10-3) para los sensores de temperatura integrados y seleccionar si la calibración se hará manual o automáticamente:

- La opción **Manual** permite introducir directamente una temperatura absoluta, tal como se la mide, por ejemplo usando un sensor de referencia externo.
- La función Eco (opcional) es necesaria para implementar el programa de conveniencia Eco. Calcula la diferencia de temperatura espacial entre la temperatura dominante en el lugar de la muestra (tal como la registra el sensor de muestras, accesorio a solicitarse por separado) y la temperatura actual de la cámara (tal como la mide el sensor integrado del horno). El controlador integrado usa el resultado para calcular un límite de temperatura que marque el final del proceso de secado. El preajuste de temperatura en el horno durante el proceso de calibración Eco debe ser igual al máximo valor teórico de temperatura para el proceso de secado. El programa de conveniencia Eco, así como su utilización, se describen con más detalle en la sección "Programa de conveniencia ECO (sólo para unidades de mesa)" en página 7-40.
- La función "Sample" (opcional) permite calibrar la temperatura de un sensor (accesorio, debe encargarse aparte) conectado al horno introduciendo directamente un valor absoluto medido con un dispositivo de medición externo.

NOTA	Requisitos previos para la calibración
Mantenga las condiciones ar especificados del horno y as completamente cerrado ante	nbientales dentro de los límites egúrese de que el regulador esté s de iniciar la calibración.
Condiciones ambientales van afectar el resultado de la ruti a un desajuste del controlado control de temperatura.	riables y/o un regulador abierto pueden na de calibración, lo cual puede conducir or y a una operación no confiable del

<b>F</b>	Presione 💭 para activar la barra del menú, luego use
	🔊 o < para seleccionar el ícono Ajustes y presione
	para confirmar.
2010-04-06 1:33рм	El ítem de menú Error del submenú Ajustes aparece en la pantalla multifunción.
2010-04-06 1:33рм	Presione ⋗ para conmutar al ítem de menú
	Calibración y confirme la selección con 🗾.
2010-04-06 1:33рм	En la pantalla de selección Calibración, presione 🗮
Calibration ∢ Manual ▶	para elegir la opción preseleccionada Manual.
2010-04-06 1:33рм	En el diálogo de ajustes emergente, fije la temperatura
Calibration	medida con el sensor de referencia externa usando
36.9°C	o v confirme sus datos con .
	El valor recientemente ingresado será almacenado y
	usado para calibrar los sensores de temperatura
	internos con el valor medido por el sensor de referencia.
	La pantalla regresa a su modo predeterminado.
	El locito Ajustes de la barra de menu se apagara.

#### Tabla 7-28 Introduzca manualmente la temperatura de referencia de calibración

#### Tabla 7-29 Calibración ECO de temperatura (sólo para unidades de mesa)

	Coloque el sensor de temperatura en el lugar donde se
	colocará la muestra. Guíe el cable del sensor a través
	del puerto de acceso en el panel posterior del horno y
	conéctelo al conector correspondiente de la pared
	posterior del equipo.
<b>F</b>	Presione 픚 para activar la barra del menú, luego use
	🔊 o < para seleccionar el ícono Ajustes y presione
	para confirmar.

2010-04-06 1:33рм	El ítem de menú Error del submenú Ajustes aparece en la pantalla multifunción.
2010-04-06 1:33рм	Presione <i>para conmutar al ítem de menú</i> Calibración y confirme la selección con <i>L</i> .
2010-04-06 1:33⊧м Calibration ♦ Есо ▶	En la pantalla de selección <b>Calibration</b> presione para seleccionar la opción Eco y confirme dicha selección con .
2010-04-06 1:33PM Calibration Eco ♦ Off ► 2010-04-06 1:33PM Calibration Eco ♦ On ►	El estado operativo actual aparece en el panel visualizador, representado por la palabra intermitente Off. Presione Derivativa para conmutar al estado On, luego - para confirmar. La opción seleccionada On permanece en el panel visualizador por unos pocos segundos permitiéndole seguir la secuencia.
2010-04-06 1:35рм Eco Calib. Running	El mensaje de confirmación Eco Calib. Running aparece en el panel visualizador multifunción para indicar que la autocalibración está en proceso.
×	El ícono Ajustes de la barra de menú se apagará. Una vez cumplimentado el proceso de calibración, la visualización retorna a su modo predeterminado.

Tabla 7-29 Calibración ECO de temperatura (sólo para unidades de mesa)

el sensor de	e muestras
	Coloque el sensor de temperatura en el lugar donde se colocará la muestra. Lleve el cable del sensor a través el puerto de acceso de conductos a un dispositivo externo de medición de temperatura.
	de muestras se haya estabilizado y suministre una lectura constante de temperatura.
<b>&gt;</b>	Presione — para activar la barra del menú, luego use o o para seleccionar el ícono Ajustes y presione — para confirmar.
2010-04-06 1:33рм ♦ Error ▶	El ítem de menú Error del submenú Ajustes aparece en la pantalla multifunción.
2010-04-06 1:33рм	Presione ⋗ para conmutar al ítem de menú
<ul> <li>♦ Calibration </li> </ul>	Calibración y confirme la selección con 픋.
2010-04-06 1:33թм Calibration ∢Sample ▶	En la pantalla de selección Calibración, presione 🛁 para elegir la opción Sample.
2010-04-06 1:33рм Calibration 36.9°С	En el diálogo de ajustes emergente, fije la temperatura medida con el sensor de referencia externa usando o  y confirme sus datos con  .
	El valor recientemente ingresado será almacenado y usado para calibrar el controlador integrado con el valor medido por el sensor de muestras. La pantalla regresa a su modo predeterminado. El ícono Ajustes de la barra de menú se apagará.

### Fecha y Hora

La opción **Ajustes -> Hora / Fecha** permite conmutar entre los formatos de pantalla de hora y fecha internacional y para configurar la hora y la fecha del reloj interno. Se puede elegir entre dos formatos de pantalla:

- El formato de fecha europeo *DD.MM.YYYY* y el formato horario de 24 horas. Ejemplo: 07.04.2010 and 15:05.
- El formato de fecha de EEUU estándar *YYYY-MM-DD* y el formato de 12 horas con el sufijo *AM/PM*. Ejemplo: 2010-04-07 and 3:05 PM.

#### Tabla 7-31 Configurar el formato de la fecha

F	Presione 🗮 para activar la barra del menú, luego use
	🔊 o < para seleccionar el ícono Ajustes y presione
	para confirmar.
2010-04-06 1:33рм	El ítem de menú Error del submenú Ajustes aparece en la pantalla multifunción.
♦ Error >	
2010-04-07 3:05рм	Presione ≥ para conmutar al ítem de menú
▲ Time / Date ▶	Hora/Fecha y confirme la selección con 🛁.
2010-04-07 3:05рм	El ítem de menú <b>Date</b> parpadea en el panel de pantalla multifuncional.
Time / Date ▲ Date ▶	Elija la opción preseleccionada Date presionando 📇.
	El campo de la fecha comenzará a parpadear en la
	esquina superior izquierda del panel multifunción.
07.04.2010 3:05рм	Presione ⋗ o < para conmutar al formato de fecha
Time / Date	deseado DD.MM.YYYY o YYYY-MM-DD y confirme la
	selección con 🗮.
	El campo de fecha en la esquina superior izquierda
	campiara su apariencia segun su selección (y dejara de parpadear).
	La pantalla regresa a su modo predeterminado.
	El ícono Ajustes de la barra de menú se apagará.

F	Presione 🗮 para activar la barra del menú, luego use
	🔊 o < para seleccionar el ícono Ajustes y presione
	para confirmar.
2010-04-06 1:33рм	El ítem de menú Error del submenú Ajustes aparece en la pantalla multifunción.
2010-04-07 3:05рм	Presione ⋗ para conmutar al ítem de menú
◀ Time / Date ▶	Hora/Fecha y confirme la selección con 🛁.
2010-04-07 3:05рм	El ítem de menú Date parpadea en el panel de pantalla
Time / Date ◀ Date ▶	multifuncional.
2010-04-07 3:05рм Time / Date	En el ítem de menú intermitente <b>Date</b> , presione <b>&gt;</b> para conmutar a la opción <b>Time</b> y confirme la selección
▲ Time ▶	
	con 🛁. El componente de menú Time destella en el panel de
◆ Time ▶	con —. El componente de menú <b>Time</b> destella en el panel de pantalla multifunción, junto al campo de hora en la esquina superior derecha.
<ul> <li>▲ Time ▶</li> <li>2010-04-07 15:05</li> <li>Time / Date</li> <li>▲ Time ▶</li> </ul>	<ul> <li>con .</li> <li>El componente de menú Time destella en el panel de pantalla multifunción, junto al campo de hora en la esquina superior derecha.</li> <li>Presione &gt;&gt; o &lt;&lt; para conmutar al formato de hora deseado hh:mm o hh:mm AM/PM y confirme la</li> </ul>
<ul> <li>▲ Time ▶</li> <li>2010-04-07 15:05</li> <li>Time / Date</li> <li>▲ Time ▶</li> </ul>	<ul> <li>con .</li> <li>El componente de menú Time destella en el panel de pantalla multifunción, junto al campo de hora en la esquina superior derecha.</li> <li>Presione &gt;&gt; o &lt;&gt; para conmutar al formato de hora deseado hh:mm o hh:mm AM/PM y confirme la selección con .</li> </ul>
<ul> <li>▼ Time ▶</li> <li>2010-04-07 15:05</li> <li>Time / Date</li> <li>▼ Time ▶</li> </ul>	<ul> <li>con .</li> <li>El componente de menú Time destella en el panel de pantalla multifunción, junto al campo de hora en la esquina superior derecha.</li> <li>Presione o o para conmutar al formato de hora deseado hh:mm o hh:mm AM/PM y confirme la selección con .</li> <li>El campo de hora en la esquina superior izquierda cambiará su apariencia según su selección (y dejará de parpadear).</li> </ul>

#### Tabla 7-32 Configurar el formato de la hora

Tabla 7-33 Configurar	la	fecha	у	hora
-----------------------	----	-------	---	------

<u></u>	Presione 🗮 para activar la barra del menú, luego use
	🔊 o < para seleccionar el ícono Ajustes y presione
	para confirmar.
2010-04-06 1:33рм ∢ Error ▶	El ítem de menú Error del submenú Ajustes aparece en la pantalla multifunción.
2010-04-07 3:05рм	Presione ⋗ para conmutar al ítem de menú
	Hora/Fecha y confirme la selección con 🛁.
2010-04-07 3:05 <sub>РМ</sub> Time / Date ↓ Date ▶	El ítem de menú <b>Date</b> parpadea en el panel de pantalla multifuncional.
2010-04-07 3:05рм	En el ítem de menú intermitente Date (Fecha), presione
Time / Date	para conmutar a la opción Adjust (ajustar) y confirme
♦ Set	la selección con 🚛.
	El ítem de menú Set (ajustar) destella en el panel
	visualizador multifunción, junto a la sección de año del
	campo de fecha en la esquina superior izquierda.
	Presione ▶ o < para fijar el año y confirme los
	ajustes con 📜.
	La selección intermitente se desplaza a la sección de
	noras y minutos usando 🥙 o 🕙 y confirme cada
	dato con 🧮.
F	Cuando usted confirma con el botón después de
	esquina superior derecha será actualizado según sus
	datos (y dejará de parpadear).
	La pantalla regresa a su modo predeterminado.
	El ícono Ajustes de la barra de menú se apagará.

### Unidad de visualización de temperatura

El ítem de menú **Settings ->°C / °F** permite conmutar el horno usado para visualizar las temperaturas en Grados Centígrados o Fahrenheit.

**Nota** Esta configuración no afecta el registro de datos vía interfaz RS-232. Todos los datos de temperatura que ingresan a la computadora para propósitos de documentación de parámetros operativos se entregan en  $\mathcal{C}$ .

Tabla 7-34 Conmutar la unidad de visualización de temperatura

F	Presione 🗮 para activar la barra del menú, luego use
	🔊 o < para seleccionar el ícono Ajustes y presione
	para confirmar.
2010-04-06 1:33рм	El ítem de menú Error del submenú Ajustes aparece en la pantalla multifunción.
2010-04-07 10:31ам	Presione ⋗ para conmutar al ítem de menú °C / °F y
<b>€</b> °C / °F <b>→</b>	confirme la selección con 💭.
2010-04-07 10:31AM °C / °F € € €	El panel del monitor multifunción cambia a la siguiente pantalla de selección, con el monitor de temperatura actualmente inactiva destellando (°C o °F) (El ajuste de fábrica es °F).
2010-04-07 10:31am °C / °F ∮°F ▶	Confirme la selección con 🚬.
F	Las temperaturas en el campo de visualización (ítem D1 en figura 7-1 en página 7-1) y panel de ajustes Set (ítem D5 en la misma figura) se visualizarán en la unidad seleccionada recientemente. La pantalla regresa a su modo predeterminado. El ícono Ajustes de la barra de menú se apagará.

### Protección de proceso

La Protección de Procesos interviene cuando un programa ingresa al modo "hold" (mantenimiento de temperatura). Cuando la temperatura medida por el sensor integrado diverge del valor teórico en más de 0,5 °C (32,9 °F), el tiempo restante en modo "hold" se suspende hasta que la diferencia de temperaturas entre valor real y el teórico vuelva ser inferior a 0,5 °C (32,9 °F).

Tabla 7	-35 Encend	iendo y apag	ando la p	rotección d	le proceso
---------	------------	--------------	-----------	-------------	------------

F	Presione 🗮 para activar la barra del menú, luego use
	🔊 o < para seleccionar el ícono Ajustes y presione
	para confirmar.
2010-04-06 1:33рм	El ítem de menú Error del submenú Ajustes aparece en
♦ Error ▶	la pantalla multifuncion.
2010-04-07 10:31ам	Presione ⋗ para conmutar al ítem de menú
Protection	Protection y confirme la selección con 🗾.
2010-04-07 12:43рм	En la pantalla de selección emergente, presione ⋗
Protection	para conmutar entre del estado On a Off.
	Para activaria, conmute del estado Off a On
	respectivamente. Confirme la selección con 💳.
2010-04-07 12:43рм	
2010-04-07 12:43рм Protection	
2010-04-07 12:43թм Protection ∢Off ▶	
2010-04-07 12:43PM Protection	La pantalla regresa a su modo predeterminado.

### Preferencias de programa

El ítem de menú **Settings ->Program** soporta las preferencias descritas a continuación para la ejecución de programas de usuario (ver "Programación" en página 7-40.) usados para controlar las operaciones con el horno. Si un programa termina después de haberse ejecutado a través de uno o más ciclos (ver sección "Prompt de ciclos de programa" a continuación), se da una señal

audible simple consistente en 5 bips si no se presenta una condición de error. Entonces el horno se apaga, salvo que el usuario haya establecido las preferencias para mantenerlo funcionando, como se explica en la sección "Modo "after End" (después de finalizar el programa)" en página 7-37.

#### Prompt de ciclos de programa

Este ítem de menú permite habilitar un prompt (consulta al usuario, función desactivada para transportar el horno de fábrica) que solicita introducir el número de veces que debe repetirse un programa (ver "Iniciando un programa" en página 7-42) cada vez que el usuario ejecuta un programa.

	Presione 💭 para activar la barra del menú, luego use
	🔊 o < para seleccionar el ícono Ajustes y presione
	para confirmar.
2010-04-06 1:33рм	El ítem de menú Error del submenú Ajustes aparece en
€ Error ►	la pantalla multifunción.
2010-04-09 13:43рм	Presione ⋗ para conmutar al ítem de menú Program
∢ Program ▶	y confirme la selección con 🗮.
2010-04-09 13:43рм	En la pantalla de selección emergente, seleccione la
Program ∢ Loops	opción preseleccionada Loops (bucles) presionando 🛁.

2010-04-09 13:43⊧м Program	En la pantalla de selección subsiguiente, habilite el prompt (desactivado en un estado "de transporte")
↓ On ▶	mediante ⋗ para pasar la selección del estado Off al
	$\mathbf{On}$ y confirme la selección con $$ .
	Para desactivar el prompt vuelva a conmutar del
2010-04-09́ 13:43⊦м Program	estado $On$ al $Off$ . Confirme la opción seleccionada
Loops	presionando 🗮.
♦ Off ▶	Si escoge Off, el horno se apaga en el próximo ciclo de
	programa. Usted puede cambiar este comportamiento
	predeterminado modificando la "Modo "after End"
	(después de finalizar el programa)" preferencia
	explicada en la siguiente sección.
	La pantalla regresa a su modo predeterminado.
	El ícono Ajustes de la barra de menú se apagará.

#### Tabla 7-36 Activación / Desactivación del prompt de ciclos de programa

#### Modo "after End" (después de finalizar el programa)

Adicionalmente, este comando permite habilitar un segundo prompt (también desactivado para el transporte del horno desde fábrica) que aparece cuando el usuario ejecuta un programa y solicita introducir el modo de operación que debe seguir el horno tras completarse el programa (ver "Iniciando un programa" en página 7-42):

- Off: Esta opción no solicita la confirmación del usuario antes de iniciar el programa. Esta opción apaga el equipo al finalizar el programa.
- On: Cuando el usuario inicia la ejecución de un programa, esta opción consulta si el horno deberá apagarse, o continuar funcionando en modo de mantenimiento (hold) con los últimos ajustes hasta que el usuario confirme la terminación del programa.

Puede suspenderse un temporizador semanal activo hasta que el mensaje **Program End** sea cancelado. Al reanudarse el conteo del temporizador, el horno buscará establecer el valor teórico que usaba antes de iniciarse el programa (incluyendo velocidad del ventilador y posición del regulador).

<b>F</b>	Presione 🗮 para activar la barra del menú, luego use
	🔊 o < para seleccionar el ícono Ajustes y presione
	para confirmar.
2010-04-06 1:33рм ♦ Error ▶	El ítem de menú Error del submenú Ajustes aparece en la pantalla multifunción.
2010-04-09 13:43թм ∢ Program ▶	Presione Deriver al selección con Confirme la selección con Confirme la selección con
2010-04-09 13:43рм Program (Loops) 2010-04-09 13:43рм Program ( Mode after End )	En la pantalla de selección emergente presione para conmutar de la opción preseleccionada <b>Loops</b> a <b>Mode after End</b> , entonces confirme la selección con $\stackrel{\blacksquare}{\longleftarrow}$ .
2010-04-09 13:43PM Program Mode after End ▲ On ▲ 2010-04-09 13:43PM Program Mode after End ▲ Off ▲	En la pantalla de selección subsiguiente, presione  o o o para conmutar entre los estados On (Mantenimiento de temperatura) y Off (ajuste de fábrica); luego confirme la selección con  .
×	La pantalla regresa a su modo predeterminado. El ícono Ajustes de la barra de menú se apagará.

#### Tabla 7-37 Cambiando el Modo after End
### Configuración

El ítem de menú **Settings -> Configuration** le permite al usuario introducir un código de cuatro dígitos que carga un juego específico de parámetros operativos para el horno – por ejemplo, para realizar la selección de tensión descrita en la sección "Conectar la Interfaz RS 232" en página 5-12.

Tabla 7-38 Introducción del código de cuatro dígitos

<b>F</b>	Presione 💭 para activar la barra del menú, luego use
	🔊 o < para seleccionar el ícono Ajustes y presione
	para confirmar.
2010-04-06 1:33рм ∢Error ▶	El ítem de menú Error del submenú Ajustes aparece en la pantalla multifunción.
2010-04-12 12:00рм	Presione ⋗ para conmutar al ítem de menú
<ul> <li>♦ Configuration &gt;</li> </ul>	Configuration y confirme la selección con 🗮.
2010-04-12 12:00PM Configuration ↓ 0208 ▶	El panel visualizador multifunción presentará el promt similar al ejemplo mostrado a la izquierda, con el primer dígito del código de configuración de cuatro dígitos intermitente. Programe el primer dígito del código de configuración
2010-04-12 12:00թм Configuration ♦ 0240 ▶	con (o) y confirme su ajuste con . Programe los tres dígitos restantes exactamente como se describe arriba.
	Confirmando la entrada del último dígito con , la nueva configuración se activará inmediatamente.
F	La pantalla regresa a su modo predeterminado. El ícono Ajustes de la barra de menú se apagará. Introduzca el código de cuatro dígitos "1234" para resetear la calibración de usuario. Introduzca el código de cuatro dígitos '4321' para resetear la calibración del sensor de muestras.



## Programación

El ítem de menú **Program** permite al usuario crear, almacenar y ejecutar hasta 10 programas para flujos de proceso automáticos. Cada uno de los 10 programas puede consistir en un máximo de 10 pasos.

Las siguientes propiedades pueden definirse para cada paso, dependiendo de la configuración de hardware actual:

- duración en hh:mm
- valor teórico de temperatura
- nivel de velocidad de ventilador (1 a 5)
- posición del regulador (de 0 a 3)

Para facilitar la creación de sus programas, los usuarios pueden escoger copiar y editar programas existentes.

### **End of Program**

El final del programa se indica mediante una señal audible (5 bips). El horno se apaga o pasa al modo Temperature Hold (mantenimiento de temperatura), según se haya preestablecido en las preferencias en "Cambiando el Modo after End" en página 7-38). El display muestra el mensaje **Program End**, así como la temperatura del calor residual.

### Programa de conveniencia ECO (sólo para unidades de mesa)

Con los hornos Heratherm de las series OGH-S y OMH-S, el programa de conveniencia Eco permite apagar automáticamente un horno en cualquier punto de un proceso de secado, tan pronto como estén secas las muestras. Esto requiere conectar un sensor de muestras (accesorio a solicitarse por separado) a un puerto dedicado sobre el panel posterior del horno, insertándolo en la muestra a secar.

Durante el proceso de secado, se compara continuamente la temperatura real de la cámara con la temperatura de muestra medida por el sensor de muestras. Al comienzo del proceso de secado, el efecto refrigerante producido por la evaporación evita que la muestra alcance el límite de temperatura específico preajustado durante la rutina de calibración Eco. Tan pronto como se haya evaporado toda la humedad de la muestra a secar, el sensor de muestras medirá un incremento de temperatura. Si la monitorización de temperatura de muestra indica que se ha sobrepasado el límite, el proceso de secado es terminado y el horno se apaga automáticamente.

El límite de temperatura se calcula a partir del diferencial entre las temperaturas actual de cámara y de muestra, lo cual significa que depende tanto de la ubicación de la muestra dentro de la cámara del equipo como del valor teórico de temperatura para el proceso de secado. Por ello, la implementación del programa de conveniencia Eco requiere de una calibración Eco preliminar del sensor de muestras. Esta rutina de calibración debe repetirse cada vez que se hayan modificado sensiblemente, ya sea la ubicación de la muestra o bien, el valor teórico de temperatura para el proceso de secado.

El resultado de la calibración Eco sólo interviene en los criterios de terminación de programa y no tiene impacto sobre la lectura de temperatura del sensor de muestras en el panel visualizador multifunción.

## Preparación del horno para el programa de conveniencia ECO - Procedimiento de calibración Eco

- 1. Conecte el sensor de muestras al puerto de la parte posterior del panel del horno e insértelo a través del puerto de acceso de conductos (que también se encuentra en el panel posterior del equipo) en la cámara del horno. El panel visualizador multifunción muestra la temperatura medida por el sensor de muestras.
- 2. Instale el sensor de muestras en el lugar previsto de la muestra a secar en la cámara del equipo.

**Nota** La muestra a secar no debe colocarse en la cámara del horno al realizar una calibración Eco.

- 3. Establezca el valor teórico de temperatura para el futuro proceso de secado con el horno.
- 4. Ajuste el regulador en el punto 3 para su apertura completa, como sería el caso durante el proceso de secado. Una vez alcanzado el valor teórico de temperatura previsto, prevea un tiempo de espera de

Una vez alcanzado el valor teórico de temperatura previsto, prevea un tiempo de espera de 2 hs. para que el horno logre estabilizarse en un estado térmico.

5. Seleccione la opción **Eco** del menú **Settings -> Calibration** y confirme con 📰 . Ahora se iniciará la calibración Eco.

#### Carga del horno con la muestra a secar

- 1. Deje enfriar el horno.
- 2. Coloque la muestra a secar en el lugar dentro de la cámara para el cual se haya realizado la calibración Eco más reciente.
- 3. Inserte el sensor de muestras dentro de la muestra.

#### Flujo del programa "ECO Convenience"

1. Programe en el horno el valor teórico de temperatura para el proceso de secado para el cual Ud. haya realizado la calibración Eco más reciente.

- 2. Ejecute el programa Eco según se explica en tabla 7-39, "Iniciando un programa," en página 42. Seleccione la opción Eco en lugar de P1...P10. Ninguno de los promts posibles aparecerá en este caso. El visualizador multifunción muestra el mensaje **Eco** mientras se ejecuta el programa de conveniencia Eco.
- 3. Una vez que el controlador integrado del horno haya llevado la temperatura al valor teórico especificado, comienza la monitorización del límite de temperatura de muestra.
- 4. El proceso de secado finaliza al excederse el límite de temperatura. La finalización del programa se indica mediante una señal audible (5 bips), y aparece el mensaje **Eco End!** en el panel visualizador multifunción, procediendo a apagar, llegado el caso.

### Instrucciones

	1
	Presione ਦ para activar la barra del menú, luego use
	🔊 o < para seleccionar el ícono Programa y
	presione 🗮 para confirmar.
2010-04-13 10:25AM	En la pantalla de selección mostrada a la izquierda,
Program	seleccione la opción preseleccionada Start (inicio)
◆ Start →	presionando 🗮.
NOTA	El prompt Stop Timer al comenzar a ejecutarse un programa Si el temporizador está corriendo, una solicitud adicional Stop Timer aparece en este momento (ver "Detener un tem- porizador de apagado antes de que expire" en página 7-14.). Al seleccionar Yes se aborta la temporización e inicia el pro- grama seleccionado. Al seleccionar No se deja continuar el temporizador. Sin embargo, si un programa de usuario y el temporizador se superponen, el programa que va a iniciarse tiene prioridad sobre el temporizador, es decir, el temporiza- dor será puesto en espera y sólo se iniciará después que el programa hava finalizado.
2040 04 42 40.25	En la lista de programas existentes (por ejemplo: P1P4)
Start	mostrada a la izquierda, conserve la selección por defecto
P1 P2 P3 P4 Eco	P1 o escoja otro programa mediante ⋗ (o < ).
	En cualquiera de los casos, presione ਦ para iniciar
	el programa.

### Tabla 7-39 Iniciando un programa

### Tabla 7-39 Iniciando un programa

2010-04-13 10:25AM Start Loops ∢ 1 ▶	En caso de habérselo activado en las preferencias para programas (ver "Activación / Desactivación del prompt de ciclos de programa" en página 7-36.), aparece el prompt Loops solicitando introducir el número deseado de ciclos de programa (= Loops). Establezca el número de ciclos mediante (>>>>>>>>>>>>>>>>>>>>>>>>>>>>>>>>>>>>
2010-04-13 10:25AM Start Mode after End (Off) 2010-04-13 10:25AM Start Mode after End (Hold Settings)	En caso de habérselo activado en las preferencias para programas (ver "Cambiando el Modo after End" en página 7-38.), aparece el prompt Mode after End. Usted puede usar este diálogo para seleccionar el modo de operación deseado después de finalizada la ejecución de un programa, o sea, si se prevé que el horno pase al modo Standby o al modo Temperature Hold al finalizar el programa. Utilice v o v para escoger el modo de operación deseado y confirme la selección con
13.04.2010 P1[L1]3:02 2 11:05 8888 8888 3 4 5	El panel visualizador multifuncional presenta una barra de progreso para el programa seleccionado, similar al ejemplo mostrado a la izquierda. Las funciones de los elementos de display se explican en tabla 7-41 en página 7-44 más abajo.
	El ícono <b>Program</b> en la barra de menú se ilumina para indicar que un programa de usuario está en marcha.

Las funciones de los elementos de display (a excepción de los campos de fecha y hora; ver tabla 7-2 en página 7-3) se explican en la tabla siguiente.

N⁰	Función
1	<ul> <li>Estos tres campos muestran la siguiente información sobre la ejecución del programa:</li> <li>ID del programa activo – por ejemplo: P1 por Programa 1</li> <li>número del ciclo de programa actual – por ejemplo, L1 para el primer ciclo (aparece sólo si el prompt para el número de ciclos de programa está activado; ver "Prompt de ciclos de programa" en página 7-36)</li> <li>tiempo restante general - por ejemplo: 3:02</li> <li>Nota El tiempo restante no incluye ninguna fase cuya duración se haya establecido en 00:00 para alcanzar un calentameinto o enfriamiento rápido.</li> </ul>
2	Este gráfico de barras utiliza una cifra variable de barras para representar el avance de la ejecución del programa: Tres barras, por ejemplo, indican paso Nº 3, cuatro barras el paso Nº 4 de un programa, etc. Pueden aparecer hasta 10 barras en este lugar, equivalentes a un máximo de 10 pasos (ver más arriba).
3	El ícono de <b>Ventilador</b> se ilumina para indicar que el ventilador fue encendido por el paso de programa actual. El nivel de velocidad del ventilador puede esta- blecerse directamente encima del gráfico de barras ubicado directamente por encima del ícono (ver tabla 7-2 en página 7-3).
4	El ícono de <b>Regulador</b> se ilumina para indicar que el regulador fue abierto por el paso de programa actual. La posición del regulador puede determinarse del gráfico de barras ubicado directamente por encima del ícono (ver tabla 7-2 en página 7-3).
5	El ícono <b>Program</b> en la barra de menú se ilumina para indicar que un programa de usuario está en marcha.

### Tabla 7-40 Elementos de display activos durante la ejecución de un programa

### Tabla 7-41 Cancelando un programa

1 1 1 1 1 1 1 1 1 1 1 1 1 1	Mientras se ejecuta un programa (como se indica en la barra intermitente, número de programa y tiempo restante en el visualizador multifuncional y el ícono de <b>Programa</b> iluminado en la barra de menú) presione $\leftarrow$ 0 $\equiv$ 0
2010-04-16 3:07 <sub>РМ</sub> Stop Program? ∢ No Yes <b>)</b>	Si el prompt mostrado a la izquierda aparece en el visualizador multifunción, presione < para escoger la opción Yes y confirme la selección con 든.
	La pantalla regresa a su modo predeterminado. El ícono <b>Program</b> de la barra de menú se apagará.

### Tabla 7-42 Crear un nuevo programa

	Presione 🗮 para activar la barra del menú, luego use
	🔊 o < para seleccionar el ícono Programa y
	presione 🗮 para confirmar.
2010-04-13 10:25AM Program	En la pantalla de selección emergente presione para conmutar de la opción preseleccionada Start a New (nuevo), y confirme la selección con .
2010-04-13 10:25AM New P4 P5 P6 P7 P8 P9 P10	En la sección siguiente, la pantalla conserva su selección por defecto (por ejemplo $P4$ ) o seleccione cualquier otro programa de la lista de slots libres de memoria de programa (el ejemplo a la izquierda muestra slots de
	En cualquiera de los casos, presione er para comenzar a crear un nuevo programa.
	Aparece una pantalla de programación para el primer paso del programa, que se representa con una barra intermitente vacía. Si el primer paso se va a prolongar más de una hora
	(se admite hasta 23 horas), use <i>para fijar el conteo</i> horario para la duración del primer paso y confirme sus
	Luego, establezca el conteo de minutos mediante (o presione :00 para aceptar cero minutos) y confirme los ajustes con

### Tabla 7-42 Crear un nuevo programa

NOTA	Programar la duración de un paso de programa
NOTA	Si deja el valor por defecto para la duración de un paso <b>00:00</b> , usted le dice al horno que caliente o refrigere al valor teórico de temperatura tan rápido como sea posible. En cambio, si introduce un lapso temporal, el horno tratará de alcanzar el valor teórico según una función de rampa ascendente o descendente. Para cada paso que implique cambios de temperatura, dicha rampa se visualiza como una figura de cuña sobre la barra de progreso.
Set	La selección se desplaza al panel de ajustes Set, el cual
	comenzará a destellar.
	Utilice ⋗ para establecer el valor teórico de temperatura
	deseado y confirme sus ajustes con 픋.
	<b>Nota</b> Si necesita crear un programa que involucre una fase siguiente de enfriamiento, debe tratar de programar el último paso con un valor teórico de temperatura de 0 °C (32 °F) para acelerar el enfriamiento. Vea en "Valor teórico de temperatura" en página 7-7 para mayores detalles e instrucciones.
	La selección se desplaza al gráfico de barra de la
	velocidad del ventilador, el cual comenzará a destellar.
	Si desea que el ventilador funcione a más del 20% de
	su velocidad máxima durante la ejecución del paso de
	programa actual, utilice 🕟 para establecer el nivel
	deseado de velocidad del ventilador y confirme sus
	parámetros con 🗮.
	La selección se desplaza al gráfico de barra del regulador,
	el cual comenzará a destellar.
	Si desea que el regulador permanezca abierto por la
	duración del paso de programa actual, utilice ⋗ para
	establecer la posición deseada para el regulador y
	confirme el dato con 🗮.

### Tabla 7-42 Crear un nuevo programa

New P4 ∢Next ▶	En el visualizador multifunción aparece el prompt Next (siguiente). Si desea crear un paso de programa adicional, presione
New P4 (Save)	el botón . El display retorna a la etapa inicial de la creación del programa. Ahora, la segunda barra destellará representando el segundo paso de programa. Si no desea crear otro paso de programa adicional,
	presione <b>&gt;</b> o <b>&lt;</b> para continuar a la opción Save (guardar), y presione una vez más para guardar el programa recientemente creado.
	La pantalla regresa a su modo predeterminado. El nuevo programa puede ejecutarse ahora, según se describe en "Iniciando un programa" en página 7-42.

### Tabla 7-43 Borrar un programa existente

	Presione 💭 para activar la barra del menú, luego use
	🔊 o < para seleccionar el ícono Programa y presione
	para confirmar.
2010-04-13 10:25AM	En la pantalla de selección mostrada a la izquierda,
Program ∢ Start ▶	presione ⋗ para conmutar a la opción Delete (borrar).
2010-04-13 10:25ам Program ↓ Delete ▶	
2010-04-13 10:25ам	En la lista de programas existentes (por ejemplo:P1P4;
Delete P1 P2 P3 P4	la selección por defecto es P1) mostrada a la izquierda,
	escoja cualquier programa presionando ⋗, y presione
	para confirmar el borrado.

### Tabla 7-43 Borrar un programa existente

	En el diálogo de confirmación de borrado emergente,
Delete P4	presione < para elegir la opción Yes (la selección por
♦ Yes No ▶	defecto es No), luego presione 픋 para llevar a cabo
	el borrado.
	La pantalla regresa a su modo predeterminado.
	La memoria de programa está en condiciones de
	aceptar un programa nuevo o copiado (ver "Crear un
	nuevo programa" en página 7-45 y el procedimiento
	descrito abajo en "Copiar un programa existente").

### Tabla 7-44 Copiar un programa existente

	Presione 🗮 para activar la barra del menú, luego use
	🔊 o < para seleccionar el ícono Programa y
	presione 🗮 para confirmar.
2010-04-13 10:25ам	En la pantalla de selección mostrada a la izquierda,
Program ∢ Start ▶	presione ≥ para conmutar a la opción Copy.
2010-04-13 10:25ам Program Сору	
2010-04-13 10:25ам	En la lista de programas existentes (por ejemplo
Copy P1 P2 P3 P4	P1P4; la selección por defecto es P1) mostrada a la
	izquierda, escoja cualquier programa presionando ⋗
	y presione 🗮 para iniciar el copiado.
	En la lista de slots libres de memoria de programa
to P5	emergente (por ejemplo: P5P10) presione ⋗ para
P6 P7 P8 P9 P10	escoger el slot destino de la copia, luego presione 🗮
	para confirmarlo.

### Tabla 7-44 Copiar un programa existente

	En el diálogo de confirmación de copiado emergente,
Сору Р1> Р5	presione < para elegir la opción Yes (la selección por
	defecto es No), luego presione $\stackrel{\scriptstyle{{\scriptstyle \rm MENU}}}{\leftarrow\!$
	copiado.
	La pantalla regresa a su modo predeterminado.
	El programa que usted acaba de copiar al slot de
	memoria escogido como destino (vea el procedimiento
	descrito más abajo en "Editar un programa existente").

### Tabla 7-45 Editar un programa existente

	Presione 💭 para activar la barra del menú, luego use
	🔊 o < para seleccionar el ícono Programa y
	presione 🗮 para confirmar.
2010-04-13 10:25ам	En la pantalla de selección mostrada a la izquierda,
Program ∢ Start ▶	presione ⋗ para conmutar a la opción Edit.
2010-04-13 10:25ам Program ↓ Edit ▶	
2010-04-13 10:25 AM	En la lista de programas existentes (por ejemplo:
Edit	P1P4; la selección por defecto es P1) mostrada a la
P1 P2 P3 P4	izquierda, escoja el programa que usted necesite
	presionando Ď y presione 🗮 para confirmar la
	selección.

### Tabla 7-45 Editar un programa existente

	En la pantalla de programación emergente, seleccione el
	primer paso de programa para edición presionando 🗮
P4	o utilice ⋗ para navegar a otro paso de programa y
	selecciónelo presionando 🗮.
	Edite la duración, valor teórico de temperatura, velocidad
	del ventilador y posición del regulador como se explica en
	Crear un nuevo programa en pagina 7-45.
	Si ya termino con la ultima edición, puede presionar
	▶ para ir al paso de programa siguiente, o bien 🗮
	para repasar los parámetros del paso de programa que
	acaba de editar. Si éste es el último paso, Ud. recibirá
	el siguiente prompt para salvar cambios.
Sava2	Al solicitársele guardar los cambios, presione < para
No Yes	elegir la opción Yes (la selección por defecto es $No$ ),
	luego presione ਦ para confirmar el guardado.
	Tras confirmar y guardar, el visualizador multifunción
	retornará a la barra intermitente para el paso de
	programa actual.

## Apagado

Este capítulo brinda instrucciones para el apagado del horno durante períodos prolongados, es decir, por lo menos durante varios días consecutivos.

### Apagado del horno

- 1. Extraiga los recipientes con las muestras y todos los accesorios de la cámara.
- 2. Encienda el horno desde el panel de control.
- 3. Desenchufe el cable y asegúrelo contra una reconexión accidental.
- 4. Hasta el cierre del horno, la cámara debe ventilarse continuamente. Deje abierta la puerta y asegúrela contra su cierre accidental.

8

Apagado Apagado del horno

### 9

# Limpieza y Desinfección

## Limpieza



### Limpieza de las superficies externas

Elimine minuciosamente los residuos y depósitos de suciedad usando una solución de agua tibia y detergente común.

Enjuague las superficies usando un paño limpio y agua limpia.

Después seque las superficies usando un paño limpio.

## Desinfección por limpieza y rociado

La desinfección por limpieza manual y rociado es el proceso siguiente:

- predesinfección,
- limpie de la manera correspondiente a la aplicación en curso.

*	iDesinfectantes alcohólicos!				
	Los desinfectantes con contenido alcohólico superior al 10 % pueden formar, en combinación con el aire, mezclas de gas explosivo y de fácil combustión.				
	¡Al usar dichos desinfectantes, evite el fuego abierto o la exposición al calor excesivo durante el proceso de desinfección completo!				
	Utilice tales desinfectantes únicamente en recintos adecuadamente ventilados.				
	Después de haber dejado actuar el desinfectante, seque por completo los componentes limpios del horno.				
	Cumpla con las normas de seguridad para evitar el fuego y/o peligro de explosión causados por los desinfectantes alcohólicos.				
	PRECAUCION ;Desinfectantes con contenido de cloruro!				
	Los desinfectantes con contenido de cloro pueden corroer las superficies de acero inoxidable y metal galvanizado.				
	¡Utilice únicamente desinfectantes que no dañen las superficies de acero inoxidable y metal galvanizado!				

### Preparación de la desinfección manual por limpieza y rociado





Choque eléctrico

El contacto con componentes eléctricos energizados puede causar un choque eléctrico mortal.

Antes de conectar el horno a la fuente de energía, controle el cable y el enchufe en busca de averías. ¡No utilice cables dañados para conectar el horno al suministro eléctrico!



## 

Riesgo para la salud

Las superficies de la cámara pueden estar contaminadas. El contacto con los líquidos de limpieza contaminados puede causar infecciones. Los desinfectantes pueden contener sustancias nocivas.

¡Cuando se limpia y desinfecta, cumpla siempre con las instrucciones de seguridad y las normas de higiene!

- Use guantes de seguridad.
- ¡Utilice gafas protectoras de seguridad!
- Use dispositivos de protección bucal y del sistema respiratorio, para proteger sus membranas mucosas.
- Cumpla con las instrucciones de seguridad de los fabricantes de desinfectantes y del supervisor de higiene.

### Predesinfección

- 1. Extraiga todas las muestras de la cámara y guárdelas en un lugar seguro.
- 2. Rocíe desinfectante sobre las superficies de la cámara y de los accesorios o limpie las superficies usando un desinfectante.
- 3. Permita que el desinfectante actúe como lo especifica el fabricante.



Limpieza y Desinfección Desinfección por limpieza y rociado

## Mantenimiento

El mantenimiento periódico es de misión crítica a la hora de evitar disfunciones por obsolecencia y desgaste. El no cumplimiento del mantenimiento a base regular puede resultar en:

- fluctuaciones en el rendimiento del calentamiento
- pérdida del control sobre la distribución de la temperatura en la cámara
- daños en las muestras

### Inspecciones y controles

Para asegurar el rendimiento operacional y la seguridad del hornos y sus funciones, se deben controlar a intervalos regulares los componentes listados más abajo.

#### **Controles regulares**

- Controlar el horno en cuanto a su limpieza integral, extrayendo cualquier residuo de procesos previos.
- Para evitar operar el horno sin un suministro adecuado de aire puro, controle en busca de contaminación el filtro de aire (accesorio a solicitarse por separado) en el tramo de entrada de aire.

#### Inspección semanal

• Controlar la integridad y la posición correcta de la junta de sellado.

**Nota** Si el horno se hace funcionar por encima de cierto límite de temperatura (ver "Reemplazo de la junta de sellado de la puerta" en página 10-4), el sello de la puerta debe remplazarse cada tres (3) meses (lleve una bitácora de la unidad).

- Cambie el cartucho de filtro de aire (opcional) en la entrada de aire.
- Realice un control funcional del panel de control y del controlador integrado del horno.
- Realice un control de seguridad eléctrica de acuerdo con las normas nacionales pertinentes.

NOTA	Control funcional
Si se extrajo o inhabilitó dis el horno no deberá operars para su funcionamiento cor	positivos de seguridad para las inspecciones, e antes de haberlos reinstalado y controlado recto.

	Rep
	par

Repuestos y modificaciones por parte del usuario

Para evitar fallas de funcionamiento importantes del horno y riesgos de seguridad asociados que puedan resultar en muerte, lesiones graves o daño a el horno y otros equipos, use solamente repuestos aprobados por Thermo Electron LED GmbH. Los repuestos de terceros sin aprobación de anulan la garantía limitada.

No modifique el horno de ninguna manera sin haber obtenido previamente la autorización escrita de Thermo Electron LED GmbH. Las modificaciones no autorizadas pueden comprometer la seguridad operativa y conllevar riesgos que pueden devenir en muerte, daños serios o daño del horno y otros equipos.

### Intervalos de servicio

Durante la operación continua, se deben realizar los siguientes servicios:

### Servicio trimestral

Realizar la medición de temperatura comparativa esquematizada en la siguiente sección.

### Servicio anual

Hacer que el horno sea inspeccionada por y reciba el mantenimiento de un agente del servicio técnico autorizado.



### Preparación para la calibración de la temperatura

Para determinar el valor medido exacto del sensor de temperatura integral del horno, se debe realizar una medición comparativa de temperatura cada tres meses. Si se encuentra una desviación de temperatura importante durante este control, se necesitará una calibración de temperatura. Durante este proceso, el controlador de temperatura del horno se ajusta para el valor medido durante la medición comparativa de temperatura.

Utilice un instrumento de medición calibrado con una precisión de <  $\pm$  1 °C (1,7 °F) para este control.

Para minimizar las variaciones de temperatura durante la medición, coloque la sensor de medición en un contenedor isotérmico (como un bol lleno de glicerol) antes de posicionarlo en el cámara. Utilice el centro de la cámara como lugar de referencia para la medición comparativa.



## Procedimiento de medición por comparación

- 1. Encienda el horno con el interruptor de encendido.
- 2. Establezca el valor teórico de temperatura y deje que el horno se estabilice. Esto puede insumir varias horas.
- 3. Ubique al dispositivo de medición en el área central del soporte de malla en el medio de la cámara. Guíe el cable de conexión a través del conducto de escape de aire en el panel posterior del horno.
- 4. Cierre la puerta.
- 5. Espere hasta que el valor mostrado en el instrumento de medición se haya estabilizado.
- 6. En caso de que las lecturas de temperatura en el dispositivo de medida y el display del horno sean divergentes, utilice la lectura de temperatura del dispositivo de medición para calibrar el control, como se explica en "Introduzca manualmente la temperatura de referencia de calibración" en página 7-28.

### Procedimiento de calibración de la temperatura

Para instrucciones detalladas de cómo hacer una calibración de temperatura manual o automática, remítase a las instrucciones en la sección "Calibración" en página 7-27.

NOTA	Temperaturas de cámara excesivas
Las temperaturas de cámara pueden reducirse dejando la segundos.	a excesivas después de la calibración ls puertas abiertas durante aprox. 30

## Reemplazo de la junta de sellado de la puerta

La junta estanca de la puerta externa está ubicada en la ranura colectora. El sello de la puerta debe inspeccionarse previendo indicios de agrietamiento cada seis meses, si el horno ha sido usado a temperaturas máximas de hasta 250 °C / 482 °F, o cada tres meses si el horno fue utilizado a temperaturas máximas que superaron los 250 °C / 482 °F.

No se necesitan herramientas para su reemplazo.



## Figura 10-1 Reemplazo de la junta de la puerta (el ejemplo muestras un horno de la serie OMH)

- 1. Extraiga la junta de la ranura guía.
- 2. Empiece por el lado de la bisagra, ubicando el encaje de la nueva junta en el lugar indicado por la flecha en figura 10-1, arriba.
- 3. Presione suavemente la junta en la ranura, contemplando la circunferencia de la puerta. Al hacerlo, tenga cuide no estirar la junta de goma.
- 4. Asegúrese de que el elemento de retención del riel esté posicionado correctamente dentro de la ranura y que la junta esté a tope con el marco de la puerta.

## Reemplazo del cable de alimentación eléctrica

Si el cable de alimentación del equipo se daña, debe ser reemplazado por un repuesto original. Se prohíbe el uso de un cable estándar con una resistencia térmica inferior.

## Devoluciones para reparación

Antes de retornar cualquier material, contacte al Departamento de Atención al Cliente para solicitar un número de "Autorización de Devolución de Material" (Return Materials Authorization, RMA).

Será rechazado todo material sin un número RMA.

	ADVERTENCIA Peligro de contaminación			
El horno puede haberse usado para el tratamiento y el procesamie sustancias infecciosas, lo cual puede causar contaminación del horr componentes. Por consiguiente, antes del embarque para la devolución, es oblig de que todos los componentes del horno sean descontaminados adecuadamente				
	<ul> <li>Limpie los componentes del horno por completo, luego desinféctelos o descontamínelos (según la aplicación).</li> </ul>			
	<ul> <li>Complete y adjunte una declaración de seguridad con detalles sobre las actividades de descontaminación realizadas a los componentes que deban ser reparados.</li> </ul>			

Mantenimiento Devoluciones para reparación

## Eliminación

	Peligro de contaminación
El horno puede haberse usad sustancias infecciosas, lo cual componentes. Por ello, antes de su elimina componentes del horno sea	o para el tratamiento y el procesamiento de puede causar contaminación del horno y sus ción, es obligatorio que todos los n adecuadamente descontaminados.
Limpie los componentes desinféctelos o desconta	del horno por completo, luego mínelos (según la aplicación).
Adjunte una declaración o actividades de descontan van a ser eliminados.	le descontaminación con detalles sobre las inación realizadas a los componentes que

## Vista general de los materiales usados

Componente	Material
Componentes de aislación térmica	Lana de vidrio
Placas de circuito impreso	Los componentes eléctricos revestidos contienen diversos materiales plásticos. Los componentes montados en placas de circuito impreso contienen adhesivos de resina epóxica.
Componentes de plástico, en general	Ver etiquetado del material
Carcasa exterior	Lámina de acero galvanizada, pintada
Panel posterior del horno	Lámina de acero galvanizada
Puerta externa	Lámina de acero galvanizada, pintada
Panel interno de puerta	OMH y OGH: acero inoxidable 1.4301
Membrana de protección del panel de control y visualizador	Polietileno
Calefactor	Alambres de resistencia calefactora recu- biertos de acero inoxidable
Recipientes de la cámara	Acero inoxidable 1.4301

11

Componente	Material
Bandejas de malla metálica	Acero, plateado al cromo
Junta estanca, sellado del marco de puerta	Silicona
Rueda de ventilador	Acero inoxidable 1.4016 (sólo series OMH y OMH-S)
Cables	Conductores multifilares de cobre con recubrimiento plástico
Embalaje	Cartón corrugado, película de polietileno y espuma de poliestireno, madera sin tra- tamiento químico

# Códigos de error

Tabla 12-1 muestra a continuación los mensajes de error que podrían aparecer en la ventana del panel de control (ver "Registro de errores (log)" en página 7-26) y da instrucciones para cancelar tales alarmas.

Tabla 12-1 Códigos	s de error de	el horno	Heratherm
--------------------	---------------	----------	-----------

Mensaje y código de error	Causa raíz	Respuesta de alarma	Instrucciones para cancelar alarmas <sup>1</sup>
Door Open Error (E001)	El interruptor de puerta (sólo series OGH-S y OMH-S) ha disparado una alarma por apertura de puerta por más de 10 minutos. Alarma activada.	Alarma audible activada, relé de alarma energi- zado, mensaje aparece en el display.	Cierre la puerta.
Display Error (E002)	Error de comunicación con el display. El controla- dor integrado no pudo res- tablecer comunicación con el panel de control.	Alarma audible activada, relé de alarma energi- zado, mensaje aparece en el display. Resetee des- pués de 30 seg.	Repita el ciclo de encen- dido, desenchufando y reconectando el cordón de alimentación. Si esto no resuelve el problema, llame al service.
Mirrored Parameter Loaded (E003)	El control no pudo leer la configuración de usuario y debió recurrir a un set de parámetros de emergen- cia almacenado en una memoria reflejada.	Alarma audible activada, relé de alarma energizado. Uso alternativo de la memoria de parámetros reflejada. El equipo conti- núa funcionando sin pér- dida de funcionalidad, incluyendo ajustes espe- cíficos del usuario.	Controle los últimos ajus- tes, por ejemplo el valor teórico.
Factory Parameter Loaded (E004)	El control no pudo leer el set de parámetros refle- jado y debió recurrir a parámetros predetermina- dos de fábrica.	Uso alternativo de pará- metros predeterminados de fábrica. Alarma audi- ble activada, relé de alarma energizado, men- saje aparece en el dis- play. Los ajustes específicos de usuario, por ejemplo, las preferen- cias de unidad de visuali- zación de temperatura o programas del usuario, pueden haberse perdido.	Confirmar presionando <sup>ESC</sup> . Vuelva a progra- mar los ajustes específi- cos de cliente.

12

Mensaje y código de error	Causa raíz	Respuesta de alarma	Instrucciones para cancelar alarmas <sup>1</sup>
Default Parameter Loaded (E005)	El control no pudo leer los parámetros preajustados en fábrica y debió recurrir a valores predeterminados.	Uso alternativo de paráme- tros predeterminados, alarma audible activada, relé de alarma energizado, mensaje aparece en el dis- play. El equipo está com- pletamente no-operativo.	Llame al service.
Power Down Error (E007)	Se interrumpió el suminis- tro eléctrico (corte de energía) mientras funcio- naba el equipo.	Alarma audible activada, relé de alarma energi- zado, mensaje aparece en el display.	Controle el suministro eléctrico. Encienda el equipo y luego cancele la alarma presionando
Program Error (008)	Error al procesar un pro- grama del usuario.	El programa es abortado. Alarma audible activada, relé de alarma energi- zado.	Confirmar presionando (ESC).
Fan Error (E009)	La velocidad del ventilador está fuera de rango. ((sólo series OMH y OMH-S)	Alarma audible activada, relé de alarma energi- zado, mensaje aparece en el display. Calefactor inha- bilitado por controlador.	Confirmar presionando <sup>ESC</sup> . Incrementar o decrementar la velocidad del ventilador desde el panel de control. Si esto no resuelve el problema, llame al service.
Damper Error (E010)	Posición del regulador indefinida.	Alarma audible activada, relé de alarma energi- zado, mensaje aparece en el display.	Confirmar presionando <sup>ESC</sup> . Ajuste posición del regulador en el panel de control. Si esto no resuelve el problema, llame al service.
Config Error (E012)	Error general de configu- ración del equipo.	Alarma audible activada, relé de alarma energizado, mensaje aparece en el dis- play. El equipo está com- pletamente no-operativo.	Llame al service.
OTP error (E013)	El contacto Klixon se abrió.	Fallo en la protección por sobretemperatura. Alarma audible activada, relé de alarma energizado, men- saje aparece en el display. El puenteo sobre el con- tacto Klixon ha fallado. (El contacto Klixon se abrió)	Llame al service.
Incorrect voltage (E014)	La tensión aplicada es excesiva o insuficiente.	Alarma audible activada, un mensaje aparece en el display.	Aplique la tensión correcta indicada en la placa de características y tome nota del error.

### Tabla 12-1 Códigos de error del horno Heratherm

#### Mensaje y código de Instrucciones para Causa raíz Respuesta de alarma cancelar alarmas <sup>1</sup> error Sensor Error (E100) Sensor de proceso ave-Alarma audible activada, Llame al service. riado. El valor real medido relé de alarma energiestá fuera de rango. zado, mensaje aparece en el display. Control transferido al sensor de referencia. Si ambos sensores están defectuosos. inhabilite todos los circuitos de control. El valor medido real Protección de proceso Temperature Too High Llame al service. (E101) excede el rango admisiactivada, el control conti-(El valor real supera el ble. El triac es defectuoso. núa sobre el valor teórico. punto teórico (circuito Alarma audible activada, de calefacción defecrelé de alarma energituoso)) zado, mensaje aparece en el display. El valor medido real cae Temperature too low Alarma audible activada. Controle las líneas de CA (E102) (únicamente en por debajo del rango relé de alarma energien cuanto a baja tensión y equipos con interrupadmisible. La tensión de zado, mensaje aparece en solucione el problema, si tor de puerta). sólo alimentación podría ser el display. Control de temes necesario. Si esto no OGH-S y OMH-S. resuelve el problema, insuficiente. peratura continúa. llame al service. Temperature Not Plau-La diferencia entre el con-El equipo usa el sensor Si el error no desaparece, sible (E103) trol y los sensores de refeque indica la mayor temllame al servicio. rencia supera la peratura para el servocontrol. Alarma audible desviación máxima admisible, imposibilitando la activada, relé de alarma medición. energizado, mensaje aparece en el display. El error admite un reset, pero no resetea. Calibration Value Too El valor de referencia de Recurre a la referencia de Controle el correcto fun-High (E104) calibración calculado calibración previa. Alarma cionamiento del sensor de sobre la base de la audible activada, relé de referencia externo y rementrada de usuario está alarma energizado, menplácelo si es necesario. Si cercano al límite superior saje aparece en el display. esto no resuelve el propara las referencias de blema, llame al service. calibración. Calibration value too El valor de referencia de Recurre a la referencia de Controle el correcto funlow (E105) calibración calculado calibración previa. Alarma cionamiento del sensor de sobre la base de la audible activada, relé de referencia externo y rementrada de usuario está alarma energizado, menplácelo si es necesario. Si superio del límite inferior saje aparece en el display. esto no resuelve el propara las referencias de blema, llame al service. calibración. Constant Sensor Sig-Ninguno de los lugares Control transferido al sen-Llame al service. nal (E106) decimales de la salida del sor de referencia. Alarma conversor A/D para el audible activada, relé de sensor de proceso ha alarma energizado, mencambiado a lo largo de un saje aparece en el disperíodo especificado. play. Si ambos sensores están averiados, se desactivarán todos los circuitos de control.

#### Tabla 12-1 Códigos de error del horno Heratherm

Mensaje y código de error	Causa raíz	Respuesta de alarma	Instrucciones para cancelar alarmas <sup>1</sup>
Constant Reference Sensor Signal (E107)	Ninguno de los lugares decimales de la salida del conversor A/D para el sensor de proceso ha cambiado a lo largo de un período especificado.	Control continúa con el sensor de proceso, alarma audible activada, relé de alarma energi- zado, display muestra un mensaje. Si ambos senso- res están averiados, se desactivarán todos los cir- cuitos de control.	Llame al service.
Constant Sample Sen- sor Signal (E108)	Ninguno de los lugares decimales de la salida del conversor A/D para el sensor de proceso ha cambiado a lo largo de un período especificado.	Alarma audible activada, relé de alarma energi- zado, mensaje aparece en el display.	Llame al service.
Heating Relay Error (E109)	La medición de tensión ha indicado una avería en el relé del circuito de cale- facción.	Equipo completamente no-operativo, alarma audi- ble activada, relé de alarma energizado, dis- play muestra un mensaje.	Asegúrese de retirar el enchufe eléctrico de las líneas de CA. Llame al service.
Heating Triac Error (E110)	La medición de tensión ha indicado una avería en el triac.	Alarma audible activada, relé de alarma energizado, mensaje aparece en el dis- play. Alarma acústica retorna tras el reseteo.	Llame al service.
Temperature Too High (E111)	El valor medido real excede el rango admisible.	Alarma audible activada, relé de alarma energi- zado, mensaje aparece en el display. Calefactor apa- gado hasta recuperar el límite superior de histére- sis. Las operaciones de servocontrol continúan. La alarma puede resetearse, apagándose cuando deje de haber diferencia entre el valor real y el teórico. Nota: Este error no indica un triac defectuoso.	Abra la puerta para acele- rar el enfriamiento. Con- trole si el equipo fue cargado con un objeto caliente; en tal caso, quí- telo. Asegúrese de que el equipo haya funcionado más de 10 minutos con al menos una bandeja perfo- rada y con la puerta sin abrir. Si esto no resuelve el problema, llame al service.
Sensor Error (E112)	El valor real medido está fuera de rango.	Alarma audible activada, relé de alarma energi- zado, mensaje aparece en el display. Control conti- núa con el sensor de pro- ceso. Si ambos sensores están defectuosos, inhabi- lite todos los circuitos de control.	Llame al service.
Sensor Error (E113)	El valor real medido está fuera de rango.	Alarma audible activada, relé de alarma energi- zado, mensaje aparece en el display.	Llame al service.

Tabla 12-1 Códigos de error del horno Heratherm

### Tabla 12-1 Códigos de error del horno Heratherm

Mensaje y código de error	Causa raíz	Respuesta de alarma	Instrucciones para cancelar alarmas <sup>1</sup>
ADC Error (E114)	El conversor A/D no entrega datos plausibles.	Alarma audible activada, relé de alarma energi- zado, mensaje aparece en el display.	Llame al service.
Watchdog error (E115)	Prueba del watchdog (vigía) fallida al encen- dido.	Alarma audible activada, relé de alarma energi- zado, mensaje aparece en el display.	Llame al service.

1.La rectificación de un error se considera exitosa cuando la alarma sonora deja de sonar, el relé de alarma suelta y desaparece el mensaje de la pantalla del panel de control.

Códigos de error

## **Datos técnicos**

Los datos técnicos son válidos sólo para un equipo vacío equipado con tres bandejas, un cuerpo exterior pintado a soplete y para una tensión de línea de 230 V/50 Hz. Otras opciones pueden afectar el rendimiento especificado.

Parámetro	Unidad	OGH 60	OGH 100	OGH 180	OGH-S 60	OGH-S 100	OGH-S 180
Proceso							
Atmósfera de la cámara Mín. <sup>1</sup>	°C/°F	50/122	50/122	50/122	50/122	50/122	50/122
Máx.	°C/°F	330/626	330/626	330/626	330/626	330/626	330/626
Desviación de temperatura del valor teórico a 150 ℃ (302 °F), espacial. Valor máx./Valor típico	К	±3 / ±2,5	±3,5 / ±3	±3 / ±2,5	±3 / ±2,5	±3,5 / ±3	±3 / ±2,5
Desviación de temperatura del valor teórico a 150 ℃ (302 °F), sostenido. Valor máx./Valor típico	К	±0,3 / ±0,3	±0,4 / ±0,3	±0,3 / ±0,3	±0,3 / ±0,3	±0,4 / ±0,3	±0,3 / ±0,3
Tiempo de calenta- miento (cámara des- ocupada, de 25 °C (77 °F) a 98% de tem- peratura de ajuste de 150 °C (302 °F)	min	22/20	25/22	25/22	22/20	25/22	25/22
Tiempo de recuperación (cámara desocupada, puerta abierta por 30 s, a temperatura de ajuste). Valor máx./Valor típico	min	9/8	10/8	9/8	9/8	10/8	9/8
Tiempo de calentamiento con refuerzo (boost) de calor activo (para temperatura de ajuste de 300 ℃ / 572 °F)	min	35	26	41	35	26	41
Disipación calórica al ambiente (a temperatura de ajuste de 150 °C (302 °F) y temperatura ambiente de 25 °C (77 °F)	W	170 ± 10%	210 ± 10%	290 ± 10%	170 ± 10%	210 ± 10%	290 ± 10%

#### Tabla 13-1 Datos técnicos - Series OGH y OGH-S

13

				0011400							
Parametro	Unidad	OGH 60	OGH 100	OGH 180	OGH-S 60	OGH-S 100	OGH-S 180				
Cambio de aire máx. a 150 °C/302 °F	h⁻¹	26	14	17	26	14	17				
Dimensiones globales											
Altura	mm/in	720/	820/	920/	720/	820/	920/				
		28,3	32,3	36,2	28,3	32,3	36,2				
Ancho	mm/in	530/	640/	640/	530/	640/	640/				
		20,8	25,2	25,2	20,8	25,2	25,2				
Profundidad	mm/in	565/ 25,2	565/ 25,2	738/ 29,1	565/ 25,2	565/ 25,2	738/ 29,1				
Peso global	kg/lbs	44/97	55/121	69/152	44/97	55/121	69/152				
Capacidad de carga											
Capacidad de carga por bandeja	kg/lbs		25/55			25/55					
Máxima capacidad de carga general por equipo	kg/lbs	50/110	50/110	75/165	50/110	50/110	75/165				
Datos eléctricos											
Potencia nominal	W	1810	3100	3100	1810	3100	3100				
Corriente máxima	А	7,9	13,5	13,5	7,9	13,5	13,5				
Sistema de puesta a tierra (ej. 1/N/PE)		1/N/PE	1/N/PE	1/N/PE	1/N/PE	1/N/PE	1/N/PE				
Frecuencia de línea eléctrica	Hz		50/60			50/60					
Tensión de línea +/- 10 %	V	230	2	30	230	230 230					
Sistema de protección IP			IP 20			IP 20					
Clase de protección			I			I					
Categoría de sobretensión según IEC 60364-4-443					II						
Protección del equipo, lado edilicio	А		16		16						
Protección fusible del equipo, en PCB	A		2 x 16		2 x 16						
Condiciones ambientale	es										
Temperatura ambiente mínima	°C/°F		18/65		18/65						
Temperatura ambiente máxima	°C/°F		32/90		32/90						
Humedad máxima en servicio, sin condensación	% h.r./ % r.H.	80, s	sin condens	ación	80, sin condensación						
Temperatura mínima de almacenamiento	℃/℉		20/68			20/68					
Temperatura máxima de almacenamiento	°C/°F		60/140		60/140						

#### Tabla 13-1 Datos técnicos - Series OGH y OGH-S

Parámetro	Unidad	OGH 60	OGH 100	OGH 180	OGH-S 60	OGH-S 100	OGH-S 180
Humedad máxima en depósito, sin condensación	% h.r./ % r.H.	90, s	in condens	ación	90, sin condensación		
Tiempo de aclimatación pos transporte	h	2			2		
Nivel de ruido	dB(A)	34			34		
Grado de contaminación según IEC EN 61010-1		2			2		
Condiciones de emplaz	amiento						
Altitud máxima sobre el nivel del mar	m/y ASL	2000/2187			2000/2187		
Espacio lateral mínimo	mm/in		50/2		50/2		
Espacio frontal mínimo	mm/in	590 / 23,2	690 / 27,2	814 / 32	590 / 23,2	690 / 27,2	814 / 32
Espacio mínimo a la pared posterior	mm/in	80/3,2			80/3,2		
Espacio inferior mínimo	mm/in	200/8			200/8		
Espacio superior mínimo	mm/in	300/12			300/12		

#### Tabla 13-1 Datos técnicos - Series OGH y OGH-S

<sup>1</sup> Al seleccionar temperaturas de una temperatura ambiente de +10 °C, la válvula de alimentación de aire debe estar abierta y no debe entrar calor adicional al aparato.

### Tabla 13-2 Datos técnicos - Series OMH-S y OMH

Parámetro	Unidad	OMH 60	OMH 100	OMH180	OMH 60-S	OMH 100-S	OMH 180-S
Proceso							
Atmósfera de la cámara Mín. <sup>1</sup>	℃/℉	50/122	50/122	50/122	50/122	50/122	50/122
Máx.	°C/°F	330/626	330/626	330/626	330/626	330/626	330/626
Desviación de temperatura del valor teórico a 150 °C (302 °F), espacial. Valor máx./Valor típico.	К	±2,2 / ±1,8	±1,5 / ±1,3	±2,2 / ±1,8	±2,2 / ±1,8	±1,5 / ±1,3	±2,2 / ±1,8
Desviación de temperatura del valor teórico a 150 °C (302 °F), sostenido. Valor máx./Valor típico.	К	±0,25/±0,2	±0,25/±0,2	±0,25/±0,2	±0,25/±0,2	±0,25/±0,2	±0,25/±0,2
Tiempo de calentamiento (cámara desocupada, de 25 ℃ (77 °F) a 98% de temperatura de ajuste de 150 °C (302 °F) Valor máx./Valor típico	min	20/18	18/16	20/18	20/18	18/16	20/18

Parámetro	Unidad	OMH 60	OMH 100	OMH180	OMH 60-S	OMH 100-S	OMH 180-S	
Tiempo de recuperación (cámara desocupada, puerta abierta por 30 s, a temperatura de ajuste). Valor máx./Valor típico.	min	5/4	5/4	5/4	5/4	5/4	5/4	
Tiempo de calentamiento con refuerzo (boost) de calor activo (para temperatura de ajuste de 300 °C / 572 °F)	min	41	25	36	41	25	36	
Disipación calórica al ambiente (a temperatura de ajuste de 150 $^{\circ}$ C (302 $^{\circ}$ ) y temperatura ambiente de 25 $^{\circ}$ C (77 $^{\circ}$ F)	W	259 ± 10%	358 ± 10%	424 ± 10%	259 ± 10%	358 ± 10%	424 ± 10%	
Cambio de aire máx. a 150 °C/302 °F	h⁻¹	33	56	32	33	56	32	
Dimensiones globales								
Altura	mm/in	720/ 28,3	820/ 32,3	920/ 36,2	720/ 28,3	820/ 32,3	920/ 36,2	
Ancho	mm/in	530/ 20,8	640/ 25,2	640/ 25,2	530/ 20,8	640/ 25,2	640/ 25,2	
Profundidad	mm/in	565/ 25,2	565/ 25,2	738/ 29,1	565/ 25,2	565/ 25,2	738/ 29,1	
Peso global	kg/lbs	44/97	55/121	69/152	44/97	55/121	69/152	
Capacidad de carga								
Capacidad de carga por bandeja	kg/lbs		25/55			25/55		
Máxima capacidad de carga general por equipo	kg/lbs	50/110	50/110	75/165	50/110	50/110	75/165	
Datos eléctricos								
Potencia nominal	W	1400	3060	3060	1400	3060	3060	
Corriente máxima	Α	6,1	13,3	13,3	6,1	13,3	13,3	
Sistema de puesta a tierra (ej. 1/N/PE)		1/N/PE	1/N/PE	1/N/PE	1/N/PE	1/N/PE	1/N/PE	
Frecuencia de línea eléctrica	Hz		50/60			50/60		
Tensión de línea +/- 10 %	V	230	23	30	230	23	30	
Sistema de protección IP			IP 20			IP 20		
Clase de protección		Ι			l			
Categoría de sobretensión según IEC 60364-4-443						II		
Protección del equipo, lado edilicio	А		16			16		

### Tabla 13-2 Datos técnicos - Series OMH-S y OMH
Parámetro	Unidad	OMH 60	OMH 100	OMH180	OMH 60-S	OMH 100-S	OMH 180-S
Protección fusible del equipo, en PCB	А	2 x 16			2 x 16		
Condiciones ambientales							
Temperatura ambiente mínima	°C/°F	18/65			18/65		
Temperatura ambiente máxima	°C/°F	32/90			32/90		
Humedad máxima en servicio, sin condensación	% h.r./ % r.H.	80, sin condensación			80, sin condensación		
Temperatura mínima de almacenamiento	°C/°F		20/68			20/68	
Temperatura máxima de almacenamiento	°C/°F	60/140				60/140	
Humedad máxima en depósito, sin condensación	% h.r./ % r.H.	90, :	90, sin condensación		90, sin condensación		
Tiempo de aclimatación pos transporte	h		2			2	
Nivel de ruido	dB(A)	45	52	52	45	52	52
Grado de contaminación según IEC EN 61010-1		2			2		
Condiciones de emplaz	amiento						

#### Tabla 13-2 Datos técnicos - Series OMH-S y OMH

Altitud máxima sobre el nivel del mar	m/y ASL	2000/2187			2000/2187		
Espacio lateral mínimo	mm/in	50/2			50/2		
Espacio frontal mínimo	mm/in	590	690	814	590	690	814
Espacio mínimo a la pared posterior	mm/in	80/3,2			80/3,2		
Espacio inferior mínimo	mm/in	200/8			200/8		
Espacio superior mínimo	mm/in	300/12			300/12		

<sup>1</sup> Al seleccionar temperaturas de una temperatura ambiente de +10 °C, la válvula de alimentación de aire debe estar abierta y no debe entrar calor adicional al aparato.

#### Tabla 13-3 Datos técnicos - Serie OMH

Parámetro	Unidad	OMH 400	OMH 750	OMH 750-3P
Proceso				
Atmósfera de la cámara Mín. <sup>1</sup>	℃/℉	50 ℃/122 ℉	50 ℃/122 ℉	50 ℃/122 ℉
Máx.	°C/°F	250 °C/482 °F	250 °C/482 °F	300 ℃/572 °F
Desviación de temperatura del valor teórico a 150 °C (302 °F), espacial. Valor máx./Valor típico.	К	±2,5/±2,1	±3,5/±3,1	±3,5/±3,1

#### Tabla 13-3 Datos técnicos - Serie OMH

Parámetro	Unidad	OMH 400	OMH 750	OMH 750-3P
Desviación de temperatura del valor teórico a 150 ℃ (302 °F), sostenido.	К	±0,4/±0,3	±0,5/±0,4	0,5/±0,4
Tiempo de calentamiento (cámara desocupada, de 25 ℃ (77 °F) a 98% de temperatura de ajuste de 150 ℃ (302 °F) Valor máx./Valor típico	min	35	50	50
Tiempo de recuperación (cámara desocu- pada, puerta abierta por 30 s, a temperatura de ajuste). Valor máx./Valor típico.	min	<7 / <5	<12 / <10	<12 / <10
Disipación calórica al ambiente (a temperatura de ajuste de 150 °C (302 °F) y temperatura ambiente de 25 °C (77 °F)	W	630 ±10%	990 ±10%	990 ±10%
Cambio de aire máx. a 150 °C (302 °F)	h⁻¹	23	17	17
Dimensiones globales				
Altura (con ruedas)	mm/in		1655/ 65,2	
Ancho	mm/in	755/ 29,7	/ 1215/ / 47,8	
Profundidad	mm/in	770/ 30,3		
Peso global	kg/lbs	135/298	185/408	185/408
Capacidad de carga por bandeja	kg/lbs	40/88		
Máxima capacidad de carga general por equipo	kg/lbs	75/165 150/330		/330
Datos eléctricos				
Potencia nominal	W	3000	3000	5750
Corriente máxima	А	10,5	13,0	12,0
Sistema de puesta a tierra (ej. 1/N/PE)		1/N/PE	1/N/PE	3/N/PE
Frecuencia de línea eléctrica	Hz		50/60	
Tensión de línea +/- 10 %	V	23	0	400
Sistema de protección IP			IP 20	
Clase de protección			I	
Categoría de sobretensión según IEC 60364-4-443		II		
Protección del equipo, lado edilicio	А	16 3 x 16		3 x 16
Protección fusible del equipo, en PCB	А	2 x 16 3 x 16		3 x 16
Condiciones ambientales				
Temperatura ambiente mínima	°C/°F	°F 18/65		
Temperatura ambiente máxima	°C/°F	32/90		
Humedad máxima en servicio, sin condensación	% h.r./ % r.H.	80	), sin condensació	'n

#### Tabla 13-3 Datos técnicos - Serie OMH

Parámetro	Unidad	OMH 400	OMH 750	OMH 750-3P	
Temperatura mínima de almacenamiento	°C/°F		20/68		
Temperatura máxima de almacenamiento	℃/℉	60/140			
Humedad máxima en depósito, sin condensación	% h.r./ % r.H.	90, sin condensación		on	
Tiempo de aclimatación pos transporte	h	ו 2			
Nivel de ruido	dB(A)	no ruido próprio			
Grado de contaminación según IEC EN 61010-1		2			
Condiciones de emplazamiento					
Altitud máxima sobre el nivel del mar	m/y ASL	2000/2187			
Espacio lateral mínimo	mm/in	n/in 120/4,7			
Espacio frontal mínimo	mm/in	810 /31,9 670 / 26,4		26,4	
Espacio mínimo a la pared posterior	mm/in		120/5,9		
Espacio superior mínimo	mm/in	n 200/8			

<sup>1</sup> Al seleccionar temperaturas de una temperatura ambiente de +10 °C, la válvula de alimentación de aire debe estar abierta y no debe entrar calor adicional al aparato.

Datos técnicos

## Piezas de repuesto y accesorios

Material Nº	Descripción
50027662	Filtro de aire puro para estufas y hornos de secado Heratherm
50073715	Fusibles 5,0 A 250V T 6, 3 X 32 mm
50126665	Adaptador de apilado Heratherm 60L
50126666	Adaptador de apilado Heratherm 100L
50126667	Adaptador de apilado Heratherm 180L
50127102	Juego de filtro de aire puro para hornos Heratherm
50127431	Puerta exterior para equipos Heratherm IGS 60, IMH 60, IMH 60-S, OMS 60, OMH 60, OMH 60-S, OGS 60, OGH 60 y OGH 60-S con un tope del lado izquierdo
50127432	Puerta exterior para equipos Heratherm IGS 100, IMH 100, IMH 100-S, OMS 100, OMH 100, OMH 100-S, OGS 100, OGH 100 y OGH 100-S con tope del lado izquierdo
50127433	Puerta exterior para equipos Heratherm IGS 180, IMH 180, IMH 180-S, OMS 180, OMH 180, OMH 180-S, OGS 180, OGH 180 y OGH 180-S con tope del lado izquierdo
50127434	Puerta exterior para equipos Heratherm IGS 60, IMH 60, IMH 60-S, OMS 60, OMH 60, OMH 60-S, OGS 60, OGH 60 y OGH 60-S con tope del lado derecho
50127435	Pies de apilado par incubadores y estufas y hornos de secado Heratherm
50127436	Juego de junta de puerta para Heratherm 60L
50127437	Juego de junta de puerta para Heratherm 100L
50127438	Juego de junta de puerta para Heratherm 180L
50127439	Juego de bloqueo de puerta para Heratherm 60L / 100L / 180 L
50127443	Pies niveladores para incubadores y estufas y hornos de secado Heratherm
50127444	Juego de manija derecha de puerta para incubadoras, así como estufas y hornos de secado Heratherm
50127445	Juego de manija izquierda de puerta para incubadoras, así como estufas y hornos de secado Heratherm
50127446	Juego de manija derecha de puerta para incubadoras, así como estufas y hornos de secado Heratherm

Material Nº	Descripción
50127447	Juego de manija izquierda de puerta para incubadoras, así como estufas y hornos de secado Heratherm
50127448	Sensor de temperatura para estufas y hornos de secado Heratherm
50127450	Bisagra superior de puerta para incubadoras, así como estufas y hornos de secado Heratherm
50127451	Bisagra inferior de puerta para incubadoras, así como estufas y hornos de secado Heratherm
50127455	Puerta exterior para equipos Heratherm IGS 100, IMH 100, IMH 100-S, OMS 100, OMH 100, OMH 100-S, OGS 100, OGH 100 y OGH 100-S con tope del lado derecho
50127456	Puerta exterior para equipos Heratherm IGS 180, IMH 180, IMH 180-S, OMS 180, OMH 180, OMH 180-S, OGS 180, OGH 180 y OGH 180-S con tope del lado derecho
50127458	Juego de panel operativo para incubadoras, así como estufas y hornos de secado Heratherm de Protocolo Avanzado y de Seguridad de Protocolo de Avanzado
50127462	Juego de inserto electrónico sin ventilador de placa madre para incuba- doras, así como estufas y hornos de secado Heratherm de Protocolo Avanzado y de Seguridad de Protocolo Avanzado
50127463	Cable de placa madre para incubadoras, así como estufas y hornos de secado Heratherm
50127469	Interruptor de puerta para el montaje en la parte derecha de los incubadores et las estufas y los hornos de secado Heratherm
50127470	Interruptor de puerta para el montaje en la parte izquierda de los incubadores et las estufas y los hornos de secado Heratherm
50127477	Limitador de temperatura para OMH, OMH-S, OGH, OGH-S
50127480	Gancho de cierre magnético de puerta para incubadoras, así como estufas y hornos de secado Heratherm con tope de puerta a la derecha
50127481	Gancho de cierre de puerta para incubadoras, así como estufas y hornos de secado Heratherm con tope de puerta a la izquierda
50127482	Gancho de cierre magnético de puerta para incubadoras, así como estufas y hornos de secado Heratherm con tope de puerta a la derecha
50127483	Gancho magnético de cierre de puerta para incubadoras, así como estufas y hornos de secado Heratherm con tope de puerta a la izquierda
50127499	Serpentines calefactores para Heratherm OGS 60, OGH 60, OGH 60-S para una tensión de 120 V
50127500	Serpentines calefactores para Heratherm OGS 60, OGH 60, OGH 60-S para una tensión de 230 V
50127502	Serpentines calefactores para Heratherm OGS 100, OGH 100, OGH 100-S para una tensión de entre 208 V y 240 V

Material Nº	Descripción
50127503	Serpentines calefactores para Heratherm OGS 180, OGH 180, OGH 180-S para una tensión de entre 208 V y 240 V
50127504	Serpentines calefactores para Heratherm OMS 60, OMH 60, OMH 60-S para una tensión de 120 V
50127509	Juego de sistema de ventilador para Heratherm OMS 60, 120 V
50127510	Juego de sistema de ventilador para Heratherm OMS 100 y OMS 180, para 208-240 V
50127511	Juego de sistema de ventilador para Heratherm IMH 60, IMH 60-S, OMH 60 y OMH 60-S, IMH 100, IMH 100-S, IMH 180 y OMH 180-S, de 120 V
50127512	Juego de sistema de ventilador para Heratherm OMH 100, OMH 100, IMH 180, IMH 100-S y OMH 180, para 208 a 240 V
50127513	Juego de sistema de ventilador para Heratherm OMS 60, 230 V
50127514	Juego de sistema de ventilador para Heratherm OMS 100 y OMS 180, para 230 V
50127515	Serpentines calefactores para Heratherm OMS 60, OMH 60, OMH 60-S para una tensión de 230 V
50127519	Serpentines calefactores para Heratherm OMS 100, OMH 100, OMH 100-S, OMS 180, OMH 180-S para un voltaje entre 208 y 240 V
50127532	Ventilador de circulación OMH / OMH-S D = 180 mm (7,1 inch), H = 28 mm (1,1 inch)
50127544	Junta de puerta ventana para estufas y hornos de secado Heratherm
50127555	Juego de sistema de ventilador para Heratherm IMH 60, IMH 60-S, OMH 60 y OMH 60-S, IMH 100, IMH 100-S, IMH 180 y OMH 180-S, de 230 V
50127556	Juego de sistema de ventilador para Heratherm OMH 100, OMH 100-S, IMH 180, IMH 100-S y OMH 180, para 230 V
50127557	Perilla de entrada de aire mecánica para Heratherm OMS 60 y OGS 60
50127558	Perilla de entrada de aire mecánica para Heratherm OMS 100, OGS 100, OMS 180 y OGS 180
50127559	Juego de deflector eléctrico de aire para hornos Heratherm de Protocolo Avanzado y de Seguridad de Protocolo Avanzado
50127566	Filtro de aire puro para OMH, OGH, OMH-S, OGH-S
50127662	Filtro de aire puro OGH / OGH-S / OMH / OMH-S
50127741	Base de soporte con ruedas para Heratherm 60L
50127742	Base de soporte con ruedas para Heratherm 100L
50127743	Base de soporte con ruedas para Heratherm 180L
50127761	Bandeja de malla de alambre OGS 60 / OGH 60 / OGH-S 60, incl. 2 soportes de las bandejas

Material Nº	Descripción
50127762	Bandeja de malla de alambre OGS 100 / OGH 100 / OGH-S 100, incl. 2 soportes de las bandejas
50127763	Bandeja de malla de alambre OGS 180 / OGH 180 / OGH-S 180, incl. 2 soportes de las bandejas
50127764	Bandeja de malla metálica OMS 60/100/180 / OMH 60/100/180 / OMH 60/100/180-S, incl. 2 soportes de bandejas
50127765	Bandeja de malla metálica OMS 100 / OMH 100 / OMH 100-S, incluyendo 2 abrazaderas de soporte de las bandejas
50127766	Bandeja de malla metálica OMS 180 / OMH 180 / OMH 180-S, incluyendo 2 abrazaderas de soporte de las bandejas
50127767	Sensor de muestras OGH 60-S, OGH 100-S, OGH 180-S, OMH 60-S, OMH 100-S, OMH 180-S
50127773	Bandeja perforada en acero inoxidable IMH 60 / IMH 60-S / OMH 60 / OMH 60-S/OMS 60/100/180, incl. 2 soportes de bandejas.
50127774	Bandeja perforada en acero inoxidable IMH 100 / IMH 100-S / OMH 100 / OMH 100-S/OMS 60/100/180, incl. 2 soportes de bandejas.
50127777	Bandeja perforada en acero inoxidable IMH 180 / IMH 180-S / OMH 180 / OMH 180-S/OMS 60/100/180, incl. 2 soportes de bandejas.
50127861	Resorte de retención para los incubadores et las estufas y los hornos de secado Heratherm
50127862	Riel de soporte para Heratherm IGS 60, IMH 60, IMH 60-S, OMS 60, OMH 60, OMH 60-S
50127863	Riel de soporte para Heratherm IGS 100, IMH 100, IMH 100-S, OMS 100, OMH 100, OMH 100-S
50127864	Riel de soporte para Heratherm IGS 180, IMH 180, IMH 180-S, OMS 180, OMH 180, OMH 180-S
50127914	Bandeja de acero inoxidable perforado OGS 60 / OGH 60 / OGH 60-S, incl. 2 soportes de las bandejas
50127925	Bandeja de acero inoxidable perforado OGS 100 / OGH 100 / OGH 100-S, incl. 2 soportes de las bandejas
50127926	Bandeja de acero inoxidable perforado OGS 180 / OGH 180 / OGH 180-S, incl. 2 soportes de las bandejas
50128182	Ventilador de refrigeración de panel operativo principal para estufas y hornos de secado Heratherm de Protocolo Avanzado y de Seguridad de Protocolo Avanzado
50128184	Conexión del sensor de muestras para incubadores y estufas y hornos de secado Heratherm
50128186	Juego de bombillas de ventana de estufas y hornos de secado Heratherm
50128237	Juego de llaves para manija de puerta con cerrojo, Heratherm

Material Nº	Descripción
50128880	Bandeja de malla metálica reforzada OGS 180 /OGH 180/ OGH 180-S
50128881	Bandeja de malla metálica reforzada OMS 180 /OMH 180/ OMH 180-S
50128887	Juego tubo flexible de aire caliente para las unidades Heratherm empotradas
50130347	Almohadillas de fibra de vidrio Ø 20 mm (0,8")
50130348	Almohadillas de fibra de vidrio Ø 50 mm (2")
50130657	Juego de junta de puerta de Viton, 60 L Heratherm
50130658	Juego de junta de puerta de Viton, 100 L Heratherm
50130659	Juego de junta de puerta de Viton, 180 L Heratherm
50134094	Juego de bobina calefactora OGS 750, 230 V
50134115	Juego de bobina calefactora OGS 400, 750, 230 V
50134118	Juego de bobina calefactora OMH 750, 230 V
50134119	Juego de bobina calefactora OMH 750, 3 fases+N
50134124	Juego de bobina calefactora OGS 750, 3 fases
50134125	Juego de bobina calefactora OGS 750, 3 fases
50134315	Juego cable de bus DS compl. 400 / 750
50134326	Junta de puerta 400 L HTM
50134327	Junta de puerta 750 L HTM
50134328	Juego de junta de perfil 750 L HTM
50134329	Juego de clips conec. puerta 400 / 750 HTM
50134333	Juego castors 400 / 750 HTM
50134334	Juego rack de bandejas compl. HTM 400 / 750
50135043	Juego inserto eléctrico compl. HTM H pie de piso
50135044	Juego inserto eléctrico compl. HTM S pie de piso
50135055	Juego inserto eléctrico compl. HTM 3 PH para piso
50135058	Juego cerrojo puerta 750 izq. compl. HTM
50135059	Juego cerrojo puerta 750 der. compl. HTM
50135060	Carcasa externa puerta izq. HTM 400
50135061	Carcasa externa puerta der. HTM 400
50135062	Carcasa externa puerta izq. HTM 750
50135063	Carcasa externa puerta der. HTM 750
50135151	Juego sensor de temperatura compl. hornos HTM p/piso
50135152	Juego corte sobretemperatura hornos HTM 350 p/piso

Material Nº	Descripción
50135153	Juego cerrojo puerta 400 der. compl. HTM
50135154	Juego cerrojo puerta 400 izq. compl. HTM

15

# Dispositivo de registro

Tipo de horno:			Número de pieza:		
Número de serie:			Número de reparaciones:		
Lugar			Nota del operador:		
Tareas realizadas		Aclaracion	nes	Fecha	Firma

Dispositivo de registro

### Contacto

Vista general de la Organización de Venta Internacional Thermo Fisher

#### Dirección postal en Alemania:

Thermo Electron LED GmbH Robert-Bosch-Straße 1 D - 63505 Langenselbold

#### Consultas desde Alemania:

 Teléfono

 Ventas
 0800 1 536376

 Service
 0800 1 112110

 Fax
 Ventas/Service 0800 1 112114

 E-Mail
 info.labequipment.de@thermofisher.com

#### Consultas desde Europa, Medio Oriente y África:

Teléfono.	+ 49 (0) 6184 / 90-6940
Fax	+ 49 (0) 6184 / 90-6772
Email	info.labequipment.de@thermofisher.com

#### Dirección postal EEUU:

Thermo Scientific 275 Aiken Road Asheville, NC 28804 EEUU

#### Consultas desde América del Norte:

Teléfono	+1 800-879 7767
Fax	+1 828-658 0363
Email	info.labequipment@thermofisher.com

#### Consultas desde América Latina:

Teléfono	+1 828-658 2711
Fax	+1 828-645 9466
Email	info.labequipment@thermofisher.com

#### Consultas desde Asia del Pacífico:

Teléfono	+852-2711 3910
Fax	+852-2711 3858
Email	info.labequipment@thermofisher.com

16

Thermo Fisher Scientific, Inc. 81 Wyman Street P.O. Box 9046 Waltham, MA 02454-9046 United States

www.thermo.com

