Thermo Scientific Orion Star serie A210

1 1 Mar 22 M

Medidores de sobremesa de electroquímica 68X001262 • de junio de 2015





Nota importante

Lea esta guía de usuario detenidamente antes de utilizar el medidor. Cualquier uso distinto al de estas instrucciones puede anular su garantía y provocar un daño permanente al medidor.

Información de contacto

Para obtener asistencia técnica para productos Thermo Scientific™ Orion™, póngase en contacto con el soporte técnico por correo electrónico en <u>wai.techservbev@thermofisher.com</u> o por teléfono. Dentro de los Estados Unidos, llame

al 1-800-225-1480 y fuera de los Estados Unidos, llame al +1 -978-232-6000 o envíe un fax al +1 -978-232-6031.

Para obtener información adicional sobre el producto, póngase en contacto con su distribuidor autorizado local, su representante de ventas técnicas local de Thermo Scientific Orion o póngase en contacto con nosotros mediante la información de Water and Laboratory Products (WLP) de la página final de este manual.

Aplicaciones y recursos técnicos

Visite <u>www.thermoscientific.com/water</u> para visualizar productos de Thermo Scientific Orion y descargar documentación sobre el producto, guías de usuario y manuales, actualizaciones de software y las últimas aplicaciones y recursos técnicos.

Índice

CAPÍTULO 1 Introducción al medidor	5
Descripción general del medidor	5
Lista de embalaje	6
Indicaciones de uso	6
CAPÍTULO 2 Principios básicos del medidor	7
Uso del adaptador de alimentación universal	7
Instalación de las pilas (fuente de alimentación opcional)	8
Conexión del soporte y portaelectrodos	g
Uso del portaelectrodos	
Conexiones del medidor	11
Teclado del medidor	
Teclas de función	14
Pantalla del medidor	15
Iconos de la pantalla de medida	17
Icono de estado del electrodo de pH	
Modelos de medidor y capacidades de medida	
Mantenimiento del medidor	
CAPÍTULO 3 Menús de configuración del medidor	20
Menú de configuración principal	
Navegación general en el menú de configuración	
Menús de método, modo y temperatura por canal	25
Menú del método	
Menú de modo y ajustes	
Menú Temperature	
Menú de configuración Instrument Settings	
Menú Log View	
Registro de datos	
Registro de calibración	
Menú Diagnostics	
Procedimiento de prueba automática del medidor	
Procedimiento de prueba de estabilidad del electrodo	
CAPITULO 4 El uso del canal pH o pH/ISE	52
Preparación del electrodo y el medidor	
Procedimiento de calibración de pH	
Edición de calibración de pH	
Procedimiento de calibración ORP (Modo mV relativo)	
Procedimiento de calibración ORP (E _H Units)	
Procedimiento de calibración ISE	
Edición de calibración ISE	
Procedimiento de medida	61

CAPÍTULO 5 El uso del canal de conductividad	62
Preparación del medidor y el sensor	62
Procedimiento de calibración de conductividad	63
Edición de calibración de conductividad	65
Procedimiento de calibración de introducción de constante de celda	
certificada alternativa	65
Estándares de conductividad frente a Tabla de temperatura	66
Procedimiento de verificación de medidor de conductividad	67
Procedimiento de medida	68
CAPÍTULO 6 El uso del canal OD/RDO	69
Preparación del medidor y el sensor	69
Procedimiento de calibración de oxígeno disuelto	70
Calibración de aire (aire saturado con agua)	71
Calibración de agua (agua saturada con aire)	71
Calibración manual (Winkler)	72
Calibración de definición cero	73
Procedimiento de medida	74
CAPÍTULO 7 Transferencia de datos y actualizaciones	
de software	75
Almacenamiento de datos y ajustes de transferencia	75
Ajuste del tipo de lectura de la medida	75
Ajustes del registro de datos	77
Ajustes de exportación de datos	77
Compatibilidad y requisitos de la impresora	80
Compatibilidad y requisitos del ordenador	81
Uso del cable de ordenador USB a serie	
Uso del cable de ordenador USB	
Software de comunicación Orion Star Com	87
Conexión del medidor con otros programas informáticos	
Protocolos de control remoto del medidor Star A200-A300	
Procedimiento de actualización de software del medidor	97
CAPÍTULO 8 Servicios de atención al cliente	102
Sugerencias de solución de problemas	103
Procedimiento de restablecimiento de fábrica del medidor	104
Procedimiento de restablecimiento del usuario del medidor	104
Declaración de conformidad	105
Conformidad RAEE	105
Declaración de conformidad	106
Especificaciones del medidor	107
Accesorios, electrodos y soluciones del medidor	118

CAPÍTULO 1 Introducción al medidor

Descripción general del medidor

Los medidores de sobremesa Thermo Scientific[™] Orion Star[™] serie A210 cuentan con una pantalla gráfica del medidor, informativa y fácil de leer, con retroiluminación, calibración en pantalla e instrucciones de configuración para un manejo intuitivo y fácil para el usuario con una formación mínima. El teclado completo con teclas de función específicas de menú y útiles teclas de acceso directo que permite la navegación y el control del medidor de forma rápida y eficiente. La interfaz de usuario multilingüe permite la personalización del medidor en una serie de idiomas locales y pueden añadirse nuevos idiomas a través de las actualizaciones del software.

Las medidas pueden realizarse rápidamente y de forma fiable con el indicador de estabilidad de lectura en pantalla y los modos de lectura seleccionables: Auto-Read (Lectura automática), Timed (Sincronizada) o Continuous (Continua) con función de duración indefinida. Data Log (Registro de datos) recopila hasta 2000 conjuntos de medidas con marca de fecha y hora, y la memoria del medidor no volátil conserva todos los datos, incluso cuando hay un fallo del suministro eléctrico. Utilice el software de Thermo Scientific[™] Orion[™] Star Com[™] para agilizar la transferencia de datos del medidor a un ordenador a través de USB o RS232; exporte los datos a una hoja de cálculo Excel o un archivo de valores separados por comas (.csv) e imprima los datos en una impresora local o en red.

Utilice la sonda de agitador Orion Star para el agitado de muestras controlado y activado por el medidor con cinco velocidades de agitado seleccionables. Utilice la sonda de agitador con el portaelectrodos incluido para simplificar la colocación dentro o fuera de soluciones y eliminar la necesidad de barras de mezcla magnéticas y placas.

La carcasa del medidor resistente a salpicaduras y polvo con calificación IP54 ofrece la flexibilidad de colocar el medidor donde sea necesario: en la mesa de trabajo o montado en una pared para ahorrar espacio. Para obtener una mayor adaptación al sitio, encienda el medidor utilizando el adaptador de alimentación universal o cuatro pilas AA opcionales. Hay seis medidores de sobremesa Orion Star serie A210 disponibles para adaptarse a sus necesidades de medidas exactas.

Medidor de sobremesa de pH Orion Star A211

Mida pH, mV, mV relativo u ORP con temperatura

Conductímetro Orion Star A212

Mida la conductividad, STD, salinidad o resistividad con temperatura

Medidor de RDO/OD Orion Star A213

Mida oxígeno disuelto como % de saturación o concentración con temperatura utilizando el sensor de OD polarográfico u óptico RDO[®].

Medidor de pH/ISE Orion Star A214

Mida la concentración de iones utilizando un electrodo de ion selectivo (ISE), pH, mV, mV relativo u ORP con temperatura

Medidor de pH/conductividad Orion Star A215

Mida pH, mV, mV relativo u ORP con temperatura en el canal uno; y mida la conductividad, STD, salinidad o resistividad con temperatura en el canal dos.

Medidor de pH/RDO/OD Orion Star A216

Mida pH, mV, mV relativo u ORP con temperatura en el canal uno y

mida el oxígeno disuelto como % de saturación o concentración con temperatura en el canal dos.

Lista de embalaje

Los medidores de sobremesa Orion Star serie A210 y los kits de medidor incluyen los siguientes elementos:

- Soporte y portaelectrodos conectables al medidor
- Adaptador de alimentación universal
- CD de documentación
- Guía de inicio rápido impresa
- Cable de interfaz de ordenador
- Certificado de prueba del medidor

Todos los medidores Orion Star serie A210 han sido probados completamente y certificados por Thermo Fisher Scientific. También se incluye un certificado de calibración con cada medidor. Para ver contenido específico del kit y el medidor, consulte el apartado Información para pedidos. Visite nuestro sitio web en

<u>http://www.thermoscientific.com/OrionMeters</u> para descargar el software complementario de ordenador para la transferencia de datos Orion Star Com y el controlador de USB Star serie A200/A300.

Indicaciones de uso

Lea esta guía de referencia detenidamente. Cualquier uso distinto al de estas instrucciones puede anular su garantía y provocar un daño permanente al medidor.



CAPÍTULO 2 Principios básicos del medidor

Uso del adaptador de alimentación universal

En los medidores de sobremesa Orion Star serie A210 se incluye un adaptador de alimentación universal (número de catálogo 1010003) con clavijas de enchufes US, EU, UK y China. El adaptador de alimentación universal está específicamente diseñado para su uso con los medidores de sobremesa Star serie A210. El uso de otros adaptadores de alimentación puede dañar el medidor y anular la garantía.

- 1. Seleccione la clavija de enchufe apropiada para la toma de corriente que se utilizará.
- 2. Retire la cubierta transparente de plástico de la ranura de la parte trasera del adaptador de alimentación.
- 3. Inserte la clavija de enchufe apropiada en la ranura de la parte trasera del adaptador de alimentación.



 Conecte el adaptador de alimentación montado a una toma de corriente y a la entrada de suministro eléctrico del medidor. Además, se recomienda un protector contra sobretensión o sistema de alimentación ininterrumpida (SAI).

Instalación de las pilas (fuente de alimentación opcional)

- 1. Seleccione cuatro pilas alcalinas AA.
- 2. Confirme que el medidor está apagado.
- 3. Dele la vuelta, con la pantalla orientada hacia abajo, en una superficie seca y limpia.
- 4. Retire la tapa del compartimento de las pilas: empuje hacia abajo la lengüeta del compartimento de las pilas y levante la tapa del compartimento de las pilas.



- 5. Coloque las nuevas pilas en el compartimento para pilas. La orientación de las mismas debe seguir las indicaciones de la carcasa del compartimento.
- 6. Sustituya la tapa del compartimento de las pilas y dé la vuelta al medidor, con la pantalla orientada hacia arriba.

Conexión del soporte y portaelectrodos

El portaelectrodos puede conectarse a cualquier lateral del medidor y pueden conectarse hasta dos soportes a cada medidor. También hay disponible una base ponderada (número de catálogo STARA-HB) para mantener el soporte sin conectar al medidor.

 Abra la caja que contiene el soporte del electrodo. Esta caja incluirá una placa base, un brazo del electrodo, perno pivotante, portaelectrodos, separador ATC y clip RDO.



- 2. Dele la vuelta, con la pantalla orientada hacia abajo, en una superficie seca y limpia.
- 3. Identifique el lateral del medidor en que se instalará el soporte y retire el tornillo entre los círculos de ese lado del medidor.
- 4. Alinee la placa base del soporte con los círculos del medidor.
- 5. Sustituya el tornillo del paso 3 para conectar la placa base al medidor.
- 6. Dé la vuelta al medidor, con la pantalla orientada hacia arriba.
- 7. Inserte el brazo del electrodo en el poste de metal de la placa base.
- 8. Conecte el portaelectrodos al brazo del electrodo mediante el perno pivotante.



Uso del portaelectrodos

Coloque los electrodos en el soporte para un fácil movimiento dentro y fuera de los contenedores durante la calibración, la medida de la muestra y el almacenamiento. Las posiciones recomendadas de los electrodos se muestran a continuación.



10 Manual de usuario del medidor portátil Orion Star serie A320

Conexiones del medidor

El siguiente diagrama describe todas las conexiones posibles del medidor en los medidores de sobremesa Orion Star serie A210. Algunos medidores de sobremesa Star serie A210 tendrán menos conexiones, dependiendo de las capacidades de medida del medidor.



Conector	Función	Modelos de medidor aplicables
Power (alimenta- ción)	Conecte el adaptador de alimentación universal (incluido con el medidor) para encender el medidor	Todos los medidores Star serie A210
Stirrer (Agitador)	Conecte una sonda de agitado Orion Star (número de catálogo 096019) para el agitado de muestras controlado y activado por el medidor con cinco velocidades seleccionables.	Todos los medidores Star serie A210
USB	Conecte el cable USB para la comunicación y la transferencia de datos bidireccional a través de USB a una impresora u ordenador	Todos los medidores Star serie A210
RS232	Conecte el cable RS232 para la comunicación y la transferencia de datos bidireccional a través de RS232 a una impresora u ordenador	Todos los medidores Star serie A210
BNC	Conecte un electrodo para pH, un electrodo redox/ORP o un electrodo de ion selectivo (ISE) con el conector BNC	Medidores Star A211, Star A214, Star A215 y Star A216
Ref.	Conecte un electrodo de referencia de media celda con un conector con punta de pines de 2,5 mm estándar	Medidores Star A211, Star A214, Star A215 y Star A216
	Conecte una sonda de temperatura ATC con el conector MiniDIN de 8 pines	Medidores Star A211, Star A214 y Star A216
ATC/CON	Conecte un sensor de conductividad o una sonda de temperatura ATC con el conector MiniDIN de 8 pines	Medidores Star A212 y Star A215
Toma de tierra	Utilícese para reducir las interferencias generadas por otros equipamientos	Todos los medidores Star serie A210
OD/RDO	Conecte un sensor de oxígeno disuelto (polarográfico u óptico RDO) con el conector MiniDIN de 9 pines.	Medidores Star A213 y Star A216

Los medidores de sobremesa Orion Star serie A210 son compatibles con los mismos electrodos y sensores utilizados con los medidores Thermo Scientific[™] Orion [™] Versa Star[™] y los anteriores medidores Thermo Scientific[™] Orion Star[™] y Star Plus. Esto incluye los electrodos de pH/ATC Thermo Scientific[™] Orion [™] ROSS Ultra[™] Triode[™], los electrodos de pH Thermo Scientific[™] Orion [™] ROSS[™] y los sensores de conductividad Thermo Scientific[™] Orion[™] DuraProbe[™]. Conexiones del medidor de pH Orion Star A211



Conexiones del conductímetro Orion Star A212



Conexiones del medidor de RDO/OD Orion Star A213



Conexiones del medidor de pH/ISE Orion Star A214



Conexiones del medidor de pH/conductividad Orion Star A215



Conexiones del medidor de pH/RDO/OD Orion Star A216



Teclado del medidor

El teclado del medidor de sobremesa Orion Star serie A210 incluye teclas de función específicas de menú que actualizan la pantalla para un funcionamiento eficiente y rápido del medidor y teclas de acceso directo para disfrutar de una navegación sencilla para realizar la medida, y menús de registros de calibración, registros de datos y configuración.



Nombre e icono de la tecla	Función
f1, f2, f3	Pulse las teclas de función f1, f2 y f3 para realizar la acción mostrada sobre cada tecla en la pantalla.
power	Pulse la tecla de power (encendido) para encender el medidor. Cuando el medidor esté encendido, pulse y suelte la tecla power para activar o desactivar la retroiluminación de la pantalla. Mantenga pulsada la tecla de power durante unos tres segundos para apagar el medidor.
measure measure (esc)	Pulse la tecla measure (medir) (esc) en el modo de medición Auto-Read para iniciar una nueva medida. Pulse la tecla measure (esc) para salir del modo o menú actual y volver al modo de medida.
log view log view	Pulse la tecla log view (ver registro) para acceder al registro de datos y al registro de calibración en el modo de medida.
stirrer	Pulse la tecla stirrer (agitador) para activar y desactivar la sonda de agitador en el modo de medida temporizado o continuo y el modo de edición de calibración.

Nombre e icono de la tecla	Función
setup / flecha arriba (▲)	Pulse la tecla setup (configuración) para acceder al menú de configuración desde el modo de medida. Pulse la tecla flecha arriba (▲) para desplazar hacia arriba una lista de elementos.
mode / flecha derecha (►)	Pulse la tecla mode (modo) para cambiar el modo de medida del canal mostrado. Pulse la tecla flecha derecha (►) para desplazar hacia la derecha una lista de elementos.
log/print / flecha abajo (▼)	Pulse la tecla log/print (registrar/imprimir) para registrar y/o imprimir manualmente una medida, dependiendo del modo de medida y los ajustes de salida de datos seleccionados. Pulse la tecla flecha abajo (▼) para desplazar hacia abajo una lista de elementos.
Hold / flecha izquierda (◄)	En el modo de medida continuo, pulse la tecla hold (bloqueo) para bloquear (congelar) la medida actualmente mostrada y púlsela de nuevo para desbloquear (descongelar) la medida. Pulse la tecla flecha izquierda (◄) para desplazar hacia la izquierda una lista de elementos.

Teclas de función

La siguiente imagen muestra las teclas **f1**, **f2** y **f3** con sus acciones correspondientes encima de cada una de ellas en la pantalla del medidor. Al pulsar la tecla **f1 (cal)** (calibración), solicitará al medidor acceder al modo de calibración. Al pulsar la tecla **f2 (sample ID)** (ID de la muestra), solicitará al medidor acceder al modo de configuración de ID de la muestra. Al pulsar la tecla **f3 (setup)**, solicitará al medidor acceder al menú de configuración principal.



Pantalla del medidor

Ejemplos de pantalla de medida

Las siguientes pantallas solo son ejemplos. Las pantallas reales del medidor variarán dependiendo de los parámetros de configuración seleccionados, los datos de calibración activos, etc.

Pantalla de medida de pH Orion Star A211



Pantalla del medidor de RDO/OD Orion Star A213





Pantalla del medidor de pH/ISE Orion Star A214

Pantalla del medidor de pH/conductividad Orion Star A215

Visualice cada canal por separado o ambos canales simultáneamente mediante la tecla **f3 (channel)** (canal).



Pantalla del medidor de pH/RDO/OD Orion Star A216

Visualice cada canal por separado o ambos canales simultáneamente mediante la tecla **f3 (channel)**.



Iconos de la pantalla de medida

lcono de la pantalla	Descripción
+	Se muestra cuando el medidor funciona con alimentación de CA
	Se muestra cuando el medidor funciona con pilas.
뼵	Indica cuándo se exportan los datos a un ordenador o impresora
60	Indica cuándo se graba una medida en el registro de datos
-	Se muestra cuando se establece una alarma y esta se activa
R\$282	Indica el puerto RS232 seleccionado para la conexión del ordenador o impresora
¥	Indica el puerto USB seleccionado para la interfaz del ordenador o impresora
25.0 °C	Muestra las medidas de temperatura activas e indica la fuente como un sensor de temperatura ATC o valor de temperatura introducido manualmente (MAN)
HOLD	Se muestra cuando se pulsa la tecla hold y las medidas mostradas se congelan; pulse la tecla hold una segunda vez para desbloquear la función hold.
N	Indica que una calibración se ha completado con éxito, parpadea cuando la alarma de la calibración está establecida y la alarma se activa
Ę,	Indica el estado del electrodo de pH como bueno (dos barras), medio (una barra) o malo (tachado), según la última calibración guardada y la estabilidad del electrodo
Σŝ	Indica el número de método de medida activo; M100 y M200 son los métodos predeterminados e indican que no se está utilizando un método protegido con contraseña
ready	El indicador de estabilidad parpadeará stabilizing (estabilizando) mientras se cambia la medida y mostrará ready (listo) cuando la medida sea estable
AR	Cuando el tipo de lectura se establece en Auto-Read, el icono parpadeará mientras se estabiliza la lectura y permanecerá estático cuando la lectura se mantenga estable y fija en la pantalla.
п	Indica que la función ID de la muestra está activa; el número asignado al operador se muestra a la derecha del icono
[†]	Indica que la función ID de usuario está activa; el nombre asignado al operador se muestra a la derecha del icono

Icono de estado del electrodo de pH

Los medidores de pH Orion Star A211, los medidores de pH/ISE Orion Star A214, los medidores de pH/conductividad Orion Star A215 y los medidores de pH/RDO/OD Orion Star A216 incluyen un icono de estado del electrodo de pH. En el modo de medida, el icono de estado del electrodo de pH indica el desempeño de este, basado en la última calibración guardada y en la estabilidad de medida del electrodo.

Icono	Estado del electrodo de pH
3	El estado del electrodo es bueno. La pendiente del electrodo está en el intervalo de 95,1 % a 104,9 %.
£	El estado del electrodo es medio. La pendiente del electrodo está en el intervalo de 85,1 % a 95 % o de 105 % a 114,9 %.
£	El estado del electrodo es malo. La pendiente del electrodo es igual o inferior a 85 % o igual o superior a 115 %. Consulte el manual del electrodo de pH para obtener instrucciones sobre cómo limpiar, mantener y reparar el electrodo.

Nota: Esto es un indicador general del estado general del electrodo. Consulte siempre el manual de usuario del electrodo de pH para obtener información específica sobre el intervalo de pendiente recomendado para cada electrodo de pH.

Modelos de medidor y capacidades de medida

La siguiente tabla muestra los modelos disponibles de los medidores de sobremesa Orion Star serie A210 y sus modos de medida disponibles. Todas las medidas incluyen temperatura.

Modelo de medidor	Canal 1 Modos de medida	Canal 2 Modos de medida	Modos de temperatura
Medidor de pH Star A211	pH mV RmV ORP	n/d	Automatic (Automático) Manual
Conductímetro Star A212	Conductivity (Conductividad) TDS (STD) Salinity (Salinidad) Resistivity (Resistividad)	n/d	Automatic Manual
Medidor de RDO/OD Star A213	DO as % Saturation (OD como % de saturación) DO as mg/L (OD como mg/L)	n/d	Automatic
Medidor de pH/ISE Star A214	pH mV RmV ORP ISE	n/d	Automatic Manual
Star A215 pH/Conductividad Medidor	pH mV RmV ORP	Conductivity TDS Salinity Resistivity	Automatic Manual
Medidor de pH/RDO/OD Star A216	pH mV RmV ORP	DO as % Saturation DO as mg/L	Automatic Manual (solo canal 1)

Mantenimiento del medidor

- Para el mantenimiento rutinario del medidor, límpielo con un paño húmedo. Si fuese necesario, puede utilizarse detergente con agua caliente o tibia.
- El mantenimiento del medidor puede realizarse de forma diaria, semanal o mensual, según lo requiera el entorno de funcionamiento.
- Retire del medidor o módulo cualquier sustancia vertida inmediatamente utilizando el procedimiento de limpieza adecuado para el tipo de vertido.



CAPÍTULO 3 Menús de configuración del medidor

Menú de configuración principal

El menú de configuración principal de los medidores Orion Star serie A210 contiene en una única ubicación los menús para los ajustes de medida, ajustes del instrumento, registros de datos y calibración y diagnósticos del medidor.

Navegación general en el menú de configuración

- 1. En el modo de medida, pulse la tecla **setup** para acceder al menú de configuración principal.
- Pulse la tecla ▲, ▼, ◀ o ▶ para desplazarse o resaltar un icono del menú de configuración y pulse la tecla f3 (select) (seleccionar) para acceder a los submenús del menú seleccionado.
- 3. Pulse la tecla ▲ o ▼ para desplazarse o resaltar una opción de submenú y pulse la tecla **f3 (select)** para acceder a los parámetros del submenú seleccionado.
- 4. Realice los pasos adecuados para establecer los parámetros y ajustes en el menú seleccionado.
 - a. Para seleccionar un valor de la lista de opciones, pulse la tecla ▲ o ▼ para resaltar el valor deseado y después pulse la tecla f3 (select) para establecer el valor.
 - b. Para introducir un valor numérico, utilice la ventana emergente de introducción de números.
 - i. Pulse la tecla **f3 (edit)** (editar) para abrir la ventana de introducción de números.



- Pulse la tecla ▲, ♥, ◀ o ▶ para resaltar un número, punto decimal o signo negativo y después pulse la tecla f3 (enter) (introducir) para seleccionar el elemento resaltado. Repita hasta que el valor deseado se muestre en la parte superior de la ventana.
- iii. Pulse la tecla **f2 (done)** (realizado) para guardar el valor y salir de la ventana de introducción de números.
- 5. Pulse la tecla **f1 (back)** (atrás) para navegar fuera de un menú y pulse la tecla **measure (esc)** para volver al modo de medida en cualquier momento.

				• /	
ICONOS V	v descrinciones	del menu	de conti	duración	principal
1001100		aci mena		guruoion	principui

Icono	Descripción	Modelos de medidor aplicables
Chaund	Utilice el menú pH Channel (Canal pH) para personalizar los ajustes de medida, calibración y alarma para pH, mV, RmV (mV relativo), ORP y la temperatura	Medidor de pH Star A211 Medidor de pH/conductividad Star A215 Medidor de pH/RDO/OD Star A216
Channel	Utilice el menú pH/ISE Channel (Canal pH/ISE) para personalizar los ajustes de medida, calibración y alarma para pH, mV, RmV (mV relativo), ORP, ISE y la temperatura	Medidor de pH/ISE Star A214
Channel	Utilice el menú COND Channel (Canal COND) para personalizar los ajustes de medida, calibración y alarma para la conductividad, la salinidad, los STD, la resistividad y la temperatura	Conductímetro Star A212 Medidor de pH/conductividad Star A215
Channel	Utilice el menú DO/RDO Channel (Canal OD/RDO) para personalizar los ajustes de medida, calibración y alarma para el oxígeno disuelto y la temperatura	Medidor Star de RDO/OD Star A213 Medidor de pH/RDO/OD A216
Settings	Utilice el menú Instrument Settings (Ajustes del instrumento) para actualizar los ajustes del medidor para la transferencia de datos, registro de datos, fecha y hora, idioma, sonidos, velocidad rotatoria, contraste de la pantalla, apagado automático, ID de usuario e ID de la muestra	Todos los medidores Star serie A210
View Log	Acceda al menú View Log (Ver registro) para ver, exportar o borrar datos guardados en el registro de datos y ver o imprimir las diez calibraciones más recientes por canal en el registro de calibración.	Todos los medidores Star serie A210
Diagnosties	Acceda al menú Diagnostics (Diagnóstico) para iniciar un restablecimiento del medidor, realizar una prueba automática del medidor, comprobar la estabilidad del electrodo o visualizar el número de serie del medidor y la revisión de software.	Todos los medidores Star serie A210

Ejemplos de pantallas del menú de configuración principal



Menú de configuración del medidor de pH/conductividad Star A215



Menú de configuración del medidor de pH/ISE Star A214



Menú de configuración del medidor de pH/RDO/OD Orion Star A216



Diagrama de flujo del menú de configuración principal

* Consulte la siguiente figura para obtener una lista detallada de los ajustes de medida, calibración y alarma en el menú de modo y ajustes por canal.



Diagrama de flujo del menú de modo y ajustes por canal

Menús de método, modo y temperatura por canal

En los menús pH Channel, pH/ISE Channel, COND Channel y DO/RDO Channel, se encuentran los submenús de Method (Método), Mode and Settings (Modo y ajustes) y Temperature (Temperatura), que pueden utilizarse para personalizar los ajustes de medida, calibración y alarma para cada canal. Los menús de los canales mostrados en cada medidor dependerán del modelo del medidor y sus capacidades de medida.

Modelo de medidor	Menús de canales disponibles	
Medidor de pH Orion Star A211	pH Channel	
Conductimetro Orion Star A212	COND Channel	
Medidor de RDO/OD Orion Star A213	RDO/DO Channel	
Medidor de pH/ISE Orion Star A214	pH/ISE Channel	
Medidor de pH/conductividad	pH Channel	
Orion Star A215	COND Channel	
Madidaraa da al URDO/OD Orian Star A216	pH Channel	
	RDO/DO Channel	

- Method: utilice el menú de configuración Método para crear, cargar, copiar, editar o eliminar métodos protegidos con contraseña.
- **Mode and Settings**: seleccione este menú para revisar y actualizar los ajustes de medida, calibración y alarma para el canal seleccionado.
- Temperature: utilice este menú para introducir manualmente un valor de temperatura de muestra, establecer las unidades de temperatura en °C o °F, realizar una calibración de temperatura en una sonda ATC o un sensor de oxígeno disuelto/conductividad con temperatura integrada y establecer la fuente de entrada de temperatura para medidores con fuentes de temperatura dobles.

Menú del método

Ahorre hasta diez métodos por canal en el menú de método para obtener una recopilación rápida y sencilla de los ajustes de medida, calibración y alarma personalizados por canal.

El método predeterminado (M100 o M200) se actualiza cuando los ajustes de medida, calibración y alarma por canal se cambian en el menú de modo y ajustes y el método predeterminado no está protegido por contraseña.

Los métodos personalizados (M101-M110 o M201-M210) pueden protegerse mediante una contraseña de tres a ocho dígitos. Al crear, cargar y activar un método protegido en el modo de medida y realizar una calibración, esa calibración se guardará en el método para que cada vez que este se cargue, también se cargue la calibración correspondiente.

Los métodos protegidos resultan útiles cuando se utilizan dos o más electrodos en un canal. Por ejemplo, un electrodo de pH y un electrodo de ion selectivo (ISE) utilizados en la misma entrada BNC del medidor de pH/ISE Star A214 o un sensor de conductividad de rango bajo y un sensor de conductividad de rango estándar utilizados en la misma entrada MiniDIN de 8 pines del conductímetro Star A212.

Creación de un nuevo método mediante los ajustes actuales del medidor

- 1. En el modo de medida, pulse la tecla setup.
- Pulse la tecla ▲, ▼, ◀ o ► para resaltar pH Channel, pH/ISE Channel, COND Channel o DO/RDO Channel y pulse la tecla f3 (select).
- Pulse la tecla ▲ o ▼ para resaltar Method y pulse la tecla f3 (select) para acceder a la lista de métodos.
- Pulse la tecla ▲ o ▼ para resaltar Current Settings (Ajustes actuales) y pulse la tecla f3 (save) (guardar).
- Pulse la tecla ▲ o ▼ para resaltar un método abierto (M101-M110 o M201-M210) y guardarlo en los ajustes actuales específicos del canal y, a continuación, pulse la tecla f2 (accept) (aceptar).
 - a. Los métodos abiertos no tienen fecha, hora o modo visibles en la lista de métodos.
- Utilice la ventana emergente de introducción de números para crear una contraseña específica para el método (de 3 a 8 dígitos).
 - Pulse la tecla f3 (edit) para acceder a la ventana de introducción de números.
 - b. Pulse la tecla ▲, ▼, ◀ o ► para resaltar un número, pulse la tecla f3 (enter) para seleccionar el número y repítalo hasta que la contraseña deseada se muestre en la parte superior de la ventana de introducción de números.



- c. Pulse la tecla **f2 (done)** para aceptar la contraseña y salir de la ventana de introducción de números.
- 7. Pulse la tecla f2 (accept) para guardar la contraseña introducida.
- 8. El medidor volverá a la lista de métodos y el número de método seleccionado tendrá la fecha, la hora y el modo visibles en la lista de métodos.

Carga de un método

Utilice la función Load para activar un método protegido para su uso en el modo de medida. Si no se ha creado ningún método protegido, el método activo será el que esté abierto. No se necesita contraseña para cargar un método protegido.

- 1. En el modo de medida, pulse la tecla setup.
- 2. Pulse la tecla ▲, ▼, ◀ o ▶ para resaltar <u>pH Channel</u>, <u>pH/ISE Channel</u>, <u>COND</u> <u>Channel</u> o <u>DO/RDO Channel</u> y pulse la tecla **f3 (select)**.
- 3. Pulse la tecla ▲ o ▼ para resaltar <u>Method</u> y pulse la tecla **f3 (select)** para acceder a la lista de métodos.
- 4. Pulse la tecla ▲ o ▼ para resaltar un método protegido que se cargará en el modo de medida y pulse la tecla **f2 (load)** (cargar).
 - a. Los métodos protegidos tienen la fecha, la hora y el modo visibles en la lista de métodos.
- 5. El medidor continuará automáticamente con el modo de medida.

Copia de un método

Utilice la función de copiar para guardar los ajustes de medida, calibración y alarma específicos del canal de un método existente a uno nuevo, de manera que los ajustes del método original se conserven y los ajustes del método nuevo puedan modificarse. Cuando se utilice la función de copiar para crear un nuevo método, debe crearse una nueva contraseña para que se guarde el nuevo método.

- 1. En el modo de medida, pulse la tecla setup.
- 2. Pulse la tecla ▲, ▼, ◀ o ▶ para resaltar <u>pH Channel</u>, <u>pH/ISE Channel</u>, <u>COND</u> <u>Channel</u> o <u>DO/RDO Channel</u> y pulse la tecla **f3 (select)**.
- Pulse la tecla ▲ o ▼ para resaltar <u>Method</u> y pulse la tecla f3 (select) para acceder a la lista de métodos.
- Pulse la tecla ▲ o ▼ para resaltar un método protegido que se copiará en un método protegido abierto y pulse la tecla f3 (options) (opciones).
 - a. Los métodos protegidos tienen la fecha, la hora y el modo visibles en la lista de métodos.
- Utilice la ventana emergente de introducción de números para introducir la contraseña específica para el método seleccionado.
 - a. Pulse la tecla f3 (edit) para acceder a la ventana de introducción de números.
 - b. Pulse la tecla ▲, ▼, ◀ o ► para resaltar un número, pulse la tecla f3 (enter) para seleccionar el número y repítalo hasta que la contraseña correcta se muestre en la parte superior de la ventana de introducción de números.
 - c. Pulse la tecla f2 (done) para aceptar la contraseña y salir de la ventana de introducción de números.
- 6. Pulse la tecla f2 (accept) para enviar la contraseña para el método seleccionado.
- 7. Pulse la tecla ▲ o ▼ para resaltar Copy (Copiar) y pulse la tecla f2 (accept).
- 8. Pulse la tecla ▲ o ▼ para resaltar un método abierto y pulse la tecla **f2 (accept)**. a. Los métodos abiertos no tienen fecha, hora o modo visibles en la lista de métodos.
- 9. Utilice la ventana emergente de introducción de números para crear una contraseña específica para el nuevo método.

- a. Pulse la tecla f3 (edit) para acceder a la ventana de introducción de números.
- b. Pulse la tecla ▲, ▼, ◀ o ▶ para resaltar un número, pulse la tecla f3 (enter) para seleccionar el número y repítalo hasta que la contraseña deseada se muestre en la parte superior de la ventana.
- c. Pulse la tecla f2 (done) para aceptar la contraseña y salir de la ventana de introducción de números.
- 10. Pulse la tecla f2 (accept) para enviar la nueva contraseña para el método nuevo.
- 11. El medidor volverá a la lista de métodos y el número de método nuevo tendrá la fecha, la hora y el modo visibles en la lista de métodos.

Edición o eliminación de un método

Utilice la función de editar para modificar los ajustes de medida, calibración y alarma específicos del canal de un método protegido existente. Utilice la función de eliminar para convertir un método protegido existente en un método abierto.

- 1. En el modo de medida, pulse la tecla setup.
- 2. Pulse la tecla ▲, ▼, ◀ o ▶ para resaltar <u>pH Channel</u>, <u>pH/ISE Channel</u>, <u>COND</u> <u>Channel</u> o <u>DO/RDO Channel</u> y pulse la tecla **f3 (select)**.
- Pulse la tecla ▲ o ▼ para resaltar Method y pulse la tecla f3 (select) para acceder a la lista de métodos.
- Pulse la tecla ▲ o ▼ para resaltar un método protegido y pulse la tecla f3 (options).
 a. Los métodos protegidos tienen la fecha, la hora y el modo visibles en la lista de métodos.
- Utilice la ventana emergente de introducción de números para introducir la contraseña específica para el método seleccionado.
 - a. Pulse la tecla f3 (edit) para acceder a la ventana de introducción de números.
 - b. Pulse la tecla ▲, ▼, ◀ o ► para resaltar un número, pulse la tecla f3 (enter) para seleccionar el número y repítalo hasta que la contraseña correcta se muestre en la parte superior de la ventana de introducción de números.
 - c. Pulse la tecla f2 (done) para aceptar la contraseña y salir de la ventana de introducción de números.
- Pulse la tecla f2 (accept) para enviar la contraseña para el método seleccionado.
- Pulse la tecla ▲ o ▼ para resaltar Edit o Delete (Eliminar) y pulse la tecla f2 (accept).
 - a. Si se ha seleccionado Edit:
 - i. Realice los pasos adecuados para establecer los ajustes de medida, calibración y alarma específicos del canal en el método seleccionado.
 - ii. Cuando haya editado los ajustes necesarios, pulse la tecla **f1 (back)** para volver a la lista de métodos.
 - b. Si se ha seleccionado Delete:
 - i. Pulse la tecla **f2 (yes)** (sí) para confirmar la eliminación del método seleccionado.
 - ii. El método se eliminará y el medidor continuará automáticamente con el modo de medida.

Menú de modo y ajustes

Opciones generales del menú de modo y ajustes

Tipos de lectura

La opción Read Type (Tipo de lectura) determina cómo muestra, registra y exporta las medidas el medidor. Asegúrese de que se activan Instrument Settings y exportación de datos adecuados en el menú de ajustes del instrumento antes de registrar o exportar medidas.

- Auto-Read:
 - Pulse la tecla measure (esc) para iniciar la medida. El icono AR parpadea a medida que el valor cambiante se estabiliza. Cuando es estable, el icono AR deja de parpadear y la medida queda fija en la pantalla hasta que se pulsa de nuevo la tecla measure (esc).
 - Cuando está activada, la medida estable se guarda automáticamente en el registro de datos y se exporta a una impresora o un ordenador.
- Continuous:
 - Los valores de medida se actualizan constantemente en la pantalla y el icono Stabilizing o Ready indica el estado de estabilidad de la medida.
 - Cuando esté activada, pulse la tecla log/print para guardar una medida en el registro de datos y exportarla a una impresora o un ordenador.
- Timed:
 - Los valores de medida se actualizan constantemente en la pantalla y el icono Stabilizing o Ready indica el estado de estabilidad de la medida.
 - Cuando está activada, las medidas se guardan en el registro de datos y se exportan a una impresora o un ordenador en el intervalo de tiempo seleccionado a partir de 3 segundos (00:00:03) hasta 24 horas (24:00:00).

Información de ajuste de estabilidad

El ajuste Stability (Estabilidad) determina cuándo se reconoce una medida como estable por el medidor, a través de un algoritmo que evalúa las fluctuaciones de medida. Cuando se selecciona Smart Stability, se utiliza la resolución de medida junto con el algoritmo. A continuación, se muestran los valores aproximados de mV por minuto, estimados en condiciones ideales de medida. Los valores son solo teóricos. Los valores reales variarán dependiendo de las condiciones concretas de medida.

	do 0 1 o 1 díaito			
resolucion de 0,1 o 1 dígito 7,6 (Estability (Estabilidad rápida)		7,6	Fast Stability (Estabilidad rápida)	7,6
Resolución de 0,01 o 2 dígitos 2,9 Medium Stability (Estabilidad media)	de 0,01 o 2 dígitos s	2,9	Medium Stability (Estabilidad media)	2,9
Resolución de 0,001 o 3 dígitos 1,0 Slow Stability (Estabilidad lenta)	de 0,001 o 3 dígitos s	1,0	Slow Stability (Estabilidad lenta)	1,0

nuto

Ajuste promedio

El ajuste promedio permite al operador seleccionar Averaging (Promedio) de Automatic Smart para obtener una estabilidad de medida más rápida o la opción de apagado para no activar ningún promedio de medida una vez cumplidos los criterios de estabilidad

Ajuste de alarma límite

El ajuste Limit Alarm (Alarma límite) permite que las medidas puedan evaluarse a través de valores límite superiores y/o inferiores. Si una medida está por encima del valor límite superior o por debajo del valor límite inferior, la alarma se activará. Por ejemplo, cuando el límite superior esté establecido en 8,50 pH, la alarma se activará cuando haya una lectura de 8,51 pH o más en el modo de medida. Active el ajuste del sonido de la alarma en el menú de ajustes del instrumento para habilitar un sonido audible cuando se active la alarma.

Ajuste de alarma de calibración vencida

El ajuste de alarma de calibración vencida permite al operador establecer un intervalo de calibración personalizado en horas. Si una calibración no se realiza dentro del intervalo de calibración especificado, la alarma se activará. Active el ajuste del sonido de la alarma en el menú Instrument Settings para habilitar un sonido audible cuando se active la alarma.

Menú Mode and Settings para el canal pH/ISE

Utilice el menú Mode and Settings para personalizar los ajustes de medida, calibración y alarma para el canal de pH de los medidores de pH Orion Star A211, los medidores de pH/conductividad Orion Star A215 y los medidores de pH/RDO/OD Orion Star A216. Los ajustes predeterminados del medidor aparecen en **negrita**.

Parámetro	Ajustes	Detalles
Measure Mode (Modo de medida)	 pH mV RmV ORP 	Establezca el modo de medida mostrado, el modo seleccionado determina qué ajustes adicionales se muestran y el tipo de calibración que realizará el medidor
Read Type (Tipo de lectura)	 Auto Continuous Timed 00:00:03 	Establezca el tipo de lectura para determinar la manera en que el medidor toma una medida y cuándo se envía esta al registro de datos y/o dispositivo externo
Resolution (Resolución) (solo pH)	1 decimal 0,1 2 decimales 0,01 3 decimales 0,001	Establezca la resolución del valor de medida de pH
Buffer Group (Grupo búfer) (solo pH)	• USA • DIN	Establezca el grupo búfer para el reconocimiento de búfer automático durante las calibraciones de pH USA: búferes pH 1,68, 4,01, 7,00, 10,01 y 12,46 DIN: búferes pH 1,68, 4,01, 6,86 y 9,18
Stability (Estabilidad)	 Smart Stability Fast Medium Slow 	Establezca cuándo una medida se reconoce como estable. Smart Stability compensa las condiciones de medida y optimiza el tiempo de respuesta del medidor
Averaging (Promedio)	Off Automatic Smart	Establezca el promedio para obtener una estabilidad de medida más rápida. Automatic Start compensa las condiciones de medida y optimiza el tiempo de respuesta del medidor
Limit Alarm (Alarma límite)	 Alarma límite Off (apagado) On (encendido) Alarm Settings (Ajustes de alarma) High (alto) High/Low (alto/bajo) Low (bajo) High Limit (Límite superior) Low Limit (Límite inferior) 	Establezca la alarma límite en apagado o encendido. Cuando la alarma límite esté encendida, establezca los ajustes de alarma como alto, alto/bajo o bajo y después introduzca los valores límites adecuados de la medida. La alarma límite se activa cuando la medida está por encima del límite superior o por debajo del límite inferior.
Cal Due Alarm (Alarma de calibración vencida) (solo pH, RmV y ORP)	 Cal Due Alarm On Off Cal Due Limit (Límite de calibración vencido) 12 Hrs (12 horas) 	Establezca la alarma de calibración vencida en apagado o encendido. Cuando la alarma de calibración vencida esté encendida, introduzca el intervalo de tiempo de calibración en horas (de 1 a 9998 horas) La alarma de calibración vencida se activa después de que pase el tiempo introducido sin haber realizado una calibración.

Ajuste de grupo búfer

Los medidores de pH Orion Star A211, los medidores de pH/conductividad Orion Star A215 y medidores de pH/RDO/OD Orion Star A216 son capaces de reconocer automáticamente valores de búfer de pH dentro del conjunto de búfer seleccionado durante una calibración de pH. Durante la calibración de pH, el medidor utiliza el conjunto de búfer de pH seleccionado y la lectura mV sin procesar del electrodo de pH en el búfer para reconocer y mostrar el valor del búfer debe estar dentro de una unidad de pH (aproximadamente ±59 mV) del valor mV teórico del búfer para que el medidor lo reconozca.

Conjunto de búfer USA			
Búfer de pH	Intervalo mV		
1,68	De +255 mV a +374 mV		
4,01	De +117 mV a +236 mV		
7,00	De -59 mV a +59 mV		
10,01	De -237 mV a -119 mV		
12,46	De -382 mV a -264 mV		

Conjunto de búfer DIN		
Búfer de pH	Intervalo mV	
1,68	De +255 mV a +374 mV	
4,01	De +117 mV a +236 mV	
6,86	De -51 mV a +67 mV	
9,18	De -189 mV a -70 mV	

Prueba de un electrodo de pH para el reconocimiento de búfer automático

Utilice el siguiente procedimiento para confirmar que la lectura mV sin procesar del electrodo de pH está dentro de una unidad de pH (±59 mV) de la lectura mV teórica del búfer de pH y, por lo tanto, verifique que el electrodo de pH en uso es capaz de realizar el reconocimiento de búfer automático.

- 1. Prepare el electrodo de pH de acuerdo con el manual de electrodos. Establezca el modo de medida del medidor en mV.
- Enjuague el electrodo de pH con agua destilada, séquelo con un paño sin pelusa y colóquelo en un búfer de pH 4,01 a aproximadamente 25 °C.
- 3. Espere a que la medida se estabilice y registre el valor mV del búfer de pH 4,01 cuando esté estable la medida.
- 4. Retire el electrodo de pH del búfer de pH 4,01.
- 5. Enjuague el electrodo de pH con agua destilada, séquelo con un paño sin pelusa y colóquelo en un búfer de pH 7,00 a aproximadamente 25 °C.
- 6. Espere a que la medida se estabilice y registre el valor mV del búfer de pH 7,00 cuando esté estable la medida.
- 7. La lectura mV del electrodo de pH debería ir de +117 mV a +236 mV en el búfer de pH 4 y de -59 mV a +59 mV en el búfer de pH 7. Si las lecturas mV están en intervalos correctos, el electrodo de pH es capaz de realizar el reconocimiento de búfer automático. Si las lecturas mV no están en intervalos correctos, introduzca manualmente los valores de búfer de pH durante la calibración de pH.

Menú Mode and Settings para el canal pH/ISE

Utilice el menú Mode and Settings para personalizar los ajustes de medida, calibración y alarma para el canal de pH/ISE de los medidores de pH/ISE Orion Star A214. Los ajustes predeterminados del medidor aparecen en negrita.

Parámetro	Ajustes	Detalles
Measure Mode	• pH • mV • RmV • ISE	Establezca el modo de medida mostrado, el modo seleccionado determina qué ajustes adicionales se muestran y el tipo de calibración que realizará el medidor
Read Type	Auto Continuous Timed 00:00:03	Establezca el tipo de lectura para determinar la manera en que el medidor toma una medida y cuándo se envía esta al registro de datos y/o dispositivo externo
Resolution (solo pH)	1 decimal 0,1 2 decimales 0,01 3 decimales 0,001	Establezca la resolución del valor de medida de pH
Buffer Group (solo pH)	• USA • DIN	Establezca el grupo búfer para el reconocimiento de búfer automático durante las calibraciones de pH USA: búferes pH 1,68, 4,01, 7,00, 10,01 y 12,46 DIN: búferes pH 1,68, 4,01, 6,86 y 9,18
Resolution (solo ISE)	1 dígito significativo 2 dígitos significativos 3 dígitos significativos	Establezca la resolución del valor de medida de concentración de iones (ISE)
Measure Unit (solo ISE)	 ppm Percentage (%) M ppb mg/L None (Ninguno) 	Establezca las unidades mostradas con el valor de medida de concentración de iones (ISE)
Blank Correct (Correc. blanco) (solo ISE)	 Yes (Sí) No 	Establezca la función de corrección con blanco automática en apagado o encendido para las medidas ISE. Cuando la función está encendida, se utiliza un algoritmo para compensar la respuesta no lineal del electrodo en estándares y muestras de bajo nivel para obtener una estabilidad de medida mejorada
Stability	 Smart Stability Fast Medium Slow 	Establezca cuándo una medida se reconoce como estable. Smart Stability compensa las condiciones de medida y optimiza el tiempo de respuesta del medidor
Averaging	Off Automatic Smart	Establezca el promedio para obtener una estabilidad de medida más rápida. Automatic Start compensa las condiciones de medida y optimiza el tiempo de respuesta del medidor
Limit Alarm	 Limit Alarm Off On Alarm Settings High High/Low Low High Limit Low Limit 	Establezca la alarma límite en apagado o encendido. Cuando la alarma límite esté encendida, establezca los ajustes de alarma como alto, alto/bajo o bajo y después introduzca los valores límites adecuados de la medida. La alarma límite se activa cuando la medida está por encima del límite superior o por debajo del límite inferior.
Cal Due Alarm (solo pH, RmV, ORP e ISE)	 Cal Due Alarm On Off Cal Due Limit 12 Hrs 	Establezca la alarma de calibración vencida en apagado o encendido. Cuando la alarma de calibración vencida esté encendida, introduzca el intervalo de tiempo de calibración en horas (de 1 a 9998 horas) La alarma de calibración vencida se activa después de que pase el tiempo introducido sin haber realizado una calibración.

Parámetro	Ajustes	Detalles
Low Lv Stability (Estabilidad de bajo nivel)	• Off • On	Establezca la función de estabilidad de bajo nivel en apagado o encendido para las calibraciones ISE. Cuando la función está encendida, el tiempo de estabilización mínimo se incrementa (~3-5 minutos) para mejorar la precisión de los estándares de calibración de bajo nivel

Ajuste de grupo búfer

Los medidores de pH/ISE Orion Star A214 son capaces de reconocer automáticamente valores de búfer de pH dentro del conjunto de búfer seleccionado durante una calibración de pH. Durante la calibración de pH, el medidor utiliza el conjunto de búfer de pH seleccionado y la lectura mV sin procesar del electrodo de pH en el búfer para reconocer y mostrar el valor del búfer en la temperatura medida. La lectura mV sin procesar del electrodo de pH en el búfer debe estar dentro de una unidad de pH (aproximadamente ±59 mV) del valor mV teórico del búfer para que el medidor lo reconozca.

Ajuste de corrección con blanco

Los medidores pH/ISE Orion Star A214 ofrecen la opción de utilizar una corrección con blanco (no lineal) para las medidas ISE cuando se realiza una calibración multipunto. La función de corrección con blanco automático utiliza un algoritmo para compensar la falta de linealidad del electrodo de ion selectivo en las muestras y estándares de bajo nivel.

Con la función de corrección con blanco activada, el medidor decide si la corrección con blanco es la mejor estrategia de medida al analizar la respuesta del electrodo durante una calibración multipunto. No debe ejecutarse un blanco por separado. Gráficamente, la corrección con blanco es equivalente al dibujo de una suave curva a través de los tres puntos más bajos de la calibración multipunto y la extrapolación a una concentración de cero basada en el supuesto del comportamiento nernstiano del electrodo.

Las calibraciones multipunto en el límite inferior son convenientes cuando la respuesta de un electrodo de ion selectivo no es lineal y no puede caracterizarse con una calibración de uno o dos puntos. Normalmente, esto se considera como una pendiente de electrodo baja. Normalmente, el electrodo tiene un comportamiento nernstiano, pero se está observando el efecto de un blanco. Consulte la siguiente figura.



Este blanco puede ser un blanco reactivo verdadero, trazas de ion de muestra en reactivos o puede ser el valor de "cieno" del electrodo. Podría ser también una interferencia de los reactivos que se hace aparente a bajos niveles de ion de muestra, o podría ser una combinación de estos efectos. La versión explicada de la ecuación de Nernst usada tradicionalmente para la corrección con blanco es la siguiente:

E = Eo + S * registro (C + b)

donde b es el blanco

En una calibración multipunto, se genera un conjunto de ecuaciones y se evalúan las relaciones entre ellas. Por ejemplo, las ecuaciones generadas en una calibración de tres puntos serían:

 $E_1 = Eo + S * registro (C_1 + b)$ $E_2 = Eo + S * registro (C_2 + b)$ $E_3 = Eo + S * registro (C_3 + b)$

El medidor evalúa las relaciones entre los tres potenciales E_1 , E_2 y E_3 , y las tres concentraciones C_1 , C_2 y C_3 . Si las relaciones indican que la corrección con blanco es deseable, se calculará un blanco automáticamente y la falta de linealidad se corregirá de forma nernstiana. Si no se cumplen las condiciones adecuadas, el blanco se establece en cero y cada segmento de la calibración multipunto se trata de forma independiente.

Cuando se dan las tres siguientes condiciones, se invoca la corrección con blanco.

- La concentración del primer estándar es cero, o la pendiente del electrodo entre el primer y el segundo estándar es menos que la pendiente entre el segundo y el tercero.
- 2. Las diferencias de potencial entre los puntos son significativas. Por ejemplo, E_3 E_1 > 10 mV
- 3. El algoritmo de corrección con blanco converge en valores de blanco y pendiente razonables. Las condiciones en los pasos 1 y 2 evitan el fallo de convergencia en la mayoría de situaciones. Sin embargo, la pendiente obtendrá cualquier valor necesario para la convergencia para un valor de blanco calculado de 3 x C₃.

Cuando se implementa una corrección con blanco, el valor de pendiente para el electrodo puede quedar fuera del intervalo de valores normalmente considerados aceptables durante una calibración ordinaria. Para obtener mejores resultados, los estándares de calibración deberían estar próximos al intervalo para las concentraciones de muestra previstas y deberían soportar la concentración de muestra prevista. Cuando no se dan las condiciones 1, 2 y 3, los datos de calibración son gestionados por el método de calibración multipunto descrito anteriormente.

En las calibraciones con más de tres puntos, se utiliza una combinación de métodos. Si los tres puntos más bajos cumplen los criterios y se utiliza la calibración multipunto para el resto de puntos, se utiliza la corrección con blanco automática. La pendiente calculada en el algoritmo de corrección con blanco automática y las pendientes para cada segmento adicional se utilizan para calcular la pendiente media.

Ajuste de estabilidad de bajo nivel

Los medidores de pH/ISE Orion Star A214 ofrecen la opción de utilizar la función de estabilidad de bajo nivel al calibrar electrodos de ion selectivo. La función de estabilidad de bajo nivel mejora la precisión de las medidas ISE de concentración baja al ajustar el tiempo para los puntos de calibración de los estándares de bajo nivel, lo que permite un mayor tiempo de estabilización para el electrodo en los estándares de calibración. El tiempo de estabilización normalmente se amplía aproximadamente de tres a cinco minutos por punto de calibración, pero esto variará según las lecturas reales del electrodo durante la calibración.

Menú Mode and Settings para el canal COND

Utilice el menú Mode and Settings para personalizar los ajustes de medida, calibración y alarma para el canal de conductividad de los medidores de conductividad Orion Star A212 y los medidores de pH/conductividad Orion Star A215. Los ajustes predeterminados del medidor aparecen en negrita.

Parámetro	Ajustes	Detalles
Measure Mode	 Conductivity Salinity TDS Resistivity 	Establezca el modo de medida mostrado, el modo seleccionado determina qué ajustes adicionales se muestran y el tipo de calibración que realizará el medidor
Read Type	Auto Continuous Timed 00:00:03	Establezca el tipo de lectura para determinar la manera en que el medidor toma una medida y cuándo se envía esta al registro de datos y/o dispositivo externo
Cell K (Celda K)	Celda K 0,4750	Introduzca el valor de la constante de celda nominal (K) del sensor de conductividad para la calibración de conductividad automática
Type (Tipo) (Solo salinidad)	 Practical Salinity (Salinidad práctica) Sea Water (Agua marina) 	Establezca el tipo de medida de salinidad como salinidad práctica (psu) o agua marina natural (ppm)
Factor TSD (Factor STD) (solo STD)	Factor STD 0,49	Introduzca el valor del factor STD para las medidas de sólidos disueltos totales
Temp. ref. (Temperatura ref.)	 +5 °C +10 °C +15 °C +20 °C +25 °C 	Establezca la temperatura de referencia para las medidas compensadas de temperatura. Las lecturas se ajustan a la temperatura de referencia seleccionada cuando el ajuste de compensación de temperatura está activado
Temp. Comp. (Comp. temperatura)	 Off Linear (Lineal) nLFn nLFu EP 	Establezca el tipo de compensación de temperatura como lineal, nLFn (agua natural no lineal), nLFu (agua ultrapura no lineal), EP (compensación desactivada con la alarma cuando los valores están fuera de los requisitos de EP para agua ultrapura) o apagado
Temp. Coeff. (Coef. temperatura)	Temp. Coefficient 2,10	Introduzca el coeficiente de temperatura utilizado con el ajuste de compensación de temperatura lineal
Stability	 Smart Stability Fast Medium Slow 	Establezca cuándo una medida se reconoce como estable. Smart Stability compensa las condiciones de medida y optimiza el tiempo de respuesta del medidor
Averaging	Off Automatic Smart	Establezca el promedio para obtener una estabilidad de medida más rápida. Automatic Start compensa las condiciones de medida y optimiza el tiempo de respuesta del medidor
Limit Alarm	 Limit Alarm Off On Alarm Settings High High/Low Low High Limit Low Limit 	Establezca la alarma límite en apagado o encendido. Cuando la alarma límite esté encendida, establezca los ajustes de alarma como alto, alto/bajo o bajo y después introduzca los valores límites adecuados de la medida. La alarma límite se activa cuando la medida está por encima del límite superior o por debajo del límite inferior.
Cal Due Alarm (Solo conductividad)	 Cal Due Alarm On Off Cal Due Limit 12 Hrs 	Establezca la alarma de calibración vencida en apagado o encendido: introduzca el intervalo de tiempo de calibración en horas (de 1 a 9998 horas) La alarma de calibración vencida se activa después de que pase el tiempo definido sin realizar una calibración
Cell Type (Tipo de celda)	 Standard (Estándar) USP (Farmacopea estadounidense) 	Establezca el tipo de sensor de conductividad como estándar (la mayoría de sensores) o USP (sensores de dos celdas en agua ultrapura)
Ajuste de celda K

Los conductímetros Orion Star A212 y los medidores de pH/conductividad Orion Star A215 son capaces de reconocer automáticamente el estándar de conductividad 100 µS, el estándar de conductividad 1413 µS y el estándar de conductividad 12,9 mS de Thermo Scientific Orion cuando el valor de la constante de celda nominal (K) del sensor de conductividad se introduce en el menú de configuración.

Número de catálogo	Descripción
011008	Estándar de Conductividad Orion 100 μ S, 5 x 60 mL
011007	Estándar de Conductividad Orion 1413 µS, 5 x 60 mL
01100710	Estándar de conductividad Orion 1413 μ S, 10 bolsas de uso individual
011006	Estándar de conductividad 12,9 mS, 5 x 60 mL
01100610	Estándar de conductividad Orion 12,9 mS, 10 bolsas de uso individual

Los valores de la constante de celda nominal (K) para los sensores de conductividad Thermo Scientific Orion compatibles con los medidores de conductividad Star serie A210 se enumeran a continuación.

Número de catálogo	Descripción	Rango de medición	Constante de celda nominal
013005MD	Sensor de temperatura/conductividad de 4 celdas Orion con cable de 1,5 metros (5 pies)	De 1 μS/cm a 200 mS/cm	0,475 cm⁻¹
013010MD	Sensor de temperatura/conductividad de 4 celdas Orion con cable de 3 metros (10 pies)	De 1 μS/cm a 200 mS/cm	0,475 cm⁻¹
013020MD	Sensor de temperatura/conductividad de 4 celdas Orion con cable de 6 metros (20 pies)	De 1 μS/cm a 200 mS/cm	0,475 cm⁻¹
013025MD	Sensor de temperatura/conductividad de 4 celdas Orion con cable de 10 metros (32 pies)	De 1 μS/cm a 200 mS/cm	0,475 cm⁻¹
013605MD	Sensor de temperatura/conductividad de 4 celdas Orion con cable de 1,5 metros (5 pies)	De 10 μS/cm a 200 mS/cm	0,55 cm ⁻¹
013610MD	Sensor de temperatura/conductividad de 4 celdas Orion con cable de 3 metros (10 pies)	De 10 μS/cm a 200 mS/cm	0,55 cm ⁻¹
013016MD	Sensor de temperatura/conductividad de 2 celdas de agua pura Orion con cable de 1,5 metros (5 pies)	De 0,01 μS/cm a 300 μS/cm	0,1 cm ⁻¹
011510MD	Sensor de temperatura/conductividad de 2 celdas Orion con cable de 3 metros (10 pies)	De 10 μS/cm a 200 mS/cm	1,0 cm ⁻¹
011050MD	Sensor de temperatura/conductividad de 2 celdas Orion con cable de 1,5 metros (5 pies)	De 1 μS/cm a 20 mS/cm	1,0 cm ⁻¹
018020MD	Sensor de conductividad de 2 celdas de gran rango Orion con cable de 1,5 metros (5 pies)	De 10 μS/cm a 2000 mS/cm	10 cm-1

Temperatura de referencia, Compensación de temperatura y Ajustes de coeficiente

Los conductímetros Orion Star A212 y los medidores de pH/conductividad Orion Star A215 ofrecen varias opciones para medidas de conductividad compensada por temperatura. La temperatura tiene un gran efecto sobre el valor de conductividad de una solución. La función de compensación de temperatura permite al medidor utilizar las lecturas de temperatura y conductividad de muestra para calcular y mostrar la conductividad prevista para la muestra en la temperatura de referencia seleccionada, como 25 °C. Para utilizar de forma precisa la función de compensación de temperatura para las muestras que se medirán, establezca la temperatura de referencia requerida como 5 °C, 10 °C, 15 °C, 20 °C o 25 °C y utilice un sensor de conductividad con una sonda de temperatura ATC integrada. Las opciones para la compensación de temperatura son:

 Lineal: utiliza el valor de coeficiente de temperatura para aplicar el factor de corrección de porcentaje constante a cada cambio gradual de temperatura. Algunos valores de coeficiente lineal comunes son:

Solución (de 25 °C a 50 °C)	Coeficiente de temperatura (%/°C)
Jarabe de azúcar	5,64
Agua ultrapura	4,55
Ácido sulfúrico al 98 %	2,84
Sal (NaCl)	2,12
Amoniaco diluido	1,88
NaOH al 5 %	1,72
HCl al 10 %	1,32
Ácido sulfúrico al 5 %	0,96

- nLFn (agua no desgasificada ultrapura no lineal): aplica un factor de corrección no constante a las muestras de agua pura que tienen tiempos de respuesta variables a los cambios de temperatura. Este modo está previsto para las lecturas compensadas de temperatura de aguas de baja conductividad en equilibrio con el dióxido de carbono en el aire, como las lecturas de agua pura cerca de 1 µS/cm a 25 °C y aguas naturales con una composición comparable a las aguas superficiales, de pozo o subterráneas naturales.
- nLFu (agua no desgasificada ultrapura no lineal): aplica un factor de corrección no constante a las muestras de agua pura que tienen tiempos de respuesta variables a los cambios de temperatura. Este modo está previsto para las lecturas compensadas de temperatura de agua ultrapura que no contiene ni aire ni dióxido de carbono, como el agua ultrapura (resistencia de 18 megaohmios o superior) directamente desde la fuente sin aireación.
- EP: no se aplica ninguna corrección de temperatura (la compensación de temperatura está desactivada) y se muestra una advertencia si el valor de conductividad medido está por encima del requisito EP para el agua pura en la temperatura de muestra medida.
- Apagado: no se aplica ninguna corrección de temperatura y el valor de conductividad real se muestra en la temperatura de muestra medida.

Ajuste del tipo de salinidad

Los conductímetros Orion Star A212 y los medidores de pH/conductividad Orion Star A215 ofrecen dos tipos de salinidad: salinidad práctica y agua marina natural. Las medidas de salinidad práctica se basan en la relación de la medida de muestras con la lectura de una solución de cloruro potásico (KCI) estándar a 15 °C. Los conductímetros Orion Star serie A210 compensarán automáticamente la temperatura e informarán del resultado de salinidad práctica previsto como unidades de salinidad práctica (psu) a 15 °C al utilizar un sensor de conductividad con temperatura integrada. Las medidas de agua marina natural utilizan una convención histórica conocida como UNESCO 1966 y los conductímetros Orion Star serie A210 informarán del resultado del agua marina natural como partes por mil (ppm).

Ajuste del factor STD

Los conductímetros Orion Star A212 y los medidores de pH/conductividad Orion Star A215 miden el STD como la cantidad total de material inorgánico disuelto en una solución. El material inorgánico disuelto lleva una corriente que se mide con el sensor de conductividad. Puesto que hay una relación directa entre la conductividad y el STD, las lecturas de conductividad se utilizan para estimar la presencia de material inorgánico a través del factor STD introducido en el menú de configuración.

El método estándar de determinación STD implica evaporar la muestra para secarla a 180 °C y pesar el residuo. El factor STD se calcula al tomar el peso del residuo y dividirlo por la conductividad de muestra. Las posteriores lecturas de conductividad se multiplican por el factor STD para determinar el valor STD de la muestra.

Ajuste del tipo de celda

Los conductímetros Orion Star A212 y los medidores de pH/conductividad Orion Star A215 aceptan sensores de conductividad de 2 y 4 celdas y permiten que el tipo de sensor de conductividad se establezca como Estándar (la mayoría de sensores de conductividad) o USP (para el sensor de conductividad de agua ultrapura, número de catálogo 013016MD, al desactivar la compensación de temperatura).

Menú Mode and Settings para el canal OD/RDO

Utilice el menú Mode and Settings para personalizar los ajustes de medida, calibración y alarma para el canal de oxígeno disuelto de los medidores de RDO/OD Orion Star A213 y los medidores de pH/RDO/OD Orion Star A216. Los ajustes predeterminados del medidor aparecen en negrita.

Parámetro	Ajustes	Detalles
Measure Mode	 DO - Polarographic (OD: polarográfico) RDO Auto 	Establezca el tipo de sensor de oxígeno disuelto conectado al medidor como polarográfico OD u óptico RDO, o seleccione Auto para que el medidor detecte automáticamente el tipo de sensor de oxígeno disuelto conectado y actualice el modo de medida mostrado.
Measure Unit	mg/L% sat	Establezca las unidades mostradas con el valor de medida del oxígeno disuelto
Resolution (% de saturación)	1 decimal 1 2 decimales 0,1	Establezca la resolución del valor de medida del oxígeno disuelto como saturación en porcentaje
Resolution (mg/L)	1 decimal 0,1 2 decimales 0,01	Establezca la resolución del valor de medida del oxígeno disuelto como mg/L
Read Type	Auto Continuous Timed 00:00:03	Establezca el tipo de lectura para determinar la manera en que el medidor toma una medida y cuándo se envía esta al registro de datos y/o dispositivo externo
Baro Press (Presión barométrica)	• Auto • Manual 760,0 mm Hg	Establezca la fuente de compensación de presión barométrica automática como el barómetro interno (Automática) o como un valor de presión barométrica introducido manualmente (Manual)
Salinity Correct (Corrección de salinidad)	• Manual 0,0	Introduzca el valor de salinidad de las muestras para la corrección de salinidad automática de las medidas de oxígeno disuelto.
Stability	 Smart Stability Fast Medium Slow 	Establezca cuándo una medida se reconoce como estable. Smart Stability compensa las condiciones de medida y optimiza el tiempo de respuesta del medidor
Averaging	Off Automatic Smart	Establezca el promedio para obtener una estabilidad de medida más rápida. Automatic Start compensa las condiciones de medida y optimiza el tiempo de respuesta del medidor
Limit Alarm	 Limit Alarm Off On Alarm Settings High High/Low Low High Limit Low Limit 	Establezca la alarma límite en apagado o encendido. Cuando la alarma límite esté encendida, seleccione los ajustes de alarma como alto, alto/bajo o bajo y después introduzca los valores límites de la medida. La alarma límite se activa cuando una medida está por encima del límite superior introducido o por debajo del límite inferior.
Cal Due Alarm	 Cal Due Alarm On Off Cal Due Limit 12 Hrs 	Establezca la alarma de calibración vencida en apagado o encendido. Cuando la alarma de calibración vencida esté encendida, introduzca el intervalo de tiempo de calibración en horas (de 1 a 9998 horas) La alarma de calibración vencida se activa después de que pase el tiempo introducido sin haber realizado una calibración.
RDO Cap Life (Vida del tapón RDO) (solo RDO)	RDO Cap Life 0.0 Serial Number (Número de serie) 0	Vea el resto de la vida del tapón del sensor de OD óptico RDO expresada en días y el número de serie

Ajuste de compensación de la presión barométrica

Los medidores de RDO/OD Orion Star A213 y los medidores de pH/RDO/OD Orion Star A216 tienen un barómetro interno que puede utilizarse para realizar lecturas automáticas de oxígeno disuelto compensado por presión. La entrada manual de la presión barométrica puede utilizarse para las lecturas automáticas de oxígeno disuelto compensado por presión, que resultan útiles cuando el oxígeno disuelto se mide con un sensor sumergido o un recipiente presurizado. El valor de presión debe introducirse como mm Hg.

1 mm Hg = 0,03937 in Hg = 1,3332 hPa (mBar) = 0,01934 PSI

Ajuste de corrección de salinidad

Los medidores de RDO/OD Orion Star A213 y los medidores de pH/RDO/OD Orion Star A216 son capaces de realizar correcciones de salinidad automáticas para lecturas de oxígeno disuelto (medidas como mg/L) cuando el valor de salinidad de la muestra se introduce manualmente como ppm (partes por mil) en el menú de configuración.

Conducti- vidad a 20 °C (mS/cm)	Valor de corrección de salinidad (ppm)	Conducti- vidad a 20 °C (mS/cm)	Valor de corrección de salinidad (ppm)	Conducti- vidad a 20 °C (mS/cm)	Valor de corrección de salinidad (ppm)
5	3	20	13	35	25
6	4	21	14	36	25
7	4	22	15	37	26
8	5	23	15	38	27
9	6	24	16	39	28
10	6	25	17	40	29
11	7	26	18	42	30
12	8	27	18	44	32
13	8	28	19	46	33
14	9	29	20	48	35
15	10	30	21	50	37
16	10	31	22	52	38
17	11	32	22	54	40
18	12	33	23	56	42
19	13	34	24		

Calculado utilizando las tablas oceanográficas internacionales, volumen 1, Instituto Nacional de Oceanografía de Gran Bretaña, Womley, Godaming, Surrey, Inglaterra y Unesco, París 1971

Menú Temperature

Los medidores de sobremesa Orion Star serie A210 ofrecen un menú de temperatura utilizado para personalizar los ajustes de temperatura de cada canal mediante la introducción manual de un valor de temperatura, la selección de las unidades de temperatura como °C o °F y la realización de una calibración de temperatura en una sonda ATC, sensor de conductividad o sensor de oxígeno disuelto con temperatura integrada. Los medidores de pH/RDO/OD Orion Star A216 también ofrecen la opción de utilizar una fuente de entrada de temperatura para ambos canales de medida.

Modelo de medidor	odelo de Canal Menú Temperature		Opciones de menú		
		 Manual Temp Value (Valor Temp manual) 	Introducir valor de temperatura		
Medidor de pH	pH Channel	 Temperature Unit (Unidad de temperatura) 	Celsius Fahrenheit		
		 Temperature Calibration (Calibración de temperatura) 	• ATC		
	0010	Manual Temp Value	Introducir valor de temperatura		
Star A212	COND Channel	Temperature Unit	Celsius Fahrenheit		
		Temperature Calibration	• ATC		
		Temperature Unit	Celsius		
Medidor de RDO/OD Star A213	DO/RDO Channel	 Temperature Calibration (solo sensor de OD polarográfico) 	 Solution Temperature (Temperatura de solución) Membrane Temperature (Temperatura de membrana) Sol & Mem Temp (Temperatura sol. y memb.) 		
Medidor de	pH/ISE Channel	Manual Temp Value	 Introducir valor de temperatura 		
pH/ISE Star A214		Temperature Unit	Celsius Fahrenheit		
		Temperature Calibration	• ATC		
		Manual Temp Value	Introducir valor de temperatura		
	pH Channel	Temperature Unit	Celsius Fahrenheit		
Medidor de pH/		Temperature Calibration	• ATC		
Star A215		Manual Temp Value	Introducir valor de temperatura		
	COND Channel	Temperature Unit	Celsius Fahrenheit		
		Temperature Calibration	• ATC		
		Manual Temp Value	Introducir valor de temperatura		
		Temperature Unit	Celsius Fahrenheit		
Medidor de pH/RDO/OD Star A216	pH Channel	Temperature Calibration	 ATC Solution Temperature Membrane Temperature Sol & Mem Temperature ATC, Sol & Mem Temp 		
		Temperature Input	ATC. DO Probe Manual		
		Temperature Unit	Celsius Fahrenheit		
	DO/RDO Channel	 Temperature Calibration (solo sensor de OD polarográfico) 	Solution TemperatureMembrane TemperatureSol & Mem Temp		

Valor de temperatura manual

- 1. En el modo de medida, pulse la tecla setup.
- Pulse la tecla ▲, ▼, ◀ o ▶ para resaltar <u>pH Channel</u>, <u>pH/ISE Channel</u> o <u>COND</u> <u>Channel</u> y pulse la tecla f3 (select).
- 3. Pulse la tecla ▲ o ▼ para resaltar <u>Temperature</u> y pulse la tecla f3 (select).
- Pulse la tecla ▲ o ▼ para resaltar <u>Manual Temp Value</u> y pulse la tecla f3 (select).
- 5. Pulse la tecla **f3 (edit)** para acceder a la ventana emergente de introducción de números e introduzca el valor de temperatura de muestra.
 - a. Pulse la tecla ▲, ▼, ◀ o ▶ para resaltar un número, punto decimal o signo negativo y después pulse la tecla
 f3 (enter) para seleccionar el elemento resaltado. Repita hasta que el valor de la temperatura de muestra se visualice en la parte superior de la ventana.



- b. Pulse la tecla f2 (done) para guardar el valor y salir de la ventana de introducción de números.
- Pulse la tecla f1 (back) para volver al menú de temperatura principal y pulse la tecla measure (esc) para volver al modo de medida.

Unidad de temperatura

- 1. En el modo de medida, pulse la tecla setup.
- 2. Pulse la tecla ▲, ▼, ◀ o ▶ para resaltar <u>pH Channel</u>, <u>pH/ISE Channel</u>, <u>COND</u> <u>Channel</u> o <u>DO/RDO Channel</u> y pulse la tecla **f3 (select)**.
- 3. Pulse la tecla ▲ o ▼ para resaltar <u>Temperature</u> y pulse la tecla f3 (select).
- Pulse la tecla ▲ o ▼ para resaltar <u>Temperature Unit</u> y pulse la tecla f3 (select).
- Pulse la tecla ▲ o ▼ para resaltar <u>Celsius</u> o <u>Fahrenheit</u> y pulse la tecla f3 (select).
- 6. Pulse la tecla measure (esc) para salir al modo de medida.

Calibración de temperatura

Al utilizar los medidores de sobremesa Orion Star serie A210, puede realizarse una calibración de temperatura en una sonda ATC, en un sensor de conductividad con temperatura integrada o en un sensor de oxígeno disuelto polarográfico con temperatura integrada cuando se conecta el sensor o la sonda al medidor. Utilice la función de calibración de temperatura solo cuando sea necesario, ya que el medidor tiene una precisión de temperatura relativa de $\pm 0,1$ °C y las sondas ATC tienen precisiones variables, normalmente entre $\pm 0,5$ °C y ± 2 °C. Puesto que la compensación de temperatura calculada durante la calibración se aplica a todas las futuras medidas de temperatura, recalibre la temperatura si utiliza una sonda ATC diferente.

- 1. En el modo de medida, pulse la tecla setup.
- Pulse la tecla ▲, ▼, ◀ o ▶ para resaltar <u>pH Channel</u>, <u>pH/ISE Channel</u>, <u>COND Channel</u> o <u>DO/RDO Channel</u> y pulse la tecla **f3 (select)**.
- 3. Pulse la tecla ▲ o ▼ para resaltar <u>Temperature</u> y pulse la tecla **f3 (select)**.
- Pulse la tecla ▲ o ▼ para resaltar <u>Temperature Calibration</u> y pulse la tecla f3 (select).
- 5. Pulse la tecla \blacktriangle o \triangledown para resaltar <u>ATC</u> y pulse la tecla **f3 (select)**.
 - a. Si utiliza el medidor Star A213, pulse la tecla ▲ o ▼ para resaltar <u>Solution</u> <u>Temperature</u>, <u>Membrane Temperature</u> o <u>Sol & Mem Temperature</u> y pulse la tecla f3 (select).
 - b. Si utiliza el medidor Star A216, pulse la tecla ▲ o ▼ para resaltar <u>ATC</u>, <u>Solution</u> <u>Temperature</u>, <u>Membrane Temperature</u>, <u>Sol & Mem Temperature</u> o <u>ATC</u>, <u>Sol & Mem Temp</u> y pulse la tecla **f3 (select)**.
- Coloque la sonda y/o sensor ATC en una solución con una temperatura estable y conocida. Deben utilizarse dos termómetros trazables según NIST para medir y verificar la temperatura.
- Espere a que el valor de temperatura del medidor se estabilice y, a continuación, pulse la tecla f3 (edit) para acceder a la ventana emergente de introducción de números e introduzca el valor de temperatura.
 - a. Pulse la tecla ▲, ▼, ◀ o ▶ para resaltar un número, punto decimal o signo negativo y después pulse la tecla f3 (enter) para seleccionar el elemento resaltado. Repita hasta que el valor deseado de temperatura se muestre en la parte superior de la pantalla.
 - b. Pulse la tecla f2 (done) para salir de la ventana de introducción de números.
- 8. Pulse la tecla f2 (accept) para guardar el valor de temperatura introducido.
- 9. Pulse la tecla **f1 (meas)** (medir) para salir al modo de medida.

Calibración de temperatura para medidores con múltiples entradas de temperatura

Cuando un sensor de oxígeno disuelto polarográfico con dos entradas de temperaturas se conecta con un medidor de RDO/OD Orion Star A213 o un medidor de pH/RDO/OD Orion Star A216, la entrada de temperatura de la solución y la entrada de temperatura de la membrana pueden calibrarse juntas o por separado.

Cuando una sonda ATC y un sensor de oxígeno disuelto polarográfico se conectan al medidor de pH/RDO/OD Orion Star A216, la calibración de temperatura para ambas entradas puede hacerse al mismo tiempo.

Entrada de temperatura

Al utilizar el medidor de pH/RDO/OD Orion Star A216, puede establecerse un sensor de oxígeno disuelto con temperatura integrada como la fuente de entrada de temperatura para todas las medidas. Por ejemplo, al utilizar un electrodo de pH sin una sonda ATC, el sensor de oxígeno disuelto con temperatura integrada puede utilizarse como la fuente de temperatura para las medidas de pH, siempre que tanto el electrodo de pH como el sensor de oxígeno disuelto estén colocados en la misma solución.

- 1. En el modo de medida, pulse la tecla setup.
- 2. Pulse la tecla ▲, ▼, ◀ o ► para resaltar <u>pH Channel</u> y pulse la tecla f3 (select).
- 3. Pulse la tecla \blacktriangle o \triangledown para resaltar <u>Temperature</u> y pulse la tecla **f3 (select)**.
- 4. Pulse la tecla ▲ o ▼ para resaltar <u>Temperature Input</u> y pulse la tecla f3 (select).
- 5. Pulse la tecla ▲ o ▼ para resaltar <u>ATC</u>, <u>DO Probe</u> o y pulse la tecla **f3 (select)**.
- 6. Pulse la tecla measure (esc) para salir al modo de medida.

Menú de configuración Instrument Settings

Utilice el menú Instrument Settings para actualizar los ajustes del medidor para la transferencia de datos, registro de datos, fecha y hora, idioma, sonidos, velocidad rotatoria, contraste de la pantalla, apagado automático, ID de usuario e ID de la muestra.

Ajustes del	Opciones	Detalles
instrumento		
Export Data (Exportar datos)	 Printing (Impresión) Off On Comm Setup (Ajuste general) USB RS232 Data Format (Formato de datos) Printer (Impresora) PC (CSV) Comm Config (Configuración general) 1200 9600 2400 19 200 4800 38 400 	Acceda al menú de exportación de datos para establecer los ajustes de transferencia para la exportación de la información de calibración y medida a una impresora u ordenador. Active la impresión para permitir la transferencia de datos del medidor a un dispositivo externo, seleccione el formato de transferencia de datos como impresora (texto estándar) o PC (texto delimitado por comas) y establezca la velocidad de baudios del medidor para que coincida con la del ordenador o la impresora. El medidor detectará automáticamente cuando un cable RS232 o USB se conecte a un puerto.
Data log (Registro de datos)	OffOn	Active el registro de datos para guardar los datos de medida en el registro de datos de acuerdo con el tipo de lectura seleccionado
Date and Time (Fecha y hora)	 Date (Fecha) DD/MM/YY (DD/MM/AA) MM/DD/AA Date: 01/01/12 Time (Hora) Reloj de 12 horas Time: 02:30 p. m. Reloj de 24 horas Time: 14:30 	Utilice el ajuste de fecha para seleccionar el formato de fecha como día/mes/año (DD/MM/AA) o mes/día/año (MM/DD/AA) y después introduzca los valores de mes, día y año Utilice el ajuste de la hora para establecer el formato de hora como 12 o 24 horas y para introducir los valores de hora y minuto
Idioma	 English Español Deutsch Français Italiano 中又(Chino) 	Establezca el idioma que se utilizará en la interfaz del medidor. Hay idiomas adicionales disponibles (incluidos portugués y coreano) junto con las actualizaciones de software en nuestro sitio web en <u>www.thermoscientific.com/OrionMeters</u>
Key Press Beep (Sonido al pulsar teclas)	OffOn	Active o desactive el sonido audible cada vez que se pulse una tecla en el medidor
Alarm Beep (Sonido de alarma)	OffOn	Active o desactive el sonido audible cuando se activa una alarma
Stirrer (Agitador)		Ajuste los parámetros de la velocidad del agitador desde uno (menor velocidad) a cinco (mayor velocidad) al utilizar la sonda de agitador de Orion Star
Contrast (Contraste)		Aumente o reduzca el contraste de la pantalla para mejorar la visualización de esta en distintas condiciones de iluminación

Ajustes del instrumento	Opciones	Detalles
Auto Shut Off (Apagado automático)	OffOn	Active esta función para apagar automáticamente el medidor cuando no se pulsa ninguna tecla en 20 minutos
User ID (ID de usuario)	ABCDE	Introduzca un valor de ID de usuario (hasta seis caracteres) utilizando la ventana de entrada alfanumérica
Sample ID (ID de la muestra)	 Off Manual Auto Increment (Incremento automático) 	Establezca la ID de la muestra en apagado, manual (hasta seis caracteres alfanuméricos) o incremento automático (hasta seis números que incrementan automáticamente para cada medida)

Menú Log View

Utilice el menú Log View para acceder al registro de datos y al registro de calibración Toda la información de los registros de calibración y datos se protegen gracias a la memoria no volátil del medidor, de manera que la información almacenada se guarda en el medidor incluso cuando no está conectado al suministro eléctrico.

Registro de datos

Los medidores de sobremesa Orion Star A210 ofrecen un registro de datos de 2000 puntos. Cada punto incluye medidas mostradas de forma activa, dependiendo del modelo del medidor y la configuración de pantalla, con marca de fecha y hora. Cuando la función de registro de datos está desactivada, el tipo de lectura seleccionado para cada canal mostrado (Auto-Read, Continuous o Timed) determina cómo se guarda el punto en el registro de datos.

- 1. En el modo de medida, pulse la tecla setup.
- 2. Pulse la tecla ▲, ▼, ◀ o ▶ para resaltar <u>View Log</u> y pulse la tecla f3 (select).
- 3. Pulse la tecla ▲ o ▼ para resaltar <u>Data Log</u> y pulse la tecla f2 (select).
- 4. El medidor mostrará una lista de los puntos de registro de datos. La lista muestra el número secuencial de puntos y la fecha y la hora en que se guardó este.
- 5. Para ver la información de la medida para un punto individual:
 - a. Pulse la tecla ▲ o ▼ para resaltar un punto y, a continuación, pulse la tecla
 f2 (select).
 - b. Pulse la tecla f1 (back) para volver a la lista de los puntos del registro de datos.
- 6. Para exportar el registro de datos a una impresora u ordenador:
 - a. Pulse la tecla **f3 (options)**, la tecla ▲ o ▼ para resaltar Log Export (Exportación de registro) y luego pulse la tecla **f2 (accept)**.
 - b. Pulse la tecla ▲ o ▼ para resaltar <u>Today</u>, <u>Last 10</u>, <u>Range</u> o <u>All</u> (Hoy, Últimos 10 o Todos) y luego pulse la tecla **f2 (accept)**.
 - Si se selecciona Range (Intervalo), pulse la tecla ▲ o ▼ para resaltar el número de los puntos primero (<u>from</u> (desde)) y último (<u>to</u> (hasta)) que desea exportar, utilice la ventana emergente de introducción de números para editar los números; a continuación, pulse la tecla f2 (accept) para guardar el rango introducido.
 - c. Pulse la tecla **f2 (yes)** para confirmar que el medidor está conectado a una impresora o un ordenador. El medidor automáticamente volverá a la lista de registro de datos cuando se complete la exportación.
- 7. Para borrar la memoria del registro de datos:
 - a. Pulse la tecla **f3 (options)**, la tecla ▲ o ▼ para resaltar Log Clear (Borrar registro) y luego pulse la tecla **f2 (accept)**.
 - b. Pulse la tecla **f2 (yes)** para confirmar la eliminación de todo el registro de datos de la memoria del medidor.
 - c. Pulse la tecla f3 (edit), utilice la ventana emergente de introducción de números para introducir la contraseña predeterminada del medidor de 111111 y, a continuación, pulse la tecla f2 (done).
 - d. Pulse la tecla f2 (accept) y espere a que se borre el registro de datos.

Registro de calibración

Los medidores Orion Star serie A210 guardan hasta diez de las calibraciones más recientes por canal.

- 1. En el modo de medida, pulse la tecla setup.
- Pulse la tecla ▲, ▼, ◀ o ▶ para resaltar View Log y pulse la tecla f3 (select).
- Pulse la tecla ▲ o ▼ para resaltar Calibration Log (Registro de calibración) y pulse la tecla f2 (select).
- Pulse la tecla ▲ o ▼ para resaltar <u>pH-Channel</u>, <u>pH/ISE-Channel</u>, Conductivity-Channel o DO/RDO-Channel y pulse la tecla **f2 (select)**.
 - a. Si se selecciona pH-Channel, pulse la tecla ▲ o ▼ para resaltar <u>pH</u>, <u>RmV</u> u <u>ORP</u> y pulse la tecla **f2 (select)**.
 - b. Si se selecciona pH/ISE-Channel, pulse la tecla ▲ o ▼ para resaltar <u>pH</u>, <u>RmV</u>, <u>ORP</u> o <u>ISE</u> y pulse la tecla **f2 (select)**.
 - c. Si se selecciona Conductivity-Channel, pulse la tecla ▲ o ▼ para resaltar <u>Conductivity</u>, <u>Resistivity</u>, <u>TDS</u> o <u>Salinity</u> y pulse la tecla **f2 (select)**.
 - d. Si se selecciona Canal OD/RDO, pulse la tecla ▲ o ▼ para resaltar <u>DO</u> o <u>RDO</u> y pulse la tecla **f2 (select)**.
- El medidor mostrará una lista de calibraciones para el canal y modo de calibración seleccionados. La lista muestra el número secuencial del registro de calibración y la fecha y la hora en que se completó la calibración.
- Para ver información detallada de una calibración individual, pulse la tecla ▲ o ▼ para resaltar una calibración y, a continuación, pulse la tecla f2 (select).
 - a. Para exportar la calibración a una impresora u ordenador pulse la tecla **f2 (print)** (imprimir).

Therne Scientific (c)	2011
ATTI TH	
Netar a/a	201036
UN DEV	1.5.1
Calibration Report	
No. of the second second	
12/2//14 10043800	
POARS 1	4 44
	2100 5
	280.0 EV
resperature	40×0 G
POINT A	4 44 44
20	174 PR
Terri a ra futa	25.0.7
penperature	0.0+11 10
reant a	7 00 -10
P4	1 . U.V. (241)
Der Mennen and Frank	2.4.V (B) V 2.6. (A) (A)
being d	dated to
FOINT N	10.01.00
and a second	-122 6
Hanna Paratura	25.0.0
Hoine A	**** *
	10:00 -0
1	200.4 mV
Table A Table 1 (1994)	75.0.2
witness 1	57.7 4
#2.000 P	90.5 8
#1 com #	97.5 8
ol-med	50.4 4
TT:	3.5 mt
N 2	0.0 10
F .)	0.0 mV
24	1.7 MV
Average Slope	98.0 %
Calibration	\$3
Operator	
Righatione	

Ejemplo de registro de calibración exportado, Formato de datos de impresora:

Menú Diagnostics

Utilice el menú Diagnostics para iniciar un restablecimiento del medidor, llevar a cabo una prueba de estabilidad del electrodo, realizar una prueba automática o ver las revisiones de software y el número de serie del módulo y medidor.

- Factory Reset (Rest. valores fábrica): todos los ajustes se restablecen a los valores predeterminados de fábrica. El registro de datos, el registro de calibración y los métodos se eliminan. Utilice la contraseña predeterminada del medidor de 111111 para finalizar el restablecimiento.
- User Reset (Rest. usuario): todos los ajustes de medida, calibración y alarma específicos del canal, así como los ajustes del instrumento se restablecen a los parámetros predeterminados de fábrica y se eliminan los métodos. El registro de datos y el registro de calibración se conservan.
- Self Test (Prueba automática): prueba la precisión del medidor. El medidor también realiza una prueba automática en cada encendido.
- Stability Test (Prueba de estabilidad): prueba la estabilidad de un electrodo al medir la lectura de entrada sin procesar por minuto de ruido.
- About Meter (Acerca del medidor): muestra el modelo del medidor, el número de serie y la revisión de software.

Procedimiento de prueba automática del medidor

- 1. En el modo de medida, pulse la tecla setup.
- 2. Pulse la tecla \blacktriangle , \triangledown , \triangleleft o \triangleright para resaltar <u>Diagnostics</u> y pulse la tecla **f3 (select)**.
- Pulse la tecla ▲ o ▼ para resaltar <u>Self Test</u> y pulse la tecla f2 (select).
- 4. Desconecte todos los electrodos y sensores del medidor, conecte el tapón de acortamiento BNC a la entrada BNC si procede y pulse la tecla **f2 (yes)**.
- 5. Pulse la tecla f2 (yes) para iniciar la prueba automática.
- El medidor realizará la prueba automática. Cuando se muestra <u>Self Test Passed</u> (Prueba automática superada), pulse la tecla **f1 (esc)** y el medidor pasará automáticamente al modo de medida.

Procedimiento de verificación del medidor (solo canales pH y pH/ISE)

- Tras completar la prueba automática, el medidor volverá al modo de medida. Mantenga todos los electrodos y sensores desconectados del medidor y el tapón de acortamiento BNC conectado a la entrada BNC del medidor.
- Establezca el modo de medida a pH. El medidor debería leer una constante de 7000 pH ±0,002 pH.
 - a. Si el medidor no lee una constante de 7000 pH ±0,002 pH, realice una calibración de pH de un punto con el tapón de acortamiento BNC conectado a la entrada BNC. Establezca el valor de pH en 7000 y el valor de pendiente en 100,0. Consulte el apartado de calibración de pH para obtener instrucciones detalladas.
- Establezca el modo de medida a mV. El medidor debería leer una constante de 0,0 mV ±0,2 mV.
 - a. Si el medidor no lee una constante de 0,0 mV ±0,2 mV, con el tapón de acortamiento BNC conectado a la entrada BNC, póngase en contacto con el soporte técnico.

Procedimiento de prueba de estabilidad del electrodo

- 1. En el modo de medida, pulse la tecla setup.
- 2. Pulse la tecla ▲, ▼, ◀ o ▶ para resaltar <u>Diagnostics</u> y pulse la tecla f3 (select).
- Pulse la tecla ▲ o ▼ para resaltar <u>Stability Test</u> y pulse la tecla f2 (select).
- Pulse la tecla ▲ o ▼ para resaltar <u>pH-Channel</u>, <u>pH/ISE-Channel</u>, <u>Conductivity-Channel</u> o <u>DO/RDO-Channel</u> y pulse la tecla **f2 (next)** (siguiente).
- 5. Asegúrese de que el electrodo o el sensor que se va a probar esté conectado al canal seleccionado en el paso anterior y pulse la tecla **f2 (yes)**.
- 6. Coloque el electrodo o el sensor en una solución adecuada a una temperatura estable y pulse la tecla **f2 (yes)**.
 - a. Se recomienda el búfer de pH 7,00 de Thermo Scientific Orion (número de catálogo 910107, 910760 o 910710) para probar un electrodo de pH.
 - b. Se recomienda el estándar ORP de Thermo Scientific Orion (número de catálogo 967901 o 967961) para probar un electrodo ORP.
 - c. Se recomienda un estándar de calibración con una concentración moderada de ion de interés para la prueba de un electrodo de ion selectivo (ISE).
 - d. Se recomienda el estándar 1413 pS de Thermo Scientific Orion (número de catálogo 011007 o 01100710) para probar la mayoría de sensores de conductividad. Se recomienda el estándar 100 pS de Thermo Scientific Orion (número de catálogo 011008) para probar la mayoría de sensores de conductividad de agua ultrapura.
 - e. Se recomienda una manga de calibración de aire saturado con agua para probar los sensores de oxígeno disuelto polarográficos y ópticos RDO.
- 7. El medidor realizará la prueba de estabilidad y mostrará la lectura de entrada sin procesar por minuto de ruido en la pantalla.
- El medidor mostrará <u>Stability Test Passed</u> (Prueba de estabilidad superada) o <u>Stability Test Failed</u> (Prueba de estabilidad no superada). Pulse la tecla **f1 (esc)** para salir al modo de medida.

4

CAPÍTULO 4 El uso del canal pH o pH/ISE

Utilice el medidor de pH Orion Star A211, el medidor de pH/conductividad Orion Star A215 o el medidor de pH/RDO/OD Orion Star A216 para medir pH, mV, mV relativo, ORP y temperatura. Utilice el medidor de pH/ISE Orion Star A214 para medir pH, concentración de iones utilizando un ISE (electrodo de ion selectivo), mV, mV relativo, ORP y temperatura. El siguiente capítulo ofrece instrucciones sobre la conexión de los electrodos, la calibración del sistema y la medida de estos parámetros.

Preparación del electrodo y el medidor

- Familiarícese con el funcionamiento general del medidor y prepare el medidor y los accesorios de conformidad con las instrucciones del <u>Capítulo 2: Principios básicos del</u> <u>medidor</u>.
 - Se recomienda colocar los electrodos en el soporte portaelectrodos conectado al medidor para un fácil movimiento dentro y fuera de los contenedores durante la calibración, la medida y el almacenamiento.
- 2. Consulte las instrucciones del <u>Capítulo 3: Menús de configuración del medidor</u> para personalizar los ajustes de medida, calibración y alarma dentro del menú de configuración del canal pH o el canal pH/ISE y los ajustes de transferencia de datos, registro de datos, fecha y hora, idioma, sonidos, velocidad rotatoria, contraste de la pantalla, apagado automático, ID de usuario e ID de la muestra dentro del menú de configuración de ajustes del instrumento.
- Conecte todos los electrodos, sondas y sensores necesarios a las entradas del medidor adecuadas.
 - a. Conecte un electrodo de ion selectivo, ORP, o pH a la entrada BNC del medidor.
 - b. Conecte una sonda de temperatura ATC a la entrada MiniDIN de 8 pines del medidor.

- c. Conecte un electrodo de media celda de referencia a la entrada de punta del medidor (para utilizar junto con un electrodo de detección de media celda con conector BNC).
- d. Conecte la sonda de agitador (número de catálogo 096019) a la entrada del medidor del agitador.
- 4. Prepare todos electrodos conectados para el uso tal como se indica en los manuales de los electrodos.

Procedimiento de calibración de pH

Los medidores de pH Orion Star A211, los medidores de pH/ISE Orion Star A214, los medidores de pH/conductividad Orion Star A215 y los medidores de pH/RDO/OD Orion Star A216 pueden llevar a cabo una calibración de pH utilizando de uno a cinco búferes con facilidad para seguir las instrucciones que aparecen en la pantalla para ayudar con el proceso de calibración.

Utilice siempre búferes de pH nuevos y seleccione los búferes que soportan el pH de muestra y están alejados de una a tres unidades de pH. Prepare el electrodo de pH de acuerdo con las instrucciones del manual de electrodos. Conecte todos los electrodos que se deben utilizar a las entradas del medidor adecuadas.

Asegúrese de que el modo de medida del medidor se haya establecido en pH. Para los medidores Star A215 y Star A216, pulse la tecla **f3 (channel)** hasta que se muestre el canal de pH en el modo de medida.

- 5. Pulse la tecla f1 (cal) para iniciar la calibración.
 - a. Si se muestra más de un canal en el modo de medida, pulse la tecla ▲ o ▼ para resaltar el <u>pH-Channel</u> y pulse la tecla **f2 (select)**.
- Enjuague el electrodo de pH y los demás electrodos en uso con agua destilada, séquelos con suavidad con un paño sin pelusas para quitar el exceso de agua y colóquelos en el búfer de pH.
- 7. Cuando el electrodo y el búfer estén listos, pulse la tecla f3 (start) (iniciar).
 - a. Si utiliza una sonda de agitador, esta empezará el agitado cuando se pulse la tecla **f3 (start)** y detendrá el agitado cuando la lectura se estabilice.
- 8. Espere hasta que el valor pH del medidor se estabilice y lleve a cabo alguna de las acciones siguientes:
 - a. Pulse la tecla **f2 (accept)** para aceptar el valor de búfer de pH determinado por la función de reconocimiento de búfer automático.

o bien

- b. Pulse la tecla **f3 (edit)** para acceder a la ventana emergente de introducción de números e introduzca manualmente el valor de búfer de pH.
 - Pulse la tecla ▲, ▼, ◀ o ▶ para resaltar un número o punto decimal, pulse la tecla f3 (enter) para seleccionar el elemento resaltado y repítalo hasta que el valor deseado se muestre en la parte superior de la pantalla.
 - ii. Pulse la tecla **f2 (done)** para guardar el valor y salir de la ventana de introducción de números.
 - iii. Pulse la tecla **f2 (accept)** para aceptar el valor introducido manualmente.

- 9. Pulse la tecla f2 (next) para continuar con el búfer de pH siguiente y repita los pasos del 2 al 4 o pulse la tecla f3 (cal done) (cal. realizada) para guardar y finalizar la calibración. Si se utilizan los cinco búferes, la calibración se guardará y finalizará una vez se haya aceptado el quinto valor.
 - a. Al realizar una calibración de un punto, pulse la tecla f2 (accept) para aceptar el valor de pendiente mostrado o pulse la tecla f3 (edit), utilice la ventana emergente de introducción de números para introducir un nuevo valor de pendiente, pulse la tecla f2 (done) y, a continuación, pulse la tecla f2 (accept).
- El medidor mostrará el resumen de la calibración, incluida la pendiente, y exportará los datos al registro de calibración. Pulse la tecla f1 (meas) para continuar con el modo de medida o pulse la tecla f2 (print) para exportar los datos de calibración a una impresora u ordenador.

Edición de calibración de pH

Cuando se muestra el resumen de calibración, se puede seleccionar la opción de edición de calibración para reparar los puntos individuales sin tener que realizar una calibración completa.

- 1. En la pantalla de resumen de calibración (después del paso 5 del procedimiento de calibración de pH), pulse la tecla **f3 (cal edit)** (edición de calibración)
- Pulse la tecla ▲ o ▼ para resaltar el punto de calibración que hay que editar y pulse la tecla f2 (select).
- 3. Pulse la tecla ▲ o ▼ para resaltar <u>Remeasure</u>, <u>Edit</u> o <u>Delete</u> (Volver a medir, Editar o Eliminar) y pulse la tecla **f2** (select).
 - a. Si se selecciona Remeasure, repita los pasos del 2 al 4 en el Procedimiento de calibración pH.
 - b. Si se selecciona Edit, pulse la tecla f3 (edit), utilice la ventana emergente de introducción de números para introducir un nuevo valor de búfer de pH, pulse la tecla f2 (done) y, a continuación, pulse la tecla f2 (accept). Seleccione otro punto de calibración para editar o pulse la tecla f1 (back).
 - c. Si se selecciona Delete, el punto de calibración se eliminará. Si la eliminación del punto lo cambia por una calibración de un punto, introduzca la pendiente con la ventana emergente de introducción de números.
- El medidor mostrará un resumen de la calibración actualizado y exportará los datos al registro de calibración. Pulse la tecla f1 (meas) para continuar con el modo de medida o pulse la tecla f2 (print) para exportar los datos de calibración a una impresora u ordenador.

Procedimiento de calibración ORP (Modo mV relativo)

Los medidores de pH Orion Star A211, los medidores de pH/ISE Orion Star A214, los medidores de pH/conductividad Orion Star A215 y los medidores de pH/RDO/OD Orion Star A216 pueden llevar a cabo una calibración ORP en el modo mV relativo (RmV) utilizando un estándar con facilidad para seguir las instrucciones que aparecen en la pantalla para ayudar con el proceso de calibración.

Utilice el estándar ORP de Thermo Scientific Orion (número de catálogo 967901 o 967961) o cualquier otro estándar ORP según lo desee. Prepare el electrodo de acuerdo con las instrucciones del manual de electrodos. Conecte todos los electrodos que se deben utilizar a las entradas del medidor adecuadas.

Asegúrese de que el modo de medida del medidor se haya establecido en RmV. Para los medidores Star A215 y Star A216, pulse la tecla **f3 (channel)** hasta que se muestre el canal de RmV en el modo de medida.

- 1. Pulse la tecla f1 (cal) para iniciar la calibración.
 - a. Si se muestra más de un canal en el modo de medida, pulse la tecla ▲ o ▼ para resaltar el <u>pH-Channel</u> y pulse la tecla **f2 (select)**.
- Enjuague el electrodo y los demás electrodos en uso con agua destilada, séquelos con suavidad con un paño sin pelusas para quitar el exceso de agua y colóquelos en el estándar.
- 3. Cuando el electrodo y el estándar estén listos, pulse la tecla f3 (start).
 - a. Si utiliza una sonda de agitador, esta empezará el agitado cuando se pulse la tecla **f3 (start)** y detendrá el agitado cuando la lectura se estabilice.
- 4. Espere a que el valor mV del medidor se estabilice (indicado por el icono intermitente **stabilizing** o el icono estático **ready**) y realice una de las siguientes acciones:
 - a. Pulse la tecla f2 (accept) para aceptar el mV, o
 - Pulse la tecla f3 (edit) para acceder a la ventana emergente de introducción de números e introduzca manualmente el valor mV previsto en la temperatura medida.
 - Pulse la tecla ▲, ▼, ◀ o ▶ para resaltar un número, punto decimal o signo negativo y pulse la tecla f3 (enter) para seleccionar el elemento resaltado y repítalo hasta que el valor deseado se muestre en la parte superior de la pantalla.
 - ii. Pulse la tecla **f2 (done)** para guardar el valor y salir de la ventana de introducción de números.
 - iii. Pulse la tecla **f2 (accept)** para aceptar el valor introducido manualmente.
- 5. El medidor mostrará el resumen de la calibración, incluida la compensación mV, y exportará los datos al registro de calibración. Pulse la tecla f1 (meas) para continuar con el modo de medida o pulse la tecla f2 (print) para exportar los datos de calibración a una impresora u ordenador.

Procedimiento de calibración ORP (E_H Units)

Los medidores de pH Orion Star A211, medidores de pH/ISE Orion Star A214, medidores de pH/conductividad Orion Star A215 y medidores de pH/RDO/OD Orion Star A216 pueden llevar a cabo una calibración ORP automática de un punto al valor EH (resistente al medio ambiente) (relativo al electrodo de hidrógeno estándar, 420 mV a 25 °C) cuando se utilizan los siguientes elementos:

- Estándar ORP de Thermo Scientific Orion (número de catálogo 967901 o 967961)
- Electrodo ORP Thermo Scientific Orion (número de catálogo 9678BNWP, 9778BNWP, 9180BNMD o 9179BNMD)
- Solución de llenado 4 M KCI Thermo Scientific Orion (número de catálogo 900011) o gel 4 M KCI (en el electrodo 9179BNMD)

El estándar ORP de Orion está clasificado como no peligroso, estable y no cambiará con el tiempo, de manera que es un estándar ideal para cualquier sistema de medida ORP. Utilice siempre un estándar de ORP nuevo. Prepare el electrodo ORP de acuerdo con las instrucciones del manual de electrodos. Conecte todos los electrodos que se deben utilizar a las entradas del medidor adecuadas.

Asegúrese de que el modo de medida del medidor se haya establecido en ORP. Para los medidores Star A215 y Star A216, pulse la tecla **f3 (channel)** hasta que se muestre el canal de ORP en el modo de medida.

- 1. Pulse la tecla f1 (cal) para iniciar la calibración.
 - a. Si se muestra más de un canal en el modo de medida, pulse la tecla ▲ o ▼ para resaltar el <u>pH-Channel</u> y pulse la tecla **f2 (select)**.
- Enjuague el electrodo y los demás electrodos en uso con agua destilada, séquelos con suavidad con un paño sin pelusas para quitar el exceso de agua y colóquelos en el estándar.
- 3. Cuando el electrodo y el estándar estén listos, pulse la tecla f3 (start).

Si utiliza una sonda de agitador, esta empezará el agitado cuando se pulse la tecla **f3 (start)** y detendrá el agitado cuando la lectura se estabilice.

- Espere a que el valor mV del medidor se estabilice (indicado por el icono intermitente stabilizing o el icono estático ready) y pulse f2 (accept) cuando el valor mV sea estable.
- El medidor mostrará el resumen de la calibración, incluida la compensación mV, y exportará los datos al registro de calibración. Pulse la tecla f1 (meas) para continuar con el modo de medida o pulse la tecla f2 (print) para exportar los datos de calibración a una impresora u ordenador.

Procedimiento de calibración ISE

Los medidores pH/ISE Orion Star A214 pueden llevar a cabo una calibración de concentración de iones utilizando de uno a cinco estándares con facilidad para seguir las instrucciones que aparecen en la pantalla para ayudar con el proceso de calibración.

Utilice siempre estándares nuevos y seleccione los estándares que sujetan la concentración de la muestra y están alejados una década (diez veces) de la concentración. Para preparar los estándares de calibración de forma precisa de una solución de existencias, se recomienda una dilución seriada mediante pipetas. Si se añade el ISA a las muestras, el mismo ISA debe añadirse a todos los estándares antes de la calibración para garantizar un factor de dilución uniforme. Si se utiliza más de un estándar para la calibración, empiece con el estándar de concentración más bajo y trabaje hasta llegar al final al estándar de calibración más alto.

Prepare el electrodo de ion selectivo de acuerdo con las instrucciones del manual de electrodos. Conecte todos los electrodos que se deben utilizar a las entradas del medidor adecuadas. Asegúrese de que el modo de medida del medidor se haya establecido en ISE.

- 1. Pulse la tecla f1 (cal) para iniciar la calibración.
- Enjuague el electrodo de ion selectivo y los demás electrodos en uso con agua destilada, séquelos con suavidad con un paño sin pelusas para quitar el exceso de agua y colóquelos en el estándar.
- 3. Cuando el electrodo y el estándar estén listos, pulse la tecla f3 (start).
 - a. Si utiliza una sonda de agitador, esta empezará el agitado cuando se pulse la tecla **f3 (start)** y detendrá el agitado cuando la lectura se estabilice.
- 4. Espere hasta que el valor del medidor se estabilice y lleve a cabo alguna de las acciones siguientes:
 - a. Pulse la tecla f2 (accept) para aceptar el valor de concentración mostrado, o
 - b. Pulse la tecla **f3 (edit)** para acceder a la ventana emergente de introducción de números e introduzca manualmente el valor de concentración.
 - i. Pulse la tecla ▲, ♥, ◀ o ▶ para resaltar un número o punto decimal, pulse la tecla f3 (enter) para seleccionar el elemento resaltado y repítalo hasta que el valor deseado se muestre en la parte superior de la pantalla.
 - ii. Pulse la tecla **f2 (done)** para guardar el valor y salir de la ventana de introducción de números.
 - iii. Pulse la tecla **f2 (accept)** para aceptar el valor introducido manualmente.
- 5. Pulse la tecla f2 (next) para continuar con el estándar siguiente y repita los pasos del 2 al 4 o pulse la tecla f3 (cal done) para guardar y finalizar la calibración. Si se utilizan los cinco estándares, la calibración se guardará y finalizará una vez se haya aceptado el quinto valor.
 - a. Al realizar una calibración de un punto, pulse la tecla f2 (accept) para aceptar el valor de pendiente mostrado o pulse la tecla f3 (edit), utilice la ventana emergente de introducción de números para introducir un nuevo valor de pendiente, pulse la tecla f2 (done) y, a continuación, pulse la tecla f2 (accept).
- 6. El medidor mostrará el resumen de la calibración, incluida la pendiente, y exportará los datos al registro de calibración. Pulse la tecla f1 (meas) para continuar con el modo de medida o pulse la tecla f2 (print) para exportar los datos de calibración a una impresora u ordenador.

Edición de calibración ISE

Cuando se muestra el resumen de calibración, se puede seleccionar la opción de edición de calibración para reparar los puntos individuales sin tener que realizar una calibración completa.

- 1. En la pantalla de resumen de calibración (después del paso 5 del Procedimiento de calibración ISE), pulse la tecla **f3 (cal edit)** (editar calibración).
- Pulse la tecla ▲ o ▼ para resaltar el punto de calibración que hay que editar y pulse la tecla f2 (select).
- Pulse la tecla ▲ o ▼ para resaltar <u>Remeasure</u>, <u>Edit</u> o <u>Delete</u> y pulse la tecla f2 (select).
 - a. Si se selecciona Remeasure, repita los pasos del 2 al 4 en el Procedimiento de calibración ISE.
 - b. Si se selecciona Editar, pulse la tecla f3 (edit), utilice la ventana emergente de introducción de números para introducir un nuevo valor estándar, pulse la tecla f2 (done) y, a continuación, pulse la tecla f2 (accept). Seleccione otro punto de calibración para editar o pulse la tecla f1 (back).
 - c. Si se selecciona Delete, el punto de calibración se eliminará. Si la eliminación del punto lo cambia por una calibración de un punto, introduzca la pendiente con la ventana emergente de introducción de números.
- El medidor mostrará un resumen de la calibración actualizado y exportará los datos al registro de calibración. Pulse la tecla **f1 (meas)** para continuar con el modo de medida o pulse la tecla **f2 (print)** para exportar los datos de calibración a una impresora u ordenador.

Procedimiento de medida

Prepare los electrodos de acuerdo con las instrucciones de los manuales de electrodos. Conecte todos los electrodos que se deben utilizar a las entradas del medidor adecuadas. Asegúrese de que los electrodos se hayan calibrado recientemente y que funcionen correctamente.

Asegúrese de que el modo de medida del medidor se haya establecido en el parámetro de medida deseado. Para los medidores Orion Star A215 y Orion Star A216, pulse la tecla **f3 (channel)** hasta que se muestre el canal o canales deseados en el modo de medida.

- Enjuague los electrodos con agua destilada o una solución adecuada, séquelos con suavidad con un paño sin pelusas para quitar el exceso de agua y colóquelos en la muestra.
- 2. Inicie la medida y espere a que se realice la lectura para estabilizar o alcance el tiempo predefinido.
 - Auto-Read: Pulse la tecla measure (esc) para iniciar la medida. Si utiliza una sonda de agitador, esta empezará el agitado cuando se pulse la tecla measure (esc) y detendrá el agitado cuando la medida se estabilice.
 - b. Continuous: La medida empezará inmediatamente cuando se encuentre en el modo de medida. Si utiliza una sonda de agitador, pulse la tecla stirrer para iniciar y detener el agitado.
 - c. Timed: La medida empezará inmediatamente cuando se encuentre en el modo de medida. Si utiliza una sonda de agitador, pulse la tecla stirrer para iniciar y detener el agitado.
- 3. Una vez la medida sea estable o alcance el tiempo definido, registre todos los parámetros aplicables.
 - a. Auto-Read: Una vez la medida sea estable, se bloqueará y se mantendrá en la pantalla y el icono AR permanecerá sólido. Si la función de registro de datos está activada, la medida se exportará al registro de datos.
 - b. Continuous: El icono intermitente Stabilizing se actualizará al icono fijo Ready cuando la medida esté estable. Si la función de registro de datos está activada, pulse la tecla log/print para exportar la medida al registro de datos.
 - c. Timed: Las medidas se registrarán en el intervalo de tiempo predefinido. Si la función de registro de datos está activada, se mostrará el icono 🔚 cuando haya transcurrido cada intervalo de tiempo y la medida se exporta al registro de datos.
- Quite los electrodos de la muestra, enjuáguelos con agua destilada o una solución adecuada, séquelos con suavidad con un paño sin pelusas para quitar el exceso de agua y colóquelos en la muestra siguiente.
- Repita los pasos del 2 al 4 en todas las muestras. Cuando se hayan medido todas las muestras, guarde los electrodos de acuerdo con las instrucciones de los manuales de electrodos.



CAPÍTULO 5 El uso del canal de conductividad

Utilice el medidor de conductividad Orion Star A212 o el medidor de pH/conductividad Orion Star A215 para medir la conductividad, la salinidad, los SDT, la resistividad y la temperatura. El capítulo siguiente ofrece instrucciones sobre la conexión de los sensores, la calibración del sistema y la medida de estos parámetros.

Preparación del medidor y el sensor

- Familiarícese con el funcionamiento general del medidor y prepare el medidor y los accesorios de conformidad con las instrucciones del <u>Capítulo 2: Principios básicos del</u> <u>medidor</u>.
 - a. Se recomienda colocar los sensores en el soporte portaelectrodos conectado al medidor para un fácil movimiento dentro y fuera de los contenedores durante la calibración, la medida y el almacenamiento.
- 2. Consulte las instrucciones del <u>Capítulo 3 Menús de configuración del medidor</u> para personalizar los ajustes de medida, calibración y alarma dentro del menú de configuración del COND Channel y los ajustes de transferencia de datos, registro de datos, fecha y hora, idioma, sonidos, velocidad rotatoria, contraste de la pantalla, apagado automático, ID de usuario e ID de muestra dentro del menú de configuración de ajustes del instrumento.
- Conecte todos los electrodos, sondas y sensores necesarios a las entradas del medidor adecuadas.
 - a. Conecte un sensor de conductividad a la entrada del medidor MiniDIN de 8 pines. Si el sensor de conductividad tiene temperatura integrada, la temperatura también se medirá cuando el sensor de conductividad esté conectado al medidor.
 - b. Conecte la sonda de agitador (número de catálogo 096019) a la entrada del medidor del agitador.
- 4. Prepare todos los sensores conectados para el uso tal como se indica en los manuales de los sensores.

Procedimiento de calibración de conductividad

Los medidores de conductividad Orion Star A212 y los medidores de pH/conductividad Orion Star A215 pueden llevar a cabo una calibración de conductividad utilizando de uno a cinco estándares de conductividad con facilidad siguiendo las instrucciones que aparecen en la pantalla para ayudar con el proceso de calibración o mediante la introducción manual del valor de la constante de la celda de conductividad certificada (K).

Utilice siempre los nuevos estándares y seleccione los estándares cercanos a la conductividad de la muestra prevista. Prepare el sensor de conductividad de acuerdo con las instrucciones del manual de sensores. Conecte todos los sensores que se deben utilizar a las entradas del medidor adecuadas.

Asegúrese de que el modo de medida del medidor se haya establecido en conductividad (Cond). Para los medidores Star A215, pulse la tecla **f3 (channel)** hasta que se muestre el canal de conductividad en el modo de medida.

Nota: Para realizar una calibración automática, introduzca la constante de la celda nominal del sensor de conductividad en el menú de configuración específico del canal y utilice los estándares Orion 100 μS, 1413 yS o 12,9 mS.

- 1. Pulse la tecla **f1 (cal)** para iniciar la calibración.
 - a. Si se muestra más de un canal en el modo de medida, pulse la tecla ▲ o ▼ para resaltar el **Conductivity-Channel** y pulse la tecla **f2 (select)**.
- Enjuague el sensor de conductividad y los demás electrodos en uso con agua destilada, séquelos con suavidad con un paño sin pelusas para quitar el exceso de agua y colóquelos en el estándar de conductividad.
- 3. Cuando el sensor y el estándar estén listos, pulse la tecla f3 (start).
 - a. Si utiliza una sonda de agitador, esta empezará el agitado cuando se pulse la tecla **f3 (start)** y detendrá el agitado cuando la lectura se estabilice.
- 4. Espere hasta que el valor del medidor se estabilice y lleve a cabo alguna de las acciones siguientes:
 - Pulse la tecla f2 (accept) para aceptar el valor de conductividad determinado por la función de reconocimiento estándar automático (solo los estándares Orion 100 μS, 1413 μS y 12,9 mS), o
 - b. Pulse la tecla f3 (edit) para acceder a la ventana emergente de introducción de números e introduzca manualmente el valor del estándar de conductividad previsto en la temperatura medida.
 - i. Pulse la tecla ▲, ▼, ◀ o ► para resaltar un número o punto decimal, pulse la tecla f3 (enter) para seleccionar el elemento resaltado y repítalo hasta que el valor deseado se muestre en la parte superior de la pantalla.
 - ii. Pulse la tecla **f2 (done)** para guardar el valor y salir de la ventana de introducción de números.
 - iii. Pulse la tecla **f2 (accept)** para aceptar el valor introducido manualmente.

- 5. Pulse la tecla f2 (next) para continuar con el estándar siguiente y repita los pasos del 2 al 4 o pulse la tecla f3 (cal done) para guardar y finalizar la calibración. Si se utilizan los cinco estándares, la calibración se guardará y finalizará una vez se haya aceptado el quinto valor.
- El medidor mostrará el resumen de la calibración y exportará los datos al registro de calibración. Pulse la tecla f1 (meas) para continuar con el modo de medida o pulse la tecla f2 (print) para exportar los datos de calibración a una impresora u ordenador.

Edición de calibración de conductividad

Cuando se muestra el resumen de calibración, se puede seleccionar la opción de edición de calibración para reparar los puntos individuales sin tener que realizar una calibración completa.

- 1. En la pantalla de resumen de calibración (después del paso 5 del Procedimiento de calibración de conductividad), pulse la tecla **f3 (cal edit)**.
- Pulse la tecla ▲ o ▼ para resaltar el punto de calibración que hay que editar y pulse la tecla f2 (select).
- 3. Pulse la tecla ▲ o ▼ para resaltar <u>Remeasure</u>, <u>Edit</u> o <u>Delete</u> y pulse la tecla **f2** (select).
 - a. Si se ha seleccionado Remeasure, repita los pasos del 2 al 4 en el Procedimiento de calibración.
 - b. Si se ha seleccionado Edit, pulse la tecla f3 (edit), utilice la ventana emergente de introducción de números para introducir un nuevo valor estándar de conductividad, pulse la tecla f2 (done) y, a continuación, pulse la tecla f2 (accept). Seleccione otro punto de calibración para editar o pulse la tecla f1 (back).
 - c. Si se selecciona Delete, el punto de calibración se eliminará.
- 4. El medidor mostrará un resumen de la calibración actualizado y exportará los datos al registro de calibración. Pulse la tecla f1 (meas) para continuar con el modo de medida o pulse la tecla f2 (print) para exportar los datos de calibración a una impresora u ordenador.

Procedimiento de calibración de introducción de constante de celda certificada alternativa

Al utilizar el procedimiento de calibración de introducción de constante de celda, introduzca el valor de constante de celda certificada (K) que está impreso en el cable del sensor de conductividad o en el certificado de calibración.

- 1. Pulse la tecla f1 (cal) para iniciar la calibración.
 - a. Si se muestra más de un canal en el modo de medida, pulse la tecla ▲ o ▼ para resaltar el <u>Conductivity-Channel</u> y pulse la tecla <u>f2 (select)</u>.
- 2. Pulse la tecla f2 (cell K).
- Espere a que el valor de conductividad del medidor se estabilice y, a continuación, pulse la tecla f3 (edit) para acceder a la ventana emergente de introducción de números e introduzca el valor de constante de celda certificada.
 - a. Pulse la tecla ▲, ▼, ◀ o ► para resaltar un número o punto decimal, pulse la tecla f3 (enter) para seleccionar el elemento resaltado y repítalo hasta que el valor deseado se muestre en la parte superior de la pantalla.
 - b. Pulse la tecla **f2 (done)** para guardar el valor y salir de la ventana de introducción de números.
 - c. Pulse la tecla f2 (accept) para aceptar el valor de constante de celda introducido.
- Pulse la tecla f2 (accept) para guardar y finalizar la calibración y exportará los datos al registro de calibración. El medidor continuará automáticamente con el modo de medida.
- 5. El sistema de medida de conductividad ahora está calibrado. Compruebe la calibración leyendo un estándar de conductividad conocido o de conformidad con su SOP.

Estándares de conductividad frente a Tabla de temperatura

Temperatura	Estándar Orion	Estándar Orion	Estándar Orion	Estándar Orion	Estándar Orion
(°C)	111,9 mS (mS/cm)	12,9 mS (mS/cm)	1413 pS (µS/cm)	147 pS (µS/cm)	100 pS (µS/cm)
0	65,10	7,135	//6	81	54
1	66,84	7,344	799	83	56
2	68,59	7,555	822	86	58
3	70,35	7,768	846	88	59
4	72,12	7,983	870	91	61
5	73,91	8,200	894	93	63
6	/5,/0	8,418	918	96	64
1	77,50	8,638	943	98	66
8	79,32	8,860	968	101	68
9	81,15	9,084	992	103	70
10	82,98	9,309	1017	106	72
11	84,83	9,535	1043	108	73
12	86,69	9,763	1068	111	/5
13	88,56	9,993	1094	114	//
14	90,45	10,22	1119	116	79
15	92,34	10,46	1145	119	81
16	94,24	10,69	1171	122	83
17	96,15	10,93	1198	125	85
18	98,08	11,16	1224	127	87
19	100,0	11,40	1251	130	88
20	102,0	11,64	1277	133	90
21	103,9	11,88	1304	136	92
22	105,9	12,12	1331	138	94
23	107,9	12,36	1358	141	96
24	109,9	12,61	1386	144	98
25	111,9	12,85	1413	147	100
26	113,9	13,10	1441	150	102
27	115,9	13,35	1468	153	104
28	117,9	13,59	1496	156	106
29	120,0	13,84	1524	159	108
30	122,0	14,09	1552	161	110
31	124,1	14,34	1580	164	112
32	126,2	14,59	1608	167	114
33	128,3	14,85	1636	170	117
34	130,4	15,10	1665	173	119
35	132,5	15,35	1693	176	121
36	134,6	15,61	1722	179	123
37	136,7	15,86	1751	182	125
38	138,9	16,12	1780	185	127
39	141,0	16,37	1808	188	129
40	143,2	16,63	1837	191	131
41	145,4	16,89	1866	194	134
42	147,6	17,15	1896	197	136
43	149,8	17,40	1925	200	138
44	152,0	17,66	1954	203	140
45	154,2	17,92	1983	206	142
46	156,4	18,18	2013	209	145
47	158,7	18,44	2042	212	147
48	160,9	18,70	2071	215	149
49	163,2	18,96	2101	219	151
50	165,4	19,22	2130	222	154

Procedimiento de verificación de medidor de conductividad

Utilice el kit de verificación del medidor de conductividad de la serie Orion Star, número de catálogo 1010001, para verificar la precisión de las medidas de conductividad en los medidores de conductividad Orion Star A212 y los medidores de pH/conductividad Orion Star A215. El valor de conductividad medido de cada reostato debería incluirse en la precisión relativa del reostato ($\pm 0,1$ % de la conductancia del reostato real) más la precisión relativa del medidor ($\pm 0,5$ % de lectura ± 1 dígito para lecturas mayores a 3 μ S/cm o $\pm 0,5$ % de lectura ± 0.01 jS/cm para lecturas menores o iguales a 3 μ S/cm).

- 1. Asegúrese de que el kit de reostato de conductividad esté certificado y tenga una fecha de calibración válida.
- 2. En el modo de medida, pulse la tecla f1 (cal).
 - a. Si se muestra más de un canal en el modo de medida, pulse la tecla ▲ o ▼ para resaltar el <u>Conductivity-Channel</u> y pulse la tecla **f2 (select)**.
- 3. Pulse la tecla f2 (cell K).
- 4. Pulse la tecla **f3 (edit)** para acceder a la ventana emergente de introducción de números e introduzca un valor de constante de celda de 1,0000.
 - a. Pulse la tecla ▲, ▼, ◀ o ▶ para resaltar un número o punto decimal, pulse la tecla f3 (enter) para seleccionar el elemento resaltado y repítalo hasta que el valor deseado se muestre en la parte superior de la pantalla.
 - b. Pulse la tecla **f2 (done)** para guardar el valor y salir de la ventana de introducción de números.
 - c. Pulse la tecla f2 (accept) para aceptar el valor de constante de celda introducido.
- 5. El medidor continuará automáticamente con el modo de medida y la <u>Constante de</u> <u>celda:</u> 1,0000 aparecerá en la pantalla.
- Conecte un reostato al medidor, pulse la tecla measure (esc) para iniciar la medida, espere a que el valor de medida se estabilice y, a continuación, registre el valor de conductividad mostrado del reostato.
- 7. Repita el paso 6 en todos los seis reostatos del kit.

Número de reostato	Resistencia nominal	Conductancia nominal	Intervalo de aceptación nominal
1010001-A	1000 ΚΩ	1 μS	de 0,984 a 1,016 μS
1010001-B	100 KΩ	10 µS	de 9,930 a 10,07 μS
1010001-C	10 KΩ	100 μS	de 99,30 a 100,7 μS
1010001-D	1 KΩ	1000 μS	de 993,0 a 1007 µS
1010001-E	100 Ω	10 mS	de 9,930 a 10,07 mS
1010001-F	10 Ω	100 mS	de 99,30 a 100,7 mS

Nota: Estos son los valores nominales únicamente de los reostatos. La resistencia y la conductancia reales se imprimen en cada reostato y el intervalo de aceptación real debe calcularse a partir del valor real en cada reostato.

Procedimiento de medida

Prepare los sensores de acuerdo con las instrucciones del manual de sensores. Conecte todos los electrodos, sensores y sondas que se deben utilizar a las entradas del medidor adecuadas. Asegúrese de que los sensores se hayan calibrado recientemente y que funcionen correctamente.

Asegúrese de que el modo de medida del medidor se haya establecido en el parámetro de medida deseado. Para los medidores Star A215, pulse la tecla **f3 (channel)** hasta que se muestre el canal o canales deseados en el modo de medida.

- Enjuague los sensores con agua destilada o una solución adecuada, séquelos con suavidad con un paño sin pelusas para quitar el exceso de agua y colóquelos en la muestra.
- 2. Inicie la medida y espere a que se realice la lectura para estabilizar o alcance el tiempo predefinido.
 - Auto-Read: Pulse la tecla measure (esc) para iniciar la medida. Si utiliza una sonda de agitador, esta empezará el agitado cuando se pulse la tecla measure (esc) y detendrá el agitado cuando la medida se estabilice.
 - b. Continuous: La medida empezará inmediatamente cuando se encuentre en el modo de medida. Si utiliza una sonda de agitador, pulse la tecla **stirrer** para iniciar y detener el agitado.
 - c. Timed: La medida empezará inmediatamente cuando se encuentre en el modo de medida. Si utiliza una sonda de agitador, pulse la tecla stirrer para iniciar y detener el agitado.
- 3. Una vez la medida sea estable o alcance el tiempo definido, registre todos los parámetros aplicables.
 - a. Auto-Read: Una vez la medida sea estable, se bloqueará y se mantendrá en la pantalla y el icono AR permanecerá sólido. Si la función de registro de datos está activada, la medida se exportará al registro de datos.
 - b. Continuous: El icono intermitente Stabilizing se actualizará al icono fijo Ready cuando la medida esté estable. Si la función de registro de datos está activada, pulse la tecla log/print para exportar la medida al registro de datos.
 - c. Timed: Las medidas se registrarán en el intervalo de tiempo predefinido. Si la función de registro de datos está activada, se mostrará el icono 🔚 cuando haya transcurrido cada intervalo de tiempo y la medida se exporta al registro de datos.
- Quite los sensores de la muestra, enjuáguelos con agua destilada o una solución adecuada, séquelos con suavidad con un paño sin pelusas para quitar el exceso de agua y colóquelos en la muestra siguiente.
- Repita los pasos del 2 al 4 en todas las muestras. Cuando se hayan medido todas las muestras, guarde los sensores de acuerdo con las instrucciones de los manuales de sensores.



CAPÍTULO 6 El uso del canal OD/RDO

Utilice el medidor de RDO/OD Orion Star A213 o el medidor de pH/RDO/OD Orion Star A216 para medir el oxígeno disuelto como saturación de porcentaje o mg/L y la temperatura mediante los sensores de oxígeno disuelto polarográfico u óptico Orion RDO. El capítulo siguiente ofrece instrucciones sobre la conexión de los sensores, la calibración del sistema y la medida de estos parámetros.

Preparación del medidor y el sensor

- Familiarícese con el funcionamiento general del medidor y prepare el medidor y los accesorios de conformidad con las instrucciones del <u>Capítulo 2: Principios básicos del</u> medidor.
 - Se recomienda colocar los sensores en el soporte portaelectrodos conectado al medidor para un fácil movimiento dentro y fuera de los contenedores durante la calibración, la medida y el almacenamiento.
- 2. Consulte las instrucciones del <u>Capítulo 3 Menús de configuración del medidor</u> para personalizar los ajustes de medida, calibración y alarma dentro del menú de configuración del canal de OD/RDO y los ajustes de transferencia de datos, registro de datos, fecha y hora, idioma, sonidos, velocidad rotatoria, contraste de la pantalla, apagado automático, ID de usuario e ID de muestra dentro del menú de configuración de ajustes del instrumento.
- Conecte todos los electrodos, sondas y sensores necesarios a las entradas del medidor adecuadas.
 - a. Conecte un sensor de oxígeno disuelto polarográfico u óptico RDO a la entrada del medidor MiniDIN de 9 pines. El medidor detectará automáticamente cuál es el sensor de oxígeno disuelto que está conectado. La temperatura también se medirá cuando el sensor de oxígeno disuelto esté conectado al medidor.
 - b. Conecte la sonda de agitador (número de catálogo 096019) a la entrada del medidor del agitador.
- 4. Prepare todos los sensores conectados para el uso tal como se indica en los manuales de los sensores.

Procedimiento de calibración de oxígeno disuelto

Nota: Un sensor polarográfico de oxígeno disuelto se debe polarizar antes de usarse. El sensor polarográfico de oxígeno disuelto se polariza de forma continuada al conectarse con el medidor. Si el sensor de oxígeno disuelto polarográfico es nuevo, se ha reparado o no se ha conectado al medidor, conecte el sensor al medidor, encienda el medidor y espere 30 minutos hasta que el sensor se polarice.

Los medidores de RDO/OD Orion Star A213 y los medidores de pH/RDO/OD Orion Star A216 pueden realizar las calibraciones siguientes:

- Air (Aire) (aire saturado con agua): es el método más sencillo y preciso y utiliza la manga de calibración incluida con los sensores de oxígeno más disuelto. Para obtener la mayor precisión, la temperatura de calibración debe coincidir con la temperatura prevista de la muestra. Humedezca la esponja de la manga de calibración con agua destilada e introduzca el sensor en la manga. De modo alternativo, se puede utilizar una botella de DBO con únicamente agua destilada suficiente para cubrir la parte inferior sin tocar el sensor.
- Water (Agua) (agua saturada con aire): este método utiliza agua que está 100 % saturada con aire. Introduzca aire en una muestra de agua durante un período prolongado, preferiblemente durante la noche.
- Manual (Winkler): este método utiliza una muestra de agua con una concentración conocida de oxígeno disuelto y se utiliza normalmente para calibrar el sensor de OD en la válvula a través de un método de titulación Winkler. Debido a posibles errores de titulación, este método es, por naturaleza, menos preciso.
- Set Zero (Definición cero): este método utiliza una solución sin oxígeno para añadir un punto cero a una calibración de aire saturado con agua o de agua saturada con aire existente. Normalmente se recomienda esta calibración cuando las medidas están por debajo de la saturación del 10 % o de 1 mg/L.

Prepare el sensor de oxígeno disuelto de acuerdo con las instrucciones del manual de sensores. Conecte todos los sensores que se deben utilizar a las entradas del medidor adecuadas.

Asegúrese de que el modo de medida del medidor se haya establecido en oxígeno disuelto (OD o RDO). Para los medidores Star A216, pulse la tecla **f3 (channel)** hasta que se muestre el canal de oxígeno disuelto en el modo de medida.

Calibración de aire (aire saturado con agua)

- 1. Pulse la tecla **f1 (cal)** para iniciar la calibración.
 - a. Si se muestra más de un canal en el modo de medida, pulse la tecla ▲ o ▼ para resaltar el <u>DO/RDO-Channel</u> y pulse la tecla f2 (select).
- 2. Pulse la tecla ▲ o ▼ para resaltar <u>Air</u> (Aire) y pulse la tecla **f3 (select)**.
- Prepare el sensor de OD y la manga de calibración y permita que alcancen su punto de equilibrio.
- Cuando el sensor de OD y la manga de calibración estén listos, pulse la tecla f3 (start).
- 5. Espere hasta que el valor de oxígeno disuelto en el medidor se estabilice.
 - Con un sensor de OD polarográfico, se mostrará el 102,3 % cuando la lectura se estabilice.
 - b. Con un sensor de OD óptico RDO, se mostrará el 100,0 % cuando la lectura se estabilice.
- 6. Pulse la tecla f2 (cal done) para guardar y finalizar la calibración.
- El medidor mostrará el resumen de la calibración y exportará los datos al registro de calibración. Pulse la tecla f1 (meas) para continuar con el modo de medida o pulse la tecla f2 (print) para exportar los datos de calibración a una impresora u ordenador.

Calibración de agua (agua saturada con aire)

- 8. Pulse la tecla f1 (cal) para iniciar la calibración.
 - a. Si se muestra más de un canal en el modo de medida, pulse la tecla ▲ o ▼ para resaltar el <u>DO/RDO-Channel</u> y pulse la tecla f2 (select).
- Pulse la tecla ▲ o ▼ para resaltar <u>Water</u> (Agua) y pulse la tecla f3 (select).
- 10. Prepare el sensor de OD y el agua saturada con aire al 100 % y permita que alcancen su punto de equilibrio.
 - a. Utilice una película de parafina de plástico para sellar la zona abierta entre el sensor de OD y el recipiente que contiene el agua saturada con aire al 100 %.
- Cuando el sensor de OD y el equipo de calibración estén listos, pulse la tecla f3 (start).
 - a. Si utiliza una sonda de agitador, esta empezará el agitado cuando se pulse la tecla **f3 (start)** y detendrá el agitado cuando la lectura se estabilice.
- 12. Espere hasta que el valor de oxígeno disuelto en el medidor se estabilice.
 - a. Se mostrará 100,0 % cuando la lectura se estabilice.
- 13. Pulse la tecla f2 (cal done) para guardar y finalizar la calibración.
- 14. El medidor mostrará el resumen de la calibración y exportará los datos al registro de calibración. Pulse la tecla f1 (meas) para continuar con el modo de medida o pulse la tecla f2 (print) para exportar los datos de calibración a una impresora u ordenador.

Calibración manual (Winkler)

- 1. Pulse la tecla f1 (cal) para iniciar la calibración.
 - a) Si se muestra más de un canal en el modo de medida, pulse la tecla ▲ o ▼ para resaltar el <u>DO/RDO-Channel</u> y pulse la tecla **f2 (select)**.
- 2. Pulse la tecla ▲ o ▼ para resaltar Manual y pulse la tecla f3 (select).
- Prepare el sensor de OD y la solución de calibración y permita que alcancen su punto de equilibrio.
 - a. Utilice una película de parafina de plástico para sellar la zona abierta entre el sensor de OD y el recipiente que contiene la solución de calibración.
- Cuando el sensor de OD y el equipo de calibración estén listos, pulse la tecla f3 (start).
 - a. Si utiliza una sonda de agitador, esta empezará el agitado cuando se pulse la tecla **f3 (start)** y detendrá el agitado cuando la lectura se estabilice.
- 5. Espere hasta que el valor de oxígeno disuelto en el medidor se estabilice y lleve a cabo alguna de las acciones siguientes:
 - a. Pulse la tecla **f2 (accept)** para aceptar el valor de oxígeno disuelto mostrado.

o bien

- b. Pulse la tecla **f3 (edit)** para acceder a la ventana emergente de introducción de números e introduzca manualmente el valor de oxígeno disuelto.
 - i. Pulse la tecla ▲, ▼, ◀ o ▶ para resaltar un número o punto decimal, pulse la tecla f3 (enter) para seleccionar el elemento resaltado y repítalo hasta que el valor deseado se muestre en la parte superior de la pantalla.
 - ii. Pulse la tecla **f2 (done)** para guardar el valor y salir de la ventana de introducción de números.
- 6. Pulse la tecla **f2 (cal done)** para guardar y finalizar la calibración.
- El medidor mostrará el resumen de la calibración y exportará los datos al registro de calibración. Pulse la tecla f1 (meas) para continuar con el modo de medida o pulse la tecla f2 (print) para exportar los datos de calibración a una impresora u ordenador.
Calibración de definición cero

Debe realizarse una calibración de aire (aire saturado con agua) o de agua (agua saturada con aire) antes de que se lleve a cabo una calibración de definición cero.

Prepare una solución de sulfito de sodio disolviendo aproximadamente 15,0 gramos de Na₂SO₃ en unos 250 mL de agua destilada. Transfiera la solución a un frasco o una botella de DBO y utilice una película de parafina de plástico para sellar la botella. Se puede añadir una pequeña cantidad de sal de cobalto a la solución de sulfito de sodio para que actúe como indicador y cambie de color cuando la solución de sulfito de sodio ya no tenga cero contenido de oxígeno.

- 1. Pulse la tecla f1 (cal) para iniciar la calibración.
 - a. Si se muestra más de un canal en el modo de medida, pulse la tecla ▲ o ▼ para resaltar el <u>DO/RDO-Channel</u> y pulse la tecla **f2 (select)**.
- Pulse la tecla ▲ o ▼ para resaltar <u>Set Zero</u> (Definición cero) y pulse la tecla f3 (select).
- 3. Prepare el sensor de OD y la solución de cero oxígeno y permita que alcancen su punto de equilibrio.
 - a. Utilice una película de parafina de plástico para sellar la zona abierta entre el sensor de OD y el recipiente que contiene la solución de calibración.
- Cuando el sensor de OD y el equipo de calibración estén listos, pulse la tecla f3 (start).
 - a. Si utiliza una sonda de agitador, esta empezará el agitado cuando se pulse la tecla **f3 (start)** y detendrá el agitado cuando la lectura se estabilice.
- 5. Espere hasta que el valor de oxígeno disuelto en el medidor se estabilice.
 - a. Se mostrará 0,0 % para P2 cuando la lectura se estabilice.
- 6. Pulse la tecla **f2 (cal done)** para guardar y finalizar la calibración.
- El medidor mostrará el resumen de la calibración y exportará los datos al registro de calibración. Pulse la tecla f1 (meas) para continuar con el modo de medida o pulse la tecla f2 (print) para exportar los datos de calibración a una impresora u ordenador.

Procedimiento de medida

Prepare los sensores de acuerdo con las instrucciones del manual de sensores. Conecte todos los electrodos, sensores y sondas que se deben utilizar a las entradas del medidor adecuadas. Asegúrese de que los sensores se hayan calibrado recientemente y que funcionen correctamente.

Asegúrese de que el modo de medida del medidor se haya establecido en el parámetro de medida deseado. Para los medidores Star A216, pulse la tecla **f3 (channel)** hasta que se muestre el canal o canales deseados en el modo de medida.

- Enjuague los sensores con agua destilada o una solución adecuada, séquelos con suavidad con un paño sin pelusas para quitar el exceso de agua y colóquelos en la muestra.
- 2. Inicie la medida y espere a que se realice la lectura para estabilizar o alcance el tiempo predefinido.
 - Auto-Read: Pulse la tecla measure (esc) para iniciar la medida. Si utiliza una sonda de agitador, esta empezará el agitado cuando se pulse la tecla measure (esc) y detendrá el agitado cuando la medida se estabilice.
 - b. Continuous: La medida empezará inmediatamente cuando se encuentre en el modo de medida. Si utiliza una sonda de agitador, pulse la tecla stirrer para iniciar y detener el agitado.
 - c. Timed: La medida empezará inmediatamente cuando se encuentre en el modo de medida. Si utiliza una sonda de agitador, pulse la tecla stirrer para iniciar y detener el agitado.
- 3. Una vez la medida sea estable o alcance el tiempo definido, registre todos los parámetros aplicables.
 - a. Auto-Read: Una vez la medida sea estable, se bloqueará y se mantendrá en la pantalla y el icono AR permanecerá sólido. Si la función de registro de datos está activada, la medida se exportará al registro de datos.
 - b. Continuous: El icono intermitente Stabilizing se actualizará al icono fijo Ready cuando la medida esté estable. Si la función de registro de datos está activada, pulse la tecla log/print para exportar la medida al registro de datos.
 - c. Timed: Las medidas se registrarán en el intervalo de tiempo predefinido. Si la función de registro de datos está activada, se mostrará el icono 🔚 cuando haya transcurrido cada intervalo de tiempo y la medida se exporta al registro de datos.
- Quite los sensores de la muestra, enjuáguelos con agua destilada o una solución adecuada, séquelos con suavidad con un paño sin pelusas para quitar el exceso de agua y colóquelos en la muestra siguiente.
- Repita los pasos del 2 al 4 en todas las muestras. Cuando se hayan medido todas las muestras, guarde los sensores de acuerdo con las instrucciones de los manuales de sensores.



CAPÍTULO 7 Transferencia de datos y actualizaciones de software

Almacenamiento de datos y ajustes de transferencia

Los ajustes del medidor que afectan al almacenamiento y a la transferencia de datos incluyen el tipo de lectura de la medida, los ajustes de los datos de exportación y el ajuste del registro de datos. Los ajustes de fecha y hora deben revisarse y actualizarse según sea necesario.

Ajuste del tipo de lectura de la medida

El tipo de lectura seleccionada determina cuándo el medidor envía las mediciones al registro de datos, si el registro de datos está activado en el menú de configuración, y cuándo el medidor exporta las medidas a una impresora o un ordenador, si hay una impresora o un ordenador correctamente conectados al medidor y los ajustes de datos de exportación están activados en el menú de configuración. Seleccione un tipo de lectura: Auto-Read, Continuous o Timed.

Auto-Read

Pulse la tecla measure (esc) para iniciar la medida. El icono **AR** parpadea a medida que el valor cambiante se estabiliza. Cuando la medida es estable, el icono **AR** deja de parpadear y la medida queda fija en la pantalla hasta que se pulsa de nuevo la tecla **measure (esc)**. Cuando los ajustes del registro de datos y de datos de exportación están activados, la medida estable se guarda automáticamente en el registro de datos y se exporta a una impresora o un ordenador.

Continuous

Los valores de medida se actualizan constantemente en la pantalla y el icono **Stabilizing** o **Ready** indica el estado de estabilidad de la medida. Cuando los ajustes del registro de datos y de datos de exportación estén activados, pulse la tecla **log/print** para guardar una medida en el registro de datos y exportarla a una impresora o un ordenador.

Timed

Los valores de medida se actualizan constantemente en la pantalla y el icono **Stabilizing** o **Ready** indica el estado de estabilidad de la medida. Cuando los ajustes del registro de datos y de datos de exportación están activados, las medidas se guardan en el registro de datos y se exportan a una impresora o un ordenador en los intervalos de tiempo seleccionados a partir de 3 segundos (00:00:03) hasta 24 horas (24:00:00).

Cómo ajustar el tipo de lectura de la medida:

- 1. En el modo de medida, pulse la tecla setup para acceder al menú de configuración principal.
- Pulse la tecla ▲, ▼, ◀ o ▶ para resaltar <u>pH Channel</u>, <u>pH/ISE Channel</u>, <u>COND Channel</u> o <u>DO/RDO Channel</u> y pulse la tecla **f3 (select)**.
- 3. Pulse la tecla \blacktriangle o \triangledown para resaltar <u>Mode and Settings</u> y pulse la tecla **f3 (select)**.
- 4. Pulse la tecla ▲ o ▼ para resaltar <u>Read Type</u> y pulse la tecla **f3 (select)**.
- Pulse la tecla ▲ o ▼ para resaltar <u>Auto</u>, <u>Continuous</u> o <u>Timed</u> y pulse la tecla
 f3 (select) para definir el modo de medida deseado.
 - a. Si se ha seleccionado Timed, pulse la tecla ◀ o ► para resaltar las horas (00:00:00), los minutos (00:00:00) o los segundos (00:00:00).
 - b. Pulse la tecla **f3 (edit)** para acceder a la ventana emergente de introducción de números.
 - c. Pulse la tecla ▲, ▼, ◀ o ▶ para resaltar un número; pulse la tecla f3 (enter) para seleccionar el número y repítalo hasta que un valor de dos dígitos se muestre en la parte superior de la ventana.
 - d. Pulse la tecla **f2 (done)** para guardar el valor y salir de la ventana de introducción de números.
 - e. Repítalo hasta que se introduzca el intervalo de tiempo correcto para las horas, los minutos y los segundos. Se pueden introducir intervalos a partir de 3 segundos (00:00:03) hasta 24 horas (24:00:00).
- 6. Pulse la tecla **f1 (back)** para navegar fuera del menú y, a continuación, pulse la tecla **measure (esc)** para volver al modo de medida.

Ajustes del registro de datos

Para guardar hasta 2000 conjuntos de medidas en el registro de datos del medidor, active el ajuste de registro de datos en el menú de configuración.

- En el modo de medida, pulse la tecla setup para acceder al menú de configuración principal.
- Pulse la tecla ▲, ▼, ◀ o ▶ para resaltar <u>Settings</u> (Ajustes) y pulse la tecla f3 (select).
- 3. Pulse la tecla \blacktriangle o \triangledown para resaltar <u>Data Log</u> y pulse la tecla **f3 (select)**.
- Pulse la tecla ▲ o ▼ para resaltar <u>On</u> y pulse la tecla f3 (select).
- Pulse la tecla f1 (back) para navegar fuera del menú y pulse la tecla measure (esc) para volver al modo de medida.

Ajustes de exportación de datos

Para transferir las medidas a un ordenador o una impresora, active la función de exportación de datos en el menú de configuración y actualice los ajustes de exportación para que coincidan con los ajustes del dispositivos externo, incluida la velocidad de baudios. La conexión de comunicación del medidor predeterminada es RS232 y el medidor detectará y ajustará automáticamente el ajuste de conexión cuando un cable USB activo esté conectado al medidor.

- En el modo de medida, pulse la tecla setup para acceder al menú de configuración principal.
- 2. Pulse la tecla \blacktriangle , \triangledown , \triangleleft o \triangleright para resaltar <u>Settings</u> y pulse la tecla f3 (select).
- Pulse la tecla ▲ o ▼ para resaltar Export Data (Exportar datos) y pulse la tecla f3 (select).
- Pulse la tecla ▲ o ▼ para resaltar <u>Printing</u> (Impresión) y pulse la tecla f3 (select).
 Pulse la tecla ▲ o ▼ para resaltar <u>On</u> y pulse la tecla f3 (select).
- Pulse la tecla ▲ o ▼ para resaltar <u>Data Format</u> (Formato de datos) y pulse la tecla
 f3 (select). Pulse la tecla ▲ o ▼ para resaltar <u>Printer</u> o <u>PC (CSV)</u> y pulse la tecla
 f3 (select).
- Pulse la tecla ▲ o ▼ para resaltar Comm Config (Configuración general) y pulse la tecla f3 (select). Pulse la tecla ▲ o ▼ para resaltar 1200, 2400, 4800, 9600, 19200 o 38400 como velocidad de baudios del medidor y pulse la tecla f3 (select).
- 7. Pulse la tecla **f1 (back)** para navegar fuera del menú y pulse la tecla **measure (esc)** para volver al modo de medida.

Ejemplos de transferencia de datos

Calibración de pH de 5 puntos, Formato CSV, Vista de HyperTerminal:

HT - HyperTerminal	6
Eile Edit View Call Iransfer Help	
ටළ මෙදී මෙපී ඒ	
A211 pH,X01036,3.04,CH-1,PH,01/03/15 12:52:00,5,1.2.00,pH,290.5,mV,25.0,C,2,4.01 ,pH,174,3,mV,25.0,C,3,7.00,pH,-0.0,mV,25.0,C,4.10.01,pH,-173.9,mV,25.0,C,5,12.00 ,pH,-289.7,mV,25.0,C,97.7,%,98.5,%,97.7,%,98.4,%,1.4,mV,-0.0,mV,-0.0,mV,1.3,mV.9 8.1,%,#1	· Trum
	+
Connected 0:00:12 Auto detect 9600 8-N-1 SCHOLL CAPS NUM Capiture Print echo	υĭ

Medida de pH, Formato CSV, Vista de HyperTerminal:

Li HT - HyperTorminal	×
Eile Edit View Call Transfer Help	
2 😹 📾 🖇 🗠 🗑 📽	
R211 pH.X01036.3.04.ABCDE.01/03/15 13:03:07.SAMPLE.CH-1.pH.7.000.pH.0.0. wV.25. .C.98.1.%.M100.#9 -	8
Connected 0:00:21 Auto detect 9600 8-N-1 SCROLL CAPS NUM Capitale Print echo	

Calibración de pH de 5 puntos, Formato de impresora:

Thermo Scientific (c) :	2011
A211 pH	100000
Meter S/N	X01036
SW Ret	3.04
Calibration Report	
PH .	
12/27/14 16:43:00	
Point 1	
pM	2.00 pH
mV	290.5 mV
Temperature	25.0 C
Point 2	
рH	4.01 pH
ΨV	174.3 mV
Temperature	25.0 C
Point 3	
5)H	7.00 pH
mv	0.0 mV
Temperature	25.0 C
Point 4	
ъH	10.01 cH
mV	-173.6 mV
Temperature	25.0 C
Point 5	
ъM	12.00 pH
m'V	-289.4 mV
Temperature	25.0 0
Slopel	97.7 %
Slope2	98.5 %
Slope3	97.5 %
Slope4	98.4 %
RI C	1.5 mW
F 2	0.0 mV
83	0.0.00
T A	1.7
	+ + 7 MAY.
Average Slope	98.0 *
Calibration	#3
Operator	
Bignature	

Medición de pH, Formato de impresora:

Thermo Scientific (c)	2011
AZ11 pH	10000000
Meter S/N	X01036
SW Rev	3.04
User ID ABCDE	
01/03/15 12:58:34	
SampleID SAMPLE	
DH	7.000 pH
νm	0.0 mV
Temperature	25.0 C
Slope	98.1 %
Method#	MICO
Calibration	\$1
Operator	2.2
Signature	

Compatibilidad y requisitos de la impresora

Los medidores de sobremesa Orion Star serie A210 pueden imprimir directamente en la impresora de inyección de tinta de la serie Orion Star, número de catálogo 1010006. Los datos de medida y calibración enviados desde el medidor a la impresora de inyección de tinta de la serie Orion Star se formatean automáticamente para ajustar la anchura del papel cuando se ha seleccionado el formato de datos de la impresora dentro del ajuste Export Data. La impresora de inyección de tinta de la serie Orion Star tiene una velocidad de 9600 baudios e incluye el cable de impresora RS232 necesario para conectar un medidor Orion Star serie A210 a la impresora.



La velocidad de baudios de los medidores de Orion Star serie A210 se puede definir en 1200, 2400, 4800, 9600, 19 200 o 38 400 para la comunicación con distintas impresoras. Los ajustes fijos del medidor son los siguientes:

Número de bits de datos:	8
Bits de parada:	1
Paridad:	Ning.
Control de flujo:	XON/XOFF

El medidor enviará las medidas y los datos de calibración a la impresora si la opción de impresora está activada en el menú de configuración. De modo alternativo, las medidas y los datos de calibración se pueden enviar al registro de datos y al registro de calibración. Desde el registro de datos y el registro de calibración, el operador puede imprimir un punto de registro de datos simple, un intervalo de puntos de registro de datos o todo el registro de datos.

Compatibilidad y requisitos del ordenador

Los medidores de la serie Orion Star A210 pueden enviar medidas y datos de calibración a un ordenador en un formato delimitado por comas que resulta fácil de analizar en programas informáticos como Excel.

La velocidad de baudios de los medidores Orion Star serie A210 se puede definir en 1200, 2400, 4800, 9600, 19 200 o 38 400. Los ajustes fijos del medidor son los siguientes:

Número de bits de datos:	8
Bits de parada:	1
Paridad:	Ning.
Control de flujo:	XON/XOFF

El medidor enviará las medidas y los datos de calibración al ordenador si la opción de impresora está activada en el menú de configuración. De modo alternativo, las medidas y los datos de calibración se pueden enviar al registro de datos y al registro de calibración. Desde el registro de datos y el registro de calibración, el operador puede imprimir un punto de registro de datos simple, un intervalo de puntos de registro de datos o todo el registro de datos.

El ordenador debe estar actualizado con las últimas actualizaciones de Windows instaladas y tener la versión 4.0 de .NET Framework con las últimas actualizaciones.

Todos los ajustes de gestión de potencia y de protectores de pantalla deben estar desactivados cuando el medidor está conectado con el ordenador. El ordenador no puede entrar en modo de suspensión o de espera mientras el medidor está transfiriendo datos. Los ordenadores portátiles deben conectarse a una fuente de alimentación; no se recomienda que funcionen con alimentación de la batería.

Se pueden utilizar tres cables de ordenador para conectar los medidores de sobremesa Orion Star serie A210 con un ordenador: el cable de ordenador USB a serie de la serie Orion Star (N.º de cat. 1010005), cable de ordenador RS232 de la serie Orion Star (N.º de cat. 1010053) o un cable de ordenador estándar Mini B USB a USB.

Cable de ordenador	Medidor Conexión	Conexión de ordenador	Controlador necesario
Cable de ordenador USB a serie, N.º de cat. 1010005	RS232	USB	Sí
Cable de ordenador RS232, N.º de cat. 1010053	RS232	RS232	No
Cable de ordenador USB (cable estándar Mini B USB a USB)	USB (Mini B)	USB	Sí

Uso del cable de ordenador USB a serie

Encienda el medidor y conecte el cable de ordenador USB a serie, en primer lugar, al medidor y, a continuación, al ordenador. El controlador de USB a serie debe estar instalado cuando el medidor se conecte a un ordenador mediante el cable de ordenador USB a serie. El controlador de USB a serie es compatible con los sistemas operativos Windows[®] XP, Windows Vista[®], Windows 7 y Windows 8 de Microsoft[®].

- 1. Conecte el medidor a una toma de pared con el adaptador de alimentación universal (número de catálogo 1010003) o instale cuatro pilas AA y encienda el medidor.
- 2. Conecte el cable de ordenador USB a serie, en primer lugar, a la entrada RS232 del medidor y, a continuación, a la entrada USB del ordenador.
- 3. Después de conectar el cable de ordenador USB a serie al ordenador, el ordenador identificará automáticamente el dispositivo e instalará el controlador. Una vez que se haya instalado el controlador del dispositivo, el cable de ordenador USB a serie se podrá utilizar para transferir datos del medidor al ordenador mediante el software informático Orion Star Com, HyperTerminal o un programa similar y enviar instrucciones remotas desde el ordenador al medidor.



- Registre la ubicación del puerto COM del cable USB a serie, por ejemplo, COM8. Para ver la ubicación del puerto COM, abra la herramienta Administrador de dispositivos del ordenador y expanda la opción Puertos.
 - a. Para Windows 7: Haga clic en el botón Inicio y, a continuación, en Panel de control. En la ventana del Panel de control, haga clic en Administrador de dispositivos si la visualización es por iconos grandes o iconos pequeños o haga clic en Sistema y seguridad y, a continuación, en Administrador de dispositivos si la visualización es por categoría. En la ventana Administrador de dispositivos, expanda la opción Puertos.
 - b. Para Windows XP: Haga clic en el botón Inicio y, a continuación, en Panel de control. En la ventana Panel de control, haga clic en el enlace Rendimiento y mantenimiento y, a continuación, en el icono Sistema o, para la vista clásica del Panel de control, haga doble clic en el icono Sistema. En la ventana Propiedades del sistema, haga clic en la pestaña Hardware y, a continuación, en el botón Administrador de dispositivos. En la ventana Administrador de dispositivos, expanda la opción Puertos.



 El medidor ahora puede transferir datos al ordenador mediante el puerto COM virtual y el software informático Orion Star Com, HyperTerminal o un programa similar.

Nota: Si el ordenador no instala automáticamente el controlador del dispositivo, abra la herramienta Administrador de dispositivos del ordenador, expanda la opción Otros dispositivos, haga doble clic en el dispositivo e instale el controlador manualmente. Es posible que los controladores de algunos RS232 de USB a adaptadores deban descargarse del sitio web del fabricante. Por ejemplo, cuando utilice un adaptador Tripp Lite, vaya a <u>http://www.tripplite.com/support/downloads/</u>, escriba el número de modelo del adaptador (p. ej. U209- 000-R) y seleccione el controlador adecuado para el sistema operativo del ordenador.



Uso del cable de ordenador USB

El controlador de USB Orion Star A200-A300 debe estar instalado cuando el medidor se conecte a un ordenador mediante el cable de ordenador USB. El controlador de USB Orion Star A200-A300 es compatible con los sistemas operativos Windows® XP, Windows Vista®, Windows 7 y Windows 8 de Microsoft®. Cuando utilice un cable de ordenador USB, encienda siempre el medidor y conecte el cable de ordenador USB, en primer lugar, al medidor y, a continuación, al ordenador.

- Vaya a <u>www.thermoscientific.com/OrionMeters</u>, descargue la carpeta del controlador de USB Orion Star A200-A300 y descomprima/extraiga los archivos al escritorio del ordenador.
- 2. Conecte el adaptador de alimentación al medidor y encienda el medidor.
- 3. Conecte el cable USB, en primer lugar, a la entrada USB Mini B del medidor y, a continuación, a la entrada USB de un ordenador.
- 4. Una vez que se haya conectado el cable USB al ordenador, el ordenador intentará buscar el software del controlador del dispositivo y, a continuación, aparecerá un mensaje indicando que el software del controlador del dispositivo no se ha instalado correctamente. Abra la herramienta Administrador de dispositivos del ordenador.
 - a. Para Windows 7: Haga clic en el botón Inicio y, a continuación, en Panel de control. En la ventana del Panel de control, haga clic en Administrador de dispositivos si la visualización es por iconos grandes o iconos pequeños o haga clic en Sistema y seguridad y, a continuación, en Administrador de dispositivos si la visualización es por categoría. En la ventana Administrador de dispositivos, expanda la opción Puertos.
 - b. Para Windows XP: Haga clic en el botón Inicio y, a continuación, en Panel de control. En la ventana Panel de control, haga clic en el enlace Rendimiento y mantenimiento y, a continuación, en el icono Sistema o, para la vista clásica del Panel de control, haga doble clic en el icono Sistema. En la ventana Propiedades del sistema, haga clic en la pestaña Hardware y, a continuación, en el botón Administrador de dispositivos. En la ventana Administrador de dispositivos, expanda la opción Puertos.
- 5. Localice y haga doble clic en el dispositivo del puerto COM virtual (CDC) en la herramienta Administrador de dispositivos.



6. Se abrirá la ventana de propiedades del puerto COM virtual (CDC). Haga clic en el botón Actualizar controlador.

General	Driver Details	perioce .	-
1	Vintual COM Port	(CDC)	
	Device type:	Other devices	
	Manufacturer:	Unknown	
	Location	Location 0	
The The Tot	drivers for this devic re is no driver select ind a driver for this di	e are not installed. (Code 28) ad for the device information set or eleme evice. click Lincate Driver	nt
		Update Drive	t

7. Haga clic en la opción "Buscar software de controlador en el equipo".

1	Update Driver Software - Virtual COM Port (CDC)	
Hov	w do you want to search for driver software?	
•	Search automatically for updated driver software Windows will search your computer and the Internet for the latest driver software for your device, unless you've disabled this feature in your device installation settings.	
•	Browse my computer for driver software Locate and install driver software manually.	
		Cance

8. Haga clic en el botón Examinar, defina el escritorio del ordenador como ubicación y haga clic en el botón Siguiente.

Browse for driver software on your	computer
Search for driver software in this location:	
C\Users\j.smith\Desktop	• Browse
Let me pick from a list of device This list will show installed driver software of the software	e drivers on my computer
driver software in the same canegory as i	the device.

- 9. Espere hasta que el software del controlador esté instalado y acepte los mensajes de advertencia.
- 10. Una vez que se haya completado la instalación, registre la ubicación del puerto COM y haga clic en el botón Close.



11. El medidor ahora puede transferir datos al ordenador mediante el puerto COM virtual y el software informático Orion Star Com, HyperTerminal o un programa similar.

Software de comunicación Orion Star Com

El software de comunicación Orion Star Com es compatible con los medidores de sobremesa Orion Star serie A210; además, se encuentra disponible una descarga gratuita en <u>www.thermoscientific.com/OrionMeters</u>. El software Star Com facilita la transferencia de los datos de calibración y medida del medidor a un ordenador y, a continuación, permite que los datos se exporten a un archivo Excel (.xls) o a un archivo de valores separados por comas (.csv). Los datos transferidos también se pueden imprimir desde el ordenador.

Los medidores de sobremesa Orion Star serie A210 deben tener la revisión de software 2.59 o superior para poder utilizarlos con el software Star Com. El software del controlador USB de Orion Star serie A200/A300 debe instalarse en el ordenador si se va a utilizar el puerto USB para conectar el medidor con un ordenador.

Star Care								0.00
	<u>.</u>	. 💐 📑	<u>.</u>				4	
	_	_	_	416233490		Mar Celan		-
Channe 1		FH Measure	ment	Method 100	Made Tarle Ma	VEHSA STAL		8
					Date (Term	100 at 1 1 1 1 1 1 1 1 1 1 1 1 1 1 1 1 1 1		0
		8 00	DH		Contraction in the			-
		0.00			Chestrate 25 Sec	Restaura at		
25.0 °C (M	ANI			-57.8 mV	Moder & Mr.	And a set		Cher.
The number of the local division of the				Damages (1) Debarriel-CT	pril to dise	ACCU per BE SHAV		1.20
	1	15			Company Ann	and comme		E
	121	100		And in the local division of the local divis	44		10	1000
Contract of the last	Tax.	1000 at	an or of	- March				100
# ROOTALIMAN	111	1000 100	-28/510	36A)				Equat
10000 4 A 20 M FBA	818	6-000 am	20.1210	1000A33				
\$5/00714 H.(28 AB/MR	. #118	5-000 (m	10.0°C	-268-ly				
\$150014 x 28-54788	111	8-008 ptr	28.0.95	(989.4)				
1 1000 H N2030HW	111	and the	100.0110	(404(4)				
1 months of the control of the		BOOK of	20.010	- Marine - M				
a substruct of the local date	104	#1001 ptr	10.010	(SPAN)				
NUMBER OF TRADE	101	\$1000 pin	28.0 10	(HORNE)				
DOUBLING NOT BEEN PM	P11	BURY at	10.1110	(9044)				
900014400044	PH	6:007 64	38.1110	BRAN				
10000714-0-08.04786	111	800001 per	10.110	(HORN)				
1000214 x 28.387M	111	6/007.01	20.010	(9044)				
80007432944298	194	ROOM with	20.0 %	SOM				
THE REAL PROPERTY OF THE PARTY	111	81001.001	10.1110	0041				
10000214-02834/4E	10	5-001 pt	- 20.0 °C	and the second s			-	
THE REPORT OF THE REPORT								

ž 🕰 📩	<u>8 🙁 </u>	2 1	_	_	4	-
Channel 1	PH Messurement	Method 100	V Modes Type Modes Type No. Schwerz Tose Charren Ha Umered Ha Umer Hi	Dary Cours Second cours and do 248 Second courses for the	e Tra	16 14
25.0 °C (MAN) General Michael	0.001	-57.8 mV	Tarraina Distance 13 Ma Moder 13 Ma Moder 13 Ma Moder 13 Ma Moder Philippe Philippe Philippe Tarraina Vivian Tarraina Vivian Tarraina Santa	Samit, s. di Ebeografi Versite Addituti 20 a we 20 a Coloma 20 a Coloma		100 B
						-
				9		-
		6.997				I

Conexión del medidor con otros programas informáticos

Si utiliza el puerto USB del medidor para conectarlo con el ordenador, el software del controlador USB Orion Star serie A200/A300 (puerto COM virtual) debe instalarse en el ordenador y el medidor debe actualizarse a la revisión de software 3.04 o superior.

- 1. Encienda el medidor y conéctelo al ordenador mediante el cable de interfaz de ordenador adecuado.
- 2. Abra un programa de comunicación informático que utilice puertos COM. Serán aplicables programas como Terminal e HyperTerminal.
- Seleccione el puerto COM virtual correcto y conéctelo. Los ejemplos siguientes son de HyperTerminal:

	11.00		
FutBolings			
Dis per second 1000		Course To-	
Data bite 8		Contrast III	111.4
Purity Notes		rtyper 5	arting.
		Torus durade for	file phone member that you want to doin
Step 54k		Quality/region	Linker (States (S
Dow tarted bioto		Argue contre	10
		Etune rumber	
	Instante Derlauffer	Cognettoing	COM7
OR. Cance	And		
OS. Cano	full Aver		
OK Cares			
Denter Manager Denter Manager Die Johan Siew Belg In die Tal I die Tal die Aufli			
Denter Manager Den gaban Sjew Bela In de amina (12 min de jar de da 1 Mantana			
Device Manager De Johan Sjew Bela In Golan (2010) (2010) As (2010) Monitors - 10 Percent adapters - 10 Percent adapters			
0. Conser Device Manager bir & John View Bris bir & Immilia (Birmil Ac) pr AcAs ■ Manatan ■ Promotion Automos ■ Promotion Autom			
On Outcome Device Manager Device Manager De Annager De Manager De Device Annager			
On Overan Device Manager Bit Bit Status Service Manager Bit Bit Status Service Manager Bit Bit Bit Bit <td></td> <td></td> <td></td>			
Ore Conserver Manager Ein Schure Verwights Manager			

- 4. Compruebe que la exportación de datos del medidor esté activada. Ajuste la velocidad de baudios del medidor con el ajuste del ordenador. Si selecciona una velocidad de baudios mayor, la transferencia de datos será más rápida.
- 5. Los datos se transferirán del medidor al ordenador, en función del ajuste del tipo de lectura del medidor y de la instrucción.

Nota: Si el medidor está apagado por algún motivo (apagado automático, pérdida de potencia, etc.), deberá restablecerse la conexión repitiendo los pasos anteriores.

Protocolos de control remoto del medidor Star A200-A300

Reglas del uso de control remoto

- CR (Retorno de carro, ASCII 13) se utiliza para finalizar una instrucción. Siempre que se reciba este carácter, el búfer interno se procesará.
- Todos los caracteres excepto NL (Nueva línea, ASCII 10) son importantes. El carácter NL se ignorará. El medidor no distingue entre mayúsculas y minúsculas, por lo que todos los caracteres en minúscula se convertirán a mayúsculas internamente.
- Solo se puede ejecutar una instrucción a la vez. No se puede emitir una nueva instrucción hasta que la instrucción actual haya acabado de procesarse. Cuando la instrucción actual haya finalizado, emitirá un aviso ">" al usuario indicando que se puede introducir una nueva instrucción.
- 4. Las instrucciones vacías (como un CR) se ignorarán y se emitirá un nuevo aviso.
- 5. Active los ajustes del medidor Export Data y Data Log. Los ajustes de comunicación predeterminados del medidor son:

Bits por segundo:	9600
Bits de datos:	8
Paridad:	Ning.
Bits de parada:	1
Control de flujo:	Ning.

Motor de control remoto

El motor de control remoto recibe la señal del puerto serie/USB y la procesa según sea necesario. Las instrucciones enviadas a la interfaz de control remoto serán en formato "OPCODE <OPERAND(s)> c_R". Los avances de línea se ignorarán. No se pueden emitir nuevas instrucciones hasta que la instrucción anterior se haya completado y se haya dado un aviso. Se muestra un aviso como el símbolo de mayor (">") seguido por un espacio.

Instrucciones de control remoto del medidor Star A200-A300

Pulse la tecla **Esc** en el teclado del ordenador para detener una instrucción que se esté ejecutando.

Instrucción	Detalles de la instrucción
	GETMEAS <cr></cr>
	Imprime inmediatamente la medida del canal actual.
	GETMEAS Data Count <cr></cr>
	Imprime la medida del canal actual para una cantidad de veces determinada.
	Ejemplo: GETMEAS 2
	GETMEASTIMED CH_Channel Combination, Time Interval <cr></cr>
GETMEAS	Define el canal o canales que se medirán y el intervalo de tiempo de la medida.
	Las medidas se envían al registro de datos del medidor:
	Combinación de canales: Números de canales que se medirán.
	Intervalo de tiempo: El intervalo de tiempo de recuperación de la medida en
	segundos.
	Ejempio: GETMEASTIMED CH_12 5 <ck></ck>
	STOP <cr></cr>
	GETCAL MODE <cr></cr>
	Imprime todos los datos de calibración del modo de canal. Si no hay disponible pinguna calibración, so devolvorá ol valor ">" para regibir la instrucción siguionte
OFTON	Datos de calibración de impresión para el MODO específico:
GETCAL	
	GEILUG «CR»
	dato, se devolverá el valor ">" para recibir la instrucción siguiente. Fiemplo:
	GETLOG <cr></cr>
GETLOG	GETLOG START, END <cr></cr>
	Imprime los datos de un intervalo específico. Si no hay datos disponibles, se
	devolverá el valor ">" para recibir la instrucción siguiente.
	Ejemplo: GETLOG 10 100 <cr></cr>
	SYSTEM <cr></cr>
SYSTEM	Imprime el modelo del medidor, el número de serie y la versión de software
	Ejemplo: 329, 12345, 2,53
	SETRTC <u>YY MM DD HH MM SS</u> <cr></cr>
SETRTC	Defina la fecha y la hora (en formato de 24 horas) del medidor
	Ejemplo: SETRTC 13 07 15 08 30 00
SETMODE	SETMODE MODE <cr></cr>
	Defina el MODO de la medida del canal actual de la siguiente manera:
	PH COND OD
	RMV RES RDO
	ORP SALT
	ISE STD

Instrucción	Detalles de la instrucción
GETMODE	GETMODE <u>CHANNEL</u> <cr></cr>
	Imprime el modo de la medida del canal.
	Ejemplo: GETMODE 1
SETCSV	SETCSV <cr></cr>
	Define el formato de impresión a CSV
SETKEYLOC	SETKEYLOCK <u>NUMBER</u> <cr></cr>
	Para deshabilitar el teclado: Número = 0
IX	Para habilitar el teclado Número= 1

Datos de medida del medidor procedentes de la instrucción GETMEAS

Modo de medida	Formato PC (CSV)
рH	Modelo de medidor, número de serie, revisión de software, id. de usuario, fecha y hora, id. de la muestra, canal, modo, valor de pH, unidad de pH, valor de mV, unidad de mV, valor de temperatura, unidad de temperatura, valor de pendiente, unidad de pendiente, n.º de método, n.º de registro
	Ejemplo : A211 pH, X01036, 3.04, ABCDE, 01/03/15 16:05:41, SAMPLE, CH-1, pH, 7,000, pH, 0,0, mV, 25,0, C, 98,1,%, M100, #1 <cr></cr>
mV	Modelo de medidor, número de serie, revisión de software, id. de usuario, fecha y hora, id. de la muestra, canal, modo, valor de mV, unidad de mV, valor de temperatura, unidad de temperatura, n.º de método, n.º de registro
	Ejemplo : A211 pH, X01036, 3.04, ABCDE, 01/03/15 16:05:41, SAMPLE, CH-1, mV, 0,0, mV, 25,0, C, M100, #2 <cr></cr>
RmV	Modelo de medidor, número de serie, revisión de software, id. de usuario, fecha y hora, id. de la muestra, canal, modo, valor de mV relativo, unidad de mV relativa, valor de mV de compensación, unidad de mV de compensación, valor de temperatura, unidad de temperatura, n.º de método, n.º de registro
	Ejemplo : A211 pH, X01036, 3,04, ABCDE, 01/03/15 16:05:41, SAMPLE, CH-1, RmV, 0,0, RmV, 0,0, mV, 25,0, C, M100, #3 <cr></cr>
ORP	Modelo de medidor, número de serie, revisión de software, id. de usuario, fecha y hora, id. de la muestra, canal, modo, valor de ORP, unidad de ORP, valor de mV, unidad de mV, valor de temperatura, unidad de temperatura, n.º de método, n.º de registro
	Ejemplo: A211 pH, X01036, 3,04, ABCDE, 01/03/15 16:05:41, SAMPLE, CH-1, ORP, 0,0, mV, 0,0, mV, 25,0, C, M100, #4 <cr></cr>
ISE	Modelo de medidor, número de serie, revisión de software, ID de usuario, fecha y hora, ID de la muestra, canal, modo, valor de ISE, unidad de ISE, valor de mV, unidad de mV, valor de temperatura, unidad de temperatura, valor de pendiente, unidad de pendiente, n.º de método, n.º de registro
	Ejemplo: A214 pH/ISE, X01037, 3,04, ABCDE, 01/03/15 16:05:41, SAMPLE, CH-1, ISE, 1, ppb, 0,0, mV, 25,0, C, 59,2, mV/dec, M100, #1 <cr></cr>

Modo de medida	Formato PC (CSV)
Conducti- vidad	Modelo de medidor, número de serie, revisión de software, id. de usuario, fecha y hora, id. de la muestra, canal, modo, valor de conductividad, unidad de conductividad, valor de conductancia, unidad de conductancia, valor de temperatura, unidad de temperatura, valor de coeficiente de temperatura, unidad de coeficiente de temperatura, valor de temperatura de referencia, unidad de constante de celda, n.º de método, n.º de registro
	Ejemplo: A212 Cond, X01038, 3,04, ABCDE, 01/03/15 16:05:41, SAMPLE, CH-1, COND, 936,41, uS/cm, 1977,59, uS, 25,0, C, 2,1,%/C, 25,0, C, 0,4750, /cm, M100, #1 <cr>.</cr>
STD	Modelo de medidor, número de serie, revisión de software, id. de usuario, fecha y hora, id. de la muestra, canal, modo, valor de STD, unidad de STD, valor de conductancia, unidad de conductancia, valor de temperatura, unidad de temperatura, valor de temperatura de referencia, unidad de temperatura de referencia, valor de constante de celda, n.º de método, n.º de registro
	Ejemplo: A212 Cond, X01038, 3,04, ABCDE, 01/03/15 16:05:41, SAMPLE, CH-1, STD, 460, ppm, 1975, uS, 25,0, C, 2,1,%/C, 25,0, C, 0,4750, /cm, M100, #2 <cr></cr>
Salinidad	Modelo de medidor, número de serie, revisión de software, id. de usuario, fecha y hora, id. de la muestra, canal, modo, valor de salinidad, unidad de salinidad, valor de conductancia, unidad de conductancia, valor de temperatura, unidad de temperatura, tipo de salinidad, valor de temperatura de referencia, unidad de temperatura de referencia, valor de constante de celda, unidad de constante de celda, n.º de método, n.º de registro
	Ejemplo: A212 Cond, X01038, 3,04, ABCDE, 01/03/15 16:05:41, SAMPLE, CH-1, SALT, 0,5112, psu, 1973,120, uS, 25,0, C, Salinidad práctica, 15,0, C, 0,4750, /cm, M100, #3 <cr></cr>
Resistividad	Modelo de medidor, número de serie, revisión de software, id. de usuario, fecha y hora, id. de la muestra, canal, modo, valor de resistividad, unidad de resistividad, valor de conductancia, unidad de conductancia, valor de temperatura, unidad de temperatura, valor de coeficiente de temperatura, unidad de coeficiente de temperatura, valor de temperatura de referencia, unidad de temperatura de referencia, unidad de constante de celda, n.º de método, n.º de registro
	Ejemplo: A212 Cond, X01038, 3,04, ABCDE, 01/03/15 16:05:41, SAMPLE, CH-1, RES, 1068, Ohm- cm, 1982, uS, 25,0, C, 2,1,%/C, 25,0, C, 0,4750, /cm, M100, #4 <cr></cr>
OD % de saturación	Modelo de medidor, número de serie, revisión de software, id. de usuario, fecha y hora, id. de la muestra, canal, modo, porcentaje de valor de saturación, porcentaje de unidad de saturación, valor de concentración, unidad de concentración, valor actual, unidad actual, valor de temperatura de solución, unidad de temperatura de solución, valor de temperatura de membrana, unidad de temperatura de membrana, valor de presión barométrica, valor de corrección de salinidad, unidad de corrección de salinidad, valor de pendiente, unidad de pendiente, n.º de método, n.º de registro
	Ejemplo: A213 DO/RDO, X01039, 3,04, ABCDE, 01/03/15 16:05:41, SAMPLE, CH-1, DO, 0,3,% Sat, 0,03, mg/L, 3,4, nA, 24,5, C, 24,6, C, 761,0, mmHg, 0,0, ppt, 11,800, Na/%Sat, M100, #1 <cr></cr>

Modo de medida	Formato PC (CSV)
OD mg/L	Modelo de medidor, número de serie, revisión de software, id. de usuario, fecha y hora, id. de la muestra, canal, modo, valor de concentración, unidad de concentración, porcentaje de valor de saturación, unidad de valor de saturación, valor actual, unidad actual, valor de temperatura de solución, unidad de temperatura de solución, valor de temperatura de membrana, unidad de temperatura de membrana, valor de presión barométrica, unidad de presión barométrica, unidad de presión de salinidad, unidad de corrección de salinidad, valor de pendiente, unidad de pendiente, n.º de método, n.º de registro
	Ejemplo: A213 DO/RDO, X01039, 3,04, ABCDE, 01/03/15 16:05:41, SAMPLE, CH-1, DO, 0,03, mg/L, 0,3,% sat, 3.4, nA, 24,5, C, 24,6, C, 761,0, mmHg, 0,0, ppt, 11,800, Na/%Sat, M100, #2 <cr></cr>
RDO % de	Modelo de medidor, número de serie, revisión de software, id. de usuario, fecha y hora, id. de la muestra, canal, modo, porcentaje de valor de saturación, porcentaje de unidad de saturación, valor de concentración, unidad de concentración, valor de presión parcial, unidad de presión parcial, valor de temperatura de solución, unidad de temperatura de solución, valor de presión barométrica, unidad de presión barométrica, unidad de presión de salinidad, unidad de valor de salinidad, valor de pendiente, unidad de pendiente, n.º de método, n.º de registro
	Ejemplo: A213 DO/RDO, X01039, 3,04, ABCDE, 01/03/15 16:05:41, SAMPLE, CH-1, RDO, 100,0,% Sat, 8,40, mg/L, 138,5, Torr, 24,0, C, 761,0, mmHg, 0,0, ppt, 1,383, Torr/%Sat, M100, #3 <cr></cr>
RDO mg/L	Modelo de medidor, número de serie, revisión de software, id. de usuario, fecha y hora, id. de la muestra, canal, modo, valor de concentración, unidad de concentración, porcentaje de valor de saturación, unidad de valor de saturación, valor de presión parcial, unidad de presión parcial, valor de temperatura de solución, unidad de temperatura de solución, valor de presión barométrica, unidad de presión barométrica, unidad de presión de salinidad, unidad de corrección de salinidad, valor de presión de presión de presión de solución, valor de presión barométrica, unidad de presión barométrica, unidad de presión de salinidad, unidad de corrección de salinidad, valor de pendiente, unidad de pendiente, n.º de método, n.º de registro
	Ejemplo: A213 DO/RDO, X01039, 3,04, ABCDE, 01/03/15 16:05:41, SAMPLE, CH-1, RDO, 8,40, mg/L, 100,0,% sat, 138,5, Torr, 24,0, C, 761,0, mmHg, 0,0, ppt, 1,383, Torr/%Sat, M100, #4 <cr></cr>
	Modelo de medidor, número de serie, revisión de software, id. de usuario, fecha y hora, id. de la muestra (seguido por los datos de medida de cada canal, como se muestra anteriormente)
Multicanal	Ejemplo: A215 pH/Cond, X01040, 3,04, ABCDE, 01/03/15 16:05:41, SAMPLE, CH-1, pH, 7,000, pH, 0,0, mV, 25,0, C, 98,1,%, M100, CH-2, COND, 936,41, uS/cm, 1977,59, uS, 25,0, C, 2,1,%/C, 25,0, C, 0,4750, /cm, M200, #1 <cr></cr>

Modo Cal	Formato PC (CSV)
	Modelo de medidor, número de serie, revisión de software, canal, modo, fecha y hora de calibración, puntos de calibración totales
рН	repetido para cada punto de calibración: Punto de calibración, valor de pH, unidad de pH, valor de mV, unidad de mV, valor de temperatura, unidad de temperatura,
	repetido para cada segmento punto a punto:Valor de pendiente, unidad de pendiente,
	repetido para cada segmento punto a punto: Valor de compensación, unidad de compensación, valor de pendiente medio, unidad de pendiente media, número de calibración
	Ejemplo (calibración de 5 puntos):
	A211 pH, X01036, 3,04, CH-1, pH, 01/03/15 16:05:41, 5, 1, 2,00, pH, 290,5, mV, 25,0, C, 2, 4,01, pH, 174,3, mV, 25,0, C, 3, 7,00, pH, 0,0, mV, 25,0, C, 4, 10,01, pH -173,9, mV, 25,0, C, 5, 12,00, pH, -289,7, mV, 25,0, C, 97,7,%, 98,5,%, 97,7,%, 98,4,%, 1,4, mV, 0,0, mV, 0,0, mV, 1,3, mV, 98,1,%, #1 <cr></cr>
RmV	Modelo de medidor, número de serie, revisión de software, canal, modo, fecha y hora de calibración, valor mV bruto, unidad mV bruta, valor de compensación mV relativa, unidad de compensación mV relativa, valor mV de referencia, unidad mV de referencia, número de calibración
	Ejemplo:
	A211 pH, X01036, 3,04, CH-1, RmV, 01/03/15 16:05:41, 0,0, mV, 0,0, mV, 0,0, mV, #1 <cr></cr>
ORP	Modelo de medidor, número de serie, revisión de software, canal, modo, fecha y hora de calibración, valor mV bruto, unidad mV bruta, valor de compensación, unidad de compensación, valor de temperatura, unidad de temperatura, número de calibración
	Ejemplo:
	A211 pH, X01036, 3,04, CH-1, ORP, 01/03/15 16:05:41, 176,9, mV, 242,6, mV, 25,0, C, #1 <cr></cr>
	Modelo de medidor, número de serie, revisión de software, canal, modo, fecha y hora de calibración, puntos de calibración totales
ISE	repetido para cada punto de calibración:Punto de calibración, valor concentración, unidad de concentración, valor de mV, unidad de mV, valor de temperatura, unidad de temperatura,
	repetido para cada segmento punto a punto:Valor de pendiente, unidad de pendiente,
	repetido para cada segmento punto a punto:Valor de compensación, unida de compensación, valor de pendiente medio, unidad de pendiente media, valo en blanco, unidad en blanco, número de calibración
	Ejemplo (calibración de 5 puntos):
	A214 pH/ISE, X01036, 3,04, CH-1, ISE, 01/03/15 16:05:41, 5, 1, -1, ppb, 0,0, mV 25,0, C, 2, 2, ppb, 19,0, mV, 25,0, C, 3, 300, ppb, 150,1, mV, 25,0, C, 4, -2000, ppb, 200,1, mV, 25,0, C, 5, 4000, ppb, 210,1, mV, 25,0, C, 63,1, mV/dec, 60,2, mV/dec, 60,7, mV/dec, 33,2, mV/dec, 0,0, mV, 0,9 mV, 0,0 mV, -0,2 mV, 90,4, mV 54,3, mV/dec, 0,0000, ppb, #1 <cr></cr>

Datos de calibración del medidor a partir de la instrucción GETCAL

Modo Cal	Formato PC (CSV)
Conductividad	Modelo de medidor, número de serie, revisión de software, canal, modo, fecha y hora de calibración, puntos de calibración totales, repetido para cada punto de calibración: Punto de calibración, valor de conductividad, unidad de conductividad, valor de conductancia, unidad de conductancia, valor de temperatura, unidad de temperatura (MTC/ATC), tipo de calibración, factor de calibración, constante de celda media (K), número de calibración
Conductividad	Ejemplo (calibración de 5 puntos): A212 Cond, X01038, 3,04, CH-1, Cond, 01/03/15 16:05:41, 5, 1, 52,8, uS/cm, 111,1, uS, 25,0, C, Manual, 0,4750, 2, 68,6, uS/cm, 143,1, uS, 25,0, C, Manual, 0,4956, 3, 100, uS/cm, 200,2, uS, 25,0, C, Manual, 0,5511, 4, 158,4, uS/cm, 333,4, uS, 25,0, C, Manual, 0,4375, 5, 475, uS/cm, 1000, uS, 25,0, C, Manual, 0,4749, 0,4868 #1 <cr></cr>
OD (Calibración de aire o agua)	Modelo de medidor, número de serie, revisión de software, canal, modo, fecha y hora de la calibración, punto de calibración, tipo de calibración, valor de saturación de calibración, unidad de saturación de calibración, valor actual de calibración de definición cero, unidad actual de calibración, valor de temperatura de solución, unidad de temperatura de solución, valor de temperatura de solución, unidad de temperatura de membrana, valor de temperatura de membrana, unidad de temperatura de membrana, valor de temperatura de membrana, unidad de calibración, valor de temperatura de membrana, unidad de temperatura de membrana, valor de presión de calibración, unidad de pendiente de calibración de calibración, número de calibración
	Ejemplo: A213 DO/RDO, X01039, 3,04, CH-1, DO, 01/03/15 16:05:41, 1, Aire automático, 102,3,%, 0,0, nA, 405,1, nA, 5,0, C, 5,0, C, 736,1, mmHg, 0,0, ppt, 4,1, nA/% sat, 1, #1 <cr></cr>
OD (Calibración manual)	Modelo de medidor, número de serie, revisión de software, canal, modo, fecha y hora de la calibración, punto de calibración, tipo de calibración, valor de concentración de calibración, unidad de concentración de calibración, valor actual de calibración de definición cero, unidad actual de calibración, valor de temperatura de solución, unidad de temperatura de solución, valor de temperatura de membrana, unidad de temperatura de membrana, valor de presión de calibración, unidad de presión de calibración, unidad de pendiente de calibración, unidad de pendiente de calibración, unidad de pendiente de calibración, número de calibración, número de calibración de pendiente de calibración de calibración de calibración de pendiente de calibración de calibración de temperatura de temperatur
	Ejemplo: A213 DO/RDO, X01039, 3,04, CH-1, DO, 01/03/15 16:05:41, 1, Manual, 12,62, mg/l, 0,0, nA, 405,1, nA, 5,0, C, 5,0, C, 736,1, mmHg, 0,0, ppt, 4,1, nA/%sat, 1, #2 <cr></cr>
RDO (Calibración de aire o agua)	Modelo de medidor, número de serie, revisión de software, canal, modo, fecha y hora de la calibración, número de serie de la sonda, punto de calibración, tipo de calibración, valor de saturación de calibración, unidad de saturación de calibración, valor de presión parcial de definición cero, unidad de presión parcial de definición cero, unidad de presión parcial de calibración, valor de temperatura de solución, unidad de temperatura de solución, valor de presión barométrica, unidad de presión barométrica, valor de parcial de calibración, unidad de pendiente de calibración, unidad de pendiente de calibración, unidad de pendiente de calibración parcial de calibración, número de calibración
	Ejemplo: A213 DO/RDO, X01039, 3,04, CH-1, RDO, 01/03/15 16:05:41, 123454, 1, Aire automático, 100,0,%, 8,2, Torr, 149,6, Torr, 24,3, C(ATC), 749,7, mmHg, 0,1, ppt, 1,52, Torr/%sat, 1, #1 <cr></cr>

Modo Cal	Formato PC (CSV)
RDO (Calibración manual)	Modelo de medidor, número de serie, revisión de software, canal, modo, fecha y hora de la calibración, número de serie de la sonda, punto de calibración, tipo de calibración, valor de concentración de calibración, unidad de concentración de calibración, valor de presión parcial de definición cero, unidad de presión parcial de definición cero, valor de presión parcial de calibración, unidad de presión parcial de calibración, unidad de temperatura de solución, valor de presión barométrica, unidad de presión barométrica, valor de salinidad, unidad de salinidad, valor de pendiente de calibración, unidad de pendiente de calibración, punto de calibración, número de calibración
	Ejemplo: A213 DO/RDO, X01039, 3,04, CH-1, RDO, 12/16/12, 09:21:00, 123454, 1, Manual, 6,9, mg/l, 8,2, Torr, 149,6, Torr, 24,3, C(ATC), 749,7, mmHg, 0,1, ppt, 1,52, Torr/% sat, 1, #1 <cr></cr>

Procedimiento de actualización de software del medidor

El programa de actualización de software para los medidores de las series Orion Star A200 y Orion A300 se ha probado únicamente con los sistemas operativos Windows 7, XP y Vista de Microsoft. Trabajamos constantemente para actualizar nuestros programas; sin embargo, el programa de actualización de software, actualmente, no se ha probado con los otros sistemas operativos.

Todos los ajustes de gestión de potencia y de protectores de pantalla deben estar desactivados cuando el medidor está conectado con el ordenador. El ordenador no puede entrar en modo de suspensión o de espera mientras se está llevando a cabo la actualización de software. Muchos ordenadores portátiles entran en modo de suspensión o de espera cuando se cierra la tapa, por lo que la tapa debe permanecer abierta durante toda la actualización de software. Los ordenadores portátiles también deben conectarse a una fuente de alimentación cuando se lleva a cabo la actualización de software; no se recomienda realizar la actualización de software con el ordenador funcionando con alimentación de la batería.

Nota: Realice una copia de seguridad de los datos del medidor antes de actualizar el software.

- Vaya a <u>www.thermoscientific.com/OrionMeters</u>, descargue la última versión de software de los medidores de la serie Orion Star A200-A300 y descomprima/extraiga los archivos de software al escritorio del ordenador.
 - a. Existen dos versiones de la actualización de software del medidor de la serie Orion Star A200-A300: para los sistemas operativos de Windows de 32 bits y para los sistemas operativos de Windows de 64 bits. Par determinar cuál debe utilizar con su ordenador:
 - i. Para Windows 7: Haga clic en el botón Inicio y, a continuación, en Panel de control. En la ventana del Panel de control, haga clic en Sistema si la visualización es por iconos grandes o iconos pequeños o haga clic en Sistema y seguridad y, a continuación, en Sistema si la visualización es por categoría. En la ventana Sistema, junto a Tipo de sistema, se mostrará 32 bits o 64 bits.
 - ii. Para Windows XP: Haga clic en el botón Inicio y, a continuación, en Panel de control. En la ventana Panel de control, haga clic en el enlace Rendimiento y mantenimiento y, a continuación, en el icono Sistema o, para la vista clásica del Panel de control, haga doble clic en el icono Sistema. En la ventana Propiedades del sistema, haga clic en la pestaña General y, debajo de Sistema, si se muestra "edición x64", utilice la versión de 64 bits y, si no se muestra "edición x64", utilice la versión de 32 bits.

 Desinstale cualquier versión anterior de la actualización de software del medidor de la serie Orion Star A200-A300 con la herramienta Programas y características del ordenador.



- Asegúrese de que el cable de interfaz del ordenador esté conectado al medidor y al ordenador y que se haya instalado completamente mediante las instrucciones que aparecen en la sección <u>Uso del cable de ordenador USB a serie</u> o <u>Uso del cable de</u> <u>ordenador USB</u>.
- Encienda el medidor. Para garantizar una alimentación sin interrupciones durante todo el proceso de actualización, utilice el adaptador de alimentación con los medidores de sobremesa y pilas completamente cargadas o el adaptador de alimentación con los medidores portátiles.
- Abra la carpeta que contiene los archivos de software descomprimidos/extraídos y haga doble clic en el archivo setup.exe para instalar el programa de instalación en el ordenador.



 Siga todos los avisos para instalar el programa de instalación en el ordenador. Si aparece algún mensaje de advertencia, seleccione la opción de ejecución del software.



 Después de la instalación correcta del programa de instalación, aparecerá un nuevo icono con el nombre StarA200_300SoftwareUpgrade en el escritorio del ordenador. Haga doble clic en el icono para iniciar el programa.

ORION A200 & A300 Series bootloader v0.21	
ORION A200 and Firmware U	A300 Series ograde
SCIENTIF	0
	Press Next to Proceed Next > Close

- 8. Siga todos los avisos, incluidas la revisión y la aceptación del contrato de licencia, para iniciar la actualización de software. Si aparece algún mensaje de advertencia, seleccione la opción de ejecución del software.
- 9. Defina los parámetros de la interfaz para ejecutar el programa de actualización de software.
 - a. Cuando utilice un cable de ordenador USB a serie, seleccione Serial (Serie) como método de conexión, introduzca la ubicación del puerto COM (visible en Administrador de dispositivos debajo de Puertos) e introduzca la velocidad de baudios del medidor actual (el ajuste predeterminado es 9600).

ORION A200 & A300 Series bootloa	ader v0.21.304		
1. Choose connection method 2. Attach meter to PC 3. Click Upgrade Firmware	 Serial 	COM8	9600
		Up	grade Firmware
			1

b. Cuando utilice un cable de ordenador USB, la actualización de software detectará la conexión USB (el controlador USB debe estar instalado antes de iniciar la actualización de software).

- 10. Haga clic en el botón Upgrade Firmware (Actualizar firmware). Espere hasta que se haya instalado la actualización en el medidor.
 - a. **Advertencia:** La actualización de software no debe interrumpirse una vez que se ha hecho clic en el botón Update Firmware.La interrupción de la actualización de software puede dañar el medidor y hacer que no se encienda correctamente.

ORION A200 & A300 Series bootic	ader v0.21.304			
1 Choose connection method 2 Atlach meter to PC 3 Click Upgrade Firmware	③ Serial	COMS	9600 pade Fernware	
Starting init communication Set BOOT FLAGS BINs loading bootstrap loading				

 La actualización tardará aproximadamente 15 minutos. Una vez que la actualización de software se haya completado, el cuadro de diálogo mostrará "Done!" (¡Realizado!)

1. Choose connecton method 2. Attach meter to PC 3. Click Upgrade Firmware	🥥 Senal	COM8	9600
		Lip	grade Firmware
ain code loading ating for reboot eanup			

12. Cierre la actualización de software y desconecte el cable del medidor y del ordenador.

Nota: Si aparece el mensaje siguiente y el cable USB se está utilizando, apague el medidor desconectando el adaptador de alimentación o quite las pilas, desconecte el cable USB, cierre la actualización de software y vuelva a intentar el procedimiento con el cable RS232.

DRDN ADD & ADD Server barrhower VL/LIDH	Trad in Ma
Passand memory hope	
Please remove batteries and power	
wait 30 seconda	
then reconnect power	
	1000
Sadry Gary Advise Town Halp down watch in Analytic Watch	



CAPÍTULO 8 Servicios de atención al cliente

Si tiene alguna pregunta o necesita ayuda, póngase en contacto con nuestros especialistas en soporte técnico:

- Correo electrónico wai.techservbev@thermofisher.com
- En los Estados Unidos, llame al 1-800-225-1480
- Fuera de los Estados Unidos, llame al +1 -978-232-6000, o mediante fax al +1 -978-232-6031

Para obtener información adicional sobre el producto, póngase en contacto con su distribuidor autorizado local, su representante de ventas técnicas local de Thermo Scientific Orion o póngase en contacto con nosotros mediante la información de Water and Laboratory Products (WLP) de la página final de este manual.

Visite <u>www.thermoscientific.com/water</u> para visualizar productos de Thermo Scientific Orion y descargar documentación sobre el producto, guías de usuario y manuales, actualizaciones de software y otras aplicaciones y recursos técnicos.

Para obtener la información de garantía más actual, consulte la tarjeta de garantía de Thermo Scientific Orion incluida en el CD de documentación del medidor Thermo Scientific Orion Star serie A210; también se encuentra disponible en línea en www.thermoscientific.com/water.

Sugerencias de solución de problemas

Si se detecta algún problema mientras utiliza el medidor Orion Star serie A210, asegúrese que se utiliza el adaptador de alimentación correcto e intente realizar un ciclo de alimentación en el medidor: desconecte el adaptador de alimentación del medidor, espere 15 segundos y vuélvalo a conectar.

Problema del medidor	Acción recomendada
El valor de medida parpadea 9999 y muestra Over Range (Encima del rango) o Under Range (Debajo del rango)	La medida se encuentra fuera del intervalo de medida permitido. Compruebe que se muestre el canal correcto en la pantalla del medidor. Asegúrese de que el electrodo o el sensor estén completamente conectados al medidor.
El medidor se bloquea cuando está conectado a un ordenador a través del cable USB.	Desconecte el cable USB del medidor y el ordenador y, a continuación, desconecte y vuelva a conectar el adaptador de alimentación del medidor. Restablezca la conexión USB del medidor y el ordenador.
El teclado del medidor no responde al utilizar el software informático Star Com	La opción de bloqueo de teclado está activada en el programa informático Star Com. Para desactivar el bloqueo de teclado: en el programa Star Com, haga clic en el icono Settings, desmarque la casilla que hay junto a Keypad Lock (Bloqueo de teclado) y pulse el icono Save (Guardar).
El valor de medida se detiene y no cambia	El tipo de lectura está definido en el modo de lectura automática (icono AR que se muestra en la pantalla). Pulse la tecla measure (esc) para tomar una nueva medida o utilice el menú de configuración para cambiar el tipo de lectura a continua.
La pantalla del medidor queda en blanco, muestra líneas aleatorias o detenciones intermitentes	Asegúrese que se utiliza la fuente de alimentación correcta para el medidor de la serie Star A. Esta fuente de alimentación es distinta de la suministrada con los medidores Thermo Scientific Orion Versa Star. Se recomienda el uso de un protector contra sobretensión o sistema de alimentación ininterrumpida (SAI). Reinicie el medidor.
El medidor no reconoce automáticamente el búfer de pH durante la calibración.	Compruebe que se haya seleccionado el conjunto de búfer correcto en el menú de configuración. El medidor utiliza lecturas mV sin procesar para reconocer el búfer. A medida que el electrodo se deteriore o se ensucie, sus lecturas mV fallarán. Compruebe los búferes y limpie el electrodo de acuerdo con las instrucciones del manual de electrodos.
El medidor no reconoce el estándar de conductividad durante la calibración	Compruebe que se haya introducido la constante de la celda nominal correcta en el menú de configuración para que se pueda calibrar el sensor de conductividad. La constante de la celda se imprime normalmente en el cable del sensor de conductividad. Compruebe que el estándar de conductividad se pueda reconocer automáticamente por el medidor. Vuelva a calibrar el sensor de conductividad con un estándar de conductividad nuevo.
El medidor no reconoce un sensor de oxígeno disuelto polarográfico u óptico RDO.	Asegúrese de que el sensor de oxígeno disuelto esté completamente conectado al medidor y compruebe que se muestre el canal correcto en la pantalla del medidor. Espere 15 segundos tras la conexión del sensor de oxígeno disuelto para que el medidor reconozca el tipo de sensor de OD. Pulse la tecla measure (esc) para iniciar una nueva medida y actualice la pantalla del medidor.

Procedimiento de restablecimiento de fábrica del medidor

Advertencia: Un restablecimiento de fábrica borrará el registro de calibración y el registro de datos del medidor y restablecerá todos los parámetros de configuración del medidor a la configuración predeterminada de fábrica.

- 1. En el modo de medida, pulse la tecla setup.
- 2. Pulse la tecla \blacktriangle , \triangledown , \triangleleft o \triangleright para resaltar <u>Diagnostics</u> y pulse la tecla **f3 (select)**.
- Pulse la tecla ▲ o ▼ para resaltar <u>Factory Reset</u> (Rest. valores fábrica) y pulse la tecla f2 (select).
- 4. Pulse la tecla f2 (yes) para iniciar el procedimiento de restablecimiento de fábrica.
- 5. Pulse la tecla **f3 (edit)** para acceder a la ventana emergente de introducción de números e introduzca la contraseña del medidor predeterminada de 111111.
 - a. Pulse la tecla ▲, ▼, ◀ o ▶ para resaltar el número <u>1</u> y pulse la tecla <u>f3 (enter)</u> seis veces hasta que el número <u>111111</u> se muestre en la parte superior de la pantalla.
 - b. Pulse la tecla **f2 (done)** para guardar el valor y salir de la ventana de introducción de números.
- 6. Pulse la tecla f2 (accept) para iniciar el restablecimiento de fábrica.
- 7. Espere hasta que se haya realizado el restablecimiento de fábrica. Una vez finalizado, el medidor realizará un ciclo de alimentación y, a continuación, seguirá con el modo de medida. Todos los ajustes del medidor se restablecerán en la configuración predeterminada de fábrica y el registro de calibración y el registro de datos se borrarán.

Procedimiento de restablecimiento del usuario del medidor

- 1. En el modo de medida, pulse la tecla setup.
- Pulse la tecla ▲, ▼, ◀ o ▶ para resaltar <u>Diagnostics</u> y pulse la tecla f3 (select).
- Pulse la tecla ▲ o ▼ para resaltar <u>User Reset</u> (Restablecimiento de usuario) y pulse la tecla f2 (select).
- 4. Pulse la tecla f2 (yes) para iniciar el procedimiento de restablecimiento de usuario.
- 5. Espere hasta que se haya realizado el restablecimiento de usuario. Una vez finalizado, el medidor seguirá con el modo de medida. Todos los ajustes del medidor se restablecerán en la configuración predeterminada de fábrica; sin embargo, el registro de calibración y el registro de datos se conservarán.

Declaración de conformidad

Este equipo genera, utiliza y puede emitir energía de radiofrecuencia y, si no se instala y utiliza de acuerdo con el manual de instrucciones, puede causar interferencias en las comunicaciones de radio. Ha sido probado y se ha determinado que cumple con los límites de los dispositivos digitales de Clase A, de conformidad con la parte 15 de las normas del FCC, que están pensadas para proporcionar protección razonable contra dicha interferencia cuando se utilizan en un entorno comercial. Es posible que el funcionamiento de este equipo en una zona residencial provoque interferencias, en cuyo caso el usuario, a sus expensas, tendrá que tomar las medidas necesarias para corregir dichas interferencias.

"Este aparato digital no excede los límites (Clase A) para emisiones de radio desde aparatos digitales estipulados en las Normas de interferencia de radio del Departamento de Comunicaciones del Canadá".

"Le present appareil numerique n' emet pas de bruits radioelectriques depassant les limites applicables aux appareils numeriques (de la class A) prescrites dans le Reglement sur le brouillage radioelectrique edicte par le ministere des Communications du Canada."

Conformidad RAEE

Este producto debe cumplir con la directiva RAEE sobre Residuos de Aparatos Eléctricos y Electrónicos de la Unión Europea (2002/96/EC). Va marcado con el siguiente símbolo.



Para obtener más información sobre nuestra conformidad con estas directivas, las empresas de reciclaje de su país y los productos de Thermo Scientific que pueden facilitar la detección de sustancias incluidas en la Directiva RoHS, póngase en contacto con nosotros mediante la información de Water and Laboratory Products (WLP) de la página final de este manual.

Declaración de conformidad

Fabricante:Thermo Fisher Scientific IncDirección:Ayer Rajah CrescentBlk 55 #04-16/24Singapur 139949SingapurSingapur

Declara, por la presente, que los productos siguientes:

Los medidores de sobremesa Thermo Scientific Orion Star serie A210 admiten una capacidad nominal de 100 a 240 VCA, 50/60 Hz, 0,5 A

Modelos de medidores portátiles:

Medidor de pH Orion Star A211 Conductímetro Orion Star A212 Medidor de RDO/OD Orion Star A213 Medidor de pH/ISE Orion Star A214 Medidor de pH/conductividad Orion Star A215 Medidor de pH/RDO/OD Orion Star A216

Clase del equipamiento:

Los medidores Orion Star serie A para medida, control y uso en laboratorio cumplen los requisitos de compatibilidad electromagnética (CEM) de Clase A

Cumplen con las siguientes directivas y estándares:

EN61326-1:2013	Directiva de compatibilidad electromagnética (EMC)
	Material eléctrico para medida, control y uso en laboratorio. Requisitos de compatibilidad electromagnética (CEM).
EN61010-1:2010	Estándares de seguridad
UL61010-1:2012	Requisitos de seguridad para equipamiento eléctrico de medida,
CAN/CSA C22.2 No. 61010-1:2012	Control y uso en laboratorio: requisitos generales.



Cheow Kwang Chan Responsable regulador y de control de calidad

Lugar y fecha de emisión: 16 de febrero de 2015 Singapur

Especificaciones del medidor

3 years	CE	0	FC
-			

Condiciones de funcionamiento del medidor de sobremesa Orion Star serie A210		
Temperatura ambiente de funcionamiento	De 5 a 45 °C	
Humedad relativa de funcionamiento	De 5 % a 85 %, sin condensación	
Temperatura de almacenamiento	De -20 a 60 °C	
Humedad relativa de almacenamiento	De 5 % a 85 %, sin condensación	
Contaminación	Grado 2	
Sobretensión	Categoría II	
Peso del medidor	0,9 kg (2 lb)	
Dimensiones del medidor (Al x An x Pr)	11 x 18 x 24 cm (4,33 x 7,09 x 9,45 pulgadas)	
Regulación y seguridad	CE, TUV 3-1, FCC Clase A	
Potencia nominal	Entrada de CC: 9 VCC, 1 A	
	Pilas: 4 pilas AA	
Choque y vibración	Choque: prueba de caída en embalaje por ISTA #1A	
	Vibración: envío/manipulación por ISTA #1A	
Caja (diseñada para cumplir con las normas)	IP54	
Garantía	3 años	

Condiciones de funcionamiento del adaptador de alimentación universal		
Temperatura ambiente de funcionamiento	De 0 a 50 °C	
Humedad relativa de funcionamiento	De 0% a 90%, sin condensación	
Temperatura de almacenamiento	De -20 a 75 °C	
Humedad relativa de almacenamiento	De 0% a 90%, sin condensación	
Contaminación	Grado 2	
Sobretensión	Categoría II	

Especificaciones del medidor de sobremesa Orion Star serie A210		
Canales de medida	De 1 a 2	
Pantalla	LCD gráfica	
Modos de medida	Lectura automática, continua, temporizada	
Memoria del registro de datos	2000 puntos de datos con registro de fecha y hora	
Funciones del registro de datos	Registro automático con modos de medida de lectura automática y temporizada; registro manual con modo de medida continua	
Transferencia del registro de datos	Transferencia simple, de rango o de todos los puntos de datos a la impresora o al ordenador	
Edición del registro de datos	Borrar todos los puntos de datos	

Especificaciones del medidor de sobremesa Orion Star serie A210			
Registro de calibración	10 calibraciones por canal con registro de fecha y hora		
Software	Software de transferencia de datos Orion Star Com disponible gratuitamente en el sitio web		
Métodos	10 por canal con protección de contraseña		
Alarmas	Alarma de límite, alarma de calibración		
ld. de la muestra	Introducción manual, de incremento automático o apagado		
ID de usuario	Introducción manual o apagado		
Puertos de comunicación	RS232, USB		
Puertos de la sonda de agitador	1		
Fuente de alimentación	Adaptador de alimentación universal CA, 90-260 VCA, 50-60 Hz o 4 pilas AA opcionales (promedio de 800 horas de vida de las pilas)		
Interfaz de usuario en varios idiomas	Inglés, español, francés, italiano, alemán y chino Portugués y coreano disponibles con la actualización de software en la web		
Especificaciones del medidor de sobremesa de pH Orion Star A211			
---	--------------------------------	--	--
Canales de medida		1 - pH, mV, mV relativo (RmV) u ORP con temperatura	
	Intervalo	de -2,000 a 20,000	
	Resolución	0,1 / 0,01 / 0,001	
	Exactitud relativa	±0,002	
н	Puntos de calibración	Hasta 5	
P	Edición de calibración	Sí	
	Impedancia de entrada	> 10 ^{^12} ohmios	
	Estado del electrodo	Indicador en la pantalla de estado del electrodo bueno, normal o malo	
	Intervalo	±2000,0 mV	
	Resolución	0,1 mV	
mV/RmV	Exactitud relativa	\pm 0,2 mV o \pm 0,05 % de la lectura, lo que sea mayor	
	Modo mV relativo	Sí	
	Modo Eh ORP	Sí	
	Intervalo	De -5 a 105 °C (de 22 a 221 °F)	
	Resolución	0,1 °C, 0,1 °F	
Temperatura	Exactitud relativa	±0,1 °C	
remperatura	Calibración de compensación	1 punto	
	Opciones de alimentación	Manual o automática con sonda ATC	
	BNC	Electrodo de pH u ORP	
Entradas del sensor	Punta	Electrodo de referencia	
	MiniDIN de 8 pines	Sonda de temperatura ATC	

Especificaciones del medidor de sobremesa de conductividad Orion Star A212			
Canales de medida		1 - conductividad, salinidad, STD o resistividad con temperatura	
	Intervalo	De 0,001 JS a 3000 mS	
	Resolución	0,001 jS mínimo, alcance automático, hasta 4 dígitos significativos	
	Exactitud relativa	0,5 % de la lectura ±1 dígito >3 jS, 0,5 % de la lectura ±0,01 jS <3 jS	
Conducti-	Temperatura de referencia	5 °C, 10 °C, 15 °C, 20 °C, 25 °C	
vidad	Compensación de temperatura	Lineal (de 0 a 10,0 %/°C), nLFn no lineal, nLFu no lineal, E (USP), apagado	
	Constantes de celda	De 0,001 a 199,9 cm-1	
	Puntos de calibración	Hasta 5	
	Edición de calibración	Sí	
	Intervalo	De 0,06 a 80,00 psu, de 0,05 a 42,00 ppt	
Salinidad	Resolución	0,01 psu o 0,01 ppt mínimo, alcance automático	
Sainnuau	Exactitud relativa	0,5 % de la lectura ±1 dígito	
	Тіро	Salinidad práctica (psu) o agua marina natural (ppt)	
	Intervalo	De 0,001 a 200,0 ppm	
STD	Resolución	0,001 ppm mínimo, alcance automático, hasta 4 dígitos significativos	
	Exactitud relativa	0,5 % de la lectura ±1 dígito	
	Factor STD	Lineal (de 0,02 a 9,99)	
	Intervalo	De 2 Q a 100 MQ	
Resistividad	Resolución	1 Q o 0,1 MQ, alcance automático	
	Exactitud relativa	0,5 % de la lectura ±1 dígito	
	Intervalo	De -5 a 105 °C (de 22 a 221 °F)	
Temperatura	Resolución	0,1 °C, 0,1 °F	
	Exactitud relativa	±0,1 °C	
	Calibración de compensación	1 punto	
	Opciones de alimentación	Manual o automática con sensor de temperatura integrado	
Entrada del sensor	MiniDIN de 8 pines	Sensor de conductividad con temperatura integrada	

Especificaciones del medidor de sobremesa de RDO/OD Orion Star A213			
Canales de medida		1 - oxígeno disuelto como % de saturación o mg/L con temperatura	
	Polarográfico	Concentración	% de saturación
	Intervalo	De 0 a 90 mg/L	De 0 a 600 %
	Resolución	0,01 / 0,1 mg/L	0,1 / 1 %
	Exactitud relativa	±0,2 mg/L o ±2 % de la lectura, lo que sea mayor	±2 % de saturación o ±2 % de la lectura, lo que sea mayor
Oxígeno	Óptico RDO	Concentración	% de saturación
disuello	Intervalo	De 0 a 50 mg/L	De 0 a 500 %
	Resolución	0,01 / 0,1 mg/L	0,1 / 1 %
	Exactitud relativa	De ±0,1 mg/L, hasta 8 mg/L, De ±0,2 mg/L, de 8 a 20 mg/L, ±10 % de lectura hasta 50 mg/L	±2 % de saturación < 200 % de saturación, ±10 % de saturación > 200 % de saturación
Corrección de la presión barométrica		De 400 a 850 mm Hg, automática con barómetro integrado (±6 mm Hg) o introducción manual	
Factor de corrección de salinidad		De 0,0 a 45,0 ppt, automático con introducción manual de la salinidad de muestra	
Tipos de calibración		Aire saturado con agua, agua saturada con aire, manual (Winkler), punto cero	
Sensores com	patibles	Polarográfico, óptico RDO	
	Intervalo	De 0 a 50 °C (de 32 a 122 °F)	
	Resolución	0,1 °C, 0,1 °F	
Temperatura	Exactitud relativa	±0,1 °C	
Temperatura	Calibración de compensación	1 punto	
	Opciones de alimentación	Automática con sensor de temperatura integrado	
Entrada del sensor	MiniDIN de 9 pines	Sensor de oxígeno disuelto con temperatura integrada	

Especificaciones del medidor de sobremesa de pH/ISE Orion Star A214			
Canales de medida		1 - pH, mV, RmV, ORP o concentración de iones con temperatura	
	Intervalo	de -2,000 a 20,000	
	Resolución	0,1 / 0,01 / 0,001	
	Exactitud relativa	±0,002	
рH	Puntos de calibración	Hasta 5	
	Edición de calibración	Sí	
	Impedancia de entrada	> 10 ^{^12} ohmios	
	Estado del electrodo	Indicador en la pantalla de estado del electrodo bueno, normal o malo	
	Intervalo	0,0001 al 19900	
	Resolución	0,0001 mínimo, de 1 a 3 dígitos significativos (seleccionable por el usuario)	
	Exactitud relativa	\pm 0,2 mV o \pm 0,05 % de la lectura, lo que sea mayor	
ISE	Unidades	ppm, molar, mg/L,%, ppb, ninguna	
(Concentración de iones)	Puntos de calibración	Hasta 5	
	Edición de calibración	Sí	
	Características avanzadas	Pendiente segmentada (punto por punto), blanco automático seleccionable no lineal, estabilidad de intervalo de concentración baja	
	Intervalo	±2000,0 mV	
	Resolución	0,1 mV	
mV/RmV	Exactitud relativa	$\pm 0,2~\text{mV}$ o $\pm 0,05~\%$ de la lectura, lo que sea mayor	
	Modo mV relativo	Sí	
	Modo Eh ORP	Sí	
	Intervalo	De -5 a 105 °C (de 22 a 221 °F)	
	Resolución	0,1 °C, 0,1 °F	
Temperatura	Exactitud relativa	±0,1 °C	
	Calibración de compensación	1 punto	
	Opciones de alimentación	Manual o automática con sonda ATC	
Entradas del	BNC	Electrodo de pH, electrodo de ORP o electrodo de ión selectivo (ISE)	
sensor	Punta	Electrodo de referencia	
	MiniDIN de 8 pines	Sonda de temperatura ATC	

Canales de medida		2 - Canal 1: pH, mV, mV relativo (RmV) u ORP con temperatura Canal 2: conductividad, salinidad, STD o resistividad con temperatura	
	Intervalo	de -2,000 a 20,000	
	Resolución	0,1 / 0,01 / 0,001	
	Exactitud relativa	±0,002	
рН	Puntos de calibración	Hasta 5	
	Edición de calibración	Sí	
	Impedancia de entrada	> 10 ^{^12} ohmios	
	Estado del electrodo	Indicador en la pantalla de estado del electrodo bueno, normal o malo	
	Intervalo	±2000,0 mV	
	Resolución	0,1 mV	
mV/RmV	Exactitud relativa	\pm 0,2 mV o \pm 0,05 % de la lectura, lo que sea mayor	
	Modo mV relativo	Sí	
	Modo Eh ORP	Sí	
	Intervalo	De 0,001 µS a 3000 mS	
	Resolución	0,001 µS mínimo, alcance automático, hasta 4 dígitos significativos	
Conducti-	Exactitud relativa	0,5 % de la lectura ±1 dígito >3 μS, 0,5 % de la lectura ±0,01 μS <3 jS	
viuau	Temp. de referencia	5 °C, 10 °C, 15 °C, 20 °C, 25 °C	
	Comp. temp.	Lineal, nLFn no lineal, nLFu no lineal, EP (USP), apagado	
	Puntos de calibración	Hasta 5 con opción de edición de calibración	
	Intervalo	De 0,06 a 80,00 psu, de 0,05 a 42,00 ppt	
Salinidad	Resolución	0,01 psu o 0,01 ppt mínimo, alcance automático	
Samuau	Exactitud relativa	0,5 % de la lectura ±1 dígito	
	Тіро	Salinidad práctica (psu) o agua marina natural (ppt)	
	Intervalo	De 0,001 a 200,0 ppm	
STD	Resolución	0,001 ppm mínimo, alcance automático, hasta 4 dígitos significativos	
	Exactitud relativa	0,5 % de la lectura ±1 dígito	
	Factor STD	Lineal (de 0,02 a 9,99)	
	Intervalo	De 2 Ω a 100 MΩ	
Resistividad	Resolución	1 Ω o 0,1 MΩ, alcance automático	
	Exactitud relativa	0,5 % de la lectura ±1 dígito	
	Intervalo	De -5 a 105 °C (de 22 a 221 °F)	
	Resolución	0,1 °C, 0,1 °F	
	Exactitud relativa	±0,1 °C	
Temperatura	Calibración de compensación	1 punto	
	Opciones de alimentación	Manual o automática con sonda ATC o sensor de temperatura integrado	
	BNC	Electrodo de pH u ORP	
Entradas del	Punta	Electrodo de referencia	
sensor	MiniDIN de 8 pines	Sensor de conductividad con temperatura integrada o sonda ATC	

Especificaciones del medidor de sobremesa de pH/RDO/OD Orion Star A216			
Canales de medida		2 - Canal 1: pH, mV, mV relativo (RmV) u ORP con temperatura Canal 2: oxígeno disuelto como % de saturación o mg/L con temperatura	
Intervalo		de -2,000 a 20,000	
	Resolución	0,1 / 0,01 / 0,001	
	Exactitud relativa	±0,002	
	Puntos de calibración	Hasta 5	
рН	Edición de calibración	Sí	
	Impedancia de entrada	> 10 ^{^12} ohmios	
	Estado del electrodo	Indicador en la pantalla de es normal o malo	stado del electrodo bueno,
	Intervalo	±2000,0 mV	
	Resolución	0,1 mV	
mV/RmV	Exactitud relativa	±0,2 mV o ±0,05 % de la lec	tura, lo que sea mayor
	Modo mV relativo	Sí	
	Modo Eh ORP	Sí	
	Polarográfico	Concentración	% de saturación
	Intervalo	De 0 a 90 mg/L	De 0 a 600 %
	Resolución	0.01 / 0.1 mg/L	0.1/1%
	Exactitud relativa	±0,2 mg/L o ±2 % de la lectura, lo que sea mayor	±2 % de saturación o ±2 % de la lectura, lo que sea mayor
	Óptico RDO	Concentración	% de saturación
	Intervalo	De 0 a 50 mg/L	De 0 a 500 %
	Resolución	0,01 / 0,1 mg/L	0,1 / 1 %
Oxígeno disuelto	Exactitud relativa	±0,1 mg/L hasta 8 mg/L, de ±0,2 mg/L 8 a 20 mg/L, ±10 % de lectura hasta 50 mg/L	± 2 % de saturación < 200 % de saturación, ± 10 % de saturación > 200 % de saturación
	Corrección de la presión barométrica	De 400 a 850 mm Hg, automática con barómetro integrado (±6 mm Hg) o introducción manual	
	Corrección de salinidad	De 0,0 a 45,0 ppt, automático con introducción manual de la salinidad de muestra	
	Tipos de calibración	Aire saturado con agua, agua saturada con aire, manual, punto cero	
	Sensores compatibles	Polarográfico, óptico RDO	
	Intervalo	De 0 a 50 °C (de 32 a 122 °F	-)
	Resolución	0,1 °C, 0,1 °F	
Temperatura	Exactitud relativa	±0,1 °C	
	Calibración de compensación	1 punto	
	Opciones de alimentación	Manual o automática con sonda ATC o sensor de temperatura integrado	
	BNC	Electrodo de pH u ORP	
Entradas del	Punta	Electrodo de referencia	
sensor	MiniDIN de 8 pines	Sonda de temperatura ATC	
	MiniDIN de 9 pines	Sensor de oxígeno disuelto con temperatura integrada	

Nota: Todas las especificaciones pueden cambiar sin previo aviso.

Información para pedidos

N.º de cat.	Descripción
STARA2110	Medidor de pH Star A211 con portaelectrodos, adaptador de alimentación universal, CD de documentación, guía de inicio rápida impresa, cable de interfaz para ordenador y certificado de prueba del medidor
STARA2115	Kit de medidor de pH Star A211 con portaelectrodos, adaptador de alimentación universal, CD de documentación, guía de inicio rápida impresa, cable de interfaz para ordenador y certificado de prueba del medidor
	 Electrodo pH/ATC con cuerpo de cristal rellenable ROSS Ultra Triode, 8302BNUMD
	Kit de solución de almacenamiento y búfer pH ROSS, 810199
	Kit de medidor de pH Star A211 con portaelectrodos, adaptador de alimentación universal, CD de documentación, guía de inicio rápida impresa, cable de interfaz para ordenador y certificado de prueba del medidor
STARAZIII6	 Electrodo pH/ATC con cuerpo de epoxi rellenable ROSS Ultra Triode, 8157BNUMD
	Kit de solución de almacenamiento y búfer pH ROSS, 810199
STARA2117	Kit de medidor de pH Star A211 con portaelectrodos, adaptador de alimentación universal, CD de documentación, guía de inicio rápida impresa, cable de interfaz para ordenador y certificado de prueba del medidor
	Electrodo pH/ATC con cuerpo de epoxi rellenable del estándar 9157BNMD Triode
	Kit de solucion de almacenamiento y buter pH, 910199
STARA2120	universal, CD de documentación, guía de inicio rápida impresa, cable de interfaz para ordenador y certificado de prueba del medidor
STARA2125	Kit de medición de conductividad Star A212 con portaelectrodos, adaptador de alimentación universal, CD de documentación, guía de inicio rápida impresa, cable de interfaz para ordenador y certificado de prueba del medidor
	 Sensor de conductividad con cuerpo de epoxi y 4 celdas DuraProbe (K=0,475) 013005MD Estécular de Osadasticidad Osias 4442 li2 044027, Esc. 60 ml.
	Estandar de Conductividad Orion 1413 JS 011007, 5 x 60 mL Kit de medición de conductividad Star A212 con partecleatrades, edentadar de
074040406	alimentación universal, CD de documentación, guía de inicio rápida impresa, cable de interfaz para ordenador y certificado de prueba del medidor
STARAZ 120	 Sensor de conductividad de 2 celdas de agua pura (K = 0,1) 013016MD con celda de flujo extraíble
	Estándar de conductividad Orion 100 jS 011008, 5 x 60 mL
STARA2130	Medidor de RDO/OD Star A213 con portaelectrodos, adaptador de alimentación universal, CD de documentación, guía de inicio rápida impresa, cable de interfaz para ordenador y certificado de prueba del medidor
	Kit de medición de RDO/OD Star A213 con portaelectrodos, adaptador de alimentación universal, CD de documentación, guía de inicio rápida impresa, cable de interfaz para ordenador y certificado de prueba del medidor
STARA2135	Sensor de OD polarográfico con manga de calibración, 083005MD
	Kit de mantenimiento de sensor OD 080513
	Adaptador BDO, embudo y agitador
	Kit de medicion de RDO/OD Star A213 con portaelectrodos, adaptador de alimentacion universal, CD de documentación, guía de inicio rápida impresa, cable de interfaz para ordenador y certificado de prueba del medidor
STARA2136	 Sensor BDO/OD polarográfico autoagitador, 086030MD
	Kit de mantenimiento de sensor OD 080513
	Solucion de electrolitos para sensor OD 080514
STARA2140	universal, CD de documentación, guía de inicio rápida impresa, cable de interfaz para ordenador y certificado de prueba del medidor

N.º de cat.	Descripción		
STARA2145	 Kit de medición de pH/ISE Star A214 con portaelectrodos, adaptador de alimentación universal, CD de documentación, guía de inicio rápida impresa, cable de interfaz para ordenador y certificado de prueba del medidor Electrodo de pH rellenable de cristal ROSS Ultra 8102BNUWP Sonda de temperatura ATC de acero inoxidable Orion 927007MD Sonda de agitador Orion Star 096019 Kit de solución de almacenamiento y búfer pH ROSS, 810199 		
STARA2146	 Kit de medición de pH/ISE Star A214 con portaelectrodos, adaptador de alimentación universal, CD de documentación, guía de inicio rápida impresa, cable de interfaz para ordenador y certificado de prueba del medidor Electrodo de pH rellenable de cristal ROSS Ultra 8102BNUWP Electrodo de ion selectivo (ISE) para amoniaco de alto rendimiento Orion 9512HPBNWP Sonda de temperatura ATC de acero inoxidable Orion 927007MD Sonda de agitador Orion Star 096019 Estándar de amoniaco de 1000 ppm Orion 951007, 475 mL Solución ISA para amoniaco de bajo nivel Orion 951210, 475 mL Solución de almacenamiento de electrodo para amoniaco Orion 951213, 475 mL 		
STARA2147	 Kit de medición de pH/ISE Star A214 con portaelectrodos, adaptador de alimentación universal, CD de documentación, guía de inicio rápida impresa, cable de interfaz para ordenador y certificado de prueba del medidor Electrodo de pH rellenable de cristal ROSS Ultra 8102BNUWP Electrodo de ion selectivo (ISE) de fluoruro Orion 9609BNWP Sonda de temperatura ATC de acero inoxidable Orion 927007MD Sonda de agitador Orion Star 096019 Estándar de fluoruro de 1 ppm con TISAB II Orion 040906, 475 mL Estándar de fluoruro de 10 ppm con TISAB II Orion 040907, 475 mL Estándar de fluoruro de 10 ppm con TISAB II Orion 040908, 475 mL Solución Orion TISAB II 940909, 1 galón 		
STARA2148	 Kit de medición de pH/ISE Star A214 con portaelectrodos, adaptador de alimentación universal, CD de documentación, guía de inicio rápida impresa, cable de interfaz para ordenador y certificado de prueba del medidor Electrodo de pH rellenable de cristal ROSS Ultra 8102BNUWP Electrodo de ion selectivo (ISE) ROSS para sodio 8611BNWP con estándares e ISA Sonda de temperatura ATC de acero inoxidable Orion 927007MD Sonda de agitador Orion Star 096019 		
STARA2150	Medidor de pH/conductividad Star A215 con portaelectrodos, adaptador de alimentación universal, CD de documentación, guía de inicio rápida impresa, cable de interfaz para ordenador y certificado de prueba del medidor		
STARA2155	 Kit de medición de pH/conductividad Star A215 con portaelectrodos, adaptador de alimentación universal, CD de documentación, guía de inicio rápida impresa, cable de interfaz para ordenador y certificado de prueba del medidor Electrodo pH/ATC con cuerpo de epoxi rellenable ROSS Ultra Triode, 8157BNUMD Sensor de conductividad con cuerpo de epoxi y 4 celdas DuraProbe (K=0,475) 013005MD Kit de solución de almacenamiento y búfer pH ROSS, 810199 Estándar de Conductividad Orion 1413 μS 011007, 5 x 60 mL 		
STARA2160	Medidor de pH/RDO/OD Star A216 con portaelectrodos, adaptador de alimentación universal, CD de documentación, guía de inicio rápida impresa, cable de interfaz para ordenador y certificado de prueba del medidor		

N.º de cat.	Descripción
STARA2165	Kit de medición de pH/RDO/OD Star A216 con portaelectrodos, adaptador de alimentación universal, CD de documentación, guía de inicio rápida impresa, cable de interfaz para ordenador y certificado de prueba del medidor
	 Electrodo pH/ATC con cuerpo de epoxi rellenable ROSS Ultra Triode, 8157BNUMD
	Sensor de OD polarográfico con manga de calibración, 083005MD
	 Kit de solución de almacenamiento y búfer pH ROSS, 810199
	Kit de mantenimiento de sensor OD 080513
	Adaptador BDO, embudo y agitador

Accesorios, electrodos y soluciones del medidor

N.º de cat.	Descripción
IQOQ-STARA	Documentación de los medidores Star A210, Star A220 y Star A320 IQ/OQ (cualificación de instalación/cualificación operativa), válida para todas las configuraciones de medidores de la lista
STARA-BEA	El portaelectrodos conectable al medidor de sobremesa Star A y Versa Star incluye brazo del electrodo, soporte y soporte para el medidor
STARA-HB	Base ponderada independiente para uso con el portaelectrodos Star A y Versa Star
810017	Manga de almacenamiento y base para los electrodos de 12 mm de diámetro
1010003	Adaptador de alimentación universal para medidores de la serie Star A
1010053	Cable de ordenador RS232 de la serie Star
1010005	Cable y adaptador de RS232 a USB de la serie Star
1010006	Impresora de inyección de tinta de la serie Star, 110 V/220 V, con cable de impresora RS232
096019	Sonda de agitador y conector en punta Orion Star
927007MD	Sonda de temperatura ATC de Orion con cuerpo de acero inoxidable, conector MiniDIN
927005MD	Sonda de temperatura ATC de Orion con cuerpo de epoxi, conector MiniDIN
928007MD	Microsonda de temperatura ATC de Orion con punta de acero inoxidable, conector MiniDIN
8102BNUWP	Electrodo de pH rellenable con cuerpo de cristal ROSS Ultra, conector BNC
8156BNUWP	Electrodo de pH rellenable con cuerpo de epoxi ROSS Ultra, conector BNC
8172BNWP	Electrodo de pH rellenable con cuerpo de cristal ROSS Sure-Flow, conector BNC
8165BNWP	Electrodo de pH rellenable con cuerpo de epoxi ROSS Sure-Flow, conector BNC
8302BNUMD	Electrodo de pH/ATC rellenable con cuerpo de cristal ROSS Ultra Triode, conectores BNC y MiniDIN
8157BNUMD	Electrodo de pH/ATC rellenable con cuerpo de epoxi ROSS Ultra Triode, conectores BNC y MiniDIN
8107BNUMD	Electrodo de pH/ATC llenado con gel con cuerpo de epoxi ROSS Ultra Triode, conectores BNC y MiniDIN
8135BNUWP	Electrodo de pH rellenable con cuerpo de epoxi ROSS Ultra con bombilla de superficie plana, conector BNC
8163BNWP	Electrodo de pH rellenable con cuerpo de cristal y punta aguda ROSS, conector BNC
8103BNUWP	Electrodo semimicro de pH rellenable con cuerpo de cristal ROSS Ultra, conector BNC
8115BNUWP	Electrodo semimicro de pH rellenable con cuerpo de epoxi ROSS Ultra, conector BNC
8220BNWP	Electrodo micro de pH rellenable con cuerpo de cristal ROSS, conector BNC
810199	Kit de solución de almacenamiento y búfer de pH ROSS All-in-One, incluye búferes de pH 4, 7, 10; solución de almacenamiento de ROSS; solución de limpieza de electrodo de pH; botella de almacenamiento de electrodo de pH
810001	Solución de almacenamiento para electrodo de pH ROSS, 475 mL
910001	Solución de almacenamiento para electrodo de pH, 475 mL
910168	Búfer de pH 1,68 Orion, 475 mL
910104	Búfer de pH 4,01 Orion, 475 mL
910105	Búfer de pH 5,00 Orion, 475 mL
910686	Búfer de pH 6,86 Orion, 475 mL
910107	Búfer de pH 7,00 Orion, 475 mL
910918	Búfer de pH 9,18 Orion, 475 mL
910110	Búfer de pH 10,01 Orion, 475 mL
910112	Búfer de pH 12,46 Orion, 475 mL

N.º de cat.	Descripción
9678BNWP	Electrodo rellenable con cuerpo de epoxi ORP Orion Sure-Flow, conector BNC
9180BNMD	Electrodo rellenable con cuerpo de epoxi ORP/ATC Orion Triode, conectores BNC y MiniDIN
967901	Solución estándar ORP Orion, 475 mL
967961	Solución estándar ORP Orion, 5 x 60 mL
9512HPBNWP	Electrodo de ion selectivo para amoniaco de alto rendimiento Orion, conector BNC
9512BNWP	Electrodo de ion selectivo para amoniaco estándar Orion, conector BNC
951007	Estándar de amoniaco de 1000 ppm Orion, 475 mL
951211	Ajustador de fuerza iónica (ISA) para amoniaco Orion con tinte de color azul indicador del pH, 475 mL
951210	ISA para amoniaco de nivel bajo Orion con tinte de color azul indicador del pH, 475 mL
9609BNWP	Electrodo de ion selectivo de fluoruro Orion, conector BNC
940907	Estándar de fluoruro de 100 ppm Orion, 475 mL
940909	Búfer de ajuste de fuerza iónica total Orion TISAB II para análisis de fluoruros, 3,8 L
040906	Premezclado estándar de fluoruro de 1 ppm Orion con TISAB II, 475 mL
040907	Premezclado estándar de fluoruro de 2 ppm Orion con TISAB II, 475 mL
040908	Premezclado estándar de fluoruro de 10 ppm Orion con TISAB II, 475 mL
9707BNWP	Electrodo de ion selectivo de nitrato Orion, conector BNC
920707	Estándar de nitrato de 1000 ppm Orion, 475 mL
930711	Ajustador de fuerza iónica (ISA) de nitrato Orion 475 mL
930710	Solución supresora de interferencia de nitratos Orion, 475 mL
8611BNWP	Electrodo de ion selectivo para sodio ROSS, conector BNC
841108	Estándar de sodio de 1000 ppm Orion, 475 mL
841111	Ajustador de fuerza iónica (ISA) de sodio Orion 475 mL
013005MD	Sensor de temperatura/conductividad Orion DuraProbe de 4 celdas (K = 0,475), conector MiniDIN
013016MD	Sensor de temperatura/conductividad Orion Pure Water de 2 celdas (K=0,1), conector MiniDIN
018020MD	Sensor de conductividad de 2 celdas de gran rango Orion (K=10), conector MiniDIN
011008	Estándar de conductividad Orion 100 jS/cm, 5 x 60 mL
011007	Estándar de conductividad Orion 1413 jS/cm, 5 x 60 mL
011006	Estándar de conductividad de 12,9 mS/cm, 5 x 60 mL
1010001	Kit de reostato de comprobación de conductividad Orion
083005MD	Sensor de OD polarográfico Orion con manga de calibración, conector MiniDIN
086030MD	Sensor de BDO polarográfico autoagitador Orion con manga de calibración, conector MiniDIN
087010MD	Sensor de OD óptico RDO con tapón óptico, manga de calibración y protección, conector MiniDIN

Visite <u>www.thermoscientific.com/water</u> para obtener una lista completa de todos los medidores, electrodos, soluciones y accesorios Thermo Scientific Orion disponibles.

thermoscientific.com/water

© 2015 Thermo Fisher Scientific Inc. Reservados todos los derechos. RDO es una marca comercial registrada de In-Situ Inc. Microsoft, Windows y Microsoft Vista son marcas comerciales registradas de Microsoft Corporation. Todas las demás marcas comerciales son propiedad de Thermo Fisher Scientific Inc. y sus filiales.

Productos de Water and Lab

Norteamérica

Tel. gratuito: 1-800-225-1480 Tel.: 1-978-232-6000 Info.water@thermofisher.com Alemania Tel.: (49) 6184-90-6000 info.water.uk@thermofisher.com

China Tel.: (86) 21-68654588 wai.asia@thermofisher.com India Tel.: (91) 22-4157-8800 wai.asia@thermofisher.com

Singapur Tel.: (65) 6778-6876 wai.asia@thermofisher.com Japón Tel.: (81) 045-453-9175 wai.asia@thermofisher.com

Australia Tel.: (613) 9757-4300 En Australia (1300) 735-295 InfoWaterAU@thermofisher.com

