

## Thermo Scientific Orion AquaMate Benutzerhandbuch

Spektrophotometer AQ7100 Vis und AQ8100 UV-Vis AQX1MAN-DE • Revision A • Juni 2022



© 2019 Thermo Fisher Scientific Inc. Alle Rechte vorbehalten.

Alle anderen Marken sind Eigentum der Thermo Fisher Scientific Inc. und ihrer Tochtergesellschaften.

Technischen Support erhalten Sie unter: www.thermofisher.com

Dieses Dokument liegt allen Produkten von Thermo Fisher Scientific Inc. beim Kauf bei und ist beim Betrieb des Produkts zu beachten. Dieses Dokument ist urheberrechtlich geschützt. Jede teilweise oder vollständige Reproduktion dieses Dokuments ist streng untersagt, sofern keine schriftliche Genehmigung von Thermo Fisher Scientific Inc. vorliegt.

Der Inhalt dieses Dokuments kann jederzeit ohne Ankündigung geändert werden. Alle in diesem Dokument enthaltenen technischen Angaben dienen lediglich zur Information. In diesem Dokument genannte Systemkonfigurationen und spezifikationen ersetzen alle dem Käufer bereits gegebenen Informationen.

Thermo Fisher Scientific Inc. erhebt keinen Anspruch auf Vollständigkeit, Genauigkeit und Fehlerfreiheit der Informationen in diesem Dokument und übernimmt keine Haftung für Fehler, Versäumnisse, Schäden oder Verluste, die aus dem Gebrauch dieses Dokuments entstehen, selbst wenn die Informationen in diesem Dokument genau befolgt werden.

Dieses Dokument ist nicht Teil eines Kaufvertrags zwischen Thermo Fisher Scientific Inc. und einem Käufer. Dieses Dokument regelt oder ändert keine Geschäftsbedingungen; bei widersprüchlichen Informationen zwischen den beiden Dokumenten gelten die Geschäftsbedingungen.

Nur für Forschungszwecke. Dieses Gerät oder Zubehör ist kein Medizinprodukt und ist nicht für die Prävention, Diagnose, Behandlung oder Heilung von Krankheiten vorgesehen.

## Inhaltsverzeichnis

| Kapitel 1  | . 7 |
|--|-----|
| Spektrophotometer – Einführung   | . 7 |
| Spektrophotometer – Übersicht  | 7   |
| Orion AquaMate 7100 Vis Spektrophotometer                              | 8   |
| Orion AquaMate 8100 UV-Vis-Spektrophotometer                           | 8   |
| Packliste  | 8   |
| Packliste für Orion AquaMate 7100 Vis Spektrophotometer                | 8   |
| Packliste für Orion AquaMate 8100 UV-Vis Spektrophotometer             | 9   |
| Orion AquaMate-Anwenderdokumentation auf USB                           | 10  |
| Bestimmungsgemäße Verwendung   | 10  |
| Vorsichtsmaßnahmen beim Betrieb  | 10  |
| Sicherheitsbezogene und besondere Hinweise                             | 11  |
| Kapitel 2  | 12  |
| Spektrophotometer – Grundlagen   | 12  |
| Komponenten des Spektrophotometers                                     | 12  |
| Touchscreen  | 13  |
| Geräteanschlüsse   | 14  |
| Optionales Zubehör   | 14  |
| Probenkammer   | 15  |
| Probenkammer für AQ7100 und AQ8100                                     | 15  |
| Einzelküvettenhalter   | 16  |
| Merkmale der Probenschale  | 16  |
| Enthehmen – den Kuvettenhalter greifen und nach oben und vorne anheben | 16  |
| Optionale Probennaiter   | 1/  |
| Kuvettennalter ersetzen  | 10  |
| Optionaler Drucker   | 19  |
| Auswahl und Positionierung von Flaschchen und Kuveiten                 | 22  |
|  | 22  |
| Kapitel 3  | 23  |
| Geräteeinrichtung, Touchscreen und Funktionen des Orion AquaMate       | 23  |
| Bildschirmnavigation   | 23  |
| Vertraute Benutzeroberfläche   | 25  |
| Inhalt der Benutzeroberflache  | 27  |
| Bildschirm 1 Startbildschirm   | 21  |
| Bildschirm 2 – Methodenentwicklung, Diagnostik und Daten               | 34  |
| Bildschirm 3 – Mentrachweilenlange und OD600                           | 44  |
|  | 45  |
|  | 45  |
| SiliariSiari   | 40  |
| nnponteren unu ⊏xponteren von versuchen<br>Datan avportieren           | 55  |
|  | 00  |
| Kapitel 4  | 5/  |
| Testmenü für Wasseranalysen  | 57  |
| Vorprogrammierte Methoden  | 57  |
| Methodenauswahl und Versuch  | 58  |

| Droplet-Ordner-Methoden  | 58   |
|--|------|
| Methoden-Optionen  | 59   |
| Probennummerierung einer Methode   | 60   |
| Laden von Testmethoden aus dem AquaMate-Gerät                                      | 61   |
| Laufende Testmethoden für die Wasseranalyse  | 62   |
| Ein-Punkt-Anpassung der Methode  | 64   |
| Verwenden der Farbumkehrfunkion  | 65   |
| Kapitel 5  | . 66 |
| Orign AOUAfact Descention Anlaitung für Orign AgusMata                             | 66   |
| Orion AQUAIast Reagenzien – Americung für Orion Aquamate                           | . 00 |
| Onon AQUAIasi Kolonimetrische Reagenzien Kompatibel mit Onon Aquaiviate-Geraten    | 00   |
| Anweisungen zum Urion AQUAtast-Keagenz   | 68   |
| Empteniungen zur Vermeidung von Messteniern  | 68   |
| AC2002 Alkalinitat-M (Alkalinitat bis pH 4,3) Tablettentest                        | 69   |
| AC3002P Alkalinitat-P (Alkalinitat bis pH 8,2) Lablettentest                       | 70   |
|  | /1   |
| AC4P27 Aluminium-Pulverpackchen- und Flussigreagenz-Test                           | 12   |
| AC2012 Ammoniak-Tabletten-Test   | /3   |
| AC4P12 Ammoniak-Pulverpackchen- und Flussigreagenz-Test                            | /4   |
| ACR012 Ammoniak-Reaktionsrohrchen-Lest für niedrigen Bereich                       | /5   |
| ACRU11 Ammoniak-Reaktionsrohrchen-Test für hohen Bereich                           | /6   |
| AC2035 Bromtabletten-Test  |      |
| AC2017 Chlorid-Tabletten-Test  | 79   |
| AC2070 Chlortabletten-Test (treies Chlor und Gesamtchlor)                          | 80   |
| AC20/1 Chlortabletten- I est (treies Chlor)  | 82   |
| AC20/2 Chlortabletten- I est (Gesamtchlor)   | 84   |
| AC4P71 Chlorpulverpäckchen-Test (freies Chlor)                                     | 86   |
| AC4P72 Chlorpulverpäckchen-Test (Gesamtchlor)                                      | 87   |
| AC3072 Chlortabletten-Test (Gesamtchlor) für hohen Bereich                         | 88   |
| AC2099 Chlordioxid-Tabletten-Test  | 89   |
| CODL00 COD-Aufschlussröhrchen-Test für niedrigen Bereich                           | 91   |
| CODH00 COD-Aufschlussröhrchen-Test für mittleren Bereich                           | 93   |
| CODHP0 COD-Aufschlussröhrchen-Test für hohen Bereich                               | 94   |
| AC2029 Kupfertabletten-Test (freies Kupfer und Gesamtkupfer)                       | 95   |
| AC4P29 Kupferpulverpäckchen-Test (freies Kupfer)                                   | 96   |
| AC2098 Cyanursäure-Tabletten-Test  | 97   |
| AC2009 SPADNS Fluorid-Flüssigreagenz-Test  | 98   |
| AC3032T Wasserhärte-Tabletten-Test (Gesamthärte)                                   | 100  |
| AC2030 Hydrazinpulver-Test   | 101  |
| AC2078 Eisentabletten-Test (Eisen II und III)                                      | 102  |
| AC4P78 Iron (Ferro) Powder Pack Test   | 103  |
| AC4P79 Eisenpulverpäckchen-Test (Gesamteisen)                                      | 104  |
| AC2055 Mangantabletten-Test  | 105  |
| AC4P54 Mangan-Pulverpäckchen- und -Flüssigreagenztest für niedrigen Bereich        | 106  |
| AC4P55 Mangan-Pulverpäckchen-Test für hohen Bereich                                | 107  |
| AC4P42 Molybdat-Pulverpäckchen   | 108  |
| ACR007 Nitrat-Reaktionsröhrchen-Test   | 109  |
| AC2046 Nitrittabletten-Test  | 110  |
| AC4P46 Nitritpulverpäckchen-Test   | 111  |
| ACD004 Stickstoff-Aufschlussröhrchen-Test (Gesamtstickstoff) für niedrigen Bereich | 112  |
| ACD007 Stickstoff-Aufschlussröhrchen-Test (Gesamtstickstoff) für hohen Bereich     | 114  |

| AC3048 Ozontabletten-Test   | 116                |
|---|--------------------|
| AC2001 pH-Tablettentest   | 118                |
| AC3001 Flüssigreagenz-pH-Test   | 119                |
| AC2095-WA Phosphattabletten-Test (Ortho) für niedrigen Bereich  | 120                |
| AC2096 Phosphattabletten-Test (Ortho) für hohen Bereich   | 121                |
| AC4P95 Phosphat-Pulverpäckchen-Test (Ortho)   | 122                |
| ACR095 Phosphat-Reaktionsröhrchen-Test (Ortho)  | 123                |
| ACD095 Phosphat-Aufschlussröhrchen-Test (Gesamtphosphat)  | 124                |
| ACD095AH Phosphat-Aufschlussröhrchen-Test (säurehydrolysierbar)   |                    |
| AC2060 Siliziumdioxid-Tabletten-Test  | 128                |
| AC2061 Siliziumdioxid mit Phosphateliminierungs-Tabletten-Test  |                    |
| AC4P60 Siliziumdioxid-Pulverpäckchen-Test   |                    |
| AC4P82 Sulfat-Pulverpäckchen-Test   |                    |
| AC2016 Sulfid-Tabletten-Test  |                    |
| AC2065 Zinktabletten-Test   |                    |
| Farbmessung aus "Application Log" (Anwendungsprotokoll) Nr. 131<br>UVA- und UV254-Farbmessungen aus "Application Log" (Anwendungsprotokoll) |                    |
| Nr. 137   |                    |
| Kapitel 6   | 142                |
| Testmenü "Standard Curve" (Standardkurve)   | 142                |
| Konzentrationsmessungen mit der Quant Standardkurven-Anwendung (benutzerdefinie   | erte               |
| Methode)  |                    |
| Aufrufen der Quant-Anwendung  | 143                |
| Optionen für Standardkurve in Quant   | 144                |
| Einrichten der Parameter für eine Standardkurve   | 144                |
| Erstellen einer Standardkurve in Quant  | 145                |
| Messstandards für eine Standardkurve  | 145                |
| Probenmessung mit Quant   | 146                |
| Versuch beenden   | 147                |
| Daten exportieren   | 148                |
| Bearbeiten einer Standardkurve  | 149                |
| Kapitel 7   | 152                |
| ,<br>Menü für den Wellenlängen-Scan-Test  | 152                |
| Finrichten der Darameter für einen Scan   | 1 <b>52</b><br>153 |
| Erfassen eines Basislinienscans und Scannen einer Prohe   |                    |
| Berechnungen mit den Scan-Daten durchführen   | 156                |
| Die Funktion 3Pt Net" (3-Pkt Netz)  | 158                |
| Funktion zur Flächenberechnung  | 159                |
| Aufrufen einer bestehenden Scan-Methode   |                    |
| Kanital 8   | 161                |
| Rapiter 6   | 101                |
| Integrierte Aquawate-Sontware   | 101                |
| Kapitel 9   | 170                |
| Messungen der Absorption, des prozentualen Transmissionsgrads und der   |                    |
| Konzentration   | 170                |
| Messungen der Absorption und prozentualen Transmission  | 170                |
| Anwendung mit fester Wellenlänge ("Fixed") für die Testmethode "Basic A-%T-C  | ' 170              |
| Verfügbare feste Gleichungen  |                    |
| Verwendung des "C-Mode" (K-Modus) zur Konzentrationsmessung   | 173                |

| Mehrfachwellenlänge                               | 174 |
|---|-----|
| 3-Punkt-Netz                                      | 175 |
| Kinetik   | 176 |
| Kapitel 10  | 177 |
| Wartung/Pflege                                    | 177 |
| Regelmäßige Pflege                                | 178 |
| Reinigung und Wartung von Fläschchen und Küvetten | 178 |
| Reinigen der Probenkammerfenster                  |     |
| Auswechseln der Wolfram-Halogenlampe              |     |
| Betriebsdauer der Xenon-Lampe                     |     |
| Auswechseln des Xenon-Blitzlichts                 |     |
| Kapitel 11  | 186 |
| Kundendienst                                      |     |
| Technischer Support                               |     |
| Gerätespezifikationen                             |     |
| Bestellinformationen                              |     |
| Anhang A  | 192 |
| Allgemeine Geräteinformationen                    |     |
| Parameter   |     |
| Berechnungen für die Software                     |     |

## **KAPITEL 1** Spektrophotometer – Einführung

## Spektrophotometer – Übersicht

Die Spektrophotometer Orion™ AquaMate™ Vis- und UV-Vis von Thermo Scientific™ bieten folgende Funktionen und Vorteile:

- Einfache Bedienung mit mehr als 260 vorprogrammierten Methoden für gängige kolorimetrische Reagenzien
- Einfacher Zugang zu behördlich anerkannten Analysemethoden für Abwasser- und Trinkwasserproben.
- Secured Smart Method: Gesicherte Methodenauswahl für häufig verwendete Methoden
- Mit Handschuhen bedienbare Touchscreen-Oberfläche.
- Flexibilität bei der Erstellung neuer Methoden für zusätzliche Reagenzien oder Proben Erstellung neuer Methoden unter Verwendung von Kalibrierstandards oder Aktualisierung von Methoden mithilfe öffentlich zugänglicher Wellenlängen und Gleichungen
- Kompatibel mit einer Vielzahl von runden und rechteckigen Fläschchen und Probenhaltersystemen.
- Tests zur Leistungsüberprüfung stellen die Wellenlängengenauigkeit und die Funktionalität des Geräts sicher, und integrierte Filter ermöglichen die Wellenlängenüberprüfung ohne zusätzliche Geräte
- Zusätzliche Funktionen umfassen Standardkurven-Konzentrationsmessungen, Wellenlängen-Scan, Messungen mit mehreren festen Wellenlängen, Absorptionsverhältnis und Differenz
- Auf das Gerät wird ein Jahr Garantie gewährt

## Orion AquaMate 7100 Vis Spektrophotometer

Das Orion AquaMate 7100 Vis Spektrophotometer misst im Wellenlängenbereich von 325 bis 1.100 nm mit einer Wolfram-Halogen-Lampe, die für einen einfachen Austausch mit der werkseitig vorjustierten Lampe und dem Sockel konzipiert ist. Die Wolfram-Halogen-Lampe hat eine durchschnittliche erwartete Betriebsdauer von mehr als 1.000 Stunden.

## Orion AquaMate 8100 UV-Vis-Spektrophotometer

Das Orion AquaMate 8100 UV-Vis-Spektrophotometer misst im Wellenlängenbereich von 190 nm bis 1.100 nm mit einem Xenon-Blitzlicht, das keine Aufwärmzeit benötigt und für eine durchschnittliche Betriebsdauer von 3–5 Jahren ausgelegt ist.

## Packliste

#### Packliste für Orion AquaMate 7100 Vis Spektrophotometer

- Vis Spektrophotometer mit 7-Zoll-Farb-Touchscreen und Wolfram-Halogen-Lampe
- 12-25 mm Rundfläschchenhalter (Art.-Nr.: AQX1LWLVH)
- 24 mm Rundfläschchen, 12er-Packung (Art.-Nr.: AC2V24)
- Externes AC/DC-Universal-Netzteil, 100–240 Volt, 50–60 Hz (AQX1PWRSUP)
- Standard-Netzkabel-Bundle (Nordamerika, EU und VK) mit AQ7100
  - Kabelversion NA (Art.-Nr.: AQX1NACBL)
  - EU-Kabel (Art.-Nr.: AQX1EUCBL)
  - Kabelversion UK (Art.-Nr.: AQX1UKCBL)
- Optionales APAC-Netzkabel-Bundle (China, Australien und Indien) mit AQ7100APAC
  - Kabelversion China (Art.-Nr.: AQX1CNCBL)
  - Kabelversion Australien (Art.-Nr.: AQX1AUCBL)
  - Kabelversion Indien (Art.-Nr.: AQX1INCBL)
- AquaMate-Benutzerhandbuch, Methodenliste und Reagenzienanweisungen auf USB (Art.-Nr.: AQX1MAN)
- AquaMate "Erste Schritte"
- Wichtige Warnhinweise (mehrsprachig)
- AquaMate Hinweise zu Aufstellungsort und Sicherheit
- CE-Konformitätserklärung
- Prüfbericht als Druckversion
- Staubabdeckung
- USB-Kabel

#### Packliste für Orion AquaMate 8100 UV-Vis Spektrophotometer

- UV-Vis-Spektrophotometer mit 7-Zoll-Farb-Touchscreen und Xenon-Blitzlicht
- 12–25 mm Rundfläschchenhalter (Art.-Nr.: AQX1LWLVH)
- 24 mm Rundfläschchen, 12er-Packung (Art.-Nr.: AC2V24)
- Externes AC/DC-Universal-Netzteil, 100–240 Volt, 50–60 Hz (AQX1PWRSUP)
- Standard-Netzkabel-Bundle (Nordamerika, EU und VK) mit AQ7100
  - Kabelversion NA (Art.-Nr.: AQX1NACBL)
  - Kabelversion EU (Art.-Nr.: AQX1EUCBL)
  - Kabelversion UK (Art.-Nr.: AQX1UKCBL)
- Optionales APAC-Netzkabel-Bundle (China, Australien und Indien) mit AQ7100APAC
  - Kabelversion China (Art.-Nr.: AQX1CNCBL)
  - Kabelversion Australien (Art.-Nr.: AQX1AUCBL)
    - Kabelversion Indien (Art.-Nr.: AQX1INCBL)
- AquaMate-Benutzerhandbuch, Methodenliste und Reagenzienanweisungen auf USB (Art.-Nr.: AQX1MAN)
- AquaMate "Erste Schritte"

0

- Wichtige Warnhinweise (mehrsprachig)
- AquaMate Hinweise zu Aufstellungsort und Sicherheit
- CE-Konformitätserklärung
- Prüfbericht als Druckversion
- Staubabdeckung
- USB-Kabel

**Hinweis:** Das APAC-Netzkabel-Bundle (China, Australien und Indien) muss bei der Bestellung entweder mit AQ7100APAC oder AQ8100APAC angegeben werden.

## Orion AquaMate-Anwenderdokumentation auf USB

Die Orion AquaMate-Anwenderdokumentation auf dem USB-Stick enthält:

- AquaMate-Benutzerhandbuch, Methodenliste und Reagenzienanweisungen auf USB (Art.-Nr.: AQX1MAN)
  - o AquaMate Hinweise zu Aufstellungsort und Sicherheit
  - Wichtige Warnhinweise (mehrsprachig)
  - o AquaMate-Benutzerhandbuch, Methodenliste und Reagenzienanweisungen auf USB (AQX1MSN)
  - o AquaMate "Erste Schritte"
  - o Garantieinformationen
  - o Informationen zur WEEE/RoHS-Konformität
  - o Produktionsprüfbericht
  - o Freigegebene Firmware und Methodenbibliotheken für Wasseranalysen

### Bestimmungsgemäße Verwendung

Bitte lesen Sie dieses Benutzerhandbuch sorgfältig durch. Die Verwendung außerhalb der in dieser Anleitung angegebenen Bestimmung kann zum Erlöschen der Garantie und zu dauerhaften Schäden am Gerät führen.

## Vorsichtsmaßnahmen beim Betrieb

**Warnung:** Betreiben Sie dieses System nur unter Beachtung der Sicherheitshinweise in diesem Handbuch und der im Lieferumfang des Systems enthaltenen Dokumentation.

Das Spektrophotometer enthält präzise optische Komponenten. Gehen Sie vorsichtig mit dem Gerät um und beachten Sie die folgenden Vorsichtshinweise.

• Lassen Sie das Gerät nach dem Auspacken auf Raumtemperatur kommen, bevor Sie es einschalten.

- Lassen Sie keine Feuchtigkeit ins Innere des Geräts eindringen
- Verschüttete Chemikalien sofort abwischen
- Das Gerät nicht fallen lassen
- · Das Gerät vor mechanische Stoßbelastung schützen
- · Das Gerät vor Staub schützen

### Sicherheitsbezogene und besondere Hinweise

Befolgen Sie alle in diesem Benutzerhandbuch aufgeführten Vorsichtsmaßnahmen. Sicherheitshinweise und spezielle Hinweise sind eingerahmt in Kästchen dargestellt.

Sicherheitshinweise und spezielle Hinweise umfassen folgende Punkte:

Hinweis: Enthält hilfreiche Zusatzinformationen

Wichtig: Anweisungen, die zu befolgen sind, um Schäden an der Systemhardware oder Datenverlust zu vermeiden

**Vorsicht:** Kennzeichnet eine gefährliche Situation, die bei Nichtvermeidung zu leichten oder mittelschweren Verletzungen führen könnte

**Warnung**: Weist auf eine Gefahrensituation hin, die bei Nichtvermeidung zu tödlichen oder schweren Verletzungen führen könnte

Dieses Dokument liegt allen Produkten von Thermo Fisher Scientific beim Kauf bei und ist beim Betrieb des Produkts zu beachten. Der Inhalt dieses Dokuments kann jederzeit ohne Ankündigung geändert werden. Sämtliche technischen Informationen in diesem Dokument dienen lediglich Referenzzwecken. In diesem Dokument genannte Systemkonfigurationen und spezifikationen ersetzen alle dem Käufer bereits gegebenen Informationen.

Thermo Fisher Scientific erhebt keinen Anspruch auf Vollständigkeit, Genauigkeit und Fehlerfreiheit der Informationen in diesem Dokument und übernimmt keine Haftung für Fehler, Versäumnisse, Schäden oder Verluste, die aus dem Gebrauch dieses Dokuments entstehen, selbst wenn die Informationen in diesem Dokument genau befolgt werden.

Dieses Dokument ist nicht Teil eines Kaufvertrags zwischen Thermo Fisher Scientific Inc. und einem Käufer. Dieses Dokument regelt oder ändert keine Geschäftsbedingungen; bei widersprüchlichen Informationen zwischen den beiden Dokumenten gelten die Geschäftsbedingungen.

Nur für Forschungszwecke. Dieses Gerät ist kein Medizinprodukt und ist nicht für die Prävention, Diagnose, Behandlung oder Heilung von Krankheiten vorgesehen.



Warnung: Explosions- und Brandgefahren sind zu vermeiden. Diese Gerät ist nicht für den Gebrauch in explosionsfähigen Bereichen vorgesehen.



## KAPITEL 2 Spektrophotometer – Grundlagen

## Komponenten des Spektrophotometers

Die nachstehende Abbildung zeigt einige an der Außenseite des Geräts sichtbare Komponenten:



#### Touchscreen

Der Touchscreen des Geräts verfügt über 3 Bildschirmansichten zur Navigation. Diese können durch Wischen von links nach rechts aufgerufen werden. Unten sehen Sie die drei Bildschirmansichten der Benutzeroberfläche, durch die Sie mit einer Wischbewegung nach links oder rechts navigieren können. Die Bildschirmansichten enthalten eine Reihe von Anwendungssymbolen, die im Folgenden kurz beschrieben werden. Nach der Auswahl einer Anwendung wird der Benutzer zum "Application Home" (Anwendungsstartbildschirm) geleitet. Hier können Methoden ausgewählt und angepasst und neue Methoden erstellt werden. Diagnose- und Versuchdaten können überprüft werden. Wenn Sie auf das Zahnradsymbol in der oberen rechten Ecke tippen, können Sie von jedem Bildschirm aus auf das Menü "Settings" (Einstellungen) zugreifen.



#### Geräteanschlüsse

#### Elektrische Anschlüsse



- · Ein/Aus-Kippschalter
- 12V DC hier wird das Kabel des Netzteils angeschlossen
- Zubehöranschluss reserviert für künftiges, optionales Zubehör
- USB-A-Anschlüsse siehe Abschnitt "Optionales Zubehör"
- Netzwerk/Ethernet-Anschluss: Verbinden Sie diesen Anschluss mithilfe eines Standard-Ethernetkabels (RJ45-RJ45) mit einem Netzwerkanschluss des lokalen Gebäudenetzwerks
- Single-USB-A unterstützt Flashspeichergeräte zur Methoden- und Datenspeicherung
- Duplex-USB-A unterstützt den Anschluss an einen Windows-Computer mit optionaler Fernsteuerungssoftware, Tastatur und Maus
- Exportieren von Daten an ein Netzwerk oder einen PC über Ethernet oder Wi-Fi-USB-Adapter (nicht abgebildet)
- Drucken über USB, Ethernet oder Wi-Fi-USB-Adapter (abgebildet)

**Warnung:** Stromschlaggefahren sind zu vermeiden. Schalten Sie das Gerät immer aus und ziehen Sie den Netzstecker aus der Steckdose, bevor Sie das Netzkabel aus dem Geräteanschluss ziehen.

#### **Optionales Zubehör**

Die USB-Anschlüsse unterstützen folgende Peripheriegeräte:

Drucker

## Probenkammer

Entfernen Sie alle Klebebänder von der Außenseite des Geräts und aus dem Inneren der Probenkammer.

### Probenkammer für AQ7100 und AQ8100





Hochfeste Scharniere fixieren den Deckel in jedem Winkel mit einem konstantem Drehmoment Magnet an der Vorderseite verschließt den Deckel lichtundurchlässig, wenn die Tür abgesenkt wird



Fenster und Linse schützen die interne Optik vor verschütteten Flüssigkeiten und Dämpfen



Die Probenhalterschale wird auf die Zapfen der Grundplatte ausgerichtet Überschüssiges Wasser läuft auf die Arbeitsfläche ab

#### Einzelküvettenhalter



Standardmäßiger 10 mm-Küvettenhalter Unterseite eines Einzelküvettenhalters

#### Merkmale der Probenschale

- Zur Aufnahme von Verschüttungen bis zu 150 ml
- Kann durch Hochziehen des Küvettenhalters entnommen werden
- Kann im Waschbecken oder in einer Spülmaschine gereinigt werden sofort trocknen!

**HINWEIS:** Reinigen Sie die Probenschale mit Wasser und einem milden Reinigungsmittel. Ethanol und Iso-Propyl-Alkohol können bei Bedarf verwendet werden, aber tränken Sie die Probenschale nicht in Alkohole. Achten Sie darauf, dass Aceton, Chlorkohlenwasserstoffe oder andere aggressive organische Lösungsmittel nicht mit der Probenschale in Berührung kommen. Der PC-ABS-Kunststoff erweicht und verfärbt sich dadurch.

## Entnehmen – den Küvettenhalter greifen und nach oben und vorne anheben



Einsetzen – den vorderen Magneten einrasten lassen. Senken Sie den Küvettenhalter ab, sodass der hintere Magnet geführt wird und einrastet

#### **Optionale Probenhalter**

Probenhalterschalen für andere Arten von Küvetten und Proben sind ebenfalls erhältlich. Das Einsetzen und Herausnehmen erfolgt auf die gleiche Weise wie beim Standard-Küvettenhalter.

## Reagenzglashalter



Halter für rechteckige Küvetten für große Schichtdicken

Filterhalter









#### Küvettenhalter ersetzen

Der Küvettenhalter und der verstellbare Filterhalter werden als Zubehörteile ohne Probenschale geliefert.



Lösen Sie die verliersichere Schraube und die Unterseite des Küvettenhalters, um diesen zu entnehmen. Montieren Sie den neuen Probenhalter auf die gleiche Weise.

#### **Optionaler Drucker**

Wenn das Gerät mit einem optionalen Drucker ausgestattet ist, gehen Sie nach der folgenden Anweisung vor und lesen Sie dann Kapitel 11 zur Druckereinrichtung über den Touchscreen.

- 1. Abdeckung vom Drucker abnehmen.
  - a. Verwenden Sie den Fingergriff
  - b. Ziehen Sie den Griff zu sich und heben Sie ihn an.



2. Legen Sie Druckerpapier in den optionalen Drucker ein.



3. Setzen Sie den Drucker von der Rückseite des AquaMate-Spektrophotometers aus in das Gerät ein



4. Richten Sie die Führungsschiene des Druckers an der Unterseite des Druckers auf die Führungsschiene des AquaMate-Spektrophotometers aus



#### Unterseite des Druckers

5. Schieben Sie den Drucker nach vorne, bis die Stecker vollständig eingerastet sind. Wenn die Stecker richtig eingerastet sind, hören Sie ein Klicken.





Nach vorne schieben

Der Drucker ist vollständig eingerastet

#### Auswahl und Positionierung von Fläschchen und Küvetten

Der kompatible Wellenlängenbereich für die verschiedenen Arten von Fläschchen und Küvetten hängt von dem verwendeten Material ab. Die Weglänge von Reagenzgläsern ist nicht so gut definiert wie die von Rechteck-Küvetten.

| Fläschchen-/Küvettentyp | Wellenlängenbereich   |
|-------------------------|-----------------------|
| Optisches Glas          | 360 nm bis > 1.100 nm |
| Borosilikatglas         | 330 nm bis > 1.100 nm |
| Quarz                   | 190 nm bis > 1.100 nm |
| Einwegprodukt:          |                       |
| Polystyrol              | > 340 nm              |
| Methacrylat             | > 300 nm              |
| Acryl                   | > 280 nm              |
| UV-durchlässig          | > 220 nm              |

**Hinweis:** Beachten Sie die Herstellerangaben und arbeiten Sie innerhalb des empfohlenen Bereichs.

Richten Sie die Fläschchen und Küvetten so aus, dass die durchsichtigen Seiten dem Lichtstrahl zugewandt sind, wobei eine durchsichtige Seite zur Vorderseite des Geräts und die andere zur Rückseite zeigt.

**Hinweis:** Stellen Sie die Fläschchen immer in der gleichen Ausrichtung in den Strahlengang des Geräts. Eine Ausrichtungsmarkierung auf dem Fläschchen unterstützt Sie bei der gleichmäßigen und korrekten Ausrichtung.

Bei Verwendung von Küvetten mit kleiner Öffnung (kleinem Volumen):

- Stets Küvetten mit schwarzer Maskierung verwenden
- Stets die gleiche Küvette für Blindwert- und Probenmessung verwenden

#### Z-Maß

Die nachstehende Abbildung veranschaulicht die Position des Lichtstrahls im Gerät. Die Spezifikationen für die Strahlengröße sind unten aufgeführt.

- Abstand vom Boden des Fläschchens/der Küvette zur Mitte des Strahls (Z-Maß): 8,5 mm
- Abmessungen des Strahlengangs: 2 mm (Breite) x 7 mm (Höhe)





# **KAPITEL 3 Geräteeinrichtung, Touchscreen und Funktionen des Orion AquaMate**

## Bildschirmnavigation

Unten sehen Sie die drei Bildschirmansichten der Benutzeroberfläche, durch die Sie mit einer Wischbewegung nach links oder rechts zu den Bildschirmen 1, 2 und 3 navigieren können. Die Bildschirmansichten enthalten eine Reihe von Anwendungssymbolen, die im Folgenden kurz beschrieben werden.

Nach der Auswahl einer Anwendung wird der Benutzer zum "Application Home" (Anwendungsstartbildschirm) geleitet. Hier können Methoden ausgewählt und angepasst und neue Methoden erstellt werden. Diagnose- und Versuchsdaten können überprüft werden, und Berichte zur Leistungsüberprüfung können angezeigt, erstellt und exportiert werden.

|  | SCREEN 1  |  |   | SCREEN 2  |  |                  | SCREEN 3           |         |
|--|---|--|---|---|--|------------------|--------------------|---------|
|  | AQ8100  | \$   |   | AQ8100  | \$   |                  | AQ8100             | \$      |
|  |   |  |   | m   | <b>.</b> →   |                  | 4                  |         |
| Millpore<br>Version 1.0                | Aquamate Orion<br>Version 1.0                                 | CHEMetrics<br>Version 1.0                                      | C-Mode  |   | Fixed  | Multi-Wavelength |                    |         |
|  |   |  | Ō   | $\odot$   |  |                  |                    |         |
| Drinking water<br>methods              | Wastewater<br>methods   | Scan   | Kinetics  | Diagnostics   | Performance<br>Verification Report   |                  |                    |         |
|  |   |  | €_  |   |  |                  |                    |         |
| Live Display                           |   |  | Data Viewer   |   |  |                  |                    |         |
|  | •••   |  |   | •••   |  |                  | •••                |         |
| METHODS: .<br>SCAN: ABS of Live Displa | AquaMate water i<br>or %T, min, max, ir<br>Y: ABS or %T at sp | method libraries<br>iterval, and speed.<br>ecified wavelength. | C-MODE: 1-<br>QUANT: Mu<br>FIXED: Dual<br>KINETICS - 1<br>DIAGNOSTI<br>PERFORMA | -point cal. @ sele<br>ulti-point cal. cur<br>I - λ and factor n<br>time based expen<br>CS: review verifie<br>NCE VERIFICATIO<br>ΈR: Logged expe | ect $\lambda$ and units.<br>rve development.<br>neasurements.<br>riments<br>cation test reports.<br>DN REPORTS<br>rimental data. | MULTI-WAVI       | ELENGTH: multi-λ s | canning |

#### Nach links oder rechts wischen, um zwischen den Bildschirmen zu wechseln

#### Vertraute Benutzeroberfläche

Die Benutzeroberfläche ist ein Touchscreen und die Funktionalität ähnelt sehr denen eines Smart-Tablets. Aktive blaue Berührungspunkte sind Bereiche, in denen eine Auswahl und Bearbeitungen vorgenommen werden können.

Wenn Sie beispielsweise, wie in der nachstehenden Abbildung gezeigt, eine SCAN-Anwendung verwenden und auf den Namen der Methode tippen, wird die Tastatur zum Bearbeiten angezeigt. Wenn Sie auf die minimale oder maximale Wellenlänge tippen, erscheint ein Tastenfeld zur Auswahl der Wellenlängen. Wenn Sie auf das Intervall- oder das Geschwindigkeitsfeld tippen, erscheinen entsprechende Popup-Tastenfelder zur Bearbeitung der Methode/des Versuchs. Schließlich können Sie mithilfe des Disketten-Symbols die Methode mit den bearbeiteten Feldern und dem Namen speichern. Andere Bereiche, auf die Sie achten sollten, sind die Ellipsen-Symbole, die einen erweiterten Zugriff auf Bearbeitungsfunktionen und mehr ermöglichen.



In diesem Abschnitt dient die Anwendung "Scan" als Beispiel. Wie gezeigt, werden durch Verwendung der Ellipse zusätzliche Felder für die gespeicherte Methode geöffnet, die folgende Funktionen enthalten:

- Überarbeitung der Methode
- Auswahl einer intelligenten Methode f
  ür den Modus "Smart Method" (Intelligente Methode).
- Methode exportieren
- Vervielfältigen der Methode (möglicherweise, um die ursprüngliche Methode intakt beizubehalten)
- Methodenlöschung (nur für Methoden der Nicht-Tropfen-Bibliothek)



Siehe Methoden importieren

## Inhalt der Benutzeroberfläche

#### **Bildschirm 1 Startbildschirm**

Bildschirm 1 wird nach dem Verschwinden der Thermo Scientific Begrüßungsanzeige beim Einschalten angezeigt. Die Touchscreen-Benutzeroberfläche basiert auf der Smart-Device-Technologie. Eine Anwendung wird durch Antippen des entsprechenden Symbols ausgewählt. Auf Bildschirm 1 befinden sich folgende Anwendungen:

- DROPLETS (TRÖPFCHEN) Fünf (5) Tröpfchen-Ordner mit entsprechenden Wasseranalysemethoden. Jede Methode, für die eine behördliche Genehmigung vorliegt, wird in den Ordnern "Drinking water" (Trinkwasser) oder "Wastewater droplet" (Abwassertröpfchen) dupliziert. Die Versionsnummern der Droplet-Ordner sind angegeben.
- SCAN ist ein Multi-Wellenlängen-Scan, der die Absorption (ABS) oder die prozentuale Transmission (%T) über einen justierbaren Wellenlängenbereich und eine justierbare Geschwindigkeit und Intervallauflösung auswählt.
- Live Display (Live-Anzeige) ist eine interaktive Anwendung f
  ür kontinuierliches Scannen in Echtzeit; es werden keine echten Daten gespeichert. "ABS" oder "%T" werden in Echtzeit angezeigt.



#### **Droplet-Ordner**

Es gibt drei Droplet-Ordner für die Haupt-Wasseranalysemethoden und zwei Ordner für "Regulatory" (behördlich genehmigte Analysen):

- 1. AquaMate Orion
- 2. Merck/Millipore
- 3. Chemetrics
- 4. Abwasser (behördlich genehmigte Analysen)
- 5. Trinkwasser (behördlich genehmigte Analysen)

Jede AquaMate-Wasseranalysemethode, die als behördlich genehmigte Abwasser- (AW) oder Trinkwassermethode (TW) qualifiziert ist, wird innerhalb der jeweiligen WW- oder DW-Ordner dupliziert. Innerhalb der einzelnen Ordner können Sie entweder nach dem numerischen Methodennamen oder nach dem Parameter suchen. Sie können zum Beispiel nach "AC2002" oder nach "Alkalinity-M" suchen. Sie können auch nach Methodenname, Erstellungsdatum, Datum der letzten Aktualisierung oder Datum der letzten Verwendung sortieren.

HINWEIS: Jede Methodenbeschreibung enthält die Wellenlänge, die geeignete Probengefäßgröße und den Messbereich. Wenn die falsche Probengefäßgröße verwendet wird, sind die Ergebnisse nicht genau.

|                           | AQ8100                        | \$                        |                        | Water Home  |                         | Water Home   |   |
|---------------------------|-------------------------------|---------------------------|------------------------|-------------|-------------------------|--|---|
|                           |                               |                           | 🔍 Search Method        |             |                         | <b>Q</b> Search Method   |   |
|                           |                               |                           |                        |             | Name 🔸                  |  | Name 🔸                                    |
|                           |                               |                           | CODL00, COD(low rang   | ge)         | 06-Mar-2019<br>06:57 PM | CODL00, COD(low range)   | 06-Mar-2019<br>07:15 PM                   |
| Millpore<br>Version 1.0   | Aquamate Orion<br>Version 1.0 | CHEMetrics<br>Version 1.0 | CODHP0, COD(High Ra    | ange)       | 06-Mar-2019<br>07:08 PM | CODH00, COD(Mid Range)   | 06-Mar-2019<br>07:15 PM                   |
|                           |                               |                           | CODH00, COD(Mid Rar    | nge)        | 06-Mar-2019<br>07:08 PM | ACR095, Phosphate, Ortho   | 05-Mar-2019<br>07:15 PM                   |
| <b>A</b>                  | <b>A</b>                      |                           | ACR095, Phosphate, O   | ortho       | 06-Mar-2019<br>07:08 PM | ACR012, Ammonia as Nitrogen(N)   | 06-Mar-2019<br>07:15 PM                   |
|                           | <b>U</b>                      |                           | ACR012, Ammonia as     | Nitrogen(N) | 06-Mar-2019<br>07:08 PM | ACD007, Nitrogen   | :   |
| Drinking water<br>methods | Wastewater<br>methods         | Scan                      | 100011 1               |             | 06-Mar-2019             | Method Description:  |   |
|                           |                               |                           | Choose Sorting Options |             |                         | Wavelength: 410nm  |   |
|                           |                               |                           | Method name            |             |                         | EPA method for Wastewater AnalysisRe<br>ACD007Vial Size: 16 mm ØMeasuremen   | agent:<br>nt Range:                       |
|                           |                               |                           | Date created           |             |                         | 5-150 mg/l NEPA-Approved Reference n<br>Approved Total Nitrogen Method NotPro<br>WWNitrogen, Total High Range, Persulf | nethod:<br>omulgated for<br>ate Digestion |
| Live Display              |                               |                           | Date last updated      |             |                         | Method, Powder & Digestion Tube  | ato Bigcotion-                            |
|                           |                               |                           |                        |             |                         |  |   |
|                           | •••                           |                           | Date last used         |             |                         | Continue   |   |
|                           |                               |                           | 4                      | 0 🗆         |                         | 4 0 🗆  |   |

**Hinweis:** Die Messmöglichkeiten für jede Methode entnehmen Sie bitte der jeweiligen Methodenbeschreibung. Das Gerät meldet Werte, die außerhalb des angegebenen Messbereichs liegen und für den speziellen Zweck des Anwenders oder behördliche Berichtsanforderungen möglicherweise nicht akzeptabel sind.

#### Wassertropfen-Methoden - Ein-Punkt-Anpassung

Wasseranalysemethoden können mit einer Ein-Punkt-Kalibrierung angepasst werden. Wählen Sie im folgenden Beispiel eine Methode aus und wählen Sie vor "Blank" (Nullwert) und "Measure" (Messung) die Option "Re-Calibrate" (Neu kalibrieren), um eine Ein-Punkt-Anpassung auf der Grundlage des vom Anwender eingegebenen Standardkonzentration-Werts durchzuführen. Dadurch wird das Feld "Calibration Factor" (Kalibrierungsfaktor) aktualisiert. Dieses Verfahren wird bei jeder Verwendung einer neuen °Reagenziencharge empfohlen, um chargenspezifische Schwankungen in der Reagenz-Zusammensetzung und andere Faktoren zu berücksichtigen, welche die Genauigkeit einer Methode mit einer festen Kalibrierungskurve beeinflussen





#### Wasseranalysemethoden mit Ausgabe in mehreren Maßeinheiten

Es gibt einige Wasseranalysemethoden, bei denen die Ergebnisse gleichzeitig in mehreren Maßeinheiten ausgeben werden können. Im folgenden Beispiel wird dargestellt, wie die Ergebnisse in drei Maßeinheiten angegeben werden können.



#### Behördlich anerkannte Wasseranalysemethoden

In den Methodenbibliotheken "Drinking Water" (Trinkwasser) und "Wastewater" (Abwasser) kann der Benutzer bei der Auswahl einer Methode die Methodenbeschreibung lesen. Darin ist auch eine Beschreibung der behördlich genehmigten Methode enthalten.

HINWEIS: Der behördlich genehmigte Methodenbereich kann auf einen anderen Bereich als den in AquaMate-Methode angegebenen Bereich beschränkt sein.

|  | Water Ho  | ome  |   |              |
|--|---|--|---|--------------|
| ${\sf Q}$ Search Method  |   |  |   |              |
|  |   |  |   | $\checkmark$ |
| CODL00, COD(low range)   |   |  |   |              |
| CODH00, COD(Mid Range  | )   |  |   |              |
| ACR095, Phosphate, Orth  | 0   |  |   |              |
| ACR012, Ammonia as Niti  | ogen(N)   |  |   |              |
| ACD007, Nitrogen   |   |  |   | •<br>•       |
| Method Description:  |   |  |   |              |
| Wavelength: 410nm<br>EPA method for Wastew<br>ACD007Vial Size: 16 mm<br>5-150 mg/l NEPA-Approv<br>Approved Total Nitrogen<br>WWNitrogen, Total, High<br>Method, Powder & Diges | ater Analy<br>n ØMeasur<br>ved Refere<br>n Method N<br>Range, Po<br>tion Tube | vsisReage<br>rement Ra<br>ence meth<br>NotPromu<br>ersulfate I | nt:<br>ange:<br>nod:<br>Ilgated fo<br>Digestior | pr<br>1      |
|  |   |  |   |              |
| C  | ontinue   |  |   |              |
| 4  | 0   |  |   |              |

#### SCAN-Anwendung

SCAN ist ein Multi-Wellenlängen-Scan, der die Absorption (ABS) oder die prozentuale Transmission (%T) über einen justierbaren Wellenlängenbereich und eine justierbare Geschwindigkeit und Intervallauflösung auswählt. Der Zweck der Funktion "SCAN" besteht darin, die Absorptions- oder Transmissionseigenschaften einer Probe oder einer mit Reagenz versetzten Probe zu bewerten. Dies ist nützlich, um die optimale Wellenlänge für die Einführung einer neuen Methode zu bestimmen.

Für die Probencharakterisierung können benutzerdefinierte Proben-Scanmethoden erstellt werden, wobei jedoch keine Konzentrationsmessungen vorgenommen werden.

|               |      |          | _         | Scanbereich:  |
|---------------|------|----------|-----------|---|
| Kethod Name   |      | New      | Accessory | Zwischen 190 nm und 1.100 nm für UV-VIS                 |
| GD1<br>Y-Axis | ABS  | %T       | SETUP     | Zwischen 325 nm und 1.100 nm für Vis-<br>Instrumente    |
| Range:        | 325  | - 1100   | nm        | Intervall:  |
| Interval      | 2 _  |          | nm        | Legt fest, wie häufig das Gerät eine Messung erfasst.   |
| Speed:        | Fast |          |           | In dieser Abbildung werden die Daten alle 2 nm erfasst. |
|               |      | Continue |           | Die Scan-Geschwindigkeiten "Fast" (Schnell) "Medium"    |

Die Scan-Geschwindigkeiten "Fast" (Schnell) "Medium" (Mittel) und "Slow" (Langsam) begrenzen die Anzahl der angebotenen Optionen für das Datenerfassungsintervall

| Speed  | Interval Options                  |
|--------|-----------------------------------|
| Fast   | 5 nm, 2 nm                        |
| Medium | 5 nm, 2 nm, 1 nm                  |
| Slow   | 5 nm, 2 nm, 1 nm, 0.5 nm, 0.2 nm, |

#### "Live-Display" (Live-Anzeige)

Im Live-Display-Modus führt das Gerät kontinuierliche Absorptions- (ABS) oder Transmissionsmessungen (%T) bei der ausgewählten Wellenlänge in Echtzeit durch.

Die Wellenlänge kann mit Hilfe der stufenlosen Skaleneinstellleiste oder durch Antippen der blauen Wellenlänge direkt bearbeitet werden. Es kann entweder "ABS" oder "%T" gewählt werden. Nach jeder Änderung sollte das System neu mit dem Blindwert angepasst werden, bevor die Live-Display-Messungen fortgesetzt werden.

Sobald das Gerät auf den Blindwert eingestellt ist, liefert das System automatisch Echtzeit- und Dauermessungen, bis das **X** in der oberen linken Ecke angetippt wird.





Für das Gerät muss zunächst der Blindwert bestimmt werden

#### Bildschirm 2 – Methodenentwicklung, Diagnostik und Daten

Wischen Sie auf Bildschirm 1 nach links, um Bildschirm 2 aufzurufen. Eine Anwendung wird durch Antippen des entsprechenden Symbols ausgewählt. Auf Bildschirm 2 befinden sich folgende Anwendungen:

- "C-Mode" (C-Modus) Wählen Sie eine Wellenlänge, geben Sie eine bekannte Standardkonzentration ein und wählen Sie die Maßeinheiten. Durch Blindwertanpassung und Messung des Standards wird ein Ein-Punkt-Kalibrierungsbericht von ABS erstellt und der "Cal Factor" (Kalibrierungsfaktor) berechnet und protokolliert.
- Quant ist eine Anwendung zur Erstellung von Kalibrierkurven. Der Benutzer kann das Verfallsdatum, die Wellenlänge, die Referenzwellenlänge, die Gleichung und die Konzentrationseinheiten auswählen und die bekannten Standards eingeben, die für die Erstellung der Kurve verwendet werden sollen. Die Methoden können unter einem Namen gespeichert und zur anschließenden Messung der Probenkonzentration verwendet werden.
- Die Funktion "Fixed" (Fest) ermöglicht die Eingabe einer Einzel- oder Multiwellenlängen-Methode auf der Grundlage von Daten, die durch Drittanbieter dokumentiert wurden. Dazu zählen "ABS" oder "%T", Wellenlänge\_1, Wellenlänge\_2, Maßeinheiten und die jeweiligen Wellenlängenfaktoren, sowie entweder eine direkte, additive, differentielle oder eine ratiometrische Gleichung. Die Methoden können unter einem Namen gespeichert und für die Messung der Konzentration nachfolgender Proben verwendet werden.
- Kinetics (Kinetik) ist ein aktiver Scan mit einer ausgewählten festen Wellenlänge und optionalen Referenzwellenlänge über einen festen Zeitraum, wobei die Daten in einem ausgewählten Intervall und Integrationszeitraum gestreamt werden. Wenn die eingestellte Versuchszeit erreicht ist, wird der Versuch beendet. Die Methoden können unter einem Namen gespeichert und zur Wiederholung eines Kinetik-Scans verwendet werden.
- Diagnostics (Diagnose) aktiviert eine Liste von Leistungstests, die am Gerät durchgeführt werden können. Ist ein geplantes Testintervall überfällig, wird ein rotes Uhrensymbol angezeigt. Diese Testintervalle sind vom Laborprotokoll abhängig. Durch Auswahl eines der verfügbaren Tests kann ein Leistungstest durchgeführt werden.
- Performance Verification Report (Leistungstestbericht) listet das Datum und die Anzahl der an diesem Tag durchgeführten Leistungstests auf. Durch Auswahl dieses Datums werden die Berichte ausgeklappt, und ein Bericht kann ausgewählt, angezeigt und gedruckt und/oder exportiert werden.
- Data Viewer (Datenanzeige) listet das Datum und die Anzahl der an diesem Tag durchgeführten Versuche auf. Durch Auswahl dieses Tages werden die Berichte ausgeklappt, und ein Bericht kann ausgewählt, angezeigt und gedruckt und/oder exportiert werden.

Die nachfolgende Abbildung "Screen 2" (Bildschirm 2) zeigt die oben beschriebenen Anwendungen. Am Ende der Seite befindet sich eine Kurzlegende.



#### C-Mode (C-Modus)

Im C-Modus kann der Benutzer eine Wellenlänge auswählen, eine bekannte Standardkonzentration eingeben und Maßeinheiten auswählen. Durch Ausblenden und Messen kann der Anwender eine Ein-Punkt-Kalibrierung entwickeln und protokollieren, wie hoch die Absorption ist und welcher Faktor erforderlich ist, um die Korrelation mit dem eingegebenen bekannten Konzentrationswert zu erreichen.

| ×           | C-M     | ē       |  |  |
|-------------|---------|---------|--|--|
| λ           |         |         |  |  |
| Std. Concen | tration | 5.000   |  |  |
| Units       |         | mg/mL   |  |  |
| Std. Absorb | ance    |         |  |  |
| Factor      |         |         |  |  |
| -           | m       | ig/mL   |  |  |
|             | Presen  | t blank |  |  |
| Blank       | Measure |         |  |  |


#### Quant

Die Anwendung Quant dient zur Entwicklung von Kalibrierkurven. Der Benutzer kann das Verfallsdatum, die Wellenlänge, die Referenzwellenlänge, die Gleichung und die Konzentrationseinheiten auswählen und die bekannten Standards eingeben, die für die Erstellung der Kurve verwendet werden sollen. Methoden können unter einem Namen gespeichert und später zur Messung einer Probenkonzentration verwendet werden



Der zur Anpassung der Standardmessdaten verwendete Gleichungstyp wird durch diese Gleichung festgelegt.

| Auswahl, ob 1 oder |  |
|--------------------|--|
| 2 Wiederholungen   |  |
| erforderlich sind  |  |

Auswahl der Maßeinheiten, in denen die Probenmessungen angezeigt und gedruckt werden



Sobald genügend eindeutige Standardkonzentrationswerte vorhanden sind, wird die Schaltfläche "Calibrate" (Kalibrieren) aktiviert.

Die Anzahl der eindeutigen Standardkonzentrationswerte wird anhand der Gleichung des Kurventyps bestimmt.







Standards hinzufügen und Kurventyp auswählen

#### **Optionen in Quant**

Wenn Sie auf die rechts neben der Wellenlänge aufgeführte Gleichung tippen, erscheint eine Auswahl von Gleichungen. Mit zunehmender Komplexität der Gleichung steigt auch die Anzahl benötigter Kurvenpunkte. Wählen Sie die für Sie am besten geeignete Gleichung. Nach Abschluss der Mehrpunktkalibrierung werden die resultierende Gleichung und die Korrelationsergebnisse (r2) angezeigt. Diese Ergebnisse beruhen auf der Qualität des Blindwertes und der Genauigkeit der erstellten und eingegebenen Standards.

Speichern Sie die Ergebnisse, indem Sie auf das blaue Disketten-Symbol oben rechts in der Ecke tippen.



#### Entwicklung voreingestellter Methoden ("Fixed")

Die Funktion "Fixed" (Fest) ermöglicht die Eingabe einer Einzel- oder Multiwellenlängen-Methode auf der Grundlage von Daten, die durch Drittanbieter dokumentiert wurden. Dazu zählen "ABS" oder "%T", Wellenlänge\_1, Wellenlänge\_2, Maßeinheiten und die jeweiligen Wellenlängenfaktoren, sowie entweder eine direkte, additive, differentielle oder eine ratiometrische Gleichung. Die Methoden können unter einem Namen gespeichert und für die Messung der Konzentration nachfolgender Proben verwendet werden



|   |   |  | 8 |
|---|---|--|---|
|   |   |  |   |
|   |   |  |   |
|   |   |  |   |
|   |   |  |   |
|   |   |  |   |
|   |   |  |   |
|   |   |  |   |
|   |   |  |   |
| Choose equatio                          |   |  |   |
| ABS(λ,)xF,                              |   |  |   |
| ABS(λ <sub>1</sub> )xF <sub>1</sub> +   | ABS(λ <sub>2</sub> )xF <sub>2</sub>           |  |   |
| ABS(λ <sub>1</sub> )xF <sub>1</sub> - A | $ABS(\lambda_1)xF_1 \cdot ABS(\lambda_2)xF_2$ |  |   |
| (ABS(λ,)xF,) /                          | / (ABS(λ <sub>2</sub> )xF <sub>2</sub> )      |  |   |

#### Messung mit der Methode "Kinetics" (Kinetik)

Die Methode "Kinetics" (Kinetik) ist ein aktiver Scan mit einer ausgewählten festen Wellenlänge und optionalen Referenzwellenlänge über einen festen Zeitraum, wobei die Daten in einem ausgewählten Intervall und Integrationszeitraum gestreamt werden. Wenn die eingestellte Versuchszeit erreicht ist, wird der Versuch beendet. Die Methoden können unter einem Namen gespeichert und zur Wiederholung eines Kinetik-Scans verwendet werden. Diese Anwendung dient dazu, die Reaktion oder den Zerfall einer Probe im Laufe der Zeit zu beobachten.



#### Diagnosemenü

Über das Menü "Diagnostics" (Diagnose) wird eine gespeicherte Liste von Leistungstest geöffnet, die mit dem Gerät durchgeführt werden können. Ist ein geplantes Testintervall überfällig, wird ein rotes Uhrensymbol angezeigt. Diese Testintervalle sind vom Laborprotokoll abhängig. Durch Auswahl eines der verfügbaren Tests kann ein Leistungstest durchgeführt werden. Tests, die kein zusätzliches Zubehör erfordern:

- Wavelength Accuracy (Wellenlängen-Genauigkeit)
- Drift at 500 nm (Drift bei 500 nm)
- Noise 0.0A at 500 nm (Rauschen 0,0 A bei 500 nm)
- Baseline Flatness (Basislinienabweichung)

|                | Performanc     | ce Verifica | tion Home    | •   |
|----------------|----------------|-------------|--------------|-----|
| Method name    |                |             |              |     |
| Wavelength A   | Accuracy Xer   | non         |              |     |
| Drift at 500nr | n              |             |              |     |
| Noise 0.0A at  | t 500nm        |             |              |     |
| Noise 1.0A at  | t 500nm        |             |              |     |
| Noise 2.0A at  | t 500nm        |             |              |     |
| Wavelengt      | h Accuracy     | / Xenon     |              |     |
| Wavelength     | ı Repeatabilit | ty test     |              | ••  |
| Wave           | elength (nm)   |             | Tolerance (r | ım) |
| 2              | 60.55          |             | +/-0.5       |     |
| 5              | 29.22          |             | +/-0.5       |     |
| 8              | 23.16          |             | +/-0.5       |     |
|                |                | Run         |              |     |
|                | Q              | 0           |              |     |

#### Leistungstestbericht

Wenn Sie "Performance Verification Reports" (Leistungstestbericht) auswählen, werden das Datum und die Anzahl der an diesem Tag durchgeführten Leistungstests aufgelistet. Wenn Sie das Datum und einen bestimmten Versuch auswählen, wird der zugehörige Bericht angezeigt. Diese Berichte werden aufgeklappt und ein Bericht kann ausgewählt, angezeigt und gedruckt und/oder exportiert werden

| Performance Verificatio                      | on Report 🔍 🔍       | ÷ | Wavele           | ength_Accuracy_X | enon_3_6_2019    | 🖶      |
|--|---------------------|---|------------------|------------------|------------------|--------|
| Time range<br>All                            |                     |   |                  | Wavelength Ac    | curacy Xenon     |        |
| <ul> <li>Wednesday, March 6, 2019</li> </ul> | 1 experiment        |   | λ (nm)           | Tolerance (nm)   | Found (nm)       | Result |
| Wavelength_Accuracy_Xenon_3                  | 8_6_2019_8:00:56_PM |   | 260.55<br>529.22 | ±0.5<br>±0.5     | 260.55<br>529.09 |        |
| <ul> <li>Sunday, March 3, 2019</li> </ul>    | 3 experiments       |   | 823.16           | ±0.5             | 822.96           | 1      |
| <ul> <li>Monday, February 5, 2018</li> </ul> | 1 experiment        |   |                  |                  |                  |        |
| Wednesday, January 24, 2018                  | 4 experiments       |   |                  |                  |                  |        |
| Tuesday, January 23, 2018                    | 1 experiment        |   |                  |                  |                  |        |
| Monday, January 22, 2018                     | 1 experiment        |   |                  |                  |                  |        |
| Friday, January 19, 2018                     | 10 experiments      |   |                  |                  |                  |        |
| Thursday, January 18, 2018                   | 3 experiments       |   |                  |                  |                  |        |
| •  |                     |   |                  |                  |                  |        |
| Generate Repo                                | rt                  |   |                  |                  |                  |        |
|  |                     |   |                  |                  |                  |        |

#### Datenanzeige

Nach der Auswahl von "Data Viewer" (Datenanzeige) werden das Datum und die Anzahl der an diesem Datum durchgeführten Versuche aufgelistet. Durch Auswahl dieses Datums werden die Berichte ausgeklappt, und ein Bericht kann ausgewählt, angezeigt und gedruckt und/oder exportiert werden. Der Versuch und alle grafischen Daten können so dargestellt werden, wie sie am Datum des Versuchs angezeigt wurden.





Nach Gleichung "adn r-Quadrat" berechneter Wert

## Bildschirm 3 – Mehrfachwellenlänge und OD600

Wischen Sie auf Bildschirm 2 nach links, um Bildschirm 3 aufzurufen. Eine Anwendung wird durch Antippen des entsprechenden Symbols ausgewählt. Auf Bildschirm 3 befinden sich folgende Anwendungen:

- Multi-wavelength (Mehrfachwellenlänge) Mit dieser Anwendung werden mehrere Messungen mit festen Wellenlängen durchgeführt. Sie ist eine schnelle Alternative zum Scannen (Anwendung "Scanning" [Scan]), wenn die zu untersuchenden Wellenlängen bekannt sind.
- OD600 reserviert

# Geräteeinstellungen

## Einstellungen

Die Geräteeinstellungen sind von jedem Hauptbildschirm aus durch Tippen auf das Zahnradsymbol in der oberen rechten Ecke zugänglich. Der Bildschirm "Settings" (Einstellungen) bietet folgende Optionen:

- Smart Start Bei Aktivierung werden nur "Smart Methods" (Intelligente Methoden) auf dem Startbildschirm angezeigt.
- Unter "Language" (Sprache) stehen die Sprachen Englisch, Deutsch, Italienisch, Spanisch, Französisch, Portugiesisch und mehrere asiatische Sprachen zur Verfügung
- Unter "Display and Sound" (Anzeige und Ton) werden die Regler in ihrer Intensität angepasst
- Unter "Network" (Netzwerk) lassen sich die Wi-Fi- und Ethernet-Einstellungen anpassen.
- Unter "Printer" (Drucker) können Netzwerk- oder optionale Thermodrucker-Einstellungen aktiviert werden
- Datum und Uhrzeit
- Festplattenspeicherplatz
- "Lamp Status" (Lampenstatus) wird im Kapitel "Wartung" behandelt



## **SmartStart**

SmartStart umfasst folgende Funktionen:

- Nur Anzeige von Methoden, die als "SmartStart" gekennzeichnet sind.
- Anzeigen oder Ausblenden des Data Viewer (Datenanzeige)
- Sperren des SmartStart-Modus mit optionalem Kennwortschutz, so dass der Anwender nur die verfügbaren Methoden nutzen kann
- Abgekürztes Einstellungsmenü, um die Einstellungs-Möglichkeiten zu begrenzen



SmartStart entsperren

Der Anwender muss über ein Passwort eines USB-Passwort-Rücksetzschlüssels verfügen, um die Geräte zu entsperren. Wenn Sie das Passwort nicht mehr wissen, wenden Sie sich an Ihren technischen Support und fordern Sie einen Schlüssel an.

## Sprache

Es stehen mehrere Sprachen zur Auswahl. Bitte wählen Sie die Sprache gemäß Ihren Anforderungen. Sollten Probleme mit der Sprachgenauigkeit auftreten, wenden Sie sich bitte an den technischen Support unter <u>wlp.techsupport@thermofisher.com.</u>

| ÷ | Language  |
|---|-----------|
| Ø | English   |
| 0 | Deutsch   |
| 0 | Italiano  |
| 0 | 日本語       |
| 0 | 한국어       |
| 0 | 中文        |
| 0 | Español   |
| 0 | Français  |
| 0 | русский   |
| 0 | ไทย       |
| 0 | Português |
|   |           |

## Anzeige, Ton und Netzwerk

Nachfolgend finden Sie die Bildschirmansichten für die Anzeige-, Ton- und Netzwerkeinstellungen.

Wenn kein Netzwerkpfad eingerichtet werden kann, wenden Sie sich bitte an Ihren Systemadministrator.



#### Druckereinstellungen

Unten sehen Sie das Fenster mit den Druckereinstellungen. Wenn Sie über einen Thermodrucker verfügen, lesen Sie bitte Abschnitt 2 dieses Handbuchs.

Für die Konfiguration kann entweder ein Thermodrucker oder ein USB-/Netzwerkdrucker ausgewählt werden. Der Berichtskopf kann so angepasst werden, dass er den Anwendungsfall des Spektrophotometers widerspiegelt.

| ÷                 | Printer                      |  |
|-------------------|------------------------------|--|
| Printer selectior | 1                            |  |
| S Thermal         | printer                      |  |
| O Network         | or USB printer               |  |
|                   |                              |  |
| Report header (Op | otional)                     |  |
| Header 1          | Thermo Fisher Scientific     |  |
| Header 2          | Water and Lab Products - WAI |  |
|                   |                              |  |
|                   |                              |  |
|                   |                              |  |
|                   | Toot                         |  |
|                   |                              |  |
|                   |                              |  |

## Datum und Uhrzeit

In diesem Einstellungsmenü können Sie das Datum, die Uhrzeit und das Zeitformat einstellen.





#### **Speicherplatz und Lampenstatus**

Auf den folgenden Bildschirmen werden zur Ansicht sowohl der verfügbare Speicher als auch die Lampenbetriebsdauer angezeigt. Nach dem Austausch der Lampe sollte die Lampenbetriebsdauer zurückgesetzt werden.

HINWEIS: Beim 7100 AquaMate sollte die Wolframlampe bei Nichtgebrauch in einen Lampenschonmodus versetzt werden. Sie schaltet sich nach 15 Minuten Nichtbenutzung automatisch ab. Bitte denken Sie daran, dass die Wolframlampe aufgewärmt werden muss, um exakte Ergebnisse zu erzielen.



#### Software-Update

Mithilfe der Funktion "Software Updates" (Software-Aktualisierungen) können zwei Aktualisierungen vorgenommen werden: Die Aktualisierung der Firmware und die Aktualisierung der Wasseranalysemethoden-Bibliothek. Wasseranalysemethoden sind nur spezifisch für AquaMate-Spektrophotometer. Es sind typische Analysemethoden für wasserhaltige Proben, in der chemische Reagenzien zur Erlangung von Analyseergebnissen eingesetzt werden. Wenn Sie aus irgendeinem Grund nicht in der Lage sind, Ihre Methodenbibliothek zu aktualisieren, wenden Sie sich bitte an den technischen Support, und geben Sie Ihr spezifisches Modell und die Seriennummer an.

#### Software-Update

Für eine Software-Aktualisierung wenden Sie sich bitte an <u>wlp.techsupport@thermofisher.com</u>, um die neueste Firmware und die neueste Methodenbibliothek zu erhalten, und speichern Sie diese auf Ihrem USB-Stick. Schließen Sie den USB-Stick an den USB-Anschluss an. Tippen Sie auf "Update Software" (Software aktualisieren). Tippen Sie auf die Version mit dem neuesten Veröffentlichungsdatum und dann auf "Update" (Aktualisieren). Das Gerät führt Sie selbständig durch den Vorgang und startet sich automatisch neu.

| ÷                | Software Updat                           | te                      | Software Updates          |                 | Current Version:          | 2.0      |
|------------------|--|-------------------------|---------------------------|-----------------|---------------------------|----------|
| Software Package | Version                                  | 2.0                     | Version<br>AquamateUpdate | R<br>eWithEng ( | elease Date<br>02-20-2019 | Eligible |
|                  |  |                         | -                         |                 |                           |          |
|                  |  |                         |                           |                 |                           |          |
| https://www      | For software updates thermofisher.com/Go | visit<br>enesysSoftware |                           |                 |                           |          |
| Upd              | ate Analyzer Me                          | ethods                  |                           |                 |                           |          |
|                  | Update Softwa                            | re                      | Cancel                    |                 | Update                    |          |
| <                | 1 0                                      |                         |                           | ⊲ 0             |                           |          |

#### Update der Wasseranalysemethoden-Bibliothek

Um die Bibliothek zu aktualisieren, wenden Sie sich an <u>wlp.techsupport@thermofisher.com</u>, um das neueste Paket mit Wasseranalysemethoden zu erhalten, und speichern Sie dieses auf Ihrem USB-Stick. Schließen Sie den USB-Stick an den USB-Anschluss an. Tippen Sie auf "Update" (Aktualisieren), um die Analyzer-Methoden zu aktualisieren. Tippen Sie auf das neueste AquaMate-Paket (z. B. aquamate 1.X) und tippen Sie auf "Activate" (Aktivieren). Das Gerät aktiviert und ersetzt die Methodenbibliotheken automatisch. Es ist kein Neustart erforderlich. Das rote X wird zu einem grünen Häkchen.



## Importieren und Exportieren von Versuchen

Wenn Ihre Messungen abgeschlossen sind, tippen Sie auf "End Experiment" (Versuch beenden). Der Versuch wird unter dem blau angezeigten Namen gespeichert.

In den unten aufgeführten Feldern können Sie folgende Aufgaben ausführen:

- Zurückgehen
- Name des Versuchs vor dem Ende bearbeiten
- Ergebnisse als Bericht an einen Drucker senden
- Ergebnisse in eine Datei exportieren (USB oder Netzwerkverbindung)



## **Daten exportieren**

Um die Daten exportieren zu können, müssen Sie einen Netzwerkpfad über Ethernet oder über WiFi eingerichtet haben. Eine typische Methode ist das direkte Speichern der Daten auf einem USB-Laufwerk. Der Bericht kann sowohl als CSV- als auch als JPEG-Datei gespeichert werden.



# Nachfolgend finden Sie ein Beispiel für das USB-Dateiverzeichnis und ein JPEG-Bild des Versuchs.

|        | Name                                | Date modified     |                    |
|--------|-------------------------------------|-------------------|--------------------|
|        | 🖀 2.1.zip                           | 4/2/2019 12:26 PM |                    |
|        | aquamatev1.5.pkg                    | 3/27/2019 1:46 PM |                    |
|        | — 👢 Aquamate Experiment Data        | 4/11/2019 1:48 PM |                    |
|        | 👢 Aquamate Log                      | 3/28/2019 6:16 PM |                    |
|        |                                     |                   |                    |
|        |                                     |                   |                    |
|        | Scan 3 6 2019 7 41 02 PM csv        | 3/28/2019 6·17 PM | Microsoft Excel Co |
|        | Aquamate_4_9_2019_4_35_06_PM.csv    | 4/11/2019 1:48 PM | Microsoft Excel Co |
|        | Scan_3_6_2019_7_41_02_PM.jpeg       | 3/28/2019 6:17 PM | JPEG image         |
| $\neg$ | Nquamate_4_9_2019_9_23_32_AM-2.jpeg | 4/11/2019 1:48 PM | JPEG image         |
|        | Nquamate_4_9_2019_9_23_32_AM-1.jpeg | 4/11/2019 1:48 PM | JPEG image         |
|        | Nquamate_4_9_2019_4_35_06_PM.jpeg   | 4/11/2019 1:48 PM | JPEG image         |

| Scan_3_6_2019_7:41:02_PM             |                                 |            | 1/1 |
|--------------------------------------|---------------------------------|------------|-----|
| 28-Mar-2019 D6:16 PM                 | instrument Serial #: 9A3V205001 |            |     |
| Method name: Quick scan              | Instrument model: AQ8100        |            |     |
| Method created: 25-Mar-2018 08:14 AM | Software Package Version: 2.0   | Signature: |     |
| Method updated: 26-Mar-2018 08:15 AM |                                 |            |     |

#### Scan: method parameters



| Sample    | ABS(550) | ABS(650) |
|-----------|----------|----------|
| Mt. Dew 1 | 0.094    | 0.065    |
| Mt. Dew 2 | 0.094    | 0.054    |
| Mt. Dew 3 | 0.094    | 0.065    |

# 

# **KAPITEL 4** Testmenü für Wasseranalysen

## Vorprogrammierte Methoden

Die Spektrophotometer Orion AquaMate 8100 UV-Vis und AquaMate 7100 Vis von Thermo Scientific enthalten über 260 vorprogrammierte Methoden zur Verwendung mit Orion<sup>™</sup> AQUAfast<sup>™</sup>, Merck und CHEMetrics-Reagenzien von Thermo Scientific<sup>™</sup>. Vorprogrammierte Methoden enthalten Werte für die Testparameter, die für die Durchführung bestimmter Reagenzien-Analysen auf dem Gerät erforderlich sind, einschließlich Wellenlänge, Probegefäß-Pfadlänge, Konzentrationsfaktoren/Kurven und Maßeinheiten. Alle behördlich anerkannten Methoden für Abwasser- oder Trinkwasseranalysen sind zwecks einfacher Zugänglichkeit in den jeweiligen Droplet-Ordnern dupliziert.

Alle vorprogrammierten Methoden und die Dokumentation sind als Methodenbibliothek auf dem Gerät gespeichert und auch auf einem USB-Speicherstick verfügbar. Die vorprogrammierten Methoden sind nur für AquaMate-Spektrophotometer geeignet. Vorprogrammierte Methoden können vom Anwender geändert oder eigene benutzerdefinierte Methoden erstellt werden, so dass jederzeit zusätzliche Parameter und Prüfmethoden hinzugefügt werden können.

AquaMate-Geräte ermöglichen eine Ein-Punkt-Justierung bei jeder vorprogrammierten Methode unter Verwendung eines bekannten Standards, um chargenpezifische Schwankungen in der chemischen Zusammensetzung der Reagenzien zu korrigieren.

Die folgenden Anweisungen gelten für die Verwendung von Orion AQUAfast-Reagenz, Merckoder CHEMetrics-Chemikalien mit dem AquaMate-Spektrophotometer. Vorprogrammierte Methoden verwenden eine bestimmte Probengefäßgröße (Pfadlänge) in der Formel. Um eine präzise Analyse zu erhalten, muss die in dieser Anleitung angegebene Probengefäßgröße verwendet werden. Die meisten AQUAfast-Reagenzienmethoden verwenden eine 24-mm-Rundküvette, Best.-Nr. AC2V24, oder eine 16-mm-Rundküvette, Best.-Nr. AC2V16. Andere Probengefäßgrößen sind in den Anleitungen zu den einzelnen Reagenzien angegeben.

# Methodenauswahl und Versuch

## Droplet-Ordner-Methoden

Wählen Sie einen beliebigen Droplet-Symbolordner, um die Methoden in diesem Ordner aufzulisten. Um die bekannten behördlichen Analysemethoden für Abwasser- und Trinkwasserproben aufzurufen, wählen Sie den entsprechenden Ordner. Innerhalb des Ordners können die Methoden nach Name, Erstellungsdatum, Datum der letzten Aktualisierung oder Datum der letzten Verwendung sortiert werden. Sie können auch innerhalb des jeweiligen Ordners nach einer Methode suchen, indem Sie entweder die Methodennummer oder den Methodenparameter eingeben.

Die Methodenbeschreibungen enthalten Angaben zu Methodennummer, Parameter, Wellenlänge, Probengefäßtyp und -größe sowie detaillierte Informationen zur Methode selbst. Befolgen Sie die Methodenanweisungen nach den Reagenzienanweisungen in Kapitel 5 oder den Herstelleranweisungen.

HINWEIS: Wenn Sie auf das Ellipsen-Symbol einer ausgewählten Methodenbeschreibung tippen, wird eine SmartStart-Option angezeigt, wenn Sie diese Methode für SmartStart aktivieren möchten.



**Hinweis:** Die Messmöglichkeiten für jede Methode entnehmen Sie bitte der jeweiligen Methodenbeschreibung. Das Gerät meldet Werte, die außerhalb des angegebenen Messbereichs liegen und für den speziellen Zweck des Anwenders oder behördliche Berichtsanforderungen möglicherweise nicht akzeptabel sind.

## **Methoden-Optionen**

Bei der ausgewählten Methode sind die blau markierten Felder normalerweise editierbar. Im folgenden Beispiel kann der "Sample Base Name (Sample)" (Name der Probenbasis [Probe]) mithilfe des beim Antippen angezeigten alphanumerischen Touchscreen-Feldes angepasst werden. Wenn das zu bearbeitende Feld vollständig ausgefüllt ist, drücken Sie die Taste "Done" (Fertig) auf der Touchscreen-Tastatur.

- Bereiten Sie die Blindlösung, den Einsatz und den Blindwert vor.
- Wenn Sie eine neue Charge oder ein neues Reagenz verwenden, können Sie einen bekannten und rückverfolgbaren Standard vorbereiten, den Standardkonzentrationswert bearbeiten und die Option "Re-Calibrate" (Neu kalibrieren) auswählen. Dadurch wird das Feld "Calibration Factor" (Kalibrierungsfaktor) aktualisiert. Andernfalls fahren Sie mit der Messung fort.
- Legen Sie die vorbereitete Probe ein und tippen Sie auf "Measure" (Messen), um die Ergebnisse zu erhalten.



#### **Probennummerierung einer Methode**

Wenn Sie die Konzentration einer Probe mit einer beliebigen Methode oder Anwendung messen, wird die Probennummer unter "Sample Base Name" (Name der Probenbasis) jedes Mal um 1 erhöht, wenn Sie die Taste "Measure" (Messen) drücken. Siehe hierzu die rot eingekreisten Bereiche.



## Laden von Testmethoden aus dem AquaMate-Gerät

- 1. Wählen Sie einen Droplet-Ordner
- 2. Suchen Sie entweder nach der Methodennummer oder nach dem Parameter
- 3. Wählen Sie die Methode aus
- 4. Überprüfen Sie die Beschreibung und stellen Sie sicher, dass Sie die angegebene Probengefäßgröße verwenden.



## Laufende Testmethoden für die Wasseranalyse

- 1. Wenn eine vorprogrammierte Testmethode geladen wurde, wird ein Versuchsfenster angezeigt.
- 2. Tippen Sie auf den Namen der Probenbasis (z. B. COD Site A), um diese mithilfe der Bildschirmtastatur zu bearbeiten.
- 3. Öffnen Sie den Probenkammerdeckel.
- 4. Fläschchen mit der Blind- oder Null-Lösung in den Probenhalter einsetzen.

**Hinweis:** Idealerweise sollte dasselbe Fläschchen oder ein auf das Fläschchen abgestimmtes Gefäß verwendet werden.

- 5. Schließen Sie den Deckel und tippen Sie auf die Funktionstaste Blank (Blind).
- 6. Öffnen Sie den Deckel und nehmen Sie das Fläschchen mit der Blind- oder Null-Lösung heraus.
- 7. Wenn eine Ein-Punkt-Kalibrierung erforderlich ist, gehen Sie wie im folgenden Abschnitt beschrieben vor.
- 8. Stellen Sie das Fläschchen mit der Probe in den Probenhalter, und schließen Sie den Deckel.
- 9. Tippen Sie auf die Funktionstaste **Measure (Messen)**. Die Ergebnisse werden angezeigt. Es können in der Regel fortlaufende Messungen für mehrere Proben durchgeführt werden.
- 10. Zum Speichern tippen Sie auf das X in der oberen linken Ecke. Damit beenden und speichern Sie den Versuch.
- 11. Die Daten werden automatisch mit Namen, Datum und Zeitstempel unter dem gewählten Namen gespeichert (z. B. "Scan\_3\_6\_2019\_7\_41\_02\_PM")

**Hinweis:** Die Blindmessung wird gespeichert und bei der Arbeit in der Probentestmethode verwendet. Die Blindmessung wird automatisch gelöscht, wenn die Einstellungen der Testmethode geändert werden, die Testmethode gespeichert oder eine neue Testmethode geladen wird.

**Hinweis:** Bei der Durchführung einer Farbumkehr-Testmethode ist nach dem Standardblindwert eine Reagenzblindmessung erforderlich. Setzen Sie das Fläschchen mit der Reagenzblindprobe ein und drücken Sie die Funktionstaste Blank (Blind). Öffnen Sie den Deckel und entnehmen Sie das Blindreagenz. Ausführliche Anweisungen finden Sie im Abschnitt <u>Verwendung der Farbumkehrfunktion</u>.

**Hinweis:** Wenn die Option "Statistics" (Statistik) auf "Off" (Aus) gesetzt ist, werden keine Statistiken angezeigt.

Nachstehend wird ein Beispiel-Bildschirm für eine COD-Methode mit typischen blau hervorgehobenen Feldern gezeigt, die möglicherweise zu beachten sind. Außerdem wird der letzte Bildschirm gezeigt, der Optionen für die Beendigung und Benennung des Versuchs sowie das Drucken und Exportieren der Ergebnisse enthält.



## Ein-Punkt-Anpassung der Methode

Jede Methode kann durch eine Ein-Punkt-Kalibrierung angepasst werden.

- Bearbeiten Sie zunächst den Standardkonzentrationswert von 1,000. Hierzu tippen Sie auf den Wert und geben den Wert der für diesen Zweck vorbereiteten Standardkonzentration ein.
- Nach der Blindwertmessung des Systems wird der vorbereitete Standard in den Probenhalter eingesetzt.
- Tippen Sie auf die Schaltfläche "Re-Calibrate" (Neu kalibrieren)
- Normalerweise ist ein Kalibrierungsfaktor von 0,7 bis 1,3 (innerhalb von ± 30 %) akzeptabel.
- Das Fläschchen kann nun in den Probenhalter gestellt und die Messung durchgeführt werden.

| ×                            | CODL00, COD(le    | ow range)             |
|------------------------------|-------------------|-----------------------|
| Sample base name<br>Test One | 8/15              | Re-calibrate          |
| Standard Con                 | centration Value  | 1.000                 |
| Calibration Fa               | ctor              | 1.000                 |
| Y = -318.719X                | 0.000<=Y<=150.000 | 0.2 0.4               |
|                              | Absort            | bance                 |
| Sample                       | ABS               | Concentration(mg/I 02 |
|                              |                   |                       |
|                              |                   |                       |
| Blank                        | Single F          | Point Calibration     |

## Verwenden der Farbumkehrfunkion

Bei der Farbumkehrmethode wird ein Reagenz verwendet, dessen Farbintensität mit zunehmender Konzentration der zu messenden Spezies in den Proben abnimmt. Die Farbumkehrmethode erfordert die Verwendung eines Reagenzblindwertes. Der Reagenzblindwert ist ein Gemisch aus dem Ausgangsreagenz und der Probe (z. B. Zink nach der Zincon-Methode) und liefert den Nullkonzentrationspunkt mit dem dunkelsten Farbwert (höchste Absorption). Die Farbe der mit dem Reagenz versetzten Proben nimmt mit steigender Konzentration ab. Die folgende Abbildung zeigt die Ergebnisse einer typischen Farbumkehrmethode. Anschließend folgt ein Überblick über die Durchführung der Farbumkehrmethode.



- Laden Sie die Testmethode im Testmenü "Water Analysis" (Wasseranalyse). Die Option "Reverse Color" (Farbumkehr) (Negative ABS) sollte für die Methode auf <u>ON (EIN)</u> gesetzt werden.
- 2. Bereiten Sie das Probenreagenz in dem für die Methode vorgesehenen Fläschchen vor.
- 3. Stellen Sie das Fläschchen in die Halterung in der Probenkammer und schließen Sie die Probenkammertür.
- 4. Drücken Sie die Funktionstaste Blank (Blind), um den Reagenzblindwert zu messen.
- 5. Öffnen Sie die Probenkammertür, und nehmen Sie das Fläschchen aus der Halterung.
- 6. Befolgen Sie die Methodenanweisung, und bereiten Sie die zu messende Probe vor.
- 7. Stellen Sie das Fläschchen in die Halterung in der Probenkammer und schließen Sie die Probenkammertür.
- 8. Drücken Sie die Funktionstaste Measure (Messen), um die Ergebnisse anzuzeigen.
- 9. Messen Sie je nach Bedarf weitere Proben.
- 10. Wenn Sie fertig sind, beenden Sie den Versuch, und exportieren oder drucken Sie die Daten.



# **KAPITEL 5** Orion AQUAfast Reagenzien – Anleitung für Orion AquaMate

# Orion AQUAfast Kolorimetrische Reagenzien Kompatibel mit Orion AquaMate-Geräten

Verwenden Sie die Angaben in der folgenden Tabelle, um den Dateinamen der Orion AQUAfast-Reagenzienmethode auf dem AQ7100 oder AQ8100 und die mit jeder Methode verbundenen Testparameter zu ermitteln. Diese Angaben finden Sie auch auf der CD mit der Benutzerdokumentation des Orion AquaMate oder auf unserer Website unter www.thermoscientific.com/water.

| Parameter   | Teile-Nr. | Methode  | Beschreibung   |
|-------------|-----------|----------|--|
| Alkalinität | AC2002    | AC2002   | Alkalinität-M Tablettenreagenz                           |
| Alkalinität | AC3002P   | AC3002P  | Alkalinität-P Tablettenreagenz                           |
| Aluminium   | AC2027    | AC2027   | Aluminium-Tablettenreagenz                               |
| Aluminium   | AC4P27    | AC4P27   | Aluminium-Pulverpäckchen und -Flüssigreagenz             |
| Ammoniak    | AC2012    | AC2012   | Ammoniak-Tablettenreagenz                                |
| Ammoniak    | AC4P12    | AC4P12   | Ammoniak-Pulverpäckchen-Reagenz                          |
| Ammoniak    | ACR012    | ACR012   | Ammoniak-Reaktionsröhrchen-Reagenz für niedrigen Bereich |
| Ammoniak    | ACR011    | ACR011   | Ammoniak-Reaktionsröhrchen-Reagenz für hohen Bereich     |
| Brom        | AC2035    | AC203524 | Brom-Tablettenreagenz                                    |
| Chlorid     | AC2017    | AC2017   | Chlorid-Tablettenreagenz                                 |
| Chlor       | AC2070    | AC207024 | Chlor-Tablettenreagenz (frei und gesamt)                 |
| Chlor       | AC2071    | AC207124 | Chlor-Tablettenreagenz (frei)                            |
| Chlor       | AC2072    | AC207224 | Chlor-Tablettenreagenz (gesamt)                          |
| Chlor       | AC4P71    | AC4P71   | Chlor-Pulverpäckchen-Reagenz (frei)                      |
| Chlor       | AC4P72    | AC4P72   | Chlor-Pulverpäckchen-Reagenz (gesamt)                    |

| Parameter             | Teile-Nr. | Methode  | Beschreibung  |
|-----------------------|-----------|----------|---|
| Chlor                 | AC3072    | AC3072   | Chlor-Tablettenreagenz (gesamt) für hohen Bereich                       |
| Chlordioxid           | AC2099    | AC209924 | Chlordioxid-Tablettenreagenz  |
| COD                   | CODL00    | CODL00   | COD-Reaktionsröhrchen-Aufschlussreagenz für niedrigen Bereich           |
| COD                   | CODH00    | CODH00   | CSB-Reaktionsröhrchen-Aufschlussreagenz für mittleren Bereich           |
| COD                   | CODHP0    | CODHP0   | COD-Reaktionsröhrchen-Aufschlussreagenz für hohen Bereich               |
| Kupfer                | AC2029    | AC202924 | Kupfer-Tablettenreagenz (frei und gesamt)                               |
| Kupfer                | AC4P29    | AC4P29   | Kupfer-Pulverpäckchen-Reagenz (frei)                                    |
| Cyanursäure           | AC2098    | AC2098   | Cyanursäure-Tablettenreagenz  |
| Fluorid               | AC2009    | AC2009   | Fluorid-SPADNS-Flüssigreagenz   |
| Härte                 | AC3032T   | AC3032TL | Wasserhärte-Tablettenreagenz (gesamt) für niedrigen Bereich             |
| Härte                 | AC3032T   | AC3032TH | Wasserhärte-Tablettenreagenz (gesamt) für hohen Bereich                 |
| Hydrazin              | AC2030    | AC2030   | Hydrazin-Pulverreagenz  |
| Eisen                 | AC2078    | AC207824 | Eisen-Tablettenreagenz (II und III)                                     |
| Eisen                 | AC4P78    | AC4P78   | Eisen-Pulverpäckchen-Reagenz (Ferro)                                    |
| Eisen                 | AC4P79    | AC4P79   | Eisen-Pulverpäckchen-Reagenz (gesamt)                                   |
| Mangan                | AC2055    | AC2055   | Mangan-Tablettenreagenz   |
| Mangan                | AC4P54    | AC4P54   | Mangan-Pulverpäckchen- und Flüssigreagenz für niedrigen<br>Bereich      |
| Mangan                | AC4P55    | AC4P55   | Mangan-Pulverpäckchen-Reagenz für hohen Bereich                         |
| Molybdat              | AC4P42    | AC4P42   | Molybdat/Molybdän-Pulverpäckchen-Reagenz                                |
| Nitrat                | ACR007    | ACR007   | Nitrat-Reaktionsröhrchen-Reagenz  |
| Nitrit                | AC2046    | AC2046   | Nitrit-Tablettenreagenz   |
| Nitrit                | AC4P46    | AC4P46   | Nitrit-Pulverpäckchen-Reagenz   |
| Stickstoff,<br>gesamt | ACD004    | ACD004   | Stickstoff-Aufschlussröhrchen-Reagenz (gesamt) für niedrigen<br>Bereich |
| Stickstoff,<br>gesamt | ACD007    | ACD007   | Stickstoff-Aufschlussröhrchen-Reagenz (gesamt) für hohen<br>Bereich     |
| Ozon                  | AC3048    | AC3048   | Ozon-Tablettenreagenz   |
| pН                    | AC2001    | AC2001   | pH-Tablettenreagenz   |
| pН                    | AC3001    | AC3001   | pH-Flüssigreagenz   |
| Phosphat              | AC2095-WA | AC2095   | Phosphat-Tablettenreagenz (Ortho) für niedrigen Bereich                 |
| Phosphat              | AC2096    | AC2096   | Phosphat-Tablettenreagenz (Ortho) für hohen Bereich                     |
| Phosphat              | AC4P95    | AC4P95   | Phosphat Pulverpäckchen-Reagenz (Ortho)                                 |
| Phosphat              | ACR095    | ACR095   | Phosphat Reaktionsröhrchen-Reagenz (Ortho)                              |
| Phosphat              | ACD095    | ACD095   | Phosphat Aufschlussröhrchen-Reagenz (gesamt)                            |
| Phosphat              | ACD095AH  | ACD095AH | Phosphat (säurehydrolysierbar) Aufschlussröhrchen-Reagenz               |
| Siliziumdioxid        | AC2060    | AC2060   | Siliziumdioxid-Tablettenreagenz   |
| Siliziumdioxid        | AC2061    | AC2061   | Siliziumdioxid mit Tablettenreagenz zur Phosphatentfernung              |
| Siliziumdioxid        | AC4P60    | AC4P60   | Siliziumdioxid-Pulverpäckchen-Reagenz                                   |
| Sulfat                | AC4P82    | AC4P82   | Sulfat-Pulverpäckchen-Reagenz   |
| Sulfid                | AC2016    | AC2016   | Sulfid-Tablettenreagenz   |
| Zink                  | AC2065    | AC2065   | Zink-Tablettenreagenz   |

# Anweisungen zum Orion AQUAfast-Reagenz

Die in den folgenden Testverfahren angegebenen Messbereiche basieren auf Standardlösungen, die unter idealen Bedingungen gemessen wurden. Diese Bereiche können je nach Art der zu messenden Probe variieren, da verschiedene Störungen einen großen Einfluss auf die Genauigkeit der Methode haben können. Da jede Probe anders ausfällt, ist die einzige Möglichkeit zur Überprüfung der Toleranz (Präzision) die Standardadditionsmethode. Bei dieser Methode wird zunächst die Originalprobe untersucht. Dann werden weitere Proben (2 bis 4) entnommen und kleine Mengen einer Standardlösung hinzugefügt, um weitere Ergebnisse zu erhalten. Die zugesetzten Mengen reichen von etwa der Hälfte bis zum Doppelten der in der Probe selbst vorhandenen Menge. Diese ergänzenden Ergebnisse ermöglichen es, die tatsächliche Konzentration der ursprünglichen Probe durch Vergleich zu schätzen.

Die Testmethoden und -bereiche können ohne vorherige Ankündigung geändert werden. Eine Liste der aktuellsten Testmethoden finden Sie unter <u>www.thermoscientific.com/water.</u>

## Empfehlungen zur Vermeidung von Messfehlern

- Reinigen Sie Fläschchen, Kappen und Rührstäbchen nach jeder Analyse gründlich, um Fehler durch Verschleppung zu vermeiden. Selbst kleinste Reagenzrückstände führen zu Fehlmessungen.
- Stellen Sie sicher, dass die Außenwände der Fläschchen trocken und sauber sind, bevor Sie die Analyse durchführen. Fingerabdrücke oder Wassertropfen auf den Lichteintrittsflächen der Fläschchen führen zu Fehlmessungen.
- Blank (Blind)- und Messverfahren sollten nach Möglichkeit mit demselben Fläschchen durchgeführt werden, da verschiedene Fläschchen leicht unterschiedliche Toleranzen aufweisen können.
- Führen Sie alle Messungen mit verschlossenen Fläschchen durch.
- Blasen an den Innenwänden des Fläschchens können zu Fehlmessungen führen. Um dies zu vermeiden, verschließen Sie das Fläschchen und entfernen Sie die Luftbläschen, indem Sie das Fläschchen vor der Durchführung des Tests schwenken.
- Geben Sie das Reagenz immer direkt aus dem Folienbeutel in die Probe. Das Reagenz darf nicht mit den Fingern oder Händen in Berührung kommen.
- Größere Temperaturunterschiede zwischen Gerät und Umgebung können zu Fehlmessungen führen – z. B. durch Kondensatbildung im Bereich der Linse oder auf der Küvette. Angegebene Toleranzen bei T = 20 °C.
- Die besten Ergebnisse erzielen Sie, wenn Sie die Proben mit einer Pipette abmessen und in die Fläschchen oder Bechergläser geben.

## AC2002 Alkalinität-M (Alkalinität bis pH 4,3) Tablettentest

#### Säure/Indikator-Methode

- 5 200 mg/l CaCO<sub>3</sub>
- 1. Laden Sie die Methode AC2002, und führen Sie sie aus.
- Befüllen Sie eine saubere AQUAfast 24-mm-Rundküvette, Bestell-Nr. AC2V24, mit 10 ml der Probe. Verschließen Sie das Fläschchen fest mit dem Deckel. Wischen Sie die Außenseite des Fläschchens ab.
- 3. Stellen Sie das Fläschchen in die Halterung in der Probenkammer und schließen Sie die Probenkammertür.
- 4. Drücken Sie die Funktionstaste Blank (Blind), um den Blindwert zu messen.
- 5. Öffnen Sie die Probenkammertür, und nehmen Sie das Fläschchen aus der Halterung.
- Geben Sie eine <u>Alka-M Tablet</u> direkt aus dem Folienbeutel in das Fläschchen. Zerdrücken Sie die Tablette mit einem sauberen Rührstab.
- Verschließen Sie das Fläschchen fest mit dem Deckel, und mischen Sie den Inhalt durch mehrmaliges Schwenken oder Umdrehen, bis sich die Tablette aufgelöst hat. Wischen Sie die Außenseite des Fläschchens ab.
- 8. Stellen Sie das Fläschchen in die Halterung in der Probenkammer und schließen Sie die Probenkammertür.
- 9. Drücken Sie die Funktionstaste Sample (Probe), um das Ergebnis in mg/l Gesamtalkalität zu erfassen.

#### Hinweise:

- Die Begriffe Gesamtalkalität, Alkalität-m, m-Wert und Alkalität bis pH 4,3 sind identisch.
- Um genaue Ergebnisse zu erzielen, müssen genau 10 ml Wasserprobe für den Test entnommen werden.

## AC3002P Alkalinität-P (Alkalinität bis pH 8,2) Tablettentest

#### Säure/Indikator-Methode

- 5 300 mg/l CaCO<sub>3</sub>
- 1. Laden Sie die Methode AC3002P, und führen Sie sie aus.
- Befüllen Sie eine saubere AQUAfast 24-mm-Rundküvette, Bestell-Nr. AC2V24, mit 10 ml der Probe. Verschließen Sie das Fläschchen fest mit dem Deckel. Wischen Sie die Außenseite des Fläschchens ab.
- 3. Stellen Sie das Fläschchen in die Halterung in der Probenkammer und schließen Sie die Probenkammertür.
- 4. Drücken Sie die Funktionstaste Blank (Blind), um den Blindwert zu messen.
- 5. Öffnen Sie die Probenkammertür, und nehmen Sie das Fläschchen aus der Halterung.
- 6. Geben Sie eine <u>Alka-P Tablet</u> direkt aus dem Folienbeutel in das Fläschchen. Zerdrücken Sie die Tablette mit einem sauberen Rührstab.
- Verschließen Sie das Fläschchen fest mit dem Deckel, und mischen Sie den Inhalt durch mehrmaliges Schwenken oder Umdrehen, bis sich die Tablette aufgelöst hat. Wischen Sie die Außenseite des Fläschchens ab.
- 8. Stellen Sie das Fläschchen in die Halterung in der Probenkammer und schließen Sie die Probenkammertür.
- 9. Drücken Sie die Funktionstaste Sample (Probe), um das Ergebnis in mg/l Gesamtalkalität zu erfassen.

#### Hinweise

- Die Begriffe Alkalinität-p, p-Wert und Alkalinität bis pH 8,2 sind identisch.
- Um genaue Testergebnisse zu erzielen, müssen genau 10 ml Wasserprobe für den Test entnommen werden.
- Diese Methode wurde aus einem volumetrischen Verfahren zur Bestimmung der Alkalinitätp entwickelt. Aufgrund nicht definierter Bedingungen können die Abweichungen von der standardisierten Methode größer sein.

## AC2027 Aluminium-Tabletten-Test

#### **Eriochrom Cyanin-R-Methode**

0,01 – 0,3 mg/l Al

- 1. Laden Sie die Methode AC2027, und führen Sie sie aus.
- Befüllen Sie eine saubere AQUAfast 24-mm-Rundküvette, Bestell-Nr. AC2V24, mit 10 ml der Probe. Verschließen Sie das Fläschchen fest mit dem Deckel. Wischen Sie die Außenseite des Fläschchens ab.
- 3. Stellen Sie das Fläschchen in die Halterung in der Probenkammer und schließen Sie die Probenkammertür.
- 4. Drücken Sie die Funktionstaste Blank (Blind), um den Blindwert zu messen.
- 5. Öffnen Sie die Probenkammertür, und nehmen Sie das Fläschchen aus der Halterung.
- 6. Geben Sie eine <u>Aluminum No. 1 Tablet</u> direkt aus dem Folienbeutel in das Fläschchen. Zerdrücken Sie die Tablette mit einem sauberen Rührstab, und mischen Sie gründlich, um die Tablette vollständig aufzulösen.
- Geben Sie eine <u>Aluminum No. 2 Tablet</u> direkt aus dem Folienbeutel in dasselbe Fläschchen. Zerdrücken Sie die Tablette mit einem sauberen Rührstab, und mischen Sie gründlich, um die Tablette vollständig aufzulösen.
- 8. Verschließen Sie das Fläschchen fest mit dem Deckel, und mischen Sie den Inhalt durch mehrmaliges Schwenken oder Umdrehen. Wischen Sie die Außenseite des Fläschchens ab.
- 9. Warten Sie, bis eine Reaktionszeit von 5 Minuten abgelaufen ist.
- 10. Stellen Sie das Fläschchen in die Halterung in der Probenkammer und schließen Sie die Probenkammertür.
- 11. Tippen Sie auf die Funktionstaste Sample (Probe), um das Ergebnis in mg/l Aluminium anzuzeigen.

#### Hinweise:

- Reinigen Sie die Fläschchen und den Messbecher vor dem Gebrauch mit Salzsäure (ca. 20 %). Spülen Sie beide gründlich mit deionisiertem Wasser.
- Um genaue Ergebnisse zu erhalten, muss die Probentemperatur zwischen 20 °C und 25 °C liegen.
- In Anwesenheit von Fluoriden und Polyphosphaten kann ein niedriges Testergebnis angezeigt werden. Die Einflüsse sind im Allgemeinen unbedeutend, es sei denn, dem Wasser wird künstlich Fluorid zugesetzt.

## AC4P27 Aluminium-Pulverpäckchen- und Flüssigreagenz-Test

#### **Eriochrom Cyanin-R-Methode**

0,01 – 0,25 mg/l Al

- 1. Laden Sie die Methode AC4P27, und führen Sie sie aus.
- 2. Verwenden Sie zwei saubere AQUAfast 24-mm-Rundfläschchen, Bestell-Nr. AC2V24, und markieren Sie eine als Blindwert.
- 3. Geben Sie 20 ml der Probe in ein 100-ml-Becherglas.
- 4. Geben Sie den Inhalt eines <u>Aluminum ECR F20 Powder Pack</u> direkt aus der Folienpackung zur Probe im Becherglas. Lösen Sie das Pulver mit einem sauberen Rührstab auf.
- 5. Geben Sie nach einer Reaktionszeit von <u>30 Minuten</u>.
- 6. den Inhalt eines <u>Hexamine F20 Powder Pack</u> direkt aus dem Folienbeutel zur selben Probe im Becherglas. Lösen Sie das Pulver mit einem sauberen Rührstab auf.
- Geben Sie 1 Tropfen <u>Aluminum ECR Masking Reagent</u> in das mit "Blindwert" markierte Gefäß. Geben Sie in das gleiche Fläschchen 10 ml der vorbereiteten Probe. (Dies ist das Blindprobengefäß).
- 8. Geben Sie die restlichen 10 ml der vorbereiteten Probe in das zweite Fläschchen (dies ist das Fläschchen).
- 9. Verschließen Sie die Fläschchen fest mit den Kappen, und mischen Sie den Inhalt durch mehrmaliges Schwenken oder Umdrehen. Wischen Sie die Außenseiten der Fläschchen ab.
- 10. Warten Sie, bis eine Reaktionszeit von 5 Minuten abgelaufen ist.
- 11. Stellen Sie das Blindprobengefäß in die Halterung in der Probenkammer, und schließen Sie die Probenkammertür.
- 12. Drücken Sie die Funktionstaste Blank (Blind), um den Blindwert zu messen.
- 13. Öffnen Sie den Probenkammerdeckel. Entnehmen Sie das Blindprobengefäß aus der Halterung.
- 14. Stellen Sie das Fläschchen in die Halterung in der Probenkammer, und schließen Sie die Probenkammertür.
- 15. Tippen Sie auf die Funktionstaste Sample (Probe), um das Ergebnis in mg/l Aluminium anzuzeigen.

#### Hinweise:

- Reinigen Sie die Fläschchen und den Messbecher vor dem Gebrauch mit Salzsäure (ca. 20 %). Spülen Sie beide gründlich mit deionisiertem Wasser.
- Um genaue Ergebnisse zu erhalten, muss die Probentemperatur zwischen 20 °C und 25 °C liegen.
- In Anwesenheit von Fluoriden und Polyphosphaten kann ein niedriges Testergebnis angezeigt werden. Die Einflüsse sind im Allgemeinen unbedeutend, es sei denn, dem Wasser wird künstlich Fluorid zugesetzt.
# AC2012 Ammoniak-Tabletten-Test

#### Indophenol-Blau-Methode

0,02 – 1 mg/I N (Ammoniak als Stickstoffträger)

- 1. Laden Sie die Methode AC2012, und führen Sie sie aus.
- Befüllen Sie eine saubere AQUAfast 24-mm-Rundküvette, Bestell-Nr. AC2V24, mit 10 ml der Probe. Verschließen Sie das Fläschchen fest mit dem Deckel. Wischen Sie die Außenseite des Fläschchens ab.
- 3. Stellen Sie das Fläschchen in die Halterung in der Probenkammer und schließen Sie die Probenkammertür.
- 4. Drücken Sie die Funktionstaste Blank (Blind), um den Blindwert zu messen.
- 5. Öffnen Sie die Probenkammertür, und nehmen Sie das Fläschchen aus der Halterung.
- 6. Geben Sie eine <u>Ammonia No. 1 Tablet</u> direkt aus dem Folienbeutel in das Fläschchen. Zerdrücken Sie die Tablette mit einem sauberen Rührstab.
- 7. Geben Sie eine <u>Ammonia No. 2 Tablet</u> direkt aus dem Folienbeutel in dasselbe Fläschchen. Zerdrücken Sie die Tablette mit einem sauberen Rührstab.
- Verschließen Sie das Fläschchen fest mit dem Deckel, und mischen Sie den Inhalt durch mehrmaliges Schwenken oder Umdrehen, bis sich die Tabletten aufgelöst haben.
   Wischen Sie die Außenseite des Fläschchens ab.
- 9. Warten Sie, bis eine Reaktionszeit von 10 Minuten abgelaufen ist.
- 10. Stellen Sie das Fläschchen in die Halterung in der Probenkammer und schließen Sie die Probenkammertür.
- 11. Tippen Sie auf die Funktionstaste Sample (Probe), um das Ergebnis in mg/l Ammoniak als Stickstoffträger anzuzeigen.

- Die Tabletten müssen in der richtigen Reihenfolge zugegeben werden.
- Die Ammoniaktablette (Ammonia No. 1) löst sich erst dann vollständig auf, wenn die Ammoniaktablette Nr. 2 (Ammonia No. 2) hinzugefügt wurde.
- Die Probentemperatur ist wichtig für eine volle Farbentwicklung. Bei einer Temperatur unter 20 °C beträgt die Reaktionszeit 15 Minuten.
- Umrechnung: mg/l NH<sub>4</sub> = mg/l N x 1,29 mg/l NH<sub>3</sub> = mg/l N x 1,22

# AC4P12 Ammoniak-Pulverpäckchen- und Flüssigreagenz-Test

#### Salicylat-Methode

0,01 - 0,8 mg/l N (Ammoniak als Stickstoffträger)

- 1. Laden Sie die Methode AC4P12, und führen Sie sie aus.
- 2. Verwenden Sie zwei saubere AQUAfast 24-mm-Rundfläschchen, Bestell-Nr. AC2V24.
- 3. Geben Sie 10 ml deionisiertes Wasser in das erste Fläschchen. (Dies ist das Blindprobengefäß.)
- 4. Geben Sie 10 ml der Probe in das zweite Fläschchen. (Dies ist das Fläschchen.)
- Geben Sie den Inhalt eines <u>Ammonia Salicylate F10 Powder Pack</u> direkt aus der Folienpackung in die Fläschchen. Verschließen Sie die Fläschchen fest mit den Kappen, und mischen Sie den Inhalt durch mehrmaliges Schwenken oder Umdrehen.
- 6. Warten Sie, bis eine Reaktionszeit von <u>3 Minuten</u> abgelaufen ist.
- Geben Sie den Inhalt eines <u>Ammonia Cyanurate F10 Powder Pack</u> direkt aus der Folienpackung zur Probe im Becherglas. Verschließen Sie die Fläschchen fest mit den Kappen, und mischen Sie den Inhalt durch mehrmaliges Schwenken oder Umdrehen. Wischen Sie die Außenseiten der Fläschchen ab.
- 8. Warten Sie, bis eine Reaktionszeit von 15 Minuten abgelaufen ist.
- 9. Stellen Sie das Blindprobengefäß in die Halterung in der Probenkammer, und schließen Sie die Probenkammertür.
- 10. Drücken Sie die Funktionstaste Blank (Blind), um den Blindwert zu messen.
- 11. Öffnen Sie den Probenkammerdeckel. Entnehmen Sie das Blindprobengefäß aus der Halterung.
- 12. Stellen Sie das Fläschchen in die Halterung in der Probenkammer, und schließen Sie die Probenkammertür.
- 13. Tippen Sie auf die Funktionstaste Sample (Probe), um das Ergebnis in mg/l Ammoniak als Stickstoffträger anzuzeigen.

#### Hinweise:

 Extrem basische oder saure Wasserproben sollten mit einer 0,5 mol/l (1 N) Schwefelsäurelösung oder 1 mol/l (1 N) Natriumhydroxidlösung auf pH 7 eingestellt werden.

| Störung                             | Störungsgrad und Korrekturen  |  |
|-------------------------------------|---|--|
| Kalzium                             | Mehr als 1.000 mg/l CaCO <sub>3</sub>   |  |
| Eisen                               | Störung in allen Konzentrationen. Bestimmen Sie zur Korrektur die Eisenkonzentration<br>in der Probe, indem Sie einen Gesamteisentest durchführen. Geben Sie die gleiche<br>Eisenkonzentration in das deionisierte Wasser (Schritt 3). Vorhandenes Eisen wird<br>erfolgreich ausgeblendet ("geblankt"). |  |
| Magnesium                           | Mehr als 6.000 mg/l CaCO <sub>3</sub>   |  |
| Nitrat                              | Mehr als 100 mg/l NO <sub>3</sub> -N  |  |
| Nitrit                              | Mehr als 12 mg/l NO <sub>2</sub> -N   |  |
| Phosphat                            | Mehr als 100 mg/l PO <sub>4</sub> -P  |  |
| Sulfat                              | Mehr als 300 mg/l SO <sub>4</sub>   |  |
| Sulfid                              | Intensiviert die Farbe  |  |
| Glycin, Hydrazin,<br>Farbe, Trübung | Weniger häufige Störungen wie Hydrazin und Glycin führen zu Farbverstärkungen in der vorbereiteten Probe. Es resultieren überhöhte Ergebnisse für Trübung und Farbe.<br>Samples (Proben) mit starken Störfaktoren müssen destilliert werden.  |  |

# ACR012 Ammoniak-Reaktionsröhrchen-Test für niedrigen Bereich

#### Salicylat-Methode

0,02 - 2,5 mg/l N (Ammoniak als Stickstoffträger)

- 1. Laden Sie die Methode ACR012, und führen Sie sie aus.
- 2. Öffnen Sie ein 16-mm-Reaktionsgefäß und geben Sie 2 ml deionisiertes Wasser dazu. (Dies ist das Blindprobengefäß.)
- 3. Öffnen Sie ein zweites 16-mm-Reaktionsgefäß und geben Sie 2 ml der Probe hinzu. (Dies ist das Fläschchen.)
- 4. Geben Sie den Inhalt eines <u>Ammonia Salicylate F5 Powder Pack</u> direkt aus der Folienpackung in die Fläschchen.
- 5. Geben Sie den Inhalt eines <u>Ammonia Cyanurate F5 Powder Pack</u> direkt aus der Folienpackung in die Fläschchen.
- 6. Verschließen Sie die Fläschchen fest mit den Kappen, und mischen Sie den Inhalt durch mehrmaliges Schwenken oder Umdrehen. Wischen Sie die Außenseiten der Fläschchen ab.
- 7. Warten Sie, bis eine Reaktionszeit von 20 Minuten abgelaufen ist.
- 8. Stellen Sie das Blindprobengefäß in die Halterung in der Probenkammer, und schließen Sie die Probenkammertür.
- 9. Drücken Sie die Funktionstaste Blank (Blind), um den Blindwert zu messen.
- 10. Öffnen Sie den Probenkammertür. Entnehmen Sie das Blindprobengefäß aus der Halterung.
- 11. Stellen Sie das Fläschchen in die Halterung in der Probenkammer, und schließen Sie die Probenkammertür.
- 12. Tippen Sie auf die Funktionstaste Sample (Probe), um das Ergebnis in mg/I Ammoniak als Stickstoffträger anzuzeigen.

- Stark alkalische oder saure Wasserproben müssen vor der Analyse auf etwa pH 7 eingestellt werden (1 mol/l Salzsäure bzw. 1 mol/l Natriumhydroxid verwenden).
- Wenn die Anwesenheit von Chlor bekannt ist, geben Sie einen Tropfen 0,1 mol/I Natriumthiosulfat f
  ür jeweils 0,3 mg/I Cl<sub>2</sub> zu einer 1-Liter-Wasserprobe hinzu.
- Eisen beeinflusst die Testergebnisse. Die Störungen können folgendermaßen eliminiert werden: Bestimmen Sie die in der Wasserprobe vorhandene Gesamteisenmenge. Zur Herstellung der Blindprobe wird anstelle von deionisiertem Wasser eine Eisen-Standardlösung mit gleicher Eisenkonzentration in das Fläschchen gegeben.
- Umrechnung:  $mg/I NH_4 = mg/I N \times 1,29$  $mg/I NH_3 = mg/I N \times 1,22$

# ACR011 Ammoniak-Reaktionsröhrchen-Test für hohen Bereich

#### Salicylat-Methode

- 1 50 mg/I N (Ammoniak als Stickstoffträger)
- 1. Laden Sie die Methode ACR011, und führen Sie sie aus.
- 2. Öffnen Sie ein 16-mm-Reaktionsgefäß und geben Sie 0,1 ml deionisiertes Wasser dazu. (Dies ist das Blindprobengefäß.)
- 3. Öffnen Sie ein zweites 16-mm-Reaktionsgefäß und geben Sie 0,1 ml der Probe hinzu. (Dies ist das Fläschchen.)
- 4. Geben Sie den Inhalt eines <u>Ammonia Salicylate F5 Powder Pack</u> direkt aus der Folienpackung in die Fläschchen.
- 5. Geben Sie den Inhalt eines <u>Ammonia Cyanurate F5 Powder Pack</u> direkt aus der Folienpackung direkt in die Reaktionsgefäße.
- 6. Verschließen Sie die Fläschchen fest mit den Kappen, und mischen Sie den Inhalt durch mehrmaliges Schwenken oder Umdrehen. Wischen Sie die Außenseiten der Fläschchen ab.
- 7. Warten Sie, bis eine Reaktionszeit von 20 Minuten abgelaufen ist.
- 8. Stellen Sie das Blindprobengefäß in die Halterung in der Probenkammer, und schließen Sie die Probenkammertür.
- 9. Drücken Sie die Funktionstaste Blank (Blind), um den Blindwert zu messen.
- 10. Öffnen Sie den Probenkammertür. Entnehmen Sie das Blindprobengefäß aus der Halterung.
- 11. Stellen Sie das Fläschchen in die Halterung in der Probenkammer, und schließen Sie die Probenkammertür.
- 12. Tippen Sie auf die Funktionstaste Sample (Probe), um das Ergebnis in mg/l Ammoniak als Stickstoffträger anzuzeigen.

- Stark alkalische oder saure Wasserproben müssen vor der Analyse auf etwa pH 7 eingestellt werden (1 mol/l Salzsäure bzw. 1 mol/l Natriumhydroxid verwenden).
- Wenn die Anwesenheit von Chlor bekannt ist, geben Sie einen Tropfen 0,1 mol/I Natriumthiosulfat f
  ür jeweils 0,3 mg/I Cl<sub>2</sub> zu einer 1-Liter-Wasserprobe hinzu.
- Eisen beeinflusst die Testergebnisse. Die Interferenzen können folgendermaßen eliminiert werden: Bestimmen Sie die in der Wasserprobe vorhandene Gesamteisenmenge. Zur Herstellung der Blindprobe wird anstelle von deionisiertem Wasser eine Eisen-Standardlösung mit gleicher Eisenkonzentration in das Fläschchen gegeben.
- Umrechnung:  $mg/I NH_4 = mg/I N \times 1,29$  $mg/I NH_3 = mg/I N \times 1,22$

# AC2035 Bromtabletten-Test

#### DPD-Methode

0,05 – 13 mg/l Br<sub>2</sub>

- 1. Laden Sie die Methode AC203524, und führen Sie sie aus.
- Befüllen Sie eine saubere AQUAfast 24-mm-Rundküvette, Bestell-Nr. AC2V24, mit 10 ml der Probe. Verschließen Sie das Fläschchen fest mit dem Deckel. Wischen Sie die Außenseite des Fläschchens ab.
- 3. Stellen Sie das Fläschchen in die Halterung in der Probenkammer und schließen Sie die Probenkammertür.
- 4. Drücken Sie die Funktionstaste Blank (Blind), um den Blindwert zu messen.
- 5. Öffnen Sie die Probenkammertür, und nehmen Sie das Fläschchen aus der Halterung.
- 6. Entleeren Sie das Fläschchen, so dass einige Tropfen der Probe im Fläschchen zurückbleiben.
- 7. Geben Sie eine <u>DPD No. 1 Tablet</u> direkt aus dem Folienbeutel in das Fläschchen. Zerdrücken Sie die Tablette mit einem sauberen Rührstab.
- 8. Füllen Sie die Probe bis zur 10-ml-Marke auf dem Fläschchen auf.
- Verschließen Sie das Fläschchen fest mit dem Deckel, und mischen Sie den Inhalt durch mehrmaliges Schwenken oder Umdrehen, bis sich die Tablette aufgelöst hat. Wischen Sie die Außenseite des Fläschchens ab.
- 10. Stellen Sie das Fläschchen in die Halterung in der Probenkammer und schließen Sie die Probenkammertür.
- 11. Tippen Sie auf die Funktionstaste Sample (Probe), um das Ergebnis in mg/l Brom anzuzeigen.

#### Hinweise:

- Alternativ kann die AC203510-Methode mit quadratischen 10-mm-Fläschchen und die AC203550-Methode mit rechteckigen 50-mm-Fläschchen verwendet werden.
   Alle Blindwert- und Probenvolumina müssen mit den in dieser Anleitung angegebenen Volumina übereinstimmen, so dass die Proben möglicherweise in separaten Behältern vorbereitet und dann in das ausgewählte Fläschchen überführt werden müssen.
- Reinigung der Fläschchen: Da viele Haushaltsreiniger (z. B. Geschirrspülmittel) reduzierende Stoffe enthalten, kann es nach deren Anwendung bei der Bestimmung von Brom zu Minderbefunden kommen. Um diesen Messfehler auszuschließen, sollten die Glasgefäße chlorzehrungsfrei sein.

- Probenvorbereitung: Bei der Vorbereitung der Probe muss das Entweichen von Bromgasen,
   z. B. beim Pipettieren oder Schütteln, vermieden werden. Die Analyse muss unmittelbar
   nach der Probenahme erfolgen.
- Die DPD-Farbentwicklung erfolgt bei einem pH-Wert von 6,2 bis 6,5. Die Reagenztablette enthält daher einen Puffer zur pH-Einstellung. Stark alkalische oder saure Wasserproben müssen vor der Analyse auf einen pH-Wert zwischen 6 und 7 gebracht werden (dazu 0,5 mol/l Schwefelsäure bzw. 1 mol/l Natronlauge verwenden).
- Überschreitung des Messbereichs: Konzentrationen von mehr als 22 mg/l Brom können zu Ergebnissen bis hin zu 0 mg/l führen. In diesem Fall muss die Wasserprobe mit bromfreiem Wasser verdünnt werden. Dazu sollten 10 ml der verdünnten Probe mit dem Reagenz versetzt und die Messung wiederholt werden.
- Oxidationsmittel wie Chlor oder Ozon sind Störfaktoren, da sie auf die gleiche Weise wie Brom reagieren.

# AC2017 Chlorid-Tabletten-Test

#### Silbernitrat/Trübungsmethode

0,5 – 25 mg/l Cl

- 1. Laden Sie die Methode AC2017, und führen Sie sie aus.
- Befüllen Sie eine saubere AQUAfast 24-mm-Rundküvette, Bestell-Nr. AC2V24, mit 10 ml der Probe. Verschließen Sie das Fläschchen fest mit dem Deckel. Wischen Sie die Außenseite des Fläschchens ab.
- 3. Stellen Sie das Fläschchen in die Halterung in der Probenkammer und schließen Sie die Probenkammertür.
- 4. Drücken Sie die Funktionstaste Blank (Blind), um den Blindwert zu messen.
- 5. Öffnen Sie die Probenkammertür, und nehmen Sie das Fläschchen aus der Halterung.
- 6. Geben Sie eine <u>Chloride T1 Tablet</u> direkt aus dem Folienbeutel in das Fläschchen. Zerdrücken Sie die Tablette mit einem sauberen Rührstab.
- 7. Geben Sie eine <u>Chloride T2 Tablet</u> direkt aus dem Folienbeutel in dasselbe Fläschchen. Zerdrücken Sie die Tablette mit einem sauberen Rührstab.
- 8. Verschließen Sie das Fläschchen fest mit dem Deckel, und mischen Sie den Inhalt durch vorsichtiges Schwenken, bis sich die Tablette aufgelöst hat. Wischen Sie die Außenseite des Fläschchens ab.
- 9. Warten Sie, bis eine Reaktionszeit von 2 Minuten abgelaufen ist.
- 10. Stellen Sie das Fläschchen in die Halterung in der Probenkammer und schließen Sie die Probenkammertür.
- 11. Tippen Sie auf die Funktionstaste Sample (Probe), um das Ergebnis in mg/l Chlorid anzuzeigen.

- Vergewissern Sie sich, dass alle Tablettenpartikel aufgelöst sind. Chlorid verursacht eine extrem fein verteilte Trübung mit milchigem Aussehen. Starkes Schütteln führt zur Bildung größerer Partikel, die zu falschen Messwerten führen können.
- Hohe Konzentrationen von Elektrolyten und organischen Verbindungen haben unterschiedliche Auswirkungen auf die Fällungsreaktionen.
- Ionen, die sich mit Silbernitrat in sauren Medien ebenfalls ablagern, wie Bromide, Iodide und Thiocyanate, beeinflussen die Analyse.
- Stark alkalisches Wasser sollte vor der Analyse erforderlichenfalls mit Salpetersäure neutralisiert werden.

## AC2070 Chlortabletten-Test (freies Chlor und Gesamtchlor) DPD-Methode

0,01 - 6 mg/l Cl<sub>2</sub>

- 1. Laden Sie die Methode AC207024, und führen Sie sie aus.
- Befüllen Sie eine saubere AQUAfast 24-mm-Rundküvette, Bestell-Nr. AC2V24, mit 10 ml der Probe. Verschließen Sie das Fläschchen fest mit dem Deckel. Wischen Sie die Außenseite des Fläschchens ab.
- 3. Stellen Sie das Fläschchen in die Halterung in der Probenkammer und schließen Sie die Probenkammertür.
- 4. Drücken Sie die Funktionstaste Blank (Blind), um den Blindwert zu messen.
- 5. Öffnen Sie die Probenkammertür, und nehmen Sie das Fläschchen aus der Halterung.
- 6. Entleeren Sie das Fläschchen, so dass einige Tropfen der Probe im Fläschchen zurückbleiben.
- 7. Geben Sie eine <u>DPD No. 1 Tablet</u> direkt aus dem Folienbeutel in das Fläschchen. Zerdrücken Sie die Tablette mit einem sauberen Rührstab.
- 8. Füllen Sie die Probe bis zur 10-ml-Marke auf dem Fläschchen auf.
- Verschließen Sie das Fläschchen fest mit dem Deckel, und mischen Sie den Inhalt durch mehrmaliges Schwenken oder Umdrehen, bis sich die Tablette aufgelöst hat. Wischen Sie die Außenseite des Fläschchens ab.
- 10. Stellen Sie das Fläschchen in die Halterung in der Probenkammer und schließen Sie die Probenkammertür.
- 11. Drücken Sie die Funktionstaste Sample (Probe), um das Ergebnis in mg/l freies Chlor anzuzeigen.
- 12. Öffnen Sie die Probenkammertür, und nehmen Sie das Fläschchen aus der Halterung.
- 13. Geben Sie eine <u>DPD No. 3 Tablet</u> direkt aus dem Folienbeutel in das Fläschchen. Zerdrücken Sie die Tablette mit einem sauberen Rührstab.
- Verschließen Sie das Fläschchen fest mit dem Deckel, und mischen Sie den Inhalt durch mehrmaliges Schwenken oder Umdrehen, bis sich die Tablette aufgelöst hat. Wischen Sie die Außenseite des Fläschchens ab.
- 15. Warten Sie, bis eine Reaktionszeit von 2 Minuten abgelaufen ist.
- 16. Stellen Sie das Fläschchen in die Halterung in der Probenkammer und schließen Sie die Probenkammertür.
- 17. Drücken Sie die Funktionstaste Sample (Probe), um das Ergebnis in mg/l Gesamtchlor zu anzuzeigen.

#### Hinweise:

- Alternativ kann die AC207010-Methode mit quadratischen 10-mm-Fläschchen und die AC207050-Methode mit rechteckigen 50-mm-Fläschchen verwendet werden. Alle Blindwertund Probenvolumina müssen mit den in dieser Anleitung angegebenen Volumina übereinstimmen, so dass die Proben möglicherweise in separaten Behältern vorbereitet und dann in das ausgewählte Fläschchen überführt werden müssen.
- Reinigung der Fläschchen: Da viele Haushaltsreiniger (z. B. Geschirrspülmittel) reduzierende Stoffe enthalten, kann es nach deren Anwendung bei der Bestimmung von Chlor zu Minderbefunden kommen. Um diesen Messfehler auszuschließen, sollten die Glasgefäße chlorzehrungsfrei sein.

- Für die Einzelbestimmung von freiem Chlor und Gesamtchlor wird empfohlen, jeweils einen eigenen Satz an Glasgefäßen zu verwenden (EN ISO 7393-2, 5.3).
- Probenvorbereitung: Bei der Vorbereitung der Probe ist das Entweichen von Chlorgasen (z. B. beim Pipettieren oder Schütteln) zu vermeiden. Die Analyse muss unmittelbar nach der Probenahme erfolgen.
- Die DPD-Farbentwicklung erfolgt bei einem pH-Wert von 6,2 bis 6,5. Die Reagenzien enthalten daher einen Puffer zur pH-Einstellung. Stark alkalische oder saure Wasserproben müssen vor der Analyse auf einen pH-Wert zwischen 6 und 7 gebracht werden (dazu 0,5 mol/I Schwefelsäure bzw. 1 mol/I Natronlauge verwenden).
- Überschreitung des Messbereichs: Konzentrationen von mehr als 10 mg/l Chlor können zu Ergebnissen bis hin zu 0 mg/l führen. In diesem Fall muss die Wasserprobe mit chlorfreiem Wasser verdünnt werden. Dazu sollten 10 ml der verdünnten Probe mit dem Reagenz versetzt und die Messung wiederholt werden.
- Trübungen können zu Fehlmessungen führen. Die Verwendung der DPD-Tablette Nr. 1 in Proben mit hohem Calciumionengehalt und/oder hoher Leitfähigkeit kann zu einer Eintrübung der Probe und damit verbundenen Fehlmessungen führen.
- Oxidationsmittel wie Brom oder Ozon sind Störfaktoren, da sie wie Chlor reagieren.

# AC2071 Chlortabletten-Test (freies Chlor)

#### DPD-Methode

0,01 - 6 mg/l Cl<sub>2</sub>

- 1. Laden Sie die Methode AC207124, und führen Sie sie aus.
- Befüllen Sie eine saubere AQUAfast 24-mm-Rundküvette, Bestell-Nr. AC2V24, mit 10 ml der Probe. Verschließen Sie das Fläschchen fest mit dem Deckel. Wischen Sie die Außenseite des Fläschchens ab.
- 3. Stellen Sie das Fläschchen in die Halterung in der Probenkammer und schließen Sie die Probenkammertür.
- 4. Drücken Sie die Funktionstaste Blank (Blind), um den Blindwert zu messen.
- 5. Öffnen Sie die Probenkammertür, und nehmen Sie das Fläschchen aus der Halterung.
- 6. Entleeren Sie das Fläschchen, so dass einige Tropfen der Probe im Fläschchen zurückbleiben.
- 7. Geben Sie eine <u>DPD No. 1 Tablet</u> direkt aus dem Folienbeutel in das Fläschchen. Zerdrücken Sie die Tablette mit einem sauberen Rührstab.
- 8. Füllen Sie die Probe bis zur 10-ml-Marke auf dem Fläschchen auf.
- Verschließen Sie das Fläschchen fest mit dem Deckel, und mischen Sie den Inhalt durch mehrmaliges Schwenken oder Umdrehen, bis sich die Tablette aufgelöst hat. Wischen Sie die Außenseite des Fläschchens ab.
- 10. Stellen Sie das Fläschchen in die Halterung in der Probenkammer und schließen Sie die Probenkammertür.
- 11. Drücken Sie die Funktionstaste Sample (Probe), um das Ergebnis in mg/l freies Chlor anzuzeigen.

#### Hinweise:

- Alternativ kann die AC207110-Methode mit quadratischen 10-mm-Fläschchen und die AC207150-Methode mit rechteckigen 50-mm-Fläschchen verwendet werden. Alle Blindwertund Probenvolumina müssen mit den in dieser Anleitung angegebenen Volumina übereinstimmen, so dass die Proben möglicherweise in separaten Behältern vorbereitet und dann in das ausgewählte Fläschchen überführt werden müssen.
- Reinigung der Fläschchen: Da viele Haushaltsreiniger (z. B. Geschirrspülmittel) reduzierende Stoffe enthalten, kann es nach deren Anwendung bei der Bestimmung von Chlor zu Minderbefunden kommen. Um diesen Messfehler auszuschließen, sollten die Glasgefäße chlorzehrungsfrei sein.

- Für die Einzelbestimmung von freiem Chlor und Gesamtchlor wird empfohlen, jeweils einen eigenen Satz an Glasgefäßen zu verwenden (EN ISO 7393-2, 5.3).
- Probenvorbereitung: Bei der Vorbereitung der Probe ist das Entweichen von Chlorgasen (z. B. beim Pipettieren oder Schütteln) zu vermeiden. Die Analyse muss unmittelbar nach der Probenahme erfolgen.
- Die DPD-Farbentwicklung erfolgt bei einem pH-Wert von 6,2 bis 6,5. Die Reagenzien enthalten daher einen Puffer zur pH-Einstellung. Stark alkalische oder saure Wasserproben müssen vor der Analyse auf einen pH-Wert zwischen 6 und 7 gebracht werden (dazu 0,5 mol/I Schwefelsäure bzw. 1 mol/I Natronlauge verwenden).
- Überschreitung des Messbereichs: Konzentrationen von mehr als 10 mg/l Chlor können zu Ergebnissen bis hin zu 0 mg/l führen. In diesem Fall muss die Wasserprobe mit chlorfreiem Wasser verdünnt werden. Dazu sollten 10 ml der verdünnten Probe mit dem Reagenz versetzt und die Messung wiederholt werden.
- Trübungen können zu Fehlmessungen führen. Die Verwendung der DPD-Tablette Nr. 1 in Proben mit hohem Calciumionengehalt und/oder hoher Leitfähigkeit kann zu einer Eintrübung der Probe und damit verbundenen Fehlmessungen führen.
- Oxidationsmittel wie Brom oder Ozon sind Störfaktoren, da sie wie Chlor reagieren.

## AC2072 Chlortabletten-Test (Gesamtchlor)

#### DPD-Methode

0,01 - 6 mg/l Cl<sub>2</sub>

- 1. Laden Sie die Methode AC207224, und führen Sie sie aus.
- Befüllen Sie eine saubere AQUAfast 24-mm-Rundküvette, Bestell-Nr. AC2V24, mit 10 ml der Probe. Verschließen Sie das Fläschchen fest mit dem Deckel. Wischen Sie die Außenseite des Fläschchens ab.
- Stellen Sie das Fläschchen in die Halterung in der Probenkammer und schließen Sie die Probenkammert
  ür.
- 4. Drücken Sie die Funktionstaste Blank (Blind), um den Blindwert zu messen.
- 5. Öffnen Sie die Probenkammertür, und nehmen Sie das Fläschchen aus der Halterung.
- 6. Entleeren Sie das Fläschchen, so dass einige Tropfen der Probe im Fläschchen zurückbleiben.
- Eine <u>DPD No. 4 Tablet</u> (oder eine <u>DPD No. 1 Tablet</u> und eine <u>DPD No. 3 Tablet</u>) direkt aus der Folienpackung in das Fläschchen geben. Zerdrücken Sie die Tablette mit einem sauberen Rührstab.
- 8. Füllen Sie die Probe bis zur 10-ml-Marke auf dem Fläschchen auf.
- Verschließen Sie das Fläschchen fest mit dem Deckel, und mischen Sie den Inhalt durch mehrmaliges Schwenken oder Umdrehen, bis sich die Tablette aufgelöst hat. Wischen Sie die Außenseite des Fläschchens ab.
- 10. Warten Sie, bis eine Reaktionszeit von 2 Minuten abgelaufen ist.
- 11. Stellen Sie das Fläschchen in die Halterung in der Probenkammer und schließen Sie die Probenkammertür.
- 12. Drücken Sie die Funktionstaste Sample (Probe), um das Ergebnis in mg/l Gesamtchlor zu anzuzeigen.

- Alternativ kann die AC207210-Methode mit quadratischen 10-mm-Fläschchen und die AC207250-Methode mit rechteckigen 50-mm-Fläschchen verwendet werden. Alle Blindwertund Probenvolumina müssen mit den in dieser Anleitung angegebenen Volumina übereinstimmen, so dass die Proben möglicherweise in separaten Behältern vorbereitet und dann in das ausgewählte Fläschchen überführt werden müssen.
- Reinigung der Fläschchen: Da viele Haushaltsreiniger (z. B. Geschirrspülmittel) reduzierende Stoffe enthalten, kann es nach deren Anwendung bei der Bestimmung von Chlor zu Minderbefunden kommen. Um diesen Messfehler auszuschließen, sollten die Glasgefäße chlorzehrungsfrei sein.

- Für die Einzelbestimmung von freiem Chlor und Gesamtchlor wird empfohlen, jeweils einen eigenen Satz an Glasgefäßen zu verwenden (EN ISO 7393-2, 5.3).
- Probenvorbereitung: Bei der Vorbereitung der Probe ist das Entweichen von Chlorgasen (z. B. beim Pipettieren oder Schütteln) zu vermeiden. Die Analyse muss unmittelbar nach der Probenahme erfolgen.
- Die DPD-Farbentwicklung erfolgt bei einem pH-Wert von 6,2 bis 6,5. Die Reagenzien enthalten daher einen Puffer zur pH-Einstellung. Stark alkalische oder saure Wasserproben müssen vor der Analyse auf einen pH-Wert zwischen 6 und 7 gebracht werden (dazu 0,5 mol/I Schwefelsäure bzw. 1 mol/I Natronlauge verwenden).
- Überschreitung des Messbereichs: Konzentrationen von mehr als 10 mg/l Chlor können zu Ergebnissen bis hin zu 0 mg/l führen. In diesem Fall muss die Wasserprobe mit chlorfreiem Wasser verdünnt werden. Dazu sollten 10 ml der verdünnten Probe mit dem Reagenz versetzt und die Messung wiederholt werden.
- Trübungen können zu Fehlmessungen führen. Die Verwendung der DPD-Tablette Nr. 1 in Proben mit hohem Calciumionengehalt und/oder hoher Leitfähigkeit kann zu einer Eintrübung der Probe und damit verbundenen Fehlmessungen führen.
- Oxidationsmittel wie Brom oder Ozon sind Störfaktoren, da sie wie Chlor reagieren.

# AC4P71 Chlorpulverpäckchen-Test (freies Chlor)

#### DPD-Methode

0,02 - 2 mg/l Cl<sub>2</sub>

- 1. Laden Sie die Methode AC4P71, und führen Sie sie aus.
- Befüllen Sie eine saubere AQUAfast 24-mm-Rundküvette, Bestell-Nr. AC2V24, mit 10 ml der Probe. Verschließen Sie das Fläschchen fest mit dem Deckel. Wischen Sie die Außenseite des Fläschchens ab.
- 3. Stellen Sie das Fläschchen in die Halterung in der Probenkammer und schließen Sie die Probenkammertür.
- 4. Drücken Sie die Funktionstaste Blank (Blind), um den Blindwert zu messen.
- 5. Öffnen Sie die Probenkammertür, und nehmen Sie das Fläschchen aus der Halterung.
- 6. Geben Sie den Inhalt eines <u>Chlorine Free-DPD / F10 Powder Pack</u> direkt aus der Folienpackung in das Fläschchen.
- Verschließen Sie das Fläschchen fest mit dem Deckel, und mischen Sie den Inhalt (für ca. 20 Sekunden) durch mehrmaliges Schwenken oder Umdrehen. Wischen Sie die Außenseite des Fläschchens ab.
- 8. Stellen Sie das Fläschchen in die Halterung in der Probenkammer und schließen Sie die Probenkammertür.
- 9. Drücken Sie die Funktionstaste Sample (Probe), um das Ergebnis in mg/l freies Chlor anzuzeigen.

#### Hinweise:

 Reinigung der Fläschchen: Da viele Haushaltsreiniger (z. B. Geschirrspülmittel) reduzierende Stoffe enthalten, kann es nach deren Anwendung bei der Bestimmung von Chlor zu Minderbefunden kommen. Um diesen Messfehler auszuschließen, sollten die Glasgefäße chlorzehrungsfrei sein.

- Für die Einzelbestimmung von freiem Chlor und Gesamtchlor wird empfohlen, jeweils einen eigenen Satz an Glasgefäßen zu verwenden (EN ISO 7393-2, 5.3).
- Probenvorbereitung: Bei der Vorbereitung der Probe ist das Entweichen von Chlorgasen (z. B. beim Pipettieren oder Schütteln) zu vermeiden. Die Analyse muss unmittelbar nach der Probenahme erfolgen.
- Die DPD-Farbentwicklung erfolgt bei einem pH-Wert von 6,2 bis 6,5. Die Reagenzien enthalten daher einen Puffer zur pH-Einstellung. Stark alkalische oder saure Wasserproben müssen vor der Analyse auf einen pH-Wert zwischen 6 und 7 gebracht werden (dazu 0,5 mol/I Schwefelsäure bzw. 1 mol/I Natronlauge verwenden).
- Überschreitung des Messbereichs: Konzentrationen von mehr als 2 mg/l Chlor (durch Pulverpäckchen) können zu Ergebnissen bis hin zu 0 mg/l führen. In diesem Fall muss die Wasserprobe mit chlorfreiem Wasser verdünnt werden. Dazu sollten 10 ml der verdünnten Probe mit dem Reagenz versetzt und die Messung wiederholt werden.
- Oxidationsmittel wie Brom oder Ozon sind Störfaktoren, da sie wie Chlor reagieren.

# AC4P72 Chlorpulverpäckchen-Test (Gesamtchlor)

#### DPD-Methode

0,02 – 2 mg/l Cl<sub>2</sub>

- 1. Laden Sie die Methode AC4P72, und führen Sie sie aus.
- Befüllen Sie eine saubere AQUAfast 24-mm-Rundküvette, Bestell-Nr. AC2V24, mit 10 ml der Probe. Verschließen Sie das Fläschchen fest mit dem Deckel. Wischen Sie die Außenseite des Fläschchens ab.
- 3. Stellen Sie das Fläschchen in die Halterung in der Probenkammer und schließen Sie die Probenkammertür.
- 4. Drücken Sie die Funktionstaste Blank (Blind), um den Blindwert zu messen.
- 5. Öffnen Sie die Probenkammertür, und nehmen Sie das Fläschchen aus der Halterung.
- 6. Geben Sie den Inhalt eines <u>Chlorine Total-DPD / F10 Powder Pack (Gesamtchlor)</u> direkt aus der Folienpackung in das Fläschchen.
- Verschließen Sie das Fläschchen fest mit dem Deckel, und mischen Sie den Inhalt (für ca. 20 Sekunden) durch mehrmaliges Schwenken oder Umdrehen. Wischen Sie die Außenseite des Fläschchens ab.
- 8. Warten Sie, bis eine Reaktionszeit von <u>3 Minuten</u> abgelaufen ist.
- 9. Stellen Sie das Fläschchen in die Halterung in der Probenkammer und schließen Sie die Probenkammertür.
- 10. Drücken Sie die Funktionstaste Sample (Probe), um das Ergebnis in mg/l Gesamtchlor zu anzuzeigen.

#### Hinweise:

 Reinigung der Fläschchen: Da viele Haushaltsreiniger (z. B. Geschirrspülmittel) reduzierende Stoffe enthalten, kann es nach deren Anwendung bei der Bestimmung von Chlor zu Minderbefunden kommen. Um diesen Messfehler auszuschließen, sollten die Glasgefäße chlorzehrungsfrei sein.

- Für die Einzelbestimmung von freiem Chlor und Gesamtchlor wird empfohlen, jeweils einen eigenen Satz an Glasgefäßen zu verwenden (EN ISO 7393-2, 5.3).
- Probenvorbereitung: Bei der Vorbereitung der Probe ist das Entweichen von Chlorgasen (z. B. beim Pipettieren oder Schütteln) zu vermeiden. Die Analyse muss unmittelbar nach der Probenahme erfolgen.
- Die DPD-Farbentwicklung erfolgt bei einem pH-Wert von 6,2 bis 6,5. Die Reagenzien enthalten daher einen Puffer zur pH-Einstellung. Stark alkalische oder saure Wasserproben müssen vor der Analyse auf einen pH-Wert zwischen 6 und 7 gebracht werden (dazu 0,5 mol/l Schwefelsäure bzw. 1 mol/l Natronlauge verwenden).
- Überschreitung des Messbereichs: Konzentrationen von mehr als 2 mg/l Chlor (durch Pulverpäckchen) können zu Ergebnissen bis hin zu 0 mg/l führen. In diesem Fall muss die Wasserprobe mit chlorfreiem Wasser verdünnt werden. Dazu sollten 10 ml der verdünnten Probe mit dem Reagenz versetzt und die Messung wiederholt werden.
- Oxidationsmittel wie Brom oder Ozon sind Störfaktoren, da sie wie Chlor reagieren.

# AC3072 Chlortabletten-Test (Gesamtchlor) für hohen Bereich

#### KI-/Säure-Methode

- 5 200 mg/l Cl<sub>2</sub>
- 1. Laden Sie die Methode AC3072, und führen Sie sie aus.
- Befüllen Sie eine saubere AQUAfast 16-mm-Rundküvette, Bestell-Nr. AC2V16, mit 10 ml der Probe. Verschließen Sie das Fläschchen fest mit dem Deckel. Wischen Sie die Außenseite des Fläschchens ab.
- 3. Stellen Sie das Fläschchen in die Halterung in der Probenkammer und schließen Sie die Probenkammertür.
- 4. Drücken Sie die Funktionstaste Blank (Blind), um den Blindwert zu messen.
- 5. Öffnen Sie die Probenkammertür, und nehmen Sie das Fläschchen aus der Halterung.
- 6. Geben Sie eine <u>Chlorine HR (KI) Tablet</u> direkt aus dem Folienbeutel in das Fläschchen. Zerdrücken Sie die Tablette mit einem sauberen Rührstab.
- 7. Geben Sie eine <u>Acidifying GP Tablet</u> direkt aus dem Folienbeutel in dasselbe Fläschchen. Zerdrücken Sie die Tablette mit einem sauberen Rührstab.
- Verschließen Sie das Fläschchen fest mit dem Deckel, und mischen Sie den Inhalt durch mehrmaliges Schwenken oder Umdrehen, bis sich die Tabletten aufgelöst haben. Wischen Sie die Außenseite des Fläschchens ab.
- Stellen Sie das Fläschchen in die Halterung in der Probenkammer und schließen Sie die Probenkammert
  ür.
- 10. Tippen Sie auf die Funktionstaste Sample (Probe), um das Ergebnis in mg/l Chlor anzuzeigen.

#### Hinweise:

• Oxidationsmittel sind Störfaktoren, da sie wie Chlor reagieren.

# AC2099 Chlordioxid-Tabletten-Test

#### DPD-Methode

0,02 - 11 mg/l ClO<sub>2</sub>

#### Chlordioxidmessung in Abwesenheit von Chlor

- 1. Laden Sie die Methode AC209924, und führen Sie sie aus.
- Befüllen Sie eine saubere AQUAfast 24-mm-Rundküvette, Bestell-Nr. AC2V24, mit 10 ml der Probe. Verschließen Sie das Fläschchen fest mit dem Deckel. Wischen Sie die Außenseite des Fläschchens ab.
- 3. Stellen Sie das Fläschchen in die Halterung in der Probenkammer und schließen Sie die Probenkammertür.
- 4. Drücken Sie die Funktionstaste Blank (Blind), um den Blindwert zu messen.
- 5. Öffnen Sie die Probenkammertür, und nehmen Sie das Fläschchen aus der Halterung.
- 6. Entleeren Sie das Fläschchen, so dass einige Tropfen der Probe im Fläschchen zurückbleiben.
- 7. Geben Sie eine <u>DPD No. 1 Tablet</u> direkt aus dem Folienbeutel in das Fläschchen. Zerdrücken Sie die Tablette mit einem sauberen Rührstab.
- 8. Füllen Sie die Probe bis zur 10-ml-Marke auf dem Fläschchen auf.
- Verschließen Sie das Fläschchen fest mit dem Deckel, und mischen Sie den Inhalt durch mehrmaliges Schwenken oder Umdrehen, bis sich die Tablette aufgelöst hat. Wischen Sie die Außenseite des Fläschchens ab.
- 10. Stellen Sie das Fläschchen in die Halterung in der Probenkammer und schließen Sie die Probenkammertür.
- 11. Tippen Sie auf die Funktionstaste Sample (Probe), um das Ergebnis in mg/l Chlordioxid anzuzeigen.

#### Chlordioxidmessung in Anwesenheit von Chlor

- 1. Laden Sie die Methode AC209924, und führen Sie sie aus.
- Befüllen Sie eine saubere AQUAfast 24-mm-Rundküvette, Bestell-Nr. AC2V24, mit 10 ml der Probe. Verschließen Sie das Fläschchen fest mit dem Deckel. Wischen Sie die Außenseite des Fläschchens ab.
- 3. Stellen Sie das Fläschchen in die Halterung in der Probenkammer und schließen Sie die Probenkammertür.
- 4. Drücken Sie die Funktionstaste Blank (Blind), um den Blindwert zu messen.
- 5. Öffnen Sie die Probenkammertür, und nehmen Sie das Fläschchen aus der Halterung.
- 6. Entleeren Sie das Fläschchen, so dass einige Tropfen der Probe im Fläschchen zurückbleiben.
- 7. Geben Sie eine <u>DPD No. 1 Tablet</u> direkt aus dem Folienbeutel in das Fläschchen. Zerdrücken Sie die Tablette mit einem sauberen Rührstab.

- Befüllen Sie eine zweite, saubere AQUAfast 24-mm-Rundküvette mit 10 ml der Probe. Geben Sie eine <u>Glycine Tablet</u> direkt aus dem Folienbeutel in das Fläschchen. Zerdrücken Sie die Tablette mit einem sauberen Rührstab. Verschließen Sie das Fläschchen fest mit dem Deckel, und mischen Sie den Inhalt durch mehrmaliges Schwenken oder Umdrehen, bis sich die Tablette aufgelöst hat.
- 9. Den Inhalt der zweiten Küvette in das erste Fläschchen umfüllen.
- Verschließen Sie das Fläschchen fest mit dem Deckel, und mischen Sie den Inhalt durch mehrmaliges Schwenken oder Umdrehen, bis sich die Tabletten aufgelöst haben. Wischen Sie die Außenseite des Fläschchens ab.
- 11. Stellen Sie das Fläschchen in die Halterung in der Probenkammer und schließen Sie die Probenkammertür.
- 12. Tippen Sie auf die Funktionstaste Sample (Probe), um das Ergebnis in mg/l Chlordioxid anzuzeigen.

#### Hinweise:

- Alternativ kann auch die AC209950-Methode mit rechteckigen 50-mm-Fläschchen verwendet werden. Alle Blindwert- und Probenvolumina müssen mit den in dieser Anleitung angegebenen Volumina übereinstimmen, so dass die Proben möglicherweise in separaten Behältern vorbereitet und dann in das ausgewählte Fläschchen überführt werden müssen.
- Reinigung der Fläschchen: Da viele Haushaltsreiniger (z. B. Geschirrspülmittel) reduzierende Stoffe enthalten, kann es nach deren Anwendung bei der Bestimmung von Chlordioxid zu Minderbefunden kommen. Um diesen Messfehler auszuschließen, sollten die Glasgefäße chlorzehrungsfrei sein.

- Probenvorbereitung: Bei der Vorbereitung der Probe ist das Entweichen von Chlordioxidgasen (z. B. beim Pipettieren oder Schütteln) zu vermeiden. Die Analyse muss unmittelbar nach der Probenahme erfolgen.
- Die DPD-Farbentwicklung erfolgt bei einem pH-Wert von 6,2 bis 6,5. Die Reagenztablette enthält daher einen Puffer zur pH-Einstellung. Stark alkalische oder saure Wasserproben müssen vor Zugabe der Tablette auf einen pH-Wert zwischen 6 und 7 gebracht werden (dazu 0,5 mol/I Schwefelsäure bzw. 1 mol/I Natronlauge verwenden).
- Überschreitung des Messbereichs: Konzentrationen von mehr als 19 mg/l Chlordioxid können zu Ergebnissen bis hin zu 0 mg/l führen. In diesem Fall muss die Wasserprobe mit chlordioxidfreiem Wasser verdünnt werden. Dazu sollten 10 ml der verdünnten Probe mit dem Reagenz versetzt und die Messung wiederholt werden.
- Oxidationsmittel wie Chlor oder Ozon sind Störfaktoren, da sie wie Chlordioxid reagieren.

## CODL00 COD-Aufschlussröhrchen-Test für niedrigen Bereich

#### Dichromataufschluss-Methode

- 0 150 mg/l O<sub>2</sub>
- 1. Ein 16-mm-COD-Reaktionsgefäß öffnen und 2 ml deionisiertes Wasser einfüllen (dies ist das Blindwert-Reagenzgefäß).
- 2. Öffnen Sie ein zweites 16-mm-Reaktionsgefäß und geben Sie 2 ml der Probe hinzu. (Dies ist das Fläschchen.)
- Die Reaktionsgefäße mit den Kappen fest verschließen und den Inhalt durch mehrmaliges Schwenken oder Umdrehen vorsichtig mischen. VORSICHT: Die Fläschchen werden beim Mischen heiß.
- 4. Die Fläschchen <u>120 Minuten lang</u> im vorgeheizten Reaktor bei einer Temperatur von 150 °C erhitzen.
- 5. VORSICHT: Die Fläschchen werden heiß.

Die Fläschchen aus dem Reaktor nehmen und auf unter 60 °C abkühlen lassen. Drehen Sie die Fläschchen mehrmals vorsichtig um, um den noch warmen Inhalt zu mischen. Lassen Sie die Fläschchen vor der Messung abkühlen. Wischen Sie die Außenseiten der Fläschchen ab.

- 6. Laden Sie die Methode CODL00, und führen Sie sie aus.
- Befüllen Sie eine saubere AQUAfast 16-mm-Rundküvette, Bestell-Nr. AC2V16, mit deionisertem Wasser (Blindprobengefäß) Verschließen Sie das Fläschchen fest mit dem Deckel. Wischen Sie die Außenseite des Fläschchens ab.
- 8. Stellen Sie das Blindprobengefäß in die Halterung in der Probenkammer und schließen Sie die Probenkammertür.
- 9. Drücken Sie die Funktionstaste Blank (Blind), um den Blindwert zu messen.
- 10. Öffnen Sie die Probenkammertür. Entnehmen Sie das Blindprobengefäß aus der Halterung.
- 11. Stellen Sie das Blindreagenzgefäß in die Halterung in der Probenkammer und schließen Sie die Probenkammertür.
- Drücken Sie die Funktionstaste Measure Rgnt Blank (Reagenzblindwert messen), um den Reagenzblindwert zu messen.
- Öffnen Sie den Probenkammert
  ür. Entnehmen Sie das Blindreagenzgef
  ä
  ß aus der Halterung.
- 14. Stellen Sie das Fläschchen in die Halterung in der Probenkammer, und schließen Sie die Probenkammertür.
- Tippen Sie auf die Funktionstaste Sample (Probe), um das Ergebnis in mg/l Sauerstoff anzuzeigen.

- Bei der Farbumkehrmethode wird ein Reagenz verwendet, dessen Farbintensität mit zunehmender Konzentration der zu messenden Spezies in den Proben abnimmt. Die Farbumkehrmethode erfordert eine Blindprobe und einen Reagenzblindwert. Die Blindprobe ist eine klare Lösung (deionisiertes Wasser) mit einer Absorption von Null. Der Reagenzblindwert ist ein Gemisch aus dem Reagenz und der Probe und liefert den Nullkonzentrationspunkt mit dem dunkelsten Farbwert (höchste Absorption). Die Farbe der mit dem Reagenz versetzten Proben nimmt mit steigender Konzentration für diese Methode ab.
- Messen Sie Proben und Blindproben mit derselben Probengefäßcharge. Die Blindprobe ist bei Lagerung im Dunkeln stabil und kann für weitere Messungen mit Fläschchen der gleichen Charge verwendet werden.
- Keine heißen Fläschchen in den Probenhalter setzen. Kühlen Sie die Fläschchen für die endgültigen Messungen auf Raumtemperatur ab.
- Schwebstoffe im Fläschchen führen zu Fehlmessungen. Aus diesem Grund ist es wichtig, die Fläschchen vorsichtig in die Probenkammer zu stellen. Der Niederschlag am Boden der Probe sollte nicht suspendiert werden.
- Reinigen Sie die Außenseiten der Fläschchen mit einem Tuch ab. Fingerabdrücke oder andere Spuren müssen entfernt werden.
- Die Samples (Proben) können gemessen werden, wenn der Chloridgehalt 1.000 mg/l nicht überschreitet.
- In Ausnahmefällen können im Wasser enthaltene Verbindungen nicht ausreichend oxidiert werden, so dass die Ergebnisse gegenüber Referenzmethoden möglicherweise niedriger ausfallen.

# CODH00 COD-Aufschlussröhrchen-Test für mittleren Bereich

#### Dichromataufschluss-Methode

0 – 1.500 mg/l O<sub>2</sub>

- 1. Öffnen Sie ein 16-mm-COD-Reaktionsgefäß und geben Sie 2 ml deionisiertes Wasser hinzu. (Dies ist das Blindprobengefäß).
- 2. Öffnen Sie ein zweites 16-mm-COD-Reaktionsgefäß und geben Sie 2 ml der Probe hinzu (dies ist das Fläschchen).
- 3. Die Reaktionsgefäße mit den Kappen fest verschließen und den Inhalt durch mehrmaliges Schwenken oder Umdrehen vorsichtig mischen. **VORSICHT:** Die Fläschchen werden beim Mischen heiß.
- 4. Die Fläschchen <u>120 Minuten lang</u> im vorgeheizten Reaktor bei einer Temperatur von 150 °C erhitzen.
- 5. VORSICHT: Die Fläschchen werden heiß. Die Fläschchen aus dem Reaktor nehmen und auf unter 60 °C abkühlen lassen. Drehen Sie die Fläschchen mehrmals vorsichtig um, um den noch warmen Inhalt zu mischen. Lassen Sie die Fläschchen vor der Messung abkühlen. Wischen Sie die Außenseiten der Fläschchen ab.
- 6. Laden Sie die Methode CODH00, und führen Sie sie aus.
- 7. Stellen Sie das Blindprobengefäß in die Halterung in der Probenkammer und schließen Sie die Probenkammertür.
- 8. Drücken Sie die Funktionstaste Blank (Blind), um den Blindwert zu messen.
- 9. Öffnen Sie den Probenkammertür. Entnehmen Sie das Blindprobengefäß aus der Halterung.
- 10. Stellen Sie das Fläschchen in die Halterung in der Probenkammer, und schließen Sie die Probenkammertür.
- 11. Tippen Sie auf die Funktionstaste Sample (Probe), um das Ergebnis in mg/l Sauerstoff anzuzeigen.

- Messen Sie Proben und Blindproben mit derselben Probengefäßcharge. Die Blindprobe ist bei Lagerung im Dunkeln stabil und kann für weitere Messungen mit Fläschchen der gleichen Charge verwendet werden.
- Keine heißen Fläschchen in den Probenhalter setzen. Kühlen Sie die Fläschchen für die endgültigen Messungen auf Raumtemperatur ab.
- Schwebstoffe im Fläschchen führen zu Fehlmessungen. Aus diesem Grund ist es wichtig, die Fläschchen vorsichtig in die Probenkammer zu stellen. Der Niederschlag am Boden der Probe sollte nicht suspendiert werden.
- Reinigen Sie die Außenseiten der Fläschchen mit einem Tuch ab. Fingerabdrücke oder andere Spuren müssen entfernt werden.
- Die Samples (Proben) können gemessen werden, wenn der Chloridgehalt 1.000 mg/l nicht überschreitet.
- In Ausnahmefällen können im Wasser enthaltene Verbindungen nicht ausreichend oxidiert werden, so dass die Ergebnisse gegenüber Referenzmethoden möglicherweise niedriger ausfallen.
- Für Proben unter 100 mg/l wird empfohlen, den Test mit dem COD-Test für den unteren Bereich (CODL00) zu wiederholen.

# CODHP0 COD-Aufschlussröhrchen-Test für hohen Bereich

#### Dichromataufschluss-Methode

- 0 15.000 mg/I O<sub>2</sub> (hoher Bereich)
- 1. Öffnen Sie ein 16-mm-COD-Reaktionsgefäß und geben Sie 0,2 ml deionisiertes Wasser hinzu. (Dies ist das Blindprobengefäß).
- 2. Öffnen Sie ein zweites 16-mm-COD-Reaktionsgefäß und geben Sie 0,2 ml der Probe hinzu (dies ist das Fläschchen).
- 3. Die Reaktionsgefäße mit den Kappen fest verschließen und den Inhalt durch mehrmaliges Schwenken oder Umdrehen vorsichtig mischen. **VORSICHT:** Die Fläschchen werden beim Mischen heiß.
- 4. Die Fläschchen <u>120 Minuten lang</u> im vorgeheizten Reaktor bei einer Temperatur von 150 °C erhitzen.
- 5. VORSICHT: Die Fläschchen werden heiß. Die Fläschchen aus dem Reaktor nehmen und auf unter 60 °C abkühlen lassen. Drehen Sie die Fläschchen mehrmals vorsichtig um, um den noch warmen Inhalt zu mischen. Lassen Sie die Fläschchen vor der Messung abkühlen. Wischen Sie die Außenseiten der Fläschchen ab.
- 6. Laden Sie die Methode CODHP0, und führen Sie sie aus.
- 7. Stellen Sie das Blindprobengefäß in die Halterung in der Probenkammer und schließen Sie die Probenkammertür.
- 8. Drücken Sie die Funktionstaste Blank (Blind), um den Blindwert zu messen.
- 9. Öffnen Sie den Probenkammertür. Entnehmen Sie das Blindprobengefäß aus der Halterung.
- 10. Stellen Sie das Fläschchen in die Halterung in der Probenkammer, und schließen Sie die Probenkammertür.
- 11. Tippen Sie auf die Funktionstaste Sample (Probe), um das Ergebnis in mg/l Sauerstoff anzuzeigen.

- Messen Sie Proben und Blindproben mit derselben Probengefäßcharge. Die Blindprobe ist bei Lagerung im Dunkeln stabil und kann für weitere Messungen mit Fläschchen der gleichen Charge verwendet werden.
- Keine heißen Fläschchen in den Probenhalter setzen. Kühlen Sie die Fläschchen für die endgültigen Messungen auf Raumtemperatur ab.
- Schwebstoffe im Fläschchen führen zu Fehlmessungen. Aus diesem Grund ist es wichtig, die Fläschchen vorsichtig in die Probenkammer zu stellen. Der Niederschlag am Boden der Probe sollte nicht suspendiert werden.
- Reinigen Sie die Außenseiten der Fläschchen mit einem Tuch ab. Fingerabdrücke oder andere Spuren müssen entfernt werden.
- Die Samples (Proben) können gemessen werden, wenn der Chloridgehalt 1.000 mg/l nicht überschreitet.
- In Ausnahmefällen können im Wasser enthaltene Verbindungen nicht ausreichend oxidiert werden, so dass die Ergebnisse gegenüber Referenzmethoden möglicherweise niedriger ausfallen.
- Für Proben unter 1.000 mg/l wird empfohlen, den Test mit dem COD-Test für den mittleren Bereich (CODH00) zu wiederholen. Für Proben unter 100 mg/l wird empfohlen, den Test mit dem COD-Test für den unteren Bereich (CODL00) zu wiederholen.

## AC2029 Kupfertabletten-Test (freies Kupfer und Gesamtkupfer)

#### Biquinolin-Methode

0,05 – 5 mg/l Cu

- 1. Laden Sie die Methode AC202924, und führen Sie sie aus.
- Befüllen Sie eine saubere AQUAfast 24-mm-Rundküvette, Bestell-Nr. AC2V24, mit 10 ml der Probe. Verschließen Sie das Fläschchen fest mit dem Deckel. Wischen Sie die Außenseite des Fläschchens ab.
- 3. Stellen Sie das Fläschchen in die Halterung in der Probenkammer und schließen Sie die Probenkammertür.
- 4. Drücken Sie die Funktionstaste Blank (Blind), um den Blindwert zu messen.
- 5. Öffnen Sie die Probenkammertür, und nehmen Sie das Fläschchen aus der Halterung.
- 6. Geben Sie eine <u>Copper No. 1 Tablet</u> direkt aus dem Folienbeutel in das Fläschchen. Zerdrücken Sie die Tablette mit einem sauberen Rührstab.
- Verschließen Sie das Fläschchen fest mit dem Deckel, und mischen Sie den Inhalt durch mehrmaliges Schwenken oder Umdrehen, bis sich die Tablette aufgelöst hat. Wischen Sie die Außenseite des Fläschchens ab.
- 8. Stellen Sie das Fläschchen in die Halterung in der Probenkammer und schließen Sie die Probenkammertür.
- Drücken Sie die Funktionstaste Sample (Probe), um das Ergebnis in mg/l freies Kupfer anzuzeigen.
- 10. Öffnen Sie die Probenkammertür, und nehmen Sie das Fläschchen aus der Halterung.
- 11. Geben Sie eine <u>Copper No. 2 Tablet</u> direkt aus dem Folienbeutel in dasselbe Fläschchen. Zerdrücken Sie die Tablette mit einem sauberen Rührstab.
- Verschließen Sie das Fläschchen fest mit dem Deckel, und mischen Sie den Inhalt durch mehrmaliges Schwenken oder Umdrehen, bis sich die Tablette aufgelöst hat. Wischen Sie die Außenseite des Fläschchens ab.
- 13. Stellen Sie das Fläschchen in die Halterung in der Probenkammer und schließen Sie die Probenkammertür.
- Drücken Sie die Funktionstaste Sample (Probe), um das Ergebnis in mg/l Gesamtkupfer anzuzeigen.

#### Hinweise:

 Alternativ kann auch die AC202950-Methode mit rechteckigen 50-mm-Fläschchen verwendet werden. Alle Blindwert- und Probenvolumina müssen mit den in dieser Anleitung angegebenen Volumina übereinstimmen, so dass die Proben möglicherweise in separaten Behältern vorbereitet und dann in das ausgewählte Fläschchen überführt werden müssen.

# AC4P29 Kupferpulverpäckchen-Test (freies Kupfer)

#### Bicinchoninat-Methode

0,05 – 5 mg/l Cu

- 1. Laden Sie die Methode AC4P29, und führen Sie sie aus.
- Befüllen Sie eine saubere AQUAfast 24-mm-Rundküvette, Bestell-Nr. AC2V24, mit 10 ml der Probe. Verschließen Sie das Fläschchen fest mit dem Deckel. Wischen Sie die Außenseite des Fläschchens ab.
- 3. Stellen Sie das Fläschchen in die Halterung in der Probenkammer und schließen Sie die Probenkammertür.
- 4. Drücken Sie die Funktionstaste Blank (Blind), um den Blindwert zu messen.
- 5. Öffnen Sie die Probenkammertür, und nehmen Sie das Fläschchen aus der Halterung.
- 6. Geben Sie den Inhalt eines <u>Cu 1 F10 Powder Pack</u> direkt aus der Folienpackung in das Fläschchen.
- 7. Verschließen Sie das Fläschchen fest mit dem Deckel, und mischen Sie den Inhalt durch mehrmaliges Schwenken oder Umdrehen. Wischen Sie die Außenseite des Fläschchens ab.
- 8. Warten Sie, bis eine Reaktionszeit von 2 Minuten abgelaufen ist.
- 9. Stellen Sie das Fläschchen in die Halterung in der Probenkammer und schließen Sie die Probenkammertür.
- 10. Drücken Sie die Funktionstaste Sample (Probe), um das Ergebnis in mg/l freies Kupfer anzuzeigen.

- Zur Bestimmung des Gesamtkupfers ist ein Aufschluss erforderlich.
- Extrem saure Wasserproben (pH 2 oder weniger) müssen vor der Zugabe des Reagenzes auf pH 4 und pH 6 gebracht werden (mit 8 mol/l Kaliumhydroxidlösung, KOH).
- Die Genauigkeit wird durch nicht aufgelöste Pulverrückstände nicht beeinträchtigt.
- Störfaktoren:

| Cyanid (CN <sup>-</sup> ) | Cyanid verhindert die volle Farbentwicklung. 0,2 ml Formaldehyd zu 10 ml Wasserprobe geben und eine Reaktionszeit von 4 Minuten abwarten (Cyanid wird maskiert). Danach den Test wie beschrieben durchführen. Das Ergebnis mit 1,02 multiplizieren, um die Probenverdünnung durch Formaldehyd zu korrigieren. |
|---------------------------|---|
| Silber (Ag+)              | Wenn die Trübung bestehen bleibt und sich schwarz färbt, ist eine<br>Silberstörung wahrscheinlich. 10 Tropfen gesättigte Kaliumchloridlösung zu<br>75 ml Wasserprobe geben. Durch einen Feinfilter filtrieren und 10 ml der<br>gefilterten Wasserprobe für die Testdurchführung verwenden.                    |

# AC2098 Cyanursäure-Tabletten-Test

#### Melamin-Methode

0 – 160 mg/l CyA

- 1. Laden Sie die Methode AC2098, und führen Sie sie aus.
- Befüllen Sie eine saubere AQUAfast 24-mm-Rundküvette, Bestell-Nr. AC2V24, mit 10 ml der Probe. Verschließen Sie das Fläschchen fest mit dem Deckel. Wischen Sie die Außenseite des Fläschchens ab.
- 3. Stellen Sie das Fläschchen in die Halterung in der Probenkammer und schließen Sie die Probenkammertür.
- 4. Drücken Sie die Funktionstaste Blank (Blind), um den Blindwert zu messen.
- 5. Öffnen Sie die Probenkammertür, und nehmen Sie das Fläschchen aus der Halterung.
- 6. Geben Sie eine <u>Cyanuric Acid Tablet</u> direkt aus dem Folienbeutel in das Fläschchen. Zerdrücken Sie die Tablette mit einem sauberen Rührstab.
- Verschließen Sie das Fläschchen fest mit dem Deckel, und mischen Sie den Inhalt durch mehrmaliges Schwenken oder Umdrehen, bis sich die Tablette aufgelöst hat (siehe Hinweise unten). Wischen Sie die Außenseite des Fläschchens ab.
- 8. Stellen Sie das Fläschchen in die Halterung in der Probenkammer und schließen Sie die Probenkammertür.
- 9. Drücken Sie die Funktionstaste Sample (Probe), um das Ergebnis in mg/l Cyanursäure anzuzeigen.

- Wenn Cyanursäure vorhanden ist, wird die Lösung eingetrübt. Kleine einzelne Partikel werden nicht unbedingt durch Cyanursäure verursacht.
- Lösen Sie die Tablette vollständig auf (schwenken Sie das Fläschchen etwa 1 Minute lang). Nicht gelöste Tablettenpartikel können zu falsch erhöhten Messergebnissen führen.
- Überschreitung des zulässigen Messbereichs: Proben mit einer Konzentration über 90 mg/l müssen mit cyanursäurefreiem Wasser verdünnt werden. 10 ml der verdünnten Probe sollten wie oben beschrieben getestet und die angezeigten Ergebnisse unter Verwendung des Verdünnungsfaktors berechnet werden.

### AC2009 SPADNS Fluorid-Flüssigreagenz-Test

#### SPADNS-Methode

0,05 – 2 mg/l F

- 1. Laden Sie die Methode AC2009, und führen Sie sie aus.
- Befüllen Sie eine saubere AQUAfast 24-mm-Rundküvette, Bestell-Nr. AC2V24, mit 10 ml deionisiertem Wasser. Verschließen Sie das Fläschchen fest mit dem Deckel. Wischen Sie die Außenseite des Fläschchens ab.
- Stellen Sie das Fläschchen in die Halterung in der Probenkammer und schließen Sie die Probenkammert
  ür.
- 4. Drücken Sie die Funktionstaste Blank (Blind), um den Blindwert zu messen.
- 5. Öffnen Sie die Probenkammertür, und nehmen Sie das Fläschchen aus der Halterung.
- Geben Sie genau 2 ml <u>SPADNS Solution</u> in das Fläschchen. VORSICHT: Das Fläschchen wird vollständig gefüllt.
- Verschließen Sie das Fläschchen fest mit dem Deckel, und mischen Sie den Inhalt durch mehrmaliges Schwenken oder Umdrehen. Wischen Sie die Außenseite des Fläschchens ab.
- Stellen Sie das Fläschchen in die Halterung in der Probenkammer und schließen Sie die Probenkammert
  ür.
- Drücken Sie die Funktionstaste Measure Rgnt Blank (Reagenzblindwert messen), um den Reagenzblindwert zu messen.
- 10. Offnen Sie die Probenkammertür, und nehmen Sie das Fläschchen aus der Halterung.
- 11. Das Fläschchen entleeren, Fläschchen und Deckel mehrmals gründlich ausspülen und dann genau 10 ml Probe in das Fläschchen füllen.
- 12. Geben Sie genau 2 ml <u>SPADNS Solution</u> in das Fläschchen. **VORSICHT:** Das Fläschchen wird vollständig gefüllt.
- Verschließen Sie das Fläschchen fest mit dem Deckel, und mischen Sie den Inhalt durch mehrmaliges Schwenken oder Umdrehen. Wischen Sie die Außenseite des Fläschchens ab.
- 14. Stellen Sie das Fläschchen in die Halterung in der Probenkammer und schließen Sie die Probenkammertür.
- 15. Tippen Sie auf die Funktionstaste Sample (Probe), um das Ergebnis in mg/l Fluorid anzuzeigen.

#### Hinweise:

 Bei der Farbumkehrmethode wird ein Reagenz verwendet, dessen Farbintensität mit zunehmender Konzentration der zu messenden Spezies in den Proben abnimmt.
 Die Farbumkehrmethode erfordert eine Blindprobe und einen Reagenzblindwert.
 Die Blindprobe ist eine klare Lösung (deionisiertes Wasser) mit einer Absorption von Null.
 Der Reagenzblindwert ist ein Gemisch aus dem Reagenz und der Probe und liefert den Nullkonzentrationspunkt mit dem dunkelsten Farbwert (höchste Absorption). Die Farbe der mit dem Reagenz versetzten Proben nimmt mit steigender Konzentration für diese Methode ab.

- Für die Tests (Reagenzblindwert- und Probenmessung) und die Ein-Punkt-Kalibrierung muss dieselbe Charge der SPADNS-Reagenzlösung verwendet werden. Die Ein-Punkt-Kalibrierung muss für jede neue Charge der SPADNS-Reagenzlösung durchgeführt werden (siehe Standard Methods 20. Aufl., 1998, APHA, AWWA, WEF 4500 F<sup>-</sup> D, 4.a).
- Bei den Tests (Messung von Blindprobe, Reagenzblindwert und Probe) und bei der Ein-Punkt-Kalibrierung sollte dasselbe Fläschchen verwendet werden, da unterschiedliche Fläschchen geringfügige Toleranzen aufweisen können.
- Die Kalibrierlösung und die Wasserproben sollten die gleiche Temperatur (+/- 1 °C) haben.
- Da das Testergebnis in hohem Maße von den exakten Proben- und Reagenzvolumina abhängt, sollten die Proben- und Reagenzvolumina immer mit einer 10 ml oder 2 ml Vollpipette (Klasse A) abgemessen werden.
- Die Genauigkeit der Testmethoden nimmt bei einer Flouridkonzentration von über 1,2 mg/l ab. Obwohl die Ergebnisse für die meisten Anwendungen ausreichend genau sind, können exaktere Ergebnisse erzielt werden, wenn die Probe vor der Verwendung im Verhältnis 1:1 verdünnt und das Ergebnis anschließend mit 2 multipliziert wird.
- Die SPADNS-Reagenzlösung enthält Arsenit. Chlorkonzentrationen bis zu 5 mg/l beeinträchtigen die Testergebnisse nicht.
- Meerwasser- und Abwasserproben müssen destilliert werden.

# AC3032T Wasserhärte-Tabletten-Test (Gesamthärte)

#### Metallphthalein-Methode

2 – 50 mg/l CaCO<sub>3</sub> (niedriger Bereich) oder 20 – 500 mg/l CaCO<sub>3</sub> (hoher Bereich)

#### Für Messungen der Gesamthärte im niedrigen Bereich:

- 1. Laden Sie die Methode AC3032TL, und führen Sie sie aus.
- Befüllen Sie eine saubere AQUAfast 24-mm-Rundküvette, Bestell-Nr. AC2V24, mit 10 ml der Probe. Verschließen Sie das Fläschchen fest mit dem Deckel. Wischen Sie die Außenseite des Fläschchens ab.
- 3. Stellen Sie das Fläschchen in die Halterung in der Probenkammer und schließen Sie die Probenkammertür.
- 4. Drücken Sie die Funktionstaste Blank (Blind), um den Blindwert zu messen.
- 5. Öffnen Sie die Probenkammertür, und nehmen Sie das Fläschchen aus der Halterung.
- 6. Geben Sie eine <u>Hardcheck P Tablet</u> direkt aus dem Folienbeutel in das Fläschchen. Zerdrücken Sie die Tablette mit einem sauberen Rührstab.
- Verschließen Sie das Fläschchen fest mit dem Deckel, und mischen Sie den Inhalt durch mehrmaliges Schwenken oder Umdrehen, bis sich die Tablette aufgelöst hat. Wischen Sie die Außenseite des Fläschchens ab.
- 8. Warten Sie, bis eine Reaktionszeit von 5 Minuten abgelaufen ist.
- 9. Stellen Sie das Fläschchen in die Halterung in der Probenkammer und schließen Sie die Probenkammertür.
- 10. Drücken Sie die Funktionstaste Sample (Probe), um das Ergebnis in mg/l Gesamthärte anzuzeigen.

#### Für Messungen der Gesamthärte im oberen Bereich:

- 1. Laden Sie die Methode AC3032TH, und führen Sie sie aus.
- Befüllen Sie eine saubere AQUAfast 24-mm-Rundküvette, Bestell-Nr. AC2V24, mit 1 ml Probe und 9 ml deionisiertem Wasser. Verschließen Sie das Fläschchen fest mit dem Deckel. Wischen Sie die Außenseite des Fläschchens ab.
- 3. Stellen Sie das Fläschchen in die Halterung in der Probenkammer und schließen Sie die Probenkammertür.
- 4. Drücken Sie die Funktionstaste Blank (Blind), um den Blindwert zu messen.
- 5. Öffnen Sie die Probenkammertür, und nehmen Sie das Fläschchen aus der Halterung.
- 6. Geben Sie eine <u>Hardcheck P Tablet</u> direkt aus dem Folienbeutel in das Fläschchen. Zerdrücken Sie die Tablette mit einem sauberen Rührstab.
- Verschließen Sie das Fläschchen fest mit dem Deckel, und mischen Sie den Inhalt durch mehrmaliges Schwenken oder Umdrehen, bis sich die Tablette aufgelöst hat. Wischen Sie die Außenseite des Fläschchens ab.
- 8. Warten Sie, bis eine Reaktionszeit von 5 Minuten abgelaufen ist.
- 9. Stellen Sie das Fläschchen in die Halterung in der Probenkammer und schließen Sie die Probenkammertür.
- 10. Drücken Sie die Funktionstaste Sample (Probe), um das Ergebnis in mg/l Gesamthärte anzuzeigen.

#### Hinweise:

• Stark alkalische oder saure Wasserproben müssen vor Zugabe der Tablette auf einen pH-Wert von 4 bis 10 gebracht werden (dazu 1 mol/l Salzsäure bzw. 1 mol/l Natronlauge verwenden).

# AC2030 Hydrazinpulver-Test

#### Dimethylaminobenzaldehyd-Methode

0,05 - 0,5 mg/l N<sub>2</sub>H<sub>4</sub>

- 1. Laden Sie die Methode AC2030, und führen Sie sie aus.
- Befüllen Sie eine saubere AQUAfast 24-mm-Rundküvette, Bestell-Nr. AC2V24, mit 10 ml der Probe. Verschließen Sie das Fläschchen fest mit dem Deckel. Wischen Sie die Außenseite des Fläschchens ab.
- 3. Stellen Sie das Fläschchen in die Halterung in der Probenkammer und schließen Sie die Probenkammertür.
- 4. Drücken Sie die Funktionstaste Blank (Blind), um den Blindwert zu messen.
- 5. Öffnen Sie die Probenkammertür, und nehmen Sie das Fläschchen aus der Halterung.
- 6. Geben Sie ein Gramm (1 g) Hydrazine Powder in das Fläschchen.
- 7. Verschließen Sie das Fläschchen fest mit dem Deckel, und mischen Sie den Inhalt durch mehrmaliges Schwenken oder Umdrehen. Wischen Sie die Außenseite des Fläschchens ab.
- 8. Warten Sie, bis eine Reaktionszeit von 10 Minuten abgelaufen ist.
- 9. Die leichte Trübung, die bei der Reagenzzugabe entsteht, muss durch Filtration entfernt werden
- 10. Stellen Sie das Fläschchen in die Halterung in der Probenkammer und schließen Sie die Probenkammertür.
- 11. Tippen Sie auf die Funktionstaste Sample (Probe), um das Ergebnis in mg/l Hydrazin anzuzeigen.

- Trübe Wasserproben müssen vor der Blindwertmessung gefiltert werden.
- Die Temperatur der Wasserprobe sollte max. 21 °C betragen.
- Verwenden Sie den Hydrazinlöffel: 1 g entspricht einem gestrichenen Löffel.
- Empfohlen werden qualitative Faltenfilterpapiere für mittlere Ausfällungen.
- Um zu pr
  üfen, ob das Reagenz gealtert ist (Lagerung 
  über einen l
  ängeren Zeitraum) f
  ühren Sie den Test wie oben beschrieben mit Leitungswasser durch. Liegt das Ergebnis 
  über der Nachweisgrenze von 0,05 mg/l, sollten Sie das Reagenz nur unter Vorbehalt verwenden, da es zu gr
  ößeren Ergebnisabweichungen kommen kann.

## AC2078 Eisentabletten-Test (Eisen II und III)

#### **PPST-Methode**

0,02 - 1 mg/l Fe

- 1. Laden Sie die Methode AC207824, und führen Sie sie aus.
- Befüllen Sie eine saubere AQUAfast 24-mm-Rundküvette, Bestell-Nr. AC2V24, mit 10 ml der Probe. Verschließen Sie das Fläschchen fest mit dem Deckel. Wischen Sie die Außenseite des Fläschchens ab.
- 3. Stellen Sie das Fläschchen in die Halterung in der Probenkammer und schließen Sie die Probenkammertür.
- 4. Drücken Sie die Funktionstaste Blank (Blind), um den Blindwert zu messen.
- 5. Öffnen Sie die Probenkammertür, und nehmen Sie das Fläschchen aus der Halterung.
- 6. Geben Sie eine <u>Iron LR Tablet</u> direkt aus dem Folienbeutel in das Fläschchen. Zerdrücken Sie die Tablette mit einem sauberen Rührstab.
- Verschließen Sie das Fläschchen fest mit dem Deckel, und mischen Sie den Inhalt durch mehrmaliges Schwenken oder Umdrehen, bis sich die Tablette aufgelöst hat. Wischen Sie die Außenseite des Fläschchens ab.
- 8. Warten Sie, bis eine Reaktionszeit von <u>5 Minuten</u> abgelaufen ist.
- 9. Stellen Sie das Fläschchen in die Halterung in der Probenkammer und schließen Sie die Probenkammertür.
- 10. Tippen Sie auf die Funktionstaste Sample (Probe), um das Ergebnis in mg/l Eisen anzuzeigen.

- Alternativ kann auch die AC207850-Methode mit rechteckigen 50-mm-Fläschchen verwendet werden. Alle Blindwert- und Probenvolumina müssen mit den in dieser Anleitung angegebenen Volumina übereinstimmen, so dass die Proben möglicherweise in separaten Behältern vorbereitet und dann in das ausgewählte Fläschchen überführt werden müssen.
- Mit dieser Methode wird das gelöste Gesamteisen als Fe<sup>2+</sup> und Fe<sup>3+</sup> bestimmt.
- Für die Bestimmung des gelösten und nicht gelösten Gesamteisens (lösliches und unlösliches Eisen) ist ein Aufschluss erforderlich. Ein Beispiel wird hier beschrieben:
  - i. 1 ml konzentrierte Schwefelsäure zu 100 ml Wasserprobe geben. Erhitzen und 10 Minuten lang kochen, bis alle Partikel aufgelöst sind. Nach dem Abkühlen die Probe mit Hilfe von Ammoniaklösung auf einen pH-Wert von 3 bis 6 einstellen. Mit deionisiertem Wasser auf das ursprüngliche Volumen von 100 ml auffüllen und gut mischen. 10 ml dieser vorbehandelten Lösung werden für die Analyse verwendet (Durchführung nach der Beschreibung der gewählten Testmethode).
  - ii. Mit organischen Verbindungen wie Korrosionsschutzmitteln behandeltes Wasser muss bei Bedarf oxidiert werden, um die Eisenkomplexe aufzuschließen. Dazu 1 ml konzentrierte Schwefelsäure und 1 ml konzentrierte Salpetersäure in 100 ml Wasserprobe geben und auf etwa die Hälfte des Volumens einkochen. Nach dem Abkühlen nach der obigen Beschreibung fortfahren.

## AC4P78 Iron (Ferro) Powder Pack Test

#### 1,10-Phenanthrolin-Methode

0,02 – 3 mg/l Fe

- 1. Laden Sie die Methode AC4P78, und führen Sie sie aus.
- Befüllen Sie eine saubere AQUAfast 24-mm-Rundküvette, Bestell-Nr. AC2V24, mit 10 ml der Probe. Verschließen Sie das Fläschchen fest mit dem Deckel. Wischen Sie die Außenseite des Fläschchens ab.
- 3. Stellen Sie das Fläschchen in die Halterung in der Probenkammer und schließen Sie die Probenkammertür.
- 4. Drücken Sie die Funktionstaste Blank (Blind), um den Blindwert zu messen.
- 5. Öffnen Sie die Probenkammertür, und nehmen Sie das Fläschchen aus der Halterung.
- 6. Geben Sie den Inhalt eines <u>Ferro F10 Powder Pack</u> direkt aus der Folienpackung in das Fläschchen.
- 7. Verschließen Sie das Fläschchen fest mit dem Deckel, und mischen Sie den Inhalt durch mehrmaliges Schwenken oder Umdrehen. Wischen Sie die Außenseite des Fläschchens ab.
- 8. Warten Sie, bis eine Reaktionszeit von <u>3 Minuten</u> abgelaufen ist.
- 9. Stellen Sie das Fläschchen in die Halterung in der Probenkammer und schließen Sie die Probenkammertür.
- 10. Tippen Sie auf die Funktionstaste Sample (Probe), um das Ergebnis in mg/l Eisen anzuzeigen.

- Das Reagenz reagiert mit allen löslichen und den meisten unlöslichen Formen von Eisen in der Wasserprobe.
- Eisenoxid erfordert einen vorherigen Aufschluss: Verwenden Sie einen milden, kräftigen oder Digesdahl-Aufschluss. Ein Beispiel für den Säureaufschluss wird hier beschrieben:
  - i. 1 ml konzentrierte Schwefelsäure zu 100 ml Wasserprobe geben. Erhitzen und 10 Minuten lang kochen, bis alle Partikel aufgelöst sind. Nach dem Abkühlen die Probe mit Hilfe von Ammoniaklösung auf einen pH-Wert von 3 bis 6 einstellen. Mit deionisiertem Wasser auf das ursprüngliche Volumen von 100 ml auffüllen und gut mischen. 10 ml dieser vorbehandelten Lösung werden für die Analyse verwendet (Durchführung nach der Beschreibung der gewählten Testmethode).
  - ii. Mit organischen Verbindungen wie Korrosionsschutzmitteln behandeltes Wasser muss bei Bedarf oxidiert werden, um die Eisenkomplexe aufzuschließen. Dazu 1 ml konzentrierte Schwefelsäure und 1 ml konzentrierte Salpetersäure in 100 ml Wasserprobe geben und auf etwa die Hälfte des Volumens einkochen. Nach dem Abkühlen nach der obigen Beschreibung fortfahren.
- Sehr stark alkalische oder saure Proben müssen vor der Analyse auf einen pH-Wert zwischen 3 und 5 eingestellt werden.
- Die Genauigkeit wird durch nicht aufgelöste Pulverrückstände nicht beeinträchtigt.
- Wasserproben mit sichtbaren Rostanteilen mindestens 5 Minuten lang reagieren lassen.

# AC4P79 Eisenpulverpäckchen-Test (Gesamteisen)

#### **TPTZ-Methode**

0,02 - 1,8 mg/l Fe

- 1. Laden Sie die Methode AC4P79, und führen Sie sie aus.
- 2. Verwenden Sie zwei saubere AQUAfast 24-mm-Rundfläschchen, Bestell-Nr. AC2V24.
- 3. Geben Sie 10 ml deionisiertes Wasser in das erste Fläschchen. (Dies ist das Blindprobengefäß.)
- 4. Geben Sie 10 ml der Probe in das zweite Fläschchen. (Dies ist das Fläschchen.)
- Geben Sie den Inhalt eines Iron TPTZ F10 Powder Pack direkt aus der Folienpackung in jedes Fläschchen. Verschließen Sie die Fläschchen fest mit den Kappen, und mischen Sie den Inhalt durch mehrmaliges Schwenken oder Umdrehen. Wischen Sie die Außenseiten der Fläschchen ab.
- 6. Warten Sie, bis eine Reaktionszeit von <u>3 Minuten</u> abgelaufen ist.
- 7. Stellen Sie das Blindprobengefäß in die Halterung in der Probenkammer und schließen Sie die Probenkammertür.
- 8. Drücken Sie die Funktionstaste Blank (Blind), um den Blindwert zu messen.
- 9. Öffnen Sie den Probenkammertür. Entnehmen Sie das Blindprobengefäß aus der Halterung.
- 10. Stellen Sie das Fläschchen in die Halterung in der Probenkammer, und schließen Sie die Probenkammertür.
- 11. Tippen Sie auf die Funktionstaste Sample (Probe), um das Ergebnis in mg/l Eisen anzuzeigen.

- Zur Bestimmung des Gesamteisens ist ein Aufschluss erforderlich. Das TPTZ-Reagenz kann die meisten unlöslichen Eisenoxide ohne Aufschluss zurückgewinnen.
- Spülen Sie alle Glasgefäße zunächst mit einer Salzsäurelösung im Verhältnis 1:1 und anschließend mit deionisiertem Wasser, um Eisenablagerungen zu entfernen, die zu leicht überhöhten Messergebnissen führen können.
- Stark alkalische oder saure Wasserproben müssen vor der Analyse auf einen pH-Wert zwischen 3 und 8 gebracht werden (dazu 0,5 mol/l Schwefelsäure bzw. 1 mol/l Natronlauge verwenden).
- Störungen: Wenn Störungen auftreten, wird die Farbentwicklung gehemmt oder es bilden sich Ausfällungen. Die folgenden Werte beziehen sich auf einen Standard mit einer Eisenkonzentration von 0,5 mg/l. Die folgenden Stoffe sind in den jeweils angegebenen Konzentrationen unbedenklich:

| Substanz                | Kein Störeinfluss auf |
|-------------------------|-----------------------|
| Kadmium                 | 4,0 mg/l              |
| Chrom ( <sup>3+</sup> ) | 0,25 mg/l             |
| Chrom (6+)              | 1,2 mg/l              |
| Kobalt                  | 0,05 mg/l             |
| Kupfer                  | 0,6 mg/l              |
| Cyanid                  | 2,8 mg/l              |

| Substanz    | Kein Störeinfluss auf |
|-------------|-----------------------|
| Mangan      | 50 mg/l               |
| Quecksilber | 0,4 mg/l              |
| Molybdän    | 4,0 mg/l              |
| Nickel      | 1,0 mg/l              |
| Nitrit-Ion  | 0,8 mg/l              |

## AC2055 Mangantabletten-Test

#### Formaldoxim-Methode

0,2 – 4 mg/l Mn

- 1. Laden Sie die Methode AC2055, und führen Sie sie aus.
- Befüllen Sie eine saubere AQUAfast 24-mm-Rundküvette, Bestell-Nr. AC2V24, mit 10 ml der Probe. Verschließen Sie das Fläschchen fest mit dem Deckel. Wischen Sie die Außenseite des Fläschchens ab.
- 3. Stellen Sie das Fläschchen in die Halterung in der Probenkammer und schließen Sie die Probenkammertür.
- 4. Drücken Sie die Funktionstaste Blank (Blind), um den Blindwert zu messen.
- 5. Öffnen Sie die Probenkammertür, und nehmen Sie das Fläschchen aus der Halterung.
- 6. Geben Sie eine <u>Manganese LR 1 Tablet</u> direkt aus dem Folienbeutel in das Fläschchen. Zerdrücken Sie die Tablette mit einem sauberen Rührstab.
- 7. Geben Sie eine <u>Manganese LR 2 Tablet</u> direkt aus dem Folienbeutel in das Fläschchen. Zerdrücken Sie die Tablette mit einem sauberen Rührstab.
- Verschließen Sie das Fläschchen fest mit dem Deckel, und mischen Sie den Inhalt durch mehrmaliges Schwenken oder Umdrehen, bis sich die Tabletten aufgelöst haben. Wischen Sie die Außenseite des Fläschchens ab.
- 9. Warten Sie, bis eine Reaktionszeit von 5 Minuten abgelaufen ist.
- 10. Stellen Sie das Fläschchen in die Halterung in der Probenkammer und schließen Sie die Probenkammertür.
- 11. Tippen Sie auf die Funktionstaste Sample (Probe), um das Ergebnis in mg/l Mangan anzuzeigen.

# AC4P54 Mangan-Pulverpäckchen- und -Flüssigreagenztest für niedrigen Bereich

#### PAN-Methode

0,01 – 0,7 mg/l Mn

- 1. Laden Sie die Methode AC4P54, und führen Sie sie aus.
- 2. Verwenden Sie zwei saubere AQUAfast 24-mm-Rundfläschchen, Bestell-Nr. AC2V24.
- Geben Sie 10 ml deionisiertes Wasser in das erste Fläschchen. (Dies ist das Blindprobengefäß.)
- 4. Geben Sie 10 ml der Probe in das zweite Fläschchen. (Dies ist das Fläschchen.)
- Geben Sie den Inhalt eines <u>Ascorbic Acid Powder Pack</u> direkt aus der Folienpackung in die Fläschchen. Verschließen Sie die Fläschchen fest mit den Kappen, und mischen Sie den Inhalt durch mehrmaliges Schwenken oder Umdrehen.
- Geben Sie in jedes Fläschchen 15 Tropfen <u>Alkaline Cyanide Reagent Solution</u>. Fügen Sie Tropfen in gleicher Größe hinzu, indem Sie die Flasche senkrecht halten und langsam zusammendrücken. Verschließen Sie die Fläschchen fest mit den Kappen, und mischen Sie den Inhalt durch mehrmaliges Schwenken oder Umdrehen.
- 7. Geben Sie in jedes Fläschchen 21 Tropfen <u>PAN Indicator Solution</u>. Fügen Sie Tropfen in gleicher Größe hinzu, indem Sie die Flasche senkrecht halten und langsam zusammendrücken. Verschließen Sie die Fläschchen fest mit den Kappen, und mischen Sie den Inhalt durch mehrmaliges Schwenken oder Umdrehen. Wischen Sie die Außenseiten der Fläschchen ab.
- 8. Warten Sie, bis eine Reaktionszeit von 2 Minuten abgelaufen ist.
- 9. Stellen Sie das Blindprobengefäß in die Halterung in der Probenkammer, und schließen Sie die Probenkammertür.
- 10. Drücken Sie die Funktionstaste Blank (Blind), um den Blindwert zu messen.
- 11. Öffnen Sie den Probenkammertür. Entnehmen Sie das Blindprobengefäß aus der Halterung.
- 12. Stellen Sie das Fläschchen in die Halterung in der Probenkammer, und schließen Sie die Probenkammertür.
- 13. Tippen Sie auf die Funktionstaste Sample (Probe), um das Ergebnis in mg/l Mangan anzuzeigen.

- Spülen Sie alle Glasgefäße zunächst mit einer 1:1-Salpetersäurelösung und dann mit deionisiertem Wasser ab.
- Wasserproben mit einer Härte von mehr als 300 mg/l CaCO<sub>3</sub>: nach Zugabe des Ascorbinsäure-Pulverpäckchens 10 Tropfen Rochelle-Salzlösung hinzufügen.
- Nach Zugabe der alkalischen Cyanid-Reagenzlösung kann sich in einigen Wasserproben eine eingetrübte oder trübe Lösung bilden. Die Trübung sollte nach Zugabe der PAN-Indikatorlösung verschwinden.
- Wasserproben mit einer Eisenkonzentration von mehr als 5 mg/l mindestens 10 Minuten lang reagieren lassen.

# AC4P55 Mangan-Pulverpäckchen-Test für hohen Bereich

#### Periodat-Oxidationsmethode

0,1 – 18 mg/l Mn (hoher Bereich)

- 1. Laden Sie die Methode AC4P55, und führen Sie sie aus.
- Befüllen Sie eine saubere AQUAfast 24-mm-Rundküvette, Bestell-Nr. AC2V24, mit 10 ml der Probe. Verschließen Sie das Fläschchen fest mit dem Deckel. Wischen Sie die Außenseite des Fläschchens ab.
- 3. Stellen Sie das Fläschchen in die Halterung in der Probenkammer und schließen Sie die Probenkammertür.
- 4. Drücken Sie die Funktionstaste Blank (Blind), um den Blindwert zu messen.
- 5. Öffnen Sie die Probenkammertür, und nehmen Sie das Fläschchen aus der Halterung.
- 6. Geben Sie den Inhalt eines <u>Manganese Citrate Buffer Powder Pack</u> direkt aus der Folienpackung in das Fläschchen.
- 7. Verschließen Sie das Fläschchen fest mit dem Deckel, und mischen Sie den Inhalt durch mehrmaliges Schwenken oder Umdrehen.
- 8. Geben Sie den Inhalt eines <u>Sodium Periodate Powder Pack</u> direkt aus der Folienpackung in das Fläschchen.
- 9. Verschließen Sie das Fläschchen fest mit dem Deckel, und mischen Sie den Inhalt durch mehrmaliges Schwenken oder Umdrehen. Wischen Sie die Außenseite des Fläschchens ab.
- 10. Warten Sie, bis eine Reaktionszeit von 2 Minuten abgelaufen ist.
- 11. Stellen Sie das Fläschchen in die Halterung in der Probenkammer und schließen Sie die Probenkammertür.
- 12. Tippen Sie auf die Funktionstaste Sample (Probe), um das Ergebnis in mg/l Mangan anzuzeigen.

- Dieser Test ist für die Bestimmung von löslichem Mangan in Wasser und Abwasser geeignet.
- Stark gepufferte Wasserproben oder extreme pH-Werte können die Pufferkapazität der Reagenzien übersteigen und erfordern eine Vorbehandlung der Probe. Wurden die Proben für die Lagerung angesäuert, so ist der pH-Wert vor dem Test mit 5 mol/l (5 N) Natronlauge auf einen Wert zwischen 4 und 5 einzustellen. Ein pH-Wert von 5 darf nicht überschritten werden, da es sonst zu Manganausfällungen kommen kann.
- Störungen:

| Störsubstanz | Störungsgrad          |
|--------------|-----------------------|
| Kalzium      | Mehr als 700 mg/l     |
| Chlorid      | Mehr als 70.000 mg/l  |
| Eisen        | Mehr als 5 mg/l       |
| Magnesium    | Mehr als 100.000 mg/l |

# AC4P42 Molybdat-Pulverpäckchen

#### Mercaptoessigsäure-Methode

0,5 – 66 mg/l MoO<sub>4</sub> (Äquivalent zu 0,3 – 40 mg/l Mo)

- 1. Laden Sie die Methode AC4P42, und führen Sie sie aus.
- Befüllen Sie eine saubere AQUAfast 24-mm-Rundküvette, Bestell-Nr. AC2V24, mit 10 ml der Probe. Verschließen Sie das Fläschchen fest mit dem Deckel. Wischen Sie die Außenseite des Fläschchens ab.
- 3. Stellen Sie das Fläschchen in die Halterung in der Probenkammer und schließen Sie die Probenkammertür.
- 4. Drücken Sie die Funktionstaste Blank (Blind), um den Blindwert zu messen.
- 5. Öffnen Sie die Probenkammertür, und nehmen Sie das Fläschchen aus der Halterung.
- Geben Sie den Inhalt eines <u>Molybdenum HR 1 F10 Powder Pack</u> direkt aus der Folienpackung in das Fläschchen. Verschließen Sie das Fläschchen fest mit dem Deckel, und mischen Sie den Inhalt durch mehrmaliges Schwenken oder Umdrehen.
- Geben Sie den Inhalt eines <u>Molybdenum HR 2 F10 Powder Pack</u> direkt aus der Folienpackung in das Fläschchen. Verschließen Sie das Fläschchen fest mit dem Deckel, und mischen Sie den Inhalt durch mehrmaliges Schwenken oder Umdrehen.
- Geben Sie den Inhalt eines Molybdenum HR 3 F10 Powder Pack direkt aus der Folienpackung in das Fläschchen. Verschließen Sie das Fläschchen fest mit dem Deckel, und mischen Sie den Inhalt durch mehrmaliges Schwenken oder Umdrehen. Wischen Sie die Außenseite des Fläschchens ab.
- 9. Warten Sie, bis eine Reaktionszeit von <u>5 Minuten</u> abgelaufen ist.
- 10. Stellen Sie das Fläschchen in die Halterung in der Probenkammer und schließen Sie die Probenkammertür.
- 11. Tippen Sie auf die Funktionstaste Sample (Probe), um das Ergebnis in mg/l Molybdat anzuzeigen.

- Trübe Wasserproben vor der Analyse mit Filterpapier und Trichter filtern.
- Stark gepufferte Wasserproben oder extreme pH-Werte sollten mit 1 mol/l Salpetersäure oder 1 mol/l Natriumhydroxid (Natronlauge) auf einen pH-Wert von annähernd 7 eingestellt werden.
- Konzentrationen über 10 mg/l Cu verursachen erhöhte Testwerte, wenn die Reaktionsdauer von 5 Minuten überschritten wird. Der Test sollte daher unbedingt wie beschrieben durchgeführt werden.
- Stoffe, die bei folgenden Konzentrationen Störfaktoren darstellen:

| Aluminium | 50 mg/l                  |
|-----------|--------------------------|
| Chrom     | 1.000 mg/l               |
| Eisen     | 50 mg/l                  |
| Nickel    | 50 mg/l                  |
| Nitrit    | In allen Konzentrationen |
# **ACR007 Nitrat-Reaktionsröhrchen-Test**

#### Chromotropsäure-Methode

- 1 30 mg/l N (Nitrat als Stickstoffträger)
- 1. Laden Sie die Methode ACR011, und führen Sie sie aus.
- 2. Öffnen Sie ein 16-mm-Reaktionsgefäß (Reagenz A) und füllen Sie 1 ml deionisiertes Wasser ein. (Dies ist das Blindprobengefäß).
- 3. Öffnen Sie ein zweites 16-mm-Reaktionsgefäß (Reagenz A) und geben Sie 1 ml der Probe hinzu. (Dies ist das Fläschchen.)
- 4. Den Inhalt eines <u>Nitrate Chromotropic Powder Pack</u> direkt aus der Folienpackung in die Fläschchen geben.
- Verschließen Sie die Fläschchen fest mit den Kappen, und drehen Sie sie vorsichtig etwa 10 Mal um, um den Inhalt zu mischen. Einige Feststoffe lösen sich möglicherweise nicht auf. Wischen Sie die Außenseite des Fläschchens ab.
- 6. Warten Sie, bis eine Reaktionszeit von <u>5 Minuten</u> abgelaufen ist.
- 7. Stellen Sie das Blindprobengefäß in die Halterung in der Probenkammer, und schließen Sie die Probenkammertür.
- 8. Drücken Sie die Funktionstaste Blank (Blind), um den Blindwert zu messen.
- 9. Öffnen Sie die Probenkammertür. Entnehmen Sie das Blindprobengefäß aus der Halterung.
- 10. Stellen Sie das Fläschchen in die Halterung in der Probenkammer, und schließen Sie die Probenkammertür.
- 11. Tippen Sie auf die Funktionstaste Sample (Probe), um das Ergebnis in mg/l Nitrat als Stickstoffträger anzuzeigen.

- Einige Feststoffe lösen sich möglicherweise nicht auf.
- Umrechnung:  $mg/I NO_3 = mg/I N \times 4,43$

# AC2046 Nitrittabletten-Test

#### N-(1-Naphthyl)-Ethylendiamin-Methode

0,01 – 0,5 mg/l N (Nitrit als Stickstoffträger)

- 1. Laden Sie die Methode AC2046, und führen Sie sie aus.
- Befüllen Sie eine saubere AQUAfast 24-mm-Rundküvette, Bestell-Nr. AC2V24, mit 10 ml der Probe. Verschließen Sie das Fläschchen fest mit dem Deckel. Wischen Sie die Außenseite des Fläschchens ab.
- 3. Stellen Sie das Fläschchen in die Halterung in der Probenkammer und schließen Sie die Probenkammertür.
- 4. Drücken Sie die Funktionstaste Blank (Blind), um den Blindwert zu messen.
- 5. Öffnen Sie die Probenkammertür, und nehmen Sie das Fläschchen aus der Halterung.
- 6. Geben Sie eine <u>Nitrite LR Tablet</u> direkt aus dem Folienbeutel in das Fläschchen. Zerdrücken Sie die Tablette mit einem sauberen Rührstab.
- Verschließen Sie das Fläschchen fest mit dem Deckel, und mischen Sie den Inhalt durch mehrmaliges Schwenken oder Umdrehen, bis sich die Tablette aufgelöst hat. Wischen Sie die Außenseite des Fläschchens ab.
- 8. Warten Sie, bis eine Reaktionszeit von 10 Minuten abgelaufen ist.
- 9. Stellen Sie das Fläschchen in die Halterung in der Probenkammer und schließen Sie die Probenkammertür.
- 10. Tippen Sie auf die Funktionstaste Sample (Probe), um das Ergebnis in mg/l Nitrit als Stickstoffträger anzuzeigen.

- Folgende Ionen können stören, da sie unter den Reaktionsbedingungen zu Ausfällungen führen: Antimon (III), Eisen (III), Blei, Quecksilber (I), Silber, Chloroplatinat, Metavanadat und Bismut. Kupfer(II)-Ionen können zu niedrigeren Messergebnissen führen, da sie die Zersetzung von Diazoniumsalz beschleunigen. Es ist in der Praxis unwahrscheinlich, dass diese störenden Ionen in so hohen Konzentrationen auftreten, dass sie signifikante Messfehler verursachen.
- Umrechnung:  $mg/I NO_2 = mg/I N \times 3,29$

# AC4P46 Nitritpulverpäckchen-Test

### **Diazotierungs-Methode (Azo)**

0,01 - 0,3 mg/I N (Nitrit als Stickstoffträger)

- 1. Laden Sie die Methode AC4P46, und führen Sie sie aus.
- Befüllen Sie eine saubere AQUAfast 24-mm-Rundküvette, Bestell-Nr. AC2V24, mit 10 ml der Probe. Verschließen Sie das Fläschchen fest mit dem Deckel. Wischen Sie die Außenseite des Fläschchens ab.
- 3. Stellen Sie das Fläschchen in die Halterung in der Probenkammer und schließen Sie die Probenkammertür.
- 4. Drücken Sie die Funktionstaste Blank (Blind), um den Blindwert zu messen.
- 5. Öffnen Sie die Probenkammertür, und nehmen Sie das Fläschchen aus der Halterung.
- 6. Geben Sie den Inhalt eines <u>Nitri 3 Powder Pack</u> direkt aus der Folienpackung in das Fläschchen.
- 7. Verschließen Sie das Fläschchen fest mit dem Deckel, und mischen Sie den Inhalt durch mehrmaliges Schwenken oder Umdrehen. Wischen Sie die Außenseite des Fläschchens ab.
- 8. Warten Sie, bis eine Reaktionszeit von 20 Minuten abgelaufen ist.
- 9. Stellen Sie das Fläschchen in die Halterung in der Probenkammer und schließen Sie die Probenkammertür.
- 10. Tippen Sie auf die Funktionstaste Sample (Probe), um das Ergebnis in mg/l Nitrit als Stickstoffträger anzuzeigen.

- Störungen:
  - o Stark oxidierende und reduzierende Substanzen stören die Messergebnisse.
  - o Kupfer- und Eisen(II)-Ionen verursachen falsch niedrige Messergebnisse.
  - Antimon-, Aurin-, Bismut-, Chloroplatinat-, Eisen-, Blei-, Quecksilber-, Metavanadatund Silberionen stören, indem sie Ausfällungen verursachen.
  - In Proben mit sehr hohen Nitratkonzentrationen (> 100 mg/l N) werden auch geringe Nitrit-Mengen anzutreffen sein. Solch hohe Nitratkonzentrationen scheinen entweder spontan oder während der Reaktionszeit des Tests in geringem Umfang zu Nitrit reduziert zu werden.

# ACD004 Stickstoff-Aufschlussröhrchen-Test (Gesamtstickstoff) für niedrigen Bereich

#### Persulfataufschluss-Methode

0,5 – 25 mg/l N (Nitrit als Stickstoffträger)

- Öffnen Sie zwei <u>TN Hydroxide LR Digestion Vials</u>, und geben Sie den Inhalt eines <u>TN</u> <u>Persulfate Reagent Powder Pack</u> direkt aus der Folie in jedes Fläschchen. Verwenden Sie einen Trichter, um das Reagenz hinzuzufügen. Wischen Sie alle Persulfat-Reagenzien ab, die auf den Deckel oder die Probengefäßgewinde gelangen können.
- 2. Geben Sie 2 ml deionisiertes Wasser in das erste Aufschlussgefäß. (Dies ist das Blindprobengefäß.)
- 3. Geben Sie 2 ml der Probe in das zweite Aufschlussgefäß. (Dies ist das Fläschchen.)
- Verschließen Sie die Fläschchen fest mit den Kappen und schütteln Sie sie mindestens 30 Sekunden lang, um den Inhalt zu mischen. Das Reagenz löst sich möglicherweise nicht vollständig auf.
- 5. Erhitzen Sie die Aufschlussgefäße <u>30 Minuten lang</u> im vorgeheizten Reaktor bei einer Temperatur von 100 °C erhitzen.
- 6. **VORSICHT:** Die Fläschchen werden heiß. Nehmen Sie die Aufschlussgefäße aus dem Reaktor, und lassen Sie sie auf Raumtemperatur abkühlen.
- Öffnen Sie die abgekühlten Aufschlussgefäße, und geben Sie den Inhalt eines <u>TN Reagent</u> <u>A Powder Pack</u> direkt aus der Folie in jedes Fläschchen. Verwenden Sie einen Trichter, um das Reagenz hinzuzufügen.
- 8. Verschließen Sie die Fläschchen fest mit den Kappen, und schütteln Sie sie mindestens 15 Sekunden lang, um den Inhalt zu mischen.
- 9. Warten Sie eine Reaktionszeit von <u>3 Minuten</u> ab.
- Öffnen Sie die Aufschlussgefäße und geben Sie den Inhalt eines <u>TN Reagent B Powder</u> <u>Pack</u> direkt aus der Folie in jedes Fläschchen. Verwenden Sie einen Trichter, um das Reagenz hinzuzufügen.
- Verschließen Sie die Fläschchen fest mit den Kappen, und schütteln Sie sie mindestens 15 Sekunden lang, um den Inhalt zu mischen. Das Reagenz wird sich nicht vollständig auflösen.
- 12. Warten Sie eine Reaktionszeit von <u>2 Minuten</u> ab.
- 13. Öffnen Sie zwei <u>TN Acid LR/HR (Reagent C) Vials</u> und geben Sie 2 ml der aufgeschlossenen, behandelten Blindprobe in das erste Fläschchen. (Dies ist die Blindprobe.)
- 14. Geben Sie in das zweite Fläschchen 2 ml der aufgeschlossenen, behandelten Probe. (Dies ist das Fläschchen.)

- 15. Verschließen Sie die Fläschchen fest mit den Kappen, und drehen Sie sie mindestens 10 Mal vorsichtig um, um den Inhalt zu mischen. Halten Sie das Fläschchen in senkrechter Position mit der Kappe nach oben. Drehen Sie das Fläschchen auf den Kopf. Warten Sie, bis die gesamte Lösung zur Kappe hinuntergelaufen ist. Bringen Sie das Fläschchen zurück in die Normalposition. Warten Sie, bis die gesamte Lösung auf den Boden des Fläschchens fließt. Dieser Vorgang ist ein einmaliges Umdrehen. Ein 10 maliges Umdrehen des Gefäßes dauert etwa 30 Sekunden. Wischen Sie die Außenseiten der Fläschchen ab. VORSICHT: Die Fläschchen werden beim Mischen warm.
- 16. Warten Sie eine Reaktionszeit von <u>5 Minuten</u> ab.
- 17. Laden Sie die Methode ACD004, und führen Sie sie aus.
- Stellen Sie das Blindprobengefäß in die Halterung in der Probenkammer, und schließen Sie die Probenkammert
  ür.
- 19. Drücken Sie die Funktionstaste Blank (Blind), um den Blindwert zu messen.
- 20. Öffnen Sie die Probenkammertür. Entnehmen Sie das Blindprobengefäß aus der Halterung.
- 21. Stellen Sie das Fläschchen in die Halterung in der Probenkammer, und schließen Sie die Probenkammertür.
- Tippen Sie auf die Funktionstaste Sample (Probe), um das Ergebnis in mg/l Stickstoff anzuzeigen.

- Während des gesamten Verfahrens sollten angemessene Sicherheitsvorkehrungen getroffen und eine gute Labortechnik angewandt werden.
- Die Volumina für Proben und Blindprobe sollten immer mit 2-ml-Vollpipetten (Klasse A) gemessen werden.
- Für jeden Probensatz ist eine Blindprobe ausreichend. Nach der Blindprobenmessung können mehrere Proben gemessen werden.
- Die Fläschchen müssen nach exakt 30 Minuten aus dem Reaktor genommen werden.
- Große Mengen an stickstofffreien, organischen Verbindungen, die in einigen Wasserproben enthalten sind, können die Wirksamkeit des Aufschlusses durch Reaktion mit dem Persulfat-Reagenz verringern. Samples (Proben), die bekanntermaßen große Mengen organischer Verbindungen enthalten, müssen verdünnt und der Aufschluss und die Messung wiederholt werden, um die Wirksamkeit des Aufschlusses zu überprüfen.
- Anwendung: für Wasser, Abwasser und Meerwasser.
- Störsubstanzen, die zu einer Konzentrationsänderung von 10 % führten: Bromid bei mehr als 60 mg/l und Chlorid bei mehr als 1.000 mg/l führen zu falsch erhöhten Ergebnissen.

# ACD007 Stickstoff-Aufschlussröhrchen-Test (Gesamtstickstoff) für hohen Bereich

#### Persulfataufschluss-Methode

- 5 150 mg/l N (Nitrit als Stickstoffträger)
- Öffnen Sie zwei <u>TN Hydroxid HR-Aufschlussgefäße</u> und geben Sie den Inhalt eines <u>TN</u> <u>Persulfat-Reagenzpulverpäckchen</u> direkt aus der Folie in jedes Fläschchen. Verwenden Sie einen Trichter, um das Reagenz hinzuzufügen. Wischen Sie alle Persulfat-Reagenzien ab, die auf den Deckel oder die Probengefäßgewinde gelangen können.
- 2. Geben Sie 0,5 ml deionisiertes Wasser in das erste Aufschlussgefäß. (Dies ist das Blindprobengefäß.)
- 3. Geben Sie 0,5 ml der Probe in das zweite Aufschlussgefäß. (Dies ist das Fläschchen.)
- Verschließen Sie die Fläschchen fest mit den Kappen, und schütteln Sie sie mindestens 30 Sekunden lang, um den Inhalt zu mischen. Das Reagenz löst sich möglicherweise nicht vollständig auf.
- 5. Erhitzen Sie die Aufschlussgefäße <u>30 Minuten lang</u> im vorgeheizten Reaktor bei einer Temperatur von 100 °C erhitzen.
- 6. **VORSICHT:** Die Fläschchen werden heiß. Nehmen Sie die Aufschlussgefäße aus dem Reaktor, und lassen Sie sie auf Raumtemperatur abkühlen.
- Öffnen Sie die abgekühlten Aufschlussgefäße, und geben Sie den Inhalt eines <u>TN Reagent</u> <u>A Powder Pack</u> direkt aus der Folie in jedes Fläschchen. Verwenden Sie einen Trichter, um das Reagenz hinzuzufügen.
- 8. Verschließen Sie die Fläschchen fest mit den Kappen, und schütteln Sie sie mindestens 15 Sekunden lang, um den Inhalt zu mischen.
- 9. Warten Sie eine Reaktionszeit von <u>3 Minuten</u> ab.
- Öffnen Sie die Aufschlussgefäße, und geben Sie den Inhalt eines <u>TN Reagent B Powder</u> <u>Pack</u> direkt aus der Folie in jedes Fläschchen. Verwenden Sie einen Trichter, um das Reagenz hinzuzufügen.
- Verschließen Sie die Fläschchen fest mit den Kappen, und schütteln Sie sie mindestens 15 Sekunden lang, um den Inhalt zu mischen. Das Reagenz wird sich nicht vollständig auflösen.
- 12. Warten Sie eine Reaktionszeit von <u>2 Minuten</u> ab.
- Öffnen Sie zwei <u>TN Acid LR/HR (Reagent C) Vials</u> und geben Sie 2 ml der aufgeschlossenen, behandelten Blindprobe in das erste Fläschchen. (Dies ist die Blindprobe.)
- 14. Geben Sie in das zweite Fläschchen 2 ml der aufgeschlossenen, behandelten Probe. (Dies ist das Fläschchen.)

- 15. Verschließen Sie die Fläschchen fest mit den Kappen, und drehen Sie sie mindestens 10 Mal vorsichtig um, um den Inhalt zu mischen. Halten Sie das Fläschchen in senkrechter Position mit der Kappe nach oben. Drehen Sie das Fläschchen auf den Kopf. Warten Sie, bis die gesamte Lösung zur Kappe hinuntergelaufen ist. Bringen Sie das Fläschchen zurück in die Normalposition. Warten Sie, bis die gesamte Lösung auf den Boden des Fläschchens fließt. Dieser Vorgang ist ein einmaliges Umdrehen. Ein 10 maliges Umdrehen des Gefäßes dauert etwa 30 Sekunden. Wischen Sie die Außenseiten der Fläschchen ab. VORSICHT: Die Fläschchen werden beim Mischen warm.
- 16. Warten Sie eine Reaktionszeit von <u>5 Minuten</u> ab.
- 17. Laden Sie die Methode ACD007, und führen Sie sie aus.
- Stellen Sie das Blindprobengefäß in die Halterung in der Probenkammer, und schließen Sie die Probenkammert
  ür.
- 19. Drücken Sie die Funktionstaste Blank (Blind), um den Blindwert zu messen.
- 20. Öffnen Sie den Probenkammerdeckel. Entnehmen Sie das Blindprobengefäß aus der Halterung.
- 21. Stellen Sie das Fläschchen in die Halterung in der Probenkammer, und schließen Sie die Probenkammertür.
- Tippen Sie auf die Funktionstaste Sample (Probe), um das Ergebnis in mg/l Stickstoff anzuzeigen.

- Während des gesamten Verfahrens sollten angemessene Sicherheitsvorkehrungen getroffen und eine gute Labortechnik angewandt werden.
- Die Volumina für Proben und Blindprobe sollten immer mit 2-ml-Vollpipetten (Klasse A) gemessen werden.
- Für jeden Probensatz ist eine Blindprobe ausreichend. Nach der Blindprobenmessung können mehrere Proben gemessen werden.
- Die Fläschchen müssen nach exakt 30 Minuten aus dem Reaktor genommen werden.
- Große Mengen an stickstofffreien, organischen Verbindungen, die in einigen Wasserproben enthalten sind, können die Wirksamkeit des Aufschlusses durch Reaktion mit dem Persulfat-Reagenz verringern. Samples (Proben), die bekanntermaßen große Mengen organischer Verbindungen enthalten, müssen verdünnt und der Aufschluss und die Messung wiederholt werden, um die Wirksamkeit des Aufschlusses zu überprüfen.
- Anwendung: für Wasser, Abwasser und Meerwasser.
- Störsubstanzen, die zu einer Konzentrationsänderung von 10 % führten: Bromid bei mehr als 60 mg/l und Chlorid bei mehr als 1.000 mg/l führen zu falsch erhöhten Ergebnissen.

# AC3048 Ozontabletten-Test

#### DPD-Methode

0,02 - 2 mg/l O<sub>3</sub>

#### Ozonmessung in Abwesenheit von Chlor

- 1. Laden Sie die Methode AC3048, und führen Sie sie aus.
- Befüllen Sie eine saubere AQUAfast 24-mm-Rundküvette, Bestell-Nr. AC2V24, mit 10 ml der Probe. Verschließen Sie das Fläschchen fest mit dem Deckel. Wischen Sie die Außenseite des Fläschchens ab.
- 3. Stellen Sie das Fläschchen in die Halterung in der Probenkammer, und schließen Sie die Probenkammertür.
- 4. Drücken Sie die Funktionstaste Blank (Blind), um den Blindwert zu messen.
- 5. Öffnen Sie die Probenkammertür, und nehmen Sie das Fläschchen aus der Halterung.
- 6. Entleeren Sie das Fläschchen, so dass einige Tropfen der Probe im Fläschchen zurückbleiben.
- Geben Sie eine <u>DPD No. 1 Tablet</u> und eine <u>DPD No. 3 Tablet</u> direkt aus der Folienverpackung in das Fläschchen. Zerdrücken Sie die Tabletten mit einem sauberen Rührstab.
- 8. Füllen Sie die Probe bis zur 10-ml-Marke auf dem Fläschchen auf.
- Verschließen Sie das Fläschchen fest mit dem Deckel, und mischen Sie den Inhalt durch mehrmaliges Schwenken oder Umdrehen, bis sich die Tabletten aufgelöst haben. Wischen Sie die Außenseite des Fläschchens ab.
- 10. Stellen Sie das Fläschchen in die Halterung in der Probenkammer, und Schließen Sie den Probenkammertür.
- 11. Warten Sie eine Reaktionszeit von <u>2 Minuten</u> ab.
- 12. Tippen Sie auf die Funktionstaste Sample (Probe), um das Ergebnis in mg/l Ozon anzuzeigen.

#### Ozonmessung in Anwesenheit von Chlor

- 1. Laden Sie die Methode AC3048, und führen Sie sie aus.
- Befüllen Sie eine saubere AQUAfast 24-mm-Rundküvette, Bestell-Nr. AC2V24, mit 10 ml der Probe. Verschließen Sie das Fläschchen fest mit dem Deckel. Wischen Sie die Außenseite des Fläschchens ab.
- 3. Stellen Sie das Fläschchen in die Halterung in der Probenkammer, und Schließen Sie den Probenkammertür.
- 4. Drücken Sie die Funktionstaste Blank (Blind), um den Blindwert zu messen.
- 5. Öffnen Sie die Probenkammertür, und nehmen Sie das Fläschchen aus der Halterung.
- 6. Entleeren Sie das Fläschchen, so dass einige Tropfen der Probe im Fläschchen zurückbleiben.

- 7. Geben Sie eine <u>DPD No. 1 Tablet</u> und eine <u>DPD No. 3 Tablet</u> direkt aus der Folienverpackung in das Fläschchen. Zerdrücken Sie die Tabletten mit einem sauberen Rührstab.
- Befüllen Sie eine zweite, saubere AQUAfast 24-mm-Rundküvette mit 10 ml der Probe. Geben Sie eine <u>Glycine Tablet</u> direkt aus dem Folienbeutel in das Fläschchen. Zerdrücken Sie die Tablette mit einem sauberen Rührstab. Verschließen Sie das Fläschchen fest mit dem Deckel, und mischen Sie den Inhalt durch mehrmaliges Schwenken, bis sich die Tablette aufgelöst hat.
- Füllen Sie den Inhalt des Fläschchens mit der Glycinlösung in das Fläschchen mit den DPD-Nr.-1- und DPD-Nr.-3-Tabletten um.
- Verschließen Sie das Fläschchen fest mit dem Deckel, und mischen Sie den Inhalt durch mehrmaliges Schwenken oder Umdrehen, bis sich die Tabletten aufgelöst haben. Wischen Sie die Außenseite des Fläschchens ab.
- 11. Stellen Sie das Fläschchen in die Halterung in der Probenkammer, und Schließen Sie den Probenkammertür.
- 12. Warten Sie eine Reaktionszeit von 2 Minuten ab.
- 13. Tippen Sie auf die Funktionstaste Sample (Probe), um das Ergebnis in mg/l Ozon anzuzeigen.

#### Hinweise:

 Reinigung der Fläschchen: Da viele Haushaltsreiniger (z. B. Geschirrspülmittel) reduzierende Stoffe enthalten, kann es nach deren Anwendung bei der Bestimmung von Ozon zu Minderbefunden kommen. Um diesen Messfehler auszuschließen, sollten die Glasgefäße chlorzehrungsfrei sein.

Vorbereitung: Legen Sie dazu alle entsprechenden Glasgefäße eine Stunde lang in Natriumhypochloritlösung (0,1 g/l) ein, und spülen Sie sie dann gründlich mit deionisiertem Wasser ab.

- Probenvorbereitung: Bei der Vorbereitung der Probe ist das Entweichen von Ozon (z. B. beim Pipettieren oder Schütteln) zu vermeiden. Die Analyse muss unmittelbar nach der Probenahme erfolgen.
- Die DPD-Farbentwicklung erfolgt bei einem pH-Wert von 6,2 bis 6,5. Die Reagenztablette enthält daher einen Puffer zur pH-Einstellung. Stark alkalische oder saure Wasserproben müssen vor Zugabe der Tablette auf einen pH-Wert zwischen 6 und 7 gebracht werden. (Dazu 0,5 mol/l Schwefelsäure bzw. 1 mol/l Natronlauge verwenden.)
- Überschreitung des Messbereichs: Konzentrationen von mehr als 6 mg/l Ozon können zu Ergebnissen bis 0 mg/l führen. In diesem Fall muss die Wasserprobe mit ozonfreiem Wasser verdünnt werden. Dazu sollten 10 ml der verdünnten Probe mit dem Reagenz versetzt und die Messung wiederholt werden.
- Oxidationsmittel wie Brom oder Chlor sind Störfaktoren, da sie wie Ozon reagieren.

# AC2001 pH-Tablettentest

#### Phenolrot-Methode

pH 6,5 – 8,4

- 1. Laden Sie die Methode AC2001, und führen Sie sie aus.
- Befüllen Sie eine saubere AQUAfast 24-mm-Rundküvette, Bestell-Nr. AC2V24, mit 10 ml der Probe. Verschließen Sie das Fläschchen fest mit dem Deckel. Wischen Sie die Außenseite des Fläschchens ab.
- 3. Stellen Sie das Fläschchen in die Halterung in der Probenkammer und schließen Sie die Probenkammertür.
- 4. Drücken Sie die Funktionstaste Blank (Blind), um den Blindwert zu messen.
- 5. Öffnen Sie die Probenkammertür, und nehmen Sie das Fläschchen aus der Halterung.
- 6. Geben Sie eine <u>Phenol Red Tablet</u> direkt aus dem Folienbeutel in das Fläschchen. Zerdrücken Sie die Tablette mit einem sauberen Rührstab.
- Verschließen Sie das Fläschchen fest mit dem Deckel, und mischen Sie den Inhalt durch mehrmaliges Schwenken oder Umdrehen, bis sich die Tablette aufgelöst hat. Wischen Sie die Außenseite des Fläschchens ab.
- 8. Stellen Sie das Fläschchen in die Halterung in der Probenkammer und schließen Sie die Probenkammertür.
- 9. Tippen Sie auf die Funktionstaste Sample (Probe), um das Ergebnis als pH-Wert anzuzeigen.

#### Hinweise:

- Wasserproben mit niedriger Alkalinität-m (unter 35 mg/l CaCO<sub>3</sub>) können falsche pH-Werte ergeben.
- pH-Werte unter 6,5 und über 8,4 können zu Ergebnissen innerhalb des Messbereichs führen. Ein Plausibilitätstest (pH-Meter) wird empfohlen.
- Die Genauigkeit der kolorimetrischen Bestimmung von pH-Werten hängt von verschiedenen Randbedingungen ab (Pufferkapazität der Probe, Salzgehalt usw.).
- Salzfehler: Korrektur der Prüfergebnisse (Durchschnittswerte) für Proben mit Salzgehalten von:

| Indikator | Salzgehalt      |                 |                 |  |
|-----------|-----------------|-----------------|-----------------|--|
| Phenolrot | 1-molare Lösung | 2-molare Lösung | 3-molare Lösung |  |
|           | – 0,21          | – 0,26          | – 0,29          |  |

 Die Werte von Parson und Douglas (1926) beruhen auf der Verwendung von Clark- und Lubs-Puffern. 1 Mol NaCl = 58,4 g/l = 5,8 %

# AC3001 Flüssigreagenz-pH-Test

#### Phenolrot-Methode

pH 6,5 – 8,4

- 1. Laden Sie die Methode AC3001, und führen Sie sie aus.
- Befüllen Sie eine saubere AQUAfast 24-mm-Rundküvette, Bestell-Nr. AC2V24, mit 10 ml der Probe. Verschließen Sie das Fläschchen fest mit dem Deckel. Wischen Sie die Außenseite des Fläschchens ab.
- Stellen Sie das Fläschchen in die Halterung in der Probenkammer und schließen Sie die Probenkammert
  ür.
- 4. Drücken Sie die Funktionstaste Blank (Blind), um den Blindwert zu messen.
- 5. Öffnen Sie die Probenkammertür, und nehmen Sie das Fläschchen aus der Halterung.
- 6. Geben Sie 6 Tropfen <u>Phenol Red Solution</u> in das Fläschchen. Fügen Sie Tropfen in gleicher Größe hinzu, indem Sie die Flasche senkrecht halten und langsam zusammendrücken.
- 7. Verschließen Sie das Fläschchen fest mit dem Deckel, und mischen Sie den Inhalt durch mehrmaliges Schwenken oder Umdrehen. Wischen Sie die Außenseite des Fläschchens ab.
- 8. Stellen Sie das Fläschchen in die Halterung in der Probenkammer und schließen Sie die Probenkammertür.
- 9. Tippen Sie auf die Funktionstaste Sample (Probe), um das Ergebnis als pH-Wert anzuzeigen.

- Bei der Analyse von gechlortem Wasser kann der Restchlorgehalt die Farbreaktion des Flüssigreagenz beeinflussen. Dies lässt sich vermeiden (ohne die pH-Messung zu beeinträchtigen), indem man der Probe vor Zugabe der Phenolrotlösung eine geringe Menge Natriumthiosulfat-Kristall (Na<sub>2</sub>S<sub>2</sub>O<sub>3</sub> • 5 H<sub>2</sub>O) zusetzt. Phenolrot-Tabletten enthalten bereits Thiosulfat.
- Aufgrund unterschiedlicher Tropfengrößen können die Ergebnisse im Vergleich zu Tabletten abweichende Genauigkeiten aufweisen. Dies kann durch die Verwendung einer Pipette (0,18 ml Phenolrotlösung entsprechen 6 Tropfen) minimiert werden.
- Verschließen Sie die Flasche nach dem Gebrauch wieder fest mit der Verschlusskappe.
- Lagern Sie die Phenolrot-Lösung an einem kühlen, trockenen Ort, am besten bei 6 bis 10 °C.

# AC2095-WA Phosphattabletten-Test (Ortho) für niedrigen Bereich

#### Phosphomolybdänsäure/Ascorbinsäure

0,05 – 4 mg/l PO<sub>4</sub>

- 1. Laden Sie die Methode AC2095, und führen Sie sie aus.
- Befüllen Sie eine saubere AQUAfast 24-mm-Rundküvette, Bestell-Nr. AC2V24, mit 10 ml der Probe. Verschließen Sie das Fläschchen fest mit dem Deckel. Wischen Sie die Außenseite des Fläschchens ab.
- 3. Stellen Sie das Fläschchen in die Halterung in der Probenkammer und schließen Sie die Probenkammertür.
- 4. Drücken Sie die Funktionstaste Blank (Blind), um den Blindwert zu messen.
- 5. Öffnen Sie die Probenkammertür, und nehmen Sie das Fläschchen aus der Halterung.
- 6. Geben Sie eine <u>Phosphate No. 1 LR Tablet</u> direkt aus dem Folienbeutel in das Fläschchen. Zerdrücken Sie die Tablette mit einem sauberen Rührstab.
- 7. Geben Sie eine <u>Phosphate No. 2 LR Tablet</u> direkt aus dem Folienbeutel in das Fläschchen. Zerdrücken Sie die Tablette mit einem sauberen Rührstab.
- Verschließen Sie das Fläschchen fest mit dem Deckel, und mischen Sie den Inhalt durch mehrmaliges Schwenken oder Umdrehen, bis sich die Tabletten aufgelöst haben. Wischen Sie die Außenseite des Fläschchens ab.
- 9. Warten Sie eine Reaktionszeit von <u>10 Minuten</u> ab.
- 10. Stellen Sie das Fläschchen in die Halterung in der Probenkammer und schließen Sie die Probenkammertür.
- 11. Tippen Sie auf die Funktionstaste Sample (Probe), um das Ergebnis in mg/l ortho-Phosphat anzuzeigen.

- Es reagieren nur ortho-Phosphat-Ionen.
- Die Tabletten müssen in der richtigen Reihenfolge zugegeben werden.
- Die Testprobe sollte einen pH-Wert zwischen 6 und 7 aufweisen.
- Störfaktoren: Höhere Konzentrationen von Cu, Ni, Cr (III), V (V) und W (VI) sind aufgrund ihrer Farbentwicklung Störfaktoren. Silikate stören die Messung nicht (sie werden durch die in den Tabletten enthaltene Zitronensäure maskiert).
- Ortho-Phosphationen reagieren mit dem Reagenz und erzeugen eine intensive blaue Färbung.
- Phosphate in organischer und kondensierter anorganischer Form (Meta-, Pyro- und Polyphosphate) müssen vor der Analyse in ortho-Phosphationen umgewandelt werden. Die Vorbehandlung der Probe mit Säure und Hitze schafft die erforderlichen Voraussetzungen für die Hydrolyse kondensierter anorganischer Formen. Organisch gebundene Phosphate werden durch Erhitzen mit Säure und Persulfat in ortho-Phosphationen umgewandelt. Die Menge organisch gebundener Phosphate kann folgendermaßen berechnet werden: mg/l Phosphat (organisch) = mg/l Phosphat (gesamt) - mg/l Phosphat (säurehydrolysierbar)
- ortho-Phosphat = reaktives Phosphor

# AC2096 Phosphattabletten-Test (Ortho) für hohen Bereich

#### Vanado-Molybdat-Verfahren

1 – 80 mg/l PO<sub>4</sub>

- 1. Laden Sie die Methode AC2096, und führen Sie sie aus.
- Befüllen Sie eine saubere AQUAfast 24-mm-Rundküvette, Bestell-Nr. AC2V24, mit 10 ml der Probe. Verschließen Sie das Fläschchen fest mit dem Deckel. Wischen Sie die Außenseite des Fläschchens ab.
- 3. Stellen Sie das Fläschchen in die Halterung in der Probenkammer und schließen Sie die Probenkammertür.
- 4. Drücken Sie die Funktionstaste Blank (Blind), um den Blindwert zu messen.
- 5. Öffnen Sie die Probenkammertür, und nehmen Sie das Fläschchen aus der Halterung.
- 6. Geben Sie eine <u>Phosphate HR P1 Tablet</u> direkt aus dem Folienbeutel in das Fläschchen. Zerdrücken Sie die Tablette mit einem sauberen Rührstab.
- 7. Geben Sie eine <u>Phosphate HR P2 Tablet</u> direkt aus dem Folienbeutel in das Fläschchen. Zerdrücken Sie die Tablette mit einem sauberen Rührstab.
- Verschließen Sie das Fläschchen fest mit dem Deckel, und mischen Sie den Inhalt durch mehrmaliges Schwenken oder Umdrehen, bis sich die Tabletten aufgelöst haben. Wischen Sie die Außenseite des Fläschchens ab.
- 9. Warten Sie eine Reaktionszeit von <u>10 Minuten</u> ab.
- 10. Stellen Sie das Fläschchen in die Halterung in der Probenkammer und schließen Sie die Probenkammertür.
- 11. Tippen Sie auf die Funktionstaste Sample (Probe), um das Ergebnis in mg/l ortho-Phosphat anzuzeigen.

- Für Proben unter 5 mg/l PO<sub>4</sub> wird empfohlen, die Probe mit dem AC2095-WA Phosphattabletten-Test (ortho) für den niedrigen Bereich zu analysieren.
- Es reagieren nur ortho-Phosphat-Ionen.
- Phosphate in organischer und kondensierter anorganischer Form (Meta-, Pyro- und Polyphosphate) müssen vor der Analyse in ortho-Phosphationen umgewandelt werden. Die Vorbehandlung der Probe mit Säure und Hitze schafft die erforderlichen Voraussetzungen für die Hydrolyse kondensierter anorganischer Formen. Organisch gebundene Phosphate werden durch Erhitzen mit Säure und Persulfat in ortho-Phosphationen umgewandelt. Die Menge organisch gebundener Phosphate kann folgendermaßen berechnet werden: mg/l Phosphat (organisch) = mg/l Phosphat (gesamt) - mg/l Phosphat (säurehydrolysierbar)
- Die ortho-Phosphationen reagieren mit dem Vanadat-Molybdat-Reagenz unter sauren Bedingungen und bilden eine gelb gefärbte Substanz.
- ortho-Phosphat = reaktives Phosphor

# AC4P95 Phosphat-Pulverpäckchen-Test (Ortho)

#### Phosphormolybdänsäure/Ascorbinsäure

0,06 – 2,5 mg/l PO<sub>4</sub>

- 1. Laden Sie die Methode AC4P95, und führen Sie sie aus.
- Befüllen Sie eine saubere AQUAfast 24-mm-Rundküvette, Bestell-Nr. AC2V24, mit 10 ml der Probe. Verschließen Sie das Fläschchen fest mit dem Deckel. Wischen Sie die Außenseite des Fläschchens ab.
- 3. Stellen Sie das Fläschchen in die Halterung in der Probenkammer und schließen Sie die Probenkammertür.
- 4. Drücken Sie die Funktionstaste Blank (Blind), um den Blindwert zu messen.
- 5. Öffnen Sie die Probenkammertür, und nehmen Sie das Fläschchen aus der Halterung.
- 6. Fügen Sie den Inhalt eines <u>Phosphat Rgt. F10 Powder Pack</u> direkt aus der Folienpackung in das Fläschchen.
- Verschließen Sie das Fläschchen fest mit der Kappe, und mischen Sie den Inhalt für 10 bis 15 Sekunden durch mehrmaliges Schwenken oder Umdrehen. Das Pulver löst sich möglicherweise nicht vollständig auf. Wischen Sie die Außenseite des Fläschchens ab.
- 8. Warten Sie eine Reaktionszeit von <u>2 Minuten</u> ab.
- 9. Stellen Sie das Fläschchen in die Halterung in der Probenkammer und schließen Sie die Probenkammertür.
- 10. Tippen Sie auf die Funktionstaste Sample (Probe), um das Ergebnis in mg/l ortho-Phosphat anzuzeigen.

- Ortho-Phosphationen reagieren mit dem Reagenz und erzeugen eine intensive blaue Färbung.
- Phosphate in organischer und kondensierter anorganischer Form (Meta-, Pyro- und Polyphosphate) müssen vor der Analyse in ortho-Phosphationen umgewandelt werden. Die Vorbehandlung der Probe mit Säure und Hitze schafft die erforderlichen Voraussetzungen für die Hydrolyse kondensierter anorganischer Formen. Organisch gebundene Phosphate werden durch Erhitzen mit Säure und Persulfat in ortho-Phosphationen umgewandelt. Die Menge organisch gebundener Phosphate kann folgendermaßen berechnet werden: mg/l Phosphat (organisch) = mg/l Phosphat (gesamt) - mg/l Phosphat (säurehydrolysierbar)
- Anwendung: für Wasser, Abwasser und Meerwasser.
- Stark gepufferte Proben oder Proben mit extremen pH-Werten sollten mit 1 mol/l Salzsäure oder 1 mol/l Natriumhydroxid (Natronlauge) auf einen pH-Wert zwischen 2 und 10 gebracht werden.
- ortho-Phosphat = reaktives Phosphor
- Störfaktoren: Starke Trübungen können zu ungleichmäßigen Ergebnissen führen.

| Störung   | Störungsgrad            |  | Störung                |
|-----------|-------------------------|--|------------------------|
| Aluminium | Mehr als 200 mg/l       |  | Nickel                 |
| Arsenat   | bei jeder Konzentration |  | Kieselsäure (Siliziumd |
| Chrom     | mehr als 100 mg/l       |  | Silikat                |
| Kupfer    | mehr als 10 mg/l        |  | Sulfid                 |
| Eisen     | mehr als 100 mg/l       |  | Zink                   |

| Störung                      | Störungsgrad            |
|------------------------------|-------------------------|
| Nickel                       | mehr als 300 mg/l       |
| Kieselsäure (Siliziumdioxid) | mehr als 50 mg/l        |
| Silikat                      | mehr als 10 mg/l        |
| Sulfid                       | bei jeder Konzentration |
| Zink                         | mehr als 80 mg/l        |

# ACR095 Phosphat-Reaktionsröhrchen-Test (Ortho)

#### Phosphormolybdänsäure/Ascorbinsäure-Methode

0,06 – 5 mg/l PO<sub>4</sub>

- 1. Laden Sie die Methode ACR095, und führen Sie sie aus.
- 2. Öffnen Sie ein16 mm <u>PO4-P Dilution Tube</u>, und geben Sie 5 ml der Probe hinzu. Wischen Sie die Außenseite des Fläschchens ab.
- 3. Stellen Sie das Fläschchen in die Halterung in der Probenkammer und schließen Sie die Probenkammertür.
- 4. Drücken Sie die Funktionstaste Blank (Blind), um den Blindwert zu messen.
- 5. Öffnen Sie den Probenkammerdeckel. Entnehmen Sie das Fläschchen aus der Halterung.
- 6. Fügen Sie den Inhalt eines <u>Phosphat-Rgt. F10 Powder Pack</u> direkt aus der Folienpackung in das Fläschchen. Verwenden Sie einen Trichter, um das Reagenz hinzuzufügen.
- Verschließen Sie das Fläschchen fest mit der Kappe, und mischen Sie den Inhalt 10 bis 15 Sekunden durch Schwenken. Das Reagenz wird sich nicht vollständig auflösen. Wischen Sie die Außenseite des Fläschchens ab.
- 8. Warten Sie eine Reaktionszeit von <u>2 Minuten</u> ab.
- 9. Stellen Sie das Fläschchen in die Halterung in der Probenkammer und schließen Sie die Probenkammertür.
- 10. Tippen Sie auf die Funktionstaste Sample (Probe), um das Ergebnis in mg/l ortho-Phosphat anzuzeigen.

- Ortho-Phosphationen reagieren mit dem Reagenz und erzeugen eine intensive blaue Färbung.
- Phosphate in organischer und kondensierter anorganischer Form (Meta-, Pyro- und Polyphosphate) müssen vor der Analyse in ortho-Phosphationen umgewandelt werden. Die Vorbehandlung der Probe mit Säure und Hitze schafft die erforderlichen Voraussetzungen für die Hydrolyse kondensierter anorganischer Formen. Organisch gebundene Phosphate werden durch Erhitzen mit Säure und Persulfat in ortho-Phosphationen umgewandelt. Die Menge organisch gebundener Phosphate kann folgendermaßen berechnet werden: mg/l Phosphat (organisch) = mg/l Phosphat (gesamt) - mg/l Phosphat (säurehydrolysierbar)
- Anwendung: für Wasser, Abwasser und Meerwasser.
- Stark gepufferte Proben oder Proben mit extremen pH-Werten sollten mit 1 mol/l Salzsäure oder 1 mol/l Natriumhydroxid (Natronlauge) auf einen pH-Wert zwischen 2 und 10 gebracht werden.
- ortho-Phosphat = reaktives Phosphor
- Störfaktoren: Starke Trübungen können zu ungleichmäßigen Ergebnissen führen.

| Störung                         | Störungsgrad      |  |
|---------------------------------|-------------------|--|
| Aluminium                       | mehr als 200 mg/l |  |
| Arsenat bei jeder Konzentration |                   |  |
| Chrom                           | mehr als 100 mg/l |  |
| Kupfer mehr als 10 mg/l         |                   |  |
| Eisen                           | mehr als 100 mg/l |  |

| Störung                      | Störungsgrad            |
|------------------------------|-------------------------|
| Nickel                       | mehr als 300 mg/l       |
| Kieselsäure (Siliziumdioxid) | mehr als 50 mg/l        |
| Silikat                      | mehr als 10 mg/l        |
| Sulfid                       | bei jeder Konzentration |
| Zink                         | mehr als 80 mg/l        |

# ACD095 Phosphat-Aufschlussröhrchen-Test (Gesamtphosphat)

#### Persulfataufschluss/Ascorbinsäure-Methode

0,02 – 1,1 mg/l P (Phosphat als Phosphorträger)

- 1. Öffnen Sie ein 16 mm <u>PO4-P Acid Reagent Digestion Tube</u>, und geben Sie 5 ml der Probe hinzu.
- Geben Sie den Inhalt eines <u>Potassium Persulfate F10 Powder Pack</u> direkt aus der Folienpackung in das Fläschchen. Verwenden Sie einen Trichter, um das Reagenz hinzuzufügen.
- 3. Verschließen Sie das Fläschchen fest mit der Kappe, und mischen Sie den Inhalt durch mehrmaliges Umdrehen.
- 4. Die Fläschchen <u>30 Minuten lang</u> bei einer Temperatur von 100 °C im vorgeheizten Reaktor erhitzen.
- 5. **VORSICHT:** Das Fläschchen wird heiß. Nehmen Sie das Fläschchen aus dem Reaktor, und lassen Sie es auf Raumtemperatur abkühlen.
- Öffnen Sie das gekühlte Aufschlussgefäß, und geben Sie 2 ml <u>1.54 N Sodium Hydroxide</u> <u>Solution</u> in das Gefäß.
- 7. Verschließen Sie das Fläschchen fest mit der Kappe, und mischen Sie den Inhalt vorsichtig durch mehrmaliges Umdrehen. Wischen Sie die Außenseite des Fläschchens ab.
- 8. Laden Sie die Methode ACD095, und führen Sie sie aus.
- 9. Stellen Sie das Fläschchen in die Halterung in der Probenkammer und schließen Sie die Probenkammertür.
- 10. Drücken Sie die Funktionstaste Blank (Blind), um den Blindwert zu messen.
- 11. Öffnen Sie den Probenkammerdeckel. Entnehmen Sie das Fläschchen aus der Halterung.
- 12. Fügen Sie den Inhalt eines <u>Phosphat-Rgt.- F10 Powder Pack</u> direkt aus der Folienpackung in das Fläschchen. Verwenden Sie einen Trichter, um das Reagenz hinzuzufügen.
- Verschließen Sie das Fläschchen fest mit der Kappe, und mischen Sie den Inhalt 10 bis 15 Sekunden durch Schwenken. Das Reagenz wird sich nicht vollständig auflösen. Wischen Sie die Außenseite des Fläschchens ab.
- 14. Warten Sie eine Reaktionszeit von <u>2 Minuten</u> ab.
- 15. Stellen Sie das Fläschchen in die Halterung in der Probenkammer und schließen Sie die Probenkammertür.
- 16. Drücken Sie die Funktionstaste Sample (Probe), um das Ergebnis in mg/l Gesamtphosphat als Phosphor (P) anzuzeigen.

#### Hinweise:

- Während des gesamten Verfahrens sollten angemessene Sicherheitsvorkehrungen getroffen und eine gute Labortechnik angewandt werden.
- Ortho-Phosphationen reagieren mit dem Reagenz und erzeugen eine intensive blaue Färbung.
- Phosphate in organischer und kondensierter anorganischer Form (Meta-, Pyro- und Polyphosphate) müssen vor der Analyse in ortho-Phosphationen umgewandelt werden. Die Vorbehandlung der Probe mit Säure und Hitze schafft die erforderlichen Voraussetzungen für die Hydrolyse kondensierter anorganischer Formen. Organisch gebundene Phosphate werden durch Erhitzen mit Säure und Persulfat in ortho-Phosphationen umgewandelt. Die Menge organisch gebundener Phosphate kann folgendermaßen berechnet werden: mg/l Phosphat (organisch) = mg/l Phosphat (gesamt) - mg/l Phosphat (säurehydrolysierbar)
- Anwendung: für Wasser, Abwasser und Meerwasser.
- Stark gepufferte Proben oder Proben mit extremen pH-Werten sollten mit 1 mol/l Salzsäure oder 1 mol/l Natriumhydroxid (Natronlauge) auf einen pH-Wert zwischen 2 und 10 gebracht werden.
- ortho-Phosphat = reaktives Phosphor

| Störsubstanz                 | Störungsgrad            |
|------------------------------|-------------------------|
| Aluminium                    | mehr als 200 mg/l       |
| Arsenat                      | bei jeder Konzentration |
| Chrom                        | mehr als 100 mg/l       |
| Kupfer                       | mehr als 10 mg/l        |
| Eisen                        | mehr als 100 mg/l       |
| Nickel                       | mehr als 300 mg/l       |
| Kieselsäure (Siliziumdioxid) | mehr als 50 mg/l        |
| Silikat                      | mehr als 10 mg/l        |
| Sulfid                       | bei jeder Konzentration |
| Zink                         | mehr als 80 mg/l        |

• Störungen: Starke Trübungen können zu ungleichmäßigen Ergebnissen führen.

• Umrechnungen:  $mg/I PO_4 = mg/I P \times 3,07$  $mg/I P_2O_5 = mg/I P \times 2,29$ 

## ACD095AH Phosphat-Aufschlussröhrchen-Test (säurehydrolysierbar)

#### Säureaufschluss/Ascorbinsäure-Methode

0,02 – 1,6 mg/l P (Phosphat als Phosphorträger)

- 1. Öffnen Sie ein 16 mm <u>PO4-P Acid Reagent Digestion Tube</u>, und geben Sie 5 ml der Probe hinzu.
- 2. Verschließen Sie das Fläschchen fest mit der Kappe, und mischen Sie den Inhalt vorsichtig durch mehrmaliges Umdrehen.
- Die Fläschchen <u>30 Minuten lang</u> bei einer Temperatur von 100 °C im vorgeheizten Reaktor erhitzen.
- 4. **VORSICHT:** Das Fläschchen wird heiß. Nehmen Sie das Fläschchen aus dem Reaktor, und lassen Sie es auf Raumtemperatur abkühlen.
- 5. Öffnen Sie das gekühlte Aufschlussgefäß und geben Sie 2 ml <u>1.00 N Sodium Hydroxide</u> <u>Solution</u> in das Gefäß.
- 6. Verschließen Sie das Fläschchen fest mit der Kappe, und mischen Sie den Inhalt vorsichtig durch mehrmaliges Umdrehen. Wischen Sie die Außenseite des Fläschchens ab.
- 7. Laden Sie die Methode ACD095AH, und führen Sie sie aus.
- 8. Stellen Sie das Fläschchen in die Halterung in der Probenkammer und schließen Sie die Probenkammertür.
- 9. Drücken Sie die Funktionstaste Blank (Blind), um den Blindwert zu messen.
- 10. Öffnen Sie den Probenkammerdeckel. Entnehmen Sie das Fläschchen aus der Halterung.
- 11. Fügen Sie den Inhalt eines <u>Phosphat-Rgt.- F10 Powder Pack</u> direkt aus der Folienpackung in das Fläschchen. Verwenden Sie einen Trichter, um das Reagenz hinzuzufügen.
- Verschließen Sie das Fläschchen fest mit der Kappe, und mischen Sie den Inhalt 10 bis 15 Sekunden durch Schwenken. Das Reagenz wird sich nicht vollständig auflösen. Wischen Sie die Außenseite des Fläschchens ab.
- 13. Warten Sie eine Reaktionszeit von 2 Minuten ab.
- 14. Stellen Sie das Fläschchen in die Halterung in der Probenkammer und schließen Sie die Probenkammertür.
- 15. Drücken Sie die Funktionstaste **Sample (Probe)**, um das Ergebnis in mg/l säurehydrolisierbarem Phosphat als Phosphor (P) anzuzeigen.

#### Hinweise:

- Während des gesamten Verfahrens sollten angemessene Sicherheitsvorkehrungen getroffen und eine gute Labortechnik angewandt werden.
- Ortho-Phosphationen reagieren mit dem Reagenz und erzeugen eine intensive blaue Färbung.
- Phosphate in organischer und kondensierter anorganischer Form (Meta-, Pyro- und Polyphosphate) müssen vor der Analyse in ortho-Phosphationen umgewandelt werden. Die Vorbehandlung der Probe mit Säure und Hitze schafft die erforderlichen Voraussetzungen für die Hydrolyse kondensierter anorganischer Formen. Organisch gebundene Phosphate werden durch Erhitzen mit Säure und Persulfat in ortho-Phosphationen umgewandelt. Die Menge organisch gebundener Phosphate kann folgendermaßen berechnet werden: mg/l Phosphat (organisch) = mg/l Phosphat (gesamt) - mg/l Phosphat (säurehydrolysierbar)
- Anwendung: für Wasser, Abwasser und Meerwasser.
- Stark gepufferte Proben oder Proben mit extremen pH-Werten sollten mit 1 mol/l Salzsäure oder 1 mol/l Natriumhydroxid (Natronlauge) auf einen pH-Wert zwischen 2 und 10 gebracht werden.
- ortho-Phosphat = reaktives Phosphor

| Störsubstanz                 | Störungsgrad            |
|------------------------------|-------------------------|
| Aluminium                    | mehr als 200 mg/l       |
| Arsenat                      | bei jeder Konzentration |
| Chrom                        | mehr als 100 mg/l       |
| Kupfer                       | mehr als 10 mg/l        |
| Eisen                        | mehr als 100 mg/l       |
| Nickel                       | mehr als 300 mg/l       |
| Kieselsäure (Siliziumdioxid) | mehr als 50 mg/l        |
| Silikat                      | mehr als 10 mg/l        |
| Sulfid                       | bei jeder Konzentration |
| Zink                         | mehr als 80 mg/l        |

• Störungen: Starke Trübungen können zu ungleichmäßigen Ergebnissen führen.

• Umrechnungen:  $mg/I PO_4 = mg/I P \times 3,07$  $mg/I P_2O_5 = mg/I P \times 2,29$ 

## AC2060 Siliziumdioxid-Tabletten-Test

#### Silicomolybdat-Methode

0,05 – 4 mg/l SiO<sub>2</sub>

- 1. Laden Sie die Methode AC2060, und führen Sie sie aus.
- Befüllen Sie eine saubere AQUAfast 24-mm-Rundküvette, Bestell-Nr. AC2V24, mit 10 ml der Probe. Verschließen Sie das Fläschchen fest mit dem Deckel. Wischen Sie die Außenseite des Fläschchens ab.
- 3. Stellen Sie das Fläschchen in die Halterung in der Probenkammer und schließen Sie die Probenkammertür.
- 4. Drücken Sie die Funktionstaste Blank (Blind), um den Blindwert zu messen.
- 5. Öffnen Sie die Probenkammertür, und nehmen Sie das Fläschchen aus der Halterung.
- 6. Geben Sie eine <u>Silica No. 1 Tablet</u> direkt aus dem Folienbeutel in das Fläschchen. Zerdrücken Sie die Tablette mit einem sauberen Rührstab.
- 7. Verschließen Sie das Fläschchen fest mit dem Deckel, und mischen Sie den Inhalt durch mehrmaliges Schwenken oder Umdrehen, bis sich die Tablette aufgelöst hat.
- 8. Warten Sie eine Reaktionszeit von <u>5 Minuten</u> ab.
- 9. Geben Sie eine <u>Silica No. 2 Tablet</u> direkt aus dem Folienbeutel in das Fläschchen. Zerdrücken Sie die Tablette mit einem sauberen Rührstab.
- Verschließen Sie das Fläschchen fest mit dem Deckel, und mischen Sie den Inhalt durch mehrmaliges Schwenken oder Umdrehen, bis sich die Tablette aufgelöst hat. Wischen Sie die Außenseite des Fläschchens ab.
- 11. Warten Sie eine Reaktionszeit von <u>1 Minute</u> ab.
- 12. Stellen Sie das Fläschchen in die Halterung in der Probenkammer und schließen Sie die Probenkammertür.
- 13. Tippen Sie auf die Funktionstaste Sample (Probe), um das Ergebnis in mg/l Siliziumdioxid anzuzeigen.

#### Hinweise:

Die Tabletten müssen in der richtigen Reihenfolge zugegeben werden.

## AC2061 Siliziumdioxid mit Phosphateliminierungs-Tabletten-Test

#### Silicomolybdat-Methode mit Phosphateliminierung

0,05 – 4 mg/l SiO<sub>2</sub>

- 1. Laden Sie die Methode AC2061, und führen Sie sie aus.
- Befüllen Sie eine saubere AQUAfast 24-mm-Rundküvette, Bestell-Nr. AC2V24, mit 10 ml der Probe. Verschließen Sie das Fläschchen fest mit dem Deckel. Wischen Sie die Außenseite des Fläschchens ab.
- Stellen Sie das Fläschchen in die Halterung in der Probenkammer und schließen Sie die Probenkammert
  ür.
- 4. Drücken Sie die Funktionstaste Blank (Blind), um den Blindwert zu messen.
- 5. Öffnen Sie die Probenkammertür, und nehmen Sie das Fläschchen aus der Halterung.
- Geben Sie eine <u>Silica No. 1 Tablet</u> direkt aus dem Folienbeutel in das Fläschchen. Zerdrücken Sie die Tablette mit einem sauberen Rührstab.
- 7. Verschließen Sie das Fläschchen fest mit dem Deckel, und mischen Sie den Inhalt durch mehrmaliges Schwenken oder Umdrehen, bis sich die Tablette aufgelöst hat.
- 8. Warten Sie eine Reaktionszeit von <u>5 Minuten</u> ab.
- Geben Sie eine <u>Silica PR Tablet</u> direkt aus dem Folienbeutel in das Fläschchen. Zerdrücken Sie die Tablette mit einem sauberen Rührstab.
- 10. Geben Sie eine <u>Silica No. 2 Tablet</u> direkt aus dem Folienbeutel in dasselbe Fläschchen. Zerdrücken Sie die Tablette mit einem sauberen Rührstab.
- Verschließen Sie das Fläschchen fest mit dem Deckel, und mischen Sie den Inhalt durch mehrmaliges Schwenken oder Umdrehen, bis sich die Tabletten aufgelöst haben. Wischen Sie die Außenseite des Fläschchens ab.
- 12. Warten Sie eine Reaktionszeit von <u>1 Minute</u> ab.
- 13. Stellen Sie das Fläschchen in die Halterung in der Probenkammer und schließen Sie die Probenkammertür.
- 14. Tippen Sie auf die Funktionstaste Sample (Probe), um das Ergebnis in mg/l Siliziumdioxid anzuzeigen.

- Die Tabletten müssen in der richtigen Reihenfolge zugegeben werden.
- Phosphationen stören unter den gegebenen Reaktionsbedingungen nicht.
- Wenn bekannt ist, dass kein Phosphat vorhanden ist, kann die Zugabe der Siliziumdioxid-PR-Tablette entfallen.

# AC4P60 Siliziumdioxid-Pulverpäckchen-Test

### Silicomolybdat-Methode

- 1 90 mg/l SiO<sub>2</sub>
- 1. Laden Sie die Methode AC4P60, und führen Sie sie aus.
- 2. Befüllen Sie eine saubere AQUAfast 24-mm-Rundküvette, Bestell-Nr. AC2V24, mit 10 ml der Probe. Die Probentemperatur sollte 15 bis 25 °C betragen. Verschließen Sie das Fläschchen fest mit dem Deckel. Wischen Sie die Außenseite des Fläschchens ab.
- 3. Stellen Sie das Fläschchen in die Halterung in der Probenkammer und schließen Sie die Probenkammertür.
- 4. Drücken Sie die Funktionstaste Blank (Blind), um den Blindwert zu messen.
- 5. Öffnen Sie die Probenkammertür, und nehmen Sie das Fläschchen aus der Halterung.
- Geben Sie den Inhalt eines <u>Silica HR Molybdate F10 Powder Pack</u> direkt aus der Folienpackung in das Fläschchen.
- 7. Verschließen Sie das Fläschchen fest mit dem Deckel, und mischen Sie den Inhalt durch mehrmaliges Schwenken oder Umdrehen.
- 8. Fügen Sie den Inhalt eines <u>Silica HR Acid Rgt. F10 Powder Pack</u> direkt aus der Folienpackung in das Fläschchen. Sind Siliziumdioxid oder Phosphat vorhanden, entsteht eine gelbe Farbe.
- 9. Verschließen Sie das Fläschchen fest mit dem Deckel, und mischen Sie den Inhalt durch mehrmaliges Schwenken oder Umdrehen.
- 10. Warten Sie eine Reaktionszeit von 10 Minuten ab.
- 11. Geben Sie den Inhalt eines <u>Silica Citric Acid F10 Powder Pack</u> direkt aus der Folienpackung in das Fläschchen. In diesem Schritt wird die durch Phosphat verursachte gelbe Farbe entfernt.
- 12. Verschließen Sie das Fläschchen fest mit dem Deckel, und mischen Sie den Inhalt durch mehrmaliges Schwenken oder Umdrehen. Wischen Sie die Außenseite des Fläschchens ab.
- 13. Warten Sie eine Reaktionszeit von <u>2 Minuten</u> ab.
- 14. Stellen Sie das Fläschchen in die Halterung in der Probenkammer und schließen Sie die Probenkammertür.
- Tippen Sie auf die Funktionstaste Sample (Probe), um das Ergebnis in mg/l Siliziumdioxid anzuzeigen.

- Gelegentlich enthalten Wasserproben Formen von Siliziumdioxid (Kieselsäure), die nur sehr langsam mit Molybdat reagieren. Die Beschaffenheit dieser Formen ist nicht bekannt.
- Eine Vorbehandlung mit Natriumhydrogencarbonat und anschließend mit Schwefelsäure macht diese Formen gegenüber Molybdat reaktiv. (Die Vorbehandlung ist in den "Standard Methods for the Examination of Water and Wastewater" [Einheitsverfahren zur Wasser- und Abwasseruntersuchung] unter "Silica Digestion with Sodium Bicarbonate" [Kieselsäureaufschluss mit Natriumbikarbonat] beschrieben.)

| Substanz | Störung  |
|----------|--|
| Eisen    | In großer Menge störend  |
| Phosphat | Keine Störung bei Konzentrationen von weniger als 50 mg/l PO <sub>4</sub><br>Bei 60 mg/l PO <sub>4</sub> beträgt der Störfaktor etwa 2 %.<br>Bei 75 mg/l PO <sub>4</sub> beträgt der Störfaktor etwa 11 %. |
| Sulfid   | Störung in allen Konzentrationen   |

# AC4P82 Sulfat-Pulverpäckchen-Test

#### Bariumsulfat/Trübungsmethode

- 5 100 mg/l SO<sub>4</sub>
- 1. Laden Sie die Methode AC4P82, und führen Sie sie aus.
- Befüllen Sie eine saubere AQUAfast 24-mm-Rundküvette, Bestell-Nr. AC2V24, mit 10 ml der Probe. Verschließen Sie das Fläschchen fest mit dem Deckel. Wischen Sie die Außenseite des Fläschchens ab.
- 3. Stellen Sie das Fläschchen in die Halterung in der Probenkammer und schließen Sie die Probenkammertür.
- 4. Drücken Sie die Funktionstaste Blank (Blind), um den Blindwert zu messen.
- 5. Öffnen Sie die Probenkammertür, und nehmen Sie das Fläschchen aus der Halterung.
- 6. Geben Sie den Inhalt eines <u>Sulpha 4/F10 Powder Pack</u> direkt aus der Folienpackung in das Fläschchen.
- 7. Verschließen Sie das Fläschchen fest mit dem Deckel, und mischen Sie den Inhalt durch mehrmaliges Schwenken oder Umdrehen. Wischen Sie die Außenseite des Fläschchens ab.
- 8. Warten Sie eine Reaktionszeit von <u>5 Minuten</u> ab.
- 9. Stellen Sie das Fläschchen in die Halterung in der Probenkammer und schließen Sie die Probenkammertür.
- 10. Tippen Sie auf die Funktionstaste Sample (Probe), um das Ergebnis in mg/l Sulfat anzuzeigen.

#### Hinweise:

• Sind Sulfationen vorhanden, wird die Lösung eingetrübt.

# AC2016 Sulfid-Tabletten-Test

### DPD-/Katalysator-Methode

0,04 – 0,5 mg/l S

- 1. Laden Sie die Methode AC2016, und führen Sie sie aus.
- Befüllen Sie eine saubere AQUAfast 24-mm-Rundküvette, Bestell-Nr. AC2V24, mit 10 ml der Probe. Verschließen Sie das Fläschchen fest mit dem Deckel. Wischen Sie die Außenseite des Fläschchens ab.
- 3. Stellen Sie das Fläschchen in die Halterung in der Probenkammer und schließen Sie die Probenkammertür.
- 4. Drücken Sie die Funktionstaste Blank (Blind), um den Blindwert zu messen.
- 5. Öffnen Sie die Probenkammertür, und nehmen Sie das Fläschchen aus der Halterung.
- 6. Geben Sie eine<u>Sulfide No. 1 Tablet</u> direkt aus dem Folienbeutel in das Fläschchen. Zerdrücken Sie die Tablette mit einem sauberen Rührstab.
- 7. Geben Sie eine <u>Sulfide No. 2 Tablet</u> direkt aus dem Folienbeutel in dasselbe Fläschchen. Zerdrücken Sie die Tablette mit einem sauberen Rührstab.
- Verschließen Sie das Fläschchen fest mit dem Deckel, und mischen Sie den Inhalt durch mehrmaliges Schwenken oder Umdrehen, bis sich die Tabletten aufgelöst haben. Wischen Sie die Außenseite des Fläschchens ab.
- 9. Warten Sie eine Reaktionszeit von 10 Minuten ab.
- 10. Stellen Sie das Fläschchen in die Halterung in der Probenkammer und schließen Sie die Probenkammertür.
- 11. Tippen Sie auf die Funktionstaste Sample (Probe), um das Ergebnis in mg/l Sulfid anzuzeigen.

- Die Tabletten müssen in der richtigen Reihenfolge zugegeben werden.
- Chlor und andere Oxidationsmittel, die mit DPD reagieren, beeinträchtigen den Test nicht.
- Um Sulfidverluste zu vermeiden, ist die Probe sorgfältig und mit minimaler Luftzufuhr zu entnehmen. Es ist wichtig, dass die Probe sofort nach der Entnahme untersucht wird.
- Die Probentemperatur sollte 20 °C betragen. Eine andere Temperatur kann zu höheren oder niedrigeren Ergebnissen führen.

# AC2065 Zinktabletten-Test

#### Zincon-Methode

0,02 – 1 mg/l Zn

- 1. Laden Sie die Methode AC2065, und führen Sie sie aus.
- Befüllen Sie eine saubere AQUAfast 24-mm-Rundküvette, Bestell-Nr. AC2V24, mit 10 ml deionisiertem Wasser. Verschließen Sie das Fläschchen fest mit dem Deckel. Wischen Sie die Außenseite des Fläschchens ab.
- 3. Stellen Sie das Fläschchen in die Halterung in der Probenkammer und schließen Sie die Probenkammertür.
- 4. Drücken Sie die Funktionstaste Blank (Blind), um den Blindwert zu messen.
- 5. Entleeren und trocken Sie das Fläschchen, und geben Sie dann 10 ml Probe in das Fläschchen.
- 6. Geben Sie eine <u>Copper / Zinc LR Tablet</u> direkt aus dem Folienbeutel in das Fläschchen. Zerdrücken Sie die Tablette mit einem sauberen Rührstab.
- Verschließen Sie das Fläschchen fest mit dem Deckel, und mischen Sie den Inhalt durch mehrmaliges Schwenken oder Umdrehen, bis sich die Tablette aufgelöst hat. Wischen Sie die Außenseite des Fläschchens ab.
- 8. Warten Sie eine Reaktionszeit von <u>5 Minuten</u> ab.
- 9. Stellen Sie das Fläschchen in die Halterung in der Probenkammer und schließen Sie die Probenkammertür.
- 10. Drücken Sie die Funktionstaste **Measure Rgnt Blank (Reagenzblindwert messen)**, um den Reagenzblindwert zu messen.
- 11. Öffnen Sie die Probenkammertür, und nehmen Sie das Fläschchen aus der Halterung.
- 12. Geben Sie eine <u>EDTA Tablet</u> direkt aus dem Folienbeutel in das Fläschchen. Zerdrücken Sie die Tablette mit einem sauberen Rührstab.
- Verschließen Sie das Fläschchen fest mit dem Deckel, und mischen Sie den Inhalt durch mehrmaliges Schwenken oder Umdrehen, bis sich die Tablette aufgelöst hat. Wischen Sie die Außenseite des Fläschchens ab.
- 14. Stellen Sie das Fläschchen in die Halterung in der Probenkammer und schließen Sie die Probenkammertür.
- 15. Tippen Sie auf die Funktionstaste Sample (Probe), um das Ergebnis in mg/l Zink anzuzeigen.

- Bei der Farbumkehrmethode wird ein Reagenz verwendet, dessen Farbintensität mit zunehmender Konzentration der zu messenden Spezies in den Proben abnimmt. Die Farbumkehrmethode erfordert eine Blindprobe und einen Reagenzblindwert. Die Blindprobe ist eine klare Lösung (deionisiertes Wasser) mit einer Absorption von Null. Der Reagenzblindwert ist ein Gemisch aus dem Reagenz und der Probe (ohne EDTA-Reagenz) und liefert den Nullkonzentrationspunkt mit dem dunkelsten Farbwert (höchste Absorption). Die Farbe der mit dem EDTA-Reagenz versetzten Proben nimmt mit steigender Konzentration für diese Methode ab.
- Die Messung des Reagenzblindwertes muss bei jeder Probenanalyse durchgeführt werden.
- Die Tabletten müssen in der richtigen Reihenfolge zugegeben werden.
- Bei einem hohen Restchlorgehalt ist die Analyse mit einer entchlorten Wasserprobe durchzuführen.

# Farbmessung aus "Application Log" (Anwendungsprotokoll) Nr. 131

#### Farbmessung – Wasserfarbe

Die Farbe unseres Wassers ist nicht nur für Trinkzwecke wichtig, sondern auch für die aquatische Ökosysteme sowie für private und industrielle Zwecke. Die Farbe im Wasser kann durch gelöste und suspendierte Stoffe hervorgerufen werden. Bei der Messung der Farbe von Wasser und Abwasser wird zwischen der scheinbaren Farbe und der echten Farbe unterschieden.<sup>1</sup> Die scheinbare Farbe ist die Farbe der Probe in der Form, wie sie erfasst wird, und umfasst die Farben, die auf die gelösten und suspendierten Stoffe im Wasser zurückzuführen sind. Die echte Farbe ist die resultierende Farbe der Probe, nachdem sie gefiltert wurde, um Einflüsse durch trübende Schwebstoffe wie Algen und Partikel auszublenden. Die echte Farbe entsteht nur durch im Wasser gelöste Stoffe, wie zum Beispiel natürliche organische Stoffe, Mineralien oder Chemikalien.

#### Methoden zur Messung der Wasserfarbe

Für die Messung der Wasserfarbe gibt es zwei Hauptansätze:

- Visuelle Verfahren Wasserproben werden visuell mit einer Reihe von Farbstandards verglichen.
- Spektrophotometrische Verfahren die Farbe wird durch die Messung bestimmt, wie viel Licht bei einer bestimmten Wellenlänge oder einer Reihe bestimmter Wellenlängen von einer Probe absorbiert oder transmittiert wird. Anschließend werden die Ergebnisse mit einem bekannten Farbstandard verglichen oder in die im Testprotokoll festgelegten Algorithmen eingespeist.

Methoden zur Farbmessung wurden von vielen wissenschaftlichen Gruppen veröffentlicht. Beispiele für spektrophotometrische Methoden sind Platin/Kobalt-Farbzahl, Tristimuluswerte und ADMI-Methoden.<sup>2</sup> Spezifische Methodenreferenzen für Wasserfarben: APHA-Standardmethoden 2120, EPA 110.1, DIN ISO 7887 und SAC 436 nm, China MEP GB 11903, NCASI Technical Bulletin 253 oder 803 und andere.

## Farbstandards und Farbeinheiten

#### Platin-Kobalt-Farbstandard und Maßeinheiten

Die gebräuchlichste Farbmessung für Wasser und Abwasser ist die Platin-Kobalt-Farbzahl (Pt-Co), auch bekannt als APHA- oder Hazen-Farbzahl. Diese Bezeichnungen werden häufig in unterschiedlichen Anwendungen verwendet, basieren aber auf identischen Verfahren. Die Messung nach dem Pt-Co/APHA/Hazen-Standard basiert auf der Hazen-Farbskala, die 1892 vom Chemiker Allen Hazen eingeführt wurde. Die Farbe ist eine hellgelbe bis braune Farbe. Der Bereich liegt zwischen 0 und 500 Platin-Farbeinheiten (PCU). Die Zwischenstandards werden aus einer 500 ppm Platin-Kobalt-Stammlösung hergestellt, die im Handel erhältlich ist<sup>3</sup> oder vom Benutzer selbst hergestellt werden kann<sup>1</sup>. Eine Farbeinheit ist die Farbe, die von 1 mg/l Platin in Form von Platinchlorid-Ionen erzeugt wird. Die durch den Vergleich mit den Platin-Kobalt-Standards gemessenen Farbwerte können je nach Verfahren als PCU-, Pt-Co-, APHA- oder Hazen-Einheiten angegeben werden. In den Vereinigten Staaten beträgt die maximal zulässige Farbzahl gemäß der *National Secondary Drinking Water Regulation* (nationale sekundäre Trinkwasserverordnung) 15 Farbeinheiten. Nach der Empfehlung der Weltgesundheitsorganisation (WHO) sollte die Farbe von Trinkwasser maximal 15 echte Farbeinheiten betragen.

#### Andere Farbstandards und Farbeinheiten

Für Proben, deren Farbe von Pt-Co-Standards abweicht, werden verschiedene Standards und Farbskalen verwendet. Dazu zählen beispielsweise: ADMI Color (American Dye Manufacturer's Institute), Gardner Color, Saybolt, Rosin, EBC (European Brewery Convention), CIE-System usw.

#### Platin-Kobalt-Farbskalen

In den Vereinigten Staaten werden in der Regel zwei Einzel-Wellenlängen-Methoden zur Farbmessung von Wasser und Abwasser verwendet. Die Farbeigenschaften ähneln denen der Pt-Co-Standards:

- Wasser/Abwasser bei 455 nm: Bei dieser Methode wird mit einem Spektrophotometer die Absorption des Lichts beim Durchgang durch die Probe bei einer Wellenlänge von 455 nm gemessen. Dieses Verfahren wird für Wasser mit natürlich vorkommenden Stoffen empfohlen, d. h. für pflanzliche Rückstände wie Blätter, Rinden, Wurzeln, Humus und Torfmaterialien. Trübe Proben sollten vor der Analyse mit einem 0,45-µm-Filter gefiltert werden, um die echte Farbe zu bestimmen. Zur Bestimmung der scheinbaren Farbe werden ungefilterte Proben gemessen. Samples (Proben) mit einer hohen Farbzahl (> 500 PCU) sollten verdünnt werden, bis die Farbe im Bereich der Standardkurve liegt. Der pH-Wert der Sample (Probe) sollte nicht angepasst werden, wenn er zwischen 4 und 10 liegt.
- Abwasser aus der Zellstoffindustrie bei 465 nm: Bei dieser Methode wird mit einem Spektrophotometer die Absorption des Lichts beim Durchgang durch die Probe bei einer Wellenlänge von 465 nm gemessen. Diese Methode ist an das NCASI Technical Bulletin 803 für Abwässer aus der Zellstoff- und Papierindustrie<sup>4</sup> (ersetzt das NCASI Technical Bulletin 253) angepasst. Die echte Farbe wird in einer Probe bestimmt, deren pH-Wert auf 7,6 +/- 0,05 pH eingestellt ist und die dann durch ein 0,8-µm-Filter geleitet wird. Die scheinbare Farbe wird anhand der ursprünglichen Probe bestimmt. Samples (Proben) mit starker Färbung sollten so verdünnt werden, dass sie in den Bereich der Standardkurve fallen.

Es werden auch Methoden mit mehreren Wellenlängen für die Farbmessung in Wasser und Abwasser empfohlen, deren Farbeigenschaften sich von denen der Platin-Kobalt-Standards unterscheiden, diese aber nicht ausschließen. Eine solche Methode ist SM 2120D-2001, die zur Berechnung von Tristimuluswerten und letztlich von Werten für dominante Wellenlänge, Farbton, Leuchtdichte und Reinheit einer Probe verwendet wird.<sup>2</sup> Bei dieser Methode prüft das Spektrophotometer eine Reihe von Punkten (z. B. 10 oder 30 Punkte) im Bereich von 400 bis 700 nm, um den Transmissionsgrad für jede Wellenlänge zu bestimmen, und verwendet die erhaltenen Werte zur Berechnung der Farbzahl nach der veröffentlichten Berechnungsmethode. Eine modifizierte Tristimulus-Methode, die in der Wasser- und Abwasserwirtschaft verwendet wird, basiert auf der Messung der prozentualen Transmission bei drei Wellenlängen (590, 540 und 438 nm).<sup>5</sup>

## Messbereich, Nachweisgrenze und Größe der Probenzelle

Genauigkeit, Präzision und Nachweisgrenzen, die mit einem bestimmten Farbmessverfahren erreichbar sind, sind von der Qualität der Geräteoptik und der Schichtdicke der Probenzelle abhängig.

- Spektrophotometer ermöglichen genauere und präzisere Ergebnisse als Kolorimeter, sind aber auch kostenaufwendiger. Niedrigste Nachweisgrenzen und eine höhere Genauigkeit bei niedrigen Konzentrationen lassen sich mit einem Spektrophotometer mit einer Probenzelle mit einer größeren Schichtdicke erreichen.
- Kolorimeter sind sinnvoll für die Durchführung von Tests außerhalb des Labors und zeigen die relative Farbintensität an. Die mit einem Kolorimeter ermittelten Farbwerte können von den mit einem Spektrophotometer ermittelten Farbwerten abweichen, insbesondere wenn die Wellenlänge des Kolorimeters nicht mit der Wellenlänge des Spektrophotometers übereinstimmt.

## Auswirkungen der Trübung auf die Farbmessung

Bei der Messung der "echten" Farbe muss die Trübung eliminiert werden, da sie Lichtstreuung verursachen und den Farbwert erhöhen kann. Die Trübung kann auch die Präzision der Messung beeinträchtigen. Diagramm 1 zeigt, dass die Filterung den Farbmesswert verringert – je trüber die Probe, desto deutlicher der Effekt. Abbildung 2 zeigt, dass die gefilterte Probe in der Regel eine bessere Präzision aufweist, insbesondere bei trüben Proben.



## Thermo Scientific Orion AquaMate Spektrophotometer für die Farbmessung

Die Orion AquaMate Spektrophotometer AQ8000 UV-Vis und AQ7000 Vis von Thermo Scientific verfügen über die erforderlichen Wellenlängen für dei Analyse jeder Farbmessung – durch vorprogrammierte Methoden, durch eine benutzerdefinierte Standardkurve oder auf Wunsch durch mehrere Wellenlängen. Die folgende Tabelle beschreibt einige der Farbmethoden, die mit den Spektrophotometern Orion AQ7000 und AQ8000 ausgeführt werden können.

| Farbmethode  | Vorprogrammierter<br>Name oder<br>benutzerdefiniert               | Probenzellen-<br>Größen               | Wellenlänge(n)  | Testprinzip  |
|--|---|---------------------------------------|---|--|
| Platin-Kobalt-<br>Farbmethode: Absorption<br>bei einer Wellenlänge<br>gemessen<br>Bereich: 0 – 500 PCU                     | CLRPT50<br>CLRPT25<br>CLRPT24<br>Benutzerdefinierte<br>Methode    | 50 mm<br>25 mm<br>24 mm<br>10 – 50 mm | 455 nm  | Wasser- und<br>Abwasserproben,<br>deren Farben<br>denen von Pt-Co-<br>Farbstandards<br>ähneln  |
| Anpassung des NCASI<br>253/803:<br>Absorptionsmessung bei<br>einer bestimmten<br>Wellenlänge.<br>Bereich: 0 – 500 PCU      | CLRPTP50<br>CLRPTP25<br>CLRPTP24<br>Benutzerdefinierte<br>Methode | 50 mm<br>25 mm<br>24 mm<br>10 – 50 mm | 465 nm  | Farbe in<br>Abwässern von<br>Zellstoffwerken   |
| ISO 7887; GB 11903:<br>Wasserqualität –<br>Untersuchung und<br>Bestimmung der Farbe  | Benutzerdefinierte<br>Methode                                     | 50 nm                                 | 436 nm<br>525 nm<br>620 nm                              | Roh- und<br>Trinkwasser,<br>Brauchwasser mit<br>geringer Färbung   |
| Tristimuluswert-Methode:<br>Der Transmissionsgrad<br>wird bei mehreren<br>Wellenlängen im Bereich<br>400 – 700 nm gemessen | Benutzerdefinierte<br>Methode                                     | 10 mm oder<br>50 mm                   | Mehrere<br>Wellenlängen<br>im Bereich 400<br>bis 700 nm | Wasser- und<br>Abwasserproben,<br>deren Farbe sich<br>von den Pt-Co-<br>Standards<br>unterscheidet,<br>diese aber nicht<br>ausschließt |

## Anwendungshinweise

Anwendungshinweise für die Farbprüfung erhalten Sie von unserer örtlichen technischen Vertriebsstelle, unserem technischen Kundendienst, unsere Webseite und andere Thermo Scientific-Webseiten. Diese Hinweise enthalten detaillierte Informationen über die Durchführung von Farbtests mit den Orion AquaMate Spektrophotometern von Thermo Scientific: AQ7000 sichtbarer Bereich und AQ8000 UV/Vis-Bereich. Folgende Anwendungshinweise sind verfügbar:

- Log 133, Farbe von Wasser und Abwasser nach der Pt-Co-Methode bei 455 nm
- Log 134, Farbzahl von Abwasser aus der Zellstoffindustrie nach der Pt-Co-Methode bei 465 nm (NCASI-Methode)

#### Referenzdokumente

- 1. Standard Methods for the Examination of Water and Wastewater, Method 2120B. www.standardmethods.org
- 2. Standard Methods for the Examination of Water and Wastewater, Method 2120. www.standardmethods.org
- 3. Available through Fisher Scientific at www.fishersci.com
- 4. NCASI, Technical Bulletin No. 253, Dec. 1971. Siehe 40CFR Abschnitt 136, Tabelle IB, Fußnote 18. <u>www.ecfr.gov</u>
- 5. EPA Color Method 110.1. http://www.umass.edu/tei/mwwp/acrobat/epa110.1colorspec.pdf

# UVA- und UV254-Farbmessungen aus "Application Log" (Anwendungsprotokoll) Nr. 137

#### UVA- und UV254-Farbmessung von Wasser

Bei dieser Methode wird mit einem Spektrophotometer die Absorption von Wasserproben, wie Trinkwasser und Quellwasser, bei einer Wellenlänge von 254 nm gemessen. Die Ergebnisse können mit organischem Kohlenstoff, Farbe und/oder Vorläufern von Desinfektions-Nebenprodukten korreliert werden. Die Ergebnisse können auch die Wirksamkeit von Aufbereitungsprozessen zur Entfernung von organischem Kohlenstoff anzeigen oder zusammen mit einem entsprechenden Ergebnis für den gesamten organischen Kohlenstoff (TOC) verwendet werden, um den spezifischen UV-Absorptionswert (SUVA) für eine Wasserprobe zu berechnen. Einzelheiten siehe Literaturhinweis 1.

## Referenzdokumente

- 1. EPA-Methode 415,3 Rev. 1.1. UV254 für SUVA. http://www.epa.gov/microbes/ordmeth.htm
- 2. Standard Methods 5910B, UV-Absorbing Organic Constituents. www.standardmethods.org
- 3. Benutzerhandbuch für Spektrophotometer Orion AquaMate 8100 UV-Vis

## **Empfohlene Ausstattung**

- Thermo Scientific Orion AquaMate AQ8000 Spektrophotometer
- Filtrationsgerät (z. B. Fisher XX1004707), 0,45-µm-Filter, 47 mm (z. B. Fisher HNWP04700)
- Vakuumquelle Sauger, Luftstrom oder Wasserstrom, handbetriebene oder elektrische Niederdruck-Vakuumpumpe
- UV-Einwegküvetten 1 cm (Fisher 14-377-009) oder Quarzprobenzellen
- Probenkarussell für 10-cm-Zelle, sofern erforderlich (Orion AQ100C)
- Orion pH-Meter und Elektrode

#### Lösungen

- 1. Deionisiertes Wasser ohne organische Bestandteile (DI).
- 2. Reagenzien für die pH-Einstellung. Natriumhydroxid, 0,1 N, Salzsäure, 0,1 N, Orion pH 4,01 und 7,00 Puffer (Orion 910104, Orion 910107).
- 3. Spektrophotometer-Kontrolllösung (SCS), optional: Organischer Kohlenstoff, KHP in pH-7-Phosphatpuffer, siehe Referenz 2 für die Zubereitung, oder kommerzieller SCS, PN 222-234700(www.unitylabservices.com).

## Lagerung und Reinigung von Probenzellen

Um reproduzierbare Ergebnisse zu erhalten, reinigen und lagern Sie die Probenzellen nach den Anweisungen im Benutzerhandbuch des Orion AquaMate-Geräts.

## Einrichtung des Messgeräts

Schalten Sie das Spektrophotometer ein. Wählen Sie eine geeignete Probenzellengröße und Methode für den zu erwartenden Gehalt an organischem Kohlenstoff in Ihren Proben. Beachten Sie dazu folgendes Diagramm. Wählen Sie die gewünschte Küvettenhalter-Position (1 oder 5 cm), indem Sie die Küvettenpositionstaste drücken. Für 10 cm installieren Sie ein Karussell mit 10 cm-Küvettenhalter. Rufen Sie über einen Computer den Inhalt des USB-Speichersticks auf. Kopieren Sie die gewünschte vorprogrammierte Methode aus dem Ordner "Orion" in den Ordner "Thermo" (siehe Hinweis 1). Nehmen Sie den USB-Speicherstick aus dem Computer, und stecken Sie ihn in den USB-Anschluss an der Vorderseite des AquaMate. Wählen Sie die Methode aus, und laden Sie sie. Drücken Sie die Funktion **Run Test (Test ausführen)**, um die Analyse zu starten. Weitere Informationen zur Auswahl der Methode entnehmen Sie bitte folgendem Diagramm.

| Probenkonzentration                    | Organischer<br>Kohlenstoff<br>> 0,5 mg/l | Organischer<br>Kohlenstoff<br>> 0,1 mg/l | Organischer<br>Kohlenstoff<br>> 0,05 mg/l |
|--|--|--|---|
| Quarzzellen-Größe<br>(Siehe Hinweis 2) | 1 cm (10 mm)                             | 5 cm (50 mm)                             | 10 cm (100 mm)                            |
| Methodenname                           | UV254_1                                  | UV254_5                                  | UV254_10                                  |

## Gerät auf Blindwert einstellen: Spektrophotometer-Test

- 1. Spülen Sie eine saubere Zelle dreimal mit VE-Wasser, wobei Sie nur die mattierten Seiten der Zelle berühren. Füllen Sie die Probenzelle mit DI-Wasser. Wischen Sie mit einem fusselfreien Tuch das Wasser von den Außenseiten ab.
- Öffnen Sie die Probenkammer und setzen Sie die mit DI-Wasser gefüllte Probenzelle (die Blindprobe) in den Probenhalter ein, wobei die klaren Seiten nach vorne und hinten zeigen müssen. Wenn sich die Probenzelle nicht im Lichtweg befindet, drücken Sie die Taste für die richtige Probenposition. Schließen Sie den Deckel, und tippen Sie auf die Funktionstaste Blank (Blind).
- 3. Testen Sie bei Bedarf ein SCS: Dieselbe Küvette mit dem vorbereiteten SCS entleeren und befüllen, abtrocknen und in den Probenhalter einsetzen. Schließen Sie den Deckel, und tippen Sie auf die Funktionstaste Sample (Probe). Notieren Sie das angezeigte Ergebnis.
- 4. Der Messwert für den SCS sollte innerhalb der gewünschten Kriterien gemäß Ihrem Qualitätssicherungsplan liegen. Siehe Abschnitt "Results" (Ergebnisse) für Beispiele.

## Probenlagerung

Die Proben werden nicht aufbewahrt. Analysieren Sie die Proben so schnell wie möglich nach der Entnahme. Proben können vor dem Test 48 Stunden lang bei < 6 °C gelagert werden. Siehe EPA-Methode 415.3 für die SUVA-Lagerung.

## Probenvorbereitung: pH-Einstellung und/oder Probenfiltration

Für Nicht-SUVA-Messungen: Sollte der pH-Wert nicht zwischen 4 und 10 liegen, stellen Sie den pH-Wert gemäß der Anleitung unter "pH Adjustment of Sample" (pH-Einstellung der Probe, S. 2) ein. Hinweis: Für eine SUVA-Bestimmung darf der pH-Wert nicht eingestellt werden.

Für UVA, UV254: Richten Sie das Filtrationsgerät mit einem 0,45-um-Filter ein. Spülen Sie den Filter mit 50 ml DI-Wasser, und verwerfen Sie das Spülwasser. Filtrieren Sie 50 ml der Probe. Messen Sie das Filtrat.

## Probenmessung

Vergewissern Sie sich, dass das Gerät ordnungsgemäß auf den Blindwert (auf "Null") eingestellt wurde. Berühren Sie nur die mattierten Seiten der Probenzelle, und spülen Sie die saubere Probenzelle mit einem Teil der gefilterten Probe ab. Füllen Sie dann die gefilterte Probe ein. Wischen Sie die Außenseiten trocken. Setzen Sie die Probenzelle in den Probenhalter ein, schließen Sie den Deckel, und tippen Sie auf die Taste **Sample (Probe)**. Notieren Sie das angezeigte Ergebnis. Die Ergebnisse können bei Bedarf auf dem USB-Stick gespeichert werden. Wenn der Absorptionsmesswert > 0,900 ist, verdünnen Sie die Probe, und messen sie erneut. Multiplizieren Sie den Messwert mit dem Verdünnungsfaktor. Wenn der Absorptionswert < 0,010 ist, sollte eine größere Probenzelle verwendet werden. Laden Sie die entsprechende Methode, und stellen Sie das Gerät auf den Blindwert ein.

# Qualitätskontrolle (QK)

Führen Sie eine SCS (Kontrolle) und Doppelproben mit jeder Charge durch, oder messen Sie QK-Proben gemäß Ihrem Qualitätssicherungsplan. Für SUVA-Tests sind die Anforderungen der EPA-Methode 415.3 zu befolgen. Siehe Referenz 1.

## pH-Einstellung der Probe

- 1. Hinweis: Verändern Sie nicht den pH-Wert von Proben, die für eine SUVA-Berechnung vorgesehen sind. Gehen Sie zum nächsten Schritt über.
- 2. Kalibrieren Sie die pH-Sonde in den Pufferlösungen pH 4,01 und 7,00.
- 3. Wärmen Sie die Probe auf Raumtemperatur auf.
- 4. Schütteln Sie die Probe, um eine Homogenität zu gewährleisten.
- 5. Messen Sie 50 ml der Probe mit einem Messzylinder in ein 100-ml-Becherglas ab.
- 6. Tauchen Sie die pH-Sonde in die Probe ein und notieren Sie den Anfangs-pH-Wert.
- Stellen Sie die Probe auf einen pH-Wert zwischen 4 und 10 ein, indem Sie tropfenweise 0,1 N Natriumhydroxid hinzufügen, um den pH-Wert zu erhöhen, oder 0,1 N Salzsäure, um den pH-Wert zu senken. Bei Bedarf kann eine anders starke Säure oder Base verwendet werden.
- 8. Beachten Sie, dass die Gesamtvolumenänderung nicht mehr als 1 % (0,5 ml) betragen sollte. Verwerfen Sie die Probe und setzen Sie sie mit einer stärkeren Säure oder Base neu an, wenn die Volumenänderung mehr als 1 % beträgt.
- 9. Notieren Sie den eingestellten pH-Wert. Fahren Sie mit dem nächsten Filtrationsschritt fort (S. 1).

# Ergebnisse der SCS-Tests mit dem Orion AQ8000-Spektrophotometer 25,0 mg/l organischer Kohlenstoff (KHP)

| Bias<br>Methode UV254_1                   | Erwartet<br>(nach SM 5910B) | Ergebnis<br>(AQ8000)   | Differenz                      | Bewertung |
|---|-----------------------------|------------------------|--------------------------------|-----------|
| Absorption                                | 0,358 cm <sup>-1</sup>      | 0,360 cm <sup>-1</sup> | 0,002 cm <sup>-1</sup> (0,6 %) | Gut       |
| Konzentration organischer<br>Kohlenstoffe | 25,0 mg/l                   | 24,9 mg/l              | 0,01 mg/l (0,4 %)              | Gut       |

**Verzerrung**: Die Messwerte eines KHP-Standards für organischen Kohlenstoff von 25,0 mg/l in Phosphatpuffer (hergestellt gemäß SM 5910B), der in einer 1-cm-Küvette getestet wurde, zeigen eine gute Genauigkeit:

- Das durchschnittliche AQ8000-Absorptionsergebnis liegt innerhalb von 0,002 Absorptionseinheiten des gemäß SM 5910B erwarteten Durchschnittswerts; 0,6 % Abweichung von der erwarteten Absorption.
- Das durchschnittliche Ergebnis der AQ8000-Konzentration an organischem Kohlenstoff (berechnet nach SM 5910B) liegt innerhalb von 0,1 mg/l des erwarteten Wertes und weist einen Unterschied von 0,4 % (99,6 % Wiederfindung) zum erwarteten Wert von 25,0 mg/l organischem Kohlenstoff auf.

| Präzision       | Zahl getesteter | Maximaler %RSD  | Ergebnis  | Bewertung |
|-----------------|-----------------|-----------------|-----------|-----------|
| Methode UV254_1 | Proben          | (nach SM 5910B) | (AQ8000)  |           |
| Absorption      | 14              | < 10,7 % RSD    | 0,3 % RSD | Gut       |

**Präzision**: Die Messwerte eines KHP-Standards für organischen Kohlenstoff von 25,0 mg/l in Phosphatpuffer (hergestellt gemäß SM 5910B), der in einer 1-cm-Küvette getestet wurde, zeigen eine gute Präzision:

 Die relative Standardabweichung (*relative standard deviation*, RSD) von 14 Testergebnissen des AQ8000 beträgt 0,3 % RSD und liegt damit deutlich unter dem nach SM 5910B erwarteten Höchstwert von 10,7 %.

- 1. Wenn sich die vorprogrammierte Methode nicht auf dem Memory Stick befindet, laden Sie die Methodendatei aus der Online-Bibliothek unter <u>www.thermoscientific.com/waterlibrary</u> herunter, oder kontaktieren Sie den technischen Kundendienst.
- 2. Alternativ können Sie für UV-Messungen entwickelte Einwegküvetten verwenden. Diese sind mit einer Küvettenlänge von 1 cm (10 mm) erhältlich.

6

# KAPITEL 6 Testmenü "Standard Curve" (Standardkurve)

# Konzentrationsmessungen mit der Quant Standardkurven-Anwendung (benutzerdefinierte Methode)

Der Standardkurventest wird verwendet, um eine quantitative Analyse mit einer Mehrpunkt-Kalibrierkurve über die Quant-Anwendung durchzuführen. Eine Kalibrierungskurve besteht aus Standards mit bekannter Konzentration. Die Konzentration der Proben wird mittels einer Kurvenangleichung gemessen.

Diese Testtechnik ist ideal, wenn kolorimetrische Reagenzien verwendet werden, für die der Testkit-Hersteller keinen Faktor oder keine Gleichung zur Ermittlung der Konzentration des Testkits angibt. Befolgen Sie bei der Erstellung einer benutzerdefinierten Methode stets die Reagenzienanweisungen des Testkit-Herstellers.

**Hinweis:** Wenn der Hersteller die Wellenlänge für das Testkit nicht angibt, sollte die Anwendung <u>Scan</u> mit einem vorbereiteten Standard zur Bestimmung der Wellenlängenparameter für das Reagenz vor der Erstellung einer Standardkurve ausgeführt werden.

## Das Standardkurven-Testverfahren eignet sich für folgende Zwecke:

- Erstellen einer Standardkurve durch Einrichten der Parameter und Messen der Standards für die Kurve
- Zum Anzeigen von Kalibrierkurvendaten
- Zum Bearbeiten einer Standardkurve, der Anzahl der Standards, Auswahl einer anderen Kurvenangleichung oder Löschen von Punkten aus der Kurve
- Messen von Proben unter Verwendung der Standardkurve und Anzeigen und Speichern von Messdaten
- Zum erneuten Aufrufen einer Standardkurven-Methode

# Aufrufen der Quant-Anwendung

Wischen Sie auf Bildschirm 1 nach links, um Bildschirm 2 aufzurufen. Die Anwendung Quant wird durch Antippen des entsprechenden Symbols ausgewählt. Die Anwendung Quant dient zur Entwicklung von Kalibrierkurven. Der Benutzer kann das Verfallsdatum, die Wellenlänge, die Referenzwellenlänge, die Gleichung und die Konzentrationseinheiten auswählen und die bekannten Standards eingeben, die für die Erstellung der Kurve verwendet werden sollen. Methoden können unter einem Namen gespeichert und später zur Messung einer Probenkonzentration verwendet werden



Standards hinzufügen und Kurventyp auswählen Der zur Anpassung der Standardmessdaten verwendete Gleichungstyp wird durch diese Gleichung festgelegt.

Auswahl, ob 1 oder 2 Wiederholungen erforderlich sind

Auswahl der Maßeinheiten, in denen die Probenmessungen angezeigt und gedruckt werden

| ÷         |          |       | 8                  | Speichern einer Methode |
|-----------|----------|-------|--------------------|-------------------------|
|           |          | tion  | Accessory<br>SETUP |                         |
|           |          |       |                    |                         |
| Ref λ     |          |       |                    |                         |
|           |          |       |                    |                         |
| Green Dy  | e Std# 1 | 1.000 |                    |                         |
| Green Dye | e Std# 2 | 2.000 |                    |                         |
| Green Dye | e Std# 3 | 4.000 |                    |                         |
| Green Dye | e Std# 4 | 5.000 |                    |                         |
| ⊕ Add S   | tandard  |       |                    |                         |
|           |          |       |                    |                         |
|           |          |       |                    |                         |

Sobald genügend eindeutige Standardkonzentrationswerte vorhanden sind, wird die Schaltfläche "Calibrate" (Kalibrieren) aktiviert.

Die Anzahl der eindeutigen Standardkonzentrationswerte wird anhand der Gleichung des Kurventyps bestimmt.



berechnetes r-Quadrat

|                               | Quant St         |       |  | 8                  |  |  |  |
|-------------------------------|------------------|-------|--|--------------------|--|--|--|
| Method Name<br>Green Dye Quan | titation         |       |  | Accessory<br>SETUP |  |  |  |
| A 62                          | <del>.</del> Y = |       |  |                    |  |  |  |
| Ref $\lambda$ –               |                  |       |  |                    |  |  |  |
|                               |                  |       |  |                    |  |  |  |
|                               |                  |       |  | •                  |  |  |  |
|                               |                  |       |  |                    |  |  |  |
|                               |                  |       |  |                    |  |  |  |
| Green Dye Std# 1              |                  | 1.000 |  | 0.257              |  |  |  |
| Green Dye Std# 2              |                  | 2.000 |  | 0.497              |  |  |  |
| Green Dye Std# 3              |                  | 4.000 |  | 0.880              |  |  |  |
| Green Dye Std∉ 4              |                  | 5.000 |  | 1.099              |  |  |  |
| 🕂 Add Standard                |                  |       |  |                    |  |  |  |
| Blank                         | Mea              | sure  |  | Sample             |  |  |  |
| J.m                           |                  |       |  |                    |  |  |  |
| Measure (Messung) wird nach d |                  |       |  |                    |  |  |  |



Kurventyp ändern, um eine neue Anpassung zu finden. Der R-Quadrat-Wert wird neu berechnet

Standards hinzufügen und Kurventyp auswählen

# Optionen für Standardkurve in Quant

# Einrichten der Parameter für eine Standardkurve

Markieren und ändern Sie die angezeigten Testparameter, einschließlich "Test Name" (Testname), "Wavelength" (Wellenlänge), "Curve Fit" (Kurvenanpassung), "Number of Standards" (Anzahl der Standards), und "Units" (Einheiten). Durch Antippen der blau markierten Parameter können Werte geändert werden. Wenn Sie zum Beispiel auf die rechts neben der Wellenlänge aufgeführte Gleichung tippen, erscheint eine Auswahl von Gleichungen. Wählen Sie die für Sie am besten geeignete Gleichung. Nach Abschluss der Mehrpunktkalibrierung werden die resultierende Gleichung und die Korrelationsergebnisse (r2) angezeigt. Diese Ergebnisse beruhen auf der Qualität des Blindwertes und der Genauigkeit der erstellten und eingegebenen Standards. Speichern Sie die Ergebnisse, indem Sie auf das blaue Disketten-Symbol oben rechts in der Ecke tippen.


## Erstellen einer Standardkurve in Quant

## Messstandards für eine Standardkurve

Sobald die Parameter für die Standardkurve festgelegt sind:

- 1. Bereiten Sie den Blindwert und die Standards vor, die für die Erstellung der Standardkurve verwendet werden sollen.
- Legen Sie die Blindprobe in die Probenkammer, schließen Sie den Deckel und tippen Sie auf die Funktionstaste Blank (Blind). Nachdem die Blindprobe gemessen wurde, nehmen Sie diese aus dem Probenkarussell.
- 3. Tippen Sie für jeden Standard auf das +, um einen Standardwert hinzuzufügen, und geben Sie die Konzentration des Standards ein.
- 4. Legen Sie die Blindprobe in die Probenkammer, schließen Sie den Deckel und tippen Sie auf die Funktionstaste Measure (Messen).
- 5. Wiederholen Sie den Vorgang für jeden Standard.
- 6. Wenn alle Standards gemessen wurden, wird die Absorption jedes Standards zusammen mit der Steigung, dem Achsenabschnitt und dem Korrelationskoeffizienten angezeigt, und es wird eine Grafik der Standardkurve angezeigt.
- 7. Geben Sie ggf. ein Verfalldatum in die Kurve ein.
- 8. Um den Test mit der Standardkurve zu speichern, tippen Sie auf das Speichern-Symbol oben rechts auf dem Bildschirm.
- 9. Alle blauen Felder können nach Bedarf ausgewählt und überarbeitet werden.
- 10. Tippen Sie auf die Funktion Sample (Probe), um die Messung der nachfolgenden Proben zu starten.

## **Probenmessung mit Quant**

Wenn Sie die Konzentration einer Probe mit einer beliebigen Methode oder Anwendung messen, wird die im Probennamen enthaltene Nummer jedes Mal um 1 erhöht, wenn Sie die Taste "Measure" (Messen) drücken. Siehe hierzu die nachfolgende Abbildung.



## Versuch beenden

Wenn Ihre Messungen abgeschlossen sind, klicken Sie auf das "X" in der oberen linken Ecke, um die Messung zu beenden. Sie müssen dann bestätigen, dass die Versuche beendet sind und die Datei gespeichert werden soll.

Die Daten können auch als Bericht ausgedruckt werden (wenn ein Netzwerkdrucker oder ein Banddrucker installiert ist) oder das Datum kann in einen Netzwerkordner oder auf ein USB-Laufwerk exportiert werden.



## **Daten exportieren**

Um die Daten exportieren zu können, müssen Sie einen Netzwerkpfad über Ethernet oder über WiFi eingerichtet haben. Eine typische Methode ist das direkte Speichern der Daten auf einem USB-Laufwerk. Der Bericht kann sowohl als CSV- als auch als JPEG-Datei gespeichert werden. Es folgt ein Beispiel für eine JPEG-Bilddatei.





## Bearbeiten einer Standardkurve

Wählen Sie auf der Startseite in Quant die aufgelistete Methodenkurve aus, die Sie überprüfen und bearbeiten möchten. Im folgenden Beispiel ist die Kurve "Chlorine Curve\_C" ausgewählt. Tippen Sie nach der Auswahl auf die Ellipse, um die Kurvenoptionen "Edit" (Bearbeiten), "Smart Method" (Intelligente Methode), "Duplicate" (Duplizieren), "Export" (Exportieren) oder "Delete" (Löschen) aufzurufen.



Sobald das Fenster mit den Kurvenoptionen erscheint, wählen Sie die Option "Edit" (Bearbeiten). Im Bearbeitungsmodus können Sie die Methode umbenennen, das Verfalldatum der Methode ändern oder eine andere Gleichung zur Darstellung der Kalibrierungskurve wählen. Tippen Sie auf die folgenden, geläufigeren blau hervorgehobenen Parameter, um sie zu überarbeiten und zu aktualisieren:

- "Method name" (Methodenname)
- "Calibration expiration" (Verfalldatum der Kalibrierung)
- "Calibration equation" (Kalibrierungsgleichung)

Wenn Sie Änderungen vornehmen, müssen Sie diese unter dem neuen Namen speichern. Standardpunkte können gelöscht oder hinzugefügt werden, allerdings muss die Kurve dann neu erstellt werden.

|                   | Quant Home   |                                 | $\leftarrow$   | Edit       | 6                  | $\leftarrow$         | Quant Stan   | dards              | 8                   |
|-------------------|--------------|---------------------------------|--|------------|--------------------|----------------------|--|--------------------|---------------------|
|                   |              | Name 🗸                          | Method name Chlorine Curve   |            | Accessory<br>SETUP | Method na<br>Chlorir | ame<br>ne Curve_C_2                                    |                    | A cessory<br>SET OF |
| 🗘 Quant 16-Jan-2  | 018_C_C_C    | 16-Jan-2018 01:33<br>PM         | Calibration expiration: N  |            |                    | Calibra              | tion expiration: Never                                 |                    | 1/30                |
| Quant 16-Jan-2    | 018_C_C      | 16-Jan-2018 01:33<br>PM         | λ 550  | Y = AX + B |                    | Y = -9.53            | 36X <sup>2</sup> + 9.936X + 1.275 r <sup>2</sup> = 1.0 |                    |                     |
| 🗘 Quant 05-Apr-20 | 019 MB3_C    | 05-Apr-2019 03:34<br>PM         | Reference λ         -           Y = 0.524X + 0.134         r² = 1.000  |            | 9                  | 4 -                  |  |                    | <b>~</b> •          |
|                   |              | 6994904919096720<br>99 <u>1</u> |  |            |                    | Absorbance           |  |                    |                     |
|                   |              |                                 | Choose Equation  |            |                    |                      |  |                    |                     |
| Chlorine Curve    |              | ×                               | Y = AX<br>Linear through zero  |            |                    | L                    |  | 0.4<br>tration     |                     |
|                   | ß            | Ē                               | Y = AX + B<br>Linear   |            | -                  | Standard<br>Didlik   | Conc<br>[m   | entration<br>g/mL] | Absorbance          |
| Edit              | Smart Method | Duplicate                       | $Y = AX^2 + BX$  |            |                    | Standar              | rd 1 0   | .120               | 2.330               |
|                   | •            |                                 | Quadratic through zero   |            |                    | Standar              | rd 2 0   | .350               | 3.585               |
|                   | Ш            |                                 | Y = AX <sup>2</sup> + BX + C<br>Quadratic                              |            |                    | Standar              | rd 3 0   | .600               | 3.804               |
|                   | Export       | Delete                          | Y = AX <sup>3</sup> + BX <sup>2</sup> + CX<br>Third order through zero |            |                    | + Add                | l standard   |                    |                     |
|                   |              |                                 | $Y = AX^3 + BX^2 + CX + D$<br>Third order                              |            |                    | Blank                | Measure  |                    | Sample              |
|                   | 0 0          | 1                               |  | 0          |                    |                      | Ø O  |                    |                     |

Sobald die Methode überarbeitet wurde, können die Probenmessungen fortgesetzt werden.



# KAPITEL 7 Menü für den Wellenlängen-Scan-Test

Mit der Anwendung "Scanning" (Scan) für Wellenlängen messen Sie die Absorption oder das prozentuale Transmissionsspektrum einer Probe. Mit Scans können Sie Bandenwellenlängen bestimmen und die Qualität eines Materials beurteilen.

Diese Testtechnik ist nützlich, wenn kolorimetrische Reagenzien verwendet werden, für die der Hersteller des Testkits keinen Wellenlängenparameter angibt. Befolgen Sie bei der Erstellung einer benutzerdefinierten Methode stets die Reagenzienanweisungen des Testkit-Herstellers.

**Hinweis:** Wenn der Hersteller keinen Faktor oder keine Gleichung zur Ermittlung der Konzentration des Testkits angibt, sollte das Verfahren <u>Standard Curve (Standardkurve)</u> mit vorbereiteten Standards verwendet werden, um den Faktor oder die Gleichung für das Reagenz zu bestimmen.

### Verwendungszweck der Scan-Tests:

- Scannen von Proben durch Einstellen von Parametern, Erstellen eines Basislinienscans und Scannen einer Probe
- Anzeige und Bearbeitung von Scandaten, einschließlich grafischer Daten, Bestimmung der Peakhöhe mithilfe einer 3-Punkt-Netzgleichung, Berechnung der Fläche unter einer Kurve und Kennzeichnung von Peaks und Kurventiefpunkte
- Aufrufen bestehender Scan-Methoden

## Einrichten der Parameter für einen Scan

SCAN ist ein Multi-Wellenlängen-Scan, der die Absorption (ABS) oder die prozentuale Transmission (%T) über einen justierbaren Wellenlängenbereich und eine justierbare Geschwindigkeit und Intervallauflösung auswählt. Der Zweck der Funktion "SCAN" besteht darin, die Absorptions- oder Transmissionseigenschaften einer Probe oder einer mit Reagenz versetzten Probe zu bewerten. Dies ist nützlich, um die optimale Wellenlänge für die Einführung einer neuen Methode zu bestimmen.

| Hethod Name |     | New           |       | Scanbereich:<br>Zwischen 190 nm und 1.100 nm für UV-VIS   |
|-------------|-----|---------------|-------|---|
| GD1         | ARS | 3/30<br>%T    | SETUP | Zwischen 325 nm und 1.100 nm für Vis-<br>Instrumente  |
| Range:      | 325 | - 1100        | nm    | Intervall:  |
| Interval    | 2   |               | nm    | Legt fest, wie häufig das Gerät eine<br>Messung erfasst.  |
| <u></u>     |     | •<br>Continue |       | In dieser Abbildung werden die Daten alle<br>2 nm erfasst.  |
|             |     | Continue      |       | Die Scan-Geschwindigkeiten "Fast" (Schnell) "Medium"<br>(Mittel) und "Slow" (Langsam) begrenzen die Anzahl der<br>angebotenen Optionen für das Datenerfassungsintervall |
|             |     |               | Sneed | Interval Ontions  |

| Speed  | Interval Options                  |
|--------|-----------------------------------|
| Fast   | 5 nm, 2 nm                        |
| Medium | 5 nm, 2 nm, 1 nm                  |
| Slow   | 5 nm, 2 nm, 1 nm, 0.5 nm, 0.2 nm, |

Tippen Sie auf die hervorgehobenen Felder und bearbeiten Sie diese:

- Überarbeiten Sie den Methodennamen für diesen Scan.
- Wählen Sie zwischen ABS oder %T für die y-Achse
- Stellen Sie den Wellenlängenbereich ein, der für die Anzeige auf der x-Achse verwendet werden soll
- Geben Sie das Intervall und die Geschwindigkeit des Scans an; beachten Sie, dass bei langsameren Scans eine höhere Auflösung der Scanintervalle verfügbar ist.
- Tippen Sie auf das hervorgehobene **Speichern-Symbol**, um die Scanmethode zu speichern

## Erfassen eines Basislinienscans und Scannen einer Probe

### Hinweis: Bitte stellen Sie sicher, dass Sie den gleichen Probengefäßhalter und die gleiche Gefäßgröße verwenden.

Sobald die Parameter für den Scan eingestellt sind, können Sie fortfahren.

- 1. Drücken Sie "Continue" (Weiter), um einen Scan-Test durchzuführen
- 2. Installieren Sie eine Blindprobe und stellen Sie den Blindwert für das Systems ein.
- 3. Installieren Sie die gewünschte Probe und tippen Sie auf **Measure (Messen)**. Dieses Verfahren kann für mehrere Proben wiederholt werden.
- 4. Nach Abschluss des Scanvorgangs können bis zu drei (3) Referenzwellenlängen ausgewählt werden, um die Ergebnisse speziell für diese Wellenlängen anzuzeigen. Im folgenden Beispiel wurden 622, 663 und 700 nm gewählt.
- 5. Tippen Sie entweder auf die Wellenlängennummer, um eine neue spezifische Referenzwellenlänge einzugeben, oder ziehen Sie die vertikale Wellenlängenlinie auf einen neuen Referenzwert.
- 6. Wenn zusätzliche Funktionen für einen Probenscan von Interesse sind, tippen Sie auf die Ellipse rechts neben der Probe, um weitere Optionen zu öffnen

Sobald der Scanvorgang abgeschlossen ist, folgen Sie den Schritten zum Speichern und Exportieren. Die grafischen Daten können mit einer Zwei-Finger-Stretch- oder Kompressions-Touchscreen-Methode manipuliert werden.



## Berechnungen mit den Scan-Daten durchführen

Tippen Sie nach Abschluss eines Scans auf das Ellipsen-Symbol, um die Optionen anzuzeigen:

- Löschen Sie die gescannte Probe
- Wählen Sie einen Peak
- Exportieren Sie die Daten

In der nachstehenden Abbildung wurde die Option "Peak Pick" (Peak auswählen) ausgewählt, und es wurden automatisch zwei Werte (654,233 und 663,910 nm) angegeben.

Mit der Stretch- und Zoomfunktion des Touchscreens kann der Peak genauer betrachtet werden, und nach der endgültigen Auswahl würde es so aussehen, als ob Peak 2 bei 663,910 (~664 nm) ausgewählt würde. Bitte beachten Sie, dass die Glättung entweder ein- oder ausgeschaltet werden kann.



Weitere verfügbare Funktionen:

- Auswahl von Peaks mithilfe von
  - Peaks only (Nur Peaks)
  - Valleys only (Nur Tiefpunkte)
  - o Peaks und Tiefpunkte
- Auswahl der maximalen Anzahl von Peaks (1 bis 20 sind zulässig)
- Auswahl einer Glättungsfunktion, die die Anzahl der Peaks reduziert.



Auswahl

## Die Funktion "3Pt Net" (3-Pkt. Netz)

Durch Umschalten der Funktionstaste **3Pt Net (3-Pkt. Netz)** wird die Peakhöhe unter Verwendung einer 3-Punkt-Netz-Gleichung aktiviert und der 3Pt Netto-Wert erscheint nun neben den drei Wellenlängen. Der Wert für die 3-Punkt-Netz-Absorption wird berechnet und mit dem Faktor 1.000 multipliziert.



## Funktion zur Flächenberechnung

Um die Fläche unter der Kurve zu berechnen, müssen nur zwei Wellenlängen angegeben werden. Wenn Sie die Funktion "3Pt Net" (3-Punkt-Netz) ausschalten, werden automatisch die Wellenlängen 1 und 2 ausgewählt, und die Fläche unter der Kurve kann berechnet werden, wenn Sie diese Funktion aktivieren.





## Aufrufen einer bestehenden Scan-Methode

Um eine bestehende Scan-Methode aufzurufen, rufen Sie Screen 1 (Bildschirm 1) auf und tippen auf das Scan-Symbol. Es wird eine Liste der erstellten Scan-Methoden angezeigt. Wählen Sie eine Option aus der Liste, tippen Sie darauf und führen Sie den Scan durch.





# **KAPITEL 8** Integrierte AquaMate-Software

## Tests und Berichte zur Leistungsüberprüfung

AquaMate-Geräte sind mit Standardtests zur Leistungsüberprüfung vorkonfiguriert. Diese Tests können nicht gelöscht oder bearbeitet werden.

### Tests zur Leistungsüberprüfung ausführen

Die Durchführung eines Tests zur Leistungsüberprüfung erfolgt ähnlich wie bei jeder anderen Versuchsmethode. Folgen Sie den Bildschirmanweisungen.

### Benutzerdefinierte Tests zur Leistungsüberprüfung

Einige Tests zur Leistungsüberprüfung können an benutzerspezifische Testanforderungen angepasst werden. Um einen Test anzupassen, muss er zunächst dupliziert werden. Häufig erfordern Standardarbeitsanweisungen eine regelmäßige Leistungsüberprüfung der Geräte. Für Tests wie Streulicht, Wellenlängengenauigkeit und photometrische Genauigkeit sind spezielle Filter und Standards erforderlich. Die Standardtests zur Leistungsüberprüfung sind für bestimmte Normen und Filter ausgelegt (siehe Tabelle 1). Diese Tests können jedoch dupliziert und an die gewünschten Standards angepasst werden.

| Tests zur Leistungsüberprüfung                                | Beschreibung  | Duplizieren<br>möglich? |
|---|---|-------------------------|
| Wavelength Accuracy Xenon<br>(Wellenlängen-Genauigkeit Xenon) | Mit diesem Test wird die Leistung der Wellenlängenachse des<br>Spektrophotometer überprüft.   | Nein                    |
|   | Scannt und lokalisiert bekannte Xenon-Emissionsspitzen und überprüft, ob sie innerhalb der Spezifikation für das Gerät genau gefunden werden.   |                         |
| Drift at 500 nm (Drift bei 500 nm)                            | Misst die Absorption bei 500 nm in 1 Minuten-Abständen über eine<br>Stunde. Measured at 0A (Messung bei 0 A) (offener Strahl)<br>Blindwerteinstellung und Meldung der maximalen Abweichung von<br>Null. Vergleicht das Ergebnis mit der Spezifikation für das Gerät.<br>Das Ergebnis sollte geringer als die Spezifikation ausfallen. | Nein                    |
| Noise 0A at 500nm<br>(Rauschen 0 A bei 500 nm)                | Zeichnet 60 ABS-Messungen im 1 Sekunden-Abstand auf. Gibt den<br>RMS-Wert (RMS = root mean square [quadratischer Mittelwert]) des   | Nein                    |
| Noise 1.0A at 500 nm<br>(Rauschen 1,0 A bei 500 nm)           | Datensatzes als Rauschwert zurück und vergleicht den Wert mit der Spezifikation für das Gerät.  |                         |
| Noise 2.0A at 500 nm<br>(Rauschen 0,0 A bei 500 nm)           | Setzen Sie für diese Messungen einen Glasfilter mit einer nominalen Absorption von 1 A oder 2 A neutraler Dichte ein.   |                         |
|   | Das Ergebnis sollte geringer als die Spezifikation ausfallen.   |                         |
|   | Hinweis: Das Rauschen fällt bei den AquaMate-Geräten so gering<br>aus, dass das Ergebnis mit mehr als den üblichen 3 Dezimalstellen<br>angezeigt wird.  |                         |

Tabelle 1. Beschreibung der einzelnen Tests und Angabe, ob die Möglichkeit zur Duplizierung besteht

| Leistungsüberprüfungstest  | Beschreibung   | Duplizieren<br>möglich? |
|--|--|-------------------------|
| Baseline Flatness 1000 to 200 nm<br>(Basislinien-Flachheit 1.000 bis 200 nm) | Misst jede systematische Abweichung vom perfekten<br>Nullpunkt beim Scannen über den gemeinsamen<br>Wellenlängenbereich. Die Daten werden geglättet, um den<br>Rauscheffekt zu eliminieren (Rauschen kann separat<br>gemessen werden). Das Ergebnis ist die maximale<br>Abweichung von Null und wird mit der Spezifikation für das<br>Gerät verglichen.  | Nein                    |
|  | Das Ergebnis sollte geringer als die Spezifikation ausfallen.  |                         |
| Stray Light SRE 220 filter<br>(Streulichtfilter SRE 220) (UV-Vis-Modelle)    | Misst Streulicht bei der angegebenen Wellenlänge.  | Ja                      |
| Stray Light SRE 400 Filter<br>(Streulichtfilter SRE 400) (Vis-Modelle)       | Der Filter ist ein Langpassfilter, der etwas oberhalb der<br>Testwellenlänge ansetzt. Bei der Testwellenlänge sollte es<br>völlig dunkel sein, d. h. 0 % T. Längere Wellenlängen<br>durchdringen den Filter, sodass jeder bei 220 nm gemessene<br>Transmissionsgrad in Wirklichkeit Photonen längerer<br>Wellenlängen sind, die "Streulicht" darstellen. Zu den<br>Streulichtquellen gehören Effekte zweiter Ordnung,<br>Unvollkommenheiten im Gitter sowie Unvollkommenheiten<br>oder Verschmutzungen auf Spiegeln. |                         |
|  | Der gemessene Transmissionsgrad wird mit der Spezifikation für das Gerät verglichen.   |                         |
|  | Der Test misst die Durchlässigkeit bei der angegebenen<br>Wellenlänge.   |                         |
|  | Das Ergebnis sollte geringer als die Spezifikation ausfallen.  |                         |
| Wavelength Accuracy<br>(Wellenlängen-Genauigkeit)                            | Mit diesem Test wird die Leistung der Wellenlängenachse des Spektrophotometer überprüft.   | Ja                      |
|  | Verwenden Sie diesen Test mit einem kalibrierten<br>Wellenlängenfilter, z. B. einem Holmium- oder<br>Didymiumglasfilter.   |                         |
|  | Der Benutzer gibt die Spitzenwellenlängen und die<br>Kalibrierungsunsicherheit aus dem Kalibrierungszertifikat ein.  |                         |
|  | Das Gerät scannt den entsprechenden Wellenlängenbereich und lokalisiert die Mitte des Peaks.   |                         |
|  | Die angegebene Wellenlänge sollte mit der im Zertifikat<br>angegebenen Wellenlänge innerhalb der Summe aus<br>Gerätespezifikation und Kalibrierungsunsicherheit<br>übereinstimmen.   |                         |

| Tests zur Leistungsüberprüfung                       | Beschreibung  | Duplizieren<br>möglich? |
|--|---|-------------------------|
| Photometric Accuracy<br>(Photometrische Genauigkeit) | Mit diesem Test wird die photometrische (Absorptions-)Leistung des Spektrophotometers überprüft.  | Ja                      |
|  | Verwenden Sie diesen Test mit einem oder mehreren kalibrierten<br>Absorptionsfiltern, wie sie im SPECTRONIC Standards 2 Kit<br>enthalten sind, um die Genauigkeit im sichtbaren Bereich zu prüfen.<br>Verwenden Sie kalibrierte Kaliumdichromat-Lösungszellen, Metall-<br>auf-Quarz-Filter oder andere anerkannte kalibrierte<br>Standardmaterialien<br>für den UV-Bereich. <sup>a</sup> Es können mehrere Filter,<br>die für dieselben Wellenlängen, aber unterschiedliche<br>Absorptionswerte kalibriert sind, in einem einzigen Test konfiguriert<br>werden. |                         |
|  | <ul> <li>Der Anwender gibt</li> <li>die Kalibrierungswellenlängen,</li> <li>die Absorptionswerte des Zertifikats</li> <li>und die Kalibrierungsunsicherheitswerte</li> </ul>  |                         |
|  | aus dem Kalibrierungszertifikat ein. Der Anwender gibt auch die<br>Leistungsspezifikation für den zu prüfenden Absorptionswert aus<br>dem technischen Datenblatt des Geräts ein. <sup>b</sup>   |                         |
|  | Das Gerät misst und protokolliert die Absorption bei jeder<br>angegebenen Wellenlänge. Es wird eine Integrationszeit von 1 s<br>verwendet.  |                         |
|  | Die angegebene Absorption sollte mit der im Zertifikat angegebenen<br>Absorption innerhalb der Summe aus Gerätespezifikation und<br>Kalibrierungsunsicherheit übereinstimmen.   |                         |

<sup>&</sup>lt;sup>a</sup> Wir empfehlen die Verwendung von Didymiumglas-Doppelstandard-Filtern, die sowohl für die Wellenlängen-Peaks als auch für die photometrische Genauigkeit kalibriert sind. Einige Kunden berichteten Probleme bei der Reproduktion von Kalibrierungswerten für diese Art von Standard, obwohl ihre Geräte die Kalibrierungswerte für andere, weithin akzeptierte und anerkannte Standards reproduzieren können. Kunden, die sich an den Kundendienst wenden, weil ein Gerät eine photometrische Genauigkeitsprüfung mit einem photometrischen Didymium-Standard nicht bestanden hat, müssen möglicherweise die photometrische Genauigkeit mit einem anderen Standard überprüfen, bevor eine Rücksendung im Rahmen der Gewährleistung genehmigt wird.

<sup>&</sup>lt;sup>b</sup> Zukünftige Softwareversionen enthalten möglicherweise eine Bezugstabelle, deren Daten auf der Grundlage des vom Benutzer eingegebenen Absorptionswerts des Zertifikats in die Gerätespezifikation eingesetzt werden.

## **Benutzerdefinierter Streulicht-Test**

|                    | PV Home   |                         |
|--------------------|-----------|-------------------------|
|                    |           |                         |
|                    |           | 01-Feb-2018<br>08:37 AM |
|                    |           | 01-Feb 2018<br>08:37 AM |
|                    |           | 01.÷eb-2018<br>08:37 AM |
|                    |           | 01-Feb-2018<br>08:37 AM |
|                    |           | 01-Feb-2018<br>08:37 AM |
|                    |           | 01.Fob.2019             |
| Stray Light SRE 22 | 20 Filter | :                       |
|                    |           | Target: %T              |
| SRE 220 Filter     | 220       | < 0.05                  |
|                    |           |                         |
|                    |           |                         |
|                    |           |                         |
|                    |           |                         |
|                    |           |                         |
|                    |           |                         |





## Benutzerdefinierter Wellenlängen-Genauigkeitstest



 ←
 Edit

 Method name

 Wavelength Accuracy-Holmium

 27/00

 Wavelength Accuracy

 Wavelength Repeatability test

 Measured in Transmittance

 ID

 Holmium Glass Filter

 10

 287.6

 1.0

 333.9

 1.0

 637.7

 1.0

 637.7

 1.0

 637.7

 1.0

 Blank

Umschalten, um einen Test der Wellenlängen-Reproduzierbarkeit durchzuführen

|                         | PV Home   | ◙                       |
|-------------------------|-----------|-------------------------|
|                         |           |                         |
| Wavelength Accuracy     | Xenon     | 01-Feb-2018<br>10:21 AM |
| Drift at 500nm          |           | 01-Feb-2018<br>10:21 AM |
| Noise 0.0A at 500nm     |           | 01-Feb-2018<br>10:21 AM |
| Noise 1.0A at 500nm     |           | 01-Feb-2018<br>10:21 AM |
| Noise 2.0A at 500nm     |           | 01-Feb-2018<br>10:21 AM |
| Baseline Flatness 1000  | 0-200 nm  | 01-Feb-2018<br>10:21 AM |
| Stray Light SRE 220 Fil | ter       | 01-Feb-2018<br>10:21 AM |
| Wavelength Accuracy     |           | 01-Feb-2018<br>10:21 AM |
| Photometric Accuracy    |           | 01-Feb-2018<br>10:21 AM |
| Custom Stray Light Te   | st        | 01-Feb-2018<br>08:37 AM |
| Wavelength Accuracy     | - Holmium | 10 195-2010<br>10 07-00 |

Umschalten zur Messung im Transmissionsmodus

## Benutzerdefinierter photometrischer Genauigkeitstest

Ein typischer Standardkit für die Prüfung der photometrischen Genauigkeit wird mit einem Kalibrierungszertifikat geliefert. In diesem Abschnitt wird erläutert, wie ein photometrischer Genauigkeitstest für benutzerdefinierte Standardkits eingerichtet wird.

## Thermo Fisher

5225 Verona Road, Bldg.1 Madison, WI 53711 USA

www.thermo.com

#### Certificate of Calibration SPECTRONIC Standards 2 Kit 840-253100

Submitted to: THERMO FISHER SCIENTIFIC 5225 VERONA ROAD MADISON, WI 53711 USA Serial Number: SA1234 Certificate Number: CC001234 Date of Calibration: 20 Apr 2015 Performed by: John Doe Test Method: 397-018500 Rev A Bench Used: MSN000 123456 Sample Temperature: 23 ± 1 °C

Certified Percent Transmittance Values and Uncertainties

| Standard ID | SA0706 -1 | SA0706 -2                                 | SA0706 -3 | SA0706 -4  |
|-------------|-----------|---|-----------|------------|
| Nominal %T  | 50 %T     | 30 %T                                     | 10 %T     | 3 %T       |
| Uncertainty | ± 0.30 %T | ± 0.18 %T                                 | ± .071 %T | ± 0.048 %T |
| 440.0 nm    | 44.51 %T  | 26.73 %T                                  | 7.78 %T   | 2.14 %T    |
| 465.0 nm    | 50.83 %T  | 33.09 %T                                  | 9.71 %T   | 2.99 %T    |
| 546.1 nm    | 51.57 %T  | 34.06 %T                                  | 9.43 %T   | 2.87 %T    |
| 590.0 nm    | 49.57 %T  | 31.26 %T                                  | 8.44 %T   | 2.42 %T    |
| 635.0 nm    | 49.78 %T  | 30.83 %T                                  | 10.02 %T  | 3.14 %T    |
|             |           | 5. A. |           |            |

| Standard ID | SA0706 -1  | SA0706 -2  | SA0706 -3  | SA0706 -4  |
|-------------|------------|------------|------------|------------|
| Nominal Abs | 0.3 A      | 0.5 A      | 1.0 A      | 1.5 A      |
| Uncertainty | ± 0.0026 A | ± 0.0026 A | ± 0.0031 A | ± 0.0070 A |
| 440.0 nm    | 0.3516 A   | 0.5730 A   | 1.1089 A   | 1.6704 A   |
| 465.0 nm    | 0.2939 A   | 0.4803 A   | 1.0127 A   | 1.5237 A   |
| 546.1 nm    | 0.2876 A   | 0.4677 A   | 1.0254 A   | 1.5419 A   |
| 590.0 nm    | 0.3048 A   | 0.5051 A   | 1.0736 A   | 1.6165 A   |
| 635.0 nm    | 0.3029 A   | 0.5110 A   | 0.9992 A   | 1.5030 A   |

8

| Thermo Fish   |                     |                                   |  |            |                 | ×                    |             |
|---|---------------------|-----------------------------------|--|------------|-----------------|----------------------|-------------|
| 5225 Verona Road, E                                       | lida.1              |                                   |  |            |                 | Method Name          |             |
| Madison, WI 53711 U                                       | ISA                 |                                   |  |            |                 | Custom Photomet      | ric Accurac |
|   |                     |                                   |  |            |                 |                      |             |
| www.thermo.com  | Cert                | ificate of Ca<br>ONIC Standards 2 | libration<br>Kit 840-253100                                      |            |                 | Photometric Accurac  | y           |
| Submitted to:   |                     | Serial N                          | lumber: SA1234   |            |                 | Certificate #: CLEV3 | 00 999003   |
| THERMO FISHER SO<br>5225 VERONA ROAL<br>MADISON, WI 53711 |                     | Certific<br>Date of<br>Perform    | ate Number: CC0012<br>Calibration: 20 Apr 20<br>and hv: John Doe | 34<br>D15  |                 | SA0706-1 SA0706-2    | Ð           |
| JSA   |                     | Test Me                           | thod: 397-018500 Re  | v A        |                 |                      |             |
|   |                     | Sample                            | Temperature: 00 _ 1  | 3          |                 | Standard ID          |             |
| Dealifier d Dearant Trans                                 |                     |                                   |  |            |                 | Nominal(Abs)         |             |
| Standard ID   | SA0706 -1           | d Uncertainties                   | SA0706 -3  | SA0706 -4  | 1               |                      |             |
| Nominal %T  | 50 %T               | 30 %T                             | 10 %T  | 3 %T       |                 | Uncertainty(Abs)     | 0.0         |
| Uncertainty   | ± 0.30 %T           | ± 0.18 %T                         | ±.071 %T   | ± 0.048 %T |                 |                      |             |
| 440.0 nm  | 44.51 %T            | 26.73 %T                          | 7.78 %T  | 2.14 %T    |                 | Specification(Abs)   | 0.0         |
| 465.0 nm  | 50.83 %T            | 33.09 %T                          | 9.71 %T  | 2.99 %T    |                 | λ (nm)               |             |
| 546.1 nm  | 51.57 %T            | 34.06 %T                          | 9.43 %T  | 2.87 %T    |                 |                      |             |
| 590.0 nm  | 49.57 %T            | 31.26 %T                          | 8.44 %T  | 2.42 %T    | Real Providence | <u>44</u> 0.0        |             |
| 635.0 nm  | 49.78 %T            | 30.83 %T                          | 10.02 %T   | 3.14 %T    |                 |                      |             |
| Certified Absorbance                                      | Values and Uncertai | nties                             |  |            |                 | 465.0                |             |
| Standard ID   | SA0706 -1           | SA0706 -2                         | SA0706 -3  | SA0706 -4  | 1               | 500.0                |             |
| Nominal Abs   | 0.3 A               | 0.5.4                             | 10.4   | 15.6       | 1               | ×25.0                |             |
| Uncertainty   | ± 0.0026 A          | ± 0.0026 A                        | ± 0.0031 A   | ± 0.0070 A |                 | 635.0                |             |
| 440.0 nm  | 0.3516 A            | 0.5730 A                          | 1.1089 A   | 1.6704 A   |                 |                      |             |
| 465.0 nm  | 0.2939 A            | 0.4803 A                          | 1.0127 A   | 1.5237 A   |                 |                      |             |
| 546.1 nm  | 0.2876 A            | 0.4677 A                          | 1.0254 A   | 1.5419 A   |                 |                      |             |
| 590.0 nm  | 0.3048 A            | 0.5051 A                          | 1.0736 A   | 1.6165 A   |                 | Blank                |             |

1.5030 A

Die Standard-ID kann verwendet werden, um den entsprechenden Filter in einem Filterset zu erkennen

Die Standard-ID kann verwendet werden, um den entsprechenden Filter in einem Filterset zu erkennen

635.0 nm

0.3029 A

0.5110 A

0.9992 A

### Zeitplan für Leistungsüberprüfung

Standardarbeitsanweisungen (SOPs) verlangen oft, dass die Leistung von Analysegeräten in regelmäßigen Abständen überprüft wird. Mit der Einstellung "Performance Verification Schedule" (Zeitplan für Leistungsüberprüfung) kann ein solches Intervall voreingestellt werden. Diese Option kann über das Menü "Settings" (Einstellungen) aufgerufen werden

| \$ | Performance Verification Schedule | >   |                      |
|----|-----------------------------------|---|----------------------|
| 0  | Lon                               |   |                      |
|    |                                   | - Performance Verification                      | Schedule             |
|    |                                   | Performance Verification Test                   |                      |
|    |                                   | Wavelength Accuracy Xenon                       |                      |
|    |                                   | Drift at 500nm                                  |                      |
|    |                                   | Noise 0.0A at 500nm                             |                      |
|    |                                   | Noise 1.0A at 500nm                             |                      |
|    |                                   | Noise 2.0A at 500nm                             |                      |
|    |                                   | Baseline Flatness 1000-200 nm                   |                      |
|    |                                   | Photometric Accuracy                            |                      |
|    |                                   | Set expiration                                  |                      |
|    |                                   | Instrument will notify when the results expire. | of this verification |
|    |                                   | Expiration interval 1                           |                      |
|    |                                   | Cancel  | ОК                   |

Das Ablaufintervall kann in Tagen angegeben werden. Wenn diese Option aktiviert ist, weist das Gerät mit einer Meldung darauf hin, dass die Ergebnisse der entsprechenden Leistungsüberprüfung abgelaufen sind. Das Zeitintervall bezieht sich auf das aktuell konfigurierte Gerätedatum. Das Gerät prüft das Ablaufdatum einmal pro Tag – entweder um 12:00 Uhr oder beim ersten Einschalten des Geräts an einem bestimmten Tag. Eine Änderung der Systemzeit nach der Konfiguration eines Verifizierungsplans kann dazu führen, dass das Gerät die Ablaufmeldung anzeigt. Es wird daher empfohlen, alle Ablaufdaten zu überprüfen, wenn die Systemzeit geändert wird.



# KAPITEL 9 Messungen der Absorption, des prozentualen Transmissionsgrads und der Konzentration

## Messungen der Absorption und prozentualen Transmission

## Anwendung mit fester Wellenlänge ("Fixed") für die Testmethode "Basic A-%T-C"

Die Anwendung mit fester Wellenlänge auf Bildschirm 2 versetzt das Gerät in den Modus "instant measurement" (Sofortmessung). Der Benutzer stellt einfach eine Probe in das Instrument und misst diese Probe. Je nach eingestelltem Modus – "Absorbance" (Absorption) (A), "%Transmittance" (proz. Transmission) (%T) oder "Concentration" (Konzentration) werden die Messergebnisse zusammen mit dem Typ der Messung, Datum und Uhrzeit, Wellenlänge und Probengefäßposition für die Messung angezeigt.

Wählen Sie ein neues festes Scan-Fenster aus und bearbeiten Sie das Fenster durch Tippen auf den Touchscreen:

- 1. Bearbeiten Sie den Methodennamen
- Wählen Sie "ABS" oder "%T"
- 3. Wählen Sie die Gleichung
- 4. Wählen Sie die Einheiten aus
- 5. Wählen Sie die Wellenlänge(n) aus
- Geben Sie den/die bekannten Faktor(en) ein 6.

| ÷                                       | New                  |         | ÷                                       | New                      |                      |
|---|----------------------|---------|---|--------------------------|----------------------|
| Method name<br>Fixed 16-Apr-2019<br>ABS | 17/30<br>%T          | Account | Method name<br>Fixed 16-Apr-2019<br>ABS | 9<br>17/30<br><b>%</b> T | Accelute             |
| Equation                                | $ABS(\lambda_1)xF_1$ |         | Equation                                | (ABS(λ1)xF1) / (ABS      | $(\lambda_2) x F_2)$ |
| Unit                                    |                      |         | Unit                                    |                          |                      |
| λ,                                      | 550                  | nm      | λ1                                      | 550                      | nm                   |
| F,                                      | 1.000                |         | F1                                      | 1.000                    |                      |
| Res                                     | ult = ABS(550)x1     |         | λ <sub>2</sub>                          | 650                      | nm                   |
|   |                      |         | F <sub>2</sub>                          | 1.000                    |                      |
|   |                      |         | Result = (A                             | BS(550)x1) / (ABS(650    | i)x1)                |
|   | Continue             |         |   | Continue                 |                      |

Je nach Komplexität der Gleichung erscheinen entweder eine oder zwei Wellenlängen und Faktoren.

Sobald die Methode eingestellt ist, folgen Sie den Schritten zur Blindwerterfassung und Messung. Das Gerät führt automatisch eine Blindwertmessung und eine Messung bei beiden Wellenlängen durch und berechnet die Ergebnisse.

## Verfügbare feste Gleichungen

Wenn Sie auf die aufgelistete Gleichung tippen, erscheint ein Pop-up-Fenster, in dem der Benutzer die gewünschte Gleichung auswählen kann. Je nach Gleichung ändern sich die verfügbaren Parameter, die bearbeitet werden können.

Es werden die folgenden Gleichungen angezeigt, die nur ABS verwenden. Ersetzen Sie "%T" nach Bedarf:

- Direkter Faktor  $ABS(\lambda_1) \times F_1$
- Additive Absorption  $[ABS(\lambda_1) \times F_1] + [ABS(\lambda_2) \times F_2]$
- Differenzielle Absorption [ABS( $\lambda_1$ ) x F<sub>1</sub>] [ABS( $\lambda_2$ ) x F<sub>2</sub>]
- Ratiometrische Absorption [  $ABS(\lambda_1) \times F_1$  ] / [  $ABS(\lambda_2) \times F_2$  ]

Im folgenden Beispiel wurde die ratiometrische Absorptionsgleichung ausgewählt, und die Ergebnisse werden nach Blindwert und Messung berechnet.



Hinweis: Die verschiedenen Gleichungen (additiv, differenziell und Verhältnis) messen die Absorption bei zwei verschiedenen Wellenlängen. Die Referenzwellenlängen-Korrektur beseitigt die Einflüsse einer Probenmatrix. Diese Methoden werden wird in der Regel in Anwendungen zur Qualitätssicherung eingesetzt und ermöglichen einen komfortablen und schnellen diagnostischen Test der Probenqualität.

## Verwendung des "C-Mode" (K-Modus) zur Konzentrationsmessung

In der Anwendung "C-Mode" (C-Modus) auf Bildschirm 2 muss eine Standardprobe mit bekannter Konzentration zur Bestimmung der Konzentration nachfolgender Proben verwendet werden. Wenn die Standardkonzentration genau bekannt ist, werden der ABS-Wert des Standards und der ABS-Wert der unbekannten Probe gemessen und die Probenkonzentration berechnet. Die Messung lässt sich mathematisch wie folgt ausdrücken:

# Standardkonzentration=ProbenkonzentrationABS des StandardsABS der unbekannten Probe

Wenn Sie "C-Mode" (C-Modus) auswählen, wird der nachstehende Bildschirm angezeigt. Durch Tippen auf die editierbaren Felder können Sie die Wellenlänge, den Standardkonzentrationswert und die Maßeinheiten ändern.

Nach dem Durchlauf des Standards wird eine Standardabsorption von 2,361 gemessen und ein Faktor von 0,424 berechnet. Alle nachfolgenden Proben können in regelmäßigen Abständen eingegeben werden, um einen direkten Messwert zu erhalten. Im folgenden Beispiel wird 1,545 mg/l gemessen.

| X C-Mode   |  |         | × C-Mode                     |                |  |  | X C-Mode                 |                |          |                                 |  |
|--|--|---------|------------------------------|----------------|--|--|--------------------------|----------------|----------|---------------------------------|--|
| λ  |  | 660     | nm                           | λ              |  |  | nm                       | λ              |          |                                 |  |
| Standard concentration 1.000   |  |         | Standard concentration 1.000 |                |  | 8  | Standard concentration 1 |                |          |                                 |  |
| Unit   |  |         |                              | Unit mg/L      |  | Unit                                     |                          |                |          |                                 |  |
| Standard absorbance  |  |         | Standard absorbance 2.361    |                | Standard abs                           | Standard absorbance 2.361                |                          |                |          |                                 |  |
| Factor   |  |         |                              | Factor         |  | 0.424                                    |                          | Factor         |          | 0.424                           |  |
| Live display does not support accessories. Ensure the single cell holder is in place |  |         |                              | Live display o | loes not support acces<br>holder is in | Mg/L<br>usories. Ensure the sir<br>place | ngle cell                | Live display o | . 545    | Mg/L<br>Isories. Ensure the sir |  |
| Present blank solution   |  |         |                              | Measuring      |  |  |                          |                | Measurin | l                               |  |
| Blank  |  | Measure |                              | Blank Measure  |  |  |                          | Blank          | Measure  |                                 |  |

# Mehrfachwellenlänge

Die Anwendung "Multi-Wavelength" (Multi-Wellenlängen) steht auf Bildschirm 3 zur Verfügung. Dadurch wird das Gerät in einen Modus versetzt, in dem für die Messung eine Blindprobe und nachfolgende Proben erforderlich sind. Mit der Anwendung "Multi-Wavelength" (Multi-Wellenlängen) werden mehrere Messungen mit festen Wellenlängen durchgeführt. Sie ist eine schnelle Alternative zum Scannen (Anwendung "Scanning" [Scan]), wenn die zu untersuchenden Wellenlängen bekannt sind. Wenn Sie auf das Symbol tippen, wird/werden folgende(n) Bildschirm(e) angezeigt. Tippen Sie auf das Symbol "+", um eine Messung hinzuzufügen.

Bis zu zwanzig (20) feste Wellenlängen können in die Wellenlängentabelle für eine Messung im ABS- oder %T-Modus eingegeben werden. In dem folgenden Beispiel wurden nur fünf (5) für die ABS-Messung ausgewählt. Tippen Sie auf das Speichern-Symbol, um die Scanmethode zu speichern. Tippen Sie auf "Continue" (Weiter), um zum Blindwert und zur Messung zu gelangen.

| <br>Multi-Wavelength Home |              | ÷                         |             | New      |       |           | X                                    | Multi-Wavel         | ength 16-Ap         | r-2019              | :                   |
|---------------------------|--------------|---------------------------|-------------|----------|-------|-----------|--------------------------------------|---------------------|---------------------|---------------------|---------------------|
|                           | $\mathbf{V}$ | Method name<br>Multi-Wave | elength 16- | Apr-2019 | 28/30 | Accessory | Sample base name                     |                     |                     |                     |                     |
|                           |              |                           | ABS         |          | %T    |           | Sample Test 1<br>ABS(400.0)<br>0.079 | ABS(500.0)<br>0.140 | ABS(600.0)<br>1.094 | ABS(700.0)<br>0.243 | ABS(800.0)<br>0.062 |
|                           |              | Wavelength                | table       |          |       |           |                                      |                     |                     |                     |                     |
|                           |              | 400.0                     | 500.0       | 600.0    | 700.0 | 800.0     |                                      |                     |                     |                     |                     |
|                           |              | λ6                        |             |          |       | λ10       |                                      |                     |                     |                     |                     |
|                           |              |                           |             |          |       | λ15       |                                      |                     |                     |                     |                     |
|                           |              | _                         |             |          |       | _         |                                      |                     |                     |                     |                     |
|                           |              | λ16                       |             |          |       | λ20       |                                      |                     |                     |                     |                     |
|                           |              | -                         |             |          |       |           |                                      |                     |                     |                     |                     |
|                           |              |                           |             |          |       |           |                                      |                     |                     |                     |                     |
|                           |              |                           |             |          |       |           |                                      |                     |                     |                     |                     |
|                           |              |                           |             |          |       |           |                                      |                     |                     |                     |                     |
|                           |              |                           |             |          |       |           |                                      |                     |                     |                     |                     |
|                           |              |                           |             |          |       |           |                                      |                     |                     |                     |                     |
|                           |              |                           |             | Continue |       |           | Blank                                |                     | Meas                | ure                 |                     |

### Erneutes Aufrufen einer bestehenden Multi-Wellenlängen-Methode

Multi-Wellenlängen-Methoden können auf der Seite "Multi-Wavelength" (Multi-Wellenlängen) aufgerufen werden, indem eine gespeicherte Methode ausgewählt wird. Wenn die Methode nicht gespeichert wurde, wird sie hier nicht angezeigt.

# 3-Punkt-Netz

Mit der Anwendung "3-Point Net" (3-Punkt-Netz) ermitteln Sie die Bandenhöhe (Peak-Höhe) anhand der geneigten Basislinie zwischen zwei Wellenlängen auf beiden Peak-Seiten. Dieser Analysetyp ist vorteilhaft, wenn die genaue Bandenhöhe für ein bestimmtes Assay erforderlich ist. Durch Multiplizieren der gemessenen Bandenhöhe mit einem Faktor ergibt sich die Konzentration des gemessenen Analyten in den zugehörigen Konzentrationseinheiten. Dieser Abschnitt umfasst:

- Einrichten von Testparametern
- Durchführen von Messungen
- Wiederaufrufen eines Tests

### Einstellung der Parameter für die 3-Punkt-Netz-Methode

### Messungen mit der 3-Punkt-Netz-Methode

Aufrufen einer bestehenden 3-Punkte-Netz-Methode

# Kinetik

Die Methode "Kinetics" (Kinetik) ist ein aktiver Scan mit einer ausgewählten festen Wellenlänge und optionalen Referenzwellenlänge über einen festen Zeitraum, wobei die Daten in einem ausgewählten Intervall und Integrationszeitraum gestreamt werden. Wenn die eingestellte Versuchszeit erreicht ist, wird der Versuch beendet. Die Methoden können unter einem Namen gespeichert und zur Wiederholung eines Kinetik-Scans verwendet werden. Diese Anwendung dient dazu, die Reaktion oder den Zerfall einer Probe im Laufe der Zeit zu beobachten.

| ÷                                 |          | 8                  | $\times$     |   |                              |  |
|-----------------------------------|----------|--------------------|--------------|---|------------------------------|--|
| Method Name<br>Green Dye Kinetics |          | Accessory<br>SETUP | Gree<br>Expe | n Dye Kinetics<br>A 628 nm<br>riment Time ( 5.0 | Ref X nm<br>Omin )           |  |
| λ                                 |          | nm                 |              |   | $00{:}00{:}00 \; \text{min}$ |  |
| Ref λ                             |          | nm                 |              |   |                              |  |
| Minutes                           |          | nds                | A contraste  |   |                              |  |
| Experiment time                   | 5.000    | min                |              |   |                              |  |
| Data interval                     | 200.000  | min                |              |   |                              |  |
| Integration time                  | 0.500    | sec                |              |   |                              |  |
|                                   |          |                    |              |   |                              |  |
|                                   |          |                    |              |   |                              |  |
|                                   |          |                    |              |   |                              |  |
|                                   |          |                    | в            | ank   |                              |  |
|                                   | Continue |                    |              | <u>ß</u>  | 2                            |  |
|                                   | Lun .    |                    |              | (")   | Z)                           |  |



Beachten Sie, dass die Messungen zu den unter "Interval" (Intervall) angegeben Zeitpunkten erfolgen.

Erstellen Sie eine neue Methode mit den dargestellten Einstellungen

> Mit der Anwendung "Kinetics" (Kinetik) messen Sie die Veränderung der Probenabsorption als Funktion der Zeit. Die lokale Steuerungssoftware ermöglicht die Ermittlung einer linearen Rate über einen bestimmten Bereich, der nach der Datenerfassung definiert werden kann. In der enzymatischen Kinetik wird zur Bestimmung der Aktivität häufig ein Faktor mit der Steigung der linearen Ratenangleichung multipliziert.

- Neuskalierung und Neuberechnung der tabellarischen Kinetik-Ergebnisse
- Anpassen der Skala für die Darstellung

Sie können wahlweise mit grafischen oder tabellarischen Daten arbeiten und dabei dieselben Funktionen nutzen. Die Position der Funktionstasten ist allerdings abhängig vom Anzeigetyp.

Hinweis: Mit der Anwendung "Kinetics" (Kinetik) kann nur jeweils ein Probe gemessen werden.



# KAPITEL 10 Wartung/Pflege

Das Spektrophotometer ist langlebig und zuverlässig, daher ist die routinemäßige Wartung minimal. Inhalt dieses Abschnitts:

- Regelmäßige Pflege
- Auswechseln der Sicherung
- Auswechseln der Wolframlampe (nur Orion AquaMate 7100 Vis-Geräte)



**Warnung:** Der Betrieb des Geräts ohne Abdeckung setzt den Bediener potenziell gefährlichen Spannungen und ultravioletter (UV-)Strahlung aus. Wir empfehlen daher, die Arbeiten, die das Entfernen der Geräteabdeckung und den Austausch der elektrischen Komponenten erfordern, nur von autorisierten Kundendienstmitarbeitern durchführen zu lassen. Zu Ihrem eigenen Schutz und zum Schutz des Geräts sollten Sie sich an einen autorisierten Kundendienst wenden, wenn Sie bei der Durchführung von Wartungsarbeiten unsicher sind.

## Regelmäßige Pflege

Die regelmäßige Pflege des Spektrophotometers erfordert nicht viel Zeit. Um den Wartungsaufwand zu minimieren und die Betriebsdauer und Leistung des Geräts zu erhöhen, beachten Sie bitte die folgenden Richtlinien:

- Bringen Sie die Staubschutzabdeckung immer an, wenn das Gerät nicht eingeschaltet ist, um zu verhindern, dass sich Staub im und auf dem Gerät ansammelt.
- Verwenden oder lagern Sie das Gerät nicht in einer korrosiven Umgebung.
- Wischen Sie die Außenseite des Geräts, einschließlich des Touchscreens, vorsichtig mit einem weichen Tuch ab, um Staub und Flecken zu entfernen. Bei Bedarf können Sie Wasser, Isopropylalkohol und andere übliche Laborreinigungsmittel verwenden.
- Wischen Sie verschüttete Flüssigkeiten immer sofort auf, um Schäden am Gerät zu vermeiden oder zu minimieren. Wenn konzentrierte Säuren oder Basen oder Kohlenwasserstoffe auf dem Gerät verschüttet werden, ist der betroffene Bereich sofort zu reinigen.

## Reinigung und Wartung von Fläschchen und Küvetten

Überprüfen Sie sorgfältig den Zustand der Fläschchen, Küvetten und anderen Probenzellen, die zur Messung der Proben verwendet werden. Wenn sie abgesplittert, gesprungen oder zerkratzt sind, müssen die beschädigten Fläschchen unbedingt entsorgt und durch neue ersetzt werden.

Die innere und äußere Sauberkeit der Fläschchen ist aus zwei Gründen für die Qualität der Ergebnisse wichtig: 1) Verunreinigendes Material kann Licht absorbieren, was zu falsch hohen Absorptionswerten führt; und 2) Verunreinigungen im Fläschchen können chemisch mit nachfolgenden Reagenzien oder Standards reagieren, die in das Fläschchen gegeben werden.

Die Reinigungsmethoden hängen in gewissem Maße von der Art des verunreinigenden Materials ab. Es ist wichtig, das zu entfernende Restmaterial im Fläschchen zu kennen. In der folgenden Tabelle finden Sie Empfehlungen zu Reinigungsmethoden, Lösungsmitteln und Materialien.

| Lösungsmittel     | Beispiele                | Empfohlene Reinigungsmethoden                 |  |  |  |  |
|-------------------|--------------------------|---|--|--|--|--|
| Wässrig           |                          | Warmes Wasser mit Reinigungsmittel            |  |  |  |  |
|                   | Proteine, Biologika, DNA | Spülung mit verdünnter Salpetersäure (< 10 %) |  |  |  |  |
|                   |                          | Mit reichlich Wasser abspülen                 |  |  |  |  |
| ) A ( in a series | Solzlögungen             | Spülung mit verdünnter Salpetersäure (< 10 %) |  |  |  |  |
| wassing           | Saiziosungen             | Mit reichlich Wasser abspülen                 |  |  |  |  |
|                   |                          | Warmes Wasser mit Reinigungsmittel            |  |  |  |  |
| Wässrig           | Basislösungen            | Spülung mit verdünnter Salpetersäure (< 10 %) |  |  |  |  |
|                   |                          | Mit reichlich Wasser abspülen                 |  |  |  |  |

| Lösungsmittel | Beispiele Empfohlene Reinigungsmethoden     |  |  |  |  |
|---------------|---|--|--|--|--|
|               |   | Spülen mit organischem Lösungsmittel                                       |  |  |  |
| Organisch     | Kohlenwasserstoffe, kleine                  | Warmes Wasser mit Reinigungsmittel   |  |  |  |
| Organisch     | Moleküle, Öle                               | Spülung mit verdünnter Salpetersäure (< 10 %)                              |  |  |  |
|               |   | Mit reichlich Wasser abspülen  |  |  |  |
| organisch     | Alkoholische Lösungen                       | Mit ähnlichem Alkohol, Aceton oder einem anderen<br>Lösungsmittel abspülen |  |  |  |
|               | _   | Mit reichlich Wasser abspülen  |  |  |  |
|               |   | Spülen mit organischem Lösungsmittel                                       |  |  |  |
| Organisch     | Säurehaltige Lösungen                       | Warmes Wasser mit Reinigungsmittel   |  |  |  |
| Organisch     |   | Spülung mit verdünnter Salpetersäure (< 10 %)                              |  |  |  |
|               |   | Mit reichlich Wasser abspülen  |  |  |  |
|               |   | Spülen mit organischem Lösungsmittel                                       |  |  |  |
| organisch     | Kohlenwasserstoffe, kleine<br>Moleküle, Öle | Warmes Wasser mit Reinigungsmittel   |  |  |  |
| organisch     |   | Spülung mit verdünnter Salpetersäure (< 10 %)                              |  |  |  |
|               |   | Mit reichlich Wasser abspülen  |  |  |  |

Wichtig: Für eine lange Lebensdauer muss das Fläschchen sauber gehalten werden.

- Lagern Sie die Fläschchen oder Küvetten zwischen den Anwendungen niemals über längere Zeit in einem Wasser- oder Lösungsmittelbad. Wenn das verwendete Lösungsmittel trocknet, können sich Verunreinigungen im Wasser oder Lösungsmittel auf der Innenseite des Fläschchens oder der Küvette ablagern und dauerhafte Schäden verursachen.
- Verwenden Sie zum Abwischen der optischen Oberflächen nur Objektivreinigungstücher/papier oder ein feines, weiches Tuch. Die meisten Papierprodukte (z. B. Kosmetiktücher, Papiertücher usw.) enthalten Holzfasern, die das Material der Fläschchen oder Küvetten beschädigen können.
- Stellen Sie sicher, dass alle Fläschchen oder Küvetten am Ende des Tages gut gereinigt sind und nach dem Trocknen in einem geeigneten Behältnis aufbewahrt werden.

| Begriff                          | Definition  |
|----------------------------------|---|
| Verdünnte Säure                  | Verdünnte Salpetersäure (< 10 %)  |
| Säure                            | Salzsäure (5 M) oder Salpetersäure (5 M) (siehe Anmerkung unten)  |
| Spülen mit<br>Lösungsmittel      | Spülen Sie mit dem Lösungsmittel, das ursprünglich zur Auflösung Ihres<br>Analyten verwendet wurde  |
| Mit reichlich Wasser<br>abspülen | Verwenden Sie reines Wasser (deionisiert, destilliert, RO) und spülen Sie mindestens 10 Mal   |
| Reinigungsmittel                 | Verwenden Sie möglichst ein pH-neutrales Reinigungsmittel (Triton X-100), um die Säure zu verdünnen, und spülen Sie mit Wasser nach, um Rückstände zu entfernen |

**Hinweis:** Verwenden Sie keine 5M-Salpetersäure auf antireflexionsbeschichteten Fläschchen oder Küvetten.

**Wichtig:** Die Verwendung eines Ultraschall-Reinigungsbades für Ihre Fläschchen oder Küvetten wird nicht empfohlen. Jedes Reinigungsbad erzeugt eine andere Frequenz; wenn Ihr Bad also mit der Resonanzfrequenz eines Fläschchens oder einer Küvette arbeitet, wird das Fläschchen oder die Küvette brechen. Wenn ein Fläschchen oder eine Küvette in einem Ultraschallbad gereinigt wurde, kann die Garantie durch den Hersteller erlöschen.

Wichtig: Trocknen Sie die Probenzellen nicht in einem Ofen.

Mikro-Durchflusszellen können folgendermaßen gereinigt werden:

- Nach Gebrauch gut mit einem Lösungsmittel ausspülen.
- Einspritzen von verdünnten Säuren, Basen, nicht filtrierenden Reinigungsmitteln oder Bleichmitteln in kurzen Stößen durch die Zelle.
- Aufbewahrung mit destilliertem Wasser in der Zelle.

## Reinigen der Probenkammerfenster

Verwenden Sie kein Aceton oder Scheuermittel, um die Probenkammerfenster zu reinigen. Verwenden Sie stattdessen eine nicht scheuernde Laborreinigungslösung (z. B. eine handelsübliche Messzellen-Reinigungslösung), destilliertes Wasser oder Alkohol.

Verwenden Sie die entsprechende Flüssigkeit und ein weiches, fusselfreies Tuch, um die Fenster zu reinigen. Üben Sie nicht zu viel Druck aus, sonst kann die Oberfläche der Fenster beschädigt werden. Entfernen Sie alle Fingerabdrücke.
#### Auswechseln der Wolfram-Halogenlampe

Dieses Verfahren gilt für das Spektrophotometer Orion AquaMate 7100 Vis. Die Betriebsdauer der Lampe beträgt etwa 1.000 Stunden.

**Warnung:** Die Lampe wird während des Betriebs sehr heiß. Schalten Sie das Gerät vor dem Austauschen einer Lampe aus und lassen Sie es 10 Minuten lang abkühlen.

#### Auswechseln der Wolframlampe:









Anschluss für das Lampengehäuse Führungsschiene für Lampengehäuse







#### Betriebsdauer der Xenon-Lampe

Der Status der Xenon-Lampenbetriebsdauer wird im Menü "Settings" (Einstellungen) angezeigt. Nach der Wartung und dem Austausch der Lampe kann die Lampenbetriebsdauer zurückgesetzt werden.



#### Auswechseln des Xenon-Blitzlichts

Dieses Verfahren gilt für das Spektrophotometer Orion AquaMate 8100 Vis. Die typische Betriebsdauer der Lampe beträgt etwa 3 bis 5 Jahre

**Warnung:** Die Lampe wird während des Betriebs sehr heiß. Schalten Sie das Gerät vor dem Austauschen einer Lampe aus und lassen Sie es 10 Minuten lang abkühlen.

#### Auswechseln des Xenon-Blitzlichts:

- 1. Schalten Sie das Gerät aus
- 2. Nehmen Sie die obere Abdeckung ab.
  - a. Lösen Sie die beiden Schrauben auf der Rückseite des Geräts mit einem Kreuzschlitzschraubendreher Nr. 1. Hinweis: Drehen Sie die beiden Schrauben nicht ganz heraus. Entfernen Sie den Drahtlos-Dongle (falls vorhanden)



b. Ziehen Sie zwei Laschen an der Unterseite in Richtung der Gerätevorderseite.



c. Heben Sie die obere Abdeckung ab, indem Sie sie nach hinten drehen. Achten Sie dabei darauf, dass Sie die Kabel des Druckers und Displays nicht zu stark belasten. Legen Sie die Abdeckung unter die Basis des Geräts, damit der Deckel nicht umkippt.



d. Entfernen Sie mit einem Kreuzschlitzschraubendreher die zwei Schrauben, die den 9-poligen Stecker fixieren.



e. Entfernen Sie mit einem Kreuzschlitzschraubendreher die beiden Schrauben, mit denen die gefräste Halterung am Gussteil befestigt ist.



- 3. Montieren Sie die neue Lampe und die Abdeckung wieder in umgekehrter Reihenfolge.
  - a. BERÜHREN Sie das Fenster der neuen Lampe NICHT.
  - b. Seien Sie vorsichtig bei der Handhabung der neuen Lampe und deren Halterung.
  - Die Lampe ist präzise auf der Halterung ausgerichtet.
  - Die Schrauben, mit denen die Lampe an der Halterung befestigt ist, dürfen nicht verstellt werden.
  - Vergewissern Sie sich, dass die Stifte und Löcher, mit denen die schwarze Halterung am Sockel ausgerichtet ist, richtig eingerastet sind, bevor Sie die Schrauben anziehen, mit denen die neue Lampe am Sockel befestigt wird.
  - Achten Sie darauf, die Schrauben nicht zu überdrehen.

## KAPITEL 11 Kundendienst

#### **Technischer Support**

Wenn Sie Fragen haben oder Unterstützung benötigen, wenden Sie sich an unseren technischen Kundendienst:

- E-Mail: wai.techservbev@thermofisher.com
- Telefon innerhalb der USA: 1-800-225-1480
- Telefon außerhalb der USA: +1-978-232-6000 oder Fax +1-978-232-6031

Bitte wenden Sie sich wegen weiterer Produktinformationen an Ihren Vertriebspartner vor Ort, Ihren Thermo Scientific Orion Vertriebsrepräsentanten oder direkt an uns. Die Kontaktdaten des Geschäftsbereichs "Water and Laboratory Products" (WLP) finden Sie auf der Rückseite dieses Benutzerhandbuchs.

Auf unserer Website unter <u>www.thermoscientific.com/water</u> können Sie sich Thermo Scientific Orion Produkte ansehen und Produktliteratur, Software-Updates, Gebrauchsanweisungen und Benutzerhandbücher sowie zusätzliche Anwendungs- und technische Ressourcen herunterladen.

Die aktuellen Garantieinformationen finden Sie auf der Thermo Scientific Orion Garantiekarte, die sich auf der Thermo Scientific Orion AquaMate-Literatur-CD befindet und online unter <u>www.thermoscientific.com/water</u> verfügbar ist.

### Gerätespezifikationen

| Spezifikationen des Orion AquaMate 8100 UV-Vis-Spektrophotometer |  |  |
|--|--|--|
| Optisches Design   | Doppelstrahl   |  |
| Spektrale Bandbreite   | 2 nm   |  |
| Lichtquelle<br>(Typische Betriebsdauer)                          | Xenon-Blitzlampe (5 Jahre)   |  |
| Detektor   | Zweifache Silizium-Fotodioden  |  |
| Wellenlängenbereich  | 190 bis 1.100 nm   |  |
| Wellenlängen-Genauigkeit   | ± 0,5 nm   |  |
| Wellenlängen-<br>Reproduzierbarkeit                              | ± 0,2 nm   |  |
| Geschwindigkeit des<br>Wellenlängen-Scans                        | Langsam, mittel und schnell (bis zu 1.600 nm/min)  |  |
| Wellenlängen-<br>Datenauflösung                                  | 0,2 nm, 0,5 nm, 1,0 nm, 2,0 nm, 3,0 nm, 5,0 nm   |  |
| Modi für die<br>photometrische Messung                           | Absorption, Transmission (%), Konzentration  |  |
| Photometrischer Bereich  | -2 A bis +3,5 A  |  |
| Photometrische<br>Genauigkeit                                    | ±0,002A bei 0,5 A, ±0,004 A bei 1,0 A, ±0,008 A bei 2,0 A  |  |
| Photometrische<br>Reproduzierbarkeit <sup>1</sup>                | ±0,001 A bis 1 A   |  |
| Photometrisches<br>Rauschen <sup>2</sup>                         | ≤ 0,00020 A bei 0 A bei 260 und 500 nm<br>≤ 0,00030 A bei 1 A bei 260 und 500 nm<br>≤ 0,00040 A bei 2 A bei 260 und 500 nm                 |  |
| Photometrischer Drift <sup>3</sup>                               | < 0,0005 A/Hr (bei 500 nm nach dem Aufwärmen)  |  |
| Photometrisches<br>Streulicht                                    | < 1,0 %T 198 nm (KCl), < 0,05 % T bei 220 nm (Nal), < 0,03 %T bei 340 nm (NaNO2)   |  |
| Display  | 7-Zoll-Farb-Touchscreen, hochauflösend, 800 × 1.280 Pixel  |  |
| Touchscreen  | Handschuhbedienbarer Touchscreen   |  |
| Konnektivität  | USB-Anschluss Typ A für USB-Stick (Vorderseite), USB-Anschluss Typ B für Computer (Rückseite), USB-Anschluss Typ A für Drucker (Rückseite) |  |
| Abmessungen  | 35,5 × 38,5 × 19,5 cm (L × B × H)  |  |
| Gewicht  | 7,5 kg   |  |
| Anforderungen an die<br>Stromversorgung                          | 100 bis 240 V; 50 bis 60 Hz  |  |

<sup>1</sup>Gemessen bei 1,0 A bei 546 nm <sup>2</sup>RMS bei 500 nm. 60 aufeinanderfolgende Messungen. <sup>3</sup>Bei 500 nm nach 1 Stunde Aufwärmzeit

| Spezifikationen des Orion AquaMate 7100 Vis-Spektrophotometer |  |  |
|---|--|--|
| Optisches Design  | Doppelstrahl   |  |
| Spektrale Bandbreite  | 5,0 nm   |  |
| Lichtquelle<br>(Typische Betriebsdauer)                       | Wolfram-Halogen-Lampe (1.000 h)  |  |
| Detektor  | Zweifache Silizium-Fotodioden  |  |
| Wellenlängenbereich   | -3 A bis +3,5 A  |  |
| Wellenlängen-<br>Genauigkeit                                  | ± 0,5 nm   |  |
| Wellenlängen-<br>Reproduzierbarkeit                           | < ± 0,2 nm   |  |
| Geschwindigkeit des<br>Wellenlängen-Scans                     | Automatisch bis zu 1.800 nm/min  |  |
| Wellenlängen-<br>Datenauflösung                               | 0,2 nm, 0,5 nm, 1 nm, 2 nm, 5 nm   |  |
| Modi für die<br>photometrische Messung                        | Absorption, Transmission (%), Konzentration  |  |
| Photometrischer Bereich                                       | -3 A bis +3,5 A  |  |
| Photometrische<br>Genauigkeit                                 | ±0,002A bei 0,5 A, ±0,004 A bei 1,0 A, ±0,008 A bei 2,0 A  |  |
| Photometrische<br>Reproduzierbarkeit <sup>1</sup>             | ±0,001 A bis 1 A   |  |
| Photometrisches<br>Rauschen <sup>2</sup>                      | ≤ 0,00020 A bei 0 A bei 260 und 500 nm<br>≤ 0,00030 A bei 1 A bei 260 und 500 nm<br>≤ 0,00040 A bei 2 A bei 260 und 500 nm                 |  |
| Photometrischer Drift <sup>3</sup>                            | < 0,0010 A/Hr (bei 500 nm nach dem Aufwärmen)  |  |
| Photometrisches<br>Streulicht                                 | < 0,05 T% bei 340 nm und 400 nm  |  |
| Display   | 7-Zoll-Farb-Touchscreen, hochauflösend, 800 × 1.280 Pixel  |  |
| Touchscreen   | Handschuhbedienbarer Touchscreen   |  |
| Konnektivität   | USB-Anschluss Typ A für USB-Stick (Vorderseite), USB-Anschluss Typ B für Computer (Rückseite), USB-Anschluss Typ A für Drucker (Rückseite) |  |
| Abmessungen   | 35,5 × 38,5 × 19,5 cm (L × B × H)  |  |
| Gewicht   | 7,5 kg (19 Pfund)  |  |
| Anforderungen an die<br>Stromversorgung                       | 100 bis 240 V; 50 bis 60 Hz  |  |

<sup>1</sup>Gemessen bei 1,0 A bei 546 nm <sup>2</sup>RMS bei 500 nm. 60 aufeinanderfolgende Messungen. <sup>3</sup>Bei 500 nm nach 1 Stunde Aufwärmzeit

*Hinweis:* Änderungen im Sinne von Produktverbesserungen und Aktualisierungen vorbehalten. Die technischen Daten können sich ohne vorherige Ankündigung ändern.

### Bestellinformationen

| Bestell-Nr. | Beschreibung   |
|-------------|--|
| AQ8100      | AquaMate 8100 UV-Vis-Spektrophotometer mit vorinstallierten Methoden, Halter für<br>Fläschchen und Reagenzgläser von 12 bis 25 mm Außendurchmesser, 10-mm-<br>Vierkantküvettenhalter, Drucker, Netzkabel für Nordamerika, Europa- und VK sowie<br>Staubschutzhülle             |
| AQ8100 APAC | AquaMate 8100 UV-Vis-Spektrophotometer mit vorinstallierten Methoden, Halter für<br>Fläschchen und Reagenzgläser von 12 bis 25 mm Außendurchmesser, 10-mm-<br>Vierkantküvettenhalter, Drucker, Netzkabel für China, Indien und<br>Australien/Neuseeland sowie Staubschutzhaube |
| AQ7100      | AquaMate 7100 Vis Spektrophotometer mit vorinstallierten Methoden, Halterung für<br>Fläschchen und Reagenzgläser von 12 bis 25 mm Außendurchmesser, Netzkabel für<br>Nordamerika, Europa und VK und Staubschutzhaube   |
| AQ7100 APAC | AquaMate 7100 Vis Spektrophotometer mit vorinstallierten Methoden, Halter für<br>Fläschchen und Reagenzgläser von 12 bis 25 mm Außendurchmesser, Netzkabel für<br>China, Indien und Australien/Neuseeland und Staubschutzhülle   |
| AQX1RNDVH   | Fläschchen-/Reagenzglashalter für Fläschchen und Reagenzgläser von 12 bis<br>25 mm Außendurchmesser für Orion™ AquaMate 7100 & 8100 Spektrophotometer  |
| AQX1SQVH    | 10 mm quadratischer Einzelzellenhalter für Orion™ AquaMate 7100 & 8100<br>Spektrophotometer  |
| AQX1LWLVH   | Rechteckiger Küvettenhalter mit langer Schichtdicke für Küvetten mit 20 bis 100 mm<br>Schichtdicke für Orion™ AquaMate 7100 & 8100 Spektrophotometer   |
| AQX1FLTRHDR | Film-/Filterhalter für Filter/Objektive bis zu 50 mm Länge x 80 mm Höhe x 10 mm<br>Dicke für Orion™ AquaMate 7100 & 8100 Spektrophotometer   |
| 840-253100  | AquaMate-Kalibrierungsstandard-Set (erfordert AQX1FLTRHDR Film-/Filterhalter)  |
| AQ71LMPTGST | Ersatz-Wolfram-Halogenlampe, vorjustiert für Orion™ AquaMate 7100  |
| AQ81LMPXEN  | Ersatz-Xenon-Lampeneinheit, vorjustiert für Orion™ AquaMate 8100   |
| AQX1PWRSUP  | Netzteil für Orion™ AquaMate 7100 & 8100 Spektrophotometer +12 V 5 A   |
| AQX1AUCBL   | Netzkabel für Australien (AS/NZS 3112) für Orion™ AquaMate 7100 & 8100<br>Spektrophotometer  |
| AQX1CNCBL   | Netzkabel für China (PRC/3) für Orion™ AquaMate 7100 & 8100 Spektrophotometer  |
| AQX1EUCBL   | EU-Netzkabel (CEE 7/7) für Orion™ AquaMate 7100 & 8100 Spektrophotometer   |
| AQX1INCBL   | Netzkabel für Indien (SABS 164) für Orion™ AquaMate 7100 & 8100<br>Spektrophotometer   |
| AQX1NACBL   | Netzkabel für Nordamerika (NEMA 5-15) für Orion™ AquaMate 7100 & 8100<br>Spektrophotometer   |
| AQX1UKCBL   | Netzkabel für VK (BS 1362) für Orion™ AquaMate 7100 & 8100 Spektrophotometer   |
| AQX1PRNTR   | Snap-on-Drucker für Orion™ AquaMate 7100 & 8100 Spektrophotometer  |
| AQX1PPRPSA  | Selbstklebendes Druckerpapier für AquaMate-Drucker-Zubehör   |
| AQX1PPRSTD  | Standard-Druckerpapier für AquaMate-Drucker-Zubehör  |
| AC2V24      | 24-mm-Rundfläschchen, 12er Pack  |
| AC2V16      | 16-mm-Rundfläschchen, 10er Pack  |
| COD165      | Thermoreaktor für Aufschlussverfahren, 100 / 120 / 150 / 160 / 165 °C-<br>Temperaturregelung   |

| Bestell-Nr. | Beschreibung   |
|-------------|--|
| CODS01      | 1.000 ppm COD-Standard, 475 ml   |
| CODS10      | 10000 ppm COD-Standard, 475 ml   |
| AC2002      | Alkalinität-M, Säure-Indikator-Methode, Reagenztablette, 100 Tests                     |
| AC3002P     | Alkalinität-P, Säure-Indikator-Methode, Reagenztablette, 100 Tests                     |
| AC2027      | Aluminium, Eriochromcyanin-R-Methode, Tablettenreagenz, 50 Tests                       |
| AC4P27      | Aluminium, Eriochromcyanin R-Methode, Pulverreagenz, 100 Tests                         |
| AC2012      | Ammoniak als Stickstoffträger, Indophenolblau-Methode, Tablettenreagenz, 50 Tests      |
| AC4P12      | Ammoniak als Stickstoffträger, Salicylatmethode, Pulverreagenz, 100 Tests              |
| ACR011      | Ammoniak als Stickstoffträger, hoher Bereich, Salicylatmethode, 50 Reaktionsgefäße     |
| ACR012      | Ammoniak als Stickstoffträger, niedriger Bereich, Salicylatmethode, 50 Reaktionsgefäße |
| AC2035      | Brom, DPD-Methode, Tablettenreagenz, 100 Tests   |
| AC2017      | Chlorid, Silbernitrat/Trübungsmethode, Tablettenreagenz, 50 Tests                      |
| AC2070      | Chlor, frei und gesamt, DPD-Methode, Reagenztablette, je 50 Tests                      |
| AC2071      | Chlor, frei, DPD-Methode, Tablettenreagenz, 100 Tests                                  |
| AC2072      | Chlor, gesamt, DPD-Methode, Tablettenreagenz, 100 Tests                                |
| AC3072      | Chlor, gesamt, hoher Bereich, KI/Säure-Methode, Tablettenreagenz, 100 Tests            |
| AC4P71      | Chlor, frei, DPD-Methode, Pulverreagenz, 100 Tests                                     |
| AC4P72      | Chlor, gesamt, DPD-Methode, Pulverreagenz, 100 Tests                                   |
| AC2099      | Chlordioxid, DPD-Methode, Tablettenreagenz, 100 Tests                                  |
| CODL00      | CSB, niedriger Bereich, Dichromat-Reaktor-Aufschlussmethode, 25 Aufschlussröhrchen     |
| CODH00      | CSB, mittlerer Bereich, Dichromat-Reaktor-Aufschlussmethode, 25 Aufschlussröhrchen     |
| CODHP0      | COD, hoher Bereich, Dichromat-Reaktor-Aufschlussmethode, 25 Aufschlussröhrchen         |
| AC2029      | Kupfer, frei und gesamt, Butinolin-Methode, Tablettenreagenz, 50 Tests                 |
| AC4P29      | Kupfer, frei, Bicinchoninat-Methode, Pulverreagenz, 100 Tests                          |
| AC2098      | Cyanursäure, Melamin-Methode, Tablettenreagenz, 100 Tests                              |
| AC2009      | Fluorid, SPADNS-Methode, Flüssigreagenz, 50 Tests                                      |
| AC3032T     | Härte, gesamt, Metallphthalein-Methode, Tablettenreagenz, 100 Tests                    |
| AC2030      | Hydrazin, 4-(Dimethyl-amino)-benzaldehyd-Methode, Pulverreagenz, 30 Tests              |
| AC2078      | Eisen, II & III, PPST-Methode, Tablettenreagenz, 100 Tests                             |
| AC4P78      | Eisen, Ferro, 1,10-Phenanthrolin-Methode, Pulverreagenz, 100 Tests                     |
| AC4P79      | Eisen, gesamt, TPTZ-Methode, Pulverreagenz, 100 Tests                                  |
| AC2055      | Mangan, Formaldoxim-Methode, Tablettenreagenz, 50 Tests                                |
| AC4P54      | Mangan, niedriger Bereich, PAN-Methode, Pulverreagenz, 100 Tests                       |

| Bestell-Nr. | Beschreibung  |
|-------------|---|
| AC4P55      | Mangan, hoher Bereich, Periodat-Oxidationsmethode, Pulverreagenz, 100 Tests                             |
| AC4P42      | Molybdat/Molybdän, Mercaptoessigsäure-Methode, Pulverreagenz, 100 Tests                                 |
| ACR007      | Nitrat als Stickstoffträger, chromotrope Säuremethode, 50 Reaktionsgefäße                               |
| AC2046      | Nitrit als Stickstoffträger, n-(1-Naphthyl)-Ethylendiamin-Methode, Tablettenreagenz, 100 Tests          |
| AC4P46      | Nitrit als Stickstoffträger, niedriger Bereich, Diazotierungsmethode (Azo),<br>Pulverreagenz, 100 Tests |
| ACD004      | Stickstoff, gesamt, niedriger Bereich, Persulfataufschlussmethode, 50 Aufschlussröhrchen                |
| ACD007      | Stickstoff, gesamt, hoher Bereich, Persulfataufschlussmethode, 50 Aufschlussröhrchen                    |
| AC3048      | Ozon, DPD/Glycin-Methode, Tablettenreagenz, 100 Tests   |
| AC2001      | pH-Wert, Phenolrot-Methode, Tablettenreagenz, 100 Tests   |
| AC3001      | pH-Wert, Phenolrot-Methode, Flüssigreagenz, 30 Tests  |
| AC2095-WA   | Phosphat, ortho, niedriger Bereich, Phosphomolybdänsäure/Ascorbinsäure, Tablettenreagenz, 50 Tests      |
| AC2096      | Phosphat, ortho, hoher Bereich, Vanadomolybdat-Methode, Tablettenreagenz, 50 Tests                      |
| AC4P95      | Phosphat, ortho, Phosphomolybdän/Ascorbinsäure-Methode, Pulverreagenz, 100 Tests                        |
| ACD095      | Phosphat als P-Träger, gesamt, Persulfataufschluss/Ascorbinsäuremethode, 50 Aufschlussröhrchen          |
| ACD095AH    | Phosphat als P, hydrolysierbar, Phosphomolybdän/Ascorbinsäure, 50 Aufschlussröhrchen                    |
| ACR095      | Phosphat, ortho, Phosphomolybdän/Ascorbinsäure-Methode, 50 Reaktionsröhrchen                            |
| AC2060      | Kieselsäure, Silicomolybdat-Methode, Tablettenreagenz, 50 Tests   |
| AC2061      | Kieselsäure, Reagenz zur Phosphatentfernung, 100 Tabletten  |
| AC4P60      | Kieselsäure, hoher Bereich, Silicomolybdat-Methode, Pulverreagenz, 100 Tests                            |
| AC4P82      | Sulfat, Bariumsulfat-Turbiditätsmethode, Pulverreagenz, 100 Tests                                       |
| AC2016      | Sulfid, DPD/Katalysator-Methode, Tablettenreagenz, 50 Tests   |
| AC2065      | Zink, Zincon-Methode, Tablettenreagenz, 50 Tests  |

Unter <u>www.thermoscientific.com/water</u> finden Sie eine vollständige Liste aller verfügbaren Thermo Scientific Orion Messgeräte, Elektroden, Lösungen und Zubehör.



# ANHANG A Allgemeine Geräteinformationen

#### Parameter

| Parameter   | Beschreibung   |
|---|--|
| + - X ÷   | Eingabe von mathematischen Operatoren im Taschenrechnermodus (Dienstprogramm)  |
| % of Lamp Life Used<br>(% der -Lampenbetriebsdauer<br>aufgebraucht) | Zeigt den geschätzten Prozentsatz der verbrauchten<br>Lampenbetriebsdauer an, basierend auf einer typischen Xenon-<br>Lampenbetriebsdauer von fünf Jahren (Dienstprogramm) |
| 3-Pt Net (3-Punkt-Netz)   | Berechnet die Peakhöhe aus der tangentialen Basislinie im Diagramm (Scan)  |
| Absorbance (Absorption)   | Eingabe des Absorptionswertes  |
| Accept Name<br>(Name akzeptieren)                                   | Übernimmt den angezeigten Namenseintrag<br>(Testname und Bearbeiten [Einheiten])   |
| Add Character<br>(Zeichen hinzufügen)                               | Fügt ein hervorgehobenes Zeichen zum Namenseintrag hinzu (Testname und Bearbeiten [Einheiten])   |
| Add nm (nm hinzufügen)  | Erweitert die Liste um eine Wellenlänge und einen Faktor bei Multi-<br>Wellenlängentests und einigen Leistungsüberprüfungstests  |
| Area (Fläche)   | Berechnet die Fläche unter dem Peak im Diagramm (Scan)   |
| AutoPrint (AutoDruck)   | Schaltet den automatischen Ausdruck ein oder aus   |
| Autoscale (Autoskalierung)  | Skaliert das Diagramm auf den ursprünglichen Bereich der X- und Y-<br>Achse zurück (Kinetik, Scan)   |
| Baseline Expiration<br>(Basislinien-Ablaufdatum)                    | Eingabe des Zeitpunkts, zu dem die Basislinie für Scantests erneut erfasst werden muss (Dienstprogramm)  |
| Beeper (Tonausgabe)   | Schaltet das akustische Signal für Tastendrucke ein und aus (Dienstprogramm)   |
| Calculation Baseline<br>(Berechnung Basislinie)                     | Wählt die Null-Basislinie oder die tangentiale Basislinie zur<br>Berechnung der Fläche unter dem Peak im Diagramm (Scan)   |

| Parameter  | Beschreibung   |
|--|--|
| Calculator (Rechner)   | Aktiviert den Taschenrechner-Modus (Dienstprogramm)  |
| Cell Position #<br>(Zellen-Position Nr.)   | Zeigt die Position des Fläschchens oder der Küvette im Strahlengang<br>an (nur mit den Einstellungen des Probenpositionierers "Auto 6" oder<br>"Auto 3") |
| Change Mode (Modus ändern)<br>Change to Abs<br>(Ändern in Abs.)<br>Change to %T<br>(Wechsel zu %T) | Schaltet die Messmodi um ("Basic A-%T-C" und einige Tests zur Leistungsüberprüfung)  |
| Collect Baseline<br>(Basislinie erfassen)  | Startet die Erfassung der Basislinie (Scan)  |
| Concentration (Konzentration)  | Legt den Konzentrationswert fest   |
| Conc. of Standard<br>(Konz. von Standard)  | Zeigt den eingegebenen Konzentrationswert an (Adv A-%T-C)  |
| Cursor   | Wechselt in den Cursorverfolgungsmodus, um Datenpunkte im<br>Diagramm anzuzeigen (Kinetik, Scan)   |
| ←Cursor<br>Cursor→   | Bewegt den Cursor im Diagramm nach rechts oder links und zeigt die Daten der einzelnen Punkte an (Kinetik, Scan)   |
| Curve Fit (Kurvenanpassung)  | Wählt die Berechnungsart der Linienanpassung aus (Standardkurve)   |
| Data File Name<br>(Datendateiname)   | Ermöglicht die Eingabe eines Namens für die Datendatei, wenn<br>"AutoSave" (Aut. Speichern) aktiviert ist  |
| Date Standards Measured<br>(Datum der<br>Standardmessung)  | Zeigt das Datum an, an dem die Standards zuletzt gemessen wurden (Standardkurve)   |
| Date/Time Setup<br>(Datum/Uhrzeit einstellen)  | Gibt die aktuellen Datums- und Zeiteinstellungen für das Gerät ein (Dienstprogramm)  |
| Delay Time<br>(Verzögerungszeit)   | Eingabe der Zeit vom Beginn des Tests bis zur ersten Messung;<br>ermöglicht die Äquilibrierung der Probe (Adv A-%T-C und Kinetik)                        |
| Delete Character<br>(Zeichen löschen)  | Löscht das letzte Zeichen des Namenseintrags (Testname und Bearbeiten [Einheiten])   |
| Delete File (Datei löschen)  | Löscht einen Test oder eine Datendatei aus dem Verzeichnis für gespeicherte Tests (Dienstprogramm)   |
| Delete Name (Name löschen)   | Löscht den gesamten Namen, um einen neuen Eintrag zu ermöglichen (Testname und Bearbeiten [Einheiten])   |
| Delete nm (nm löschen)   | Entfernt eine Wellenlänge und einen Faktor aus der Liste (Multiwellenlängen- und Leistungsüberprüfung)   |
| Diluent Volume (Volumen des  | Eingabe des Volumens des vor der Messung hinzugefügten   |
| Dilution Multiplier  | Zeigt den Faktor an, der zur Korrektur der Probenverdünnung  |
| (Verdünnungsmultiplikator)   | verwendet wird   |
| Display Activity<br>(Aktivität anzeigen)   | Gibt an, ob die Ergebnisse die Proteinkonzentration enthalten sollen   |
| DNA ε(260)   | Berechnet den Extinktionskoeffizienten   |
| DNA-Faktor   | Eingabe des Faktors zur Berechnung der DNA-Konzentration (DNA-Biotests)  |
| Edit (Bearbeiten)  | Änderung einer Wellenlänge und eines Faktors aus der Liste (Multiwellenlängen- und Leistungsüberprüfung)   |

| Parameter  | Beschreibung  |
|--|---|
| Edit Curve (Kurve bearbeiten)                    | Dient zum Modifizieren des Graphen (Kinetik)  |
| Edit Data (Daten bearbeiten)                     | Wählt einen Teil der Daten in einer Tabelle zur Neuberechnung des Ergebnisses aus (Kinetik, Scan)   |
| Edit Graph (Grafik bearbeiten)                   | Dient zum Modifizieren des Graphen (Scan)   |
| Edit Scale (Skala bearbeiten)                    | Ändern der Achsenskalen des Diagramms und Anzeigen einzelner Datenpunkte (Scan)   |
| Factor (Faktor)                                  | Eingabe eines Faktors zur Umrechnung eines Bezugspunkts in ein<br>Ergebnis<br>Abs. x Faktor 1 = Konzentrationsergebnis<br>Abs./min x Faktor 2 = Ergebnis der Kinetik<br>Kann eingegeben oder aus Konzentration und Absorption in "Adv A-<br>%T-C" berechnet werden  |
| Factor 1 (Faktor 1)                              | Eingabe eines Faktors zur Umrechnung eines Bezugspunkts in ein<br>Ergebnis<br>Abs(WL1) x Faktor = Ergebnis (Abs-Verhältnis, Abs-Differenz,<br>Multiwellenlänge)   |
| Factor 2 (Faktor 2)                              | Eingabe eines Faktors zur Umrechnung eines Bezugspunkts in ein<br>Ergebnis<br>Abs(WL2) x Faktor = Ergebnis (Abs-Verhältnis, Abs-Differenz,<br>Multiwellenlänge)   |
| Factor 3-31 (Faktor 3-31)                        | Eingabe eines Faktors zur Umrechnung eines Bezugspunkts in ein<br>Ergebnis<br>Abs (WL3-31) x Faktor = Ergebnis (Multiwellenlänge)   |
| Graph (Diagramm)                                 | Zeigt ein Diagramm der erfassten Daten an (Kinetik, Scan)   |
| ID # (ID-Nr.)                                    | Eingabe der numerischen Kennung für die Messung; automatische Erhöhung während des Tests bis zum Ausschalten (auf 0 gesetzt)  |
| Instrument Serial Number<br>(Geräteseriennummer) | Anzeige der Seriennummer des Geräts (Dienstprogramm)  |
| Intercept (Schnittpunkt)                         | Eingabe, wo die Linie die Y-Achse kreuzt (Abs. bei Konzentration = 0)   |
| Interval (Intervall)                             | Gibt den Wellenlängenbereich zwischen den Datenpunkten ein (Scan)   |
| Interval Time (Intervallzeit)                    | Eingabe der Zeit zwischen wiederholten Messungen (Kinetik)  |
| Linearity Value<br>(Linearitätswert)             | Gibt einen Linearitätswert ein (Kinetik)<br>Um die Linearität der Reaktion während der Messung zu bestimmen,<br>bietet das Gerät einen Linearitätsparameter. Dies ist die Differenz<br>zwischen den Absorptionsänderungen zweier Messungen, wie im<br>folgenden Beispiel gezeigt:<br>Zeit Abs. ?A Linearität<br>1 .1<br>2 .2 .1<br>3 .29 .09 P<br>4 .38 .09 P<br>5 .46 .08 P<br>6 .52 .06 F<br>Die Linearität ist das "?A" zwischen den "?A"-Berechnungen; P = Pass<br>(Bestanden) und F = Fail (Nicht bestanden) |
| Load Test (Belastungstest)                       | Lagt den markierten Test aus dem Verzeichnis der gespeicherten<br>Tests in den aktiven Speicher und stellt das Gerät auf die<br>Testparameter ein (Dienstprogramm)  |

| Parameter  | Beschreibung  |
|--|---|
| Lock/Unlock<br>(Sperren/Entsperren)                            | Wird verwendet, um gespeicherte Tests vor versehentlichem Löschen<br>oder Ändern zu schützen; fragt nach einem Passwort, damit der<br>Benutzer die Datei sperren oder entsperren kann (Dienstprogramm)                                    |
| Low/High Limits<br>(Untere/obere Grenzen)                      | Zur Eingabe der niedrigsten und höchsten akzeptablen Ergebnisse,<br>außerhalb derer das Ergebnis als "Niedrig" oder "Hoch"<br>gekennzeichnet wird (Adv A-%T-C, Standardkurve, AbsVerh., Abs<br>Diff., Kinetik, 3-Pt Net, einige Biotests) |
| Math   | Zugriff auf Bearbeitungsfunktionen des Graphen (Scan)   |
| Measure Blank (Blindprobe messen; als Funktionstaste)          | Startet die Messung der Blindprobe  |
| Measure Blank (Blindprobe messen; als Testparameter)           | Wählt die Häufigkeit des Nullabgleichs (Blindwert) des Geräts als<br>einmalig oder bei jeder Messung (Kinetik)  |
| Measurement Mode<br>(Messmodus)                                | Wählt die Art der für eine Messung gemeldeten photometrischen Daten (Abs, %T, Konz.) in A-%T-C, Kinetik, Scan, Multiwellenlänge   |
| Measure Samples<br>(Proben messen)                             | Startet die Messung von Proben  |
| Max, X   | Eingabe eines maximalen X-Wertes zur manuellen Skalierung des<br>Diagramms (Kinetik, Scan)  |
| Max, Y   | Eingabe eines maximalen Y-Wertes zur manuellen Skalierung des Diagramms (Kinetik, Scan)   |
| Min, X   | Eingabe eines minimalen X-Wertes zur manuellen Neuskalierung des Diagramms (Kinetik, Scan)  |
| Min, Y   | Eingabe eines minimalen Y-Wertes zur manuellen Skalierung des Diagramms (Kinetik, Scan)   |
| Next Cursor<br>(Nächster Cursor)                               | Auswahl eines Cursorpunktes in Funktionen, die mehr als eine<br>Cursoreinstellung verwenden: Scan-Fläche und Scan-3-Punkt-Netz-<br>Berechnungen in der Grafik (Scan)  |
| Number of Matched Cuvettes<br>(Anzahl angepasster<br>Küvetten) | Angabe der Anzahl der Fläschchen oder Küvetten, die im Korrekturprogramm ausgeführt werden sollen (maximal 5)   |
| Number of Samples<br>(Anzahl der Proben)                       | Anzahl der zu messenden Proben im Test (nicht verfügbar in Kinetik oder Scan)   |
| Number of Standards<br>(Anzahl der Standards)                  | Anzahl der zu messenden Standards für die Standardkurve   |
| Printer (Drucker)  | Auswahl des Ausgabemodus als RS-232 oder Parallel (Dienstprogramm)  |
| Protein Factor<br>(Protein-Faktor)                             | Eingabe des Faktors zur Berechnung der Proteinkonzentration (DNA-Biotests)  |
| Ref. Wavelength<br>(RefWellenlänge)                            | Eingabe eines Referenzwellenlängenwerts. Für jede gemeldete<br>Messung wird die analytische Wellenlänge und die<br>Referenzwellenlänge gemessen<br>Gemeldeter Messwert = Abs. bei Analytische WL - Abs. bei Referenz-<br>WL               |
| Ref. Wavelength Correction (RefWellenlängen-Korrektur)         | Ein-/Ausschalten der Referenz-Wellenlängenkorrektur   |
| Run Standard<br>(Standarddurchlauf)                            | Wechselt zum Bildschirm für die Eingabe von Standards   |
| Run Test (Testdurchlauf)                                       | Wechselt zum Bildschirm für die Datenerfassung  |

| Parameter  | Beschreibung   |
|--|--|
| Sample Positioner<br>(Probenpositionierer)                                 | Auswahl des Typs des Probenpositionierers:<br>1 Cell (1 Zelle) = keine Bewegung; Blindwertmessung und Messung<br>der Probe in derselben Position<br>Manual 6 (Manuell 6) = Probengefäßhalter-Karussell wird über den<br>Touchscreen des Probenpositionierers bewegt; Nullstellung immer auf<br>Position B, dann Rückkehr zur eingestellten Position zum Start der<br>Messung<br>Auto 3 = Probengefäßhalter-Karussell fährt automatisch auf B, 2, 4<br>(Nullstellung immer auf Position B, dann auf Position 2, um die<br>Messung zu starten)<br>Auto 6 = Probengefäßhalter-Karussell fährt automatisch auf die<br>Positionen B, 1, 2, 3, 4, 5 (Nullstellung immer auf Position B, dann auf<br>Position 1, um die Messung zu starten) |
| Sample Volume<br>(Probenvolumen)   | Eingabe des Gesamtvolumens der Probe (bei einigen Biotests mit Verdünnungsmultiplikator)   |
| Save Test (Test speichern)   | Speichert alle Parameter des aktuellen Tests im internen Speicher für<br>einen späteren Abruf  |
| Scan Speed<br>(Scan-Geschwindigkeit)                                       | Auswahl der Geschwindigkeit (nm/min) eines Scan als "Slow"<br>(Langsam), "Medium" (Mittel) oder "Fast" (Schnell) (Scan)  |
| Screen Contrast<br>(Bildschirmkontrast)                                    | Verbesserung der Anzeigesichtbarkeit durch Veränderung des<br>Kontrasts zwischen Hintergrund und Text (Dienstprogramm)   |
| Select Test (Test auswählen)   | Kennzeichnung eines markierten Testnamens mit ">", um den Test in<br>das SmartStart-Menü aufzunehmen (Verzeichnis für gespeicherte<br>Tests des Dienstprogramms)   |
| Set Max. X<br>(Einstellung Max. X)<br>Set Min. X<br>((Einstellung Min. X)) | Legt die Cursorposition im Diagramm als minimalen X- und maximalen<br>X-Wert für die Neuberechnung der Rate fest (Kinetik)   |
| Set nms (nms einstellen)   | Eingabe und Bearbeitung von Wellenlänge und Faktorwerten   |
| Set Options<br>(Optionen einstellen)                                       | Auswahl der Faktor-Eingabe oder der Basislinie zur Berechnung der Fläche unter dem Peak im Diagramm (Scan)   |
| Setup Correction<br>(Setup-Korrektur)                                      | Startet das Verfahren zur Erfassung der erforderlichen Daten zur Korrektur von Absorptions-Unterschieden zwischen Fläschchen oder Küvetten   |
| Slope (Steigung)   | Eingabe des Abs/Konzentrationswertes (Standardkurve)   |
| Smoothing (Glättung)   | Ein-/Ausschalten der Datenglättung (Scan)  |
| Software Revision  | Zeigt die Version der Firmware im Gerät an (Dienstprogramm)  |
| SRE Tolerance<br>(SRE-Toleranz)  | Acceptable minimum stray light (Zulässiges Mindeststreulicht)  |
| Standard Concentrations<br>(Standard-Konzentrationen)                      | Eingabe der Konzentration der Standards, die zur Erstellung der Standardkurve für den Test verwendet werden  |
| Standby  | Auswahl der seit dem letzten Tastendruck oder der letzten<br>Geräteaktivität vergangenen Zeit und Herunterfahren des Geräts, um<br>die Betriebsdauer der Lampe zu verlängern (Dienstprogramm)  |
| Start Wavelength<br>(Start-Wellenlänge)                                    | Eingabe der Anfangswellenlänge für einen Scan (Scan)   |

| Parameter  | Beschreibung   |
|--|--|
| Statistics (Statistik)   | Schaltet die Statistik ein oder aus; berechnet den Durchschnitt und die<br>Standardabweichung der Ergebnisse, wenn sie eingeschaltet ist; die<br>Statistikregister werden gelöscht, wenn sie ausgeschaltet ist, wenn das<br>Gerät ausgeschaltet wird, wenn die Testparameter geändert werden<br>und wenn der Test gespeichert oder erneut gespeichert wird (alle<br>Testtypen außer Kinetik, Scan, Multiwellenlänge) |
| Std Concentration<br>(StdKonzentration)                        | Eingabe der Konzentration des Analyten in der Standardlösung   |
| Stop Wavelength<br>(Stopp-Wellenlänge)                         | Eingabe der Abschluss-Wellenlänge für einen Scan (Scan)  |
| Stored Tests Directory<br>(Verzeichnis gespeicherter<br>Tests) | Zeigt eine Liste der im Gerät gespeicherten Tests an (Dienstprogramm)  |
| Tabular (Tabellarisch)   | Zeigt eine Liste der erfassten Daten an (Kinetik, Scan)  |
| Testname   | Der Bediener gibt einen alphanumerischen Namen (maximal<br>16 Zeichen) für den Test ein; der Name wird auf dem Datenausdruck<br>erscheinen und, wenn der Test gespeichert wird, auf dem Bildschirm<br>"Utility Test Directory" (Dienstprogramm-Testverzeichnis) (verfügbar in<br>allen Tests) angezeigt  |
| Total Run Time<br>(Gesamtlaufzeit)                             | Gibt die Zeit vom Start des Laufs bis zum Ende des Tests ein;<br>entspricht der Verzögerungszeit + Intervallzeiten + Messzeiten<br>(Kinetik)   |
| Units (Einheiten)  | Wählt oder erstellt Einheitenetiketten für Ergebnisse [alle<br>gespeicherten Tests außer "Abs Ratio" (AbsVerh.), "Scanning"<br>(Scan), "Cell Growth" (Zell-Anstieg)]   |
| Unselect Test<br>(Test deaktivieren)                           | Entfernt die Markierung ">" des hervorgehobenen Testnamens, um<br>den Test aus dem SmartStart-Menü zu entfernen (Verzeichnis der<br>gespeicherten Tests des Dienstprogramms)   |
| Wavelength (Wellenlänge)                                       | Eingabe von Werten für analytische Wellenlängen  |

## Berechnungen für die Software

| Berechnung                               | Berechnungsformel   |
|--|---|
| Standardkurven                           |   |
| Teil-Summen                              | $SX = \sum x_i$<br>$SY = \sum y_i$<br>$SXX = \sum x_i^2$<br>$SYY = \sum y_i^2$<br>$SXY = \sum x_i y_i$<br>$SQX = \sum (x_i - \bar{x})^2 = N * SXX * SX^2$<br>$SQY = \sum (y_i - \bar{y})^2 = N * SYY * SY^2$<br>$SSXY = \sum (x_i - \bar{x})^2 (y_i - \bar{y})^2 = N * SXY - SX * SY$<br>Dabei gilt:<br>xi = Konzentration des i-ten Standards<br>yi = Absorption des i-ten Standards<br>N = Anzahl der Standards   |
| Lineare Regression<br>(allgemeiner Fall) | $\begin{array}{l} A = A(c) \\\\ Dabei gilt: \\\\ A = Absorption \\\\ c = Konzentration \\\\ A(c) \text{ ist durch folgende Gleichungsform definiert:} \\\\ A(c) = a_4c^4 + a_3c^3 + a_2c^2 + a_1c + a_0 \\\\\\ Dabei gilt: \\\\ a_0 = Schnittpunkt der Y \cdot Achse \\\\ a_{1a_4} = Koeffizienten \\\\ (Berechnung der Koeffizienten erfolgt nach der Methode der kleinsten \\\\ Quadrate) \end{array}$  |
| Lineare Regression<br>durch Null         | <ul> <li>A = a1 *(c)</li> <li>Dabei gilt:</li> <li>A = Absorption</li> <li>c = Konzentration</li> <li>a1 = Steigung</li> <li>Die Berechnung der Steigung erfolgt als: a1 = SXY/SXX</li> <li>Voraussetzungen für dieses Modell:</li> <li>Steigung ist nicht gleich "Null" oder "unendlich"</li> <li>Mindestens ein Standarddatenpunkt mit einer Konzentration &gt; 0 vorhanden</li> <li>Die Absorption des Blindwertes der Konzentration 0 = 0A</li> </ul> |

| Berechnung                       | Berechnungsformel   |
|----------------------------------|---|
| Segmentiertes Modell             | <ul> <li>Voraussetzungen f ür das segmentierte Modell:</li> <li>Es sind Daten f ür mindestens zwei Standarddatenpunkte mit<br/>unterschiedlicher Konzentration und Absorption vorhanden</li> <li>Die Steigungen aller Segmente m üssen ansteigend (positiv) oder<br/>absteigend (negativ) sein</li> </ul>   |
| Gültigkeit von<br>Standardkurven | $A(c_1) > A(c_2)$ für alle $c_1 > c_2$  |
|                                  | oder  |
|                                  | $A(c_1) < A(c_2)$ für alle $c_1 > c_2$  |
|                                  | Dabei gilt:<br>A = Absorption<br>c <sub>1</sub> , c <sub>2</sub> = Konzentration  |
|                                  | Graph einer gültigen nichtlinearen Standardkurve:   |
|                                  | A3SORBANCE<br>RANGE<br>WEASURED<br>ABSORBANCE<br>UNDERNED<br>CONCEN-  |
|                                  | Wenn dies nicht der Fall ist, gibt es mehr als eine Lösung innerhalb des<br>angegebenen Bereichs und die Meldung "Curve cannot be used to determine<br>sample concentrations – it may produce ambiguous results" (Kurve kann nicht<br>zur Bestimmung der Probenkonzentrationen verwendet werden – sie kann zu<br>nicht eindeutigen Ergebnissen führen) wird angezeigt, wenn die Kurve<br>betrachtet wird. |
|                                  | Graph einer ungültigen nichtlinearen Standardkurve:   |
|                                  | MEASURED ABSORBANCE   |

| Berechnung   | Berechnungsformel  |
|--|--|
| Statistik (allgemeiner<br>Fall der linearen<br>Regression) | $\sigma = \sqrt{\frac{\sum (y_i - \overline{y})^2}{N - n - 1}}$ Dabei gilt:<br>N = Grad des Polynoms<br>$r = \frac{ SSXY }{\sqrt{SQX * SQY}}$ Die Berechnung des Korrelationskoeffizienten gilt nur für lineare<br>Regressionskurven erster Ordnung (Polynome ersten Grades)                                     |
| Lineare Regression<br>durch Nullmodell                     | $\sigma = \sqrt{\frac{SYY - (a_1 * SXY)}{N - 1}}$  |
| Absorptionsverhältnis                                      | $\frac{Abs\lambda_1}{Abs\lambda_2}  \text{oder}  \frac{Abs\lambda_1 - Abs_{ref}}{Abs\lambda_2 - Abs_{ref}}$  |
| Absorptionsdifferenz                                       | Ergebnis = Absλ <sub>1</sub> * Faktor <sub>1</sub> – Absλ <sub>2</sub> * Faktor <sub>2</sub><br>oder<br>Ergebnis = (Absλ <sub>1</sub> – Absλ <sub>ref</sub> ) * Faktor <sub>1</sub> – (Absλ <sub>2</sub> – Absλ <sub>ref</sub> ) * Faktor <sub>2</sub>   |
| 3-Punkt-Netz   | Basislinienkorrektur für Absorption =<br>$A_2 - \left(A_3 + \left([A_1 - A_2] * \frac{\lambda_3 - \lambda_2}{\lambda_3 - \lambda_1}\right)\right)$<br>3-Punkt-Netz-Absorptionskurve der Sample (Probe):  |
| 3-Punkt-Netz<br>(ASTM E16904)                              | Basislinienkorrektur für Absorption =<br>$A_{2} - \left(A_{3} + \left([A_{2} - A_{3}] * \frac{\lambda_{3} - \lambda_{1}}{\lambda_{3} - \lambda_{2}}\right)\right)$ $\downarrow 0 \qquad $ |

#### thermoscientific.com/water

© 2019 Thermo Fisher Scientific Inc. Alle Rechte vorbehalten. Microsoft und Windows sind eingetragene Warenzeichen der Microsoft Corporation. Alle übrigen Marken sind Eigentum von Thermo Fisher Scientific Inc. und seinen Tochtergesellschaften.

Wasser- und Laborprodukte

Nordamerika Gebührenfrei: 1-800-225-1480 Tel.: 1-978-232-6000 Info.water@thermofisher.com

Deutschland Tel.: (49) 6184-90-6000 info.water.uk@thermofisher.com

China Tel.: (86) 21-68654588 wai.asia@thermofisher.com

Indien Tel.: (91) 22-4157-8800 wai.asia@thermofisher.com

Singapur Tel.: (65) 6778-6876 wai.asia@thermofisher.com

Japan Tel.: (81) 045-453-9175 wai.asia@thermofisher.com

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

