



www.thermofisher.com

Potassium Tellurite 3.5%

REF SR0030J

* This instructions for use (IFU) document is intended to be read in conjunction with the IFU for Hoyle Medium Base (CM0083B).

Intended Use

SR0003J is a selective supplement to be used with Hoyle Medium Base (Dehydrated) (CM0083B) for the isolation and differentiation of *Corynebacterium diphtheriae* from wound and upper respiratory samples, including throat and nasal swabs.

The device is for professional use only, is not automated, nor is it a companion diagnostic

Summary and Explanation

Corynebacterium species are Gram-positive rods with clubbed end appearance and exist in single or pairs. The medically relevant corynebacteria include *Corynebacterium pseudotuberculosis*, *Corynebacterium ulcerans* and the most toxigenic form *Corynebacterium diphtheriae*, with four biovars: *intermedius*, *gravis*, *mitis* and *belfanti*¹. Diphtheria is an illness caused by the toxigenic strains of *C. diphtheriae* and less often by *C. ulcerans*². The only known reservoir of *C. diphtheriae* is humans³.

The clinical manifestation of diphtheria can be described as respiratory diphtheria which occur as invasive or noninvasive in the respiratory tract or cutaneous diphtheria associated with skin lesions. The mode of transmission of disease can occur either through droplets from symptomatic or asymptomatic patients or from contaminated milk or dairy products. The onset of symptoms starts with a mild fever and sore throat in two to four days. During the infection exotoxin induces the formation of a white glossy necrotic coagulum in the walls of the respiratory tract which then turns into grey to black patches in appearance leading to suffocation and aspiration of membrane which ultimately leads to death in patients of all age groups⁴. However, with a timely administration of antitoxin and antimicrobial therapy, the illness can be treated completely⁵.

Principle of Method

Potassium tellurite acts as a selective agent and has inhibitory activity against most gram-positive and gram-negative bacteria except *Corynebacterium* species. *C. diphtheriae* reduces potassium tellurite to tellurium and produce grey black-coloured colonies

Typical Formula

Potassium tellurite 2mL

Materials Provided

SR0030J: 10x 2ml vials. Yield will depend upon formulation.

Materials Required but Not Supplied

- Hoyle Medium Base (dehydrated) product code CM0083B
- Laked Horse blood SR0048C

Storage

- Store product in its original packaging between 2°C and 8°C.
- Keep container tightly closed.
- The product may be used until the expiry date stated on the label.
- Protect from moisture.
- Store away from light.
- Allow reconstituted product to equilibrate to room temperature before use.

Warnings and Precautions

- DO NOT USE SUPPLEMENTS IF YOU ARE PREGNANT OR CONTEMPLATING PREGNANCY.
- Each vial is for single use. Do not re-use.
- For in vitro diagnostic use only.
- For professional use only.
- Inspect the product packaging before first use.
- Do not use the product if there is any visible damage to the packaging or vials.
- Do not use the product beyond the stated expiry date.
- Do not use the device if signs of contamination are present.
- Do not use the device if the colour has changed or there are other signs of deterioration.
- It is the responsibility of each laboratory to manage waste produced according to their nature and degree of hazard and to have them treated or disposed of in accordance with any federal, state and local applicable regulations. Directions should be read and followed carefully. This includes the disposal of used or unused reagents as well as any other contaminated disposable material following procedures for infectious or potentially infectious products.

Refer to the Safety Data Sheet (SDS) for safe handling and disposal of the product (www.thermofisher.com).

Serious Incidents

Any serious incident that has occurred in relation to the device shall be reported to the manufacturer and the relevant regulatory authority in which the user and/or the patient is established.

Specimen Collection, Handling and Storage

Specimen should be collected and handled following local recommended guidelines, such as the UK Standards for Microbiology Investigations (UK SMI) ID 2 and Q 5.

Procedure

- Suspend 4g of Hoyle Medium Base (CM0083B) in 100ml of distilled water and bring to the boil to dissolve completely.
- Sterilise by autoclaving at 121°C for 15 minutes.
- Cool to 50°C.
- Add 5ml of Laked Horse Blood (SR0048C) and 1ml Potassium Tellurite (SR0030J).
- Mix well and pour into sterile containers

Refer to the IFU for CM0083B. IFUs can be found at www.thermofisher.com

Bibliography

1. Public Health England. 2014 'Identification of *Corynebacterium* species' ID_23i4.1 <https://www.gov.uk/government/publications/smi-id-2-identification-of-corynebacterium-species>
2. Mokrousov, Igor. "Corynebacterium diphtheriae: genome diversity, population structure and

- genotyping perspectives." *Infection, Genetics and Evolution* 9, no. 1 (2009): 1-15.
- Rajamani Sekar, S. K., B. Veeraraghavan, S. Anandan, N. K. Devanga Ragupathi, L. Sangal, and S. Joshi. "Strengthening the laboratory diagnosis of pathogenic *Corynebacterium* species in the Vaccine era." *Letters in applied microbiology* 65, no. 5 (2017): 354-365.
 - Zakikhany, Katherina, and Androulla Efstratiou. "Diphtheria in Europe: current problems and new challenges." *Future microbiology* 7, no. 5 (2012): 595-607.
 - Sharma, Naresh Chand, Androulla Efstratiou, Igor Mokrousov, Ankur Mutreja, Bhabatosh Das, and Thandavarayan Ramamurthy. "Diphtheria (Primer)." *Nature Reviews: Disease Primers* (2019).

Made in the United Kingdom	Made in the United Kingdom
----------------------------	----------------------------

©2022 Thermo Fisher Scientific Inc. All rights reserved. All other trademarks are the property of Thermo Fisher Scientific Inc. and its subsidiaries.



Oxoid Limited, Wade Road, Basingstoke, Hampshire, RG24 8PW, UK

For technical assistance please contact your local distributor.

Symbol Legend

Symbol	Definition
	Catalogue number
	In Vitro Diagnostic Medical Device
	Batch code
	Temperature limit
	Use-by date
	Keep away from sunlight
	Do not re-use
	Consult instructions for use or consult electronic instructions for use
	Contains sufficient for <n> tests
	Do not use if packaging damaged and consult instructions for use
	Manufacturer
	Authorized representative in the European Community/ European Union
	European Conformity Assessment
	UK Conformity Assessment
	Unique device identifier

Revision information

Version	Modifications Introduced
1.0	2022-08-04 Original Document



www.thermofisher.com

Potassium Tellurite 3.5%

REF SR0030J

* Le présent mode d'emploi est destiné à être lu conjointement avec le mode d'emploi relatif au milieu base Hoyle (CM0083B).

Domaine d'application

SR0003J est un complément sélectif à utiliser avec le milieu base Hoyle (déshydraté) (CM0083B) pour l'isolement et la différenciation de *Corynebacterium diphtheriae* à partir d'échantillons de plaies et de voies respiratoires supérieures, notamment les écouvillons de gorge et de nez.

Le produit est destiné à un usage professionnel uniquement, n'est pas automatisé et n'est pas un diagnostic compagnon.

Résumé et description

Les espèces *Corynebacterium* sont des bâtonnets à Gram positif avec un aspect d'extrémité en massue et existent individuellement ou en paires. Les corynebactéries médicalement pertinentes comprennent *Corynebacterium pseudotuberculosis*, *Corynebacterium ulcerans* et la forme la plus toxigène *Corynebacterium diphtheriae*, avec quatre biovars : *intermedius*, *gravis*, *mitis* et *belfanti*¹. La diphtérie est une maladie causée par les souches toxigènes de *C. diphtheriae* et moins souvent par *C. ulcerans*². Le seul réservoir connu de *C. diphtheriae* est l'homme³.

La manifestation clinique de la diphtérie peut être décrite comme une diphtérie respiratoire qui se manifeste de manière invasive ou non invasive dans les voies respiratoires ou une diphtérie cutanée associée à des lésions cutanées. Le mode de transmission de la maladie peut se faire soit par des gouttelettes provenant de patients symptomatiques ou asymptomatiques, soit par du lait ou des produits laitiers contaminés. L'apparition des symptômes commence par une légère fièvre et un mal de gorge en deux à quatre jours. Au cours de l'infection, l'exotoxine induit la formation d'un coagulum nécrotique blanc brillant dans les parois des voies respiratoires qui se transforme ensuite en taches grises à noires d'aspect conduisant à la suffocation et à l'aspiration de la membrane qui conduit finalement à la mort chez les patients de tous les groupes d'âge⁴. Cependant, avec une administration opportune d'antitoxine et une thérapie antimicrobienne, la maladie peut être traitée complètement⁵.

Principe de la méthode

Le tellurite de potassium agit comme un agent sélectif et possède une activité inhibitrice contre la plupart des bactéries à Gram positif et à Gram négatif, à l'exception de l'espèce *Corynebacterium*. *C. diphtheriae* réduit le tellurite de potassium en tellurium et produit des colonies de couleur gris noir.

Formule typique

Tellurite de potassium

2 mL

Matériel fourni

SR0030J : 10 flacons de 2 ml. Le rendement dépendra de la formulation.

Thermo

SCIENTIFIC

Matériel requis, mais non fourni

- Base de milieu Hoyle (déshydratée) code produit CM0083B
- Sang de cheval floconné SR0048C

Conservation

- Conserver le produit dans son emballage d'origine entre 2 et 8 °C.
- Garder le récipient hermétiquement fermé.
- Le produit peut être utilisé jusqu'à la date de péremption indiquée sur l'étiquette.
- Protéger de l'humidité.
- Conserver à l'abri de la lumière.
- Laisser le produit se reconstituer à température ambiante avant utilisation.

Avertissements et précautions

- N'UTILISEZ PAS LES SUPPLÉMENTS SI VOUS ÊTES ENCEINTE OU ENVISAGEZ UNE GROSSESSE.
- Chaque flacon est à usage unique. Ne pas réutiliser.
- Pour usage diagnostique in vitro uniquement.
- Usage exclusivement réservé à des professionnels.
- Inspecter l'emballage du produit avant la première utilisation.
- Ne pas utiliser le produit si l'emballage ou les flacons présentent des dommages visibles.
- Ne pas utiliser au-delà de la date de péremption indiquée.
- Ne pas utiliser le produit s'il présente des signes de contamination.
- Ne pas utiliser le produit si sa couleur a changé ou s'il présente d'autres signes de détérioration.
- Il relève de la responsabilité de chaque laboratoire de gérer les déchets produits conformément à leur nature et à leur degré de danger et de les traiter ou de les éliminer conformément aux réglementations fédérales, nationales et locales applicables. Les instructions doivent être lues et respectées scrupuleusement. Cela inclut l'élimination des réactifs utilisés ou inutilisés ainsi que de tout autre matériel jetable contaminé après les procédures impliquant des produits infectieux ou potentiellement infectieux.

Consulter la fiche de données de sécurité du matériel pour savoir comment manipuler et éliminer le produit en toute sécurité à l'adresse (www.thermofisher.com).

Incidents graves

Tout incident grave survenu en relation avec le dispositif doit être signalé au fabricant et à l'autorité réglementaire compétente dont dépendent l'utilisateur et/ou le patient.

Prélèvement, manipulation et stockage des échantillons

L'échantillon doit être collecté et manipulé conformément aux directives locales recommandées, telles que les UK Standards for Microbiology Investigations (UK SMI) ID 2 et Q 5.

Procédure

- Suspender 4 g de milieu base Hoyle (CM0083B) dans 100 ml d'eau distillée et porter à ébullition pour dissoudre complètement.
- Stériliser à l'autoclave à 121 °C pendant 15 minutes.
- Refroidir jusqu'à 50 °C.










- Ajouter 5 ml de sang de cheval fumé (SR0048C) et 1 ml de tellurite de potassium (SR0030J).
- Bien mélanger et verser dans des récipients stériles.







Se reporter au mode d'emploi de CM0083B. Les modes d'emploi sont disponibles sur www.thermofisher.com

Bibliographie

1. Public Health England. 2014 'Identification of Corynebacterium species' ID_2314.1 <https://www.gov.uk/government/publications/smi-id-2-identification-of-corynebacterium-species>
2. Mokrousov, Igor. "Corynebacterium diphtheriae: genome diversity, population structure and genotyping perspectives." Infection, Genetics and Evolution 9, no. 1 (2009): 1-15.
3. Rajamani Sekar, S. K., B. Veeraraghavan, S. Anandan, N. K. Devanga Ragupathi, L. Sangal, and S. Joshi. "Strengthening the laboratory diagnosis of pathogenic Corynebacterium species in the Vaccine era." Letters in applied microbiology 65, no. 5 (2017): 354-365.
4. Zakikhany, Katherina, and Androulla Efstratiou. "Diphtheria in Europe: current problems and new challenges." Future microbiology 7, no. 5 (2012): 595-607.
5. Sharma, Naresh Chand, Androulla Efstratiou, Igor Mokrousov, Ankur Mutreja, Bhabatosh Das, and Thandavarayan Ramamurthy. "Diphtheria (Primer)." Nature Reviews: Disease Primers (2019).

Symboles

Symbole	Définition
	Référence catalogue
	Dispositif médical de diagnostic in vitro
	Code de lot
	Limite de température
	Date limite d'utilisation
	Tenir à l'abri de la lumière directe du soleil
	Ne pas réutiliser
	Se référer aux instructions d'utilisation ou consulter les instructions d'utilisation électroniques
	Contenu suffisant pour <n> tests

	Ne pas utiliser si l'emballage est endommagé et consulter les instructions d'utilisation
	Fabricant
	Représentant agréé pour la Communauté européenne/ Union européenne
	Évaluation de la conformité européenne
	Évaluation de la conformité pour le Royaume-Uni
	Identifiant unique du dispositif
Fabriqué au Royaume-Uni	Fabriqué au Royaume-Uni

© 2022 Thermo Fisher Scientific Inc. Tous droits réservés. Toutes les autres marques sont la propriété de Thermo Fisher Scientific Inc. et de ses filiales.



Pour une assistance technique, contacter le distributeur local.

Informations de révision

Version	Modifications introduites
1.0	2022-08-04 Document original



www.thermofisher.com

Potassium Tellurite 3.5%

REF SR0030J

* Diese Gebrauchsanweisung sollte in Verbindung mit der Gebrauchsanweisung für Hoyle-Medium-Basis (CM0083B) gelesen werden.

Verwendungszweck

SR0003J ist ein Selektiv-Supplement für die Verwendung mit Hoyle-Medium-Basis (dehydriert) (CM0083B) zur Isolierung und Differenzierung von *Corynebacterium diphtheriae* aus Wundproben und Proben der oberen Atemwege, einschließlich Rachen- und Nasenabstrichen.

Das Produkt ist nur für den professionellen Gebrauch bestimmt, es ist nicht automatisiert und dient auch nicht als Begleitdiagnose.

Zusammenfassung und Erläuterung

Corynebacterium-Spezies sind grampositive Stäbchen mit keulenförmigem Aussehen und kommen einzeln oder paarweise vor. Zu den medizinisch relevanten Corynebakterien gehören *Corynebacterium pseudotuberculosis*, *Corynebacterium ulcerans* und die höchst toxische Form *Corynebacterium diphtheriae*, mit vier Biovarien: *intermedius*, *gravis*, *mitis* und *belfanti*.¹ Diphtherie ist eine Krankheit, die durch die toxigenen Stämme von *C. diphtheriae* und seltener durch *C. ulcerans* verursacht wird.² Das einzige bekannte Reservoir von *C. diphtheriae* ist der Mensch.³

Die klinische Manifestation der Diphtherie kann als respiratorische Diphtherie beschrieben werden, die invasiv oder nicht-invasiv in den Atemwegen auftritt, oder als kutane Diphtherie, die mit Hautläsionen einhergeht. Die Übertragung der Krankheit kann entweder durch Tröpfchen von symptomatischen oder asymptomatischen Patienten oder durch kontaminierte Milch oder Milchprodukte erfolgen. Der Beginn der Symptome beginnt mit leichtem Fieber und Halsschmerzen innerhalb von zwei bis vier Tagen. Während der Infektion induziert das Exotoxin die Bildung eines weißen, glänzenden, nekrotischen Koagulums in den Wänden der Atemwege, das sich dann in graue bis schwarze Flecken verwandelt, was zum Ersticken und zur Aspiration von Membranen führt, was schließlich bei Patienten aller Altersgruppen zum Tod führt.⁴ Mit einer rechtzeitigen Verabreichung von Antitoxin und einer antimikrobiellen Therapie kann die Krankheit jedoch vollständig behandelt werden.⁵

Das Prinzip der Methode

Kaliumtellurit wirkt Selektiv und hat eine hemmende Wirkung auf die meisten gram-positiven und gram-negativen Bakterien mit Ausnahme von *Corynebacterium*-Spezies. *C. diphtheriae* reduziert Kaliumtellurit zu Tellur und bildet graue, schwarz gefärbte Kolonien.

Typische Formel

Kaliumtellurit 2 ml

Mitgeliefertes Material

SR0030J: 10x 2-ml-Fläschchen. Die Ausbeute hängt von der Formulierung ab.

Thermo
SCIENTIFIC

Erforderliche, aber nicht mitgelieferte Materialien

- Hoyle-Medium-Basis (dehydriert) Produktcode CM0083B
- Getränktes Pferdeblut SR0048C

Lagerung

- Lagern Sie das Produkt in der Originalverpackung zwischen 2 °C und 8 °C.
- Behälter dicht geschlossen halten.
- Das Produkt kann bis zu dem auf dem Etikett angegebenen Verfallsdatum verwendet werden.
- Vor Feuchtigkeit schützen.
- Vor Licht geschützt aufbewahren.
- Lassen Sie das rekonstituierte Produkt vor der Verwendung auf Raumtemperatur kommen.

Warnungen und Vorsichtsmaßnahmen

- VERWENDEN SIE KEINE SUPPLEMENTS, WENN SIE SCHWANGER SIND ODER EINE SCHWANGERSCHAFT IN BETRACHT ZIEHEN.
- Jedes Fläschchen ist zum einmaligen Gebrauch bestimmt. Nicht wiederverwenden.
- Nur für die In-vitro-Diagnostik geeignet.
- Nur für den professionellen Gebrauch.
- Überprüfen Sie die Produktverpackung vor dem ersten Gebrauch.
- Verwenden Sie das Produkt nicht, wenn es sichtbare Schäden an der Verpackung oder den Fläschchen aufweist.
- Verwenden Sie das Produkt nicht nach Ablauf des angegebenen Verfallsdatums.
- Verwenden Sie das Produkt nicht, wenn es Anzeichen von Verschmutzung aufweist.
- Verwenden Sie das Produkt nicht, wenn sich die Farbe verändert hat oder andere Anzeichen einer Verschlechterung vorliegen.
- Es liegt in der Verantwortung jedes Labors, die anfallenden Abfälle entsprechend ihrer Art und ihres Gefährdungsgrades zu behandeln und sie in Übereinstimmung mit den auf Bundes-, Landes- und lokaler Ebene geltenden Vorschriften zu behandeln oder zu entsorgen. Die Gebrauchsanweisung sollte sorgfältig gelesen und befolgt werden. Dazu gehört auch die Entsorgung gebrauchter oder unbenutzter Reagenzien sowie aller anderen kontaminierten Einwegmaterialien gemäß den Verfahren für infektiöse oder potenziell infektiöse Produkte.

Beachten Sie das Sicherheitsdatenblatt (SDB) für die sichere Handhabung und Entsorgung des Produkts (www.thermofisher.com).

Schwere Zwischenfälle

Jeder schwerwiegende Zwischenfall im Zusammenhang mit dem Produkt ist dem Hersteller und der zuständigen Aufsichtsbehörde, in deren Zuständigkeitsbereich der Anwender und/oder der Patient niedergelassen ist/sind, zu melden.

Entnahme, Handhabung und Lagerung von Proben

Die Probenentnahme und -behandlung sollte gemäß den vor Ort empfohlenen Richtlinien erfolgen, wie z. B. den UK Standards for Microbiology Investigations (UK SMI) ID 2 und Q 5.

Verfahren

- Suspendieren Sie 4 g Hoyle-Medium-Basis (CM0083B) in 100 ml destilliertem Wasser und

bringen Sie es zum Kochen, um es vollständig aufzulösen.









- Sterilisieren Sie es bei 121 °C für 15 Minuten im Autoklaven.
- Auf 50 °C abkühlen.
- Geben Sie 5 ml getränktes Pferdeblut (SR0048C) und 1 ml Kaliumtellurit (SR0030J) hinzu.
- Gut mischen und in sterile Behälter füllen.




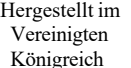
Siehe die Gebrauchsanweisung für CM0083B. Gebrauchsanweisungen finden Sie [unter www.thermofisher.com](http://www.thermofisher.com).

Bibliographie

1. Public Health England. 2014 'Identification of Corynebacterium species' ID_2314.1 <https://www.gov.uk/government/publications/smi-id-2-identification-of-corynebacterium-species>
2. Mokrousov, Igor. "Corynebacterium diphtheriae: genome diversity, population structure and genotyping perspectives." Infection, Genetics and Evolution 9, no. 1 (2009): 1-15.
3. Rajamani Sekar, S. K., B. Veeraraghavan, S. Anandan, N. K. Devanga Ragupathi, L. Sangal, and S. Joshi. "Strengthening the laboratory diagnosis of pathogenic Corynebacterium species in the Vaccine era." Letters in applied microbiology 65, no. 5 (2017): 354-365.
4. Zakikhany, Katherina, and Androulla Efstratiou. "Diphtheria in Europe: current problems and new challenges." Future microbiology 7, no. 5 (2012): 595-607.
5. Sharma, Naresh Chand, Androulla Efstratiou, Igor Mokrousov, Ankur Mutreja, Bhabatosh Das, and Thandavarayan Ramamurthy. "Diphtheria (Primer)." Nature Reviews: Disease Primers (2019).

Symbollegende

Symbol	Definition
	Katalognummer
	Medizinprodukt zum In-vitro-Diagnostikum
	Chargencode
	Temperaturgrenze
	Haltbarkeitsdatum
	Vom Sonnenlicht fernhalten
	Nicht wiederverwenden
	Gebrauchsanweisung oder elektronische Gebrauchsanweisung konsultieren

	Enthält ausreichend für <n> Tests
	Nicht verwenden, wenn die Verpackung beschädigt ist und die Gebrauchsanweisung konsultieren
	Hersteller
	Bevollmächtigter Vertreter in der Europäischen Gemeinschaft/Europäischen Union
	Europäische Konformitätsbewertung
	Konformitätsbewertung des Vereinigten Königreichs
	Eindeutige Kennung des Geräts
	Hergestellt im Vereinigten Königreich

© 2022 Thermo Fisher Scientific Inc. Alle Rechte vorbehalten.

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



Für technische Unterstützung wenden Sie sich bitte an Ihren örtlichen Händler.

Informationen zur Revision

Version	Eingeführte Modifikationen
1.0	2022-08-04 Original-Dokument



www.thermofisher.com

Potassium Tellurite 3.5%

[REF] SR0030J

* Aceste instrucțiuni de utilizare trebuie citite împreună cu instrucțiunile de utilizare pentru Hoyle Medium Base (CM0083B).

Utilizare prevăzută

SR0030J este un supliment selectiv conceput pentru a fi utilizat împreună cu Hoyle Medium Base (Dehydrated) (CM0083B) pentru izolarea și diferențierea *Corynebacterium diphtheriae* din probe prelevate din plăgi și din sistemul respirator, inclusiv de pe tamponane nazale și orofaringiene.

Dispozitivul este exclusiv de uz profesional, nu este automatizat și nici nu constituie un diagnostic complementar.

Rezumat și explicație

Speciile *Corynebacterium* sunt bacili gram-pozitivi cu capete umflate, care pot apărea individual sau în perechi. Corinebacteriile relevante din punct de vedere medical includ *Corynebacterium pseudotuberculosis*, *Corynebacterium ulcerans* și cea mai toxigenă formă, *Corynebacterium diphtheriae*, cu patru biovari: *intermedius*, *gravis*, *mitis* și *belfanti*.¹ Difteria este o boală cauzată de tulpinile toxice ale *C. diphtheriae* și mai rar de *C. ulcerans*.² Singurul rezervor cunoscut de *C. diphtheriae* sunt oamenii.³

Manifestările clinice ale difteriei sunt reprezentate de difterie respiratorie, care apare în tractul respirator și poate fi invazivă sau neinvazivă, sau difterie cutanată, asociată cu leziuni cutanate. Boala se poate transmite prin picături, de la pacienți simptomatici sau asimptomatici ori din lapte sau produse lactate contaminate. Debutul simptomelor este caracterizat de febră ușoară și dureri în gât, în termen de două-patru zile. În timpul infecției, exotoxina induce formarea unui coagul necrotic alb lucios pe pereții căilor respiratorii, care apoi se transformă în pete cenușii până la negre, care provoacă sufocarea și aspirarea membranei, și, în cele din urmă, duce la deces la pacienții din toate grupele de vârstă.⁴ Cu toate acestea, boala poate fi tratată complet prin administrarea la timp a antitoxinelor și a terapiei antimicrobiene.⁵

Principiul metodei

Teluritul de potasiu acționează ca agent selectiv și are activitate inhibitoare împotriva majorității bacteriilor gram-pozitive și gram-negative, cu excepția speciei *Corynebacterium*. *C. diphtheriae* reduce teluritul de potasiu la telurii și produce colonii de culoare gri negru.

Formula tipică

Telurit de potasiu 2 ml

Materiale furnizate

SR0030J: 10 flacoane de 2 ml. Randamentul depinde de formulă.

Materiale necesare, dar nefurnizate

- Hoyle Medium Base (dehydrated) cod produs CM0083B
- Sânge de cal hemolizat SR0048C

Depozitare

- Depozitați produsul în ambalajul original, la temperaturi între 2 °C – 8 °C.
- Păstrați recipientul închis etanș.
- Produsul poate fi utilizat până la data de expirare înscrisă pe etichetă.
- A se proteja de umiditate.
- A se păstra departe de surse de lumină.
- Lăsați produsul reconstituit să ajungă la temperatura camerei înainte de utilizare.

Avertismente și mijloace de precauție

- NU UTILIZAȚI SUPLIMENTE DACĂ SUNTEȚI SAU DORIȚI SĂ RĂMÂNEȚI ÎNSĂRCINATĂ.
- Fiecare flacon este de unică folosință. A nu se reutiliza.
- Exclusiv pentru diagnosticarea in vitro.
- Exclusiv de uz profesional.
- Inspectați ambalajul produsului înainte de prima utilizare.
- Nu utilizați produsul dacă ambalajul sau flacoanele sunt deteriorate vizibil.
- A nu se utiliza produsul după data de expirare specificată.
- Nu utilizați dispozitivul dacă există semne de contaminare.
- Nu utilizați dispozitivul dacă culoarea este modificată sau dacă există alte semne de deteriorare.
- Este responsabilitatea fiecărui laborator să gestioneze deșeurile produse, în funcție de natura și gradul de pericol, și de a le trata sau elimina în conformitate cu reglementările aplicabile federale, statale și locale. Instrucțiunile trebuie citite și urmate cu atenție. Aceasta include eliminarea reactivilor utilizați sau neutilizați, precum și a oricărui alt material contaminat de unică folosință, urmând procedurile pentru produsele infecțioase sau potențial infecțioase.

Consultați Fișa cu date de securitate a materialelor (FDSM) pentru manipularea și eliminarea în siguranță a produsului (www.thermofisher.com).

Incidente grave

Orice incident grav survenit în legătură cu dispozitivul va fi raportat producătorului și autorității de reglementare relevante a Statului Membru în care utilizatorul și/sau pacientul își are reședința.

Recoltarea, manipularea și depozitarea probelor

Probele trebuie recoltate și manipulate cu respectarea orientărilor locale recomandate, precum UK Standards for Microbiology Investigations (UK SMI) ID 2 și Q 5.

Procedură













- Suspendedați 4 g de Hoyle Medium Base (CM0083B) în 100 ml de apă distilată și aduceți la temperatura de fierbere pentru a se dizolva complet.
- Se sterilizează prin autoclavare la 121 °C timp de 15 minute.
- Se răcește la 50 °C.
- Adăugați 5 ml de sânge de cal hemolizat (SR0048C) și 1 ml telurit de potasiu (SR0030J).
- Se amestecă bine și se toarnă în recipiente sterile.




Consultați instrucțiunile de utilizare pentru CM0083B. Instrucțiunile de utilizare pot fi găsite la www.thermofisher.com

Bibliografie

1. Public Health England. 2014 'Identification of *Corynebacterium* species' ID_2314.1 <https://www.gov.uk/government/publications/smi-id-2-identification-of-corynebacterium-species>
2. Mokrousov, Igor. "Corynebacterium diphtheriae: genome diversity, population structure and genotyping perspectives." *Infection, Genetics and Evolution* 9, no. 1 (2009): 1-15.
3. Rajamani Sekar, S. K., B. Veeraraghavan, S. Anandan, N. K. Devanga Ragupathi, L. Sangal, and S. Joshi. "Strengthening the laboratory diagnosis of pathogenic *Corynebacterium* species in the Vaccine era." *Letters in applied microbiology* 65, no. 5 (2017): 354-365.
4. Zakikhany, Katherina, and Androulla Efstratiou. "Diphtheria in Europe: current problems and new challenges." *Future microbiology* 7, no. 5 (2012): 595-607.
5. Sharma, Naresh Chand, Androulla Efstratiou, Igor Mokrousov, Ankur Mutreja, Bhabatosh Das, and Thandavarayan Ramamurthy. "Diphtheria (Primer)." *Nature Reviews: Disease Primers* (2019).

Legenda simbolurilor

Simbol	Definiție
	Număr de catalog
	Dispozitiv medical pentru diagnosticarea in vitro
	Codul lotului
	Limita de temperatură
	Data expirării
	A se păstra ferit de expunere la soare
	A nu se reutiliza
	Consultați instrucțiunile de utilizare sau consultați instrucțiunile de utilizare electronice
	Conține o cantitate suficientă pentru <n> teste
	A nu se utiliza dacă ambalajul este deteriorat și consultați instrucțiunile de utilizare
	Producător
	Reprezentant autorizat în Comunitatea Europeană/ Uniunea Europeană

	Marcajul de conformitate europeană
	Marcajul de conformitate pentru Regatul Unit
	Identificatorul unic al dispozitivului
Made in the United Kingdom	Fabricat în Regatul Unit

©2022 Thermo Fisher Scientific Inc. Toate drepturile rezervate.

Toate celelalte mărci comerciale aparțin Thermo Fisher Scientific Inc. și subsidiarelor acesteia.



Oxoid Limited, Wade Road, Basingstoke, Hampshire, RG24 8PW, Marea Britanie

Pentru asistență tehnică, vă rugăm să contactați distribuitorul local.

Informații privind reviziile

Versiunea	Modificări introduse
1.0	2022-08-04 Document original



www.thermofisher.com

Potassium Tellurite 3.5%

[REF] SR0030J

* Este documento de instrucciones de uso está diseñado para utilizarlo junto con las instrucciones de uso de Hoyle Medium Base (CM0083B).

Uso previsto

SR0003J es un suplemento selectivo para utilizarse con Hoyle Medium Base (Dehydrated) (CM0083B) para el aislamiento y la diferenciación de *Corynebacterium diphtheriae* a partir de muestras de heridas y de las vías respiratorias altas, incluidos hisopos nasales y de garganta.

El dispositivo es exclusivamente para uso profesional, no está automatizado y no es una prueba diagnóstica acompañante.

Resumen y explicación

Las especies de *Corynebacterium* son bacilos grampositivos con extremos en forma de acropaquia y existen en solitario o en parejas. Las corinebacterias médicamente relevantes incluyen *Corynebacterium pseudotuberculosis*, *Corynebacterium ulcerans* y la forma más toxígena, *Corynebacterium diphtheriae*, con cuatro biovariedades: *intermedius*, *gravis*, *mitis* y *belfanti*¹. La difteria es una enfermedad causada por las cepas toxígenas de *C. diphtheriae* y, con menor frecuencia, por *C. ulcerans*². El único reservorio conocido de *C. diphtheriae* son los humanos³.

La manifestación clínica de la difteria se puede describir como difteria respiratoria que se presenta como invasiva o no invasiva en las vías respiratorias, o difteria cutánea asociada a lesiones en la piel. El modo de transmisión de la enfermedad puede ser a través de gotículas de pacientes sintomáticos o asintomáticos o de leche o productos lácteos contaminados. La aparición de los síntomas comienza con fiebre leve y garganta irritada en dos a cuatro días. Durante la infección, la exotoxina induce la formación de un coágulo necrótico blanco y brillante en las paredes de las vías respiratorias que luego se convierte en parches de color gris a negro que conducen a la sofocación y la aspiración de la membrana, lo que finalmente conduce a la muerte en pacientes de todos los grupos de edad⁴. Sin embargo, con la administración oportuna de antitoxina y terapia antimicrobiana, la enfermedad se puede tratar por completo⁵.

Principio del método

El telurito de potasio actúa como agente selectivo y tiene actividad inhibitoria contra la mayoría de las bacterias grampositivas y gramnegativas, excepto las especies de *Corynebacterium*. *C. diphtheriae* reduce el telurito de potasio a telurio y produce colonias de color negro grisáceo.

Fórmula típica

Telurito de potasio 2 ml

Materiales suministrados

SR0030J: 10 viales de 2 ml. El rendimiento dependerá de la formulación.

Thermo
SCIENTIFIC

Materiales necesarios pero no suministrados

- Hoyle Medium Base (Dehydrated), código de producto CM0083B
- Laked Horse blood SR0048C

Almacenamiento

- Almacenar el producto en su envase original entre 2 °C y 8 °C.
- Mantener el recipiente cerrado herméticamente.
- El producto se puede utilizar hasta la fecha de caducidad indicada en la etiqueta.
- Protegerlo de la humedad.
- Almacenar protegido de la luz.
- Dejar que el producto reconstituido se temple a temperatura ambiente antes de usarlo.

Advertencias y precauciones

- NO UTILIZAR SUPLEMENTOS SI ESTÁ EMBARAZADA O PREVÉ QUEDARSE EMBARAZADA.
- Cada vial es de un solo uso. No reutilizar.
- Para uso diagnóstico in vitro exclusivamente.
- Para uso profesional exclusivamente.
- Inspeccionar el envase del producto antes del primer uso.
- No utilizar el producto si hay daños visibles en el envase o los viales.
- No utilizar el producto más allá de la fecha de caducidad indicada.
- No utilizar el dispositivo si presenta signos de contaminación.
- No utilizar el dispositivo si el color ha cambiado o hay otros signos de deterioro.
- Es responsabilidad de cada laboratorio manejar los residuos generados de acuerdo con su naturaleza y grado de peligrosidad y tratarlos o eliminarlos según los reglamentos federales, estatales y locales aplicables. Es necesario leer las instrucciones y seguirlas atentamente. Esto incluye la eliminación de reactivos usados o sin usar, así como cualquier otro material desechable contaminado según los procedimientos para productos infecciosos o potencialmente infecciosos.

Consulte las instrucciones de manipulación y eliminación segura del producto en la Hoja de datos de seguridad del material (SDS) (www.thermofisher.com).

Incidentes graves

Cualquier incidente grave que se produzca en relación con el producto se debe notificar al fabricante y a la autoridad reguladora pertinente donde residan el usuario o el paciente.

Recogida, manipulación y almacenamiento de muestras

Es necesario recoger y manipular las muestras según las directrices locales recomendadas, como los Estándares para investigaciones de microbiología del Reino Unido (UK SMI) ID 2 y Q 5.

Procedimiento

- Suspensa 4 g de Hoyle Medium Base (CM0083B) en 100 ml de agua destilada y lleve a ebullición para que se disuelva por completo.
- Esterilice en autoclave a 121 °C durante 15 minutos.
- Enfríe a 50 °C.

- Añada 5 ml de Laked Horse Blood (SR0048C) y 1 ml de Potassium Tellurite (SR0030J).
- Mezcle bien y vierta en recipientes estériles.

Consulte las instrucciones de uso de CM0083B. Puede encontrar las instrucciones de uso en www.thermofisher.com

Bibliografía

1. Public Health England. 2014 'Identification of Corynebacterium species' ID_2314.1 <https://www.gov.uk/government/publications/smi-id-2-identification-of-corynebacterium-species>
2. Mokrousov, Igor. "Corynebacterium diphtheriae: genome diversity, population structure and genotyping perspectives." Infection, Genetics and Evolution 9, no. 1 (2009): 1-15.
3. Rajamani Sekar, S. K., B. Veeraraghavan, S. Anandan, N. K. Devanga Ragupathi, L. Sangal, and S. Joshi. "Strengthening the laboratory diagnosis of pathogenic Corynebacterium species in the Vaccine era." Letters in applied microbiology 65, no. 5 (2017): 354-365.
4. Zakikhany, Katherina, and Androulla Efstratiou. "Diphtheria in Europe: current problems and new challenges." Future microbiology 7, no. 5 (2012): 595-607.
5. Sharma, Naresh Chand, Androulla Efstratiou, Igor Mokrousov, Ankur Mutreja, Bhabatosh Das, and Thandavarayan Ramamurthy. "Diphtheria (Primer)." Nature Reviews: Disease Primers (2019).

Leyenda de símbolos

Símbolo	Definición
	Número de catálogo
	Producto sanitario para diagnóstico in vitro
	Código de lote
	Límite de temperatura
	Fecha de caducidad
	Mantener alejado de la luz solar
	No reutilizar
	Consultar las instrucciones de uso o las instrucciones de uso electrónicas
	Contiene la cantidad suficiente para <n> pruebas
	No utilizar si el paquete está dañado y consultar las instrucciones de uso

	Fabricante
	Representante autorizado en la Comunidad Europea/ Unión Europea
	Evaluación de conformidad europea
	Evaluación de conformidad para el Reino Unido
	Identificador único de dispositivo
Hecho en el Reino Unido	Hecho en el Reino Unido

©2022 Thermo Fisher Scientific Inc. Reservados todos los derechos.

Todas las demás marcas comerciales son propiedad de Thermo Fisher Scientific Inc. y sus filiales.



Oxoid Limited, Wade Road, Basingstoke, Hampshire, RG24 8PW, Reino Unido

Para obtener asistencia técnica, póngase en contacto con su distribuidor local.

Información de revisiones

Versión	Modificaciones introducidas
1.0	2022-08-04 Documento original