



www.thermofisher.com

Agar
Physical Appearance

Colour	Orange
Clarity	Clear
Fill weight	3.0 – 3.3g
pH	6.8 ± 0.2

Materials Provided

BO0337B: 24 x Urea Agar Slope bottles
EB0337B: 200 x Urea Agar Slope Bottles

Each bottle should only be used once.

Materials Required but Not Supplied

- Inoculating loops
- Swabs
- Collection containers
- Incubators
- Quality control organisms

Storage

- Store product in its original packaging at 2–10°C until used.
- The product may be used until the expiry date stated on the label.
- Store away from light.
- Allow product to equilibrate to room temperature before use.
- Do not incubate prior to use.

Warnings and Precautions

- 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 plates.
- 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

Specimens should be collected and handled following local recommended guidelines, such as the UK Standards for Microbiology Investigations (UK SMI) ID 16, ID 20, ID 21, ID 24,

Procedure

- Allow product to equilibrate to room temperature.

Urea Agar Slope



REF BO0337B & EB0337B

Intended Use

Urea Agar Slope (BO0337B and EB0037B) devices are non-selective media intended for use in culture media. Urea Agar Slope (BO0337B and EB0037B) devices are intended to be used to detect rapid urease activity of the *Proteae* and non-rapid urease activity of some *Enterobacteriaceae* in faecal and urine samples.

The devices are intended to be used in a diagnostic workflow to aid clinicians in determining potential treatment options for patients suspected of having *Enterobacteriaceae* infections.

The devices are for professional use only, are not automated and nor are they companion diagnostics."

Summary and Explanation

Urinary tract infections, often referred to as UTIs, are common infections that occur when bacteria invade the urinary tract by entering via the urethra. These infections can impact different parts of the urinary tract, but the most prevalent is cystitis (a bladder infection). UTIs occur more frequently in women, as the urethra is shorter, and therefore closer to the rectum. This increases the likelihood of infection as it is easier for bacteria to enter the urinary tract. Kidney infections are another form of UTI. These are more serious than cystitis, but far less common. The symptoms of cystitis can include frequent urination, or needing to urinate whilst having an empty bladder, a burning sensation or pain whilst urinating, blood in the urine and pressure or cramps in the lower abdominal area. The symptoms of kidney infections are very different. These can include chills, a fever, nausea and vomiting and pain in the lower back¹. There are many species of bacteria that cause UTIs, but the three covered in this file are *Escherichia coli*, *Staphylococcus aureus* and *Proteus mirabilis*.

Due to the severity of UTIs in a vast number of patients, it is therefore highly important to be able to isolate and identify common UTI-causing microorganisms from urine samples. Early diagnosis is vital in the prevention and reduction of urinary tract infections.

Principle of Method

Urea Agar is a lightly buffered medium containing urea and phenol red, a pH indicator. When organisms utilize urea, ammonia is formed which turns the medium alkaline. The indicator, phenol red, changes the medium colour from pale-yellow to pink red in an alkaline environment. Gelatin peptone promotes rapid growth of enteric gram-negative bacilli, permitting a decrease in incubation time. Dextrose stimulates urease activity in those organisms which hydrolyze urea slowly.

Typical Formula

	<u>grams per litre</u>
Peptone	1.0
Glucose	1.0
Sodium chloride	5.0
Disodium phosphate	1.2
Potassium dihydrogen phosphate	0.8
Phenol red	0.012

- Inoculate and streak the specimen onto the medium.
- or alternatively using aseptic technique pick a single well-isolated colony with a sterile inoculating stab needle and stab the needle several times through the agar to the bottom of the vial or tube
- Replace the cap and tighten loosely when incubating to allow gas exchange
- Incubate the stabbed agar / slant 18 - 24 hours at 37 ± 2°C.
- Visually inspect bottles to assess colony growth and colour under good lighting.

Interpretation

Urease negative, no colour change/ yellow medium indicates *Escherichia coli*.

Urease positive with pink slopes indicates *Proteus mirabilis* or *Proteus hauseri*.

Weak urease positive with pale pink slopes indicates *Klebsiella pneumoniae*.

Quality Control

It is the responsibility of the user to perform Quality Control testing taking into account the intended use of the medium, and in accordance with any local applicable regulations (frequency, number of strains, incubation temperature etc.).

The performance of this medium can be verified by testing the following reference strains.

Incubation Conditions: 18 - 24 h @ 37° ± 2°C

Positive Controls	
Inoculum level: 10,000 – 100,000 cfu	
<i>Escherichia coli</i> ATCC® 25922™	Urease negative, no colour change/ yellow medium
<i>Proteus mirabilis</i> NCTC 10975	Urease positive, pink slope
<i>Proteus hauseri</i> ATCC® 13315™	Urease positive, pink slope
<i>Klebsiella pneumoniae</i> ATCC® 13883™	Weak urease positive, pale pink slope

Limitations

The alkaline reaction produced in this medium after prolonged incubation may not be caused by urease activity. Check using medium without urea.

Do not heat or reheat the medium because urea decomposes very easily.

For the detection of urease-positive *Proteae* the reaction must be read within the first 2-6 hours of incubation.

Performance Characteristics

Accuracy has been demonstrated through review of the QC data. Correct detection of Enterobacteriaceae is confirmed by the inclusion of a well-characterised isolate in the QC processes performed as part of the manufacture of each batch of the devices. The precision of Urea Agar Slope (BO0337B) was demonstrated by an overall pass rate of 80% obtained for the product over ten months of testing (18.SEP.2021– 25.JUL.2022; 10 batches). The precision of Urea Agar Slope (EB0337B) was demonstrated by an overall pass rate of 90% obtained for the product over nine months of testing (13.SEP.2021– 13.JUN.2022; 10 batches). An overall pass-rate of 80% (based on data from the most recent 10 batches) provides sufficient data on accuracy considering the low risk of the device and it is not reflective of the product that is released for clinical use. With the normal variation in the manufacturing process, a small
















number of failures may occur, however, there is no risk to the end user as batches will be captured in QC if out of specification. This shows that the performance is reproducible, thus meeting the specifications set out in the Performance Evaluation Plan.


Urea Agar Slope (BO0337B & EB0337B) devices are tested in-house as part of the QC process since 1997 (BO0337B) and 1998 (EB0337B). For target organisms, when using 10⁴ - 10⁵ cfu inoculum of *Escherichia coli*, *Proteus mirabilis* or *Klebsiella pneumoniae* and incubating the device at 37° ± 2°C for 18-24 hours, the user can expect urease reaction as listed in this document.

Bibliography

1. Centers for Disease Control and Prevention. 2021. 'Urinary Tract Infection'. <https://www.cdc.gov/antibiotic-use/uti.html>.

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

	Importer - To indicate the entity importing the medical device into the locale. Applicable to the European Union
Made in the United Kingdom	Made in the United Kingdom



©2022 Thermo Fisher Scientific Inc. All rights reserved. ATCC and ATCC catalogue marks are a trademark of American Type Culture Collection. NCTC and NCTC catalogue marks are a trademark of National Collection of Type Cultures. All other trademarks are the property of Thermo Fisher Scientific Inc. and its subsidiaries.



Oxoid Limited, Wade Road, Basingstoke, Hampshire, RG24 8PW, United Kingdom



Peptone	1.0
Glucose	1.0
Sodium chloride	5.0
Disodium phosphate	1.2
Potassium dihydrogen phosphate	0.8
Phenol red	0.012

Revision Information

Version	Date of modifications introduced
2.0	2023-11-21



www.thermofisher.com

Urea Agar Slope



REF BO0337B & EB0337B

Intended Use

Urea Agar Slope (BO0337B and EB0037B) devices are non-selective media intended for use in culture media. Urea Agar Slope (BO0337B and EB0037B) devices are intended to be used to detect rapid urease activity of the *Proteae* and non-rapid urease activity of some *Enterobacteriaceae* in faecal and urine samples.

The devices are intended to be used in a diagnostic workflow to aid clinicians in determining potential treatment options for patients suspected of having *Enterobacteriaceae* infections.

The devices are for professional use only, are not automated and nor are they companion diagnostics."

Summary and Explanation

Urinary tract infections, often referred to as UTIs, are common infections that occur when bacteria invade the urinary tract by entering via the urethra. These infections can impact different parts of the urinary tract, but the most prevalent is cystitis (a bladder infection). UTIs occur more frequently in women, as the urethra is shorter, and therefore closer to the rectum. This increases the likelihood of infection as it is easier for bacteria to enter the urinary tract. Kidney infections are another form of UTI. These are more serious than cystitis, but far less common. The symptoms of cystitis can include frequent urination, or needing to urinate whilst having an empty bladder, a burning sensation or pain whilst urinating, blood in the urine and pressure or

cramps in the lower abdominal area. The symptoms of kidney infections are very different. These can include chills, a fever, nausea and vomiting and pain in the lower back¹. There are many species of bacteria that cause UTIs, but the three covered in this file are *Escherichia coli*, *Staphylococcus aureus* and *Proteus mirabilis*.

Due to the severity of UTIs in a vast number of patients, it is therefore highly important to be able to isolate and identify common UTI-causing microorganisms from urine samples. Early diagnosis is vital in the prevention and reduction of urinary tract infections.

Principle of Method

Urea Agar is a lightly buffered medium containing urea and phenol red, a pH indicator. When organisms utilize urea, ammonia is formed which turns the medium alkaline. The indicator, phenol red, changes the medium colour from pale-yellow to pink red in an alkaline environment. Gelatin peptone promotes rapid growth of enteric gram-negative bacilli, permitting a decrease in incubation time. Dextrose stimulates urease activity in those organisms which hydrolyze urea slowly.

Typical Formula

grams per litre

Agar

15.0

Physical Appearance

Colour	Orange
Clarity	Clear
Fill weight	3.0 – 3.3g
pH	6.8 ± 0.2

Materials Provided

BO0337B: 24 x Urea Agar Slope bottles
EB0337B: 200 x Urea Agar Slope Bottles

Each bottle should only be used once.

Materials Required but Not Supplied

- Inoculating loops
- Swabs
- Collection containers
- Incubators
- Quality control organisms

Storage

- Store product in its original packaging at 2–10°C until used.
- The product may be used until the expiry date stated on the label.
- Store away from light.
- Allow product to equilibrate to room temperature before use.
- Do not incubate prior to use.

Warnings and Precautions

- 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 plates.
- 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

Specimens should be collected and handled following local recommended guidelines, such as the UK Standards for Microbiology Investigations (UK SMI) ID 16, ID 20, ID 21, ID 24,

Procedure

- Allow product to equilibrate to room temperature.
- Inoculate and streak the specimen onto the

medium.

- or alternatively using aseptic technique pick a single well-isolated colony with a sterile inoculating stab needle and stab the needle several times through the agar to the bottom of the vial or tube
- Replace the cap and tighten loosely when incubating to allow gas exchange
- Incubate the stabbed agar / slant 18 - 24 hours at 37 ± 2°C.
- Visually inspect bottles to assess colony growth and colour under good lighting.

Interpretation

Urease negative, no colour change/ yellow medium indicates *Escherichia coli*.

Urease positive with pink slopes indicates *Proteus mirabilis* or *Proteus hauseri*.

Weak urease positive with pale pink slopes indicates *Klebsiella pneumoniae*.

Quality Control

It is the responsibility of the user to perform Quality Control testing taking into account the intended use of the medium, and in accordance with any local applicable regulations (frequency, number of strains, incubation temperature etc.).

The performance of this medium can be verified by testing the following reference strains.

Incubation Conditions: 18 - 24 h @ 37° ± 2°C

Positive Controls	
Inoculum level: 10,000 – 100,000 cfu	
<i>Escherichia coli</i> ATCC® 25922™	Urease negative, no colour change/ yellow medium
<i>Proteus mirabilis</i> NCTC 10975	Urease positive, pink slope
<i>Proteus hauseri</i> ATCC® 13315™	Urease positive, pink slope
<i>Klebsiella pneumoniae</i> ATCC® 13883™	Weak urease positive, pale pink slope

Limitations

The alkaline reaction produced in this medium after prolonged incubation may not be caused by urease activity. Check using medium without urea.

Do not heat or reheat the medium because urea decomposes very easily.

For the detection of urease-positive *Proteae* the reaction must be read within the first 2-6 hours of incubation.

Performance Characteristics

Accuracy has been demonstrated through review of the QC data. Correct detection of Enterobacteriaceae is confirmed by the inclusion of a well-characterised isolate in the QC processes performed as part of the manufacture of each batch of the devices. The precision of Urea Agar Slope (BO0337B) was demonstrated by an overall pass rate of 80% obtained for the product over ten months of testing (18.SEP.2021–25.JUL.2022; 10 batches). The precision of Urea Agar Slope (EB0337B) was demonstrated by an overall pass rate of 90% obtained for the product over nine months of testing (13.SEP.2021–13.JUN.2022; 10 batches). An overall pass-rate of 80% (based on data from the most recent 10 batches) provides sufficient data on accuracy considering the low risk of the device and it is not reflective of the product that is released for clinical use. With the normal variation in the manufacturing process, a small number of failures may occur, however, there is no risk to the end user as batches will be captured in QC if out of specification. This shows that the performance is reproducible, thus meeting the specifications set out in the
















Performance Evaluation Plan.


Urea Agar Slope (BO0337B & EB0337B) devices are tested in-house as part of the QC process since 1997 (BO0337B) and 1998 (EB0337B). For target organisms, when using 10^4 - 10^5 cfu inoculum of *Escherichia coli*, *Proteus mirabilis* or *Klebsiella pneumoniae* and incubating the device at $37^\circ \pm 2^\circ\text{C}$ for 18-24 hours, the user can expect urease reaction as listed in this document.

Bibliography

- Centers for Disease Control and Prevention. 2021. 'Urinary Tract Infection'. <https://www.cdc.gov/antibiotic-use/uti.html>.

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

	Importer - To indicate the entity importing the medical device into the locale. Applicable to the European Union
Made in the United Kingdom	Made in the United Kingdom



©2022 Thermo Fisher Scientific Inc. All rights reserved. ATCC and ATCC catalogue marks are a trademark of American Type Culture Collection.

NCTC and NCTC catalogue marks are a trademark of National Collection of Type Cultures.

All other trademarks are the property of Thermo Fisher Scientific Inc. and its subsidiaries.



Oxoid Limited, Wade Road, Basingstoke, Hampshire, RG24 8PW, United Kingdom



For technical assistance please contact your local distributor.

Revision Information

Version	Date of modifications introduced
2.0	2023-11-21



www.thermofisher.com

Thermo
SCIENTIFIC

Urea Agar Slope

HR

REF **BO0337B & EB0337B**

Namjena

Proizvodi urea agar (kosina) (BO0337B i EB0337B) neselektivni su mediji namijenjeni za uporabu u uzgojnim medijima. Proizvodi urea agar (kosina) (BO0337B i EB0337B) namijenjeni su za otkrivanje brze aktivnosti ureaze *Proteae* i spore aktivnosti ureaze nekih bakterija roda *Enterobacteriaceae* u uzorcima stolice i urina.

Proizvodi su namijenjeni za uporabu u dijagnostičkom tijeku rada kao pomoć liječnicima u određivanju potencijalnih mogućnosti liječenja bolesnika kod kojih postoji sumnja na infekcije bakterijama roda *Enterobacteriaceae*.

Proizvodi su namijenjeni samo za profesionalnu uporabu, nisu automatizirani niti su nadopuna dijagnostičkim postupcima.

Sažetak i objašnjenje

Infekcije mokraćnog sustava uobičajene su infekcije koje se javljaju kada bakterije napadnu mokraćni sustav ulaskom kroz uretru. Te infekcije mogu zahvatiti različite dijelove mokraćnog sustava, ali najčešća je cistitis (infekcija mokraćnog mjehura). Infekcije mokraćnog sustava češće se javljaju kod žena jer im je uretra kraća i stoga bliže rektumu. Time se povećava vjerojatnost infekcije jer je bakterijama lakše ući u mokraćni sustav. Infekcije bubrega još su jedan oblik infekcije mokraćnog sustava. One su ozbiljnije od cistitisa, ali daleko rjeđe. Simptomi cistitisa mogu uključivati učestalo mokrenje ili potrebu za mokrenjem dok imate prazan mjehur, osjećaj pečenja ili bol tijekom mokrenja, krv u mokraći i pritisak ili grčeve u donjem dijelu trbuha. Simptomi infekcija bubrega vrlo su različiti. Oni mogu uključivati zimicu, groznicu, mučninu i povraćanje te bol u donjem dijelu leđa¹. Postoje mnoge vrste bakterija koje uzrokuju infekcije mokraćnog sustava, ali tri vrste koje su obuhvaćene ovom datotekom jesu *Escherichia coli*, *Staphylococcus aureus* i *Proteus mirabilis*.

Zbog težine infekcija mokraćnog sustava kod velikog broja bolesnika, vrlo je važno moći izolirati i identificirati uobičajene mikroorganizme koji ih uzrokuju iz uzoraka urina. Rana dijagnoza ključna je u prevenciji i smanjenju infekcija mokraćnog sustava.

Načelo metode

Urea agar je blago puferirani medij koji sadrži ureu i fenol crveno, pH indikator. Kada organizmi iskoriste ureu, nastaje amonijak koji medij pretvara u alkalni. Indikator, fenol crveno, mijenja boju medija od blijedožute do ružičasto-crvene u alkalnoj sredini. Želatina pepton potiče brzi rast crijevnih gram-negativnih bacila, što omogućuje smanjenje vremena inkubacije. Dekstroza stimulira aktivnost ureaze u organizmima koji sporo hidroliziraju ureu.

Uobičajena formula

	grama po litri
Pepton	1,0
Glukoza	1,0
Natrijev klorid	5,0
Dinatrijev fosfat	1,2
Kalijev dihidrogenfosfat	0,8
Fenol crveno	0,012
Agar	15,0

Fizički izgled

Boja	Narančasta
Bistina	Bistro
Težina punjenja	3,0 – 3,3 g
pH	6,8 ± 0,2

Priloženi materijali

BO0337B: 24 x bočice s urea agarom (kosina)
EB0337B: 200 x bočica s urea agarom (kosina)

Svaka se bočica smije upotrijebiti samo jednom.

Potrebni materijali koji nisu isporučeni

- Inokulacijske petlje
- Brisovi
- Spremnici za prikupljanje
- Inkubatori
- Organizmi za kontrolu kvalitete

Skладиštenje

- Čuvajte proizvod u originalnom pakiranju na 2 – 10 °C do uporabe.
- Proizvod se može koristiti do isteka roka valjanosti navedenog na naljepnici.
- Čuvati podalje od svjetlosti.
- Prije uporabe pustite da proizvod postigne sobnu temperaturu.
- Nemojte inkubirati prije uporabe.

Upozorenja i mjere opreza

- Samo za in vitro dijagnostičku uporabu.
- Samo za profesionalnu uporabu.
- Pregledajte pakiranje proizvoda prije prve uporabe.
- Nemojte upotrebljavati proizvod ako ima vidljivih oštećenja na pakiranju ili pločicama.
- Nemojte upotrebljavati proizvod nakon isteka navedenog roka valjanosti.
- Nemojte upotrebljavati proizvod ako su prisutni znakovi kontaminacije.
- Nemojte upotrebljavati proizvod ako je došlo do promjene boje ili su prisutni drugi znakovi narušenja kvalitete.
- Svaki je laboratorij odgovaran za upravljanje proizvedenim otpadom u skladu s prirodom i stupnjem opasnosti otpada te za njegovu obradu ili zbrinjavanje u skladu s primjenjivim saveznom, državnim i lokalnim propisima. Potrebno je pročitati upute i pažljivo ih se pridržavati. To uključuje odlaganje iskorištenih ili neiskorištenih reagensa kao i bilo kojeg drugog kontaminiranog jednokratnog materijala pridržavajući se postupaka za zarazne ili potencijalno zarazne proizvode.

Proučite Sigurnosno-tehnički list za sigurno rukovanje i odlaganje proizvoda (www.thermofisher.com).

Ozbiljni štetni događaji

Svaki ozbiljan štetni događaj do kojeg je došlo vezano uz proizvod treba prijaviti proizvođaču i relevantnom regulatornom tijelu države u kojoj se korisnik i/ili bolesnik nalazi.

Prikupljanje uzoraka, rukovanje i skladištenje

Uzorke treba prikupiti i s njim postupiti u skladu s lokalnim i preporučenim smjernicama, kao što su Standardi za mikrobiološka istraživanja u Ujedinjenom Kraljevstvu (UK SMI) ID 16, ID 20, ID 21, ID 24.

Postupak

- Pustite da proizvod postigne sobnu temperaturu.
- Inokulirajte i razmažite uzorak na medij.
- Alternativno, aseptičkom tehnikom odaberite jednu dobro izoliranu koloniju sterilnom iglom za inokulaciju i nekoliko puta ubodite iglu kroz agar do dna bočice ili epruvete.
- Vratite poklopac i labavo ga zategnite tijekom inkubacije kako biste omogućili izmjenu plinova.
- Inkubirajte ubodeni agar / kosinu 18 do 24 sata na 37 ± 2 °C.
- Vizualno pregledajte bočice kako biste procijenili rast i boju kolonije pod dobrim osvjetljenjem.

Tumačenje

Negativno na ureazu, nema promjene boje / žuti medij ukazuje na *Escherichia coli*.

Pozitivno na ureazu s ružičastim kosinama ukazuje na *Proteus mirabilis* ili *Proteus hauseri*.

Slabo pozitivno na ureazu s blijedoružičastim kosinama ukazuje na *Klebsiella pneumoniae*.

Kontrola kvalitete

Korisnik je odgovoran za provedbu testiranja kontrole kvalitete uzimajući u obzir namjenu medija te u skladu s primjenjivim lokalnim propisima (učestalost, broj sojeva, temperatura inkubacije itd.).

Učinkovitost ovog medija može se provjeriti testiranjem sljedećih referentnih sojeva.

Uvjeti inkubacije: 18 – 24 sata na $37^\circ \pm 2$ °C

Pozitivne kontrole	
Razina inokuluma: 10.000 – 100.000 cfu	
<i>Escherichia coli</i> ATCC® 25922™	Negativno na ureazu, nema promjene boje / žuti medij
<i>Proteus mirabilis</i> NCTC 10975	Pozitivno na ureazu, ružičasta kosina
<i>Proteus hauseri</i> ATCC® 13315™	Pozitivno na ureazu, ružičasta kosina
<i>Klebsiella pneumoniae</i> ATCC® 13883™	Slabo pozitivno na ureazu, blijedoružičasta kosina

Ograničenja

Alkalna reakcija nastala u ovom mediju nakon produljene inkubacije ne mora biti uzrokovana aktivnošću ureaze. Provjerite korištenjem medija bez uree.

Nemojte zagrijavati niti ponovno zagrijavati medij jer se urea vrlo lako razgrađuje.

Za otkrivanje *Proteae* pozitivne na ureazu, reakcija se mora očitati unutar prvih 2 – 6 sati inkubacije.

Karakteristike učinkovitosti

Točnost je dokazana pregledom podataka kontrole kvalitete. Ispravno otkrivanje bakterija Enterobacteriaceae potvrđeno je uključivanjem dobro karakteriziranog izolata u postupke kontrole kvalitete koji se izvode kao dio proizvodnje svake serije proizvoda. Preciznost proizvoda urea agar (kosina) (BO0337B) dokazana je ukupnom stopom prolaznosti od 80 % dobivenom za proizvod tijekom deset mjeseci testiranja (18. rujna 2021. – 25. srpnja 2022.; 10 serija). Preciznost proizvoda urea agar (kosina) (EB0337B) dokazana je ukupnom stopom prolaznosti od 90 % dobivenom za proizvod tijekom devet mjeseci testiranja (13. rujna 2021. – 13. lipnja 2022.; 10 serija). Ukupna stopa prolaznosti od 80 % (na temelju podataka iz posljednjih 10 serija) daje dovoljno podataka o točnosti s obzirom na nizak rizik proizvoda i ne odražava proizvod koji je stavljen na tržište za kliničku uporabu. Uz uobičajene varijacije u

procesu proizvodnje, može se dogoditi mali broj neispravnosti, međutim, nema rizika za krajnjeg korisnika jer će serije biti pronađene tijekom kontrole kvalitete ako su izvan specifikacije. To pokazuje da je učinkovitost obnovljiva, čime se ispunjavaju specifikacije navedene u Planu ocjenjivanja učinkovitosti.

Proizvodi urea agar (kosina) (BO0337B i EB0337B) testirani su unutar tvrtke kao dio procesa kontrole kvalitete od 1997. (BO0337B) odnosno 1998. (EB0337B). Za ciljne organizme, kada se rabi 10^4 – 10^5 cfu inokuluma bakterija *Escherichia coli*, *Proteus mirabilis* ili *Klebsiella pneumoniae* i proizvod se inkubira na $37^\circ \pm 2$ °C tijekom 18 – 24 sata, korisnik može očekivati reakciju ureaze kako je navedeno u ovom dokumentu.

Bibliografija

1. Centers for Disease Control and Prevention. 2021. 'Urinary Tract Infection'. <https://www.cdc.gov/antibiotic-use/uti.html>.

Kazalo simbola

Simbol	Definicija
	Kataloški broj
	In vitro dijagnostički medicinski proizvod
	Broj serije
	Granica temperature
	Rok valjanosti
	Čuvati podalje od sunčeve svjetlosti
	Ne upotrebljavati višekratno
	Proučite upute za upotrebu ili elektroničke upute za upotrebu
	Sadrži dovoljno za <n> testova
	Ne upotrebljavati ako je pakiranje oštećeno; proučite upute za uporabu
	Proizvođač
	Ovlašteni zastupnik u Europskoj zajednici/ Europskoj uniji

	Europska ocjena sukladnosti
	Ocjena sukladnosti u Ujedinjenoj Kraljevini
	Jedinstvena identifikacija proizvoda
	Uvoznik – za označavanje subjekta koji uvozi medicinski proizvod u pojedinu zemlju. Primjenjivo u Europskoj uniji
	Proizvedeno u Ujedinjenoj Kraljevini



© 2022. Thermo Fisher Scientific Inc. Sva prava pridržana. ATCC i ATCC kataloške oznake zaštitni su znak Američke zbirke tipskih kultura. NCTC i NCTC kataloške oznake zaštitni su znak Nacionalne zbirke tipskih kultura. Svi ostali zaštitni znakovi vlasništvo su društva Thermo Fisher Scientific Inc. i njegovih podružnica.



Oxoid Limited, Wade Road, Basingstoke, RG24 8PW, Engleska



Za tehničku pomoć obratite se svom lokalnom distributeru.

Informacije o reviziji

Verzija	Datum uvedenih izmjena
2.0	2023-11-21



www.thermofisher.com

Thermo
SCIENTIFIC

Kaliumdihydrogenphosphat 0,8
Fenolrødt 0,012
Agar 15,0

Urea Agar Slope

DA

REF BO0337B & EB0337B

Tilsigtet anvendelse

Urea-agar-hældningsanordninger (BO0337B og EB0037B) er ikke-selektive medier beregnet til brug i dyrkningsmedier. Urea-agar-hældningsanordninger (BO0337B og EB0037B) er beregnet til at blive brugt til at detektere hurtig ureaseaktivitet af *Proteae* og ikke-hurtig ureaseaktivitet hos nogle *Enterobacteriaceae* i fæces- og urinprøver.

Enhederne er beregnet til at blive brugt i en diagnostisk arbejdsdag for at hjælpe klinikere med at bestemme potentielle behandlingsmuligheder for patienter, der mistænkes for at have *Enterobacteriaceae* infektioner.

"Enhederne er kun til professionel brug, er ikke automatiserede og er heller ikke ledsagende diagnostik."

Oversigt og forklaring

Urinvejsinfektioner, ofte omtalt som UTI'er, er almindelige infektioner, der opstår, når bakterier invaderer urinvejene ved at trænge ind via urinrøret. Disse infektioner kan påvirke forskellige dele af urinvejene, men den mest udbredte er blærebetændelse (en blæreinfektion). UTI'er forekommer hyppigere hos kvinder, da urinrøret er kortere og derfor tættere på endetarmen. Dette øger sandsynligheden for infektion, da det er lettere for bakterier at trænge ind i urinvejene. Nyreinfektioner er en anden form for UTI. Disse er mere alvorlige end blærebetændelse, men langt mindre almindelige. Symptomerne på blærebetændelse kan omfatte hyppig vandladning eller behov for at tisse, mens du har en tom blære, en brændende fornemmelse eller smerte under vandladning, blod i urinen og tryk eller kramper i den nedre del af maven. Symptomerne på nyreinfektioner er meget forskellige. Disse kan omfatte kulderystelser, feber, kvalme og opkastning og smerter i lænden¹. Der er mange arter af bakterier, der forårsager UTI'er, men de tre, der er omfattet af denne fil, er *Escherichia coli*, *Staphylococcus aureus* og *Proteus mirabilis*.

På grund af sværhedsgraden af UTI'er hos et stort antal patienter, er det derfor meget vigtigt at være i stand til at isolere og identificere almindelige UTI-fremkaldende mikroorganismer fra urinprøver. Tidlig diagnose er afgørende for forebyggelse og reduktion af urinvejsinfektioner.

Metodens principper

Urea-agar er et let bufret medium, der indeholder urinstof og phenolrødt, en pH-indikator. Når organismer udnytter urinstof, dannes der ammoniak, som gør mediet alkalisk. Indikatoren, phenolrød, ændrer mellemfarven fra bleggul til pinkrød i et alkalisk miljø. Gelatinpepton fremmer hurtig vækst af enteriske gramnegative baciller, hvilket tillader et fald i inkubationstid. Dextrose stimulerer ureaseaktivitet i de organismer, som langsomt hydrolyserer urea.

Typisk formel

	gram per liter
Pepton	1,0
Glukose	1,0
Natriumklorid	5,0
Dinatriumfosfat	1,2

Fysisk fremtoning

Farve	Orange
Klarhed	Ryd
Fill vægt	3,0 -3,3 g
pH	6,8 ± 0,2

Leverede materialer

BO0337B: 24 x urea agar hældningsflasker
EB0337B: 200 x urea agar hældningsflasker

Hver flaske bør kun bruges én gang.

Nødvendige materialer, som ikke medfølger

- Inokuleringsløkker
- Pødepindsprøve
- Indsamlingsbeholdere
- Inkubatorer
- Kvalitetskontrolorganismer

Opbevaring

- Opbevar produktet i dets originale emballage ved 2-10 °C, indtil det bruges.
- Produktet kan bruges indtil den udløbsdato, der står på etiketten.
- Opbevares væk fra lys.
- Lad produktet opnå stuetemperatur før brug.
- Må ikke inkuberes før brug.

Advarsler og forholdsregler

- Kun til in vitro-diagnostisk brug.
- Kun til professionel brug.
- Efterse produktets emballage, før det bruges første gang.
- Brug ikke produktet, hvis der er synlige skader på emballagen eller pladerne.
- Brug ikke produktet efter den anførte udløbsdato.
- Brug ikke enheden, hvis der er tegn på kontaminering.
- Brug ikke enheden, hvis farven er ændret, eller der er andre tegn på nedbrydning.
- Det er hvert laboratoriums ansvar at håndtere produceret affald i overensstemmelse med dets art og grad af fare og at få det behandlet eller bortskaffet i overensstemmelse med alle gældende føderale, statslige og lokale regler. Vejledninger skal læses og følges omhyggeligt. Dette omfatter bortskaffelse af brugte eller ubrugte reagenser samt ethvert andet kontamineret engangsmateriale i henhold til procedurer for infektiøse eller potentielt infektiøse produkter.

Se sikkerhedsdatabladet (SDS) for sikker håndtering og bortskaffelse af produktet (www.thermofisher.com).

Alvorlige hændelser

Enhver alvorlig hændelse, der er opstået i forbindelse med udstyret, skal rapporteres til producenten og den relevante tilsynsmyndighed, hvor brugeren og/eller patienten er etableret.

Prøveindsamling, håndtering og opbevaring

Prøver skal indsamles og håndteres efter lokale anbefalede retningslinjer, såsom UK Standards for Microbiology Investigations (UK SMI) ID 16, ID 20, ID 21, ID 24,

Procedure

- Lad produktet ækvilibrere til stuetemperatur.
- Inokulér og strøg prøven på mediet.
- eller alternativt ved brug af aseptisk teknik, vælg en enkelt velisoleret koloni med en steril podestiknål og stik nålen flere gange gennem agaren til bunden af hætteglasset eller røret
- Sæt hættten på igen og spænd løst til under inkubering for at tillade gasudveksling
- Inkuber den stukkede agar/skråning i 18-24 timer ved 37 ± 2 °C.
- Inspicer flaskerne visuelt for at vurdere kolonivækst og farve under god belysning.

Fortolkning

Urease negativ, ingen farveændring/gult medium indikerer *Escherichia coli*.

Urease positiv med lyserøde skråninger indikerer *Proteus mirabilis* eller *Proteus hauseri*.

Svag urease positiv med lyserøde skråninger indikerer *Klebsiella pneumoniae*.

Kvalitetskontrol

Det er brugerens ansvar at udføre kvalitetskontroltest under hensyntagen til den tilsigtede brug af mediet og i overensstemmelse med lokale gældende regler (hyppighed, antal stammer, inkubationstemperatur osv.).

Ydeevnen af dette medie kan verificeres ved at teste følgende referencestammer.

Inkubationsbetingelser: 18-24 timer @ 37 ± 2 °C

Positiv kontroller	
Inokulum niveau: 10.000-100.000 cfu	
<i>Escherichia coli</i> ATCC® 25922™	Urease negativ, ingen farveændring/ gul medium
<i>Proteus mirabilis</i> NCTC 10975	Urease positiv, lyserød hældning
<i>Proteus hauseri</i> ATCC® 13315™	Urease positiv, lyserød hældning
<i>Klebsiella pneumoniae</i> ATCC® 13883™	Svag ureasepositiv, lyserød hældning

Begrænsninger

Den alkaliske reaktion produceret i dette medium efter langvarig inkubation er muligvis ikke forårsaget af ureaseaktivitet. Tjek med medium uden urinstof.

Opvarm eller genopvarm ikke mediet, da urinstof nedbrydes meget let.

For påvisning af urease-positive *Proteae* skal reaktionen aflæses inden for de første 2-6 timer efter inkubation.

Ydelseskarakteristik

Nøjagtighed er blevet demonstreret gennem gennemgang af kvalitetskontrol-dataene. Korrekt påvisning af Enterobacteriaceae bekræftes ved at inkludere et velkarakteriseret isolat i kvalitetskontrol-processerne, der udføres som en del af fremstillingen af hver batch af enhederne. Præcisionen af Urea Agar Slope (BO0337B) blev demonstreret ved en samlet beståelsesrate på 80 % opnået for produktet over ti måneders test (18.SEP.2021-25.JUL.2022; 10 batches). Præcisionen af Urea Agar Slope (EB0337B) blev demonstreret ved en samlet beståelsesrate på 90 % opnået for produktet over ni måneders test (13.SEP.2021-13.JUN.2022; 10 batches). En samlet beståelsesrate på 80 % (baseret på data fra de seneste 10 batches) giver tilstrækkelige data om nøjagtighed i betragtning af enhedens lave risiko, og denafspejler ikke

det produkt, der er frigivet til klinisk brug. Med den normale variation i fremstillingsprocessen kan der forekomme et lille antal fejl, dog er der ingen risiko for slutbrugeren, da batches vil blive fanget i kvalitetskontrol, hvis de ikke er i specifikation. Dette viser, at ydeevnen er reproducerbar og dermed opfylder specifikationerne i Performance Evaluation Plan.





Urea-agar-hældningsenheder (BO0337B & EB0337B) er testet internt som en del af kvalitetskontrol-processen siden 1997 (BO0337B) og 1998 (EB0337B). Til målorganismer, ved brug af 10^4 - 10^5 cfu inokulum af *Escherichia coli*, *Proteus mirabilis* eller *Klebsiella pneumoniae* og inkubering af enheden ved 37 ± 2 °C i 18-24 timer kan brugeren forvente en ureasereaktion som angivet i dette dokument.

Bibliografi

1. Centers for Disease Control and Prevention. 2021. 'Urinary Tract Infection'. <https://www.cdc.gov/antibiotic-use/uti.html>.

Symbolforklaring

Symbol	Definition
	Katalognummer
	In vitro-diagnostisk medicinsk udstyr
	Batchkode
	Temperaturgrænse
	Sidste anvendelsesdato
	Holdes væk fra sollys
	Må ikke genbruges
	Se brugsanvisningen eller den elektroniske brugsanvisning
	Tilstrækkeligt indhold til <n> tests
	Må ikke bruges, hvis emballagen er beskadiget, og se brugsanvisningen
	Fabrikant
	Autoriseret repræsentant i Det Europæiske Fællesskab/Den Europæiske Union

	Europæisk overensstemmelsesvurdering
	Britisk konformitetsvurdering
	Unik udstyrsidentifikation
	Importør – Angiver den enhed, der importerer det medicinske udstyr til regionen/området. Gælder for EU
Made in the United Kingdom	Lavet i Storbritannien



© 2022 Thermo Fisher Scientific Inc. Alle rettigheder forbeholdes.

ATCC og ATCC-katalogmærker er varemærker tilhørende American Type Culture Collection.

NCTC- og NCTC-katalogmærker er varemærker tilhørende National Collection of Type Cultures.

Alle andre varemærker tilhører Thermo Fisher Scientific Inc. og dets datterselskaber.



Oxoid Limited, Wade Road, Basingstoke, RG24 8PW, England



Kontakt din lokale distributør i forbindelse med hjælp til tekniske spørgsmål.

Revisionsoplysninger

Version	Dato for indførte ændringer
2.0	2023-11-21



www.thermofisher.com

Thermo
SCIENTIFIC

Glukose	1,0
Natriumchlorid	5,0
Dinatriumphosphat	1,2
Kaliumdihydrogenphosphat	0,8
Phenolrot	0,012
Agar	15,0

Urea Agar Slope

DE

REF **BO0337B & EB0337B**

Verwendungszweck

Harnstoff-Agar-Neigung-Produkte (BO0337B und EB0337B) sind nicht-selektive Medien, die für die Verwendung in Kulturmedien bestimmt sind. Harnstoff-Agar-Neigung-Produkte (BO0337B und EB0337B) sind zur Bestimmung der schnellen Urease-Aktivität von *Proteae* und der nicht-schnellen Urease-Aktivität von einigen *Enterobacteriaceae* in Fäkal- und Urinproben vorgesehen.

Diese Produkte sind für die Verwendung in einem diagnostischen Arbeitsablauf vorgesehen, um Kliniker bei der Bestimmung möglicher Behandlungsoptionen für Patienten mit Verdacht auf *Enterobacteriaceae*-Infektionen zu unterstützen.

Die Produkte sind nur für den professionellen Gebrauch bestimmt, sie sind nicht automatisiert und sie sind auch keine Begleitdiagnostik.

Zusammenfassung und Erläuterung

Harnwegsinfektionen, oft auch als UTI bezeichnet, sind häufige Infektionen, die auftreten, wenn Bakterien durch die Harnröhre in die Harnwege eindringen. Diese Infektionen können verschiedene Teile der Harnwege betreffen, am häufigsten ist allerdings die Zystitis (eine Blasenentzündung). UTI treten häufiger bei Frauen auf, da die Harnröhre kürzer ist und somit näher am Enddarm liegt. Dies erhöht die Wahrscheinlichkeit einer Infektion, da Bakterien leichter in die Harnwege gelangen können. Niereninfektionen sind eine weitere Form von UTI. Diese sind schwerwiegender als eine Zystitis, aber weitaus seltener. Zu den Symptomen einer Zystitis gehören häufiges Wasserlassen oder Harndrang bei leerer Blase, Brennen oder Schmerzen beim Wasserlassen, Blut im Urin und Druck oder Krämpfe im Unterleibsbereich. Die Symptome von Niereninfektionen sind sehr unterschiedlich. Diese können Schüttelfrost, Fieber, Übelkeit und Erbrechen sowie Schmerzen im unteren Rückenbereich umfassen¹. Es gibt viele Spezies von Bakterien, die UTI verursachen. Die drei Spezies, die in diesem Dokument behandelt werden, sind *Escherichia coli*, *Staphylococcus aureus* und *Proteus mirabilis*.

Aufgrund des Schweregrads von UTI bei einer großen Anzahl von Patienten ist es daher äußerst wichtig, häufige UTI-Verursachende Mikroorganismen aus Urinproben isolieren und identifizieren zu können. Eine frühzeitige Diagnose ist entscheidend für die Vorbeugung und Reduzierung von Harnwegsinfektionen.

Das Prinzip der Methode

Harnstoff-Agar ist ein leicht gepuffertes Medium, das Harnstoff und den pH-Indikator Phenolrot enthält. Wenn Organismen Harnstoff verwerten, wird Ammoniak gebildet, welches das Medium alkalisch macht. Der Indikator Phenolrot verändert die Farbe des Mediums in alkalischem Milieu von blassgelb zu rosarot. Gelatine-Pepton fördert das schnelle Wachstum enterischer gram-negativer Bazillen und kann die Inkubationszeit verkürzen. Dextrose stimuliert die Urease-Aktivität in denjenigen Organismen, die Harnstoff langsam hydrolysieren.

Typische Formel

	<u>Gramm pro Liter</u>
Pepton	1,0

Physische Erscheinung

Farbe	Orange
Klarheit	Klar
Gewicht	der 3,0–3,3 g
Füllung	
pH	6,8 ± 0,2

Mitgeliefertes Material

BO0337B: 24 x Harnstoff-Agar-Neigung-Flaschen
EB0337B: 200 x Harnstoff-Agar-Neigung-Flaschen

Jede Flasche sollte nur einmal verwendet werden.

Erforderliche, aber nicht mitgelieferte Materialien

- Beimpfen von Schleifen
- Tupfer
- Entnahmebehälter
- Inkubatoren
- Organismen für die Qualitätskontrolle

Lagerung

- Lagern Sie das Produkt bis zur Verwendung in der Originalverpackung bei 2–10 °C.
- Das Produkt kann bis zu dem auf dem Etikett angegebenen Verfallsdatum verwendet werden.
- Vor Licht geschützt aufbewahren.
- Lassen Sie das Produkt vor der Verwendung auf Raumtemperatur kommen.
- Vor der Verwendung nicht inkubieren.

Warnungen und Vorsichtsmaßnahmen

- 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 Platten 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, zu melden.

Reaktion innerhalb der ersten 2–6 Stunden nach der Inkubation abgelesen werden.

Entnahme, Handhabung und Lagerung von Proben

Die Probenentnahme und -behandlung sollte gemäß den empfohlenen lokalen Richtlinien erfolgen, wie z. B. den UK Standards for Microbiology Investigations (UK SMI) ID 16, ID 20, ID 21 und ID 24.

Verfahren

- Lassen Sie das Produkt auf Raumtemperatur kommen.
- Inokulieren Sie die Probe und streichen Sie sie auf das Medium.
- Alternativ können Sie unter Anwendung einer aseptischen Technik eine einzelne, gut isolierte Kolonie mit einer sterilen Inokulationsnadel entnehmen und die Nadel mehrmals durch den Agar bis zum Boden des Fläschchens oder Röhrchens stechen.
- Setzen Sie die Kappe wieder auf und ziehen Sie sie zum Inkubieren locker fest, um einen Gasaustausch zu ermöglichen.
- Inkubieren Sie den Stichagar/Schrägagar 18–24 Stunden lang bei $37 \pm 2 \text{ }^\circ\text{C}$.
- Untersuchen Sie die Flaschen visuell, um das Wachstum und die Farbe der Kolonien bei guter Beleuchtung zu beurteilen.

Interpretation

Eine negative Urease ohne Farbveränderung/gelbes Medium zeigt *Escherichia coli an*.

Eine positive Urease mit rosa Neigungen zeigt *Proteus mirabilis* oder *Proteus hauseri an*.

Eine schwach positive Urease mit blassrosa Neigungen zeigt *Klebsiella pneumoniae an*.

Qualitätskontrolle

Es liegt in der Verantwortung des Anwenders, Qualitätskontrolltests unter Berücksichtigung der beabsichtigten Verwendung des Mediums und in Übereinstimmung mit allen vor Ort geltenden Vorschriften (Häufigkeit, Anzahl der Stämme, Inkubationstemperatur usw.) durchzuführen.

Die Leistungsfähigkeit dieses Mediums kann durch Testen der folgenden Referenzstämme überprüft werden.

Inkubationsbedingungen: 18–24 Stunden bei $37 \pm 2 \text{ }^\circ\text{C}$

Positiv-Kontrollen	
Inokulumkonzentration: 10.000–100.000 KBE	
<i>Escherichia coli</i> ATCC® 25922™	Urease negativ, keine Farbveränderung/gelbes Medium
<i>Proteus mirabilis</i> NCTC 10975	Urease positiv, rosa Neigung
<i>Proteus hauseri</i> ATCC® 13315™	Urease positiv, rosa Neigung
<i>Klebsiella pneumoniae</i> ATCC® 13883™	Schwach positive Urease, blassrosa Neigung

Beschränkungen

Die in diesem Medium nach längerer Inkubation hervorgerufene alkalische Reaktion darf nicht durch Urease-Aktivität verursacht werden. Mit einem Medium ohne Harnstoff überprüfen.

Das Medium nicht (erneut) erhitzen, da sich Harnstoff sehr leicht zersetzt.

Für den Nachweis von urease-positiven *Proteae* muss die

Leistungsmerkmale







Die Genauigkeit wurde durch die Überprüfung der QC-Daten nachgewiesen. Der korrekte Nachweis von Enterobacteriaceae wird durch die Aufnahme eines gut charakterisierten Isolats in die QK-Prozesse bestätigt, die im Rahmen der Herstellung jeder Charge der Produkte durchgeführt werden. Die Präzision von Harnstoff-Agar-Neigung (BO0337B) wurde durch eine Gesamterfolgsquote von 80 % nachgewiesen, die über einen Testzeitraum von zehn Monaten (18. SEP. 2021 bis 25. JUL. 2022; 10 Chargen) für das Produkt erzielt wurde. Die Präzision von Harnstoff-Agar-Neigung (EB0337B) wurde durch eine Gesamterfolgsquote von 90 % nachgewiesen, die über einen Testzeitraum von neun Monaten (13. SEP. 2021 bis 13. JUN. 2022; 10 Chargen) für das Produkt erzielt wurde. Eine Gesamterfolgsquote von 80 % (basierend auf Daten der letzten 10 Chargen) liefert ausreichende Daten zur Genauigkeit unter Berücksichtigung des geringen Risikos des Produkts und spiegelt nicht das Produkt wider, das für die klinische Verwendung freigegeben ist. Bei den normalen Schwankungen im Herstellungsprozess kann es zu einer kleinen Anzahl von Fehlern kommen; allerdings besteht kein Risiko für den Endverbraucher, da Chargen im Rahmen der QK erfasst werden, wenn sie außerhalb der Spezifikationen liegen. Dies zeigt, dass die Leistung reproduzierbar ist und somit die im Leistungsbewertungsplan festgelegten Spezifikationen erfüllt.

Die Harnstoff-Agar-Neigung-Produkte (BO0337B und EB0337B) werden seit 1997 (BO0337B) und 1998 (EB0337B) im Rahmen der QK-Prozesse hausintern getestet. Für Zielorganismen bei Verwendung von 10^4 – 10^5 KBE Inokulum von *Escherichia coli*, *Proteus mirabilis* oder *Klebsiella pneumoniae* und einer Inkubation des Produkts bei $37 \pm 2 \text{ }^\circ\text{C}$ für 18–24 Stunden kann der Anwender eine Urease-Reaktion wie in diesem Dokument beschrieben erwarten.

Bibliographie

1. Centers for Disease Control and Prevention. 2021. 'Urinary Tract Infection'. <https://www.cdc.gov/antibiotic-use/uti.html>.

Symbollegende

Symbol	Definition
	Katalognummer
	Medizinprodukt zum In-vitro-Diagnostikum
	Chargencode
	Temperaturgrenze
	Haltbarkeitsdatum
	Vom Sonnenlicht fernhalten

	Nicht wiederverwenden
	Konsultieren Sie die Gebrauchsanweisung oder konsultieren Sie die elektronische Gebrauchsanweisung
	Enthält ausreichend für <n> Tests
	Nicht verwenden, wenn die Verpackung beschädigt ist und die Gebrauchsanweisung beachten
	Hersteller
	Bevollmächtigter Vertreter in der Europäischen Gemeinschaft/Europäischen Union
	Europäische Konformitätsbewertung
	Konformitätsbewertung des Vereinigten Königreichs
	Eindeutige Kennung des Produkts
	Importeur – Angabe der juristischen Person, die das Medizinprodukt in die Region importiert. Gilt für die Europäische Union.
	Hergestellt im Vereinigten Königreich

ATCC Licensed
Derivative[®]

© 2022 Thermo Fisher Scientific Inc. Alle Rechte vorbehalten.
ATCC und ATCC-Katalogmarken sind eine Marke der American Type Culture Collection.
NCTC und NCTC-Katalogmarken sind eine Marke der National Collection of Type Cultures.
Alle anderen Marken sind Eigentum der Thermo Fisher Scientific Inc. und ihrer Tochtergesellschaften.



Oxoid Limited, Wade Road, Basingstoke, RG24 8PW, England



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

Informationen zur Revision

Version	Datum der eingeführten Änderungen
2.0	2023-11-21



Urea Agar Slope



REF BO0337B & EB0337B

Προβλεπόμενη χρήση

Τα ιατροτεχνολογικά προϊόντα κεκλιμένης επιφάνειας με άγαρ ουρίας Urea Agar Slope (BO0337B και EB0037B) είναι μη εκλεκτικά μέσα που προορίζονται για χρήση σε μέσα καλλιέργειας. Τα ιατροτεχνολογικά προϊόντα Urea Agar Slope (BO0337B και EB0037B) προορίζονται για χρήση για την ανίχνευση της ταχείας δράσης ουρεάσης του *Proteae* και την μη ταχεία δράση ουρεάσης ορισμένων *Enterobacteriaceae* σε δείγματα κοπράνων και ούρων.

Αυτά τα ιατροτεχνολογικά προϊόντα προορίζονται για χρήση σε μια διαγνωστική ροή εργασιών προκειμένου να βοηθηθούν οι κλινικοί ιατροί στον καθορισμό πιθανών θεραπευτικών επιλογών για ασθενείς όπου υπάρχει υποψία ότι πάσχουν από βακτηριακή λοίμωξη που οφείλεται σε *Enterobacteriaceae*.

Τα ιατροτεχνολογικά προϊόντα προορίζονται αποκλειστικά για επαγγελματική χρήση, δεν είναι αυτοματοποιημένα και δεν αποτελούν συνοδευτικά διαγνωστικά μέσα.

Περίληψη και Επεξήγηση

Οι λοιμώξεις του ουροποιητικού συστήματος, που συχνά αναφέρονται ως ουρολοιμώξεις, είναι συχνές λοιμώξεις που προκύπτουν όταν βακτήρια εισβάλλουν στο ουροποιητικό σύστημα εισερχόμενοι μέσω της ουρήθρας. Αυτές οι λοιμώξεις μπορεί να επηρεάσουν διαφορετικά τμήματα του ουροποιητικού συστήματος, αλλά η πιο διαδεδομένη είναι η κυστίτιδα (λοίμωξη της ουροδόχου κύστης). Οι ουρολοιμώξεις εμφανίζονται συχνότερα στις γυναίκες, καθώς η ουρήθρα είναι πιο κοντή και επομένως βρίσκεται πιο κοντά στο ορθό. Αυτό αυξάνει την πιθανότητα λοίμωξης καθώς είναι ευκολότερο για τα βακτήρια να εισέλθουν στο ουροποιητικό σύστημα. Οι νεφρικές λοιμώξεις αποτελούν μια άλλη μορφή ουρολοίμωξης. Αυτές οι λοιμώξεις είναι πιο σοβαρές από την κυστίτιδα, αλλά πολύ λιγότερο συχνές. Τα συμπτώματα της κυστίτιδας μπορεί να περιλαμβάνουν συχνή ούρηση ή ανάγκη για ούρηση ενώ η κύστη είναι άδεια, αίσθημα καύσου ή πόνο κατά την ούρηση, αίμα στα ούρα και πίεση ή επώδυνες μυικές συσπάσεις στην κατώτερη περιοχή της κοιλιακής χώρας. Τα συμπτώματα των νεφρικών λοιμώξεων είναι πολύ διαφορετικά. Αυτά μπορεί να περιλαμβάνουν ρίγη, πυρετό, ναυτία και έμετο και πόνο στην οσφυϊκή χώρα¹. Υπάρχουν πολλά είδη βακτηρίων που προκαλούν ουρολοίμωξη, αλλά τα τρία είδη που καλύπτονται σε αυτό το αρχείο είναι τα *Escherichia coli*, *Staphylococcus aureus* και *Proteus mirabilis*.

Λόγω της σοβαρότητας των ουρολοιμώξεων σε ένα τεράστιο αριθμό ασθενών, είναι επομένως πολύ σημαντικό να υπάρχει η δυνατότητα απομόνωσης και ταυτοποίησης των κοινών μικροοργανισμών που προκαλούν ουρολοίμωξη από δείγματα ούρων. Η έγκαιρη διάγνωση είναι ζωτικής σημασίας για την πρόληψη και τη μείωση των λοιμώξεων του ουροποιητικού συστήματος.

Αρχή της μεθόδου

Το άγαρ ουρίας είναι ένα ελαφρώς ρυθμισμένο μέσο που περιέχει ουρία και κόκκινο της φαινόλης που είναι δείκτης του pH. Όταν οι οργανισμοί χρησιμοποιούν ουρία,

σχηματίζεται αμμωνία η οποία μετατρέπεται το μέσο σε αλκαλικό. Ο δείκτης, το κόκκινο της φαινόλης, αλλάζει το χρώμα του μέσου από ανοιχτό κίτρινο σε ροζ-κόκκινο σε αλκαλικό περιβάλλον. Η ζελατινή πεπτόνης προάγει την ταχεία ανάπτυξη των εντερικών gram-αρνητικών βακίλων, επιτρέποντας μείωση του χρόνου επώασης. Η δεξτρώζη ενεργοποιεί τη δράση της ουρεάσης σε εκείνους τους μικροοργανισμούς που υδρολύουν την ουρία αργά.

Τυπική σύνθεση

	γραμμάρια ανά λίτρο
Πεπτόνη	1,0
Γλυκόζη	1,0
Χλωριούχο νάτριο	5,0
Φωσφορικό δινάτριο	1,2
Διόξινο φωσφορικό κάλιο	0,8
Κόκκινο φαινόλης	0,012
Άγαρ	15,0

Εξωτερική εμφάνιση

Χρώμα	Πορτοκαλί
Διαύγεια	Διαφάνεια
Συμπλήρωση βάρους	30,0-3,3 g
pH	6,8 ± 0,2

Υλικά που Παρέχονται

BO0337B: 24 x φιάλες Urea Agar Slope
EB0337B: 200 x φιάλες Urea Agar Slope

Κάθε φιάλη πρέπει να χρησιμοποιείται μόνο μία φορά.

Υλικά που απαιτούνται αλλά δεν παρέχονται

- Κρίκοι ενοφθαλμισμού
- Στυλεοί
- Δοχεία συλλογής
- Επωαστήρες
- Μικροοργανισμοί ποιοτικού ελέγχου

Αποθήκευση

- Αποθηκεύστε το προϊόν στην αρχική του συσκευασία στους 2–10 °C μέχρι τη χρήση του.
- Το προϊόν μπορεί να χρησιμοποιηθεί μέχρι την ημερομηνία λήξης που αναγράφεται στην ετικέτα.
- Φυλάσσετε μακριά από το φως.
- Αφήστε το προϊόν να ισορροπήσει σε θερμοκρασία δωματίου πριν από τη χρήση.
- Μην επωάζετε πριν από τη χρήση.

Προειδοποιήσεις και προφυλάξεις

- Μόνο για in vitro διαγνωστική χρήση.
- Μόνο για επαγγελματική χρήση.
- Επιθεωρήστε τη συσκευασία του προϊόντος πριν από την πρώτη χρήση.
- Μην χρησιμοποιείτε το προϊόν εάν υπάρχει ορατή ζημιά στη συσκευασία ή στα τρυβλία.
- Μη χρησιμοποιείτε το προϊόν πέρα από την αναγραφόμενη ημερομηνία λήξης.
- Μη χρησιμοποιείτε το ιατροτεχνολογικό προϊόν εάν υπάρχουν σημάδια επιμόλυνσης.
- Μη χρησιμοποιείτε το ιατροτεχνολογικό προϊόν εάν το χρώμα έχει αλλάξει ή υπάρχουν άλλα σημάδια φθοράς.
- Είναι ευθύνη κάθε εργαστηρίου να διαχειρίζεται τα απόβλητα που παράγονται σύμφωνα με τη φύση και τον βαθμό επικινδυνότητάς τους και να τα αντιμετωπίζει ή να τα απορρίπτει σύμφωνα με τους ομοσπονδιακούς πολιτειακούς και τοπικούς ισχύοντες κανονισμούς. Οι οδηγίες πρέπει να διαβάζονται και να ακολουθούνται προσεκτικά. Αυτό περιλαμβάνει την απόρριψη χρησιμοποιημένων ή αχρησιμοποίητων αντιδραστηρίων καθώς και οποιουδήποτε άλλου

μολυσμένου υλικού μιας χρήσης, ακολουθώντας διαδικασίες για μολυσματικά ή δυνητικά μολυσματικά προϊόντα.

Ανατρέξτε στο Δελτίο Δεδομένων Ασφάλειας Υλικού (SDS) για ασφαλή χειρισμό και απόρριψη του προϊόντος στη διεύθυνση (www.thermofisher.com).

Σοβαρά Συμβάντα

Κάθε σοβαρό συμβάν που έχει προκύψει σε σχέση με το ιατροτεχνολογικό προϊόν πρέπει να αναφέρεται στον κατασκευαστή και στην σχετική ρυθμιστική αρχή του κράτους στο οποίο είναι εγκατεστημένος ο χρήστης ή/και ο ασθενής.

Συλλογή, χειρισμός και αποθήκευση δειγμάτων

Τα δείγματα θα πρέπει να συλλέγονται και να χειρίζονται σύμφωνα με τις τοπικές συνιστώμενες οδηγίες, όπως τα Πρότυπα του HB για Μικροβιολογικές Έρευνες (UK SMI) ID 16, ID 20, ID 21, ID 24,

Διαδικασία

- Αφήστε το προϊόν να ισορροπήσει σε θερμοκρασία δωματίου πριν από τη χρήση.
- Ενοφθαλμίστε και απλώστε το δείγμα πάνω στο μέσο.
- ή εναλλακτικά χρησιμοποιώντας άσηπτη τεχνική, επιλέξτε μια μεμονωμένη και καλώς απομονωμένη αποικία με μια αποστειρωμένη βελόνα ενοφθαλμισμού και τρυπήστε το άγαρ με τη βελόνα αρκετές φορές έως το κατώτατο μέρος του φιαλιδίου ή του σωλήνα
- Επανατοποθετήστε το καπάκι και σφίξτε το χαλαρά κατά την επώαση για να επιτρέψετε την ανταλλαγή αερίων
- Επώαστε το τρυπημένο άγαρ / κεκλιμένη επιφάνεια για 18 - 24 ώρες στους $37 \pm 2^\circ\text{C}$.
- Επιθεωρήστε οπτικά τις φιάλες για να αξιολογήσετε την ανάπτυξη και το χρώμα της αποικίας κάτω από επαρκή φωτισμό.

Ερμηνεία

Αρνητική ουρεάση, καμία αλλαγή χρώματος/ κίτρινο μέσο υποδηλώνει παρουσία *Escherichia coli*.

Θετική ουρεάση με ροζ κεκλιμένη επιφάνεια υποδηλώνει παρουσία *Proteus mirabilis* ή *Proteus hauseri*.

Ασθενώς θετική ουρεάση με ανοιχτό ροζ χρώμα κεκλιμένης επιφάνειας υποδηλώνει παρουσία *Klebsiella pneumoniae*.

Έλεγχος ποιότητας

Είναι ευθύνη του χρήστη να πραγματοποιήσει δοκιμές Ποιοτικού Ελέγχου λαμβάνοντας υπόψη την προβλεπόμενη χρήση του μέσου και σύμφωνα με τυχόν τοπικούς ισχύοντες κανονισμούς (συχνότητα, αριθμός στελεχών, θερμοκρασία επώασης κ.λπ.).

Η επίδοση αυτού του μέσου μπορεί να επαληθευτεί δοκιμάζοντας τα ακόλουθα στελέχη αναφοράς.

Συνθήκες επώασης: 18 - 24 ώρες στους $37 \pm 2^\circ\text{C}$

Θετικοί μάρτυρες	
Επίπεδο ενοφθαλμισμού: 10.000-100.000 cfu	
<i>Escherichia coli</i> ATCC® 25922™	Αρνητική ουρεάση, καμία χρωματική αλλαγή/ κίτρινο μέσο
<i>Proteus mirabilis</i> NCTC 10975	Θετική ουρεάση, ροζ κεκλιμένη επιφάνεια
<i>Proteus hauseri</i> ATCC® 13315™	Θετική ουρεάση, ροζ κεκλιμένη επιφάνεια

<i>Klebsiella pneumoniae</i> ATCC® 13883™	Ασθενώς θετική ουρεάση, ανοιχτό ροζ χρώμα κεκλιμένης επιφάνειας
--	---

Περιορισμοί

Η αλκαλική αντίδραση που παράγεται σε αυτό το μέσο μετά από παρατεταμένη επώαση μπορεί να μην προκαλείται από τη δράση της ουρεάσης. Ελέγξτε χρησιμοποιώντας μέσο χωρίς ουρία.

Μην θερμαίνετε ή μην θερμαίνετε εκ νέου το μέσο διότι η ουρία αποσυντίθεται πολύ εύκολα.

Για την ανίχνευση θετικού στην ουρεάση *Proteae* η αντίδραση πρέπει να διαβαστεί εντός των 2-6 ωρών από την επώαση.

Χαρακτηριστικά απόδοσης




Η ακρίβεια έχει αποδειχθεί μέσω της ανασκόπησης των δεδομένων ποιοτικού ελέγχου. Η σωστή ανίχνευση των στελεχών *Enterobacteriaceae* επιβεβαιώνεται με τη συμπερίληψη καλά χαρακτηρισμένων απομονωθέντων στελεχών στις διαδικασίες ποιοτικού ελέγχου (QC) που εκτελούνται ως μέρος της κατασκευής κάθε παρτίδας των ιατροτεχνολογικών προϊόντων. Η ακρίβεια του Urea Agar Slope (BO0337B) αποδείχθηκε από ένα συνολικό ποσοστό επιτυχίας 80% που επιτεύχθηκε για το προϊόν σε διάστημα δέκα μηνών δοκιμής (18.ΣΕΠΤ.2021–25.ΙΟΥΛ.2022, 10 παρτίδες). Η ακρίβεια του Urea Agar Slope (EB0337B) αποδείχθηκε από ένα συνολικό ποσοστό επιτυχίας 90% που επιτεύχθηκε για το προϊόν σε διάστημα εννέα μηνών δοκιμής (13.ΣΕΠΤ.2021–13.ΙΟΥΝ.2022, 10 παρτίδες). Ένα συνολικό ποσοστό επιτυχίας 80% (βάσει δεδομένων από τις πιο πρόσφατες 10 παρτίδες) παρέχει επαρκή δεδομένα για την ακρίβεια λαμβάνοντας υπόψη τον χαμηλό κίνδυνο του ιατροτεχνολογικού προϊόντος και δεν αντικατοπτρίζει το προϊόν που βρίσκεται σε κυκλοφορία για κλινική χρήση. Στα πλαίσια της φυσιολογικής διακίμανσης στη διαδικασία κατασκευής, ενδέχεται να προκύψει ένας μικρός αριθμός αστοχιών, ωστόσο, δεν συντρέχει κανένας κίνδυνος για τον τελικό χρήστη, καθώς οι παρτίδες θα υποβληθούν σε έλεγχο QC και οι αστοχίες θα εντοπιστούν, στην περίπτωση που είναι εκτός προδιαγραφών. Αυτό δείχνει ότι η απόδοση είναι αναπαραγωγίμη, πληρώνοντας έτσι τις προδιαγραφές που ορίζονται στο Σχέδιο Αξιολόγησης Απόδοσης.

Τα ιατροτεχνολογικά προϊόντα Urea Agar Slope (BO0337B και EB0337B) υποβάλλονται σε εσωτερικές δοκιμές ως μέρος της διαδικασίας ποιοτικού ελέγχου από το 1997 (BO0337B) και το 1998 (EB0337B). Για οργανισμούς στόχους, όταν χρησιμοποιείται ενοφθαλμισμό $10^4 - 10^5$ cfu από *Escherichia coli*, *Proteus mirabilis* ή *Klebsiella pneumoniae* και το ιατροτεχνολογικό προϊόν επωαστεί στους $37 \pm 2^\circ\text{C}$ για 18-24 ώρες, ο χρήστης μπορεί να αναμένει αντίδραση ουρεάσης όπως αναφέρεται σε αυτό το έγγραφο.

Βιβλιογραφία

1. Centers for Disease Control and Prevention. 2021. 'Urinary Tract Infection'. <https://www.cdc.gov/antibiotic-use/uti.html>.

Υπόμνημα συμβόλων

Σύμβολο	Ορισμός
	Αριθμός Καταλόγου
	In Vitro Διαγνωστικό Ιατροτεχνολογικό Προϊόν
	Κωδικός Παρτίδας



Για τεχνική βοήθεια, επικοινωνήστε με τον τοπικό διανομέα σας.

Πληροφορίες Αναθεώρησης

Έκδοση	Ημερομηνία τροποποιήσεων που εισήχθησαν
2.0	21-11-2023

	Όριο θερμοκρασίας
	Ημερομηνία λήξης
	Φυλάσσετε μακριά από το ηλιακό φως
	Να μην επαναχρησιμοποιείται
	Συμβουλευτείτε τις οδηγίες χρήσης ή συμβουλευτείτε τις ηλεκτρονικές οδηγίες χρήσης
	Περιέχει επαρκή αριθμό για <n> δοκιμές
	Μην το χρησιμοποιείτε εάν η συσκευασία είναι κατεστραμμένη και συμβουλευτείτε τις οδηγίες χρήσης
	Κατασκευαστής
	Εξουσιοδοτημένος αντιπρόσωπος στην Ευρωπαϊκή Κοινότητα/ Ευρωπαϊκή Ένωση
	Ευρωπαϊκή Αξιολόγηση Συμμόρφωσης
	Αξιολογήθηκε η Συμμόρφωση του Ηνωμένου Βασιλείου
	Μοναδικό αναγνωριστικό ιατροτεχνολογικού προϊόντος
	Εισαγωγέας - Υποδεικνύει την οντότητα που εισάγει το ιατροτεχνολογικό προϊόν στη συγκεκριμένη τοποθεσία. Ισχύει για την Ευρωπαϊκή Ένωση
Made in the United Kingdom	Κατασκευάζεται στο Ηνωμένο Βασίλειο



©2022 Thermo Fisher Scientific Inc. Με την επιφύλαξη παντός δικαιώματος.

Τα σήματα καταλόγου ATCC και ATCC αποτελούν εμπορικό σήμα της American Type Culture Collection.

Τα σήματα καταλόγου NCTC και NCTC αποτελούν εμπορικό σήμα του National Collection of Type Cultures.

Όλα τα άλλα εμπορικά σήματα αποτελούν ιδιοκτησία της Thermo Fisher Scientific Inc. και των θυγατρικών της.





www.thermofisher.com

Thermo
SCIENTIFIC

Kaliumdihydrogenfosfat	0,8
Fenol rød	0,012
Agar	15,0

Urea Agar Slope

NO

REF

BO0337B OG EB0337B

Tiltenkt bruk

Urea Agar Slope (BO0337B og EB0337B)-enheter er ikke-selektive medier beregnet for bruk i dyrkningsmedier. Urea Agar Slope (BO0337B og EB0337B) enheter er tiltenkt brukt til å påvise rask ureaseaktivitet til *Proteae* og ikke-rask ureaseaktivitet hos noen *Enterobacteriaceae* i avførings- og urinprøver.

Enheterne er tiltenkt brukt i en diagnostisk arbeidsflyt for å hjelpe klinikere med å bestemme potensielle behandlingsalternativer for pasienter som mistenkes å ha *Enterobacteriaceae*-infeksjoner.

Enheterne er kun for profesjonell bruk, er ikke automatiserte og er heller ikke ledsagende diagnostikk.»

Sammendrag og forklaring

Urinveisinfeksjoner er vanlige infeksjoner som oppstår når bakterier invaderer urinveiene ved å komme inn via urinrøret. Disse infeksjonene kan påvirke forskjellige deler av urinveiene, men den mest utbredte er blærebetennelse (en blæreinfeksjon). Urinveisinfeksjoner forekommer hyppigere hos kvinner, da urinrøret er kortere, og derfor nærmere endetarmen. Dette øker sannsynligheten for infeksjon da det er lettere for bakterier å komme inn i urinveiene. Nyreinfeksjoner er en annen form for urinveisinfeksjon. Disse er mer alvorlige enn blærebetennelse, men langt mindre vanlige. Symptomene på blærebetennelse kan inkludere hyppig vannlating, eller behov for å urinere mens du har en tom blære, en brennende følelse eller smerte under vannlating, blod i urinen og trykk eller kramper i nedre del av magen. Symptomene på nyreinfeksjoner er svært forskjellige. Disse kan omfatte frysninger, feber, kvalme og oppkast og smerter i korsryggen¹. Det er mange arter av bakterier som forårsaker UVI, men de tre som dekkes i denne filen er *Escherichia coli*, *Staphylococcus aureus* og *Proteus mirabilis*.

På grunn av alvorlighetsgraden av urinveisinfeksjon hos et stort antall pasienter, er det derfor svært viktig å kunne isolere og identifisere vanlige mikroorganismer som forårsaker urinveisinfeksjon fra urinprøver. Tidlig diagnose er avgjørende for å forebygge og redusere urinveisinfeksjoner.

Metodeprinsipp

Urea-agar er et lett bufret medium som inneholder urea og fenolrødt, en pH-indikator. Når organismer bruker urea, dannes ammoniakk som gjør mediet alkalisk. Indikatoren, fenolrød, endrer mediefargen fra blekgul til rosarød i et alkalisk miljø. Gelatin pepton fremmer rask vekst av enteriske gramnegative basiller, noe som tillater en reduksjon i inkubasjonstiden. Dekstrose stimulerer ureaseaktivitet i organismer som hydrolyserer urea sakte.

Typisk formel

	gram per liter
Pepton	1,0
Glukose	1,0
Natriumklorid	5,0
Dinatriumfosfat	1,2

Fysisk utseende

Farge	Oransje
Klarhet	Klar
Fyllvekt	3,0 – 3,3 g
pH	6,8 ± 0,2

Materialer som følger med

BO0337B: 24 x Urea Agar Slope-flasker
EB0337B: 200 x Urea Agar Slope-flasker

Hver flaske skal kun brukes én gang.

Materialer som er nødvendige, men som ikke følger med

- Inokuleringsøser
- Bomullspinne
- Oppsamlingsbeholdere
- Inkubatorer
- Kvalitetskontrollorganismer

Oppbevaring

- Oppbevar produktet i originalemballasjen ved 2–10 °C til det skal brukes.
- Produktet kan brukes til utløpsdatoen som er angitt på etiketten.
- Må ikke utsettes for lys.
- La produktet nå likevekt i romtemperatur før bruk.
- Ikke inkuber før bruk.

Advarsler og forholdsregler

- Kun for in vitro-diagnostisk bruk.
- Kun til profesjonell bruk.
- Inspiser produktets emballasje før første gangs bruk.
- Ikke bruk produktet hvis det er synlige skader på emballasjen eller produktet.
- Produktet må ikke brukes etter den angitte utløpsdatoen.
- Ikke bruk enheten hvis det er tegn på kontaminering.
- Ikke bruk enheten hvis fargen har endret seg eller det er andre tegn på forringelse.
- Det er hvert laboratoriums ansvar å håndtere avfall som produseres i henhold til deres natur og grad av fare, og å få det behandlet eller kastet i samsvar med eventuelle føderale, statlige og lokale gjeldende forskrifter. Instruksjonene skal leses og følges nøye. Dette inkluderer kassering av brukte eller ubrukte reagenser samt alle andre kontaminerte engangsmaterialer etter prosedyrer for smittefarlige eller potensielt smittsomme produkter.

Se sikkerhetsdatabladet (SDS) for sikker håndtering og kassering av produktet (www.thermofisher.com).

Alvorlige hendelser

Enhver alvorlig hendelse som har oppstått i forbindelse med bruk av enheten, skal rapporteres til produsenten og den relevante tilsynsmyndigheten der brukeren og/eller pasienten er etablert.

Prøveinnsamling, håndtering og oppbevaring

Prøver skal samles og håndteres i henhold til lokale

anbefalte retningslinjer, slik som UK Standards for Microbiology Investigations (UK SMI) ID 16, ID 20, ID 21, ID 24,

Prosedyre

- La produktet komme i likevekt i romtemperatur.
- Inokuler og stryk ut prøven på mediet.
- eller alternativt ved bruk av aseptisk teknikk velg en enkelt godt isolert koloni med en steril inokulerende stikk nål og stikk nålen flere ganger gjennom agaren til bunnen av hetteglasset eller røret
- Sett på hetten og stram løst til under inkubering for å tillate gassutveksling
- Inkuber den stukkede agaren /skråstilt i 18 - 24 timer ved 37 ± 2 °C.
- Inspiser flaskene visuelt for å vurdere kolonivekst og farge under god belysning.

Tolkning

Urease negativ, ingen fargeendring/ gult medium indikerer *Escherichia coli*.

Urease positiv med rosa hevinger indikerer *Proteus mirabilis* eller *Proteus hauseri*.

Svak ureasepositiv med blekrosa hevinger indikerer *Klebsiella pneumoniae*.

Kvalitetskontroll

Det er brukerens ansvar å utføre kvalitetskontrolltesting under hensyntagen til tiltenkt bruk av mediet, og i samsvar med lokale gjeldende forskrifter (frekvens, antall stammer, inkubasjonstemperatur osv.).

Ytelsen til dette mediet kan verifiseres ved å teste følgende referansestammer.

Inkubasjonsbetingelser: 18 - 24 timer ved 37 ± 2 °C

Positive kontroller	
Inokulurnivå: 10 000 – 100 000 cfu	
<i>Escherichia coli</i> ATCC® 25922™	Urease negativ, ingen fargeendring/ gul medium
<i>Proteus mirabilis</i> NCTC 10975	Urease positiv, rosa heving
<i>Proteus hauseri</i> ATCC® 13315™	Urease positiv, rosa heving
<i>Klebsiella pneumoniae</i> ATCC® 13883™	Svak ureasepositiv, blekrosa heving

Begrensninger

Den alkaliske reaksjonen som produseres i dette mediet etter langvarig inkubasjon er kanskje ikke forårsaket av ureaseaktivitet. Kontroller med medium uten urea.

Ikke varm opp eller varm opp mediet på nytt fordi urea brytes ned veldig lett.

For påvisning av urease-positiv *Proteae* må reaksjonen avleses innen de første 2-6 timene av inkubasjonen.

Ytelseegenskaper

Nøyaktighet har blitt demonstrert gjennom gjennomgang av QC-dataene. Riktig påvisning av Enterobacteriaceae bekreftes ved inkludering av et godt karakterisert isolat i kvalitetskontrollprosessene som utføres som en del av produksjonen av hvert sett av enhetene. Presisjonen til Urea Agar Slope (BO0337B) ble demonstrert ved en samlet bestått rate på 80 % oppnådd for produktet over ti måneders testing (18.SEP.2021–25.JUL.2022; 10 sett). Presisjonen












til Urea Agar Slope (EB0337B) ble demonstrert ved en total bestått rate på 90 % oppnådd for produktet i løpet av ni måneders testing (13.SEP.2021–13.JUN.2022; 10 sett). En samlet bestått-rate på 80 % (basert på data fra de siste 10 settene) gir tilstrekkelige data om nøyaktighet med tanke på den lave risikoen for enheten, og den reflekterer ikke produktet som er utgitt for klinisk bruk. Med normal variasjon i produksjonsprosessen kan det oppstå et lite antall feil, men det er ingen risiko for sluttbrukeren da sett vil bli fanget opp i kvalitetskontrollen hvis de ikke er spesifisert. Dette viser at ytelsen er reproduserbar, og oppfyller dermed spesifikasjonene i ytelseevalueringen.


Urea Agar Slope (BO0337B & EB0337B)-enheter er testet internt som en del av kvalitetskontrollprosessen siden 1997 (BO0337B) og 1998 (EB0337B). For målorganismer, ved bruk av 10^4 - 10^5 cfu inokulum av *Escherichia coli*, *Proteus mirabilis* eller *Klebsiella pneumoniae* og inkubering av enheten ved 37 ± 2 °C i 18-24 timer, kan brukeren forvente ureasereaksjon som oppført i dette dokumentet.

Bibliografi

1. Centers for Disease Control and Prevention. 2021. 'Urinary Tract Infection'. <https://www.cdc.gov/antibiotic-use/uti.html>.

Symbolforklaring

Symbol	Definisjon
	Katalognummer
	In vitro-diagnostisk medisinsk utstyr
	Settkode
	Temperaturgrense
	Utløpsdato
	Holdes unna sollys
	Må ikke brukes flere ganger
	Se bruksanvisning eller konsulter elektronisk bruksanvisning
	Inneholder tilstrekkelig mengde til <n> tester
	Ikke bruk hvis emballasjen er skadet, se bruksanvisningen
	Produsent

EC REP	Autorisert representant innen det europeiske fellesskap / EU
CE	Europeisk samsvarsvurdering
UK CA	Samsvarsvurdering for Storbritannia
UDI	Unik enhetsidentifikator
	Importør - for å angi foretaket som importerer det medisinske utstyret til stedet. Gjelder for EU
Made in the United Kingdom	Produsert i Storbritannia



©2022 Thermo Fisher Scientific Inc. Med enerett. ATCC- og ATCC-katalogmerkene er et varemerke for American Type Culture Collection. NCTC- og NCTC-katalogmerkene er et varemerke for National Collection of Type Cultures. Alle andre varemerker tilhører Thermo Fisher Scientific Inc. og dets datterselskaper.



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



For teknisk assistanse, vennligst kontakt din lokale distributør.

Revisjonsinformasjon

Versjon	Dato for innførte endringer
2,0	2023-11-21



www.thermofisher.com

Urea Agar Slope

REF **BO0337B Y EB0337B**

Uso previsto

Los dispositivos Urea Agar Slope (BO0337B y EB0337B) son medios no selectivos destinados a su uso en medios de cultivo. Los dispositivos Urea Agar Slope (BO0337B y EB0337B) están destinados al uso en la detección rápida de actividad de ureasa de la *Proteae* y la actividad no rápida de ureasa de ciertas *Enterobacteriaceae* en muestras de heces y orina.

Los dispositivos están destinados al uso en un flujo de trabajo de diagnóstico para ayudar a los médicos a determinar posibles opciones de tratamiento para pacientes con presuntas infecciones por *Enterobacteriaceae*.

Los dispositivos son exclusivamente para uso profesional, no están automatizados y no son pruebas diagnósticas complementarias.

Resumen y explicación

Las infecciones urinarias, a menudo denominadas IU, son infecciones comunes que se producen cuando las bacterias invaden las vías urinarias al ingresar a través de la uretra. Estas infecciones pueden afectar diferentes partes de las vías urinarias, pero la más frecuente es la cistitis (infección de la vejiga). Las infecciones urinarias (IU) ocurren con mayor frecuencia en las mujeres, ya que la uretra es más corta y, por lo tanto, está más cerca del recto. Esto aumenta la probabilidad de infección, ya que es más fácil que las bacterias entren en las vías urinarias. Las infecciones renales son otra forma de IU. Son más graves que la cistitis, pero mucho menos comunes. Los síntomas de la cistitis pueden incluir micción frecuente o necesidad de orinar con la vejiga vacía, sensación de ardor o dolor al orinar, sangre en la orina y presión o calambres en el área abdominal inferior. Los síntomas de las infecciones renales son muy diferentes. Estos pueden incluir escalofríos, fiebre, náuseas, vómitos y dolor en la zona lumbar¹. Hay muchas especies de bacterias que causan IU, pero las tres que se abordan en este documento son *Escherichia coli*, *Staphylococcus aureus* y *Proteus mirabilis*.

Debido a la gravedad de las IU en una gran cantidad de pacientes, es muy importante poder aislar e identificar los microorganismos comunes causantes de las IU a partir de muestras de orina. El diagnóstico precoz es fundamental para la prevención y reducción de las infecciones urinarias.

Principio del método

Urea Agar es un medio ligeramente tamponado que contiene urea y rojo fenol, un indicador de pH. Cuando los microorganismos utilizan urea, se forma amoníaco que convierte el medio en alcalino. El indicador, rojo fenol, cambia el color del medio de amarillo pálido a rojo rosado en un entorno alcalino. La peptona de gelatina favorece el crecimiento rápido de bacilos gramnegativos entéricos, lo que permite una disminución del tiempo de incubación. La dextrosa estimula la actividad de la ureasa en aquellos microorganismos que hidrolizan la urea lentamente.

Fórmula típica

	gramos por litro
Peptona	1,0
Glucosa	1,0
Cloruro de sodio	5,0
Fosfato disódico	1,2
Dihidrogenofosfato de potasio	0,8
Rojo fenólico	0,012
Agar	15,0

Apariencia física

Color	Naranja
Claridad	Transparente
Peso de relleno	3,0-3,3 g
pH	6,8 ± 0,2

Materiales suministrados

BO0337B: 24 frascos de Urea Agar Slope
EB0337B: 200 frascos de Urea Agar Slope

Cada frasco es de un solo uso exclusivamente.

Materiales necesarios pero no suministrados

- Asas de inoculación
- Hisopos
- Recipientes de recogida
- Incubadoras
- Microorganismos de control de calidad

Almacenamiento

- Almacenar el producto en su envase original a 2 °C-10 °C hasta que se vaya a utilizar.
- El producto se puede utilizar hasta la fecha de caducidad indicada en la etiqueta.
- Almacenar protegido de la luz.
- Dejar que el producto se temple a temperatura ambiente antes de usarlo.
- No incubar antes de usar.

Advertencias y precauciones

- 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 embalaje o las placas.
- 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 esté establecido 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 16, ID 20, ID 21 e ID 24.

Procedimiento

- Deje que el producto se temple a temperatura ambiente.
- Inocule y siembre la muestra en el medio.
- O bien, como alternativa, con una técnica aséptica, recoja una sola colonia bien aislada con una aguja punzante de inoculación estéril y pinche con la aguja varias veces a través del agar hasta el fondo del vial o tubo.
- Vuelva a colocar la tapa y ajústela sin apretar mientras se esté incubando para permitir el intercambio de gases.
- Incube el agar pinchado/en pendiente durante 18-24 horas a 37 ± 2 °C.
- Inspeccione visualmente los frascos para evaluar el crecimiento y el color de las colonias con una iluminación adecuada.

Interpretación

Un resultado de ureasa negativo, sin cambio de color/medio amarillo indica *Escherichia coli*.

Un resultado de ureasa positivo con pendientes rosas indica *Proteus mirabilis* o *Proteo hauseri*.

Un resultado de ureasa positivo débil con pendientes de color rosa pálido indica *Klebsiella pneumoniae*.

Control de calidad

Es responsabilidad del usuario realizar las pruebas de control de calidad teniendo en cuenta el uso previsto del medio y de acuerdo con las normativas locales aplicables (frecuencia, número de cepas, temperatura de incubación, etc.).

Es posible verificar el rendimiento de este medio probando las cepas de referencia siguientes.

Condiciones de incubación: 18-24 h a 37 °C \pm 2 °C

Controles positivos	
Nivel de inóculo: 10.000-100.000 ufc	
<i>Escherichia coli</i> ATCC® 25922™	Resultado de ureasa negativo, sin cambio de color/medio amarillo
<i>Proteus mirabilis</i> NCTC 10975	Resultado de ureasa positivo, con pendiente rosa
<i>Proteus hauseri</i> ATCC® 13315™	Resultado de ureasa positivo, con pendiente rosa
<i>Klebsiella pneumoniae</i> ATCC® 13883™	Resultado de ureasa positivo débil, con pendiente rosa pálido

Limitaciones

La reacción alcalina que se produce en este medio tras una incubación prolongada puede no estar provocada por la actividad de la ureasa. Realice una comprobación utilizando medio sin urea.

No caliente ni recaliente el medio, ya que la urea se descompone muy fácilmente.

Para la detección de *Proteae* con resultado de ureasa positivo, la reacción debe leerse dentro de las primeras 2-6 horas de incubación.

Características de rendimiento

Se ha demostrado la precisión mediante la revisión de los


datos de control de calidad. La detección correcta de cepas de Enterobacteriaceae se confirma mediante la inclusión de un aislado bien caracterizado en los procesos de control de calidad realizados como parte de la fabricación de cada lote de los dispositivos. La precisión de Urea Agar Slope (BO0337B) quedó demostrada por una tasa general de aprobación del 80 % obtenida para el producto durante diez meses de prueba (18.SEP.2021-25.JUL.2022; 10 lotes). La precisión de Urea Agar Slope (EB0337B) quedó demostrada por una tasa general de aprobación del 90 % obtenida para el producto durante nueve meses de prueba (13.SEP.2021-13.JUN.2022; 10 lotes). Una tasa general de aprobación del 80 % (basada en los datos de los 10 lotes más recientes) proporciona datos suficientes sobre la precisión teniendo en cuenta el bajo riesgo del dispositivo y no refleja el producto que se lanza para uso clínico. Con la variación normal en el proceso de fabricación, puede ocurrir un pequeño número de fallos; sin embargo, no hay riesgo para el usuario final, ya que los lotes se detectarán en el control de calidad si no cumplen con las especificaciones. Esto demuestra que el rendimiento es reproducible, cumpliendo así con las especificaciones establecidas en el plan de evaluación del rendimiento.

Los dispositivos Urea Agar Slope (BO0337B y EB0337B) se han probado internamente como parte del proceso de control de calidad desde 1997 (BO0337B) y 1998 (EB0337B). En el caso de los microorganismos diana, al usar un inóculo de 10^4 - 10^5 ufc de *Escherichia coli*, *Proteus mirabilis* o *Klebsiella pneumoniae* e incubar el dispositivo a 37 °C \pm 2 °C durante 18-24 horas, el usuario puede esperar una reacción a la ureasa como la que se indica en este documento.

Bibliografía

1. Centers for Disease Control and Prevention. 2021. 'Urinary Tract Infection'. <https://www.cdc.gov/antibiotic-use/uti.html>.

Legenda de símbolos

Símbolo	Definición
	Numero de catalogo
	Producto sanitario para diagnóstico in vitro
	Código de lote
	Límite de temperatura
	Fecha de caducidad
	Mantener alejado de la luz solar
	No reutilizar
	Consulte las instrucciones de uso o consulte 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 la conformidad para el Reino Unido
	Identificador único de dispositivo
	Importador: Indicación de la entidad que importa el dispositivo médico a la ubicación local. Aplicable a la Unión Europea.
	Hecho en el Reino Unido



© 2022 Thermo Fisher Scientific Inc. Todos los derechos reservados.

ATCC y las marcas del catálogo de ATCC son marcas registradas de American Type Culture Collection.

NCTC y las marcas del catálogo de NCTC son marcas comerciales de National Collection of Type Cultures.

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



Oxoid Limited, Wade Road, Basingstoke,
RG24 8PW, Inglaterra



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

Información de revisiones

Versión	Fecha de las modificaciones introducidas
2.0	2023-11-21