


www.thermofisher.com

Nutrient Agar Slope

REF BO0336B, BO0336E, EB0336B

EN

Intended Use

Nutrient Agar Slope (BO0336B, BO0336E and EB0336B) devices are non-selective media intended to be used for the maintenance of microorganisms isolated from other culture media products previously inoculated using clinical samples (e.g. urine and wound).

Nutrient Agar Slope (BO0336B, BO0336E and EB0336B) devices are intended to be used in a diagnostic workflow to support the clinician for the maintenance of bacteria from clinical samples. The devices are for professional use only, are not automated and nor are they companion diagnostics.

Summary and Explanation

Nonfastidious microorganisms are organisms that grow and replicate without special nutritional supplements or conditions. They can synthesize all organic molecules required for their growth. Nonfastidious microorganisms can grow in nutrient sparse environments and can survive in a wider temperature range. This groups of bacteria are involved in causing illnesses and infections in humans and includes *E. coli* and *S. aureus*^{1,2,3,4,5}.

The first step in processing a clinical specimen in order to isolate organisms in the laboratory is the inoculation of culture medium. Cell culture is the most commonly used method for cultivating and isolating intracellular bacteria from clinical samples. There are two possible approaches. First, culture conditions can be standardised to allow for the efficient growth of a diverse range of bacteria. The second strategy consists isolating a previously known bacterium and exposing it to specially adapted culture conditions in order to optimise bacterial growth. Non-selective culture media contain no inhibitors and should allow the growth of the majority of microorganisms found in clinical samples⁶.

Principle of Method

Nutrient Agar Slope (BO0336B, BO0336E and EB0336B) devices are non-selective media intended to be used for the maintenance of microorganisms isolated from other culture media products previously inoculated using clinical samples. Nutrient Agar Slope (BO0336B, BO0336E and EB0336B) devices contain peptone and yeast extract to provide a source of nitrogen, essential vitamins and amino acids necessary for the growth of bacteria. Starch is added to absorb any toxic metabolites and sodium chloride maintains osmotic balance. Agar slopes allow bacteria to grow on a larger surface area in a tube and are suitable for the maintenance of bacterial cultures.

Typical Formula

	g/L
Special peptone	23.0
Starch	1.0
Sodium chloride	5.0
Agar	10.0

* Adjusted as required to meet performance standards

Physical Appearance

Colour	Straw
Clarity	Clear
pH	7.3 ± 0.2
Fill weight: BO0336B	3.0g - 3.3g
BO0336E	10.0 – 10.3g
EB0336B	3.0 – 3.3g

Materials Provided

BO0336B – 24 x 3ml nutrient agar in bijous.....	24	Σ
BO0336E – 24 x 10ml nutrient agar in universals.....	24	Σ
EB0336B – 200 x 3ml nutrient agar in bijous.....	200	Σ

Materials Required but Not Supplied

- 1) Inoculating loops
- 2) Swabs
- 3) Collection containers
- 4) Incubator
- 5) Quality control organisms

Storage

- Store product in its original packaging at 2-25°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 bottles.
- 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) Q5.

Procedure

- (1) Allow product to equilibrate to room temperature.
- (2) 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
- (3) Replace the cap and tighten loosely when incubating to allow gas exchange
- (4) Incubate the stabbed agar / slant 18 hours at 37°C.
- (5) Visually inspect plates to assess colony growth and colour under good lighting.

Interpretation

Agar slopes allow bacteria to grow on a larger surface area in a tube and are suitable for the maintenance of bacterial cultures.

1-2 mm cream colonies indicate the growth of *Staphylococcus aureus* ATCC® 25923, and 1-2 mm cream colonies indicate the growth of *Escherichia coli* ATCC® 25922.

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 hours @ 37°C

This media can be tested with the following strains:

Positive Controls	
Colony count is ≥ 50% of the control medium count. Inoculum level 10-100 cfu	
<i>Staphylococcus aureus</i> ATCC® 25923	1-2 mm cream colonies
<i>Escherichia coli</i> ATCC® 25922	1-2 mm cream colonies

Performance Characteristics

Accuracy has been demonstrated through review of the QC data. Maintenance of nonfastidious microorganisms 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 Nutrient Agar Slope (BO0336B) devices was demonstrated by an overall pass rate of 100% obtained for the product over 11 months of testing (17.04.2021 – 02.03.2022; 10 batches), the precision of Nutrient Agar Slope (BO0336E) devices was demonstrated by an overall pass rate of 100% obtained for the product over 29 months of testing (23.07.2019 – 07.01.2022; 10 batches) and the precision of Nutrient Agar Slope (EB0336B) devices was demonstrated by an overall pass rate of 100% obtained for the product over two months of testing (03.11.2021 – 05.01.2022; 10 batches).

These media devices have been tested in-house as part of the QC process since they were first manufactured; Nutrient Agar Slope (BO0336B, BO0336E and EB0336B) devices have been on the market for at least 20 years. When using 10 - 100 cfu inoculum of *Escherichia coli* and *Staphylococcus aureus* and incubating the device at 37°C for 18 hours, the user can maintain organisms with colony size and morphology as listed in this document.

Limitations

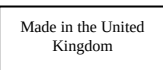
Nutrient Agar Slope (BO0336B, BO0336E and EB0336B) devices are non-selective media intended to be used for the maintenance of microorganisms isolated from other culture media products previously inoculated using clinical samples, and there are no expectations of specimens negative for the assay but maintain nonfastidious microorganisms in a condition that can mimic the test condition.

Bibliography

1. World Health Organization (WHO). 2017. 'Microbial Fact Sheets: Guidelines for drinking-water quality'. https://www.who.int/water_sanitation_health/publications/gdwq4-with-add1-chap11.pdf.
2. Public Health England. 2015. 'Identification of *Enterobacteriaceae*'. UK Standards for Microbiology Investigations ID 16 (4). https://assets.publishing.service.gov.uk/government/uploads/system/uploads/attachment_data/file/423601/ID_16i4.pdf
3. Public Health England. 2015. 'Identification of Vero cytotoxin-producing *Escherichia coli* including *Escherichia coli* O157'. UK Standards for Microbiology Investigations ID 22 (4). https://assets.publishing.service.gov.uk/government/uploads/system/uploads/attachment_data/file/436870/ID_22i4.pdf
4. Public Health England. 2020. 'Identification of *Staphylococcus* species, *Micrococcus* species and *Rothia* species'. UK Standards for Microbiology Investigations ID 07 (4). https://assets.publishing.service.gov.uk/government/uploads/system/uploads/attachment_data/file/887570/UK_SMI_ID_07i4.pdf
5. Barton, Michelle, Michael Hawkes, Dorothy Moore, John Conly, Lindsay Nicolle, Upton Allen, Nora Boyd, et al. 2006. 'Guidelines for the Prevention and Management of Community-Associated Methicillin- Resistant *Staphylococcus aureus*: A Perspective for Canadian Health Care Practitioners.' *Canadian Journal of Infectious Diseases and Medical Microbiology* 17 (c): 4C-24C. <https://doi.org/10.1155/2006/971352>
6. The Australian Society for Microbiology. 2012. "Guidelines for Assuring Quality of Medical Microbiological Culture Media: Culture Media Special Interest Group: The Australian Society for Microbiology: bringing Microbiologists together: 1–32. <https://static1.squarespace.com/static/5c184125b98a781241e53ab6/t/5db2996892f5c03a7ff21a0f/1571985772520/Guidelines+for+the+Quality+Assurance+of+Medical+Microbiological+culture+media+2nd+edition+July+2012.pdf>

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



©2025 Thermo Fisher Scientific Inc. All rights reserved. ATCC and ATCC catalogue marks are a trademark of American Type Culture Collection. 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.

Revision Information

Version	Date of issue and modifications introduced
2.0	10 June 2025 Minor formatting changes


www.thermofisher.com

Nutrient Agar Slope

REF BO0336B, BO0336E, EB0336B

ES

Uso previsto

Los dispositivos Nutrient Agar Slope (BO0336B, BO0336E and EB0336B) son medios no selectivos que se utilizan para el mantenimiento de microorganismos aislados de otros productos de medios de cultivo inoculados previamente con muestras clínicas (p. ej., orina y heridas).

Los dispositivos Nutrient Agar Slope (BO0336B, BO0336E and EB0336B) están diseñados para el uso en un flujo de trabajo de diagnóstico para ayudar al médico en el mantenimiento de bacterias procedentes de muestras clínicas. Los dispositivos son exclusivamente para uso profesional, no están automatizados y no son un diagnóstico complementario.

Resumen y explicación

Los microorganismos no exigentes son organismos que crecen y se multiplican sin suplementos nutritivos ni condiciones especiales. Pueden sintetizar todas las moléculas orgánicas necesarias para su crecimiento. Los microorganismos no exigentes pueden crecer en ambientes con escasez de nutrientes y sobrevivir en un intervalo de temperaturas más amplio. Este grupo de bacterias están involucrados en enfermedades e infecciones en humanos e incluyen *E. coli* y *S. aureus*^{1,2,3,4,5}.

El primer paso en el procesamiento de una muestra clínica para aislar organismos en el laboratorio es inocular el medio de cultivo. El cultivo celular es el método más habitual para cultivar y aislar bacterias intracelulares procedentes de muestras clínicas. Hay dos enfoques posibles. El primero consiste en estandarizar las condiciones de cultivo para permitir el crecimiento eficiente de una amplia gama de bacterias. La segunda estrategia consiste en aislar una bacteria conocida previamente y exponerla a condiciones de cultivo adaptadas especialmente para optimizar el crecimiento bacteriano. Los medios de cultivo no selectivos no contienen ningún inhibidor y deben permitir el crecimiento de la mayoría de los microorganismos que se encuentran en las muestras clínicas⁶.

Principio del método

Los dispositivos Nutrient Agar Slope (BO0336B, BO0336E and EB0336B) son medios no selectivos que se utilizan para el mantenimiento de microorganismos aislados de otros productos de medios de cultivo inoculados previamente con muestras clínicas. Los dispositivos Nutrient Agar Slope (BO0336B, BO0336E and EB0336B) contienen peptona y extracto de levadura para proporcionar una fuente de nitrógeno, vitaminas esenciales y aminoácidos necesarios para el crecimiento de bacterias. Se añade almidón para absorber cualquier metabolito tóxico y el cloruro de sodio mantiene el equilibrio osmótico. El agar en pendiente permite que las bacterias crezcan en una superficie mayor en un tubo y resulta adecuado para mantener cultivos bacterianos.

Fórmula típica

	g/l
Peptona especial	23,0
Almidón	1,0
Cloruro de sodio	5,0
Agar	10,0

* Ajustado según sea necesario para cumplir con los estándares de rendimiento

Apariencia física

Color	Pajizo
Claridad	Transparente
pH	7,3 ± 0,2
Peso de relleno: BO0336B	3,0 g-3,3 g
BO0336E	10,0-10,3 g
EB0336B	3,0-3,3 g

Materiales suministrados

BO0336B – 24 x 3 ml agar nutritivo en bijou.....	24	▽
BO0336E – 24 x 10 ml agar nutritivo en universal.....	24	▽
EB0336B – 200 x 3 ml agar nutritivo en bijou.....	200	▽

Materiales necesarios pero no suministrados

- 1) Asas de inoculación
- 2) Hisopos

- 3) Recipientes de recogida
- 4) Incubadora
- 5) Organismos de control de calidad

Almacenamiento

- Almacenar el producto en su envase original a 2 °C-25 °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.
- Deje 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 los frascos.
- 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) Q5.

Procedimiento

- (1) Deje que el producto se temple a temperatura ambiente.
- (2) 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.
- (3) Vuelva a colocar la tapa y ajústela sin apretar mientras se esté incubando para permitir el intercambio de gases.
- (4) Incubar el agar pinchado/en pendiente durante 18 horas a 37 °C ± 2 °C.
- (5) Inspeccione visualmente las placas para evaluar el crecimiento y el color de las colonias con una iluminación adecuada.

Interpretación

El agar en pendiente permite que las bacterias crezcan en una superficie mayor en un tubo y resulta adecuado para mantener cultivos bacterianos.

La presencia de colonias de color crema de 1-2 mm indica el crecimiento de *Staphylococcus aureus* ATCC® 25923 y la de colonias de color crema de 1-2 mm indica el crecimiento de *Escherichia coli* ATCC® 25922.

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 h a 37 °C ± 2 °C, aeróbica. Este medio se puede probar con las cepas siguientes:

Controles positivos	
El recuento de colonias es ≥50 % del recuento del medio de control.	
Nivel de inóculo: de 10 a 100 UFC	
<i>Staphylococcus aureus</i> ATCC® 25923	Colonias de 1-2 mm de color crema
<i>Escherichia coli</i> ATCC® 25922	Colonias de 1-2 mm de color crema

Características de rendimiento

Se ha demostrado la precisión mediante la revisión de los datos de control de calidad. EL mantenimiento de microorganismos no exigentes se confirma mediante la inclusión de aislados bien caracterizados en los procesos de control de calidad realizados como parte de la fabricación de cada lote de dispositivos.

La precisión de los dispositivos Nutrient Agar Slope (BO0336B) se demostró mediante una tasa global de corrección del 100 %

obtenida para el producto durante 11 mes de prueba (17.04.2021 – 02.03.2022; 10 lotes); la precisión de los dispositivos Nutrient Agar Slope (BO0336E) se demostró mediante una tasa global de corrección del 100 % obtenida para el producto durante 29 mes de prueba (23.07.2019 – 07.01.2022; 10 lotes) y la precisión de los dispositivos de Nutrient Agar Slope (EB0336B) se demostró mediante una tasa global de corrección del 100 % obtenida para el producto durante dos meses de prueba (03.11.2021 – 05.01.2022; 10 lotes).

Los dispositivos se prueban internamente como parte del proceso de control de calidad. En el caso de los organismos objetivo, al utilizar un inóculo de 10-100 ufc de *Escherichia coli* o *Staphylococcus aureus* e incubar el dispositivo a 37 °C ± 2 °C durante 18 horas, es posible recuperar organismos con la morfología adecuada tal como se describe en este documento.

Limitaciones

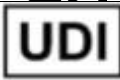
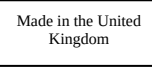
Los dispositivos de Nutrient Agar Slope (BO0336B, BO0336E and EB0336B) son medios no selectivos que se utilizan para mantener microorganismos aislados en otros productos de medios de cultivo previamente inoculados utilizando muestras clínicas, y no hay expectativas de que las muestras sean negativas para la prueba, pero mantienen microorganismos no exigentes en condiciones que pueden imitar las de la prueba.

Bibliografía

1. World Health Organization (WHO). 2017. 'Microbial Fact Sheets: Guidelines for drinking-water quality'. https://www.who.int/water_sanitation_health/publications/gdwg4-with-add1-chap11.pdf.
2. Public Health England. 2015a. 'Identification of *Enterobacteriaceae*'. UK Standards for Microbiology Investigations ID 16 (4). https://assets.publishing.service.gov.uk/government/uploads/system/uploads/attachment_data/file/423601/ID_16i4.pdf
3. Public Health England. 2015b. 'Identification of Vero cytotoxin-producing *Escherichia coli* including *Escherichia coli* O157'. UK Standards for Microbiology Investigations ID 22 (4). https://assets.publishing.service.gov.uk/government/uploads/system/uploads/attachment_data/file/436870/ID_22i4.pdf
4. Public Health England. 2020a. 'Identification of *Staphylococcus* species, *Micrococcus* species and *Rothia* species'. UK Standards for Microbiology Investigations ID 07 (4). https://assets.publishing.service.gov.uk/government/uploads/system/uploads/attachment_data/file/887570/UK_SMI_ID_07i4.pdf
5. Barton, Michelle, Michael Hawkes, Dorothy Moore, John Conly, Lindsay Nicolle, Upton Allen, Nora Boyd, et al. 2006. 'Guidelines for the Prevention and Management of Community-Associated Methicillin-Resistant *Staphylococcus aureus*: A Perspective for Canadian Health Care Practitioners.' *Canadian Journal of Infectious Diseases and Medical Microbiology* 17 (c): 4C-24C. <https://doi.org/10.1155/2006/971352>
6. The Australian Society for Microbiology. 2012. "Guidelines for Assuring Quality of Medical Microbiological Culture Media: Culture Media Special Interest Group: The Australian Society for Microbiology: bringing Microbiologists together: 1–32. <https://static1.squarespace.com/static/5c184125b98a781241e53ab6/t/5db2996892f5c03a7ff21a0f/1571985772520/Guidelines+for+the+Quality+Assurance+of+Medical+Microbiological+culture+media+2nd+edition+July+2012.pdf>

Leyenda de símbolos

Symbol	Definition
	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
	Consultar las instrucciones de uso o consultar 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



©2025 Thermo Fisher Scientific Inc. Reservados todos los derechos. ATCC y las marcas del catálogo de ATCC son marcas registradas de American Type Culture Collection.



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



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

Información de revisiones

Versión	Fecha de publicación y modificaciones introducidas
2.0	10 Junio 2025 Pequeños cambios de formato