



MicroTest™ M6™ Multi-Microbe Media

EN

REF R12530 M6 1.5mL	Σ 72
REF R12533 M6 3mL	Σ 72
REF R12557 M6 Traditional Swab Kit 1.5mL	Σ 100
REF R12560 M6 Traditional Flocked Swab Kit 1.5mL	Σ 100
REF R12561 M6 MicroTip Flock Swab Kit 1.5mL	Σ 100
REF R12562 M6 Universal Flocked Swab Kit 1.5mL	Σ 100
REF R12563 M6 Universal Flocked Swab Kit 3mL	Σ 100
REF R12568 M6 Traditional Flocked Swab Kit 3mL	Σ 100
REF R12569 M6 MicroTip Flocked Swab Kit 3mL	Σ 100
REF R12582 M6 Universal Swab Kit 1.5mL	Σ 100

INTENDED USE

Remel MicroTest™ M6™ is a liquid medium recommended for the transport of clinical specimens to the laboratory for microbiological procedures for viruses, chlamydiae, and mycoplasmas, which include *Mycoplasma* and *Ureaplasma* species. The device is used in a diagnostic workflow to aid clinicians in the treatment options for patients suspected of having viral and bacterial infections

The device is not automated, is for professional use only and is not a companion diagnostic.

SUMMARY AND EXPLANATION

Viruses, chlamydiae, and mycoplasmas are susceptible to adverse environmental conditions and require specific transport media to maintain viability (and infectivity) while in transit to the laboratory.^{1,2} M6™ is ideal for this purpose because it contains protein for stabilization, antibiotics to minimize bacterial and fungal contamination, and a buffer to maintain a neutral pH.³

PRINCIPLE

M6™ medium consists of modified Hank's balanced salt solution supplemented with bovine serum albumin, gelatin, sucrose, and glutamic acid. The pH is buffered with HEPES buffer and phenol red is the indicator. Vancomycin, amphotericin B, and colistin are selective agents which inhibit growth of competing bacteria and yeast. The medium is isotonic and non-toxic to mammalian host cells enabling it to be used in shell vial procedures. Cryoprotectants are added to ensure organism viability during freeze-thaw cycles. M6™ does not contain whole bovine serum which has been shown to inhibit attachment and growth of myxoviruses and paramyxoviruses.

REAGENTS (CLASSICAL FORMULA)*

Reactive Ingredients:	
Hank's Balanced Salts	HEPES Buffer
Bovine Serum Albumin	Phenol Red
Gelatin	Vancomycin
Sucrose	Amphotericin B
L-Glutamic Acid	Colistin

pH 7.3 ± 0.2 @ 25°C

*Adjusted as required to meet performance standards.

PRODUCT VARIANTS AND PACKAGING CONFIGURATIONS

MicroTest™ M6™ media tubes are supplied in screw-cap tubes containing either 1.5mL or 3mL of transport medium plus glass beads (unless stated differently). Swabs with several applicator-shaft options are available to facilitate collection of specimens from various patient sites. Refer to individual product descriptions in the Remel catalog or visit www.thermofisher.com for specific information about the materials supplied.

Tubes may be supplied alone or with the following specimen collection options:

Swab Type	Product Code
Polyester Tip with Plastic Shaft*	R12557
Regular Flocked Swab	R12560, R12562, R12563, R12568
Micro-Tip Flocked Swab	R12561, R12562, R12563, R12569
Polyester Swab with Plastic Applicator	R12582

PRECAUTIONS

This product is for *in vitro* diagnostic use and should be used by properly trained individuals. Precautions should be taken against the dangers of microbiological hazards by properly sterilizing specimens, containers, and media after use. Directions should be read and followed carefully. Please refer to the Safety Data Sheet (SDS) on company website and product labeling for information on potentially hazardous components.

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

In the event of malfunction, do not use the device.

STORAGE

This product is ready for use and no further preparation is necessary. Store product in its original container at 2-30°C until used. Do not overheat. Do not incubate prior to use. Improper storage will result in a loss of antimicrobial activity.

PRODUCT DETERIORATION

This product should not be used if (1) there is evidence of contamination, (2) there is evidence of leakage, (3) the color has changed from light pink, (4) the expiration date has passed, or (5) there are other signs of deterioration.

SPECIMEN COLLECTION, STORAGE, AND TRANSPORT

Specimens should be collected and handled following recommended guidelines.³⁻⁵ To maintain optimum viability, transport specimens to the laboratory as soon as possible. Although M6™ can maintain even fragile organisms for relatively long periods of time at room temperature, specimens should be refrigerated at 2-8°C or kept on wet ice (or equivalent) following collection and while in transit.

If there will be a delay in processing, freeze specimens at or below -70°C and transport on dry ice. Ship specimens in compliance with federal, state, and hospital guidelines. Specimens should be processed as soon as possible after being received in the laboratory. Dispose of all parts in accordance with national, federal and state laws and hospital and laboratory guidelines.

MATERIALS REQUIRED BUT NOT SUPPLIED

(1) Syringes, needles, (2) Sterile forceps, (3) Supplemental media, (4) Inoculating loop, loop sterilization device, (5) Incubators, alternative environmental systems, (6) Quality control organisms, (7) Materials required for organism identification. Refer to appropriate references for necessary equipment required in specimen collection.³⁻⁵

PROCEDURE

Proper specimen collection is critical for successful isolation and identification of infectious organisms. Specimens should be collected as soon as possible after onset of disease.

1. Aseptically remove cap from vial.
2. Insert swab into medium.
3. Break swab shaft evenly at the scored line. Use sterile pair of scissors if additional trimming is needed
4. Replace vial cap and close tightly.
5. Label with appropriate patient information.
6. Send to the laboratory for processing with minimal delay.

INTERPRETATION

This transport medium serves as a vehicle for maintaining organism viability while transporting the specimen to the laboratory.

QUALITY CONTROL

All lot numbers of MicroTest™ M6™ have been tested for microbial contamination, toxicity to host cells, ability to maintain viability of desired agents (listed below) and have been found to be acceptable. Testing of control organisms should be performed in accordance with established laboratory quality control procedures. If aberrant quality control results are noted, patient results should not be reported.

CONTROL

Herpes simplex type 2 ATCC® VR-734
 Respiratory syncytial virus ATCC® VR-1540
 Cytomegalovirus ATCC® VR-977
Chlamydia trachomatis ATCC® VR-880
Mycoplasma pneumoniae ATCC® 15531
Mycoplasma hominis ATCC® 23114
Ureaplasma urealyticum ATCC® 27618

LIMITATIONS

- Specimens should be handled aseptically.
- Condition, timing, and volume of specimen collected for culture can significantly impact culture and detection results. Follow recommended guidelines for specimen collection.³⁻⁵
- M6™ is intended for use as a collection and transport medium for viral, *Chlamydia*, *Ureaplasma*, and *Mycoplasma* agents only. This medium can serve as a cryoprotectant for clinical viruses, including cytomegalovirus and varicella-zoster virus.
- Because calcium alginate swabs are toxic for many enveloped viruses and may interfere with fluorescent antibody tests, they should not be used for specimen collection. Wooden shaft swabs may contain toxins and formaldehydes and should not be used. Cotton- or polyester-tipped swabs are suitable when specimen collection by a swab is appropriate.
- Specimens for isolation of mycoplasmas should be processed within 48 hours of specimen collection.
- The bag we provide with the kit in is not intended to be used as a containment device for patient samples. Please refer to and follow hospital or clinic guidelines for transporting samples.

EXPECTED VALUES

Results obtained will largely depend on proper and adequate specimen collection, as well as on timely transport and processing in the laboratory.

PERFORMANCE CHARACTERISTICS













M6™ transport medium was compared to commercial and standard transport media routinely used for the transport and maintenance of viral, *Chlamydia*, *Ureaplasma*, and *Mycoplasma* agents. The results were equivalent or superior to the media in the comparison. The percentage of recovery for each agent when stored at 2-8°C was as follows:

ORGANISM	8 HOURS	24 HOURS	48 HOURS
Respiratory Syncytial Virus	75%	64%	58%
Herpes Simplex Type 1 Mcntyre		88%	49%
Cytomegalovirus (AD 169)		58%	27%
Influenza A H3N2 Shanghai 87		87%	74%
<i>Chlamydia trachomatis</i>	82%	43%	33%
<i>Mycoplasma pneumoniae</i>	Recovered	Recovered	Recovered
<i>Mycoplasma hominis</i>	Recovered	Recovered	Recovered
<i>Ureaplasma urealyticum</i>	Recovered	Recovered	Recovered

BIBLIOGRAPHY

- Isenberg, H.D. 2004. Clinical Microbiology Procedures Handbook. 2nd ed., Vol. 3. ASM Press, Washington, D.C.
- Mahony, J.B. and M.A. Chernesky. 1985. J. Clin. Microbiol. 22:865-867.
- Miller, J.M. 1999. A Guide to Specimen Management in Clinical Microbiology. 2nd ed. ASM Press, Washington, D.C.
- Murray, P.R., E.J. Baron, J.H. Jorgensen, M.L. Landry, and M.A. Tenover. 2007. Manual of Clinical Microbiology. 9th ed. ASM Press, Washington, D.C.
- Forbes, B.A., D.F. Sahm, and A.S. Weissfeld. 2007. Bailey and Scott's Diagnostic Microbiology. 12th ed. Mosby Elsevier, St. Louis, MO.

Symbol Legend

	Catalog Number
	In Vitro Diagnostic Medical Device
	Batch Code (Lot Number)
	Consult Instructions for Use (IFU)
	Temperature Limitation (Storage Temp.)
	Use By (Expiration Date)
	Contains sufficient for <n> tests
	Do not use if packaging damaged
	Authorized European Representative
	Manufacturer
	UK Conformity Assessment
	European Conformity Assessment



Remel Inc.
 12076 Santa Fe Trail Drive
 Lenexa, KS 66215, USA
www.thermofisher.com/microbiology
 (800) 255-6730
 International: (913) 888-0939



For technical information contact your local distributor

©2022 Thermo Fisher Scientific Inc. All rights reserved.

MicroTest™ and M6™ are trademarks of Thermo Fisher Scientific and its subsidiaries
 ATCC and ATCC catalogue marks are a trademark of American Type Culture Collection.

IFU 12530, Revised 2022-06

Printed in U.S.A.