

## Technical Data Sheet

# **Tryptic Soy Agar + LTHThio Contact -ICR**

Ordering number: 1.46797.0020 / 1.46797.0200

Tryptic Soy Agar + LTHThio Contact- ICR is designed for the determination of the total microbial count on dry, sanitized surfaces and personnel in Isolators and Clean Rooms.

Ten contact plates each with a diameter of 55 mm are triple-bagged in transparent, hydrogen peroxide impermeable sleeves. The product is gamma-irradiated in the final packaging at a dose of 9-20 kGy. The sleeves consist of polypropylene with a barrier of PE-EVOH-PE.

The formulation of the basic medium (Soybean-Casein Digest Agar) is prepared according to the recommendations of the current European, Japanese and United States Pharmacopoeia (EP, 2.6.12.; JP, 4.05 and USP, 61) and supplemented with neutralizers.

Further plate designs are available:

- TSA LTHThio Contact – ICR+ (article number 146783): 55 mm lockable contact plates, triple-bagged, gamma-irradiated; intended for microbial monitoring of dry, sanitized surfaces and personnel in cleanrooms and isolators. The plate design allows aerobic, microaerophilic and anaerobic incubation.
- TSA + LTHThio Sedi. - ICR (article number 146786): 90 mm settle plates, triple-bagged, gamma-irradiated; intended for microbial monitoring of air (passive and active) and personnel in cleanrooms and isolators. The plate design allows aerobic incubation only.
- TSA + LTHThio Sedi.- ICR+ (article number 146787): 90 mm lockable settle plates, triple-bagged, gamma-irradiated; intended for viable air monitoring (passive and active) and personnel testing in cleanrooms and isolators. The plate design allows aerobic, microaerophilic and anaerobic incubation.

## **Mode of Action**

Tryptic Soy Agar (TSA, Soybean Casein Digest Agar) is a complex medium for cultivation and isolation of a wide range of bacteria, yeasts and molds. The medium is supplemented with pyruvate in order to provide an efficient neutralization of hydrogen peroxide for use in isolators. Internal studies confirmed the neutralization efficiency of the neutralizers lecithin, polysorbate

(Tween®) 80, histidine and sodium thiosulfate for disinfectants containing the following active agents:

- Alcohol (70 % ethanol or isopropyl alcohol)
- Aldehyde
- Dichloroisocyanurate
- Glucoprotamine
- Hydrogen Peroxide
- Peracetic acid
- Phenols (low pH value)
- Low concentrated quaternary ammonium compounds

The neutralizing efficiency towards residues of disinfectants in use should be validated at the application site. For neutralization of high concentrated quaternary ammonium compounds and/or polyhexamethylene biguanides the use of Neutralizer A Contact Plate is recommended (article number 146697).

### Typical Composition

Ingredient	Amount per liter
Casein Peptone	15 g/l
Soy Peptone	5 g/l
NaCl	5 g/l
Polysorbate (Tween®) 80	5 ml/l
Lecithin	0.7 g/l
Histidine	0,5 g/l
Sodium thiosulfate	0.05 g/l
Agar	15 g/l

The appearance of the medium is clear and yellowish. The pH value is in the range of 7.1-7.5. The medium can be adjusted and/or supplemented according to the performance criteria required.

### Application and Interpretation

The plates are introduced into cleanrooms grade A or B by removing one bag in each material lock. For use in isolators the inner bag has a hole in the sealing to hang up the bag during decontamination. Do not leave plates which are unprotected (unwrapped) in an isolator during decontamination.

Each plate is provided with a label including a data matrix code for paperless plate identification. The code consists of a two-dimensional 20-digit serial number, which harbors the following information: Digits 1-3: here code 802 (corresponds to article 146797); digits 4-9: lot number; digits 10-14: batch specific individual number; digits 15-20: expiration date (YY/MM/DD).

Please check each agar plate before using it on sterility and pay attention to aseptic handling in order to avoid false positive results.

According to ISO 14698 the plates are opened and the agar surface is pressed on the dry surface to be tested for some seconds with a steady pressure. Similar recommendations are included in the PDA technical report No.13. Afterwards the plates are closed and transferred to an incubator. To protect the plates from secondary contamination during transport and incubation outside of the cleanroom zone, sterile transport bags (article number 146509) may be used. Residues of culture medium should be removed from the surface after sampling.

Merck, Millipore, the Vibrant M, Sigma-Aldrich are trademarks of Merck KGaA, Darmstadt, Germany or its affiliates. . All other trademarks are the property of their respective owners. [merckgroup.com/life-science](http://merckgroup.com/life-science)  
We provide information and advice to our customers on application technologies and regulatory matters to the best of our knowledge and liability, but without obligation or liability. Existing laws and regulations are to be observed in all cases by our customers. This also applies in respect to any right of third parties. Our information and advice do not relieve our customers of their own responsibility for checking the

© 2018 Merck KGaA, Darmstadt, Germany and/or its affiliates. All Rights Reserved.

35 °C for 48 to 72 hours, while the plates for determination of the total yeast and mold count should be incubated at 20 to 25 °C for 5 to 7 days. Individual incubation conditions can be chosen and should be validated at the application side.

Several recommendations are given by different guidelines for incubation: according to USP <1116> the plates used for environmental monitoring should be incubated between 20 and 35 °C for not less than 72 hours. According to the FDA Aseptic Guide the plates for determination of the total aerobic bacterial count should be incubated at 30 to

Finally, the number of CFU per plate is examined.

Grown colonies are recommended to be identified.

### Storage and Shelf Life

The product can be used for sampling until the expiry date if stored upright, protected from light and properly sealed at +15 °C to +25 °C.

Condensation can be prevented by avoiding quick temperature shifts and mechanical stress.

The testing procedures as described on the CoA can be started up to the expiry date printed on the label.

### Disposal

Please mind the respective regulations for the disposal of used culture medium (e.g. autoclave for 20 min at 121 °C, disinfect, incinerate etc.).

### Quality Control

Control Strains	ATCC #	Inoculum CFU	Incubation	Expected Result Recovery in %
<i>Staphylococcus aureus</i>	6538	10-100	20-24 h at 30-35 °C	50-200
<i>Staphylococcus aureus</i> in the presence of 50µl Cutasept F	6538	10-100	20-24 h at 30-35°C	50-200
<i>Pseudomonas aeruginosa</i>	9027	10-100	20-24 h at 30-35 °C	50-200
<i>Bacillus subtilis</i>	6633	10-100	20-24 h at 30-35 °C	50-200

Merck, Millipore, the Vibrant M, Sigma-Aldrich are trademarks of Merck KGaA, Darmstadt, Germany or its affiliates. . All other trademarks are the property of their respective owners. [merckgroup.com/life-science](http://merckgroup.com/life-science)  
We provide information and advice to our customers on application technologies and regulatory matters to the best of our knowledge and liability, but without obligation or liability. Existing laws and regulations are to be observed in all cases by our customers. This also applies in respect to any right of third parties. Our information and advice do not relieve our customers of their own responsibility for checking the

© 2018 Merck KGaA, Darmstadt, Germany and/or its affiliates. All Rights Reserved.

<i>Candida albicans</i>	10231	10-100	44-48 h at 30-35 °C	50-200
<i>Aspergillus brasiliensis</i>	16404	10-100	44-48 h at 30-35°C	50-200

Please refer to the actual batch related Certificate of Analysis.

## Literature

EU GMP Medicinal Products for Human and Veterinary use (2008): Annex1 Manufacture of Sterile Medicinal Products.

European Pharmacopoeia 9.0 (2016): 2.6.12. Microbial examination of non-sterile products (total viable aerobic count).

Guidance for Industry (2004): Sterile Drug Products Produced by Aseptic Processing - Current Good Manufacturing Practice.

ISO 14698-1:2003: Cleanrooms and associated controlled environments - Biocontamination control - Part 1: General principles and methods.

Japanese Pharmacopoeia 16th edition (2011): 4.05 Microbial Limit Test.

PDA Technical Report No. 13 (2014 Revised): Fundamentals of an Environmental Monitoring Program.

United States Pharmacopoeia 41 NF 36 (2018): <61> Microbiological Examination of Non-Sterile Products: Microbial Enumeration Tests; <1116> Microbiological Control and Monitoring of Aseptic Processing Environments.

## Ordering Information

Product	Cat. No.	Pack size
Tryptic Soy Agar + LTHThio Contact - ICR+	1.46783.0020	20 x 55 mm plates
Tryptic Soy Agar + LTHThio Contact - ICR+	1.46783.0200	200 x 55 mm plates
Tryptic Soy Agar + LTHThio Sedi.- ICR	1.46786.0020	20 x 90 mm plates
Tryptic Soy Agar + LTHThio Sedi. - ICR	1.46786.0120	120 x 90 mm plates
Transport Bags, Sterile	1.46509.0125	25 x 5 bags

Merck, Millipore, the Vibrant M, Sigma-Aldrich are trademarks of Merck KGaA, Darmstadt, Germany or its affiliates. . All other trademarks are the property of their respective owners. [merckgroup.com/life-science](http://merckgroup.com/life-science)  
We provide information and advice to our customers on application technologies and regulatory matters to the best of our knowledge and liability, but without obligation or liability. Existing laws and regulations are to be observed in all cases by our customers. This also applies in respect to any right of third parties. Our information and advice do not relieve our customers of their own responsibility for checking the

© 2018 Merck KGaA, Darmstadt, Germany and/or its affiliates. All Rights Reserved.