

BOVIGAM™ Avian Tuberculin PPD Stimulating Ag

For the laboratory diagnosis of tuberculosis in animals by the Bovine Gamma Interferon Test

Catalog Number 7600065

Pub. No. MAN0017038 Rev. A.0

NOT FOR SKIN TESTING



WARNING! Read the Safety Data Sheets (SDSs) and follow the handling instructions. Wear appropriate protective eyewear, clothing, and gloves. Safety Data Sheets (SDSs) are available from thermofisher.com/support.



WARNING! POTENTIAL BIOHAZARD. Read the biological hazard safety information at this product's page at thermofisher.com. Wear appropriate protective eyewear, clothing, and gloves.

Introduction

Applied Biosystems™ BOVIGAM™ Avian Tuberculin PPD Stimulating Ag is a sterile, preservative free, liquid containing purified protein derivatives prepared from the filtrate of a heat-killed *Mycobacterium avium* (strain D4ER) grown on synthetic medium. It has been standardized by Guinea pig potency testing as described in European Pharmacopoeia and is issued at strength of 25,000 IU/mL.

General instructions

It is essential that all blood samples collected for BOVIGAM™ TB Kit and BOVIGAM™ 2G TB Kit (bovine gamma interferon test) testing are taken from a cleaned site into a tube containing heparin as anticoagulant. Gently invert tubes several times to ensure complete mixing of blood and heparin. Blood samples should be transported to the testing laboratory at ambient temperature (22±3°C, avoid extremes) and used in the BOVIGAM™ TB Kit and BOVIGAM™ 2G TB Kit assay within 30 hours of collection.

IMPORTANT! Blood must not be frozen or stored in a refrigerator.

BOVIGAM™ (bovine gamma interferon) tests

BOVIGAM™ TB Kit and BOVIGAM™ 2G TB Kit are rapid, blood-based assays of cell mediated immunity (CMI) for the diagnosis of bovine tuberculosis in cattle and other ruminants. Tuberculin PPD is mixed with lymphocytes in whole blood and incubated overnight at 37°C. Blood plasma is then collected from above the cultures and assayed for interferon-γ (IFN-γ) production using the BOVIGAM™ TB Kit BOVIGAM™ 2G TB Kit enzyme immunoassay (EIA).

General precautions

Laboratory safety

National Safety Regulations must be strictly followed.

Usage

For stimulation in 96-well cell culture plates

Dilute 110 µL of BOVIGAM™ Avian Tuberculin PPD Stimulating Ag in 890 µL of RPMI medium for final assay concentration of 250 IU/mL. Mix blood samples evenly immediately before use. For comparative testing, dispense **three** 250 µL aliquots of blood into culture vessels. To one aliquot, add 25 µL of diluted stock and mix thoroughly. To the other aliquots add 25 µL of sterile PBS to one and 25 µL of BOVIGAM™ Bovine Tuberculin PPD Stimulating Ag (Cat. No. 7600060) to the other. Aseptic procedures should be followed using a sterile pipetting apparatus.

For stimulation in 24-well cell culture plates

Dilute 160 µL of BOVIGAM™ Avian Tuberculin PPD Stimulating Ag in 840 µL of RPMI medium for final assay concentration of 250 IU/mL. Mix blood samples evenly immediately before use. For comparative testing, dispense **three** 1.5 mL aliquots of blood into culture vessels. To one aliquot, add 100 µL of diluted stock and mix thoroughly. To the other aliquots add 100 µL of nil control (sterile PBS) to one and 100 µL of BOVIGAM™ Bovine Tuberculin PPD Stimulating Ag (Cat. No. 7600060) to the other. Aseptic procedures should be followed using a sterile pipetting apparatus.

Interpretation

(See also the respective Package Inserts of BOVIGAM™ TB Kit and BOVIGAM™ 2G TB Kit for further details.)

Compare the absorbance values of the nil antigen, avian and bovine PPD samples for each animal.

Positive = OD Bovine PPD – Nil Antigen ≥ 0.1; and
OD Bovine PPD – Avian PPD ≥ 0.1

Negative = OD Bovine PPD – Nil Antigen < 0.1; or
OD Bovine PPD – Avian PPD < 0.1

Retesting

Retesting may be performed at any time after the last test.

Storage

Store, protected from light, at 2–8°C. Do not freeze.

Presentation

BOVIGAM™ Avian Tuberculin PPD Stimulating Ag is available in 5 mL vials.

Customer and technical support

Technical support: visit thermofisher.com/askaquestion

Visit thermofisher.com/support for the latest in services and support, including:

- Worldwide contact telephone numbers
- Order and web support
- User guides, manuals, and protocols
- Certificates of Analysis
- Safety Data Sheets (SDSs; also known as MSDSs)

NOTE: For SDSs for reagents and chemicals from other manufacturers, contact the manufacturer.

Limited product warranty

Life Technologies Corporation and/or its affiliate(s) warrant their products as set forth in the Life Technologies' General Terms and Conditions of Sale found on Life Technologies' website at www.thermofisher.com/us/en/home/global/terms-and-conditions.html. If you have any questions, please contact Life Technologies at thermofisher.com/support.



Prionics Lelystad B.V. | Platinastraat 33 | 8211 AR Lelystad | The Netherlands

The information in this guide is subject to change without notice.

DISCLAIMER: TO THE EXTENT ALLOWED BY LAW, LIFE TECHNOLOGIES AND/OR ITS AFFILIATE(S) WILL NOT BE LIABLE FOR SPECIAL, INCIDENTAL, INDIRECT, PUNITIVE, MULTIPLE, OR CONSEQUENTIAL DAMAGES IN CONNECTION WITH OR ARISING FROM THIS DOCUMENT, INCLUDING YOUR USE OF IT.

Revision history of Pub. No. MAN0017038 (English)

Rev.	Date	Description
A.0	7 June 2017	New document. Converted the legacy document [Bovigam_Lelystad_Tuberculin_PPD_Avian_v1.4_e.pdf] to the current document template, with associated updates to the publication number, limited license information, warranty, trademarks, and logos.

Important Licensing Information: These products may be covered by one or more Limited Use Label Licenses. By use of these products, you accept the terms and conditions of all applicable Limited Use Label Licenses.

©2017 Thermo Fisher Scientific Inc. All rights reserved. All trademarks are the property of Thermo Fisher Scientific and its subsidiaries unless otherwise specified.

For Veterinary Use Only. For In Vitro Use Only.

thermofisher.com/support | thermofisher.com/askaquestion

thermofisher.com

7 June 2017