

# A 20/20<sup>n</sup> Luminometer Method for Cambrex MycoAlert<sup>®</sup> Mycoplasma Detection Assay



## 1. INTRODUCTION

The Turner BioSystems' 20/20<sup>n</sup> Luminometer in combination with the MycoAlert<sup>®</sup> Assay from Cambrex Bio Science provides a convenient and rapid method for mycoplasma detection. Mycoplasma are common contaminants of cells grown in culture. With the MycoAlert Assay System, mycoplasma are detected within minutes, making it easy and reliable to use at every cell passage.

The MycoAlert assay follows a 2-step format. MycoAlert Reagent is added to the supernatant from cultured cells lysing any mycoplasma present and releasing mycoplasma enzymes. An initial luminescent measurement provides a baseline level of ATP for a particular sample. The addition of MycoAlert Substrate allows the mycoplasma enzymes to catalyze the conversion of ADP to ATP. The final luminescent measurement checks for an increase in ATP due to the presence of these enzymes.

The combination of the 20/20<sup>n</sup> Luminometer with the MycoAlert Assay is capable of detecting very low levels of mycoplasma. (Figure 1).

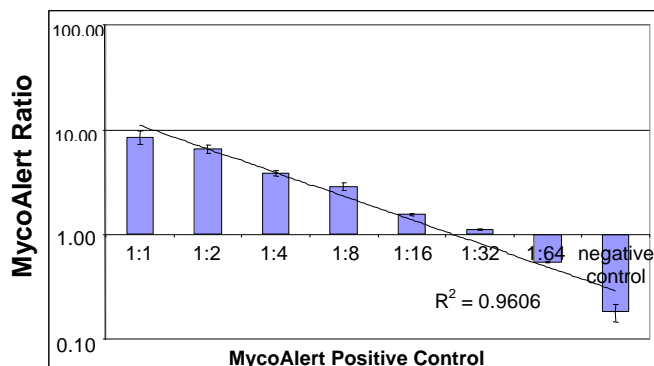


Figure 1. MycoAlert Assay performed on the 20/20<sup>n</sup> Luminometer using MycoAlert Positive Control.

## 2. MATERIALS REQUIRED

- 20/20<sup>n</sup> Luminometer (P/N 2030-000)
- 1.5 mL microfuge tube holder
- Centrifuge
- 1.5 mL microfuge tubes
- Luminometer tubes/cuvettes
- Pipettes (50–200  $\mu$ L and 200–1000  $\mu$ L) and the appropriate pipette tips
- MycoAlert Mycoplasma Detection Assay (LT07-118); which contains:
  - LT27-217 MycoAlert Reagent. Lyophilized. 2 x 600  $\mu$ L vials.
  - LT27-218 MycoAlert Assay Buffer. 1 x 10 ml bottle.
  - LT27-221 MycoAlert Substrate. Lyophilized. 2 x 600  $\mu$ L vials.

## 3. EXPERIMENT PROTOCOL

**NOTE:** Wear gloves to prevent ATP contamination from your hands during reagent preparation and sample analysis.

### 3.1 Reagent Preparation

3.1.1 Add 600  $\mu$ L of MycoAlert Assay Buffer into one vial containing the lyophilized MycoAlert Reagent. Replace white screw cap and mix gently. Allow the reagent to equilibrate for 15 minutes at room temperature.

3.1.2 Add 600  $\mu$ L of MycoAlert Assay Buffer into one vial containing the lyophilized MycoAlert Substrate. Replace the green screw cap and mix gently. Allow the reagent to equilibrate for 15 minutes at room temperature.

## 3.2 Instrument Setup

3.2.1 Turn ON the 20/20<sup>n</sup>. A 10-minute warm up period is recommended but not necessary. warm up period is recommended but not necessary.

3.2.2 Choose “Protocols” at the bottom of the touchscreen. Select “Default Protocol.” The Default Protocol measures samples for 1 second.

3.2.3 Touch “OK” to go to the Home Screen.

## 3.3 Sample Preparation and Analysis

3.3.1 Transfer 2 mL of cell culture or culture supernatant into a centrifuge tube and pellet any cells at 1500 rpm (200 x g) for 5 minutes.

3.3.2 Transfer 100 µL of cleared supernatant into a 1.5 mL clear microfuge tube.

3.3.3 Add 100 µL of MycoAlert<sup>®</sup> Reagent (white screw cap) to the 1.5 mL clear microfuge tube and incubate at room temperature (for 5 minutes).

3.3.4 Insert the microfuge tube into the tube holder of the 20/20<sup>n</sup>. Close the lid and touch “Measure Luminescence.” Record the result as Reading A.

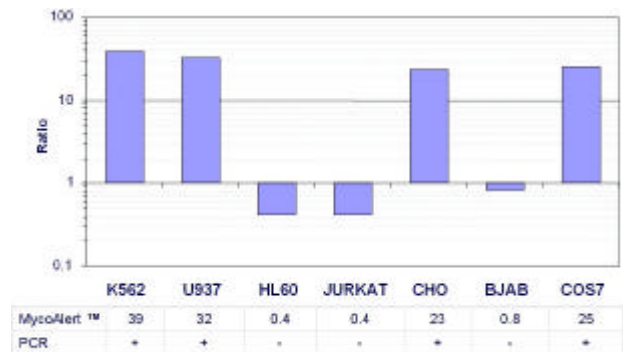
3.3.5 Add 100 µL of MycoAlert Substrate to the same microfuge tube and incubate at room temperature (for 10 minutes).

3.3.6 Repeat step 3.3.4 and record this result as Reading B.

3.3.7 Calculate ratio =  $\frac{\text{Reading B}}{\text{Reading A}}$

## 4. RESULTS

4.1 A ratio < 1 indicates an uninfected cell line. Mycoplasma infected cells routinely produce ratios greater than 1.



MycoAlert™ compared to Stratagene's Mycoplasma Plus™ PCR kit. Positive cell lines infected with M. hyorhina  
**Table 1. Example of cells analyzed for mycoplasma infection using the MycoAlert system.**

4.2 Borderline ratios (e.g. 1.0-1.3) should be re-tested after 24 hours in quarantine. Any cultures maintained in quarantine may be re-tested in 24 to 48 hours to see if the ratios increase.

## 5. ABOUT CAMBREX

MycoAlert is a registered trademark of CBM Intellectual Properties, Inc. MycoAlert Assay covered by U.S. patent and U.S. patent pending. Orders for Cambrex products may be placed by:

Phone: (800) 638-8174

Fax: (301) 845-8338

E-mail: [biotechserv@cambrex.com](mailto:biotechserv@cambrex.com)

Internet: <http://www.cambrex.com/bioproductions>

Mailing Address:

Cambrex Bio Science Rockland, Inc.  
 191 Thomaston Street  
 Rockland, ME 04841  
 USA

## **6. ABOUT TURNER BIOSYSTEMS**

20/20<sup>n</sup> is a trademark of Turner BioSystems.  
Orders for Turner BioSystems' products may be placed by:

Phone: (408) 636-2400 or  
Toll Free: (888) 636-2401 (US and  
Canada)  
Fax: (408) 737-7919

Web Site: [www.turnerbiosystems.com](http://www.turnerbiosystems.com)

E-Mail: [sales@turnerbiosystems.com](mailto:sales@turnerbiosystems.com)

Mailing Address:

Turner BioSystems, Inc.  
645 N. Mary Avenue  
Sunnyvale, CA 94085 USA