



## Egg Yolk with Potassium Tellurite

355-4201  
355-4205

### DEFINITION

Egg yolk with potassium tellurite is included in the complete Baird Parker medium formula. It is to be added to both of our presentations (dehydrated and ready to use).

### PRESENTATION

#### Ready to use

5 ml x 1 vial

code 355-4201

25 ml x bottle

code 355-4205

### STORAGE

- Ready to use: + 2°C to 8°C
- Expiration date and batch number are shown on the package.

### QUALITY CONTROL OF MANUFACTURER

Every product manufactured and marketed by Bio-Rad is subject to a quality-assurance procedure at all stages, from the reception of raw materials to the marketing of the end-product. Each batch of finished product undergoes quality control and is marketed only if it satisfies the acceptability criteria.

Documentation relative to the production and control of each batch is kept on file.