

## Yeast Extract/Additive

(Purified powdered yeast autolysate for bacteriology)

## ode 356-4343

### DEFINITION

This hydrosoluble yeast extract, specially purified and neutralised, is suitable for the preparation of most culture media for bacteriology. It offers numerous advantages compared to paste yeast extracts.

### PRESENTATION

#### Dehydrated

500 g

code 356-4343

### AVERAGE CHARACTERISTICS

Appearance Brown powder  
Density 0.6 to 0.7  
Odour Weak, non-putrid  
Final pH = 7.2 ± 0.2

Humidity Under 7%  
Sulfated ash 15%  
Sodium chloride Under 10%  
Total nitrogen 10%  
Amino nitrogen 4.5%  
Phosphorus (P<sub>2</sub>O<sub>5</sub>) 4%  
Calcium 0.1 to 0.2%  
Magnesium 0.25%  
Copper Less than 10 p.p.m.  
Iron Less than 100 p.p.m.  
Insoluble particles Negative (none)  
Precipitation after autoclaving Negative  
Culture tests Satisfactory

### STORAGE

- +15°C to 25°C, in carefully-sealed cask in a cool, dry place.
- Expiration date and batch number are shown on the package.

### UTILISATION

1 g of standard paste yeast extract can be replaced by 0.8 g of this powdered extract. Media prepared with this are pale, clear and do not precipitate on autoclaving. Used at the dose level of 0.3% to 0.5%, it is conducive to an excellent growth of all bacteria.

### PERFORMANCES / QUALITY CONTROL OF THE TEST

The growth performances of the media are verified with the following strains:

STRAINS	Results after 24h culture at 37°C
<i>Salmonella Enteridis</i> ATCC 13076	Good growth
<i>Staphylococcus aureus</i> ATCC 25923	Good growth
<i>Staphylococcus epidermidis</i> ATCC 12228	Good growth
<i>Streptococcus faecalis</i> var. <i>zymogenes</i> ATCC 29212	Good growth
<i>Streptococcus pyogenes</i> ATCC 19615	Good growth
<i>Listeria Monocytogenes</i>	Good growth

### 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.

### KEY WORDS

Yeast extract / Food products / Water / Enumeration / Isolation / Medium.