

Fast-Media® Blas TB

Microwaveable media for selection and propagation of blasticidin resistant *E. coli*

Catalog # fas-bl-l

For research use only

Version # 13F26-MM

PRODUCT INFORMATION

Contents:

- 20 individual sealed pouches of Fast-Media® Blas TB. Each pouch contains the necessary amount of powder for the preparation of **200 ml of Terrific Broth (TB) based liquid medium** supplemented with **blasticidin S**.

Effective concentration: Blasticidin 100 µg/ml

Storage and stability:

- Fast-Media® Blas TB are shipped at room temperature, and must be stored in a dry and cool place. They are stable for 2 years at room temperature.

- When properly prepared, Fast-Media® Blas TB broths are stable for 4 weeks at 4°C, retaining sterility and selective properties.

Quality control:

The high quality and performance of each formulation are tested with *E. coli* K12 derived strains. *E. coli* transformed with a plasmid carrying the blasticidin resistance gene are used as positive controls for Fast-Media® Blas TB.

METHOD

For customer convenience, the following procedure is directly printed on each pouch.

1. Pour the pouch contents into a clean borosilicate glass bottle or flask.
2. Add 200 ml of distilled or deionized water.
3. Mix thoroughly by swirling the glass bottle or flask.
4. Heat in a microwave oven on MEDIUM power setting (about 450W) until bubbles start to appear (about 3 minutes).

Do not heat in a closed container.

5. Swirl gently to mix the preparation and re-heat for 30 seconds. Swirl gently again.
6. Repeat step 4 if necessary until the medium is completely dissolved. Do not overboil.
7. Allow the medium to cool to 50-55 °C before use.

Caution: Any solution heated in a microwave oven may become superheated and suddenly boil when moved or touched. Handle with extreme care. Wear heat-proof gloves.

Note: Do not repeat this above procedure once the medium is prepared because the antibiotic will be adversely affected.

SPECIAL HANDLING

Caution should be exercised during handling of Fast-Media® due to potential allergenic properties of antibiotics. Wear protective gloves, do not breathe the dust.

FAST-MEDIA® FEATURES

Fast-Media® offer researchers a quick and convenient way to prepare 200 ml of sterile *E. coli* growth medium in about five minutes using a **microwave** instead of an autoclave.

Fast-Media® is supplied with a choice of antibiotics for selection (see table below), and chromogenic substrates, for the growth or selection of *E. coli* transformant colonies, as well as detection of blue/white colonies. Fast-Media® Base is supplied without selective antibiotics.

Several Fast-Media® are available:

- **Fast-Media® TB**, Terrific Broth based liquid medium
- **Fast-Media® LB**, Lysogeny Broth (LB) based liquid medium
- **Fast-Media® Agar**, LB based solid medium
- **Fast-Media® Agar X-Gal**, LB based solid medium containing IPTG and X-Gal
- **Fast-Media® Agar X-Gluc** LB based solid medium containing X-Gluc.

Fast-Media®	Agar	Agar X-Gal	Agar X-Gluc	LB	TB
Base	X			X	
Ampicillin	X	X		X	X
Blasticidin	X	X			X
Hygromycin	X	X			X
Kanamycin	X	X		X	X
Puromycin	X				X
Zeocin™	X	X	X		X

RELATED PRODUCTS

Product	Catalog Code
Fast-Media® Blas Agar	fas-bl-s
Fast-Media® Blas Agar X-Gal	fas-bl-x

TECHNICAL SUPPORT

Toll free (US): 888-457-5873

Outside US: (+1) 858-457-5873

Europe: +33 562-71-69-39

E-mail: info@invivogen.com

Website: www.invivogen.com



3950 Sorrento Valley Blvd. Suite 100
San Diego, CA 92121 - USA