

PlasmoTest™ Reagent Kit

PlasmoTest™ Reagents for 500 samples

Catalog code: rep-ptrk

<https://www.invivogen.com/plasmotest-reagent>

For research use only

Version 22C28-MM

PRODUCT INFORMATION

Contents

- 2 pouches HEK-Blue™ Detection
- 8 x 1 ml HEK-Blue™ Selection (250X concentrated solution)
- 2 x 60 ml HEK-Blue™ Water (sterile endotoxin-free water)
- 4 x 1 ml Normocin™ (50 mg/ml)
- 1 vial Negative Control
- 1 vial Positive Control

Storage and stability

- PlasmoTest™ Reagent Kit is shipped at room temperature.
 - Upon receipt, store HEK-Blue™ Selection at 4 °C or at -20 °C.*
 - Upon receipt, store Normocin™ at 4 °C or at -20 °C.*
- * The expiry date is specified on the product label.
- Upon receipt, store sealed HEK-Blue™ Detection pouches at 4 °C. Reconstituted HEK-Blue™ Detection is stable for 2 weeks at 4 °C and for 2 months at -20 °C. Protect from light.
 - Upon receipt, store lyophilized Positive and Negative Controls at 4 °C. Upon resuspension, prepare aliquots of Positive Control and store at 4 °C or at -20 °C. Store resuspended Negative Control at 4 °C.
- Note: Avoid repeated freeze-thaw cycles. Data sheets for all components are available on our website.

DESCRIPTION

HEK-Blue™ Selection

HEK-Blue™ Selection is a solution that combines several selective antibiotics. These antibiotics guarantee the persistent expression of the various transgenes introduced in HEK-Blue™-2 Cells. For cell culture maintenance, supplement the growth medium with 1X HEK-Blue™ Selection, which represents a 1:250 dilution of stock solution. Dilute HEK-Blue™ Selection directly in the culture medium.

Normocin™

Normocin™ combines three antibiotics active against mycoplasma, bacteria and fungi. For cell culture maintenance, Normocin™ is used at a concentration of 100 µg/ml, which represents a 1:500 dilution of stock solution. Dilute Normocin™ directly in the culture medium.

HEK-Blue™ Detection

HEK-Blue™ Detection is a medium specifically designed for the detection of secreted alkaline phosphatase (SEAP). In the presence of mycoplasma-contaminated samples, HEK-Blue™-2 cells secrete SEAP in the HEK-Blue™ Detection medium resulting in a color change from pink to purple/blue. HEK-Blue™ Detection is a powdered medium provided in individually sealed pouches. Each pouch allows the preparation of 50 ml of detection medium with HEK-Blue™ Water.

Positive and Negative Controls

The Positive and Negative Controls are provided lyophilized. After resuspension with 1 ml of HEK-Blue™ Water, 10X solutions of Positive and Negative Controls are obtained. They have to be diluted at 1/10 before use. Each vial allows to perform 200 tests.

METHODS

All reagents should be prepared under sterile conditions according to good laboratory practices.

HEK-Blue™ Detection Medium preparation

1. Pour the contents of one pouch of HEK-Blue™ Detection into a sterile vial/bottle.
 2. Solubilize the powder with 50 ml of endotoxin-free water.
 3. Swirl gently until powder is completely dissolved.
 4. Warm reconstituted HEK-Blue™ Detection to 37 °C for 30 minutes to 1 hour.
 5. Filter the medium through a 0.2 µm membrane into a sterile vial/bottle.
- Note: We recommend using filter units providing a large filter area to facilitate filtration.
6. Keep the HEK-Blue™ Detection medium at 37 °C before use or store at 4 °C for up to 2 weeks or at -20 °C for up to 2 months.

Positive and Negative Controls preparation

1. To obtain 10X solutions of Positive and Negative Controls, add 1 ml of HEK-Blue™ Water to each tube.
2. Mix vigorously by vortexing as the 10X Positive and Negative Controls may stick to the tube wall.
3. Prepare 50 µl aliquots of 10X Positive Controls for storage at -20 °C.
4. Prepare a 1/10 dilution of 10X Positive and Negative Controls with HEK-Blue™ Water.

10X and 1X solutions of Positive Control are stable for 1 month at 4 °C and for 6 months at -20 °C, while 10X and 1X solutions of Negative Control are stable for 6 months at 4 °C when stored properly.

RELATED PRODUCTS

Product	Description	Cat.Code
HEK-Blue™ Detection	SEAP detection medium	hb-det2
HEK-Blue™ Selection	Antibiotic mix	hb-sel
Normocin™	Antimicrobial reagent	ant-nr-1
PlasmoTest™	Mycoplasma detection kit	rep-pt1
PlasmoTest™ Controls	Controls for PlasmoTest™	pt-ctr2

TECHNICAL SUPPORT

InvivoGen USA (Toll-Free): 888-457-5873

InvivoGen USA (International): +1 (858) 457-5873

InvivoGen Europe: +33 (0) 5-62-71-69-39

InvivoGen Asia: +852 3622-3480

E-mail: info@invivogen.com

