

HEK-Blue™ Selection

Antibiotic mixture for the maintenance of certain HEK-Blue™ cells; Cell culture tested

Catalog code: hb-sel

<https://www.invivogen.com/hek-blue-selection>

For research use only

Version 24J10-MM

PRODUCT INFORMATION

Contents

10 x 1 ml tubes of HEK-Blue™ Selection 250X concentrate

Storage and stability

- HEK-Blue™ Selection is shipped at room temperature and should be stored at 4 °C or -20 °C. Avoid repeated freeze-thaw cycles.

- The expiry date is specified on the product label.

Note: HEK-Blue™ Selection is stable for 2 weeks at room temperature.

Quality control

Each lot is thoroughly tested to ensure the absence of lot-to-lot variation.

- Physicochemical characterization (pH, appearance)

- The activity of each antibiotic in HEK-Blue™ Selection has been individually confirmed in mammalian cells.

- HEK-Blue™ Selection has been functionally validated.

DESCRIPTION

HEK-Blue™ Selection is a solution that combines several selection antibiotics. These antibiotics guarantee the persistent expression of the various transgenes introduced in the HEK-Blue™ cells listed in table 2. The concentration of antibiotics has been optimized to maximize their effectiveness for the maintenance of these cells. HEK-Blue™ Selection is a simple and convenient solution as all the antibiotics required are provided in a single tube and it can be simply added to the culture at a 1:250 dilution.

IMPORTANT HEK-Blue™ Selection is only for use with a defined set of HEK-Blue™ cells as listed in table 2.

SAFETY CONSIDERATIONS

HEK-Blue™ Selection is a harmful compound. Refer to safety data sheet for handling instructions.

METHOD

Growth Medium for cell culture maintenance of the HEK-Blue™ cells listed in table 2:

DMEM 4.5 g/l glucose, 2-4 mM L-glutamine, supplemented with 10% FBS, Penicillin-Streptomycin, and 1X HEK-Blue™ Selection.

1. Warm at 37 °C before use.
2. Dilute 250-fold (v/v) directly in the culture medium.

Reagent	T25 with 5 ml medium	T75 with 15 ml medium	T175 with 30 ml medium
HEK-Blue™ Selection	20 µl	60 µl	120 µl

Table 1. Recommended volumes for HEK-Blue™ Selection

APPLICATION

Use HEK-Blue™ Selection for the maintenance of the cells listed below.

Cell lines	Cat. code
HEK-Blue™-2 cells (part of Plasmotest™ Kit)	rep-pt1
HEK-Blue™-4 cells (part of EndotoxDetect™ Kit)	rep-lps2
HEK-Blue™-hTLR2 cells	hkb-htlr2
HEK-Blue™ hTLR2-TLR1 cells	hkb-htlr21
HEK-Blue™ hTLR2-TLR6 cells	hkb-htlr26
HEK-Blue™ hTLR2 KO-TLR1/6 cells	hkb-htlr2k16
HEK-Blue™ mTLR2 cells	hkb-mtlr2
HEK-Blue™ hTLR4 cells	hkb-htlr4
HEK-Blue™ mTLR4 cells	hkb-mtlr4
HEK-Blue™ IL-1R cells	hkb-il1r
HEK-Blue™ IL-5 cells	hkb-il5
HEK-Blue™ IL-6 cells	hkb-hil6
HEK-Blue™ IL-7 cells	hkb-il7
HEK-Blue™ IL-9 cells	hkb-il9
HEK-Blue™ IL-10 cells	hkb-il10
HEK-Blue™ IL-11 cells	hkb-hil11r
HEK-Blue™ IL-12 cells	hkb-il12
HEK-Blue™ IL-17 cells	hkb-il17
HEK-Blue™ IL-17C cells	hkb-il17c
HEK-Blue™ IL-18 cells	hkb-hmil18
HEK-Blue™ IL-21 cells	hkb-il21
HEK-Blue™ IL-23 cells	hkb-il23
HEK-Blue™ IL-31 cells	hkb-il31
HEK-Blue™ IL-33 cells	hkb-hil33
HEK-Blue™ CD122/CD132 cells	hkb-il2bg
HEK-Blue™ G-CSF cells	hkb-gcsf
HEK-Blue™ GM-CSF cells	hkb-hgmcsfr
HEK-Blue™ STAT6-hSTING-R232 cells	hkb-st6r232
HEK-Blue™ TPO cells	hkb-tpo
HEK-Blue™ TSLP cells	hkb-tslp

Table 2. Cell lines requiring HEK-Blue™ Selection

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



Any questions about our selective antibiotics? Visit our FAQ page.

 **InvivoGen**
www.invivogen.com

RELATED PRODUCTS

Product	Description	Cat. Code
EndotoxDetect™	HEK-Blue™ LPS detection kit	rep-lps2
Fungin™	Antifungal agent	ant-fn-1
MycStrip®	Mycoplasma detection kit	rep-mys-10
Normocin®	Antimicrobial agent	ant-nr-05
Normocure™	Antibacterial agent	ant-noc
Plasmocin® Prophylactic	Anti-mycoplasma agent	ant-mpp
Plasmocin® Treatment	Mycoplasma removal agent	ant-mpt-1
PlasmoTest™	Mycoplasma detection kit	rep-pt1

InvivoGen provides an expanding collection of cell lines in which specific genes are over-expressed or knocked out. For more information, visit www.invivogen.com/cell-lines.

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



Any questions about our selective antibiotics? Visit our FAQ page.