

HEK-Blue™ CLR Selection

Antibiotics for maintenance of HEK-Blue™ Dectin, HEK-Blue™ Mincle and HEK-Blue™ IL-2 cells

Cell culture tested

Catalog code: hb-csm

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

For research use only

Version 21J21-MM

PRODUCT INFORMATION

Contents

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

Storage and stability:

- HEK-Blue™ CLR Selection is shipped at room temperature. Upon receipt it should be stored at 4 °C or -20 °C. Avoid repeated freeze-thaw cycles.
- The expiry date is specified on the product label.

Quality control

- Physicochemical characterization (pH, appearance)
- The activity of each antibiotic in HEK-Blue™ CLR Selection has been individually confirmed in mammalian cells.
- HEK-Blue™ CLR Selection has been functionally validated.

PRODUCT DESCRIPTION

HEK-Blue™ CLR Selection is a solution that combines several selective antibiotics. These antibiotics guarantee the persistent expression of the various transgenes introduced in HEK-Blue™ cells such as the C-type Lectin (CLR) genes, namely Dectin-1, Dectin-2, and Mincle. The use of HEK-Blue™ CLR Selection is also recommended to guarantee persistent transgene expression in HEK-Blue™ IL-2 cells.

The concentration of antibiotics has been optimized to maximize their effectiveness for the maintenance of these cells. HEK-Blue™ CLR 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™ CLR Selection is only for use with a defined set of HEK-Blue™ cells as listed in table 2.

SAFETY CONSIDERATIONS

HEK-Blue™ CLR 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™ CLR Selection.

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

Table 1. Recommended volumes for HEK-Blue™ CLR Selection

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

APPLICATION

HEK-Blue™ CLR Selection should be used for the maintenance of the cell lines listed in the table below.

Table 2. Cell lines requiring HEK-Blue™ CLR Selection

Cell lines	Cat. code
HEK-Blue™ hDectin-1a cells	hkb-hdect1a
HEK-Blue™ hDectin-1b cells	hkb-hdect1b
HEK-Blue™ mDectin-1b cells	hkb-mdect1b
HEK-Blue™ mDectin-2 cells	hkb-mdect2
HEK-Blue™ mMincle cells	hkb-mmcl
HEK-Blue™ IL-2 cells	hkb-hil2

RELATED PRODUCTS

Product	Description	Cat. Code
Fungin™	Antifungal agent	ant-fr-1
MycoStrip™	Mycoplasma detection kit	rep-mys-10
Normocin™	Antimicrobial agent	ant-nr-1
Normocure™	Antibacterial agent	ant-noc
Plasmocin™ Prophylactic	Anti-mycoplasma agent	ant-mpp
Plasmocin™ Treatment	Mycoplasma removal agent	ant-mpt
PlasmoTest™	Mycoplasma detection kit	rep-pt1

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 3622-34-80

E-mail: info@invivogen.com



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

 **InvivoGen**
www.invivogen.com