

# HEK-Blue™ Selection

Antibiotics for maintenance of HEK-Blue™ Cells

Catalog # hb-sel

For research use only

Version # 10B26-MM

## PRODUCT INFORMATION

### **Contents:**

HEK-Blue™ Selection is provided as 5 x 2 ml tubes of 250X solution.

### **Storage and stability:**

Store HEK-Blue™ Selection at 4°C for immediate use or -20°C for long term storage. Antibiotics are stable for 6 months at 4°C and 12 months when stored at -20°C. Avoid repeated freeze-thaw cycles.

## BACKGROUND

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™ cells.

### **HEK-Blue™ Growth Medium for cell culture maintenance:**

DMEM high glucose supplemented with 10% FBS, Penicillin-Streptomycin, and 1X HEK-Blue™ Selection Warm at 37°C before use and store at 4°C.

---

### 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](mailto:info@invivogen.com)

Website: [www.invivogen.com](http://www.invivogen.com)



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