

PlasmoTest™ Controls

Positive and Negative Controls allowing 200 tests

Catalog pt-ctr2

For research use only

Version # 15C04-MM

PRODUCT INFORMATION

Content:

- 1 vial of Positive Control
- 1 vial of Negative Control
- 10 ml endotoxin-free water

Storage and stability:

- PlasmoTest™ Controls (Positive and Negative Controls) are provided lyophilized and shipped at room temperature.
- Store lyophilized Positive and Negative controls at 4 °C for up to 12 months. Resuspended Negative control is stable at 4 °C for 6 months when properly stored. Upon resuspension, prepare aliquots of the Positive control and store at 4 °C for short term storage (1 month) or -20 °C for long term storage (6 months).

Note: Avoid repeated freeze-thaw cycles.

Preparation of stock solutions

Solubilize the Positive and Negative Controls by adding 1 ml of endotoxin-free water to each tube. 10X solutions of Positive and Negative Controls are obtained.

- Mix by vortexing as the 10X Positive and Negative Controls may stick to the tube wall.
- Prepare 50 µl aliquots of 10X Positive Controls for storage at -20 °C.
- Prepare a 1/10 dilution of 10X Positive and Negative Controls with endotoxin-free 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.

Stimulation of HEK-Blue™-2 cells

- Add 50 µl of 1X Negative Control in a well of a flat-bottom 96-well plate.

- Add 50 µl of 1X Positive Control in another well.

Note: Use new sterile filter-tips for each well to avoid cross-contamination.

- Add 200 µl (50,000 cells) of HEK-Blue™-2 cells suspension, prepared following the PlasmoTest™ protocol, to each well.

Note: Plasmotest™ protocol is available online at
<http://www.invivogen.com/plasmotest>

- Incubate the plate at 37 °C in a CO₂ incubator for 16-24 hours.

- Determine secreted alkaline phosphate levels with the naked eye or a spectrophotometer at 620-655 nm.

Note: Blue color indicates a positive result and pink color indicates a negative result.

RELATED PRODUCTS

Product	Catalog Code
HEK-Blue™ Detection	hb-det2
HEK-Blue™ Selection	hb-sel
Normocin™	ant-nr-1
PlasmoTest™	rep-pt2
PlasmoTest™ Reagents Kit	rep-ptrk

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 Hong Kong : +852 3-622-34-80
E-mail: info@invivogen.com

