

# HEK-Blue™ Endotoxin Standard

Standardized *E. coli* O55:B5 LPS

Catalog code: rep-hbes-10

<https://www.invivogen.com/endotoxin-standard>

For research use only

Version 22J11-MM

## PRODUCT INFORMATION

### Contents

- 10 tubes of HEK-Blue™ Endotoxin Standard. Each vial contains approximately 40-100 EU lyophilized endotoxin. The actual value of X is stated on the Certificate of Analysis.
- 1 bottle of endotoxin-free water (25 ml)

### Storage and stability

- HEK-Blue™ Endotoxin Standard is provided lyophilized and shipped at room temperature.
- Store lyophilized product at -20°C for up to 12 months.
- Resuspended product is stable at 2-8°C for 1 week when properly stored.

## DESCRIPTION

HEK-Blue™ Endotoxin Standard is a preparation of *E. coli* O55:B5 LPS standardized against FDA approved control standard endotoxin (CSE). HEK-Blue™ Endotoxin Standard is designed for use with the HEK-Blue™ LPS Detection Kit 2, a cell-based colorimetric assay for the detection of biologically active endotoxin that offers a sustainable alternative to the LAL test.

## METHOD

### Preparation of stock solution

- Reconstitute by adding 1 ml of endotoxin-free water (provided) to the content of the tube to obtain an initial stock solution at X EU/ml, where X is the amount of EU per vial.

*Note: Each vial contains approximately 40-100 EU lyophilized endotoxin. The actual value of X is stated on the Certificate of Analysis. For example, if the X value of the vial is 55 EU, when reconstituted with 1 ml endotoxin-free water it will yield a concentration of 55 EU/ml.*

- Mix vigorously by vortexing as endotoxin tends to attach to glass.
- Prepare a 1.0 EU/ml HEK-Blue™ Endotoxin Standard stock solution by diluting 1/X. For example, if the initial stock solution is at 55 EU/ml, the dilution is 1/55 or 0.1 ml of initial stock solution into 5.4 ml of endotoxin-free water.
- Store at 2-8°C. Initial and 1.0 EU/ml HEK-Blue™ Endotoxin Standard stock solutions are stable for 1 week.

## RELATED PRODUCTS

Product	Description	Catalog Code
HEK-Blue™ LPS Detection Kit 2	LPS detection assay	rep-lps2
HEK-Blue™ Selection	Mix of selective antibiotics	hb-sel
Normocin™	Antimicrobial agent	ant-nr-1
QUANTI-Blue™ Solution	SEAP detection medium	rep-qbs

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

