

# HKEB

## Heat Killed *E.coli* 0111:B4 - TLR2 & TLR4 agonist

Catalog # tlr1-hkeb2

For research use only

Version # 12F07-MM

### PRODUCT INFORMATION

#### Content:

- 10<sup>10</sup> freeze-dried cells of Heat Killed *E.coli* 0111:B4 (HKEB)
- 1.5 ml sterile endotoxin-free water

Endotoxin level: >1 EU/10<sup>9</sup> cells

#### Storage:

- HKEB is provided lyophilized and shipped at room temperature. Store at 4°C.
- Upon resuspension, prepare aliquots of HKEB and store at 4°C for short term storage or -20°C for long storage.
- Resuspended product is stable 1 month at 4°C and at least 1 year at -20°C when properly stored.

### DESCRIPTION

HKEB is a heat killed preparation of the gram negative bacterium, *E.coli* 0111:B4. Cell wall components from this bacterium, such as peptidoglycan (PGN) and lipopolysaccharide (LPS), are recognized by the toll-like receptor (TLR) 2 and TLR4<sup>1</sup>. It has been demonstrated that HKEB can stimulate TLR2 and induce the production of NF-κB and pro-inflammatory cytokines, such as IL-8<sup>2</sup>. HKEB is a potent stimulator of TLR2 and TLR4.

1. Takeuchi O. *et al.*, 1999. Differential roles of TLR2 and TLR4 in recognition of gram negative and gram-positive bacterial cell wall components. *Immunity*, 11(4):443-51.

2. van Riet E. *et al.*, 2009. Combined TLR2 and TLR4 ligation in the context of bacterial or helminth extracts in human monocyte derived dendritic cells: molecular correlates for TH1/TH2 polarization. *BMC Immunology*, 10:9.

### METHODS

#### Preparation of stock solution (10<sup>10</sup> HKEB cells/ml)

Stimulation of TLR2 and TLR4 can be achieved with 10<sup>5</sup> - 10<sup>7</sup> HKEB cells/ml.

- Add 1 ml sterile endotoxin-free water (provided) to rehydrate the pellet.
- Vortex for 10 seconds or until homogenized.

*Note: Rehydrated HKEB results in a milky suspension.*

#### Example TLR2 and TLR4 stimulation using HKEB

HKEB can be used to stimulate TLR2 or TLR4 in HEK-Blue™ TLR2 and HEK-Blue™ TLR4 cells, respectively. HEK-Blue™-TLR cells stably express an NF-κB-inducible secreted embryonic alkaline phosphatase (SEAP) and overexpress a TLR gene. For more information visit: [www.invivogen.com/hek-blue-tlr-cells](http://www.invivogen.com/hek-blue-tlr-cells)

- Add 10<sup>5</sup> - 10<sup>7</sup> HKEB cells/ml to HEK-Blue™ TLR cells (prepare cell suspension according to data sheet).

- Incubate cells and HKEB for 6 - 24 h at 37°C, 5% CO<sub>2</sub>.

- Determine TLR2 or TLR4 stimulation with HKEB by assessing cytokine expression using an ELISA, or SEAP expression using a SEAP detection medium, such as QUANTI-Blue™.

### RELATED PRODUCTS

Product	Catalog Code
HEK-Blue™ hTLR2 cells	hkb-htlr2
HEK-Blue™ hTLR4 cells	hkb-htlr4
QUANTI-Blue™	rep-qb1
<b>Other TLR2 ligands:</b>	
FSL-1 (synthetic diacylated lipoprotein)	tlr1-fsl
HKLM (heat killed <i>L.monocytogenes</i> )	tlr1-hklm
LM-MS (Lipomannan from <i>M.smegmatis</i> )	tlr1-lmm2
Pam3CSK4 (synthetic triacylated lipoprotein)	tlr1-pms

#### 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 3622-3480  
E-mail: [info@invivogen.com](mailto:info@invivogen.com)