

HKST

Heat Killed *Salmonella typhimurium*; TLR2 and TLR4 agonist

Catalog # tlr1-hkst2

<http://www.invivogen.com/hkst>

For research use only

Version # 17C15-MM

PRODUCT INFORMATION

Content:

- 10¹⁰ freeze-dried cells of Heat Killed *Salmonella typhimurium* strain CDC 6516-60 (HKST)
- 1.5 ml endotoxin-free water

Storage:

- HKST is provided lyophilized and shipped at room temperature. Store at 4°C.
- Upon resuspension, prepare aliquots of HKST and store at -20°C.
- Resuspended product is stable for 1 year at -20°C when properly stored.

DESCRIPTION

HKST is a heat killed preparation of the Gram negative bacterium, *Salmonella typhimurium*. Recognition of HKST is mediated by TLR2 and TLR4^{1,2}. TLR2 and TLR4 recognize cell wall components from HKST, such as peptidoglycan (PGN) and lipopolysaccharide (LPS) resulting in the production of pro-inflammatory cytokines, such as IL-6 and TNF- α ¹. InvivoGen's HKST is a potent agonist of TLR2 and TLR4.

1. Lembo A. *et al.*, 2003. Differential Contribution of Toll-Like Receptors 4 and 2 to the Cytokine Response to *Salmonella enterica* Serovar Typhimurium and *Staphylococcus aureus* in Mice. *Infect Immun.* 71(10):6058-62. 2. Arpaia N. *et al.*, 2011. TLR signaling is required for virulence of an intracellular pathogen Cell, 144(5):675-688.

METHODS

Preparation of stock suspension

To prepare a stock suspension at 10¹⁰ cells/ml:

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

Note: Rehydrated HKST results in a cloudy suspension.

TLR2 and TLR4 activation using HKST

Activation of TLR2 and TLR4 by HKST can be determined using HEK-Blue™ TLR2 and HEK-Blue™ TLR4, respectively. HEK-Blue™ TLR cells stably express an NF- κ B-inducible secreted embryonic alkaline phosphatase (SEAP) and overexpress a TLR gene.

For more information visit: www.invivogen.com/hek-blue-tlr-cells

In the HEK-Blue™-TLR cells, the optimal concentration of HKST is 10⁴ cells/ml to activate TLR2 and 10⁶ cells/ml to activate TLR4.

- Add 10⁴-10⁶ HKST cells/ml to HEK-Blue™ TLR cells (prepare cell suspension according to data sheet).
- Incubate cells and HKST for 6-24 h at 37°C, 5% CO₂.
- Determine TLR2 or TLR4 stimulation with HKST by assessing cytokine expression using an ELISA, or SEAP expression using a SEAP detection medium, such as HEK-Blue™ Detection.

RELATED PRODUCTS

Product	Catalog Code
HEK-Blue™ hTLR2 cells	hkb-htlr2
HEK-Blue™ hTLR4 cells	hkb-htlr4
HEK-Blue™ Detection	hb-det2
Other TLR2 ligands:	
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
Other TLR4 ligands:	
LPS-EB Ultrapure (LPS from <i>E. coli</i> O111:B4)	tlr1-3pelps
LPS-EK Ultrapure (LPS from <i>E. coli</i> K12)	tlr1-peklps
MPLA (monophosphoryl lipid A from <i>S. minnesota</i>)	tlr1-mpla

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