

# FLA-BS

## Flagellin from *B. subtilis*; TLR5 ligand

Catalog # tlr1-bsfla

<http://www.invivogen.com/fla-bs>

**For research use only**

Version # 17L04-MM

### PRODUCT INFORMATION

#### **Content:**

- 100 µg FLA-BS (flagellin from *Bacillus subtilis*)
- 1.5 ml endotoxin-free water

#### **Storage:**

- FLA-BS is shipped at room temperature and should be stored at -20 °C.
- Upon resuspension, prepare aliquots of FLA-BS and store at -20 °C. Resuspended product is stable for 6 months at -20 °C when properly stored. Avoid repeated freeze-thaw cycles.

#### **Quality Control:**

- Endotoxin levels: <0.05 EU/µg
- Biological activity has been confirmed using HEK-Blue™ hTLR5 cells.

### DESCRIPTION

FLA-BS, a ~32 kDa protein, is isolated from the Gram-positive bacteria *Bacillus subtilis*. Flagellin, the principal component of the flagella present on many Gram-negative and Gram-positive bacteria, is a proinflammatory molecule recognized by distinct types of pattern recognition receptors (PRRs); the surface localized Toll-like receptor (TLR5)<sup>1</sup> and the cytosolic NOD-like receptors (NLRs), NLRC4 and NAIP5<sup>2</sup>. Extracellular flagellin is detected by TLR5 resulting in MyD88-mediated NF-κB activation, cytokine and nitric oxide production depending on the nature of the TLR5 signaling complex<sup>3</sup>. Intracellular flagellin is detected by NLRC4 (also known as IPAF) and NAIP5. Recognition by NLRC4 and NAIP5, leads to inflammasome assembly, triggering caspase-1 activation of IL-1β and IL-18.

**1. Hayashi F. et al., 2001.** The innate immune response to bacterial flagellin is mediated by Toll-like receptor 5. *Nature* 410(6832):1099-103. **2. Zhao et al., 2011.** The NLRC4 inflammasome receptors for bacterial flagellin and type III secretion apparatus. *Nature*. 2011 Sep 14;477(7366):596-600. **3. Mizel SB. et al., 2003.** Induction of macrophage nitric oxide production by Gram-negative flagellin involves signaling via heteromeric Toll-like receptor 5/Toll-like receptor 4 complexes. *J Immunol.* 170(12):6217-23.

### METHODS

#### **Preparation of stock solution (500 µg/ml):**

- Stimulation of TLR5 can be achieved with FLA-BS at a concentration of 10 ng -10 µg/ml.
- Open vial lid carefully to avoid any loss of product.
  - Add 200 µl of the endotoxin-free water provided and mix by pipetting. Do not vortex.

#### **TLR5 stimulation using FLA-BS**

FLA-BS can be used to stimulate TLR5 in HEK-Blue™ TLR5 cells. HEK-Blue™ TLR5 cells stably overexpress the TLR5 gene and an NF-κB-inducible secreted embryonic alkaline phosphatase (SEAP). For more information, visit: [www.invivogen.com/hek-blue-tnfr5](http://www.invivogen.com/hek-blue-tnfr5)

- Dispense 20 µl of FLA-BS (10 ng -10 µg/ml final concentration) per well of a 96-well plate.
- Prepare a cell suspension of HEK-Blue™ TLR5 cells according to the data sheet.
- Add 180 µl of HEK-Blue™ TLR5 cell suspension per well.
- Incubate the plate for 6-24 h at 37 °C, 5% CO<sub>2</sub>.
- Collect 20 µl of supernatant and add to a well of a 96-well plate containing 180 µl of QUANTI-Blue™.
- Incubate the plate at 37 °C for 1-3 h.
- Determine SEAP levels using a spectrophotometer at 620-655 nm.

### RELATED PRODUCTS

| Product   | Catalog Code |
|---|--------------|
| HEK-Blue™ hTLR5 cells (human TLR5)                            | hkb-htlr5    |
| HEK-Blue™ mTLR5 cells (mouse TLR5)                            | hkb-mtlr5    |
| QUANTI-Blue™  | rep-qb1      |
| <b>Other TLR5 ligands:</b>                                    |              |
| FLA-BS Ultrapure (flagellin from <i>B. subtilis</i> )         | tlr1-pbsfla  |
| FLA-ST Ultrapure (flagellin from <i>S. typhimurium</i> )      | tlr1-epstfla |
| RecFLA-ST (recombinant flagellin from <i>S. typhimurium</i> ) | tlr1-flic    |

#### 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)