

Validation data for HKBF

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

For research use only

Version # 17L15-MM

HKBF is a pasteurized preparation of nontoxigenic *Bacteroides fragilis* (NTBF), a bile-resistant, Gram-negative, obligate anaerobe present in normal intestinal microbiota. Recent research has focused on the role of this bacteria in immunity and inflammation. Indeed, the presence of *B. fragilis* exerts a protective effect in animal models of colitis. Specifically, it is believed that the capsular polysaccharide A (PSA) of *B. fragilis* directly acts on regulatory T cells (Tregs) and also induces anti-inflammatory cytokine production primarily in a TLR2-dependent manner.

Stimulation of InvivoGen's HEK-Blue™ hTLR2 reporter cells with HKBF results in a dose-dependent induction of the NF-κB signaling pathway (Figure 1). HKBF does not activate InvivoGen's HEK-Blue™ hTLR4 reporter cells confirming the absence of endotoxin contamination.

Evaluation of HKBF in HEK-Blue™ hTLR2 and HEK-Blue™ hTLR4 cells

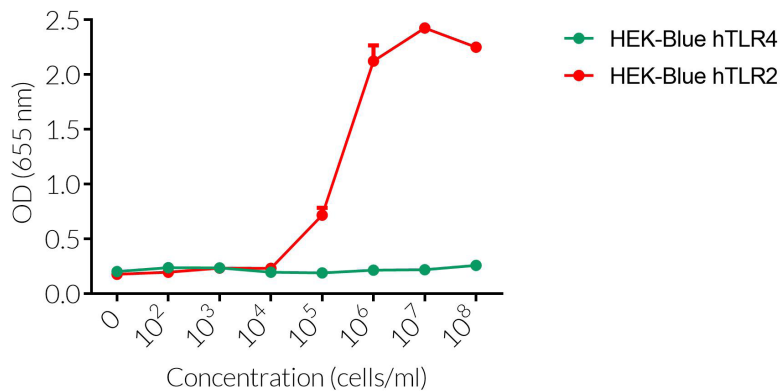


Figure 1: HKBF induces a dose-dependent response in HEK-Blue™ hTLR2 cells.

HEK-Blue™ hTLR2 and HEK-Blue™ hTLR4 cells were stimulated with increasing concentrations of HKBF. After overnight incubation, the NF-κB response was determined using QUANTI-Blue™, a SEAP detection reagent, and by reading the optical density (OD) at 655 nm.

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