

Validation data for HEK-Blue™ mMincle cells

For research use only

Version # 15B09-MM

HEK-Blue™ mMincle cells stably express the murine Mincle (mMincle) gene, as well as the genes of the Mincle-NF-κB signaling pathway. They also stably express an NF-κB-inducible secreted embryonic alkaline phosphatase (SEAP) reporter gene. These reporter cells are activated by Mincle ligands, such as trehalose-6,6-dibehenate (TDB) and heat-killed *M. tuberculosis* (HKMT). They do not respond to other CLR ligands such as zymosan, a Dectin-1 and TLR2 ligand.

Response of HEK-Blue™ mMincle cells to CLR agonists

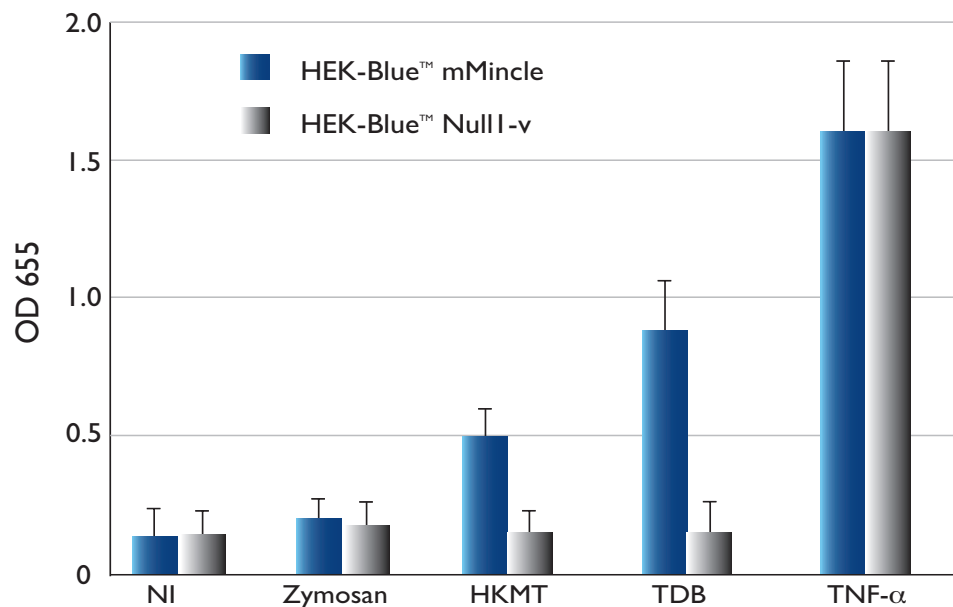


Figure 1: HEK-Blue™ mMincle and HEK-Blue™ Null1-v cells (parental cell line) were stimulated with various CLR agonists: zymosan (100 µg/ml), HKMT (100 µg/ml), TDB (10 µg/ml), and TNF-α (10 ng/ml). TNF-α, an NF-κB activator, was used as a positive control. NF-κB-induced SEAP activity was assessed using HEK-Blue™ Detection 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