

Validation data for HEK-Blue™ mDectin-2 cells

For research use only

Version # 15B18-MM

HEK-Blue™ mDectin-2 cells are engineered HEK293 cells that stably express the mouse Dectin-2, as well as the genes of the Dectin-2-NF-κB signaling pathway. In addition these cells express an NF-κB-inducible SEAP (secreted embryonic alkaline phosphatase) reporter gene. These reporter cells are activated by Dectin-2 ligands. They do not respond to other CLR ligands such as trehalose-6,6-dibehenate (TDB), a Mincle ligand (see figure 1).

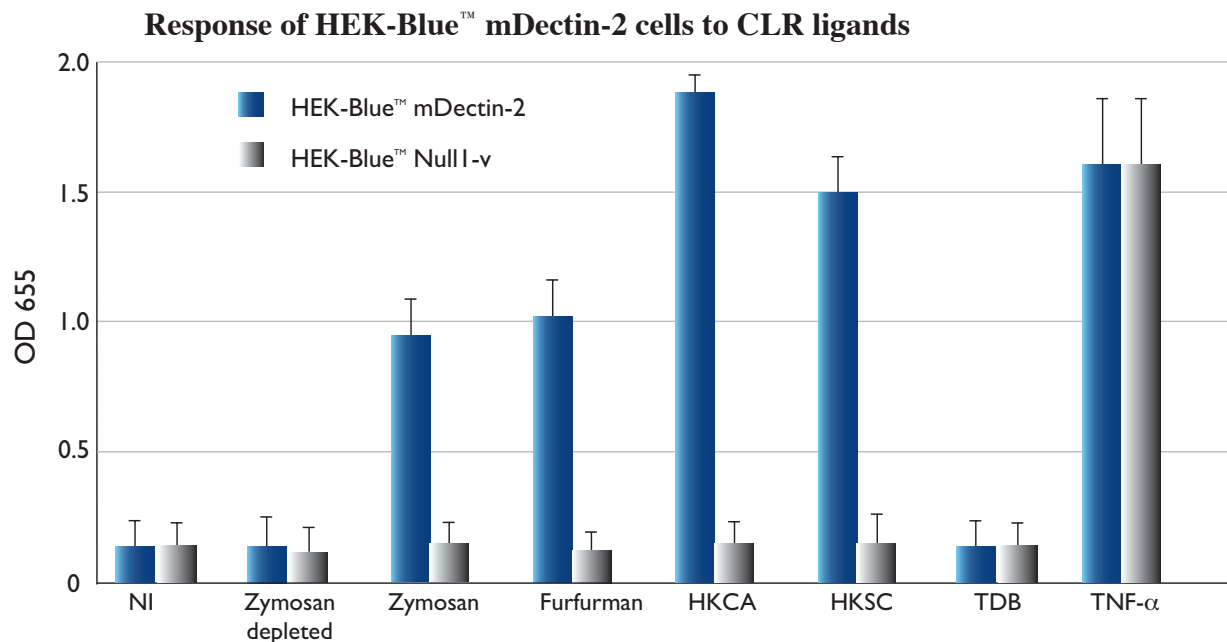


Figure 1: HEK-Blue™ mDectin-2 and HEK-Blue™ Null1-v cells (parental cell line) were stimulated with various CLR agonists: zymosan depleted (100 µg/ml), zymosan (10 µg/ml), furfurman (10 µg/ml), HKCA (1x10⁸ cells), HKSC (1x10⁸ cells), TDB (10 µg/ml), and TNF-α (100 ng/ml). TNF-α, an NF-κB activator, was 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. Non-induced cells (NI) served as negative controls.

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