

# Validation data for HEK-Blue™ mDectin-1b cells

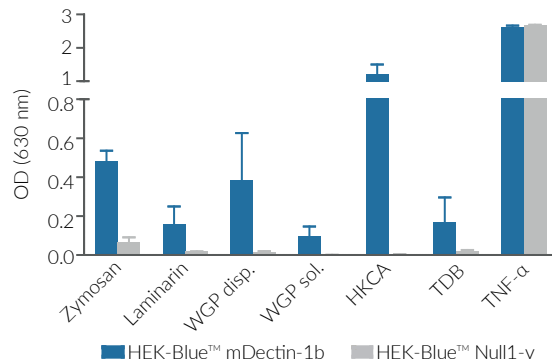
<http://www.invivogen.com/hek-blue-mdectin1b>

For research use only

Version 18C08-NJ

Dectin-1 is alternatively spliced into 2 major isoforms: a full-length A isoform and a 'stalkless' B isoform, which do not induce the same response to soluble and particulate  $\beta$ -glucans. We have engineered HEK-Blue cells that stably express high levels of murine Dectin-1b isoform and an NF- $\kappa$ B-inducible secreted alkaline phosphatase (SEAP) reporter gene. These cells also express genes involved in the Dectin-1 signaling pathway leading to NF- $\kappa$ B activation. HEK-Blue™ mDectin-1b reporter cells are activated by Dectin-1 ligands, however, their responses to soluble and particulate  $\beta$ -glucans differ. These cells respond better to particulate ligands such as Heat Killed *Candida albicans* (HKCA), Zymosan, or WGP dispersible, than to soluble ligands, such as Laminarin and WGP soluble. HEK-Blue™ mDectin-1b cells do not respond to other CLR ligands such as trehalose-6,6-dibehenate (TDB), a Mincle ligand. HEK-Blue™ mDectin-1b reporter cells allow to determine the biological activity of soluble and particulate compounds in a specific manner.

## Evaluation of NF- $\kappa$ B responses to Dectin-1 ligands



**NF- $\kappa$ B responses of HEK-Blue™ mDectin-1b and HEK-Blue™ Null 1-v cells (control cell line) to Dectin-1 ligands.** Cells were incubated with particulate ligands such as Zymosan (10  $\mu$ g/ml), WGP dispersible (100  $\mu$ g/ml) and HKCA (3 x 10<sup>6</sup> cells/ml), or soluble ligands such as Laminarin (100  $\mu$ g/ml), WGP soluble (10  $\mu$ g/ml) or TDB (10  $\mu$ g/ml). TNF- $\alpha$  (10 ng/ml) was used as a positive control. After 24h, SEAP activity was assessed in the supernatant using QUANTI-Blue™, by reading the optical density (OD) at 630 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 3622-3480  
E-mail: [info@invivogen.com](mailto:info@invivogen.com)