

Validation data for WGP Dispersible

<https://www.invivogen.com/wgp>

For research use only

Version 22L13-AK

Whole glucan particles (WGP) are β -glucans notable for their ability to modulate the immune response. InvivoGen provides WGP in two forms, WGP Dispersible (WGP® Dispersible), and WGP Soluble (WGP® Soluble). WGP Dispersible is a particulate *Saccharomyces cerevisiae* β -glucan preparation that triggers Dectin-1-dependent responses, including phagocytosis and the induction of pro-inflammatory cytokines. Unlike its soluble control (WGP Soluble), WGP Dispersible acts as a Dectin-1 agonist and efficiently activates human (h)Dectin-1a and hDectin-1b (Figure 1).

WGP® Dispersible and WGP® Soluble are registered trademarks of Biothera - Wellmune®.

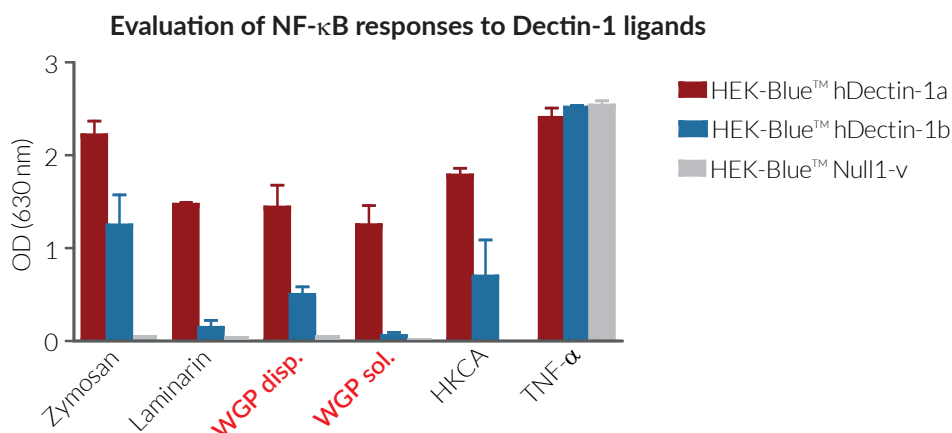


Figure 1. NF- κ B responses to Dectin-1 ligands in HEK-Blue™-derived cell lines. HEK-Blue™ hDectin-1a, HEK-Blue™ hDectin-1b and HEK-Blue™ Null 1-v cells (control cell line) were incubated with particulate ligands such as Zymosan (10 μ g/ml), WGP Dispersible (100 μ g/ml) and HKCA (3×10^6 cells/ml), or soluble ligands such as Laminarin (100 μ g/ml), WGP Soluble (10 μ g/ml) or TDB (10 μ g/ml). TNF- α (10 ng/ml) was used as a positive control. After overnight incubation, the activation of NF- κ B was assessed by measuring the activity of SEAP in the supernatant using QUANTI-Blue™ Solution. Data are shown as optical density (OD) at 630 nm (mean \pm SEM).

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 Asia: +852 3622-3480

E-mail: info@invivogen.com