

Validation data for WGP Soluble

<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 Soluble is a soluble *Saccharomyces cerevisiae* β -glucan preparation that inhibits human (h)Dectin-1b-dependent responses. At the same time, it efficiently binds and activates the hDectin-1a (Figure 1). Furthermore, this soluble ligand displays an inhibitory activity when cells are incubated with particulate agonists such as Zymosan or Heat Killed *Candida albicans* (HKCA) and thus, can be used as a control for Dectin-1b-dependent signaling (Figure 2).

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

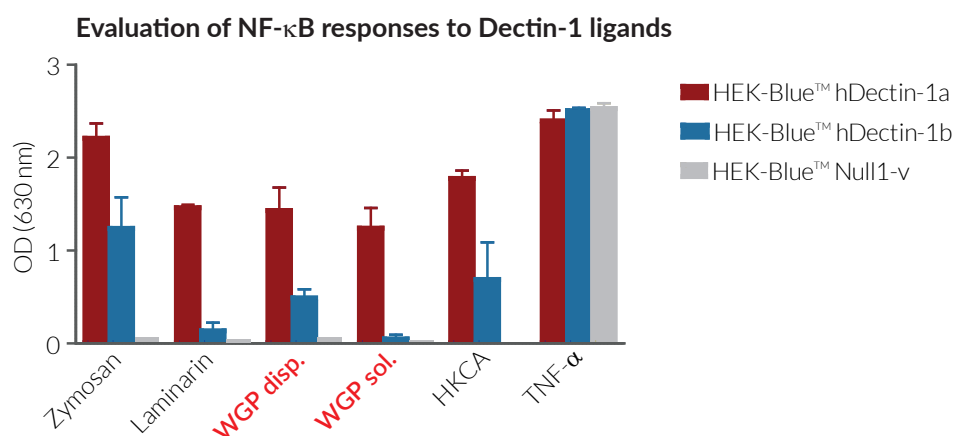


Figure 1. NF- κ B responses to Dectin-1 ligands in HEK-Blue™-derived cell lines. HEK-Blue™ hDectin-1a, HEK-Blue™ hDectin-1b and HEK-Blue™ Null 1-v cells (control cell line) were incubated with particulate ligands such as Zymosan (10 μ g/ml), WGP Dispersible (100 μ g/ml) and HKCA (3 $\times 10^6$ cells/ml), 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).

Dose-dependent inhibition HEK-Blue™ hDectin-1b responses by WGP Soluble

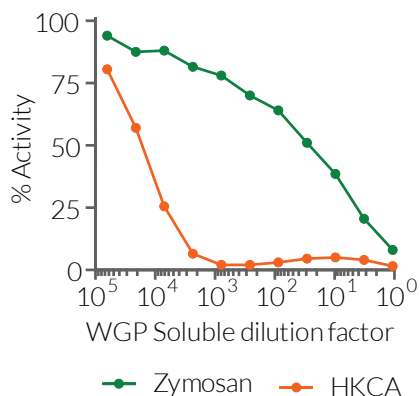


Figure 2. Dose-dependent inhibition HEK-Blue™ hDectin-1b responses by WGP Soluble. Cells were incubated with particulate ligands such as Zymosan (30 μ g/ml) or HKCA (3 $\times 10^6$ cells/ml) together with the soluble ligands WGP Soluble (highest concentration 1 mg/ml). After 24h, SEAP activity was assessed in the supernatant using QUANTI-Blue™ Solution. Data are presented as the percentage of SEAP activity over non-induced cells.

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