

# Validation data for $\beta$ -GlcCer

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

For research use only

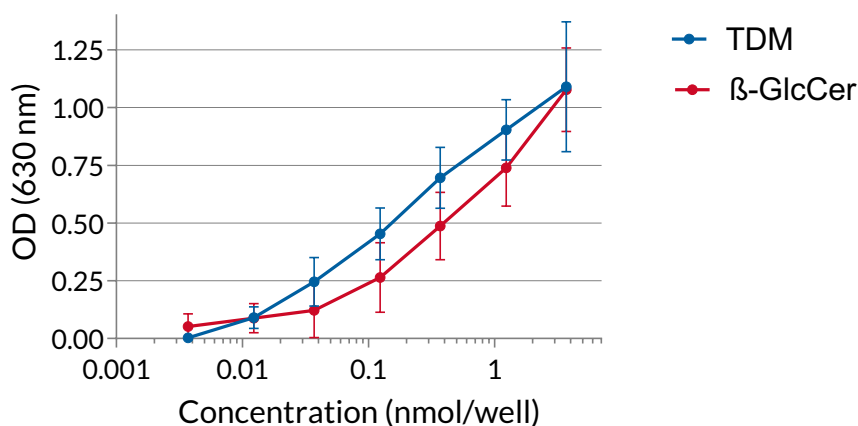
Version 19H05-MM

Beta-glucosylceramide ( $\beta$ -GlcCer) is a glycolipid that activates the macrophage-inducible C-type lectin (Mincle) receptor. Mincle is a member of the Dectin-2 family, that recognizes a variety of exogenous and endogenous stimuli, such as mycobacteria, certain fungi, and necrotic cells.  $\beta$ -GlcCer, an intracellular metabolite, is recognized as a damage-associated molecular pattern (DAMP) and associated with cell death.

Using InvivoGen's HEK-Blue™ Mincle reporter cells, which co-express an NF- $\kappa$ B-inducible SEAP (secreted embryonic alkaline phosphatase) reporter protein and Mincle,  $\beta$ -GlcCer was identified as an endogenous Mincle agonist and compared to the mycobacteria-derived trehalose-6,6-dimycolate (TDM), an extensively studied exogenous Mincle agonist<sup>1</sup>. Stimulation of InvivoGen's HEK-Blue™ hMincle reporter cells with  $\beta$ -GlcCer results in a dose-dependent induction of the NF- $\kappa$ B signaling pathway (data shown below). Upon  $\beta$ -GlcCer recognition, Mincle interacts with the Fc receptor common  $\gamma$ -chain (FcR $\gamma$ ) triggering Syk-dependent signaling resulting in NF- $\kappa$ B, NFAT, and AP-1 activation<sup>1</sup>.

1. Nagata M. *et al.*, 2017. Intracellular metabolite  $\beta$ -glucosylceramide is an endogenous Mincle ligand possessing immunostimulatory activity. PNAS. 114(16): E3285-E3294.

## Evaluation of $\beta$ -GlcCer in HEK-Blue™ hMincle cells



### Response of HEK-Blue™ hMincle cells to $\beta$ -GlcCer and TDM.

HEK-Blue™ hMincle cells were stimulated with increasing concentrations of  $\beta$ -GlcCer and TDM. After overnight incubation, the NF- $\kappa$ B response was determined using QUANTI-Blue™ Solution, a SEAP detection reagent, and 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)