

# Validation data for GlcC<sub>14</sub>C<sub>18</sub>

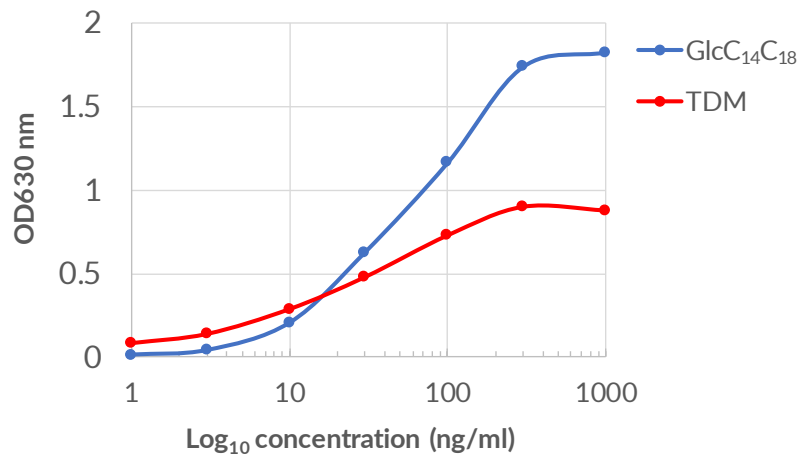
<https://www.invivogen.com/glc-c14c18>

For research use only

Version 18J23-MM

Glucosyl-6-tetradecyloctadecanoate (GlcC<sub>14</sub>C<sub>18</sub>) is a C<sub>6</sub>-branched glycolipid that activates the macrophage-inducible C-type lectin (Mincle) receptor. This glucose monoester was developed using a rational design to obtain a structurally simple molecule that binds and activates Mincle. Upon GlcC<sub>14</sub>C<sub>18</sub> recognition, Mincle interacts with the Fc receptor common  $\gamma$ -chain (FcR $\gamma$ ) triggering Syk-CARD9-dependent signaling and NF- $\kappa$ B activation. Stimulation of InvivoGen's HEK-Blue™ hMincle reporter cells with GlcC<sub>14</sub>C<sub>18</sub> results in a dose-dependent induction of the NF- $\kappa$ B signaling pathway.

## Evaluation of GlcC<sub>14</sub>C<sub>18</sub> in HEK-Blue™ hMincle cells



### GlcC<sub>14</sub>C<sub>18</sub> induces a dose-dependent response in HEK-Blue™ hMincle cells.

HEK-Blue™ hMincle cells were stimulated with increasing concentrations of GlcC<sub>14</sub>C<sub>18</sub> and trehalose-6,6-dimycolate (TDM, also known as cord factor), an extensively studied Mincle agonist. After overnight incubation, the NF- $\kappa$ B response was determined using QUANTI-Blue™, 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 3622-34-80  
E-mail: [info@invivogen.com](mailto:info@invivogen.com)