

Validation data for Hemozoin

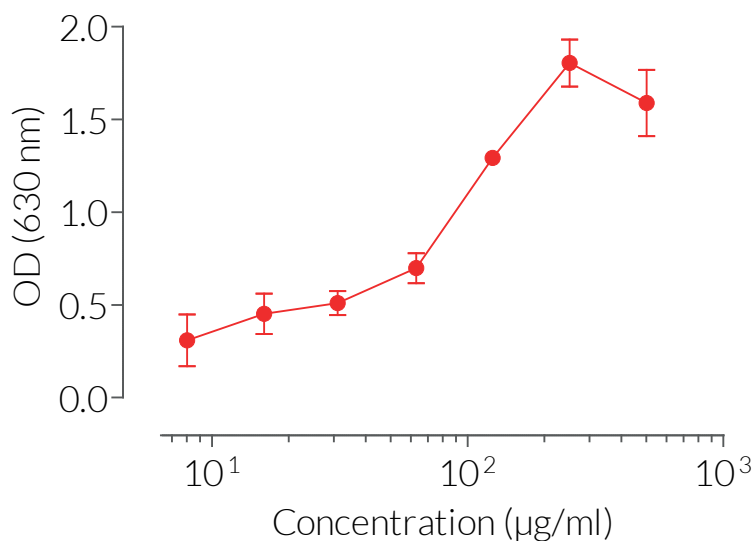
<https://www.invivogen.com/hemozoin>

For research use only

Version 22I28-AK

Hemozoin (HZ) is a metabolically crystallized byproduct generated by the malaria parasite during infection. HZ-induced production of IL-1 β and IL-18 is linked to the activation of the NLRP3 inflammasome. The biological activity of the synthetic hemozoin has been validated using InvivoGen's THP1-Null2 cells. The production of IL-1 β by THP1-Null2 cells was measured using HEK-Blue™ IL-1 β cells. Treatment with hemozoin induced IL-1 β secretion, an indicator of NLRP3 inflammasome activation, in a dose-dependent manner.

Dose-dependent IL-1 β induction



Mature IL-1 β production in THP1-Null2 cells. Human THP-1 monocytes were primed for 3 hours with LPS-EK(10 μg/ml) prior to the incubation with increasing concentrations of Hemozoin. The next day, the supernatant was incubated with HEK-Blue™ IL-1 β cells for 24 hours and IL-1 β was assessed using QUANTI-Blue™ Solution. The optical density (OD) was read 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 Asia: +852 3622-3480

E-mail: info@invivogen.com