

Validation data for Indirubin

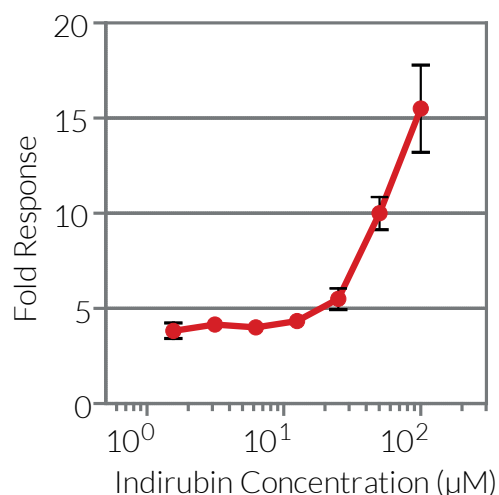
<https://www.invivogen.com/indirubin>

For research use only

Version 19E07-NJ

Indirubin (2-(2-oxo-1H-indol-3-ylidene)-1H-indol-3-one) is a dietary-derived indole-based ligand of the cytosolic aryl hydrocarbon receptor (AhR). Indirubin is a more potent agonist of human AhR than murine AhR. Stimulation of InvivoGen's HepG2-Lucia™ AhR cells with indirubin results in a dose-dependent induction of the AhR genomic signaling pathway, evidenced by the Cyp1a1-induced expression of the Lucia luciferase reporter.

Evaluation of cellular AhR activation by indirubin



Indirubin induces a dose-dependent response in HepG2-Lucia™ AhR cells. The cells were incubated with increasing concentrations of indirubin. After overnight incubation, the AhR activation was assessed by determining Lucia luciferase activity in the supernatant using QUANTI-Luc™. Data are expressed as fold responses, as compared to non-induced cells. DMSO final concentration impacts cell viability and AhR activation. We recommend to include a control dilution of DMSO in your assay. Using our protocol, a 100 µM indirubin concentration includes 2% DMSO.

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