

# Validation data for FICZ

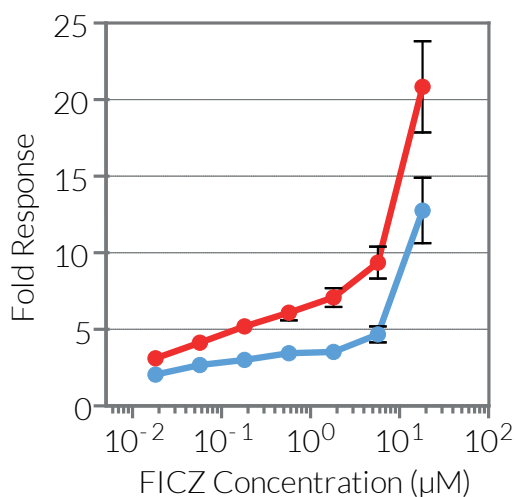
<https://www.invivogen.com/ficz>

For research use only

Version 23A19-MM

FICZ (6-formylindolol[3,2-b]carbazole) is a highly potent agonist for the cytosolic aryl hydrocarbon receptor (AhR) that results from tryptophan conversion upon UV-dependent photo-oxidation or H<sub>2</sub>O<sub>2</sub>-mediated oxidative stress. Stimulation of InvivoGen's HepG2-Lucia™-AhR and HT29-Lucia™-AhR cells with FICZ results in a dose-dependent induction of the AhR genomic signaling pathway, evidenced by the Cyp1a1-induced expression of Lucia luciferase reporter.

## Evaluation of cellular AhR activation by FICZ



—●— HepG2-Lucia™ AhR Cells    —●— HT29-Lucia™ AhR Cells

FICZ induces a dose-dependent response in HepG2-Lucia™ AhR and HT29-Lucia™ AhR cells. The cells were incubated with increasing concentrations of FICZ. After overnight incubation, the AhR activation was assessed by determining Lucia luciferase activity in the supernatant using QUANTI-Luc™ 4 Lucia/Gaussia. Data are expressed as fold responses, as compared to 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](mailto:info@invivogen.com)