

Validation data for CH-223191

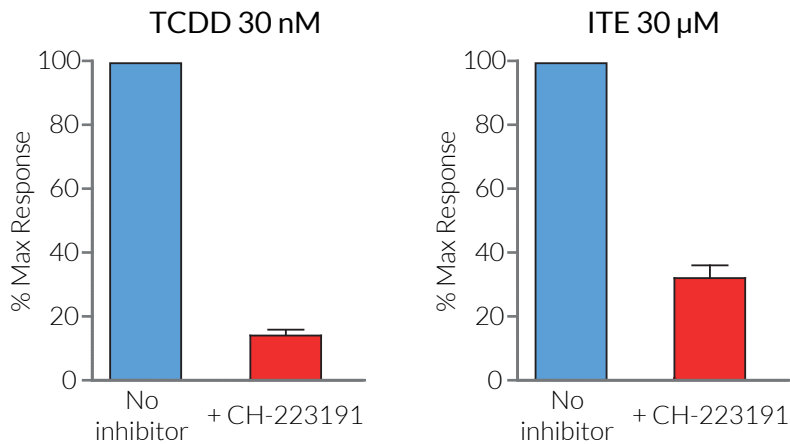
<https://www.invivogen.com/ch223191>

For research use only

Version 23A19-MM

CH-223191 is a synthetic antagonist for the cytosolic aryl hydrocarbon receptor (AhR) that was first described as a competitive ligand of the xenobiotic 2,3,7,8-tetrachlorodibenzo-p-dioxin (TCDD). CH-223191 blocks TCDD-mediated nuclear translocation and DNA binding of AhR. This AhR inhibitor exerts a ligand-selective antagonism and appears to be more effective on halogenated aromatic hydrocarbons such as TCDD, than on polycyclic aromatic hydrocarbons and non halogenated aromatic hydrocarbons such as FICZ (6-formylindolo[3,2-b]carbazole) and ITE (2-(1'H-indole-3'-carbonyl)-thiazole-4-carboxylic acid methyl ester), respectively. Co-incubation of InvivoGen's HepG2-Lucia™-AhR cells with CH-223191, and TCDD or ITE, results in a significant inhibition (70% to 85%) of the AhR genomic signaling pathway, evidenced by a decreased Cyp1a1-induced expression of Lucia luciferase reporter.

Evaluation of cellular AhR inhibition by CH-223191



Inhibition of AhR activity by CH-223191 in HepG2-Lucia™ AhR cells. The cells were stimulated with 30 nM TCDD or 30 μM ITE in the presence of 10 μM CH-223191. After overnight incubation, the AhR response was assessed by determining Lucia luciferase activity in the supernatant using QUANTI-Luc™ 4 Lucia/Gaussia. Percentages of the maximal response for the ligand with no inhibitor are shown.

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