

# PRODUCT INFORMATION

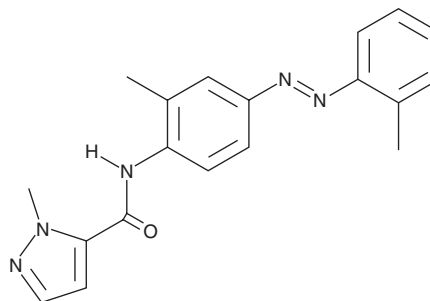


CH 223191

Item No. 16154

CAS Registry No.: 301326-22-7  
Formal Name: 1-methyl-N-[2-methyl-4-[2-(2-methylphenyl)diazenyl]phenyl]-1H-pyrazole-5-carboxamide

MF: C<sub>19</sub>H<sub>19</sub>N<sub>5</sub>O  
FW: 333.4  
Purity: ≥95%  
UV/Vis.: λ<sub>max</sub>: 233, 342 nm  
Supplied as: A crystalline solid  
Storage: -20°C  
Stability: ≥4 years



Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

## Laboratory Procedures

CH 223191 is supplied as a crystalline solid. A stock solution may be made by dissolving the CH 223191 in the solvent of choice, which should be purged with an inert gas. CH 223191 is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide. The solubility of CH 223191 in these solvents is approximately 0.1, 25, and 30 mg/ml, respectively.

## Description

The aryl hydrocarbon receptor (AhR) is a ligand-activated transcription factor that up-regulates many xenobiotic metabolizing enzymes, particularly in response to planar aromatic hydrocarbons.<sup>1</sup> CH 223191 is a potent and specific antagonist of AhR that blocks activation by 2,3,7,8-tetrachlorodibenzo-*p*-dioxin (TCDD, IC<sub>50</sub> = 0.03 μM).<sup>2</sup> It prevents the induction of cytochrome P450 1A1 by TCDD in HepG2 cells and in livers of mice.<sup>2</sup> Through its effects on AhR, CH 223191 suppresses Th17 cell differentiation both *in vitro* and *in vivo* and alters the expression of homeobox transcription factors necessary for the differentiation of embryonic stem cells to cardiomyocytes.<sup>3,4</sup>

## References

1. Gu, Y.Z., Hogenesch, J.B., and Bradfield, C.A. The PAS superfamily: Sensors of environmental and developmental signals. *Annu. Rev. Pharmacol. Toxicol.* **40**, 519-561 (2000).
2. Kim, S.-H., Henry, E.C., Kim, D.-K., *et al.* Novel compound 2-methyl-2H-pyrazole-3-carboxylic acid (2-methyl-4-*o*-tolylazo-phenyl)-amide (CH-223191) prevents 2,3,7,8-TCDD-induced toxicity by antagonizing the aryl hydrocarbon receptor. *Mol. Pharmacol.* **69**(6), 1871-1878 (2006).
3. Veldhoen, M., Hirota, K., Christensen, J., *et al.* Natural agonists for aryl hydrocarbon receptor in culture medium are essential for optimal differentiation of Th17 T cells. *J. Exp. Med.* **206**(1), 43-49 (2009).
4. Wang, Q., Chen, J., Ko, C.I., *et al.* Disruption of aryl hydrocarbon receptor homeostatic levels during embryonic stem cell differentiation alters expression of homeobox transcription factors that control cardiomyogenesis. *Environ. Health Perspect.* **121**(11-12), 1334-1343 (2013).

### WARNING

THIS PRODUCT IS FOR RESEARCH ONLY - NOT FOR HUMAN OR VETERINARY DIAGNOSTIC OR THERAPEUTIC USE.

### SAFETY DATA

This material should be considered hazardous until further information becomes available. Do not ingest, inhale, get in eyes, on skin, or on clothing. Wash thoroughly after handling. Before use, the user must review the complete Safety Data Sheet, which has been sent via email to your institution.

### WARRANTY AND LIMITATION OF REMEDY

Buyer agrees to purchase the material subject to Cayman's Terms and Conditions. Complete Terms and Conditions including Warranty and Limitation of Liability information can be found on our website.

Copyright Cayman Chemical Company, 09/28/2022

## CAYMAN CHEMICAL

1180 EAST ELLSWORTH RD  
ANN ARBOR, MI 48108 · USA

PHONE: [800] 364-9897  
[734] 971-3335

FAX: [734] 971-3640

CUSTSERV@CAYMANCHEM.COM  
WWW.CAYMANCHEM.COM