

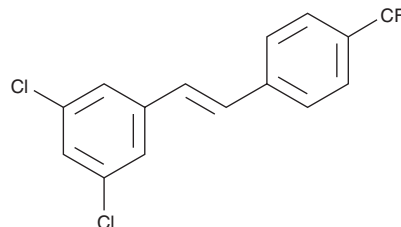
# PRODUCT INFORMATION



## CAY10465

Item No. 10006546

**CAS Registry No.:** 688348-33-6  
**Formal Name:** 1,3-dichloro-5-[(1E)-2-[4-(trifluoromethyl)phenyl]ethenyl]-benzene  
**MF:** C<sub>15</sub>H<sub>9</sub>Cl<sub>2</sub>F<sub>3</sub>  
**FW:** 317.1  
**Purity:** ≥98%  
**UV/Vis.:** λ<sub>max</sub>: 211, 229, 245, 298, 309 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

CAY10465 is supplied as a crystalline solid. A stock solution may be made by dissolving the CAY10465 in the solvent of choice, which should be purged with an inert gas. CAY10465 is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF). The solubility of CAY10465 in ethanol is approximately 20 mg/ml and approximately 30 mg/ml in DMSO and DMF.

CAY10465 is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, CAY10465 should first be dissolved in DMSO and then diluted with the aqueous buffer of choice. CAY10465 has a solubility of approximately 0.5 mg/ml in a 1:1 solution of DMSO:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

### Description

The aryl hydrocarbon receptor (AhR) is a ligand-dependent intracellular transcription factor whose ligands include some of the most infamous xenobiotics, including dioxin, benzo[a]pyrene, and numerous polyaromatics from soot and coal tar.<sup>1</sup> CAY10465 is an analog of resveratrol acting as a potent and selective AhR agonist, with a K<sub>i</sub> of 0.2 nM.<sup>2</sup> CAY10465 is inactive as a ligand for the estrogen receptor even at 100 μM.

### References

1. Denison, M.S. and Nagy, S.R. Activation of the aryl hydrocarbon receptor by structurally diverse exogenous and endogenous chemicals. *Annu. Rev. Pharmacol. Toxicol.* **43**, 309-334 (2003).
2. de Medina, P., Casper, R., Savouret, J.-F., *et al.* Synthesis and biological properties of new stilbene derivatives of resveratrol as new selective aryl hydrocarbon modulators. *J. Med. Chem.* **48(1)**, 287-291 (2005).

#### 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, 11/13/2023

#### 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