

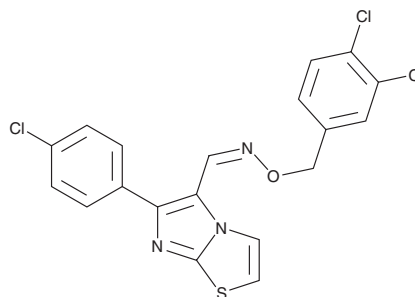
# PRODUCT INFORMATION



## Z-CITCO

Item No. 39172

**CAS Registry No.:** 2583138-04-7  
**Formal Name:** [C(Z)]-6-(4-chlorophenyl)-imidazo[2,1-b]thiazole-5-carboxaldehyde O-[(3,4-dichlorophenyl)methyl]oxime  
**MF:** C<sub>19</sub>H<sub>12</sub>Cl<sub>3</sub>N<sub>3</sub>OS  
**FW:** 436.7  
**Purity:** ≥95%  
**UV/Vis.:** λ<sub>max</sub>: 203, 257, 315 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

Z-CITCO is supplied as a crystalline solid. A stock solution may be made by dissolving the Z-CITCO in the solvent of choice, which should be purged with an inert gas. Z-CITCO is soluble in organic solvents such as DMSO and dimethyl formamide. The solubility of Z-CITCO in these solvents is approximately 10 and 20 mg/ml, respectively.

### Description

Z-CITCO is an agonist of the constitutive androstane receptor (CAR; EC<sub>50</sub> = 3.9 μM for the human receptor in a reporter assay) and the *cis* isomer of the CAR agonist CITCO (Item No. 16027).<sup>1</sup>

### Reference

1. Diethelm-Varela, B., Kumar, A., Lynch, C., *et al.* Stereoisomerization of human constitutive androstane receptor agonist CITCO. *Tetrahedron* **79**, 131886 (2021).

#### 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, 06/30/2023

#### 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](http://WWW.CAYMANCHEM.COM)