

PRODUCT INFORMATION



Chloroguanabenz (acetate)

Item No. 26423

Formal Name: (E)-2-(2,3,6-trichlorobenzylidene)hydrazine-1-carboximidamide, monoacetate

MF: C₈H₇Cl₃N₄ • C₂H₄O₂

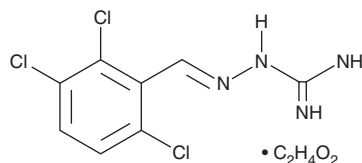
FW: 325.6

Purity: ≥95%

Supplied as: A 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

Chloroguanabenz (acetate) is supplied as a solid. A stock solution may be made by dissolving the chloroguanabenz (acetate) in the solvent of choice, which should be purged with an inert gas. Chloroguanabenz (acetate) is soluble in organic solvents such as ethanol and DMSO.

Description

Chloroguanabenz is an antiprion agent and a derivative of the α_2 -adrenergic receptor agonist guanabenz (Item No. 10851).¹ It inhibits prion formation in yeast and mammalian cell-based assays. Chloroguanabenz reduces the levels of both soluble and aggregated forms of the truncated Huntingtin derivative Htt48 in an HEK293T cellular model of Huntington's disease.²

References

1. Tribouillard-Tanvier, D., Béringue, V., Desban, N., *et al.* Antihypertensive drug guanabenz is active in vivo against both yeast and mammalian prions. *PLoS One* **3(4)**, e1981 (2008).
2. Bertolotti, A. and Blondel, M. Use of chlorine guanabenz derivatives for treating polyglutamine expansion associated diseases. *Centre National de la Recherche Scientifique*. **WO2008/041133** (2011).

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, 12/20/2022

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