

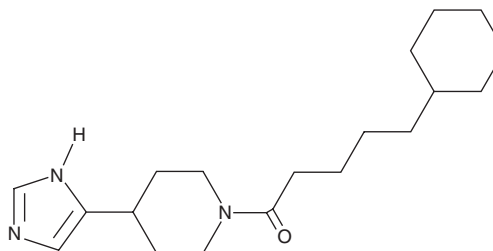
PRODUCT INFORMATION



GT 2016

Item No. 11996

CAS Registry No.: 152241-24-2
Formal Name: 5-cyclohexyl-1-[4-(1H-imidazol-5-yl)-1-piperidinyl]-1-pentanone
MF: $C_{19}H_{31}N_3O$
FW: 317.5
Purity: $\geq 98\%$
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

GT 2016 is supplied as a crystalline solid. A stock solution may be made by dissolving the GT 2016 in the solvent of choice, which should be purged with an inert gas. GT 2016 is soluble in organic solvents such as ethanol and DMSO.

Description

GT 2016 is an antagonist of histamine H_3 receptors ($K_i = 43.8$ nM).¹ It is selective for H_3 over H_1 and H_2 receptors ($IC_{50} = >10,000$ nM for both) and lacks activity against kidney and brain histamine methyltransferase at concentrations up to $3 \mu\text{M}$. *In vivo*, GT 2016 (10 and 30 mg/kg) increases histamine release in the rat parietal cortex.

Reference

1. Tedford, C.E., Yates, S.L., Pawlowski, G.P., *et al.* Pharmacological characterization of GT-2016, a non-thiourea-containing histamine H_3 receptor antagonist: *In vitro* and *in vivo* studies. *J. Pharmacol. Exp. Ther.* **275**(2), 598-604 (1995).

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, 11/08/2022

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