

PRODUCT INFORMATION

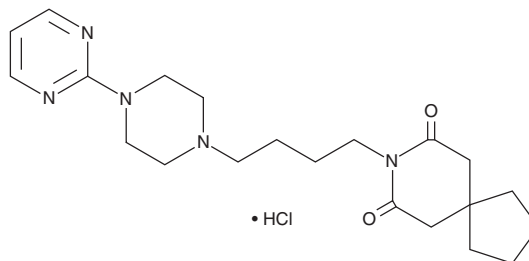


Buspirone (hydrochloride)

Item No. 22028

CAS Registry No.: 33386-08-2
Formal Name: 8-[4-[4-(2-pyrimidinyl)-1-piperazinyl]butyl]-8-azaspiro[4.5]decane-7,9-dione, monohydrochloride

Synonym: MJ9022-1
MF: C₂₁H₃₁N₅O₂ • HCl
FW: 422.0
Purity: ≥98%
UV/Vis.: λ_{max}: 237, 299 nm
Supplied as: A crystalline solid
Storage: -20°C
Stability: ≥5 years



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

Description

Buspirone (hydrochloride) (Item No. 22028) is an analytical reference standard that is categorized as an anxiolytic.¹ It modulates serotonin and dopamine receptors, undergoes rapid first-pass metabolism, and is metabolized to phenylpiperazine (Item No. 11203), which has high affinity for α-adrenergic receptors.² This product is intended for research and forensic applications.

References

1. Gale, C. and Oakley-Brown, M. Anxiety disorder. *B.M.J.* **321(7270)**, 1204-1207 (2000).
2. Rënyi, L., Evenden, J.L., Fowler, C.J., et al. The pharmacological profile of (R)-3,4-dihydro-N-isopropyl-3-(N-isopropyl-N-propylamino)-2H-1-benzopyran-5-carboxamide, a selective 5-hydroxytryptamine_{1A} receptor agonist. *J. Pharmacol. Exp. Ther.* **299(3)**, 883-893 (2001).

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, 06/13/2023

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