

PRODUCT INFORMATION

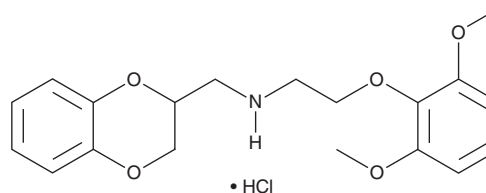


WB 4101 (hydrochloride)

Item No. 29528

CAS Registry No.: 2170-58-3
Formal Name: N-[2-(2,6-dimethoxyphenoxy)ethyl]-2,3-dihydro-1,4-benzodioxin-2-methanamine, monohydrochloride

MF: $C_{19}H_{23}NO_5 \cdot HCl$
FW: 381.9
Purity: $\geq 95\%$
Supplied as: A crystalline solid
Storage: $-20^{\circ}C$
Stability: ≥ 4 years



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

Laboratory Procedures

WB 4101 (hydrochloride) is supplied as a crystalline solid. A stock solution may be made by dissolving the WB 4101 (hydrochloride) in the solvent of choice, which should be purged with an inert gas. WB 4101 (hydrochloride) is soluble in organic solvents such as DMSO and dimethyl formamide. The solubility of WB 4101 (hydrochloride) in these solvents is approximately 1 and 5 mg/ml, respectively.

Description

WB 4101 is an antagonist of α -adrenergic receptors (α -ARs).¹ It inhibits norepinephrine-induced contraction of isolated rat vas deferens ($pA_2 = 9.8$). WB 4101 also inhibits contractions induced by carbachol (carbamoylcholine; Item No. 14486) in isolated guinea pig taenia caeci ($pA_2 = 8.9$).²

References

1. Mottram, D.R. and Kapur, H. The α -adrenoceptor blocking effects of a new benzodioxane. *J. Pharm. Pharmacol.* **27**(4), 295-296 (1975).
2. Butler, M. and Jenkinson, D.H. Blockade by WB 4101 of α -adrenoceptors in the rat vas deferens and guinea-pig taenia caeci. *Eur. J. Pharmacol.* **52**(3-4), 303-311 (1978).

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, 10/06/2022

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