

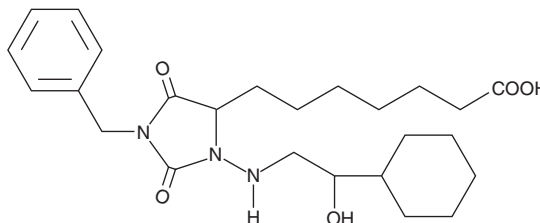
PRODUCT INFORMATION



BW A868C

Item No. 12060

CAS Registry No.: 118675-50-6
Formal Name: 3-[(2-cyclohexyl-2-hydroxyethyl)amino]-2,5-dioxo-1-(phenylmethyl)-4-imidazolidineheptanoic acid
MF: C₂₅H₃₇N₃O₅
FW: 459.6
Purity: ≥98%
Supplied as: A solution in ethanol
Storage: -20°C
Stability: ≥2 years



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

Laboratory Procedures

BW A868C is supplied as a solution in ethanol. To change the solvent, simply evaporate the ethanol under a gentle stream of nitrogen and immediately add the solvent of choice. Solvents such as DMSO and dimethyl formamide purged with an inert gas can be used. The solubility of BW A868C in these solvents is approximately 20 and 30 mg/ml, respectively.

BW A868C is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, the ethanolic solution of BW A868C should be diluted with the aqueous buffer of choice. The solubility of BW A868C in PBS (pH 7.2) is approximately 0.25 mg/ml. We do not recommend storing the aqueous solution for more than one day.

Description

BW A868C is a hydantoin compound that is structurally related to the DP receptor agonist BW 245C.^{1,2} BW A868C antagonizes the prostaglandin D₂ and BW 245C-induced activation of human platelet adenylate cyclase. It reverses the anti-aggregatory effects of DP receptor agonists in a dose-dependent manner. Likewise, BW A868C antagonizes the accumulation of cAMP in rabbit non-pigmented ciliary epithelial cells.³ It has virtually no effect on human TP, IP, EP₁, EP₂, and FP receptors. The K_i for BW A868C is approximately 1.7 nM.⁴

References

1. Caldwell, A.G., Harris, C.J., Stepney, R., *et al.* Hydantoin prostaglandin analogues, potent and selective inhibitors of platelet aggregation. *J. Chem. Soc. Chem. Comm.* 561-562 (1979).
2. Trist, D.G., Collins, B.A., Wood, J., *et al.* The antagonism by BW A868C of PGD₂ and BW245C activation of human platelet adenylate cyclase. *Br. J. Pharmacol.* **96**(2), 301-306 (1989).
3. Bhattacharjee, P., Jacobs, N., Coca-Prados, M., *et al.* Identification of prostanoid receptors in rabbit non-pigmented ciliary epithelial cells. *Exp. Eye Res.* **62**(5), 491-498 (1996).
4. Giles, H., Leff, P., Bolofo, M.L., *et al.* The classification of prostaglandin DP-receptors in platelets and vasculature using BW A868C, a novel, selective and potent competitive antagonist. *Br. J. Pharmacol.* **96**(2), 291-300 (1989).

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, 02/27/2024

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