

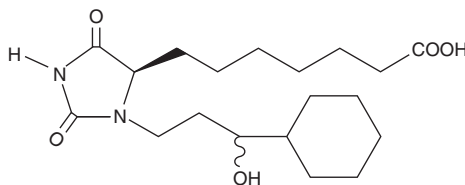
PRODUCT INFORMATION



BW 246C

Item No. 12055

CAS Registry No.: 65705-83-1
Formal Name: (4R)-(3-[(3R,S)-3-cyclohexyl-3-hydroxypropyl]-2,5-dioxo)-4-imidazolidineheptanoic acid
Synonym: 8-*epi* BW 245C
MF: C₁₉H₃₂N₂O₅
FW: 368.5
Purity: ≥98%
Stability: ≥1 year at -20°C
Supplied as: A crystalline solid



Laboratory Procedures

For long term storage, we suggest that BW 246C be stored as supplied at -20°C. It should be stable for at least one year.

BW 246C is supplied as a crystalline solid. A stock solution may be made by dissolving the BW 246C in an organic solvent purged with an inert gas. BW 246C is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide. The solubility of BW 246C in these solvents is approximately 50 mg/ml.

Further dilutions of the stock solution into aqueous buffers or isotonic saline should be made prior to performing biological experiments. Ensure that the residual amount of organic solvent is insignificant, since organic solvents may have physiological effects at low concentrations. Organic solvent-free aqueous solutions of BW 246C can be prepared by directly dissolving the crystalline compound in aqueous buffers. The solubility of BW 246C in PBS (pH 7.2) is approximately 2.2 mg/ml. We do not recommend storing the aqueous solution for more than one day.

Description

BW 245C (Item No. 12050) is the less active C-8 diastereomer DP₁ receptor agonist BW 246C.^{1,2} The activity of BW 245C is 70-fold greater than that of BW 246C.³

References

1. Boie, Y., Sawyer, N., Slipetz, D.M., *et al.* Molecular cloning and characterization of the human prostanoid DP receptor. *J. Biol. Chem.* **270**, 18910-18916 (1995).
2. Town, M., Casals-Stenzel, J., Schillinger, E. Pharmacological and cardiovascular properties of a hydantoin derivative, BW 245 C, with high affinity and selectivity for PGD₂ receptors. *Prostaglandins* **25**, 13-28 (1983).
3. Caldwell, A.G., Harris, C.J., Stepney, R., *et al.* Heterocyclic prostaglandin analogues. Part 2. Hydantoins and other imidazole analogues. *J. Chem. Soc. Perkin Trans. I* 495-505 (1980).

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, 08/11/2015

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