

PRODUCT INFORMATION

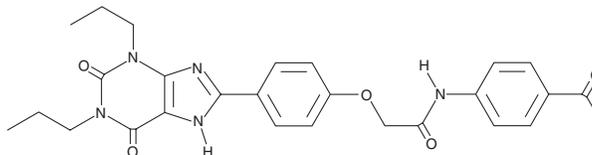


MRS1706

Item No. 17417

CAS Registry No.: 264622-53-9
Formal Name: N-(4-acetylphenyl)-2-[4-(2,3,6,9-tetrahydro-2,6-dioxo-1,3-dipropyl-1H-purin-8-yl)phenoxy]-acetamide

MF: C₂₇H₂₉N₅O₅
FW: 503.6
Purity: ≥95%
UV/Vis.: λ_{max}: 252, 307 nm
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

MRS1706 is supplied as a crystalline solid. A stock solution may be made by dissolving the MRS1706 in the solvent of choice, which should be purged with an inert gas. MRS1706 is soluble in DMSO at a concentration of approximately 0.3 mg/ml.

Description

MRS1706 is a selective adenosine A_{2B} receptor inverse agonist with K_i values of 1.39, 157, 112, and 230 nM for human A_{2B}, A₁, A_{2A}, and A₃ receptors, respectively.^{1,2}

References

1. Kim, Y.-C., Ji, X., Melman, N., *et al.* Anilide derivatives of an 8-phenylxanthine carboxylic congener are highly potent and selective antagonists at human A(2B) adenosine receptors. *J. Med. Chem.* **43(6)**, 1165-1172 (2000).
2. Li, Q., Ye, K., Blad, C.C., *et al.* ZM241385, DPCPX, MRS1706 are inverse agonists with different relative intrinsic efficacies on constitutively active mutants of the human adenosine A_{2B} receptor. *J. Pharmacol. Exp. Ther.* **320(2)**, 637-645 (2007).

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, 10/12/2022

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