

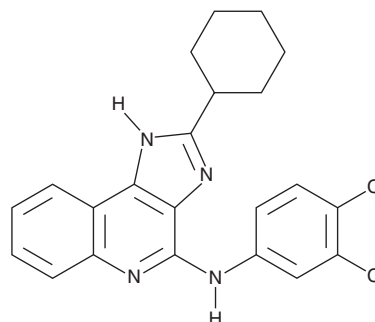
PRODUCT INFORMATION



LUF6000

Item No. 28437

CAS Registry No.: 890087-21-5
Formal Name: 2-cyclohexyl-N-(3,4-dichlorophenyl)-3H-imidazo[4,5-c]quinolin-4-amine
MF: C₂₂H₂₀Cl₂N₄
FW: 411.3
Purity: ≥98%
UV/Vis.: λ_{max}: 228, 264, 324, 337 nm
Supplied as: A 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

LUF6000 is supplied as a solid. A stock solution may be made by dissolving the LUF6000 in the solvent of choice, which should be purged with an inert gas. LUF6000 is soluble in the organic solvent DMSO.

Description

LUF6000 is a positive allosteric modulator of adenosine A₃ receptors.¹ It reduces dissociation of the agonist I-AB-MECA from A₃ receptors and increases the efficacy, but not potency, of 2-Cl-IB-MECA-induced inhibition of forskolin-stimulated cAMP accumulation by 45% in CHO cells expressing the human receptor when used at a concentration of 10 μM. LUF6000 (100 μg/kg, three times per day) reduces joint edema and decreases levels of PI3K, IKK, IκB, and NF-κB in peripheral blood mononuclear cells (PBMCs) in a rat model of adjuvant-induced arthritis.² It reduces knee edema in a rat model of osteoarthritis induced by monosodium iodoacetate (MIA) when administered at a dose of 100 μg/kg twice per day.

References

1. Göblyös, A., Gao, Z.-G., Brussee, J., *et al.* Structure-activity relationships of new 1H-imidazo[4,5-c]quinolin-4-amine derivatives as allosteric enhancers of the A₃ adenosine receptor. *J. Med. Chem.* **49**(11), 3354-3361 (2006).
2. Cohen, S., Barer, F., Bar-Yehuda, S., *et al.* A₃ adenosine receptor allosteric modulator induces an anti-inflammatory effect: *In vivo* studies and molecular mechanism of action. *Mediators Inflamm.* **2014:708746** (2014).

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, 12/05/2022

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