

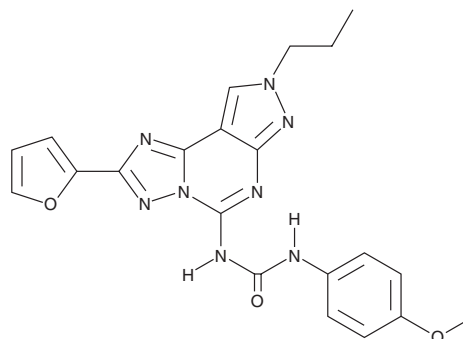
PRODUCT INFORMATION



MRE 3008F20

Item No. 43169

CAS Registry No.: 252979-43-4
Formal Name: N-[2-(2-furanyl)-8-propyl-8H-pyrazolo[4,3-e][1,2,4]triazolo[1,5-c]pyrimidin-5-yl]-N'-(4-methoxyphenyl)-urea
MF: C₂₁H₂₀N₈O₃
FW: 432.4
Purity: ≥95%
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

MRE 3008F20 is supplied as a solid. A stock solution may be made by dissolving the MRE 3008F20 in the solvent of choice, which should be purged with an inert gas. MRE 3008F20 is sparingly soluble (1-10 mg/ml) in DMSO.

Description

MRE 3008F20 is an adenosine A₃ receptor antagonist (K_i = 0.8 nM for the human receptor).¹ It is selective for the adenosine A₃ receptor over adenosine A₁, -A_{2A}, and -A_{2B} receptors (K_is = 1,197, 140, and 2,056 nM, respectively, for the human receptors). MRE 3008F20 inhibits cAMP accumulation induced by IB-MECA (piclidenosan; Item No. 28445) in CHO cells expressing the human adenosine A₃ receptor (IC₅₀ = 5.2 nM). It reverses decreases in NF-κB activation and TNF-α and IL-8 levels induced by NECA (5'-N-ethylcarboxamidoadenosine; Item No. 21420) or 2-Cl-IB-MECA (Item No. 27336) in LPS-stimulated primary human osteoarthritic synoviocytes when used at a concentration of 1 μM.²

References

1. Baraldi, P.G., Cacciari, B., Romagnoli, R., *et al.* Pyrazolo[4,3-e]1,2,4-triazolo[1,5-c]pyrimidine derivatives as highly potent and selective human A₃ adenosine receptor antagonists: influence of the chain at the N8 pyrazole nitrogen. *J. Med. Chem.* **43**, 4768-4780 (2000).
2. Varani, K., Vincenzi, F., Tosi, A., *et al.* Expression and functional role of adenosine receptors in regulating inflammatory responses in human synoviocytes. *Br. J. Pharmacol.* **160**(1), 101-115 (2010).

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, 05/12/2025

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