

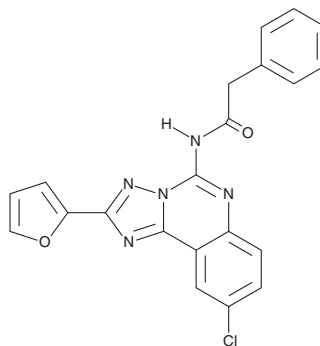
PRODUCT INFORMATION



MRS1220

Item No. 41015

CAS Registry No.: 183721-15-5
Formal Name: N-[9-chloro-2-(2-furanyl)[1,2,4]triazolo[1,5-c]quinazolin-5-yl]-benzeneacetamide
Synonym: NSC 695674
MF: C₂₁H₁₄ClN₅O₂
FW: 403.8
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

MRS1220 is supplied as a solid. A stock solution may be made by dissolving the MRS1220 in the solvent of choice, which should be purged with an inert gas. MRS1220 is slightly soluble (0.1-1 mg/ml) in acetonitrile and sparingly soluble (1-10 mg/ml) in chloroform.

Description

MRS1220 is an adenosine A₃ receptor antagonist (K_i = 0.65 nM for the human receptor).¹ It is selective for human adenosine A₃ receptors over rat adenosine A₁ and A₂ receptors (K_is = 305 and 52 nM, respectively). MRS1220 inhibits [³⁵S]GTPγS binding in HEK293 cells expressing human adenosine A₃ receptors (IC₅₀ = 7.2 nM). *In vivo*, MRS1220 (0.15 mg/kg) decreases tumor growth and angiogenesis in a rat C6 glioma model.²

References

1. Jacobson, K.A., Park, K.S., Jiang, J.L., *et al.* Pharmacological characterization of novel A₃ adenosine receptor-selective antagonists. *Neuropharmacology* **36**(9), 1157-1165 (1997).
2. Rocha, R., Torres, A., Ojeda, K., *et al.* The adenosine A₃ receptor regulates differentiation of glioblastoma stem-like cells to endothelial cells under hypoxia. *Int. J. Mol. Sci.* **19**(4), 1228 (2018).

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, 06/21/2024

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