

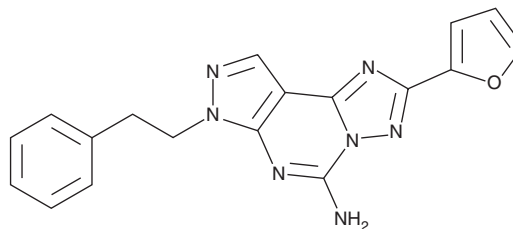
PRODUCT INFORMATION



SCH 58261

Item No. 19676

CAS Registry No.: 160098-96-4
Formal Name: 2-(2-furanyl)-7-(2-phenylethyl)-7H-pyrazolo[4,3-e][1,2,4]triazolo[1,5-c]pyrimidin-5-amine
MF: C₁₈H₁₅N₇O
FW: 345.4
Purity: ≥98%
UV/Vis.: λ_{max}: 232, 262 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

SCH 58261 is supplied as a crystalline solid. A stock solution may be made by dissolving the SCH 58261 in the solvent of choice, which should be purged with an inert gas. SCH 58261 is soluble in organic solvents such as DMSO and dimethyl formamide (DMF). The solubility of SCH 58261 in these solvents is approximately 10 and 15 mg/ml, respectively.

SCH 58261 is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, SCH 58261 should first be dissolved in DMF and then diluted with the aqueous buffer of choice. SCH 58261 has a solubility of approximately 0.33 mg/ml in a 1:2 solution of DMF:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

SCH 58261 is a competitive antagonist of the adenosine A_{2A} receptor (K_is = 2.3 and 2.0 nM in rat and bovine brain, respectively).¹ It is selective for the A_{2A} subtype, displaying an A₁/A_{2A} selectivity ratio of 50-100-fold and does not interact with A_{2B} or A₃ receptors.¹ SCH 58261 has been used to improve 6-hydroxy dopamine-induced motor deficits in a rat model of Parkinson's disease.²

References

1. Zocchi, C., Ongini, E., Ferrara, S., *et al.* Binding of the radioligand [³H]-SCH 58261, a new non-xanthine A_{2A} adenosine receptor antagonist, to rat striatal membranes. *Br. J. Pharmacol.* **117(7)**, 1381-1386 (1996).
2. Reyhani-Rad, S. and Mahmoudi, J. Effect of adenosine A_{2A} receptor antagonists on motor disorders induced by 6-hydroxydopamine in rat. *Acta Cir. Bras.* **31(2)**, 133-137 (2016).

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, 11/16/2022

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