

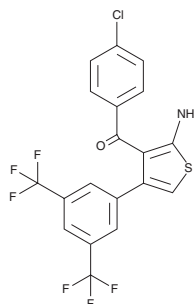
PRODUCT INFORMATION



MIPS521

Item No. 37776

CAS Registry No.: 1146188-19-3
Formal Name: [2-amino-4-[3,5-bis(trifluoromethyl)phenyl]-3-thienyl](4-chlorophenyl)-methanone
MF: $C_{19}H_{10}ClF_6NOS$
FW: 449.8
Purity: $\geq 98\%$
Supplied as: A solid
Storage: $-20^{\circ}C$
Stability: ≥ 4 years



Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

Laboratory Procedures

MIPS521 is supplied as a solid. A stock solution may be made by dissolving the MIPS521 in the solvent of choice, which should be purged with an inert gas. MIPS521 is soluble in methanol.

Description

MIPS521 is an adenosine A_1 receptor positive allosteric modulator.¹ It binds to adenosine A_1 receptors ($K_b = 11 \mu M$) and enhances adenosine-induced inhibition of forskolin-induced cAMP production in CHO cells expressing recombinant human adenosine A_1 receptors when used at concentrations ranging from 0.3 to 30 μM . *In vivo*, MIPS521 (10 and 30 μg /animal) increases the mechanical paw withdrawal threshold in a rat model of neuropathic pain induced by partial sciatic nerve ligation.

Reference

1. Draper-Joyce, C.J., R., B., J., W., *et al.* Positive allosteric mechanisms of adenosine A_1 receptor-mediated analgesia. *Nature* **597**(7877), 571-576 (2021).

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, 11/012023

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