

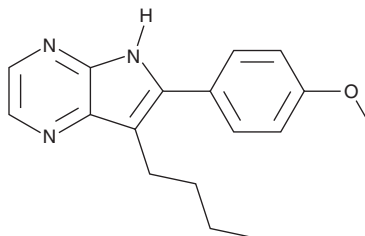
PRODUCT INFORMATION



Aloisine RP106

Item No. 21873

CAS Registry No.: 496864-15-4
Formal Name: 7-butyl-6-(4-methoxyphenyl)-5H-pyrrolo[2,3-b]pyrazine
MF: C₁₇H₁₉N₃O
FW: 281.4
Purity: ≥98%
UV/Vis.: λ_{max}: 260, 354 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

Aloisine RP106 is supplied as a crystalline solid. A stock solution may be made by dissolving the aloisine RP106 in the solvent of choice, which should be purged with an inert gas. Aloisine RP106 is soluble in organic solvents such as DMSO and dimethyl formamide (DMF). The solubility of aloisine RP106 in these solvents is approximately 12 and 25 mg/ml, respectively.

Aloisine RP106 is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, aloisine RP106 should first be dissolved in DMF and then diluted with the aqueous buffer of choice. Aloisine RP106 has a solubility of approximately 0.2 mg/ml in a 1:4 solution of DMF:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

Aloisine is an inhibitor of Cdk1/cyclin B, Cdk5/p25, and GSK3 (IC₅₀s = 0.70, 1.5, and 0.92 μM, respectively).¹ It is a derivative of the aloisines A and B, which competitively inhibit ATP binding to the catalytic subunit of CDKs and GSKs.

Reference

1. Mettey, Y., Gompel, M., Thomas, V., *et al.* Aloisines, a new family of CDK/GSK-3 inhibitors. SAR study, crystal structure in complex with CDK2, enzyme selectivity, and cellular effects. *J. Med. Chem.* **46**(2), 222-236 (2003).

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, 10/27/2022

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