

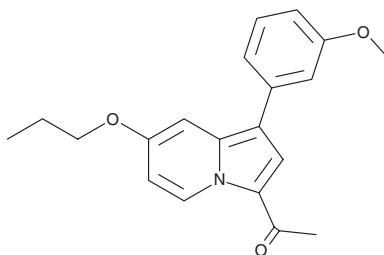
PRODUCT INFORMATION



GSK8573

Item No. 43560

CAS Registry No.: 1693766-04-9
Formal Name: 1-[1-(3-methoxyphenyl)-7-propoxy-3-indoliziny]-ethanone
MF: $C_{20}H_{21}NO_3$
FW: 323.4
Purity: $\geq 98\%$
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

GSK8573 is supplied as a solid. A stock solution may be made by dissolving the GSK8573 in the solvent of choice, which should be purged with an inert gas. GSK8573 is sparingly soluble (1-10 mg/ml) in DMSO.

Description

GSK8573 is an inactive derivative of GSK2801 (Item No. 14120), an inhibitor of bromodomain adjacent to zinc finger domain protein 2A (BAZ2A) and BAZ2B, and has been used as a negative control for binding studies for bromodomain-containing proteins.¹

Reference

1. Chen, P., Chaikuad, A., Bamborough, P., *et al.* Discovery and characterization of GSK2801, a selective chemical probe for the bromodomains BAZ2A and BAZ2B. *J. Med. Chem.* **59**(4), 1410-1424 (2015).

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, 09/04/2025

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