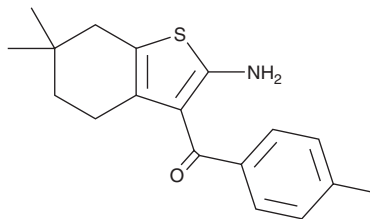


PRODUCT INFORMATION

6,6-Dimethyl-3-(4-methylbenzoyl)-4,5,6,7-tetrahydro-1-benzothiophen-2-amine Item No. 42175

CAS Registry No.: 1461715-47-8
Formal Name: (2-amino-4,5,6,7-tetrahydro-6,6-dimethylbenzo[b]thien-3-yl) (4-methylphenyl)-methanone
MF: C₁₈H₂₁NOS
FW: 299.4
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

6,6-Dimethyl-3-(4-methylbenzoyl)-4,5,6,7-tetrahydro-1-benzothiophen-2-amine is supplied as a solid. A stock solution may be made by dissolving the 6,6-dimethyl-3-(4-methylbenzoyl)-4,5,6,7-tetrahydro-1-benzothiophen-2-amine in the solvent of choice, which should be purged with an inert gas. 6,6-Dimethyl-3-(4-methylbenzoyl)-4,5,6,7-tetrahydro-1-benzothiophen-2-amine is sparingly soluble (1-10 mg/ml) in DMSO and slightly soluble (0.1-1 mg/ml) in ethanol.

Description

6,6-Dimethyl-3-(4-methylbenzoyl)-4,5,6,7-tetrahydro-1-benzothiophen-2-amine is a glucagon-like peptide 1 receptor (GLP-1R) positive allosteric modulator (PAM).¹ It increases GLP-1-induced cAMP production by 268.35% in a reporter assay using HEK293 cells expressing human GLP-1R. 6,6-Dimethyl-3-(4-methylbenzoyl)-4,5,6,7-tetrahydro-1-benzothiophen-2-amine (5 μM) induces insulin secretion in glucose-stimulated GLP-1 (7-36)-induced INS-1 832/13 cells. It decreases blood glucose levels in a glucose tolerance test in fasted mice when administered at doses of 10 or 30 mg/kg.

Reference

1. Campbell, J.A., Do, P., Li, Z., *et al.* Synthesis and biological studies of 2-aminothiophene derivatives as positive allosteric modulators of glucagon-like peptide 1 receptor. *Bioorg. Med. Chem.* **111**, 117864 (2024).

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, 10/28/2024

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