

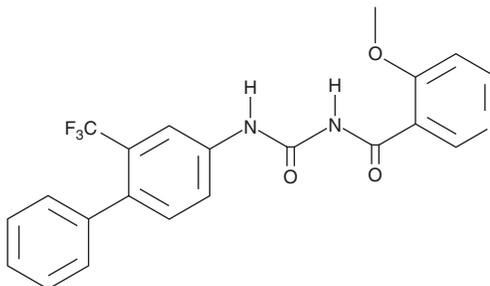
PRODUCT INFORMATION



TC-G 1006

Item No. 17721

CAS Registry No.: 1324003-64-6
Formal Name: 4-methoxy-N-[[[2-(trifluoromethyl)
[1,1'-biphenyl]-4-yl]amino]
carbonyl]-3-pyridinecarboxamide
MF: C₂₁H₁₆F₃N₃O₃
FW: 415.4
Purity: ≥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

TC-G 1006 is supplied as a solid. A stock solution may be made by dissolving the TC-G 1006 in the solvent of choice, which should be purged with an inert gas. TC-G 1006 is sparingly soluble (1-10 mg/ml) in DMSO.

Description

TC-G 1006 is a sphingosine-1-phosphate receptor 1 (S1P₁) agonist.¹ It selectively induces calcium mobilization in cells expressing S1P₁ over cells expressing S1P₂, S1P₃, S1P₄, or S1P₅ (EC₅₀s = 0.035, >25, >25, >50, and 4.3 μM, respectively). TC-G 1006 (0.3, 1, and 3 mg/kg) reduces the number of circulating lymphocytes in rats.

Reference

1. Pennington, L.D., Sham, K.K., Pickrell, A.J., *et al.* 4-Methoxy-N-[2-(trifluoromethyl)biphenyl-4-ylcarbonyl]nicotinamide: A potent and selective agonist of S1P₁. *ACS Med. Chem. Lett.* **2**(10), 752-757 (2011).

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, 02/02/2026

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