

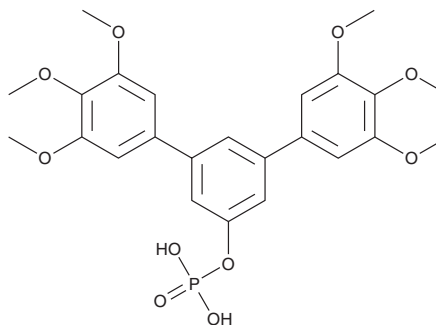
PRODUCT INFORMATION



Stafia-1

Item No. 43191

CAS Registry No.: 2582757-90-0
Formal Name: 3,3'',4,4'',5,5''-hexamethoxy-[1,1':3,1''-terphenyl]-5'-yl dihydrogen phosphate
MF: $C_{24}H_{27}O_{10}P$
FW: 506.4
Purity: $\geq 95\%$
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

Stafia-1 is supplied as a solid. A stock solution may be made by dissolving the stafia-1 in the solvent of choice, which should be purged with an inert gas. Stafia-1 is sparingly soluble (1-10 mg/ml) in ethanol and slightly soluble (0.1-1 mg/ml) in DMSO.

Description

Stafia-1 is an inhibitor of STAT5a ($K_i = 10.9 \mu M$).¹ It is selective for STAT5a over STAT5b and STAT3 (K_i s = $>200 \mu M$ for both).

Reference

1. Natarajan, K., Müller-Klieser, D., Rubner, S., *et al.* Stafia-1: A STAT5a-selective inhibitor developed via docking-based screening of in silico o-phosphorylated fragments. *Chemistry* **26**(1), 148-154 (2020).

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, 04/24/2025

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