

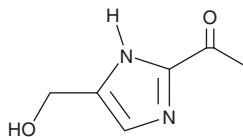
PRODUCT INFORMATION



A6770

Item No. 29972

CAS Registry No.: 1331754-16-5
Formal Name: 1-[5-(hydroxymethyl)-1H-imidazol-2-yl]-ethanone
MF: C₆H₈N₂O₂
FW: 140.1
Purity: ≥98%
UV/Vis.: λ_{max}: 280 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

A6770 is supplied as a crystalline solid. A stock solution may be made by dissolving the A6770 in the solvent of choice, which should be purged with an inert gas. A6770 is soluble in organic solvents such as DMSO and dimethyl formamide. The solubility of A6770 in these solvents is approximately 30 mg/ml.

Further dilutions of the stock solution into aqueous buffers or isotonic saline should be made prior to performing biological experiments. Ensure that the residual amount of organic solvent is insignificant, since organic solvents may have physiological effects at low concentrations. Organic solvent-free aqueous solutions of A6770 can be prepared by directly dissolving the crystalline solid in aqueous buffers. The solubility of A6770 in PBS (pH 7.2) is approximately 10 mg/ml. We do not recommend storing the aqueous solution for more than one day.

Description

A6770 is an inhibitor of sphingosine-1-phosphate (S1P) lyase.¹ It induces accumulation of [³H] sphinganine-1-phosphate ([³H]dhS1P), an S1P lyase substrate, in IT-79MTNC3 cells that endogenously express high levels of S1P lyase (EC₅₀ = <0.01 μM), an effect that is reduced in the presence of vitamin B₆ (EC₅₀ = <100 μM).

Reference

1. Ohtoyo, M., Tamura, M., Machinaga, N., *et al.* Scintillation proximity assay to detect the changes in cellular dihydrosphingosine 1-phosphate levels. *Lipids* **51**(10), 1207-1216 (2016).

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, 11/14/2022

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