

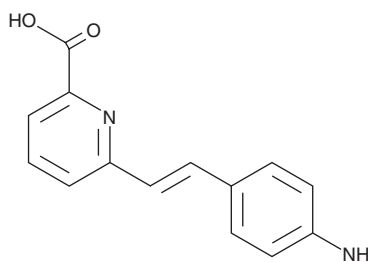
PRODUCT INFORMATION



CB-7921220

Item No. 23332

CAS Registry No.: 115453-99-1
Formal Name: 6-[2-(4-aminophenyl)ethenyl]-2-pyridinecarboxylic acid
MF: C₁₄H₁₂N₂O₂
FW: 240.3
Purity: ≥98%
UV/Vis.: λ_{max}: 353 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

CB-7921220 is supplied as a crystalline solid. A stock solution may be made by dissolving the CB-7921220 in the solvent of choice, which should be purged with an inert gas. CB-7921220 is soluble in organic solvents such as DMSO and dimethyl formamide. The solubility of CB-7921220 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 CB-7921220 can be prepared by directly dissolving the crystalline solid in aqueous buffers. The solubility of CB-7921220 in PBS, pH 7.2, is approximately 0.3 mg/ml. We do not recommend storing the aqueous solution for more than one day.

Description

CB-7921220 is an inhibitor of the adenylyl cyclase (AC) isoforms AC1 and AC6.¹ It is selective, reducing AC1 and AC6 activity by 60% while having no effect on AC2 and AC5 in Sf9 membranes stimulated by forskolin (Item No. 11018) when used at a concentration of 100 μM.

Reference

1. Brand, C.S., Hocker, H.J., Gorfe, A.A., *et al.* Isoform selectivity of adenylyl cyclase inhibitors: Characterization of known and novel compounds. *J. Pharmacol. Exp. Ther.* **347**(2), 265-275 (2013).

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, 12/15/2022

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